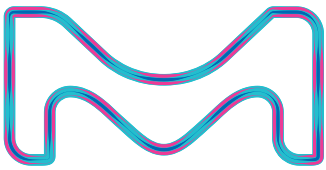


Protein Blotting Handbook

Tips and Tricks



The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

Millipore®

Preparation, Separation,
Filtration & Monitoring Products

About the Seventh Edition

With the publication of the seventh edition of the Protein Blotting Handbook, MilliporeSigma continues to keep researchers up to date on innovations in protein detection. We've added more comprehensive guidance on optimizing antibody concentrations and reducing background, along with expanded data and protocols for fluorescence detection.

This handbook represents the collective experience of our application scientists, who are actively engaged in advancing the science of protein blotting and detection. It also includes many of the most common recommendations provided by our technical service specialists who are contacted by scientists worldwide for assistance.

Better membranes, better blots.

MilliporeSigma has been a leading supplier of transfer membranes for nearly four decades. E.M. Southern used these membranes to develop the first nucleic acid transfer from an agarose gel in 1975¹. The first 0.45 µm PVDF substrate for Western blotting, Immobilon®-P membrane, was introduced in 1985, and the first 0.2 µm PVDF substrate for protein blotting and sequencing, Immobilon®-P^{5Q} membrane, was introduced in 1988.

In addition to Immobilon® membranes and reagents, MilliporeSigma provides a wide selection of other tools for protein research, including gentle protein extraction kits, rapid protein isolation with PureProteome™ magnetic beads, and fast, effective concentration with Amicon® Ultra centrifugal filters.

Where to find additional information

If you have questions or need assistance, please contact a MilliporeSigma technical service specialist, or pose your question online at: **[SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice)**

You'll also find answers to frequently asked questions (FAQs) concerning Western blotting and other related methods at: **[SigmaAldrich.com/WesternHelp](https://www.sigmaaldrich.com/WesternHelp)**

Table of Contents

INTRODUCTION	3	Mass Spectrometry	24
1 MEMBRANE SELECTION	4	Mass Spectrometry with Immobilon® PVDF Transfer Membranes	24
Immobilon® PVDF Transfer Membranes	5		
2 PROTEIN BINDING	8	5 PROTOCOLS	26
Immobilon®-P vs. Immobilon®-P ^{sq} Transfer Membranes	8	1 Protein Transfer Protocols	27
Factors Affecting Protein Binding	8	1.1 Electrotransfer: Tank (Wet) Transfer	27
		1.2 Electrotransfer: Semi-dry Transfer	28
3 BLOTTING METHODS: PRINCIPLES AND OPTIMIZATION	9	1.3 Dot Blotting/Slot Blotting: Vacuum Filtration Method	30
Filtration	9	1.4 Dot Blotting: Manual Spotting Method	31
Western Blotting	10	1.5 Optimization of Blocking Reagents	31
Separation of Complex Protein Mixtures on 1-D or 2-D Gels	10	1.6 Double Blotting to Eliminate NSB	32
Molecular Weight Markers	11	1.7 Membrane Drying Methods	32
Polyacrylamide Concentration	11	2 Protein Visualization Protocols	33
Gel Running Buffers	11	2.1 Visualization by Transillumination	33
Transfer of Proteins from Gel to Membrane	12	2.2 Visualization by Reversible Staining	34
Electrotransfer Techniques	12	Ponceau-S Red	34
Transfer Buffers	13	CPTS (Copper Phthalocyanine Tetrasulfonic Acid)	34
Functions of Methanol in Transfer Buffer	14	2.3 Visualization by Irreversible Staining	35
Factors Affecting Successful Protein Transfer	14	Coomassie® Brilliant Blue R Dye	35
Presence of SDS	14	Amido Black	35
Current and Transfer Time	15	3 Immunodetection Protocols	36
Transfer Buffer pH	15	3.1 Standard Immunodetection Method	36
Equilibration Time	15	3.2 Rapid Immunodetection Method	38
Developing a New Transfer Protocol	16	3.3 Simple Immunodetection Using the Immobilon® GO Device	39
Preparing Membrane for Protein Identification	16	3.4 Rapid Immunodetection Using the SNAP i.d.® 2.0 System	40
For Staining and Immunodetection	16	3.5 High Salt Wash to Remove Persistent Background	41
For Rapid Immunodetection and Visualization by Transillumination	16	4 Membrane Stripping Protocols	42
Storage	16	4.1 Stripping by Heat and Detergent	42
Protein Visualization	16	4.2 Stripping by Low pH	42
Transillumination	16	4.3 Stripping with the ReBlot™ Plus Western Blot Recycling Kit	42
Staining	17	5 Protein Digestion Protocol	43
Reversible Stains	17	5.1 On-Membrane Protein Digestion for Mass Spectrometry	43
Irreversible Stains	17	6 Blot Storage Protocol	44
		6.1 Preparation of Blots for Long-Term Storage	44
4 PROTEIN IDENTIFICATION	18	6 APPENDICES	45
Immunodetection	18	Standard Recipes for Western Blotting	45
Standard vs. Rapid Immunodetection Procedures	18	Recipes for mPAGE™ Precast Gel System	46
Factors Influencing Immunodetection	19	Troubleshooting	47
Buffers	19	Glossary	53
Blocking	19	References	55
Antibodies	20	Patents	56
Washing	21	7 ORDERING INFORMATION	57
Double Blotting	21		
Detection Substrates	21		
Chromogenic Detection	22		
Chemiluminescent Detection	22		
Fluorescent Detection	23		
Reprobing Immobilon® PVDF Transfer Membranes	23		

Introduction

Since its introduction in 1979, protein blotting¹⁷ has become a routine tool used to detect low amounts of proteins in complex samples or to monitor protein expression and purification. The simplest protein blotting procedure, known as dot blot or slot blot, uses vacuum filtration to transfer protein onto a microporous membrane. While this method may provide qualitative information about total protein expression levels and can be performed on multiple samples in parallel, it lacks information on protein molecular weight. Also, specificity can be compromised as protein degradation products or post-translationally modified isoforms may be detected along with the intact protein.

A more complex procedure, Western blotting, involves the separation of a protein mixture by gel electrophoresis with subsequent electrotransfer to a suitable membrane (e.g., PVDF). After proteins have been transferred onto a membrane, they can be stained for visualization and directly identified by N-terminal sequencing, mass spectrometry or immunodetection. Immunodetection involves the identification of a protein through its reaction with a specific antibody. Through spatial resolution, Western blotting provides molecular weight information on individual proteins and distinguishes isoforms, alternate processing products, and other post-translationally modified forms.

In the clinical laboratory, immunoblotting has proven useful in fields such as infectious and autoimmune diseases, allergy and others.^{48,49} Western blotting is considered to be a reliable confirmatory diagnostic test following a repeatedly reactive ELISA over the course of viral infection, and is a sensitive, unequivocal and simple assay, providing high complexity of information.⁵⁰⁻⁵²

Examples of Western blotting applications include analysis of protein expression in yeast by quantitative Western analysis⁵³, determination of protein copy number and compartmentalization⁵⁴, study of competitive protein kinase inhibition by ATP⁵⁵, and detection of genetically modified organisms in crops and foods⁵⁶.

Clearly, protein blotting remains the platform of choice for exploratory research, and is still the standard by which new antibodies and other protein detection assays (such as ELISA, bead-based assays, flow cytometry and immunohistochemistry) are evaluated. However, the need to analyze more proteins simultaneously to characterize complex networks and the associated need to conserve valuable samples has driven ongoing research into improving the sensitivity and speed of blotting techniques. Double blotting^{35,36} eliminates false positives due to strong nonspecific interactions between the blotted proteins and unrelated secondary antibodies. Far-Western blotting enables the detection of specific protein-protein interactions⁵⁷ and Southwestern blotting is used to identify proteins that interact with specific DNA sequences⁵⁸. Multistrip Western blotting has proved to increase throughput while minimizing inter-blot variability⁵⁹. A new generation of blotting technologies features reductions in the amounts of protein required to produce a signal⁶⁰ and methods to improve the quantitative power of Western blotting^{61, 62}.

1. Membrane Selection

Membrane type can influence the following factors in immunoblotting:

- Protein binding capacity
- Requirement for prewetting with alcohol
- Ability to perform multiple stripping and reprobing experiments
- Protein visualization
- Long-term blot storage
- Signal-to-noise ratio

Polyvinylidene fluoride (PVDF) and nitrocellulose membranes are most commonly used in Western blotting.

In comparison to nitrocellulose, PVDF membranes offer better protein retention, physical strength and

broad chemical compatibility² (Table 1). The higher mechanical strength and superior chemical resistance of PVDF membranes make them ideal for multiple immunodetection staining applications and reprobing. Another advantage of using PVDF membranes is that replicate lanes from a single gel can be used for various applications, such as staining with Coomassie® Blue dye followed by band excision and N-terminal sequencing, proteolysis/peptide separation/internal sequencing, and immunodetection³. Typical binding capacity of commercially available nitrocellulose membranes is 80 – 100 µg/cm², while PVDF membranes offer a binding capacity of 150 – 200 µg/cm².

In comparison to nitrocellulose membranes, PVDF membranes were shown to have better retention of total HIV antigens and improved detection of antibodies to glycosylated envelope antigens⁴.

Table 1. Membrane attributes and applications

Attributes/Applications	Nitrocellulose	PVDF
Physical strength	Poor	Good
Protein binding capacity	80 – 100 µg/cm ²	150 – 200 µg/cm ²
Solvent resistance	No	Yes
Western transfer	Yes	Yes
Total protein stain	Colloidal gold	Colloidal gold
	Ponceau-S red	Ponceau-S red
	Amido black	Amido black
	India ink	India ink
	Sypro® blot stains	Coomassie® Blue dye
Detection	Chromogenic	Chromogenic
	Chemiluminescent	Chemiluminescent
	Fluorescent	Fluorescent
	Radioactive	Chemifluorescent
		Radioactive
Double-blotting method	No	Yes
Rapid immunodetection	No	Yes
Western reprobing	Yes	Yes
Edman sequencing	No	Yes
Amino acid analysis	Yes	Yes
Binding in the presence of SDS	Poor	Good
On-membrane digestion or mass spectrometry	No	Yes
Direct MALDI-TOF MS analysis	No	Yes
Blot can be archived	No	Yes

Immobilon® PVDF Transfer Membranes

MilliporeSigma offers five PVDF membranes:

- **Immobilon®-P membrane** (0.45 μm) is a versatile substrate that is well suited for common immunoblotting applications.
- **Immobilon®-P^{sq} membrane** (0.2 μm) is ideal for protein sequencing and immunoblotting of low molecular weight proteins. It has a higher protein binding capacity and a higher retention than 0.45 μm membranes.
- **Immobilon®-FL membrane** (0.45 μm) was developed for fluorescence-based immunodetection. It has very low background fluorescence across a wide range of excitation and emission wavelengths.
- **Immobilon®-E membrane** (0.45 μm) wets out in water or buffer, eliminating the alcohol pre-wet step while retaining similar performance to Immobilon®-P membranes.

Immobilon® transfer membranes are available in rolls, cut sheets, and the Immobilon® NOW format. Precut membranes are compatible with all precast gels and most commercially available gel running systems. Yielding midi or mini sized sheets with a single cut, the Immobilon® NOW roll format provides users convenience and flexibility.

The tables on the following pages cover the properties of each type of Immobilon® transfer membrane (Table 2), as well as match precut Immobilon membranes to commercially available electrophoresis systems (Table 3) and precast gels (Table 4).



Table 2. Properties and applications of Immobilon® PVDF transfer membranes.

	Immobilon®-P transfer membrane	Immobilon®-P ^{5Q} transfer membrane	Immobilon®-FL transfer membrane	Immobilon®-E transfer membrane
Description	Optimized to bind proteins transferred from a variety of gel matrices	Uniform pore structure results in superior binding of proteins with MW <20 kDa	Optimized for fluorescence immunodetection applications	First PVDF transfer membrane for Western blotting that wets out in transfer buffer or water
Composition	PVDF	PVDF	PVDF	PVDF
Pore size	0.45 µm	0.2 µm	0.45 µm	0.45 µm
Phobicity	Hydrophobic	Hydrophobic	Hydrophobic	Hydrophobic
Applications	Western blotting	Low molecular weight	Western blotting	Western blotting
	Binding assays	Western blotting	Dot/slot blotting	
	Amino acid analysis	Amino acid analysis	Fluorescence immunodetection	
	N-terminal protein sequencing	Mass spectrometry		
	Dot/slot blotting	N-terminal protein sequencing		
	Glycoprotein visualization			
	Lipopolysaccharide analysis			
	Mass spectrometry			
Detection methods	Chromogenic	Chromogenic	Fluorescent	Chemiluminescence
	Chemifluorescent	Fluorescent	Chromogenic	Chromogenic
	Chemiluminescent	Chemifluorescent	Chemifluorescent	Radioactive
	Radioactive	Chemiluminescent	Chemiluminescent	
		Radioactive		
Protein binding capacity	Insulin: 160 µg/cm ²	Insulin: 262 µg/cm ²	Insulin: 155 µg/cm ²	Insulin: 160 µg/cm ²
	BSA: 215 µg/cm ²	BSA: 340 µg/cm ²	BSA: 205 µg/cm ²	BSA: 215 µg/cm ²
	Goat IgG: 294 µg/cm ²	Goat IgG: 448 µg/cm ²	Goat IgG: 300 µg/cm ²	Goat IgG: 294 µg/cm ²
Compatible stains reversible	Transillumination	Transillumination	Transillumination	Transillumination
	Ponceau-S	Ponceau-S	Ponceau-S	Ponceau-S red
	CPTS	CPTS	CPTS	Sypro® ruby
	Toluidine blue	Toluidine blue	Sypro® blot stains	
	Sypro® blot stains	Sypro® blot stains		
Compatible stains irreversible	Coomassie® Brilliant Blue dye	Coomassie® Brilliant Blue dye	Coomassie® Brilliant Blue dye	Coomassie® brilliant blue
	Amido black	Amido black	Amido black	Amido black
	India ink	India ink		India ink
	Colloidal gold	Colloidal gold		Colloidal gold

Immobilon® NOW membrane format

Providing the ultimate convenience in Western blotting, our Immobilon® NOW membrane format yields both midi and mini-sized blots in a single cut. All four Immobilon® membrane types (P, P^{5Q}, FL, and E) are available in the Immobilon® NOW format. In tables 3 and 4 below, Immobilon® cut sheets are paired with matching electrophoresis systems and precast gels. Alternatively, the Immobilon® NOW format can be used in the place of cut sheets, when one dimension is 8.5 cm. A column has been added to the table to indicate when the Immobilon® NOW roll format is a possible alternative to cut sheets.

Description	Cat. No.
Immobilon®-P Transfer Membrane	IPVH85R
Immobilon®-P ^{5Q} Transfer Membrane	ISEQ85R
Immobilon®-FL Transfer Membrane	IPFL85R
Immobilon®-E Transfer Membrane	IEVH85R

Table 3. Immobilon® PVDF transfer membrane cut sheets and matching electrophoresis systems

Manufacturer	Vertical Gel Box	Gel size (cm)	Immobilon® Size (cm)	Immobilon® NOW format compatible	Immobilon®-P 0.45 µm	Immobilon®-P ^{SQ} 0.2 µm	Immobilon®-FL 0.45 µm*	Immobilon®-E 0.45 µm
GE Healthcare	SE 250 Mighty Small™	8 x 7	8.4 x 7	■	IPVH07850	ISEQ07850	IPFL07810	IEVH07850
	SE 260 Mighty Small™	8 x 9.5	8 x 10	■	IPVH08100	ISEQ08100		IEVH08100
	miniVE	8 x 9.5	8 x 10	■	IPVH08100	ISEQ08100		IEVH08100
		8.5 x 10	10 x 10	■	IPVH10100	ISEQ10100	IPFL10100	
	SE 400	14 x 16	15 x 15	■	IPVH15150	ISEQ15150		
SE 600	14 x 16	15 x 15	■	IPVH15150	ISEQ15150			
	14 x 8	13.5 x 8.5	■	IPVH08130	ISEQ08130			
Bio-Rad	Mini-PROTEAN® 3 Dodeca™	8.3 x 7.3	8.4 x 7	■	IPVH07850	ISEQ07850	IPFL07810	IEVH07850
	Criterion™, Criterion™ Dodeca™	13.3 x 8.7	13.5 x 8.5	■	IPVH08130	ISEQ08130		
	PROTEAN® II xi	16 x 16	15 x 15	■	IPVH15150	ISEQ15150		
		16 x 20	26.5 x 37.5	■	IPVH00010	ISEQ00010	IPFL00010	
	PROTEAN® II XL	19.3 x 18.3	20 x 20	■	IPVH20200	ISEQ20200	IPFL20200	
	PROTEAN® Plus Dodeca™	20 x 20.5	26 x 26	■	IPVH304F0	ISEQ304F0		
Mini-PROTEAN® Tetra	8.3 x 7.3	8.4 x 7	■	IPVH07850	ISEQ07850	IPFL07810	IEVH07850	
Invitrogen	XCell SureLock™ Mini-Cell	8 x 8	8.4 x 7	■	IPVH07850	ISEQ07850	IPFL07810	IEVH07850
	XCell4 SureLock™ Midi	8 x 13	8.5 x 13.5	■	IPVH08130	ISEQ08130		
Thermo Scientific	Owl® P81, P82, P8DS	10 x 10	10 x 10	■	IPVH10100	ISEQ10100	IPFL10100	
	Owl® P82, P8DS	8 x 10	8 x 10	■	IPVH08100	ISEQ08100		IEVH08100
	Owl® P9DS	16 x 16	15 x 15	■	IPVH15150	ISEQ15150		
	Owl® P10PS	20 x 20	20 x 20	■	IPVH20200	ISEQ202 00	IPFL20200	

Table 4. Immobilon® PVDF transfer membrane cut sheets and matching pre-cast gels

Manufacturer	Vertical Gel Box	Gel size (cm)	Immobilon® Size (cm)	Immobilon® NOW format compatible	Immobilon®-P 0.45 µm	Immobilon®-P ^{SQ} 0.2 µm	Immobilon®-FL 0.45 µm*	Immobilon®-E 0.45 µm
Bio-Rad	Mini-PROTEAN™ TGX, Mini-PROTEAN™ TGX Stain-Free	8.6 x 6.8	8.4 x 7	■	IPVH07850	ISEQ07850		IPFL07810
	Criterion™ TGX, Criterion™ TGX Stain-Free, Criterion™ Tris-HCl, Criterion™ Stain-Free Tris-HCl, Criterion™ XT Bis-Tris	13.3 x 8.7	13.5 x 8.5	■	IPVH08130	ISEQ08130		
Lonza	PAGEr® Gold	9 x 10	8 x 10	■	IPVH08100	ISEQ08100		IEVH08100
		10 x 10	10 x 10	■	IPVH10100	ISEQ10100	IPFL10100	
NuSep	nUView Tris-Glycine, Stain Free	8.6 x 6.8	8.4 x 7	■	IPVH08100	ISEQ08100	IPFL07810	IEVH07850
		8 x 8	8 x 10	■	IPVH08100	ISEQ08100		IEVH08100
Invitrogen	NuPAGE®, Bolt®	8 x 8	8.4 x 7	■	IPVH07850	ISEQ07850	IPFL07810	IEVH07850
	Novex®	8 x 8	8.4 x 7	■	IPVH07850	ISEQ07850	IPFL07810	IEVH07850
	E-PAGE	13.5 x 10.8	13.5 x 8.5	■	IPVH08130	ISEQ08130		
Expedeon	RunBlue™ TEO-Tricine SDS	8 x 10	8 x 10	■	IPVH08100	ISEQ08100		IEVH08100
	RunBlue™ Tris-Bis	10 x 10	10 x 10	■	IPVH10100	ISEQ10100	IPFL10100	
Genscript	ExpressPlus™	8 x 10	8 x 10	■	IPVH08100	ISEQ08100	IPFL10100	IEVH08100
	SurePAGE™	8.6 x 6.8	8.4 x 7	■	IPVH07850	ISEQ07850	IPFL07810	IEVH07850
MilliporeSigma	mPAGE™	8.6 x 6.8	8.4 x 7	■	IPVH07850	ISEQ07850	IPFL07810	IEVH07850

2. Protein Binding

Due to the hydrophobicity of PVDF, traditional PVDF membranes will not wet out in aqueous solutions and must first be wet in $\geq 50\%$ (v/v) alcohol solution. Methanol, ethanol, and isopropanol solutions are all suitable to wet PVDF membranes. Complete wetting is evident by a change in the membrane's appearance from opaque to semi-transparent. The alcohol must be removed from the membrane prior to equilibration in transfer buffers by extensive rinsing in water. In contrast to traditional PVDF membranes, our Immobilon®-E membrane wets out directly in water or buffer, eliminating the need for alcohol pre-wetting and removal.

Binding Differences between Immobilon®-P and Immobilon®-P^{SQ} Transfer Membranes

After wetting, protein binding can be achieved by simply bringing the protein into contact with the membrane. Because binding occurs throughout the depth of the membrane, the binding capacity is determined by the internal surface area of the pores⁵. Immobilon®-P^{SQ} transfer membrane has approximately three times the internal surface area of Immobilon®-P transfer membrane, resulting in higher adsorptive capacity (Table 2, Page 6). The values listed in Table 2 represent upper limits for protein binding after saturation of the membrane surface in a non-denaturing buffer. However, the maximum binding that can be achieved will depend on the specific protocols employed, variations in the structural conformation of the proteins, the chemical nature of the buffers used, and the limitations of the methods used to apply the sample.

An example of the binding difference between Immobilon®-P and Immobilon®-P^{SQ} transfer membranes is shown in Figure 1, where protein samples were electrotransferred from a polyacrylamide gel. A fraction of the proteins passed through the Immobilon®-P transfer membrane and were captured on a second membrane placed behind the first one. In contrast, all of the proteins were bound to the Immobilon®-P^{SQ} membrane without passing through it. In this case, the tighter pore structure and higher internal membrane surface area facilitated complete adsorption of all of the transferred protein. However, immunodetection on Immobilon®-P^{SQ} transfer membrane can result in a higher background and can require more stringent washing conditions. Thus, membrane choice is dictated by the goal of the experiment: use Immobilon®-P transfer membrane for high-sensitivity detection of > 20 kDa proteins, but switch to Immobilon®-P^{SQ} transfer membrane if smaller proteins are being analyzed or 100% protein capture is necessary for peptide sequencing.

Factors Affecting Protein Binding

At the molecular level, protein adsorption results, at least in part, from the interaction of hydrophobic amino acid side chains and hydrophobic domains with the polymer surface. Matsudaira observed an 80% decline in the sequencing efficiency of small peptides after hydrophobic residues were cleaved, presumably due to washout of the peptide remnants⁶. Also, in peptide digestions, it has been observed that peptides characterized as hydrophobic often do not elute from the membrane as efficiently as more hydrophilic peptides^{7, 8}. McKeon and Lyman demonstrated that addition of Ca^{2+} ions to the transfer buffer enhanced the binding of calmodulin to Immobilon®-P transfer membrane⁹. Binding of the calcium to the protein causes formation of a hydrophobic pocket in the molecule's structure, leading to enhanced binding to the hydrophobic PVDF surface.

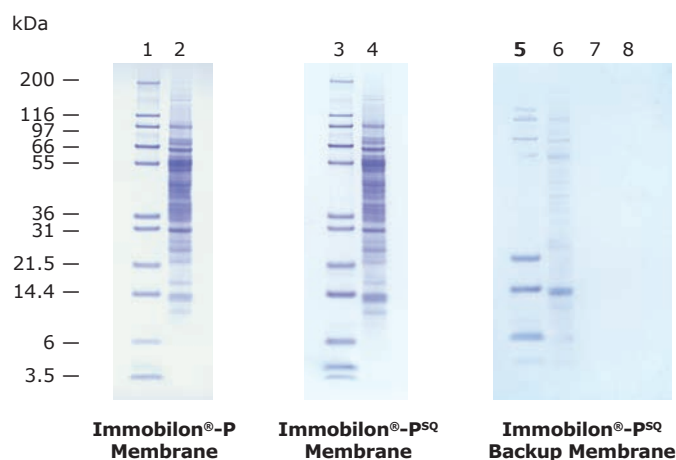


Figure 1.

Prolonged electrotransfer of proteins using Immobilon®-P and Immobilon®-P^{SQ} transfer membranes. Molecular weight standards (lanes 1,3,5,7) and calf liver lysate (lanes 2,4,6,8) were transferred to Immobilon®-P or Immobilon®-P^{SQ} membranes by the tank transfer method and stained with Coomassie® Blue dye. A sheet of Immobilon®-P^{SQ} transfer membrane was placed behind the primary membranes to capture proteins that passed through them. (Lanes 5 and 6 behind Immobilon®-P; lanes 7 and 8 behind Immobilon®-P^{SQ}.)

3. Blotting Methods: Principles and Optimization

Filtration

Proteins can be directly applied to membranes by vacuum-driven filtration (Figure 2). Proteins adsorb onto the membrane, and the other sample components are pulled through by the vacuum (Figure 2). Alternatively, the sample can be spotted directly onto the surface and allowed to dry. The proteins immobilized on the membrane can then be analyzed.

Dot blotting (Figure 3) and slot blotting are two variations of the filtration method, employing manifolds that permit sample application in dot or slot patterns. These techniques can be used as qualitative method for rapid screening or as a quantitative technique for analysis of similar samples. It is especially useful for testing the suitability of experimental design parameters to be used in complex analyses.

Another variation of the filtration method is grid immunoblotting, a technique useful for highly parallel sample analysis when the sample quantity is extremely limited and analysis cannot be performed by conventional techniques such as ELISA. For example, grid

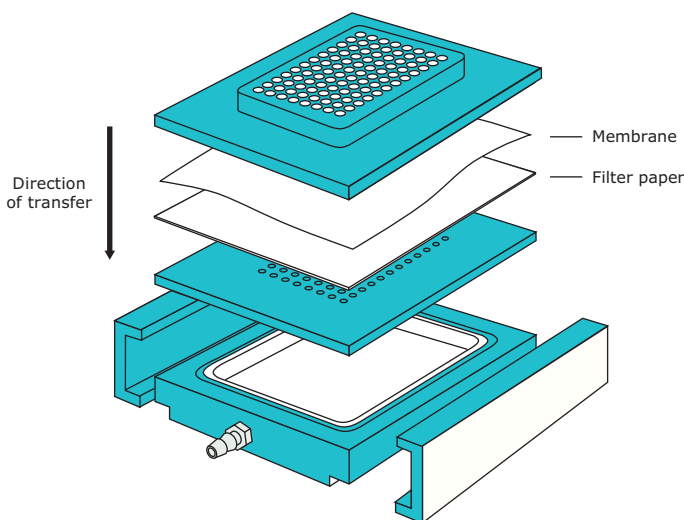


Figure 2.
A blotting system using filtration.

immunoblotting has been used in the characterization of allergen-specific antibody response with minimal amounts of patient serum¹⁰.

When preparing blots by filtration, consider the following:

- Detergents can inhibit the adsorption of proteins to the membrane. Buffers used for sample dissolution and washing should contain no more than 0.05% detergent, and only if required.
- The sample volume should be large enough to cover the exposed membrane in each well, but should not contain protein in excess of the binding capacity of the membrane.
- Samples with high particulate loads may clog the membrane, while those with high viscosity will reduce the flow rate. Particles should be removed by prefiltration or centrifugation, and only the supernatant should be applied to the membrane. Viscous samples should be diluted in buffer.

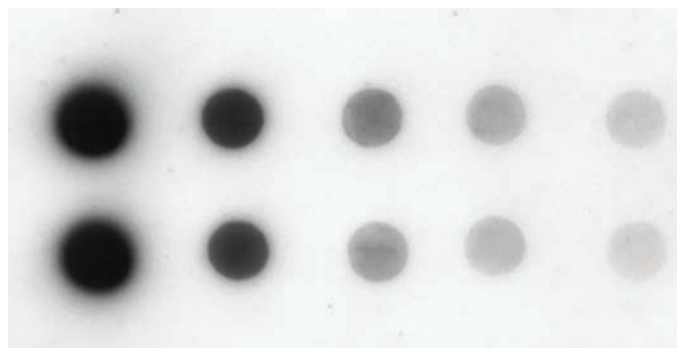


Figure 3.
Detection of transferrin on a dot blot of human serum (See Protocol 1.4. Dot Blotting, page 31) using Immobilon® Western Chemiluminescent HRP Substrate. Transferrin was detected in duplicate serial dilutions of serum with goat anti-transferrin antibody (dilution 1:10,000) and rabbit anti-goat HRP-conjugated secondary antibody (dilution 1:50,000) on Immobilon®-P membrane.

Western Blotting

Western blotting comprises the following steps:

1. Resolution of a complex protein sample in a polyacrylamide gel
2. Transfer of the resolved proteins onto a membrane
3. Identification of a specific protein on the membrane

For a successful Western blot, four requirements must be met:

- Elution from the gel — the protein must elute from the gel during transfer. If it is retained in the gel, it will not be available for analysis on the blot.
- Adsorption to the membrane — the protein must adsorb to the membrane during the transfer process. If the protein is not adsorbed, it will not be available for analysis on the blot.

- Retention during processing — the protein must remain adsorbed to the membrane during post-transfer processing of the blot.
- Accessibility during processing — the adsorbed protein must be available for antibody binding. If the protein is masked, it cannot be detected.

The sections that follow discuss theoretical and practical considerations of the protocols involved in Western blotting.

Separation of Complex Protein Mixtures in 1-D or 2-D Gels

One-dimensional (1-D) sodium dodecyl sulfate-polyacrylamide gel electrophoresis (SDS-PAGE) is commonly used to separate proteins by molecular weight prior to blotting (Figure 4). In some cases, non-denaturing conditions are used to separate native proteins. Although this method usually lacks the resolution of denaturing electrophoresis, it may be useful when the primary antibody only recognizes non-denatured proteins or when the protein's biological activity must be retained.

Two-dimensional (2-D) gel electrophoresis is the technique of choice for analyzing protein composition of cell types, tissues and fluids, and is a key technology in proteomics (Figure 5). Immunoblotting of 2-D gels provides information on molecular weight and isoelectric point and can be useful to discriminate protein isoforms generated by post-translational modifications¹¹. In some cases, protein phenotyping can be achieved by immunoblotting after only a 1-D separation by isoelectrofocusing^{12,13}.

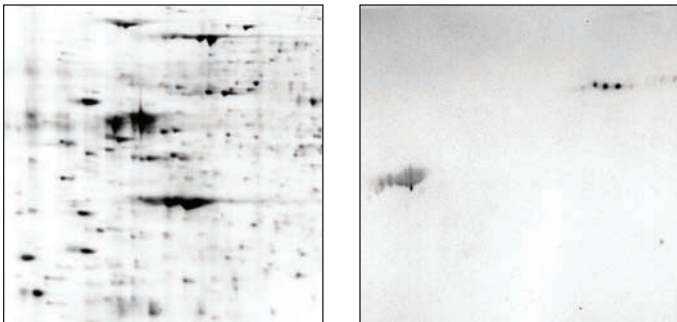


Figure 5.

Chemiluminescent detection of proteins separated by two-dimensional electrophoresis. Silver-stained 2-D gel of rat fibroblast cell line (left) and blot of the same gel (right), probed with a mouse monoclonal antibody and visualized using chemiluminescence on Immobilon[®]-P membrane.

Molecular Weight Markers

The inclusion of molecular weight (MW) standards, or markers, facilitates the estimation of the sizes of the proteins of interest after resolution by electrophoresis. Two types are available, unstained and pre-stained.

Unstained MW markers usually consist of a mixture of purified native or recombinant proteins of defined molecular weights. Visualizing their location on a gel or membrane requires a staining step.

Pre-stained markers allow monitoring of protein separation in the gel during electrophoresis and indicate transfer efficiency in subsequent blotting steps. However, they can be relatively expensive and the addition of dyes may affect protein mobility. Pre-stained markers may be less accurate for molecular weight determination, as dyes attached to the proteins may alter their ability to adsorb to the membrane during blotting.

Polyacrylamide Concentration

Polyacrylamide concentration can be homogenous throughout the gel or a gradient. The most common polyacrylamide concentration, 10%, is best suited for the separation of proteins in the range of 10 – 150 kDa. If unknown proteins are being analyzed or a broader range of separation is desired, gradient gels are recommended. For example, 4 – 12% Tris-glycine gels are suitable for proteins in the range of 30 to 200 kDa, while 10 – 20% gels will successfully separate proteins from 6 to 150 kDa. SDS-PAGE gels are usually 1.0 and 1.5 mm thick; however, for blotting and proteins transfer best out of thinner gels (= 1 mm).

Gel Running Buffers

Gel running buffers are typically composed of tris-glycine or tris-tricine and may contain 0.1% detergent, usually SDS. Tris-glycine buffer systems are useful for separating a wide range of protein molecular weights (6 – 200 kDa) and are compatible with denaturing or non-denaturing conditions. Tris-tricine systems are best for the separation of smaller proteins (< 10 kDa) that need to be reduced and denatured prior to loading. Both buffer systems are compatible with protein transfer to PVDF membranes.

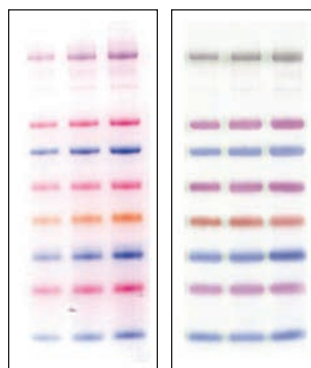


Figure 4.

ColorBurst molecular weight markers transferred to nitrocellulose after resolution by SDS-PAGE. A 10-20% Tris-Glycine gel was loaded (left to right) with 3, 5 and 7 μ L of the ColorBurst Marker. The markers were run using standard conditions on 10 x 10 cm, 1 mm thick, 10-well precast gels.

Bands transferred to nitrocellulose membranes from the gels in Figure 1 (Tris-Glycine on left and Tris-Tricine on right). Transfers were completed in 90 minutes at 70 volts with Towbin's buffer (Tris-Glycine in 20% methanol.)

Transfer of Proteins from Gel to Membrane

The process of transferring proteins from a gel to a membrane while maintaining their relative position and resolution is known as blotting. Blotting can be achieved in three different ways:

Simple diffusion¹⁴ is accomplished by laying a membrane on top of the gel with a stack of dry filter paper on top of the membrane, and placing a weight on top of the filter paper to facilitate the diffusion process³. This method can be used to transfer proteins from one gel to multiple membranes, obtaining several imprints of the same gel. The major disadvantage of the diffusion method is that transfer is not quantitative and only transfers 25 – 50% of the proteins, compared to electroblotting¹⁵.

Vacuum-assisted solvent flow¹⁶ uses the suction power of a pump to draw separated proteins from the gel onto the membrane. Both high and low molecular weight proteins can be transferred by this method; however, a smaller pore size membrane (0.2 μm) may be needed for proteins with MW < 20 kDa, since they are less readily adsorbed by the 0.45 μm membrane³. Vacuum blotting of proteins out of polyacrylamide gels is uncommon and is mostly used for nucleic acid transfer from agarose gels.

Electrotransfer Techniques

Electrophoretic elution, or electrotransfer¹⁷ is the most commonly used transfer method in protein blotting. The principal advantages are the speed and completeness of transfer compared to diffusion or vacuum blotting.

The two commonly used electrotransfer techniques are tank transfer and semi-dry transfer. Both are based on the same principles and differ only in the mechanical devices used to hold the gel/membrane stack and application of the electrical field.

In tank transfer, (Figure 6) the gel/membrane stack is completely immersed in a buffer reservoir and current is applied. Typically run at constant voltage, mixing the buffer during tank transfer typically keeps the current relatively constant. While this method is effective, tank transfer is a slow technique that requires large volumes of buffer.

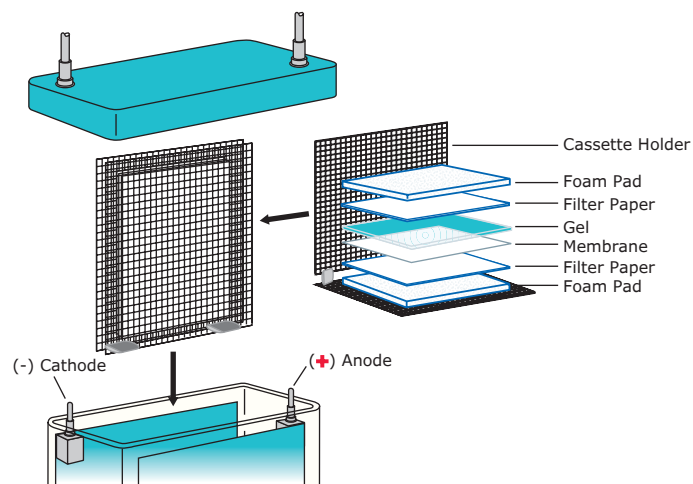


Figure 6.
Tank (wet) transfer system.

Semi-dry transfer (Figure 7) replaces the buffer reservoir with layers of filter paper soaked in buffer and is the preferred method for blotting large 2-D gels. Because the plate electrodes are in direct contact with the filter papers, the field strength across the gel is maximized offering efficient and fast (15-45 minutes) transfer. Most semi-dry transfer methods use more than one buffer system to achieve efficient transfer but have a lower buffering capacity, making this method inappropriate for prolonged transfers. Semi-dry transfer is typically run at constant current; voltage normally increases during the transfer period.

For semi-dry transfer systems, the filter papers and membrane must be cut to the same size as the gel so the current is forced to flow through the gel. Otherwise, the current will short-circuit through overlapping filter paper around the edges of the gel. In both wet and semi-dry transfer, extra caution should be taken to prevent the introduction of air bubbles between the filter paper, gel, and membrane. Bubbles prevent transfer and cause “bald spots” (i.e., areas of non-transfer) on the blot.

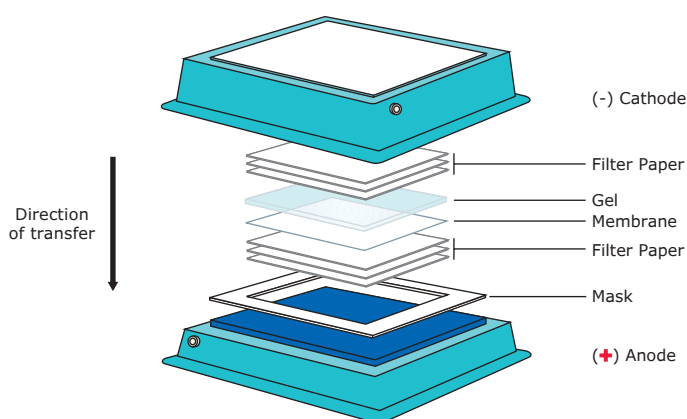


Figure 7.
Semi-dry transfer system.

Transfer Buffers

Transfer buffers must be conductive to provide electrical continuity between the electrodes. The transfer buffer must also provide a chemical environment that maintains protein solubility without preventing protein adsorption to the membrane during transfer. Additionally, most buffers undergo Joule heating during transfer. For this reason, many tank transfer systems are equipped with built-in cooling coils. The tanks can also be placed in a cold room, and the buffer can be chilled before use. In semi-dry transfer systems, the electrode plates serve as heat sinks. Their heat dissipation capacity is limited, and semi-dry systems are not normally used for prolonged transfers.

Traditional transfer buffers consist of a buffering system and methanol. Towbin buffer¹⁷, a Tris-glycine buffer, is commonly used in tank transfer because its pH (8.3) is higher than the isoelectric point (pI) of most proteins. Separated proteins (on a gel) must retain their net negative charge to migrate towards the anode and transfer to the membrane. As the buffer is mixed in the tank during transfer, ion distribution remains relatively constant.

Semi-dry systems can be run using either a single- or a three-buffer system.¹⁸ Three buffers are used because semi-dry transfer is an isotachophoretic process, in which proteins are mobilized between a leading ion and a trailing ion.¹⁹ In some cases, a three-buffer system provides better quantitative transfer. The three buffers are:

- Anode buffer I: 0.3 M Tris at pH 10.4
- Anode buffer II: 25 mM Tris at pH 10.4
- Cathode buffer: 25 mM Tris and 40 mM ϵ -aminocaproic acid at pH 9.4

Anode buffer I neutralizes excess protons generated on the surface of the anode plate. Anode buffer II contains Tris at the same pH as anode buffer I, but at a reduced concentration of 25 mM. The cathode buffer contains ϵ -aminocaproic acid, which serves as the trailing ion during transfer and is depleted from the cathode buffer as it migrates through the gel toward the anode. Review manufacturer’s recommendations for single buffers in your semi-dry system.

Although the buffer systems defined above are suitable for the majority of protein transfers, the literature contains many variations suited to specific applications. One of the most significant variations was the recommendation of 10 mM CAPS buffer at pH 11 for protein sequencing applications.⁶ The glycine in Towbin buffer caused high backgrounds in automated protein sequencers employing Edman chemistry. Changing the transfer buffer composition significantly reduced this artifact. Modifications to the buffer strength or composition should be made with care to ensure the transfer unit does not experience excessive heating.

Functions of Methanol in Transfer Buffer

Adding methanol to transfer buffers has two major functions:

- Stabilizes the dimensions of the gel
- Strips complexed SDS from the protein molecules

As a hydrogel, polyacrylamide has the capacity to absorb water. When placed in pure water, the gel's size increases (swells) in all dimensions. The degree of swelling depends on the concentration of acrylamide used in the gel. High-concentration gels expand more than low-concentration gels. Gradient gels highlight this effect quite dramatically with the more concentrated zone expanding, leading to a trapezoidal-shaped gel. Methanol concentrations of 10 to 20% minimize gel swelling, and achieve dimensional stability fairly rapidly. Lower methanol concentrations, require additional equilibration time for dimensional stability. If dimensional changes occur during transfer, protein resolution may be lost. For high MW proteins with limited solubility in methanol, elimination of the methanol can result in a significant increase in protein transfer efficiency, but this may necessitate a longer equilibration time to ensure dimensional stability.

Methanol is also critical for the protein transfer from gels containing SDS, as methanol helps to strip complexed SDS from protein molecules.²⁰ Although SDS is necessary for the resolution of individual proteins on the gel, it can be extremely detrimental to effective blotting. By imparting a high negative charge density to the protein molecule, SDS causes the protein molecule to move very rapidly through the membrane, reducing residence within the membrane pore structure and minimizing the opportunity for molecular interaction. Additionally, by coating the protein molecule, the SDS limits the ability of the protein to make molecular contact with the membrane. These effects increase as the MW of the protein decreases. By stripping off SDS, methanol reduces both effects and increases the probability that a protein molecule will bind to the membrane.

Factors Affecting Successful Protein Transfer

Presence of SDS

Some studies have suggested that proteins, in the context of an SDS-PAGE gel, bind to SDS molecules and are therefore unable to bind efficiently to the transfer membrane. Specifically, when the transfer of BSA was monitored over two hours in a standard tank transfer system, the data suggested that within a single protein band there is more than one population of molecules transferring from the gel (Figure 8). About 90% of the BSA eluted from the gel during the first 60 minutes, with an additional 7% eluting in the last 60 minutes. During the first 15 minutes, part of the eluted BSA adsorbed to the Immobilon®-P transfer membrane, while the remainder passed through and adsorbed to the backup Immobilon®-P^{SQ} transfer membrane. BSA that eluted after 15 minutes adsorbed almost exclusively to the sheet of Immobilon®-P transfer membrane.

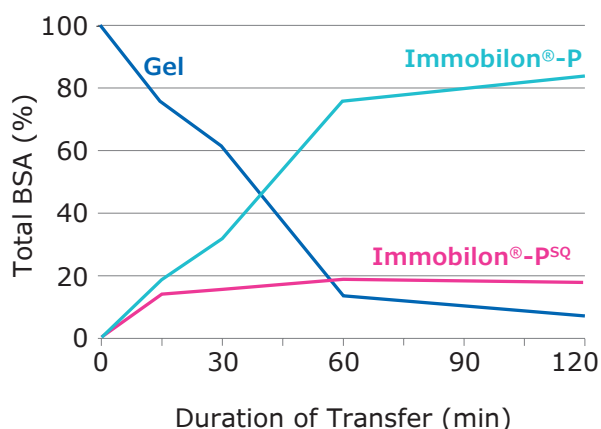


Figure 8.

Electrotransfer of BSA. 25 picomoles of ¹²⁵I-labeled BSA were resolved by SDS-PAGE on a 10–20% gradient gel. After equilibration for 5 minutes, the BSA was transferred to Immobilon®-P transfer membrane, backed up with Immobilon®-P^{SQ} transfer membrane, in a tank transfer system using 25 mM Tris, 192 mM glycine, and 10% methanol, as the transfer buffer. The system was run at 8 V/cm interelectrode distance. At 15, 30, 60, and 120 minutes, a gel/membrane cassette was removed and stained. The BSA bands were excised and counted.

The simplest interpretation is that BSA bound to residual SDS eluted from the gel to rapidly adsorb to Immobilon®-P transfer membrane. BSA that eluted more slowly was able to adsorb to the Immobilon®-P transfer membrane.

Although omitting SDS is typical, there are instances where addition of low amounts of SDS to the transfer buffer may be advantageous. One instance is when the proteins to be transferred have low solubility in the absence of SDS. Proteins associated with or integral to cellular membranes may be very hydrophobic and can precipitate within the polyacrylamide gel if SDS is removed. High MW proteins also may exhibit solubility

problems in the absence of SDS, especially after being exposed to the denaturing conditions of the gel sample buffer and the methanol used in the transfer buffer. Supplementation of the transfer buffer with SDS can help maintain sufficient solubility to permit transfer from the gel.²¹⁻²³ The SDS concentration in the transfer buffer should not exceed 0.05%, and sufficient equilibration time should be allowed to remove excess SDS from the gel.

Other methods employed to improve the transfer efficiency of high molecular weight proteins include prolonged blotting time, up to 21 hours²⁴, or the use of a composite agarose-polyacrylamide gel containing SDS and urea.²⁵

Current and Transfer Time

Appropriate current and transfer times are critical for successful blotting. Insufficient current and/or time will result in incomplete transfer. Conversely, if the current is too high, the protein molecules may migrate through the membrane too fast to be adsorbed. This can be a significant problem for smaller proteins. Blotting systems typically come with recommendations for current and transfer time.

Optimization may still be required depending on the percentage of acrylamide, buffer composition, and the MW of the protein of interest. Generally, long transfer times are best suited for tank systems, which normally require cooling of the unit and internal recirculation of the transfer buffer. In semi-dry transfer, however, prolonged blotting may result in buffer depletion, overheating and gel drying. If too much drying occurs, the unit can be damaged by electrical arcing between the electrode plates.

Transfer Buffer pH

The pH of the transfer buffer is another important factor. If a protein has an isoelectric point equal to the buffer pH, transfer will not occur. To alleviate this problem, higher pH buffers such as CAPS or lower pH buffers such as acetic acid solutions can be used.

Equilibration Time

In the early days of protein blotting (late 1970s, early 1980s), most protocols called for equilibration of the gel for 30 minutes prior to blotting. Gel sizes of 5 inches or more and thicknesses >1 mm required extended equilibration time to stabilize the gel size. Dimensional equilibrium can be reached in standard mini-gels within 5 to 10 minutes, but the kinetics of SDS stripping are significantly slower, so a minimum equilibration time of 15 minutes is recommended for most mini-gels.

Note: For samples containing small peptides, the rapid migration of peptides can occur without electrical force. In this instance, equilibration of the gel in transfer buffer should be limited to less than 10 minutes.

The running buffer in SDS-PAGE systems is supplemented with SDS, which runs into the gel behind the bromophenol blue tracking dye. Since most gels are run until the tracking dye is at the bottom of the gel, all the excess SDS remains in the gel. If the SDS isn't allowed to diffuse out of the gel prior to blotting, it will interfere with protein adsorption. Equilibration times can be extended up to 30 minutes, and a sufficient volume of transfer buffer should be used to dilute the SDS to a minimal level.

The effect of equilibration time on electrotransfer of BSA is shown in Figure 9. In this study, radioactively labeled BSA was resolved by SDS-PAGE, and the gels were equilibrated in transfer buffer for periods ranging from 0 to 30 minutes. Protein was transferred to Immobilon®-P transfer membrane backed up with a piece of Immobilon®-P^{SQ} transfer membrane to adsorb any BSA not retained by the Immobilon®-P transfer membrane. At the end of the transfer period, the BSA in the gel, on the primary blot (Immobilon®-P transfer membrane) and on the backup blot (Immobilon®-P^{SQ} transfer membrane) was quantified. Retention improved to 90% when the duration of the equilibration period was increased to 30 minutes. Other proteins have been found to behave similarly.

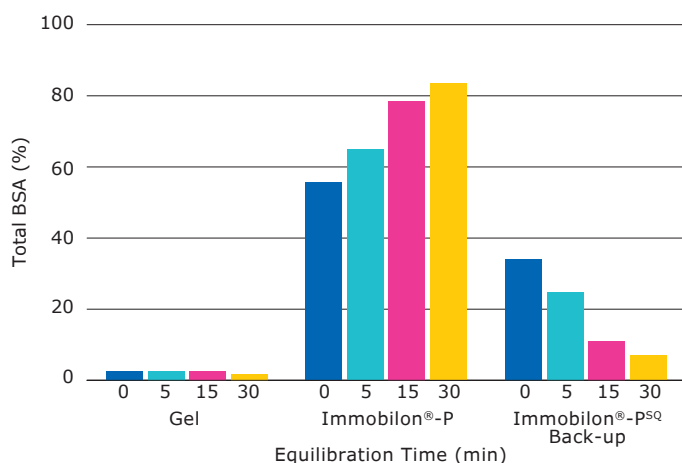


Figure 9.

Effect of equilibration time on electrotransfer of BSA to Immobilon®-P transfer membrane. ¹²⁵I-labeled BSA was resolved by SDS-PAGE on a 10–20% gradient gel. After equilibration for the times noted, the BSA was transferred to Immobilon®-P transfer membrane, backed up with Immobilon®-P^{SQ} transfer membrane, in a tank transfer system using 25 mM Tris, 192 mM glycine, and 10% methanol, as the transfer buffer. The system was run at 8 V/cm interelectrode distance. At the end of the 2-hour transfer period, the gel and membranes were stained. The BSA bands were excised and counted.

Developing a New Transfer Protocol

Although the previous section suggests that the selection of buffers and transfer conditions can be very complex, the tank transfer system defined by Towbin et al.¹⁷ and the semi-dry transfer system defined by Kyhse-Anderson¹⁸ work well for most protein samples. Both represent excellent starting points. If they prove less than optimal for a particular protein transfer conditions can be tailored to fit the biochemical peculiarities of the protein. An interesting optimization strategy for the efficient transfer of proteins over an MW range from 8,000 to > 400,000 kDa was demonstrated by Otter et al.²². The transfer buffer was supplemented with 0.01% SDS to maintain the solubility of high MW proteins and 20% methanol to enhance adsorption. The electrical field was applied in two phases. The first hour of transfer was at a low current density to slow the migration rate of low MW proteins and increase their residence time in the membrane. This was followed by a prolonged period at high current density to elute high MW proteins.

When developing a new transfer protocol or working with a new sample type, the gel should be stained to verify that all of the proteins have completely eluted from the gel. It is also highly recommended to have a lane with prestained markers in each gel to monitor the transfer efficiency. Some proteins have limited solubility in typical transfer buffers, requiring modification of the buffer chemistry to prevent precipitation. Other proteins, such as histones and ribosomal proteins, are positively charged in standard transfer buffers and will migrate toward the cathode. These proteins can be successfully transferred by placing a sheet of Immobilon®-P membrane on the cathode side of the gel. Staining the membrane after the transfer can also be helpful to ensure that the target protein is on the blot.

Another method to monitor protein transfer is to stain SDS-PAGE gels prior to electroblotting.²⁶ In this method, the gels are stained with Coomassie® Brilliant Blue dye either after, or during electrophoresis. The transferred proteins remain stained during immunodetection, providing a set of background markers for protein location and size determination.

Preparing the Membrane for Protein Identification

For Staining and Immunodetection

PVDF membranes should be washed with deionized water to remove any gel debris. The blot can then be incubated with the blocking solution.

For Rapid Immunodetection and Visualization by Transillumination

After the transfer is complete, PVDF membranes should be completely dried before continuing on to staining or immunodetection procedures. Drying enhances the adsorption of the proteins to the PVDF polymer, helping to minimize desorption during subsequent analyses. As the blotted membrane dries, it becomes opaque. This optical change is a surface phenomenon that can mask retention of water within the depth of the pores. The membrane should be dried for the recommended period to ensure that all liquid has evaporated from within the membrane's pore structure (see Protocol 1.7. Membrane Drying Methods).

Storage

PVDF membranes can be stored dry for long periods of time after protein transfer with no ill effects to the membrane or the proteins (up to two weeks at 4°C; up to two months at -20°C; for longer periods at -70°C). Some proteins, however, may be sensitive to chemical changes (e.g., oxidation, deamidation, hydrolysis) upon prolonged storage in uncontrolled environments. Long-term storage at low temperature is recommended. Prior to further analysis, dried membranes must be wetted by soaking in 100% methanol.

Protein Visualization

Transillumination

Transillumination (Figure 10, Page 17) is a visualization technique unique to PVDF membranes and was first described for Immobilon®-P transfer membran.²⁷ This technique takes advantage of a characteristic unique to PVDF membranes: areas of PVDF coated with transferred protein are capable of

wetting out in 20% methanol while the surrounding areas of PVDF are not. In the areas where the PVDF wets, it becomes optically transparent, allowing visualization of protein bands using backlighting and photographic archiving. The process is fully reversible by evaporation. Further denaturation of the proteins is unlikely as the proteins had been previously exposed to methanol during blotting. Even though this technique

does not allow for the visualization of low-abundance proteins, it can be used to assess overall transfer efficiency and the suitability of the blot for further analysis.

Staining

Staining (Figure 10) is a simple technique to make proteins visible on a blot. Staining can be used to:

- Verify proteins have transferred onto the membrane
- Determine if lanes were loaded equally
- Evaluate overall transfer efficiency, especially for a new buffer system or protein
- Identify and excise bands for peptide sequencing

Many types of stains are available, (Table 5) including organic dyes (Ponceau-S red, amido black, fast green and Coomassie® Brilliant Blue dye), fluorescent dyes (fluorescamine, coumarin) and colloidal particles (gold, silver, copper, iron and India ink).³ Dyes can be separated into two groups: reversible and irreversible stains.

Reversible Stains

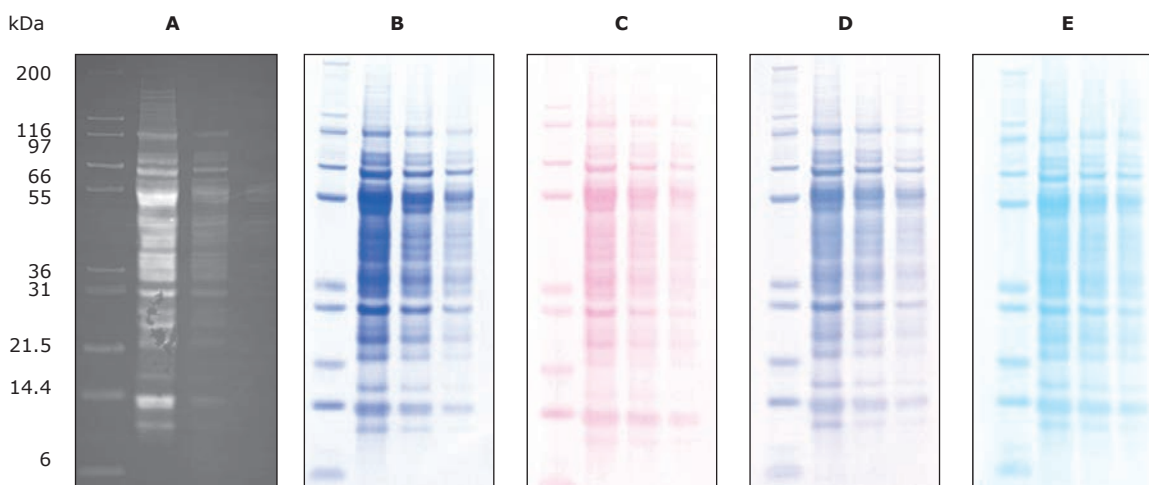


Figure 10.

Calf liver proteins are visualized after electroblotting to Immobilon®-P membranes:

(A) Transillumination, (B) Coomassie® Brilliant Blue dye, (C) Ponceau-S red, (D) Amido black and (E) CPTS total protein stains. Left to right, molecular weight standards and 12.2 µg, 6.1 µg, 3.1 µg of the lysate per lane.

Table 5. Common stains used in Western blotting

	Detection Reagent (protein per spot)	Approximate Sensitivity	Reference
Reversible	Ponceau-S red	5 µg	30
	Fast green FC	5 µg	30
	CPTS	1 µg	31
	Sypro® Ruby	1–2 ng	29
	Sypro® Rose	1–2 ng	29
Irreversible	Amido black 10B	1 µg	30
	Coomassie® Brilliant Blue R-250	500 ng	30
	India ink	100 ng	30
	Colloidal gold	4 ng	30

Reversible stains allow assessment of the blot and then can be washed from the membrane. These will not interfere with subsequent immunodetection or other analysis of the proteins on the blot. The most commonly used reversible protein stain is Ponceau-S red.

The major drawback of reversible stains is that they are less sensitive than irreversible stains. Since the staining pattern of more abundant proteins in a blot is generally a good indicator of how well low-abundance proteins transferred, this drawback can be minimized in most cases.

Alternatively, fluorescent blot stains are highly sensitive and compatible with downstream immunodetection, Edman-based sequencing, and mass spectrometry²⁸. Sypro® Ruby and Sypro® Rose protein blot stains (Thermo Fisher Scientific) can be used prior to chromogenic, fluorogenic, or chemiluminescent immunostaining procedures and provide sensitivity of about 1–2 ng/band.²⁹

Irreversible Stains

Irreversible stains generally exhibit ideal sensitivity, but can interfere with or prevent further protein analysis. Examples of irreversible stains are amido black and Coomassie® Brilliant Blue dye.

4. Protein Identification

Immunodetection

Immunodetection uses a specific antibody to detect and localize a protein blotted onto the membrane (Figure 11). The specificity of antigen-antibody binding permits the identification of a single protein in a complex sample.

When developing new immunodetection protocols, all components and their interactions must be considered. Antibody concentrations, buffer compositions, blocking agents, and incubation times must be tested empirically to determine optimal conditions. Water quality is important in all steps—small impurities can cause big problems. For instance, the enzyme activity of horseradish peroxidase is inhibited by pyrogens, a common contaminant of even high-purity water, and azide, a common preservative in antibody solutions. The quality of the blocking agents must also be considered relative to consistency and contaminants.

Standard vs. Rapid Immunodetection Procedures

There are two types of protocols for immunodetection: standard and rapid.

Standard immunodetection methods include the following steps:

1. Block unoccupied membrane sites to prevent nonspecific antibody binding.
2. Incubate the membrane with primary antibody, which binds to the protein of interest.
3. Wash to remove any unbound primary antibody.
4. Incubate the membrane with a conjugated secondary antibody, which binds the first antibody.
5. Wash to remove any unbound secondary antibody.
6. Incubate the membrane with a substrate that reacts with the conjugated secondary antibody to reveal the protein location.

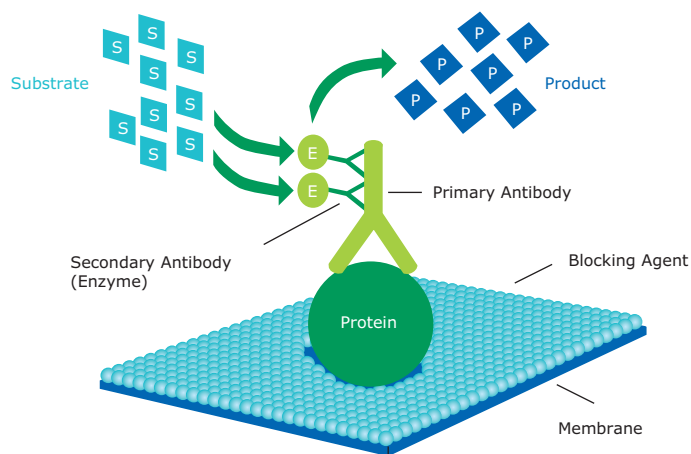


Figure 11.
Membrane-based immunodetection.

Rapid immunodetection eliminates the blocking step and reduces the time necessary for washing and incubation steps. The rapid immunodetection method works well to quickly visualize higher-abundance proteins. Standard immunodetection, however, offers higher sensitivity and requires less optimization for new sample types. Procedures for both standard and rapid immunodetection methods are outlined in Protocol 3. Table 6 compares the times required to perform the two protocols.

Table 6. Standard vs. rapid immunodetection

Step	Standard Immunodetection	Rapid Immunodetection
1. Block the membrane	1 hr	None
2. Incubate with primary antibody	1 hr	1 hr
3. Wash the membrane	3 x 10 min	3 x 5 min
4. Incubate with secondary antibody	1 hr	30 min
5. Wash the membrane	3 x 10 min	3 x 5 min
6. Add substrate	5 min	5 min
Total time	4 hr 5 min	2 hr 5 min

Factors Influencing Immunodetection

Understanding the concepts presented in the following sections on immunodetection will help in optimizing protocols for specific samples.

Buffers

The two most commonly used buffers are phosphate-buffered saline (PBS) and Tris-buffered saline (TBS). Many variations on the compositions of these buffers have been published. The key requirement of the buffer is that it should help preserve the biological activity of the antibodies. Thus, the ionic strength and pH should be fairly close to physiological conditions. PBS formulations with 10 mM total phosphate concentration work well with a wide array of antibodies and detection substrates.

Blocking

For meaningful results, the antibodies must bind only to the protein of interest and not to the membrane. Nonspecific binding (NSB) of antibodies can be reduced by blocking the unoccupied membrane sites with an inert protein, non-ionic detergent, or a protein-free blocker such as Millipore® Block noise-canceling reagents. The blocking agent should have a greater affinity for the membrane than the antibodies and should fill all unoccupied membrane binding sites without displacing the target protein. The most common blocking agents are bovine serum albumin (BSA, 0.2–5.0%), non-fat milk, casein, gelatin, dilute solutions of Tween®-20 detergent (0.05–0.1%), and Millipore® Block reagents. Tween®-20 detergent was also shown to have a renaturing effect on antigens, resulting in improved recognition by specific antibodies.^{32,33} Other detergents, such as Triton® X-100 detergent, SDS, and NP-40, are sometimes used, but can be too harsh and disrupt interaction between proteins. The blocking agent is usually dissolved in PBS or TBS buffers.

There are risks associated with blocking: a poorly selected blocking agent or excessive blocking can displace or obscure the protein of interest. Therefore, the blocking agent can be critical to successful immunodetection. For example, dry milk cannot be used with biotinylated or concanavalin-labeled antibodies, as milk contains both glycoproteins and biotin. Detection of phosphorylated proteins can be compromised if a

During incubation, the container holding the membrane should be gently agitated. A sufficient volume of buffer should be used to cover the membrane completely so that it is floating freely in the buffer. If more than one blot is placed in a container, insufficient buffer volume will cause the blots to stick together. This will limit the accessibility of the incubation solutions and can cause a variety of artifacts including high backgrounds, weak signals, and uneven sensitivity.

blocking agent containing phosphatases is used. The phosphorylated proteins on the blot could become dephosphorylated by these enzymes. It has been shown that addition of phosphatase inhibitors to the blocking solution increases the signal with phospho-specific antibody.³⁴ Finally, blocking agents are not universally suitable for all antigen-antibody combinations.

Compatibility between the blocking agent and detection reagent can be determined easily using the Dot Blotting Method described in Protocol 1.5. The blocking solution gets spotted onto a blank PVDF membrane that has been wet in methanol and equilibrated in TBS. Detection reagents are then added to the blot, the blot is incubated for 5 minutes and then exposed to X-ray film or imaged using a digital imager. Appearance of a dark spot indicates that the blocking reagent is incompatible with the detection reagent. It is important to remember that Immobilon®-P^{SQ} transfer membrane, with its higher surface area and smaller pore size than Immobilon®-P transfer membrane, binds more protein. If Immobilon®-P^{SQ} transfer membrane is substituted directly for Immobilon®-P transfer membrane in a standard Western blotting procedure, there may be insufficient blocking reagent to saturate the membrane surface. Additional washing steps may be required to reduce the background. Blocking can be conveniently optimized using the Dot Blotting Method.

Antibodies

After blocking, the blot is incubated with one or more antibodies. The first antibody binds to the target protein, and a secondary antibody binds to the first. The secondary antibody is conjugated to an enzyme or dye that is used to indicate the location of the protein.

Although the primary antibody may be labeled directly, using a secondary antibody has distinct advantages. First, more than one molecule of the secondary antibody may be able to bind to a single molecule of the primary antibody, resulting in signal amplification. Second, a labeled secondary antibody (enzyme-antibody conjugate) can be used for a large number of primary antibodies of different specificities, thereby eliminating the need to label numerous primary antibodies.

Either polyclonal or monoclonal primary antibodies are used. Polyclonal antibodies usually come in the form of antiserum or affinity-purified antibody. Monoclonal antibodies are expressed in ascites fluid or tissue culture fluid and can be used directly or as an affinity-purified preparation. It is important to remember that a denatured protein may not be recognized by an antibody raised to the native antigen. In some cases, a nondenaturing gel may be required for production of the blot. Antibodies are diluted in buffer and blocking solution to prevent nonspecific membrane binding. The antibody diluent also normally contains trace amounts of Tween®-20 or another detergent to prevent nonspecific aggregation of the antibodies. Many published protocols for chemiluminescence require 0.1% (v/v) Tween®-20 in the blocking solution and antibody diluent. It is important to recognize that concentrations above 0.05% (v/v) have the potential to wash blotted proteins from the membrane.

Elevating the concentration of Tween®-20 detergent can often reduce the background. Often, a simpler and more cost-effective strategy is to reduce the concentration of the antibodies, especially the secondary antibody (Figure 12). In addition to being specific for the protein of interest, the antibodies must not cross-react with components of the blocking buffer and should be relatively pure. Impurities in the form of other proteins or aggregates can result in nonspecific binding and increased background.

Immunodetection is an extremely sensitive method. In order to achieve a high signal-to-noise ratio and maximum sensitivity, the concentrations of primary and secondary antibodies should be optimized for each case. Generally, nonspecific signal can be reduced by higher dilution of the primary antibody or decreased protein load on the original gel. High overall background can be minimized by higher dilution of the enzyme-conjugated secondary antibody.

Optimal concentration of both primary and secondary antibodies also depends on the sensitivity of the detection reagents. Up to twenty times less antibody is required for high-sensitivity reagents (detection at femtogram level) compared to low-sensitivity reagents (detection at the picogram level). An example of antibody concentration optimization based on the detection substrate is shown in Figure 13. When using the most sensitive detection reagent (Luminata™ Forte reagent), excessive amounts of antibody led to high nonspecific binding and increased overall background (top left). In this case, increasing the primary antibody dilution from 1:1,000 to 1:10,000 improved the signal-to-noise ratio. With progressively less-sensitive detection reagents (middle and bottom rows, respectively), antibodies must be more concentrated to yield the same signal.

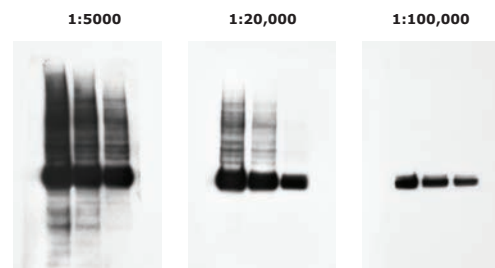


Figure 12.

Optimization of secondary antibody dilution for immunodetection of ERK 1 with Immobilon® Western Chemiluminescent HRP Substrate. Two-fold dilutions of rat liver lysate samples were applied to each gel. Electroblotted proteins were probed with rabbit anti-ERK 1 antibody (1:1,000 dilution) and HRP-conjugated goat anti-rabbit IgG (1:5,000, 1:20,000 and 1:100,000 dilutions, left to right).

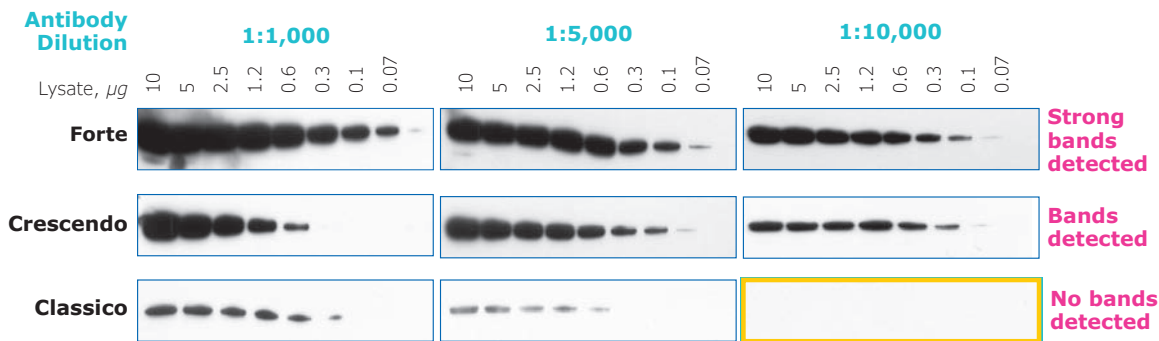


Figure 13.

The more sensitive the detection reagent, the less antibody is needed. Immunoblots containing the indicated amounts of A431 lysate were probed with different concentrations of anti-GAPDH antibody (Cat. No. MAB374) indicated in the top row, followed by an appropriate secondary antibody. Bands were visualized using the indicated Luminata™ HRP substrate and exposed to x-ray film for 5 minutes. Of the three detection reagents shown, Luminata™ Forte reagent is the most sensitive, followed by Luminata™ Crescendo and then Luminata™ Classico reagents.

Using higher-sensitivity HRP substrates was advantageous in three respects:

- Better results:** It produced stronger bands for a more quantitative blot (compare the increase in band intensities for Luminata™ Crescendo & Forte substrates at 1:10,000 dilution)
- Faster:** It took only 10 minutes to wash blot and add a new substrate relative to the 2.5 hours required to repeat antibody incubations
- Cheaper:** The HRP substrates are much cheaper than the price of antibodies

Washing

Washing the blot removes any unbound antibodies from the membrane that could cause high background and poor detection. A dilute solution of Tween®-20 (0.05% v/v) in PBS or TBS buffer is commonly used, especially when the antibody preparations are comparatively crude or used at high concentrations. As mentioned previously, concentrated detergent solutions may cause unwanted protein elution from the membrane. For highly purified antibodies, buffer alone is often sufficient for washing.

The amount of washing required is best determined experimentally. Too little washing will yield excessive background, while overwashing may elute the antibodies and reduce the signal. It is recommended that washing be performed a minimum of three times for 5 minutes each.

Persistent background can be reduced by adding up to 0.5 M sodium chloride and up to 0.2% SDS to the TBS wash buffer and extending wash time to 2 hours.

Double Blotting

An innovative method to eliminate nonspecific binding in Western blots was developed by Dr. Françoise Lasne of the Laboratoire National de Dépistage du Dopage (National Anti-Doping Laboratory).^{35,36} While studying recombinant human erythropoietin (EPO), Dr. Lasne found that recombinant and naturally occurring EPO have different isoelectric points (pI). Recombinant EPO has a pI of 4.42–5.11, while natural EPO has a more acidic pI of 3.92–4.42. However, when urine samples are blotted, the very high nonspecific binding (NSB) of the secondary antibodies makes it difficult to distinguish between recombinant and natural EPO. To eliminate the NSB, Dr. Lasne used “double blotting”. After the primary antibody is bound to the blotted protein, the antibodies are transferred to a second Immobilon®-P membrane under acidic conditions. The primary antibody desorbs from its corresponding antigen and transfers through an intermediate onto the second (double-blot) membrane. When the double-blot membrane is probed with the secondary antibody, no other proteins are present to bind nonspecifically, thus eliminating the background problem (see Protocol 1.6. Double Blotting Procedure). Double blotting has also been used in the detection of transthyretin and may be a useful method in detecting other low-concentration serum or urine proteins.

Detection Substrates

Modern immunodetection methods are based on enzyme-linked detection, utilizing secondary antibodies covalently bound to enzymes such as horseradish peroxidase (HRP) or alkaline phosphatase (AP). The conjugated enzyme catalyzes the degradation

of specific substrates, resulting in signal generation. Three types of substrates are commonly used: chromogenic, chemiluminescent, and chemifluorescent, as well as detection with fluorophore-labeled secondary antibodies. Immobilon® PVDF membranes have been tested for compatibility with all commercially available chromogenic and chemiluminescent substrates.

Chromogenic Detection

Chromogenic detection (Figure 14) uses a conjugated enzyme to catalyze a reaction resulting in the deposition of an insoluble colored precipitate, for example, the insoluble blue compound obtained through the interaction of 5-bromo-4-chloro-3-indolylphosphate (BCIP) and nitroblue tetrazolium salt (NBT).³⁷ This technique is easy to perform and requires no special equipment for analysis. However, the following should be kept in mind:

- Sensitivity of chromogenic detection is typically at least one order of magnitude lower than chemiluminescent reagents
- Production of the precipitate can interfere with enzyme activity and limit sensitivity
- The precipitate is difficult to strip from membrane, limiting reuse of the blot for detection of other proteins

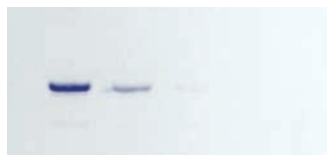


Figure 14.

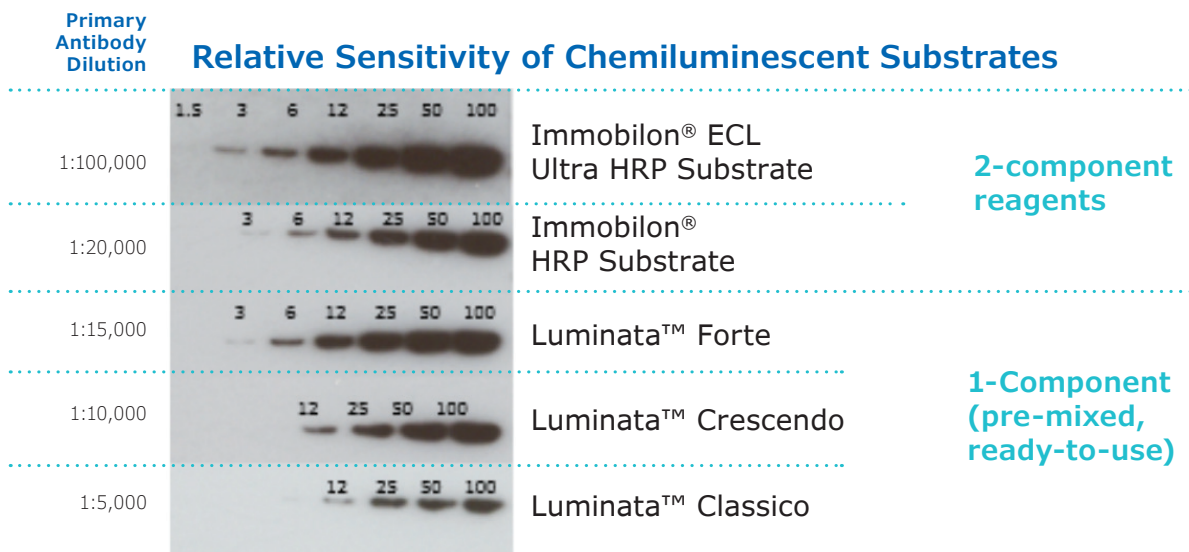
Immunodetection of transferrin in human serum with chromogenic substrate BCIP/NBT (KPL). Left to right, 5 µL human serum dilutions 1:1,000, 1:5,000, 1:25,000, 1:125,000. Electroblotted proteins were probed with goat anti-human transferrin (1:10,000 dilution) and AP-conjugated rabbit anti-goat IgG (1:30,000 dilution).

Chemiluminescent Detection

Chemiluminescent detection uses a conjugated enzyme to catalyze a reaction that results in the production of visible light. Some chemiluminescent systems are based on the formation of peroxides by horseradish peroxidase; other systems use 1,2-dioxetane substrates and the enzyme alkaline phosphatase.³⁸ This technique has the speed and safety of chromogenic detection with sensitivity levels comparable to radioisotopic detection. Detection is achieved by either exposing the blot to X-ray film or acquiring the image directly in a chemiluminescence-compatible digital imaging system, usually equipped with super cooled CCD cameras to avoid electronic noise. Reprobing is possible with chemiluminescent substrates.

Researchers can obtain varying levels of detection sensitivity with chemiluminescent substrates. Traditional or low-sensitivity substrates allow protein detection at the picogram level. While these substrates may be appropriate for routine applications, they cannot detect low abundance proteins. High-sensitivity substrates allow visualization of proteins at the femtogram level. However, use of these powerful substrates often requires primary and secondary antibody concentration optimization (Figure 13, Page 22). When switching from a low-sensitivity to a high-sensitivity substrate, it is recommended that antibody dilutions be increased to avoid excessive background and appearance of nonspecific bands.

Reagents for ECL immunodetection can be prepared using p-iodophenol (PIP) and luminol.³⁹ PIP is needed to enhance the visible light reaction by acting as a co-factor for peroxidase activity toward luminol. When phenolic enhancers are used in combination with HRP, the level of light increases about 100-fold.⁴⁰ These homemade reagents are cited to produce excellent results; however, the purity of the luminol and PIP is critical.³⁹



Fluorescent Detection

Fluorescent detection employs either a fluorophore-conjugated antibody or fluorogenic substrates that fluoresce at the site of enzyme activity (chemifluorescence). One advantage of this method is that the fluorescent signal is stable for long periods of time, and blots can be archived and re-imaged. In addition, the wide variety of fluorophores makes it possible to simultaneously detect multiple protein targets in a single sample (multiplex detection). Until recently, fluorescent detection in Western blotting was limited by the high-fluorescent background of most blotting membranes. Immobilon®-FL transfer membrane exhibits low background fluorescence across a wide range of excitation/emission wavelengths compared to other blotting membranes (Figure 15). The membrane is ideal for any application involving fluorescence-based immunodetection, including chemifluorescent substrates and multiplexing (Figure 16). In addition, Immobilon®-FL membrane can be used for standard chemiluminescent or chromogenic detection.

Reprobing Immobilon® PVDF Transfer Membranes

A single blot can be sequentially analyzed with multiple antibodies by stripping the first antibody from the blot and incubating with another. This may be especially useful for co-localization experiments and method optimization or when sample amount is limited (see Protocol 4.1 Membrane Stripping Protocols).

The stripping process disrupts the antigen antibody bonds and dissolves the antibody in the surrounding buffer. This is usually achieved either by a combination of detergent and heat or by exposure to low pH. Neither method removes the colored precipitates generated from chromogenic detection systems (e.g., BCIP, 4CN, DAB, and TMB). However, it is still possible to analyze the blot with an antibody specific for a different target protein.

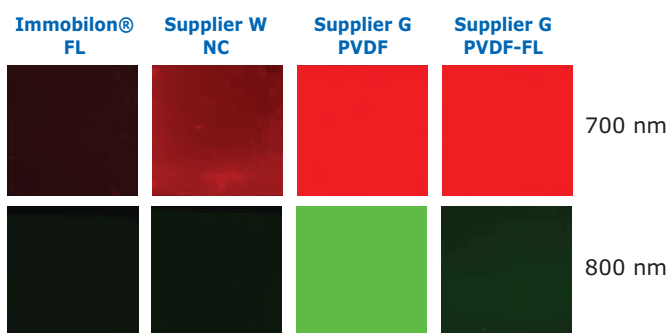


Figure 15. Comparison of background fluorescence observed in PVDF membranes from different manufacturers. Only Immobilon®-FL membrane offers low background fluorescence at both 700 and 800 nm, increasing its ability to be used in multiple excitation and emission wavelengths.

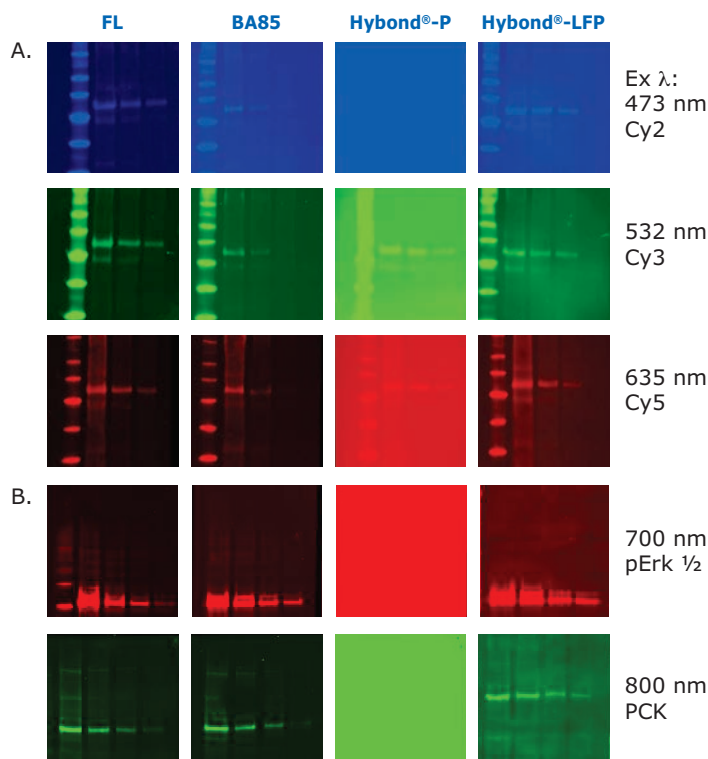


Figure 16. **A.** Two-fold dilution series of (10-2.5 µg/lane) of human brain lysate (CL302-250UG, MilliporeSigma) were resolved by SDS-PAGE and transferred onto either an Immobilon®-FL membrane, Protran® BA85 membrane, Hybond®-P membrane, or Hybond®-LFP membrane. The blots were probed with Anti-GSK3-β antibody (MAB8687, MilliporeSigma) followed by anti-rabbit IgG antibody either Cy2, Cy3, or Cy5 conjugated (AP132J, AP187C, AP187S, MilliporeSigma). The blot was scanned on Fuji FLA-5100 after drying the blot for 1 hr under vacuum. **B.** Two-fold dilution series of (10-2.5 µg/lane) of human brain lysate (CL302-250UG, MilliporeSigma) were resolved by SDS-PAGE and transferred onto either an Immobilon®-FL membrane, Protran® BA85 membrane, Hybond®-P membrane, or Hybond®-LFP membrane. The blots were probed using the SNAP i.d.® system with Anti-GSK3-β antibody (MAB8687, MilliporeSigma) followed by pErk ½ or PCK. The blot was imaged using the Odyssey® Infrared Imaging System.

Mass Spectrometry

Mass Spectrometry with Immobilon® PVDF Transfer Membranes

Mass spectrometry (MS) is an analytical method used to identify proteins on a blot. After staining the proteins on a PVDF membrane with MS-compatible dye (Coomassie® Brilliant Blue dye, Amido black or Sypro® stains), the band of interest is cut out of the membrane and subjected to tryptic digestion. The peptide fragments generated are then analyzed by mass spectrometry.⁴¹⁻⁴³ Figure 17 demonstrates a MALDI-TOF spectrum obtained for an on-membrane-digested bovine protein successfully identified as catalase.

Another method to consider is protein mass spectrometry directly from the blotted Immobilon® PVDF transfer membrane. This method usually is applied for 2-D gel separated proteins. Using a parallel process, all proteins on a gel are simultaneously digested proteolytically and electrotransferred onto an Immobilon® PVDF transfer

membrane, which is then scanned for the presence of the peptides.^{42, 44, 45} Alternatively, in a method called “chemical printing”, two-dimensionally separated proteins are first transferred to PVDF membrane and visualized, and then digested by dispensing small amounts of trypsin directly onto the spots⁴⁶. In both methods, the membrane is sprayed with matrix and directly scanned by MALDI-TOF MS. Protein identification is obtained by peptide mass fingerprinting. Figure 18 shows 2-D-separated human plasma proteins transferred to Immobilon®-P^{SQ} transfer membrane and two MALDI-TOF spectra obtained for identification of membrane-immobilized proteins.

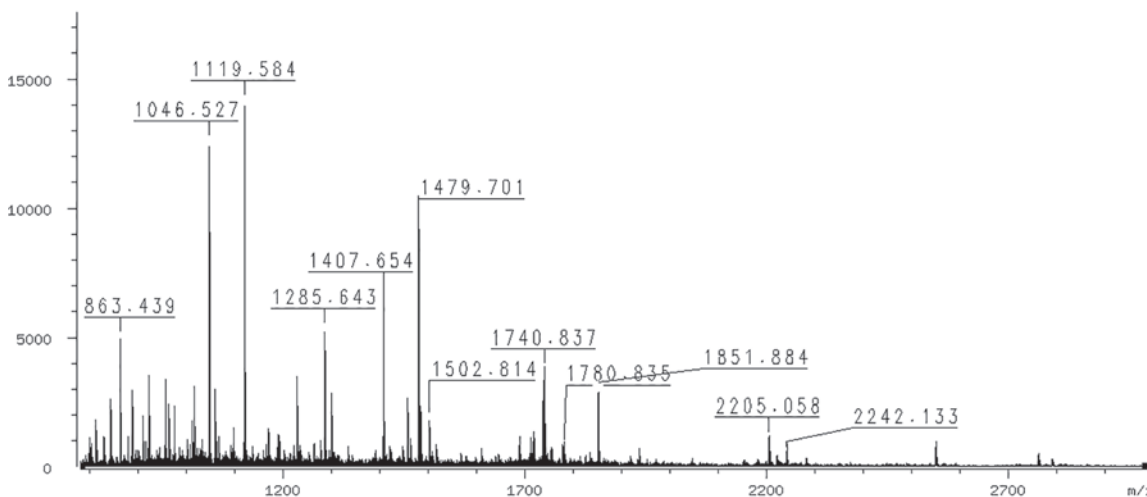


Figure 17.

MALDI-TOF spectrum of a band from Coomassie® Brilliant Blue-stained blot of liver proteins, using Protocol 5.1. On-membrane protein digestion.⁴² The protein was identified as bovine catalase, with 6.4e+06 MOWSE score and 34% coverage. Data were obtained on a Bruker® Autoflex™ mass spectrometer. The search was done using Protein Prospector.

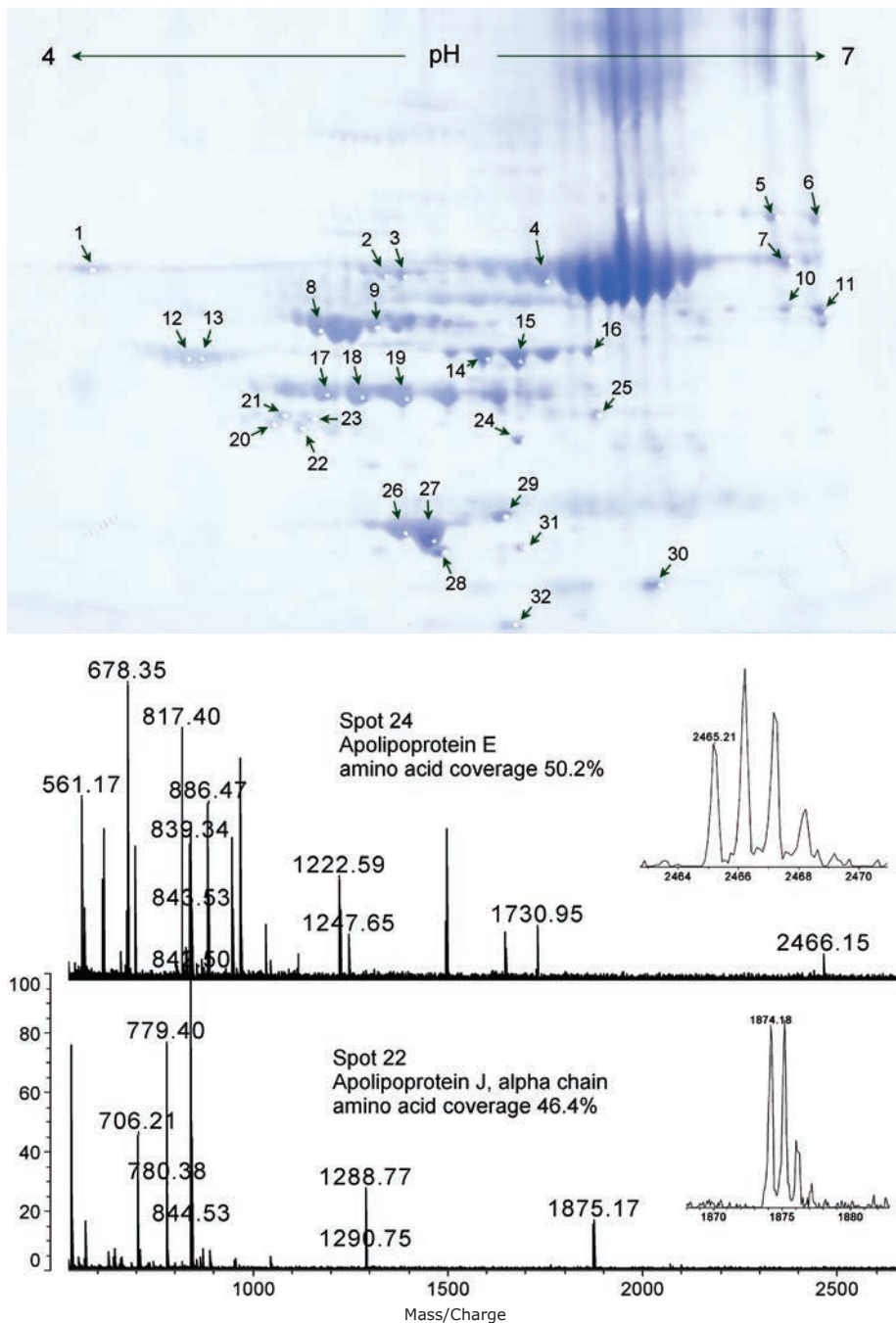


Figure 18.

Human plasma separated by 2-D electrophoresis and transferred to (left) Immobilon®-P^{5Q} transfer membrane and (right) MALDI-TOF mass spectra collected directly from the membrane surface of tryptic digested proteins. Human plasma was separated by 2-D electrophoresis, electro-blotted onto Immobilon®-P^{5Q} transfer membrane, and adhered to a MALDI target. Digestion of proteins with the endoproteinase trypsin was performed directly on the membrane surface. Nanoliters of enzyme were required for digestion and were microdispensed with drop-on-demand piezoelectric ink-jet devices onto the membrane surface using a ChIP™ (Proteome Systems and Shimadzu Corp.) instrument. The resultant peptides were analyzed with an AXIMA™-CFR (Shimadzu Corp.) MALDI-TOF MS and identified with peptide mass fingerprinting. The peptides were analyzed directly from the membrane surface, where MALDI matrix was microdispensed on top of the digested protein prior to analysis. Data courtesy of Drs. J.L. Duff, F.G. Hopwood, C.J. Hill, A.A. Gooley (Proteome Systems, Ltd., Sydney, Australia).

5. Protocols

This section of the protein blotting handbook is a compendium of the most frequently used protein blotting protocols. The methods are general enough that they can be used with all commercially available detection reagents and can be optimized using the numerous tips included in this section. These optimization tips are a product of our years of experience with transfer membranes and blotting, as well as referenced literature and feedback from our customers.

- ▶ **Tip:** Ethanol or isopropanol can be substituted for methanol in the transfer buffer.
- ▶ **Tip:** SDS in the transfer buffer (up to 0.05%) can improve transfer efficiency but may also reduce the membrane's protein retention.
- ▶ **Tip:** Immobilon® transfer membrane pre-cut sheets fit all standard mini-gel sizes and common electrophoresis systems. (Tables 3 and 4, Page 7)
- ▶ **Recommendation** – Chill transfer buffers prior to tank transfer.
- ▶ **Tip:** Thicker gels or larger proteins may require longer transfer times or increased field strength.

- ▶ **Tip:** Methanol in the transfer buffer (8 – 20%) decreases efficiency of protein elution from the gel but improves absorption to PVDF transfer membrane.
- ▶ **Tip:** Tris-glycine buffer will produce higher background in N-terminal sequencing. Use CAPS or TBE buffer for transfer if protein bonds are to be sequenced by Edman degradation.
- ▶ **Tip:** Replace 0.45 µm Immobilon®-P transfer membrane with 0.2 µm Immobilon®-P^{SQ} transfer membrane when analyzing lower molecular weight proteins.
- ▶ **Tip:** Use 0.45 µm Immobilon®-FL transfer membrane for fluorescent detection methods.

1. Protein Transfer

Protocol 1.1 Electrotransfers: Tank (Wet) Transfer

The following protocol describes the standard procedure for transferring proteins from a polyacrylamide gel (SDS-PAGE) onto an Immobilon® PVDF transfer membrane using a tank transfer system. Please review the instructions supplied with your tank transfer system for additional information.

Required Equipment and Solutions

- Polyacrylamide gel containing resolved proteins
- Immobilon® PVDF transfer membrane, cut to the same dimensions as the gel (including notched corner for orientation purposes)
- Two sheets of Immobilon® blotting filter paper or equivalent, cut to the same dimensions as the gel
- Two foam pads (compatible with the transfer tank system being used)
- Tank transfer system large enough to accommodate gel
- Methanol, Ethanol or Isopropanol
- Milli-Q® water
- Tris/glycine transfer buffer: 25 mM Tris base, 192 mM glycine, 10% (v/v) methanol, pH 8.3; or CAPS buffer: 10 mM 3-[Cyclohexylamino]-1-propanesulfonic acid (CAPS), 10% (v/v) methanol, pH 11 (adjust with NaOH)

Note: Both buffers can be prepared as 10X stock solutions and mixed with methanol prior to use.

Setup

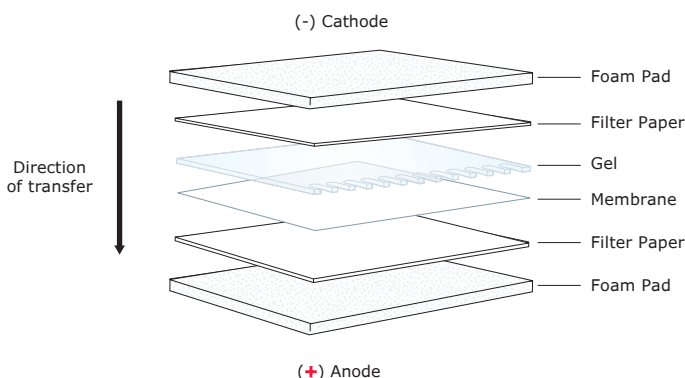
1. Prepare sufficient transfer buffer to fill the transfer tank, plus an additional 200 mL to equilibrate the gel and membrane, and wet the filter paper.
2. Remove the gel from its cassette; trim away any stacking gel and wells.
3. Immerse the gel in transfer buffer for 10 to 30 minutes (follow gel manufacturers guideline for gel incubation).
4. Soak filter papers in transfer buffer for at least 30 seconds.
5. Prepare the membrane:
 - a. Wet the membrane in methanol for 15 seconds. Membrane should uniformly change from opaque to semi-transparent.
 - b. Carefully place the membrane in Milli-Q® water and soak for 2 minutes.
 - c. Carefully place the membrane in transfer buffer and let equilibrate for at least 5 minutes.

Transfer Stack Assembly

1. Open the transfer cassette.

Important: To ensure an even transfer, remove air bubbles between layers by carefully rolling a pipette or a stirring rod over the surface of each layer in the stack. Do not apply excessive pressure to prevent damaging the membrane and gel.
2. Place a foam (fiber) pad on one side of the cassette.
3. Place one sheet of filter paper on top of the pad.
4. Place the gel on top of the filter paper.
5. Place the membrane on top of the gel.
6. Place a second sheet of filter paper on top of the stack.
7. Place a second foam pad on top of the filter paper.
8. Close the transfer cassette.

The completed transfer stack should look like this:



Protein Transfer

1. Place the transfer cassette into the electrophoresis tank so that the gel side of the cassette holder is facing the cathode (-) and the membrane side is facing the anode (+).
2. Add adequate buffer to the tank to cover the cassette holder.
3. Insert the black cathode lead (-) into the cathode jack and the red anode lead (+) into the anode jack on the transfer unit.
4. Connect the tanks anode and cathode leads to the corresponding power supply inputs.
5. If available, set up the cooling unit on the tank transfer unit according to the manufacturer's instructions.
6. Turn on the system for 1 to 2 hours at 6 to 8 V/cm inter-electrode distance. Follow the tank manufacturer's guidelines, optimizing as needed.
7. After transfer is complete, remove the transfer cassette from the tank.
8. Using forceps, carefully disassemble the transfer stack.

Protocol 1.2 Electrotransfers: Semi-dry Transfer

The following protocol describes the standard procedure for transferring proteins from a polyacrylamide gel (SDS-PAGE) onto an Immobilon® PVDF transfer membrane using a semi-dry transfer system. It is specific for semi-dry transfer devices with the anode plate serving as the base. For devices having the cathode plate as the base, consult the instruction manual for recommended buffers and transfer stack assembly. This protocol describes a three-buffer system. Single-buffer systems can also be used. Please consult manufacturer's recommendations.

- ▶ **Recommended** – Transfer proteins at constant current. If transferring at constant voltage, monitor current to make sure it doesn't exceed 0.4 amp. Start from 100 V and reduce voltage if current is too high. (Follow transfer system and gel manufactures recommendations)
- ▶ **Tip:** For semi-dry transfer systems, it is important that the filter papers and membrane are cut to the same size as the gel so that the current is forced to flow through the gel.
- ▶ **Tip:** In both types of transfer systems (tank and semi-dry), extra caution should be taken to prevent introduction of air bubbles anywhere between the filter paper, gel or membrane.
- ▶ **Tip:** Both sides of Immobilon® transfer membranes work equally well. The appearance of either side (shiny or dull) has no effect on the transfer and detection efficiency of the membrane.
- ▶ **Tip:** For samples containing small peptides, equilibration of the gel in transfer buffer should be limited to less than 10 minutes.
- ▶ **Tip:** Gels can be transferred individually or multiple gels can be transferred in a single stack.

Required Equipment and Solutions

For single transfers:

- Polyacrylamide gel containing resolved proteins
- Immobilon® PVDF transfer membrane, cut to the same dimensions as the gel (including notched corner)
- Six pieces of Whatman® 3MM filter paper or equivalent, cut to the same dimensions as the gel
- Semi-dry transfer system large enough to accommodate gel
- Anode buffer I: 0.3 M Tris, pH 10.4, 10% (v/v) methanol
- Anode buffer II: 25 mM Tris, pH 10.4, 10% (v/v) methanol
- Cathode buffer: 25 mM Tris, 40 mM 6-amino-n-caproic acid (glycine may be substituted), 10% (v/v) methanol, pH 9.4
- Methanol, 100%
- Milli-Q® water

For multiple transfers, all of the above plus the following:

- Dialysis membrane, cut to the same dimensions as the gel and wet with Milli-Q® water (The membrane should have a molecular weight exclusion small enough to retain the lowest molecular weight protein in the gel)
- Additional pieces of filter paper

Setup

1. Prepare 200 mL of each anode buffer and 400 mL of cathode buffer.
2. Remove the gel from its cassette; trim away any stacking gel.
3. Immerse the gel in 200 mL of cathode buffer for 15 minutes.
4. Soak two pieces of filter paper in anode buffer I for at least 30 seconds.
5. Soak one piece of filter paper in anode buffer II for at least 30 seconds.
6. Soak three pieces of filter paper in cathode buffer for at least 30 seconds.
7. Prepare the membrane:
 - a. Wet the membrane in methanol for 15 seconds. The membrane should uniformly change from opaque to semi-transparent.
 - b. Carefully place the membrane in Milli-Q® water and soak for 2 minutes.
 - c. Carefully place the membrane in anode buffer II and let equilibrate for at least 5 minutes.

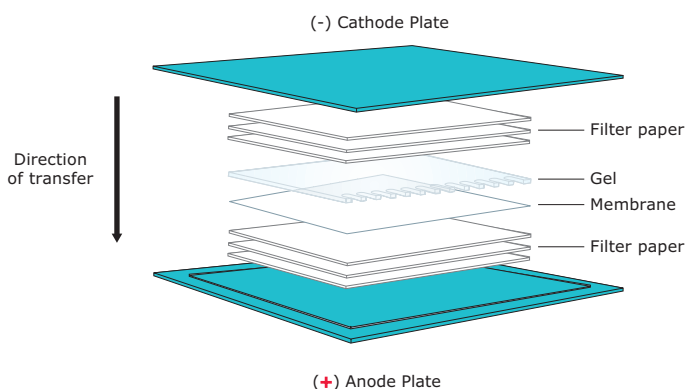
Transfer Stack Assembly

Refer to the manufacturer's specific operating instructions for the semi-dry transfer system being used.

Important: To ensure an even transfer, remove air bubbles between layers by carefully rolling a pipette or stirring rod over the surface of each layer in the stack. Do not apply excessive pressure to prevent damaging the membrane and gel.

For single transfers:

1. Place the anode electrode plate on a level bench top.
2. Place two pieces of filter paper soaked in anode buffer I in the center of the plate.
3. Place the filter paper soaked in anode buffer II on top of the first two sheets.
4. Place the membrane on top of the filter papers.
5. Place the gel on top of the membrane.
6. Place the three pieces of filter paper soaked in cathode buffer on top of the membrane.
7. Place the cathode electrode plate on top of the stack.



For multiple transfers:

1. Place the anode electrode plate on a level bench top.
2. Place two pieces of filter paper soaked in anode buffer I in the center of the plate.
3. Place the filter paper soaked in anode buffer II on top of the first two sheets.
4. Place the membrane on top of the filter papers.
5. Place the gel on top of the membrane. For the last gel, go to step 10.
6. Place a piece of filter paper soaked in cathode buffer on top of the gel.
7. Place a piece of dialysis membrane on top of the filter paper.
8. Place a piece of filter paper soaked in anode buffer II on top of the dialysis membrane.
9. Repeat steps 4 through 8 until all gels (up to the maximum for the unit) have been incorporated into the stack.
10. Place three pieces of filter paper soaked in cathode buffer on top of the last gel.
11. Place the cathode electrode plate on top of the stack.

Important: Do not bump the cathode plate cover during the run since it could disturb the alignment of the transfer stack and cause inaccurate results.

Protein Transfer

1. Insert the black cathode lead (-) into the cathode plate jack.
2. Insert the red anode lead (+) into the anode plate jack.
3. Connect the anode lead and cathode lead to their corresponding power supply inputs.
4. Turn on the power supply.
5. Set the current and let it run for the time indicated in the following chart:
6. Turn off the power supply when the transfer is complete.
7. Disconnect the system leads.
8. Remove the cover.
9. Remove and discard the filter papers.
Note: When using graphite plates, graphite particles from the anode electrode plate occasionally appear on the filter paper. These particles do not affect operation.
10. Remove the gel.
11. Remove the blotted membrane with a pair of forceps.

Current Density	Time Limit
0.8 mA/cm ² *	1–2 hours
1.2 mA/cm ²	1 hour
2.5 mA/cm ²	30–45 minutes
4.0 mA/cm ²	10–30 minutes

*The surface area (cm²) is calculated from the dimensions of the footprint of the stack on the anode plate. This value is independent of the number of gels in the stack.

Protocol 1.3 Dot Blotting/Slot Blotting: Vacuum Filtration Method

The following protocol describes a typical procedure for filtering proteins onto an Immobilon® PVDF transfer membrane. Please review the instructions supplied with your blotting unit for additional information.

- ▶ **Tip:** Detergents may inhibit the binding of proteins to the Immobilon® PVDF transfer membrane.
- ▶ **Tip:** Place a second sheet of membrane between the filter paper and primary membrane to ensure the membrane microporous structure is not compressed
- ▶ **Recommended** – Viscous samples must be diluted with buffer to reduce viscosity.

Required Equipment and Solutions

- Two sheets of Immobilon® PVDF transfer membrane, cut to size for blotting unit
- Two pieces of Whatman® 3MM filter paper or equivalent, cut to size for blotting unit
- Methanol, 100%
- Milli-Q® water
- Buffer, for sample loading and washes
- Blotting unit, dot blot or slot blot format

Setup

1. Prepare the membrane:
 - a. Wet the membrane in methanol for 15 seconds. The membrane should uniformly change from opaque to semi-transparent.
 - b. Carefully place the membrane in Milli-Q® water and soak for 2 minutes.
 - c. Carefully place the membrane in buffer and let equilibrate for at least 5 minutes.

2. Dissolve the sample in buffer. If the sample solution is cloudy, centrifuge to remove particles. If the sample is viscous, dilute with additional buffer.

Blotting Unit Assembly

See manufacturer's instructions for detailed assembly instructions.

1. Place one sheet of moistened filter paper on unit. Some units may require more than one sheet.
2. Place two sheets of the membrane on top of the filter paper.
3. Close the unit.
4. Connect to vacuum line.

Protein Transfer

1. Briefly apply the vacuum to remove excess buffer.
2. With the vacuum off, carefully pipette samples into the wells.
3. Apply vacuum to the blotting unit.
4. When all of the samples have filtered through the membrane, turn off the vacuum.
5. Add buffer to each well to wash down the sides. Apply the vacuum to filter through the wash buffer.
6. When all of the wash buffer has filtered through the membrane, turn off the vacuum.
7. To remove the blot, open the blotting unit.
8. Using forceps, carefully remove the membrane.

Protocol 1.4 Dot Blotting: Manual Spotting Method

Required Equipment and Solutions

- Immobilon® PVDF transfer membrane
- Two pieces of Whatman® 3MM filter paper or equivalent, cut to size of transfer membrane
- Paper towels
- Methanol, 100%
- Buffer

Setup

1. Prepare the membrane:
 - a. Wet the membrane in methanol for 15 seconds. The membrane should uniformly change from opaque to semi-transparent.
 - b. Carefully place the membrane in Milli-Q® water and soak for 2 minutes.
 - c. Carefully place the membrane in buffer and let equilibrate for at least 5 minutes.
2. If the sample solution is cloudy, centrifuge to remove particles. If the sample is viscous, dilute with additional buffer.
3. Wet one piece of filter paper with buffer

Transfer Stack Assembly

Assemble stack as follows (from the bottom up):

1. Place paper towels on work surface.
Note: Bottom towels should remain dry throughout blotting procedure.
2. Place dry filter paper on paper towels.
3. Place pre-wet filter paper on the dry filter paper.
4. Place the pre-wet membrane on wet filter paper.

Protein Transfer

1. Spot 1 – 5 μL of sample onto the membrane. Sample should wick into membrane.
Note: Membrane should be wet enough to absorb sample, but not so wet that sample spreads across membrane.
2. After sample is absorbed, place membrane on clean filter paper to dry.

Protocol 1.5 Optimization of Blocking Reagents

Required Equipment and Solutions

- Small piece of Immobilon® PVDF transfer membrane
- One piece of Whatman® 3MM filter paper or equivalent, cut to the membrane dimensions
- Methanol, 100%
- Wash buffer: Phosphate-buffered saline (PBS) or Tris-buffered saline (TBS) containing 0.05 – 0.1% Tween®-20 surfactant
- Blocking solution(s) to be tested
- Detection reagents to be tested
- X-ray film. CCD based digital imaging system cooled.

Setup

1. Prepare the membrane:
 - a. Wet PVDF membrane in methanol for 15 seconds. The membrane should uniformly change from opaque to semi-transparent.
 - b. Carefully place the membrane in Milli-Q® water and soak for 2 minutes.
 - c. Carefully place the membrane in buffer and let equilibrate 5 minutes with gentle shaking.
2. Wet filter paper in wash buffer.

Procedure

1. Place the membrane on wet filter paper.
2. Spot 5 μL of blocking solution onto the membrane and allow it to absorb completely.
3. Add detection reagents to the blot and incubate for 5 minutes.
4. Drain the excess reagents.
5. Cover the blot with a clean plastic wrap or sheet protector and remove any air bubbles.
6. Expose the blot to a suitable X-ray film for a few minutes. Appearance of a black spot on the film indicates cross-reactivity between the blocking reagent and the detection substrate.

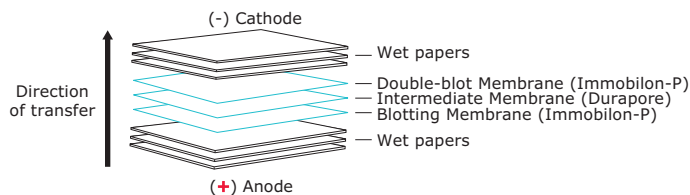
Protocol 1.6 Double Blotting: An Innovative Procedure for Eliminating the Nonspecific Binding of Secondary Antibodies

Required Equipment and Solutions

- Semi-dry transfer apparatus
- Two sheets Immobilon®-P membrane (blot and double-blot membranes)
- Durapore® hydrophilic PVDF membrane, 0.65 µm pore size (Cat. No. DVPP 000 10)
- Six pieces of Whatman® 3MM filter paper or equivalent, cut to the membrane dimensions
- 0.7% (v/v) acetic acid solution
- Phosphate-buffered saline (PBS), pH 7.4

Procedure

1. Blot resolved proteins onto Immobilon®-P membrane; block; apply primary antibody; and wash as described in Protocol 3.1.
2. Cut a sheet of Durapore® membrane to the membrane dimensions.
3. Immerse the Durapore® membrane in the acid solution for at least 10 minutes. Pre-wet the unblotted Immobilon®-P membrane in methanol for 3 seconds; rinse in water for 2 minutes; and equilibrate in acid solution for 10 minutes. Wet the stacks of filter paper in the acetic acid solution by capillary wicking.
4. Place the blotted membrane onto a stack of filter paper and cover it with the Durapore® membrane and the double-blot membrane. Quickly put a second stack of filter paper onto the double-blot membrane to prevent drying. Position this sandwich in the blotting instrument with the blotted membrane facing the anode and the double-blot membrane facing the cathode.



5. Apply a constant intensity of 0.8 mA/cm² for 10 minutes.
6. Disassemble the sandwich.
7. Rinse the double-blot membrane twice in PBS using fresh buffer each time.
8. Block the double-blot membrane.
9. If the original blotting membrane (on which the blotted proteins are retained) will be reprobbed with the same or another primary antibody, it can be kept in PBS after the acidic transfer.
10. Proceed with probing of the double-blot membrane with the appropriate secondary antibody and protein detection as described in Protocol 3.1, Page 37. See Pages 22-23 for an explanation of double blotting.

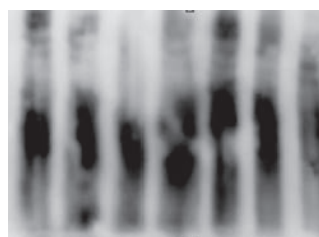
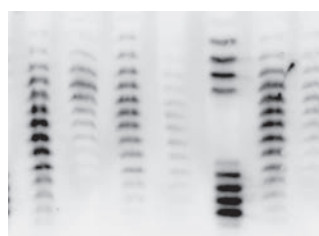


Figure 20.

Isoelectric patterns of urinary endogenous erythropoietin (EPO) without (top) and with (bottom) double blotting.
Data courtesy of Dr. Françoise Lasne, Laboratoire National de Dépistage du Dopage.



Protocol 1.7 Membrane Drying Methods

Membranes must be dried before continuing on to transillumination or rapid immunodetection procedures. After proteins have been blotted onto the membrane, rinse the membrane with Milli-Q® water. As the blotted membrane dries, it becomes opaque. The table below details four drying methods and required times. Always wait the full length of drying time to ensure that all liquid has evaporated from within the membrane's pore structure.

Drying Method	Required Time*
Soak in 100% methanol for 10 seconds. Remove from methanol and place on piece of filter paper until dry.	15 minutes
Secure between two sheets of filter paper and place in a vacuum chamber.	30 minutes
Incubate at 37°C for 1 hour	1 hour
Place on lab bench and let dry at room temperature.	2 hours

*Longer times required in higher humidity environments.

2. Protein Visualization

The following protocols describe typical procedures for staining proteins immobilized on an Immobilon® PVDF transfer membrane.

- ▶ **Tip:** Only reversible stains or transillumination should be used prior to immunodetection.
- ▶ **Tip:** To visualize fainter bands, increase the methanol concentration to 40%.
- ▶ **Tip:** When examining a stained blot, the degree of contrast is best while the membrane is still wet. For photographic purposes, use a wet blot and light transmitted through the membrane.

Protocol 2.1 Visualization by Transillumination

Transillumination is a nondestructive, reversible technique²⁷ based on the change of PVDF from opaque to semi-transparent when wet in methanol. 20% methanol will wet out regions of the membrane coated with protein, but not areas without protein.

- ▶ **Tip:** When developing a new transfer protocol or working with a new sample type, stain the gel to verify that all of the proteins have eluted from the gel.

The protein bands appear as clear areas when placed on a light box. Detection sensitivity is comparable to Coomassie® Brilliant Blue R stain when used in conjunction with photography.

Required Equipment and Solutions

- 20% (v/v) methanol
- Shallow tray, large enough to hold membrane
- White light box
- Black paper

Procedure

1. Dry the blot completely using one of the drying methods in Protocol 1.7, Page 33.
2. Fill the tray with enough 20% methanol to cover the blot.
3. Place the blot in 20% methanol for 5 minutes.
4. Place the blot on a light box and mask the areas around the blot with a sheet of black paper.
5. The bands appear as clear areas against an opaque background (Figure 10, blot A, Page 17).
6. If a permanent record is desired, photograph the wet blot.
7. When all the methanol has evaporated, the blot will return to its original appearance.

Protocol 2.2 Visualization by Reversible Staining

Ponceau-S Red

This stain produces pink bands on a light background.

Required Equipment and Solutions

- Stain: 0.2% Ponceau-S red, 1% acetic acid. Prepare by diluting stock solution (2% dye in 30% (w/v) trichloroacetic acid and 30% (w/v) sulfosalicylic acid) 1:10 in 1% (v/v) acetic acid
- Methanol, 100%
- Destain solution: 0.1 N NaOH
- Milli-Q® water
- Shallow tray, large enough to hold membrane

Procedure

1. If the blot is dry, wet in 100% methanol.
2. Fill the tray with enough stain to cover the blot.
3. Place the blot in stain and agitate for 1 minute.
4. Remove the blot and rinse thoroughly with Milli-Q® water until the desired contrast has been achieved.
5. To remove the stain completely, wash the blot with 0.1 N NaOH.

CPTS (Copper Phthalocyanine Tetrasulfonic Acid)

This stain produces turquoise blue bands on a light background.

Required Equipment and Solutions

- Stain: 0.05% (w/v) CPTS in 12 mM HCl
- Wash solution I: 12 mM HCl
- Wash solution II: 20% (v/v) methanol
- Methanol, 100%
- Milli-Q® water
- Destain solution: 0.5 M NaHCO₃
- Shallow tray, large enough to hold membrane

Procedure

1. If the blot is dry, wet in 100% methanol.
2. Fill the tray with enough stain to cover the blot.
3. Place the blot in stain and agitate for 1 minute.
4. Place the blot in wash solution I to remove excess stain and achieve the desired contrast.
5. To remove stain from the background completely, wash the blot with destain wash solution II.
6. To completely destain the proteins, agitate the blot in the destain solution until the stain has been removed.

Protocol 2.3 Visualization by Irreversible Staining

Coomassie® Brilliant Blue R Dye

This stain produces dark bands on a light background.

Important: This stain will interfere with immunodetection.

Required Equipment and Solutions

- Stain: 0.1% (w/v) Coomassie® Brilliant Blue R in 50% (v/v) methanol, 7% (v/v) acetic acid
- Wash solution I: 50% (v/v) methanol, 7% (v/v) acetic acid
- Wash solution II: 90% (v/v) methanol, 10% (v/v) acetic acid
- Methanol, 100%
- Milli-Q® water
- Shallow tray, large enough to hold membrane

Procedure

1. If the blot is dry, wet in 100% methanol.
2. Fill the tray with enough stain to cover the blot.
3. Place the blot in stain and agitate for 2 minutes.
4. Remove the blot and rinse briefly with Milli-Q® water.
5. Place the blot in wash solution I and agitate for 10 minutes to remove excess stain.
6. Place the blot in wash solution II and agitate for 10 minutes to completely destain the background.

Amido Black

This stain produces dark bands on a light background.

Important: This stain will interfere with immunodetection.

Required Equipment and Solutions

- Stain: 0.1% (w/v) amido black in 25% (v/v) isopropanol, 10% (v/v) acetic acid
- Wash solution: 25% (v/v) isopropanol, 10% (v/v) acetic acid
- Methanol, 100%
- Milli-Q® water
- Shallow tray, large enough to hold membrane

Procedure

1. If the blot is dry, wet in 100% methanol.
2. Fill the tray with enough stain to cover the blot.
3. Place the blot in stain and agitate for 2 minutes.
4. Remove the blot and rinse briefly with Milli-Q® water.
5. Place the blot in wash solution and agitate for 5 to 10 minutes to remove excess stain.

3. Immunodetection

Protocol 3.1 Standard Immunodetection Method

Standard immunodetection is performed on blotted proteins directly after electrotransfer. (If the membrane was dried after transfer, thoroughly wet the blot in methanol for 5 minutes before proceeding to immunodetection.) The unoccupied membrane binding sites on the wet blot are blocked with optimized reagents (see Protocol 1.5). While standard immunodetection requires blocking and a lengthy procedure (4 hours), this method typically requires less optimization for new sample types.

- ▶ **Tip:** Immobilon®-PSQ transfer membrane has a smaller pore size (0.2 µm) and higher surface area than Immobilon®-P transfer membrane (0.45 µm). Increased background can be expected on Immobilon®-PSQ and the blocking and wash steps will need to be adjusted accordingly.
- ▶ **Tip:** Phosphatases in the blocking solution may dephosphorylate blotted proteins.
- ▶ **Tip:** Do not use sodium azide in buffers, as it inhibits HRP activity.
- ▶ **Tip:** Do not let the blot dry out at any time during and after blocking.
- ▶ **Tip:** If more than one blot is placed in a container, insufficient buffer volume will cause the blots to stick together.
- ▶ **Tip:** Non-fat dry milk cannot be used as a blocking reagent with biotin-avidin systems.
- ▶ **Tip:** Sensitivity of chromogenic detection is typically at least an order of magnitude lower than of chemiluminescent detection.
- ▶ **Tip:** High nonspecific signal can be alleviated by higher dilution of the primary antibody or reduced protein load on the gel.
- ▶ **Tip:** Immobilon®-FL membrane can be scanned dry or wet.
- ▶ **Tip:** High overall background can be minimized by higher dilution of the enzyme-conjugated secondary antibody.

Required Solutions

- Primary antibody (specific for protein of interest)
- Secondary antibody (specific for primary antibody), labeled with alkaline phosphatase or horseradish peroxidase
- Detection reagent appropriate to the enzyme conjugate
- Phosphate-buffered saline (PBST): 10 mM sodium phosphate, pH 7.2, 0.9% (w/v) NaCl, up to 0.1% Tween®-20 detergent
- TBST: 10 mM Tris, pH 7.4, 0.9% (w/v) NaCl, up to 0.1% Tween®-20 detergent
- Blocking solution: 1% (w/v) BSA (bovine serum albumin), 0.05% Tween®-20 detergent
- Milli-Q® water

Required Equipment

- Shallow trays, large enough to hold blot
- Glass plates
- Plastic wrap (e.g., Saran™ film), freezer bag, or sheet protector

Setup

1. Dilute the primary antibody in the blocking solution to the desired working concentration.
2. Dilute the secondary antibody in the blocking solution to the desired working concentration.

Note: Enough solution should be prepared to allow for 0.1 mL of antibody solution (primary and secondary) per cm² of membrane.

Antibody Incubations

1. Place the blot in the blocking solution and incubate with agitation for 1 hour.
2. Place the blot in the primary antibody solution and incubate with agitation for 1 hour. The solution should move freely across the surface of the membrane.
3. Place the blot in PBST and wash for 10 minutes. Repeat twice with fresh buffer.
4. Place the blot in the secondary antibody solution and incubate with agitation for 1 hour at RT or 37 °C.
5. Place the blot in PBST and wash for 10 minutes. Repeat twice with fresh buffer.
6. Proceed with either chromogenic, chemiluminescent or fluorescent detection.

Chromogenic Detection

1. Prepare the substrate according to manufacturer's instructions.
2. Place the blot in a clean container and add substrate to completely cover the surface of the membrane. Incubate for 10 minutes with mild agitation or until signal reaches desired contrast.
3. Rinse the blot with Milli-Q® water to stop the reaction.
4. Store the blot out of direct light to minimize fading. Blot may be stored dry.

Chemiluminescent Detection

Follow manufacturer's instructions.

1. Prepare the substrate according to manufacturer's instructions.
2. Place the blot in a container and add substrate to completely cover the membrane. Incubate for 1 minute.
3. Drain excess substrate.
4. Place the blot on a piece of clean glass and wrap in plastic wrap or place directly into CCD based digital imaging system.
Note: A cut-to-size sheet protector or a freezer bag can also be used.
5. Gently smooth out any air bubbles.
6. In a dark room, place the wrapped membrane in a film cassette.
7. In the dark room, place a sheet of autoradiography film on top and close the cassette.
8. Expose film. Multiple exposures of 15 seconds to 30 minutes should be done to determine the optimum exposure time; 1 to 5 minutes is common. Multiple images should also be acquired. If using CCD based imaging system.

Fluorescent Detection

Required Equipment

- Autoradiography film and cassette
- Dark room
- Autoradiography film processing equipment
- Proteins blotted onto Immobilon®-FL transfer membrane and probed with antibodies
- Mylar® wrap
- Fluorescent imaging equipment

The following is a general protocol for fluorescent immunodetection. For optimal results, refer to manufacturer's protocol provided with the reagents.

Note: If using chemifluorescent reagents, follow reagent manufacturer's directions.

1. Place the blot in diluted fluorescent dye-labeled secondary antibody solution and incubate for 1 hour with gentle agitation.
2. Wash the blot with wash buffer 3–5 times for 5 minutes each.
3. Place the blot onto a piece of clean filter paper to dry.
4. If using a wrap, use Mylar®. Do not use Saran™ wrap because it permits light to shine through and quench fluorescence.
5. Image the blot using an appropriate fluorescence scanner.

Protocol 3.2 Rapid Immunodetection Method

Rapid immunodetection takes advantage of the fact that antibodies cannot bind to the hydrophobic (non-wetted) surface of the Immobilon®-P transfer membrane, but will bind to a protein immobilized on the membrane. Rapid immunodetection is compatible with both chromogenic and chemiluminescent substrates.

► **Tip:** Use the Rapid Immunodetection Method to quickly visualize higher-abundance proteins. For higher sensitivity, use Protocol 3.1.

The major advantage of rapid immunodetection is that blocking is not required, saving time and eliminating the risks involved.⁴⁷ Also, because excess antibody won't bind to a dry membrane, the amount of washing required is reduced. As a result, the total time for analysis is under two hours, as opposed to over four hours for the standard method.

Important: The blot must be thoroughly dry before beginning rapid immunodetection. Refer to Protocol 1.7.

Required Solutions

- Primary antibody (specific for protein of interest)
- Secondary antibody (specific for primary antibody), labeled with alkaline phosphatase or horseradish peroxidase
- Substrate appropriate to the enzyme conjugate.
- Phosphate buffered saline (PBS): 10 mM sodium phosphate, pH 7.2, 0.9% (w/v) NaCl
- Antibody dilution buffer: 1% (w/v) BSA (bovine serum albumin), 0.05% Tween®-20 detergent
- Methanol, 100%
- Milli-Q® water

Required Equipment

- Shallow trays, large enough to hold blot
- Plastic wrap (e.g., Saran™), freezer bag, or sheet protector
- Autoradiography film and cassette
- Dark room and autoradiography film processing equipment

Setup

1. Dry the blot completely using one of the drying methods in Protocol 1.7. Do not re-wet blot in methanol.
2. Dilute the primary antibody in dilution buffer to the desired working concentration.
3. Dilute the secondary antibody in dilution buffer to the desired working concentration.

Note: Enough solution should be prepared to allow for 0.1 mL of antibody solution (primary and secondary) per cm² of membrane.

Antibody Incubations

1. Place the blot in the primary antibody solution and incubate with agitation for 1 hour. The solution should move freely across the surface of the membrane.
2. Place the blot in PBS and wash for 5 minutes. Repeat twice with fresh buffer.
3. Place the blot in the secondary antibody solution and incubate with agitation for 30 minutes.
4. Place the blot in PBS and wash for 5 minutes. Repeat twice with fresh buffer.
5. Proceed with chromogenic, chemiluminescent or fluorescent protein detection as described in Protocol 3.1.

Protocol 3.3 Simple Immunodetection Using the Immobilon® GO Device

Immobilon® GO is a hands-free, immunodetection device designed for mini Western blot processing. The unique design uses continuous, sequential fluid convection across a flow matrix to enable autonomous antibody binding and membrane wash steps during immunodetection. Immobilon® GO is intended for one-time use and does not require an external power or vacuum source for operation. Standard non-fat dry milk TBST buffer is used to allow for easy transition from traditional Western blot protocols. Western blots can be processed in as little as 3 hours or may remain in the device overnight for processing.

General guidelines for Immobilon® GO immunodetection device

- Membrane should not be larger than 8.5 cm x 8 cm
- Immobilon® GO device requires a flat level surface for proper operation
- Avoid indenting membrane when labeling
- Once reagents have been loaded, do not move or open Immobilon® GO device until immunodetection is completed (3 hours)

Materials required for immunodetection, but not supplied:

- Non-fat dry milk (NFDM)
- Tris-buffered saline with Tween® 20 (TBS-T) (20 mM Tris-HCl, pH 7.4, 150 mM NaCl, 0.1% Tween® 20)
- Blot Roller
- Forceps
- Incubation Tray
- Primary and secondary antibodies

Setup

1. Prepare 25 mL of 2% non-fat dry milk (NFDM) blot buffer by completely dissolving 0.5 g of NFDM in 25 mL TBS-T.
2. If blotting membrane was previously dried, wet in deionized water when using nitrocellulose or with methanol when using PVDF, washing thoroughly with deionized water to remove residual alcohol.
3. Place wet membrane (protein side facing up) into a dish. Add 10 mL of 2% NFDM blot buffer on top of the membrane. Incubate for 5-10 minutes.
4. Prepare primary and secondary antibody dilution in 2% NFDM blot buffer. A total volume of 1 mL is required for each antibody.

Blot Assembly & General Protocol

1. Open Immobilon® GO device and fold back both parts of the flow matrix.
2. Position blotting membrane (protein side up), 1 cm from stage edge. Low molecular weight proteins should be facing the notched flow matrix.
3. Fold flow matrix over the membrane and add 4 mL of 2% NFDM blot buffer evenly across the flow matrix.
4. To assure even contact between the membrane and flow matrix, pull the blot roller from the loading port end toward the flow matrix tab.
5. Using the blot roller fold over the flow matrix tab. Assure all air bubbles have been removed and close contact between the flow matrix and membrane has been established.
6. Replace the lid.
7. Seal lid by gently push down in area specified by loading ports; a clicking sound indicates that lid is closed.
8. To activate Immobilon® GO device, add 2% NFDM blot buffer to the reagent loading ports in the order and volumes shown below (i.e., fill port 1A, followed by 1B).

Loading Port	Volume (mL)	
	Side A	Side B
1	0.5	0.5
2	0.5	0.5
3	0.5	0.5
4	3.0	3.0

9. Add the primary and secondary antibodies in the order and volumes as outlined in the table below.

Reagent	Loading Port	Volume (mL)	
		Side A	Side B
Primary Antibody	1	0.5	0.5
Secondary Antibody	2	0.5	0.5

10. Incubate Immobilon® GO device for a minimum of 3 hours at room temperature.
11. Remove blot from Immobilon® GO device and wash twice for 5 minutes with 50 mL TBS-T.
12. Continue with desired detection method, as described in traditional Western blotting protocols.

Protocol 3.4 Rapid Immunodetection Using the SNAP i.d.[®] 2.0 System

The SNAP i.d.[®] 2.0 Protein Detection System is a unique, vacuum-driven system, that pulls reagents through the membrane, increasing membrane contact while eliminating dependence on diffusion or agitation. This improvement in three-dimensional reagent distribution decreases the length of time required for immunodetection. While traditional Western blotting requires 4 to 24 hours for the blocking, antibody incubation and washing steps, the SNAP i.d.[®] 2.0 protocol takes only 30 minutes with no loss of signal intensity or reduction in blot quality. All immunodetection steps after protein transfer to a membrane (i.e., blocking, washing, and primary and secondary antibody incubations) can be performed with the SNAP i.d.[®] 2.0 System.

Required Solutions

- Blocking reagent such as non-fat/low fat dry milk (0.5% or less), casein, bovine serum albumin (BSA) or other commercially available blocking agents such as Block-CH reagent
- Antibodies (monoclonal and/or polyclonal)
- Detection reagents
- Wash buffer: Tris- or phosphate-buffered saline solution, pH 7.4, supplemented with Tween[®]-20 surfactant (TBST or PBST)

Required Equipment

- Vacuum source: Pump or other uniform vacuum source that can deliver a sustained minimum pressure of 410 millibar (12 in. Hg) and 34 L/min
- **Note:** Millipore[®] WP61 series chemical duty pumps can be used, but may require longer processing times
- One-liter or larger vacuum flask with stopper (for waste collection). A Millex[®]-FA50 filter (or equivalent) is recommended between the vacuum flask and the vacuum source to protect the vacuum source from contamination
- Vacuum tubing to connect vacuum flask to vacuum source
- Forceps
- Blot with transferred proteins

Setup

1. Place the SNAP i.d.[®] 2.0 base on a level bench top.
2. Attach the vacuum tubing to the back of the system by pushing the coupling insert on the end of the tubing into the quick disconnect fitting at the back of the system base until it clicks.
3. Connect the other end of the tubing to a vacuum source. Use a one-liter vacuum flask as a trap and a Millex[®]-FA50 filter (Cat. No. SLFA05010) to protect the vacuum source from contamination.

Note: Any vacuum source that can deliver 410 millibar (12 in. Hg) and 34 L/min is sufficient. If the vacuum source operates at higher than 410 millibar, the SNAP i.d.[®] 2.0 system will automatically regulate the vacuum pressure.

Blot Assembly & General Protocol

1. Hold the blot holder by the support layer (blue edges) and wet the membrane layer (white) with distilled water in the wetting tray provided. Place the wetted blot holder on the rolling pad.
2. If required, pre-wet the blot in methanol and water, then place it in the center of the blot holder with the protein side down.
3. Roll the blot gently to remove air bubbles, then close the blot holder and roll one more time.
4. Open the blot holding frame, flip the blot holder so that it is protein side up, then place it inside the frame. A notch in the blot holder ensures correct placement in the frame.
5. Close and lock the frame. Add 30 mL of blocking solution. Press the frame down and turn the system knob to apply vacuum. When frame is completely empty, TURN VACUUM OFF.
6. Apply 5 mL (Mini blot) or 10 mL (Midi blot) of primary antibody across the surface of the blot holder.
7. Incubate for 10 minutes at room temperature. Solution will be absorbed into the blot holder and surface may appear dry.
8. Press the frame down and apply vacuum. Wait 5–8 seconds until the frame is completely empty.
9. With vacuum running continuously, add 30 mL of wash buffer. Repeat the washing step 3 more times (total of 4 washes). When frame is completely empty, TURN VACUUM OFF.
10. Apply 5 mL (Mini blot) or 10 mL (Midi blot) of secondary antibody evenly across the blot holder surface. Incubate for 10 minutes at room temperature with vacuum off. Again, solution will be absorbed into the blot holder and surface may appear dry.
11. Press frame down and apply vacuum. Wait 5–8 seconds until frame is completely empty. With vacuum running continuously, add 30 mL of wash buffer. Repeat the washing step 3 more times (total of 4 washes).
12. Turn vacuum off and remove blot holder from frame. Remove blot from blot holder and incubate with the appropriate detection reagent.

Extended Primary Antibody

Incubation: One Hour to Overnight

1. Perform steps 1 through 5 of the general protocol.
2. Add 5 mL (Mini blot) or 10 mL (Midi blot) of 1X primary antibody (concentration normally used during standard immunodetection).
3. Cover the blot holding frame with the lid. If desired, remove it from the base and incubate with constant shaking. If overnight incubation is required, refrigeration is recommended. Although the surface of the blot holder may look partially dry, the blot will not dry out, since the solution is contained in the frame.

Note: If processing more than one blot for an extended period of time, the blot holding frames can be stacked to reduce the working space.

Optional: If recovering primary antibody, place antibody recovery trays into base before proceeding to step 4.

4. After the extended incubation period, place the frame back on the base, remove the lid, and follow steps 8 through 12 of the general protocol.

Note: Extended incubation periods for the secondary antibody are not required for the SNAP i.d.[®] 2.0 System. A 5X concentration of secondary antibody for 10 minutes is recommended.

Antibody Recovery

1. Perform steps 1 through 5 of the general protocol, but increase the vacuum time to 2 minutes to ensure that all the blocking solution has been removed from the grooves and channels of the frame.
 2. Remove the blot holding frame from the base and wipe any residual liquid from the bottom of the frame with a paper towel.
 3. Place the antibody collection tray into the base, making sure that the positioning holes in the antibody collection tray line up with the positioning pins in the base.
 4. Place the blot holding frame back into position on the base.
 5. Apply primary antibody and incubate as indicated in steps 6 and 7 of the general protocol.
 6. After the 10-minute incubation, turn the vacuum on and wait one minute to ensure that all of the antibody has been collected.
- Note:** When processing two frames at the same time, apply vacuum first to one side, then to the other. This ensures full vacuum force on each frame and improves volume recovery.
7. Turn the vacuum off and remove the frame. Remove the antibody collection tray.
 8. Transfer the antibody to a suitable container for storage or analysis.
 9. Place the blot holding frame back into position and continue with step 9 of the General Protocol.

Protocol 3.5 High Salt Wash to Remove Persistent Background

Required Solutions

- High salt buffer: PBS or TBS buffer supplemented with 0.5M NaCl and 0.2% SDS.

Procedure

1. Follow Protocol 3.1, through primary and secondary antibody incubations.
2. After the secondary antibody incubation, place the blot in the high salt buffer and incubate for 30 minutes with gentle shaking.
3. Rinse blot with Milli-Q[®] water and proceed with chromogenic, chemiluminescent or fluorescent protein detection as described in the Protocol 3.1.

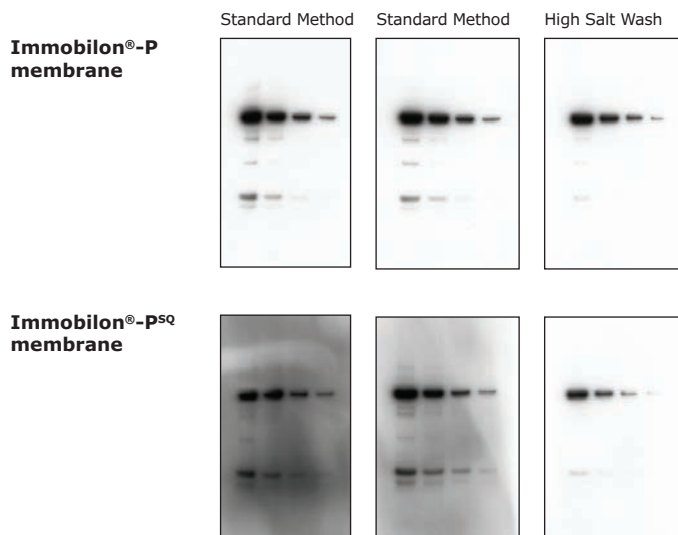


Figure 21.

Adding 0.5 M NaCl and 0.2% SDS to the wash buffer can reduce background noise on membranes with high protein binding capacity, such as Immobilon[®]-P^{SQ} membrane.

4. Membrane Stripping

Two protocols for removing antibodies from blots are presented below. The first is applicable to any chemiluminescent substrate system and uses a combination of detergent and heat to release the antibodies. The second is commonly used for applications where antibodies have to be separated from an antigen and employs low pH to deactivate the antibody binding site.

Neither method will remove the colored precipitates generated from chromogenic detection systems (e.g., BCIP, 4CN, DAB and TMB). However, it is still possible to analyze the blot with another antibody specific to a different target protein.

Important: The blot should not be allowed to dry between rounds of immunodetection. Any residual antibody molecules will bind permanently to the membrane if it is allowed to dry.

Protocol 4.1 Stripping by Heat and Detergent

Required Equipment and Solutions

- Stripping solution: 100 mM 2-mercaptoethanol, 2% (w/v) SDS, 62.5 mM Tris-HCl, pH 6.7
- Phosphate buffered saline (PBS): 10 mM sodium phosphate, pH 7.2, 0.9% (w/v) NaCl
- Shallow tray, large enough to hold the membrane

Procedure

1. In a fume hood, place the blot in stripping solution and incubate with agitation for 30 minutes at 50 °C.
2. Place the blot in buffer and agitate for 10 minutes. Repeat with fresh buffer.
3. **Optional:** Repeat the initial detection protocol (omitting the primary antibody step) to make sure that the antibodies have been inactivated or stripped from the membrane.
4. Place the blot in buffer and agitate for 10 minutes.
5. Proceed to the blocking step for the next round of immunodetection.

Protocol 4.2 Stripping by Low pH

Required Equipment and Solutions

- Stripping solution: 25 mM glycine-HCl, pH 2, 1% (w/v) SDS.
- Phosphate buffered saline (PBS): 10 mM sodium phosphate, pH 7.2, 0.9% (w/v) NaCl.
- Shallow tray, large enough to hold the membrane.

Procedure

1. Place the blot in stripping solution and agitate for 30 minutes.
2. Place the blot in buffer and agitate for 10 minutes. Repeat with fresh buffer.
3. Proceed to the blocking step for the next round of immunodetection.

Protocol 4.3 Stripping with the ReBlot™ Plus Western Blot Recycling Kit

Required Equipment and Solutions

- Standard blot or blot strips, on nitrocellulose or PVDF/nylon membrane
- Blocking solutions
- Plastic wrap for storage of blots that will not be reprobbed immediately
- Distilled water, for reagent dilution
- Plastic trays for incubation of blots or blot strips in stripping, washing and blocking solutions
- Positive and negative stripping controls

Procedure

Note: The blots or individual strips that are to be reused should be prepared for stripping immediately after their first usage. If stripping cannot be performed right away, membranes can be wrapped in plastic wrap and stored moist in PBS at 4°C. DO NOT STORE BLOTS IN DRY FORM.

1. Fill a plastic tray with appropriate amount of 1X Antibody Stripping Solution (supplied in kit).
2. Using tweezers or forceps, submerge blot or blot strips in stripping solution. Incubate with gentle mixing for 15 minutes at room temperature.
3. Fill a clean plastic tray with an equal amount of blocking buffer. Conventional blocking buffers such as 20 mM Tris HCl, pH 8.0; 150 mM NaCl; 0.1% Tween®-20 detergent; 5% non-fat dry milk or similar are appropriate.
4. Wash blots two times 5 minutes each with blocking buffer.
5. The blot is now ready for reprobing with antibodies.

5. Protein Digestion

Protocol 5.1 On-Membrane Protein Digestion for Mass Spectrometry

This method describes the preparation of proteins immobilized on Immobilon® PVDF transfer membrane for analysis by mass spectrometry.

► **Tip:** Mass spectrometry-compatible membrane stains include Coomassie® Brilliant Blue dye, amido black, and Sypro® blot stains.

Required Equipment and Solutions

- 50% methanol in Milli-Q® water
- 30% acetonitrile, 50 mM ammonium bicarbonate
- Trypsin, dissolved in Milli-Q® water at 0.1 mg/mL
- 80% acetonitrile in Milli-Q® water
- Vacuum centrifuge
- Microcentrifuge tubes
- MALDI-TOF matrix

Procedure

1. Transfer protein from the gel to the membrane (Protein Transfer Protocols).
2. Stain the blotted membrane with Coomassie® Brilliant Blue dye or amido black (Protocol 2.3).
3. Wash the stained membrane with water and dry.
4. Excise pieces of membrane containing proteins of interest with a clean scalpel and place in separate microcentrifuge tubes.
5. Destain membrane pieces with 500 µL of 50% methanol for 2 hours at room temperature.
6. Remove supernatant and dry the membrane.
7. Add 10 µL of 30% acetonitrile, 50 mM ammonium bicarbonate, and 4 µL of 0.1 mg/mL trypsin.
8. Incubate overnight at room temperature.
9. Transfer supernatants to clean microcentrifuge tubes.
10. Extract peptides with 20 µL of 80% acetonitrile for 15 minutes with sonication.
11. Pool the extracts with the previous supernatants.
12. Dry the digests in the vacuum centrifuge.
Note: Alternatively, ZipTip® SCX pipette tips can be used to purify and concentrate the peptide digest instead of drying. To do this, acidify the digest with 0.5% TFA (make sure the pH is < 3) and follow the instructions in the ZipTip® SCX Pipette Tip User Guide (Lit. No. P36444).
13. Resuspend each peptide extract in a small volume of 30% acetonitrile, 0.1% TFA.
14. Load 1 – 2 µL of the digest onto a MALDI plate, and mix with 1 µL of matrix.

6. Blot Storage

Protocol 6.1 Preparation of Protein Blots for Long-Term Storage

PVDF is a chemically resistant polymer with excellent long-term stability. For blots that need to be stored for use at a later date, storage conditions are determined by the stability of the proteins bound to the membrane. We recommend storage at low temperature, but room temperature may be adequate for some proteins.

Required Materials

- Dry blotted Immobilon® PVDF transfer membrane
- Two sheets of Immobilon® blotting filter paper or equivalent
- Two sheets of card stock or thin cardboard
- Paper clips
- Plastic bag

Note: Blots stored in a freezer should not be subjected to mechanical shock, which can cause breakage of the membrane. The blot should be allowed to come to room temperature before removal from the plastic bag. Blots may also be stored wet at 4°C in a plastic bag, but a bactericide such as sodium azide should be added to prevent bacterial growth. The azide must be thoroughly washed out of the blot prior to use as it inhibits HRP activity.

Procedure

1. Place the dry blot between two clean sheets of Whatman® 3MM paper.
2. Place the blot-filter paper sandwich between two sheets of card stock.
3. Clip the stack together along the edges. The clips should not overlap the blot.
4. Place the stack into a sealable plastic bag.
5. Close or seal the bag.
6. Store the blot at the desired temperature:
 - 4°C For up to 2 weeks
 - 20°C For up to 2 months
 - 70°C For longer-term storage

6. Appendices

Timeless Recipes and Solutions for Blotting Success

Standard recipes for Western blotting

TBS 10x (concentrated Tris-buffered saline)

[Tris]: 200 mM
[NaCl]: 1500 mM

For 1 L:

24 g Tris base (Product 93362 – FW: 121.1 g/mol)
88 g NaCl (Product 71376)
Dissolve in 900 mL Milli-Q® water. Adjust pH to 7.6 with 12 N HCl (Product H1758). Add Milli-Q® water to a final volume of 1 L.

1X solution: Mix 1 part of the 10X solution with 9 parts Milli-Q® water and adjust pH to 7.6.

TBST (Tris-buffered saline, 0.1% Tween® 20)

[Tris]: 20 mM
[NaCl]: 150 mM
[Tween® 20]: 0.1% (w/v)

For 1 L:

100 mL TBS 10X
900 mL Milli-Q® water
1 mL Tween®-20 detergent (Product P9416)

4X SDS-PAGE sample loading buffer

[Tris-HCl]: 0.2 M
[DTT]: 0.4 M
[SDS]: 277 mM, 8.0% (w/v)
[Bromophenol blue]: 6 mM
[Glycerol]: 4.3 M

For 7.5 mL:

1.5 mL 1 M Tris-HCl pH 6.5 (Product 20-160)
3 mL 1 M dithiothreitol (DTT, Product 43816)
0.6 g sodium dodecyl sulfate (SDS, Product L3771)
30 mg bromophenol blue (Product B0126)
2.4 mL glycerol (Product G5516)

Bring final volume to 7.5 mL with Milli-Q® water
If solution is orange/yellow in color, add 1 drop of 5 M NaOH (Product S8263) to adjust pH
Make 500 µL aliquots and store at -20°C

SDS-PAGE separating gel making buffer

[Tris-HCl]: 1.5 M

For 500 mL:

118.2 g Tris-HCl (Product 93363 - FW:157.60 g/mol)
Add 450 mL Milli-Q® water and adjust pH to 8.8
Add Milli-Q® water to final volume of 500 mL
Filter and degas

SDS-PAGE stacking gel making buffer

[Tris-HCl]: 1 M

For 500 mL:

78.8 g Tris-HCl (Product 93363 - FW:157.60 g/mol)
Add 450 mL Milli-Q® water and adjust pH to 6.8
Add Milli-Q® water to final volume 500 mL
Filter and degas

SDS-PAGE 10X gel running buffer

[Tris]: 248 mM
[Glycine]: 1.92 M
[SDS]: 35 mM, 1% (w/v)

For 2 L:

60 g Tris base (Product 93362 – FW: 121.1 g/mol)
288 g glycine (Product 50046)
20 g sodium dodecyl sulfate (SDS, Product L3771)
Add Milli-Q® water to yield solution with final volume of 2 L
No need to pH, filter, or degas
Dilute to 1X for running SDS-PAGE gels

10X PBS

[NaCl]: 1.37 M
[KCl]: 27 mM
[Na₂HPO₄]: 100 mM
[KH₂PO₄]: 18 mM

For 1 L:

80 g NaCl (Product 71376)
2 g KCl (Product P9541)
17.8 g Na₂HPO₄•2 H₂O (Product 71643 – FW: 177.99 g/mol)
2.45 g KH₂PO₄ (Product P9791 – FW: 136.09 g/mol)
Add to 900 mL of Milli-Q® water and adjust pH to 7.4.
Add Milli-Q® water to final volume of 1 L.

Transfer buffer 10X

[Tris]: 250 mM
[Glycine]: 1.9 M
[SDS]: 17 mM, 0.5% (w/v)

For 500 mL:

15.2 g Tris base (Product 93362 – FW: 121.1 g/mol)
72.1 g glycine (Product 50046)
2.5 g sodium dodecyl sulfate (SDS, Product L3771)
Add Milli-Q® water to final volume of 500 mL

Transfer buffer 1X

Make fresh each time

For 500 mL:

50 mL 10X transfer buffer
100 mL methanol (Product 494437)
Add Milli-Q® water to final volume of 500 mL

SDS-PAGE Coomassie® staining solution

[Coomassie® Blue R-250]: 3.0 mM
[Methanol]: 50% (v/v)
[Acetic acid]: 1.74 M, 10% (v/v)

For 500 mL:

1.25 g Coomassie® Blue R-250 (Product B7920)
225 mL methanol (Product 494437)
225 mL Milli-Q® water
50 mL glacial acetic acid (Product 45726)

SDS-PAGE destaining solution

[Methanol]: 30% (v/v)
[Acetic acid]: 1.74 M, 10% (v/v)

For 1 L:

300 mL methanol (Product 494437)
100 mL glacial acetic acid (Product 45726)
600 mL Milli-Q® water

Recipes for mPAGE™ Precast Gel System

mPAGE™ MES SDS Running Buffer

[Tris]: 50 mM
[MES]: 50 mM
[SDS]: 0.1%, 3.5 mM
[EDTA]: 1 mM

For 1000 mL:

6.06 g Tris base (Product 93362 – FW: 121.1 g/mol)
9.76 g 2-(N-Morpholino)ethanesulfonic acid
(MES, Product M3671)
1.0 g sodium dodecyl sulfate (SDS, Product L3771)
0.3 g Ethylenediaminetetraacetic acid (EDTA – Product 03609)

Dissolve in 900 mL Milli-Q® water and mix well.
Add Milli-Q® water to a final volume of 1 L. Do not adjust pH.

mPAGE™ MOPS SDS Running Buffer

[Tris]: 50 mM
[MOPS]: 50 mM
[SDS]: 0.1%, 3.5 mM
[EDTA]: 1 mM

For 1000 mL:

6.06 g Tris base (Product 93362 – FW: 121.1 g/mol)
1.046 g 3-(N-morpholino)propanesulfonic acid (MOPS,
Product 69947)
1.0 g sodium dodecyl sulfate (SDS, Product L3771)
0.3 g Ethylenediaminetetraacetic acid (EDTA – Product 03609)

Dissolve in 900 mL Milli-Q® water and mix well.
Add Milli-Q® water to a final volume of 1 L. Do not adjust pH.

mPAGE™ 4X LDS Sample Buffer

[Tris-HCl]: 424 mM
[Tris]: 564 mM
[LDS]: 8%, 294 mM
[EDTA]: 2.07 mM
[Glycerol]: 40%, 4.34 M
[Coomassie® Blue G250]: 0.88 mM
[Phenol Red]: 0.7 mM

For 10 mL:

0.666 g Tris-HCl (Product 93363 - FW:157.60 g/mol)
0.682 g Tris base (Product 93362 – FW: 121.1 g/mol)
0.800 g Lithium dodecyl sulfate (LDS, Product L9781)
0.006 g Ethylenediaminetetraacetic acid (EDTA, Product
03609)
3.2 mL Glycerol (Product G5516)
0.75 mL 1% Coomassie® Blue G250 solution
0.25 mL 1% Phenol Red solution

Combine and add Milli-Q® water to final volume 10 mL.
Store at 2-8 °C. The pH of the 1X solution is 8.5. Do not adjust
pH with acid or base

mPAGE™ 1X Transfer Buffer, pH 8.2, for Wet Transfer

[Tris]: 25 mM
[Bicine]: 25 mM
[Methanol]: 10%

For 1000 mL:

3.0 g Tris base (Product 93362 – FW: 121.1 g/mol)
4.08 g N,N-Bis(2-hydroxyethyl)glycine (Bicine, Product B8660)
100 mL methanol (Product 494437)

Combine and mix well. Add Milli-Q® water to final volume
1000 mL. Do not adjust pH.

mPAGE™ Transfer Buffer (with Methanol), pH 8.2, for Semi-Dry Transfer

[Tris]: 50 mM
[Bicine]: 50 mM
[Methanol]: 10%

For 500 mL:

3.0 g Tris base (Product 93362 – FW: 121.1 g/mol)
4.08 g N,N-Bis(2-hydroxyethyl)glycine (Bicine, Product B8660)
50 mL methanol (Product 494437)

Combine and mix well. Add Milli-Q® water to final volume
500 mL. Do not adjust pH.

mPAGE™ Gel Equilibration Buffer, pH 8.2, for Semi-Dry Transfer

[Tris]: 50 mM
[Bicine]: 50 mM

For 500 mL:

3.0 g Tris base (Product 93362 – FW: 121.1 g/mol)
4.08 g N,N-Bis(2-hydroxyethyl)glycine (Bicine, Product B8660)

Dissolve in 450 mL of Milli-Q® water and mix well. Add Milli-Q®
water to final volume 500 mL. Do not adjust pH.

Troubleshooting Blotting Problems

1. Dot/Slot (Filtration) Blotting

Symptom	Possible Cause	Remedy
Slow or no filtration of the sample through the membrane	Inadequate vacuum	Make sure the blotting unit is closed properly and the seal is intact. Make sure the vacuum source (e.g. pump) is operating properly. Seal off any open wells with a high-quality laboratory tape. Increase vacuum level.
	Membrane pores clogged	Centrifuge or filter samples to remove particulates. Dilute viscous samples with buffer.
Little or no protein observed on the blot	Not enough protein applied to the membrane	Minimize sample dilution and filter more sample through the membrane.
	Detergents (e.g., SDS) may inhibit lower molecular weight proteins from binding to the membrane	Eliminate detergents if possible.
	Stain not sensitive enough	Use a more sensitive stain.
Stained blot is not uniform	Membrane structure was compressed by filter paper	Place a second membrane in the blotting unit to protect the membrane receiving the samples.
	Air bubbles trapped in the interior of the membrane	Pre-wet membrane by laying it on the surface of the methanol. Immersing the membrane can entrap air.
	Membrane not pre-wet in methanol	Membrane must be pre-wet with methanol; entire membrane should change uniformly from opaque to semi-transparent.
	Air bubbles in the sample	Carefully pipette samples into well to avoid the formation of air bubbles.
	Not enough sample volume loaded	Sample must cover the entire exposed membrane area.
Protein smeared across the top of the membrane	Sample leaked across the wells	Make sure the blotting unit is properly assembled, closed and sealed prior to filtration.
Protein smeared across the back of the membrane	Membrane capacity was exceeded	Reduce the amount of protein loaded into the well.

2. Semi-dry or Tank Electrotransfer

Symptom	Possible Cause	Remedy
Band smeared/distorted	Membrane not uniformly wetted with methanol	The entire membrane must be pre-wet with methanol; the entire membrane should change uniformly from opaque to semi-transparent.
	Air bubbles under membrane and between other layers in the stack	Using a pipette or stirring rod, gently roll out any trapped air bubbles while assembling the stack.
	Uneven contact between gel and membrane	Make sure entire gel and membrane surfaces are in good contact.
	Too much heat generated during the transfer	The temperature of the run should not exceed 20°C. For a tank transfer, pre-chill the buffer or carry out the transfer in a cold room. For a semi-dry transfer, either shorten the run time, increase the number of filter papers, or reduce the current.
	Filter paper dried out during semi-dry transfer	Make sure filter paper is thoroughly drenched prior to transfer or use additional sheets. Be sure the stack is assembled in less than 15 minutes.
	Proteins transferred too rapidly; protein buildup on the membrane surface	Reduce the strength of the electrical field.

2. Semi-dry or Tank Electrotransfer (continued)

Symptom	Possible Cause	Remedy
Weak signal	Proteins passing through the membrane	<p>Increase the time the proteins have to interact with membrane by reducing the voltage by as much as 50%.</p> <p>Highly negatively charged proteins (due to high aspartic acid and glutamic acid content) tend to move very fast in an electric field. Decrease the voltage to slow down migration of these proteins.</p> <p>Presence of SDS in the gel may inhibit protein-binding. Equilibrate the gel in the transfer buffer for at least 15 minutes.</p> <p>Methanol concentration in transfer buffer is too low to facilitate removal of SDS. Increase the methanol to 15 – 20%, especially for smaller molecular weight proteins.</p> <p>The membrane must be pre-wet with methanol; the entire membrane should change uniformly from opaque to semi-transparent.</p> <p>Switch to Immobilon®-P^{5Q} transfer membrane.</p>
	Proteins retained in the gel	If the methanol concentration in the transfer buffer is too high, it can remove SDS from proteins and lead to protein precipitation in the gel. This would reduce the transfer of large molecular weight proteins out of the gel. If protein precipitation is an issue, the transfer buffer can be supplemented with SDS (0.01% – 0.05%) to aid in solubility. In addition, excess methanol can tend to shrink or tighten a gel, thus inhibiting transfer of large molecular weight proteins.
	Isoelectric point of the protein is at or close to the pH of the transfer buffer	A protein that has the same isoelectric point as the pH of the transfer buffer will have no net charge and thus will not migrate in an electric field. To facilitate transfer, try a higher pH buffer such as 10 mM CAPS buffer at pH 11, including 10% methanol or a lower pH buffer such as an acetic acid buffer.
	Poor detection when urea is used in the gel and/or transfer buffer	Reduce the temperature by using a circulating buffer setup or run your transfer in a cold room. Urea in the presence of heat can cause carbamylation of proteins, which can change the charge of amino acids in a protein. This could affect the epitopes essential for antibody recognition and binding.
	Incomplete transfer of proteins	Stain the gel to check for residual proteins. If transfer was not complete, review your transfer technique.
	Poor protein retention	Once transfer is complete, be sure to dry the membrane completely to obtain optimal binding and fixation of the proteins. This should be done prior to any downstream detection method.
No signal	No transfer of proteins	Check for the gel and membrane orientation during the transfer process. Use pre-stained molecular weight standards to monitor transfer.
Poor transfer of small molecular weight proteins	Insufficient protein retention SDS interferes with binding of small molecular weight proteins	Switch to Immobilon®-P ^{5Q} transfer membrane. Remove SDS from the transfer solution.
	Low methanol concentration in the transfer buffer	Use higher percentage of methanol (15% – 20%) in the transfer buffer.
	Insufficient protein binding time	A lower voltage may optimize binding of small proteins to the membrane.
	Current doesn't pass through the membrane	Cut membrane and blotting paper exactly to the gel size; do not allow overhangs.
Poor transfer of large molecular weight proteins (~ > 80 kDa)	Methanol concentration is too high	Reducing the methanol concentration to 10% (v/v) or less should help in the transfer of large molecular weight proteins by allowing the gel to swell. Moreover, a lower methanol percentage would also reduce SDS loss from the proteins and reduce protein precipitation in the gel. Proteins > 200 kDa are not as sensitive to interference from the SDS in binding to membrane as are proteins < 100 kDa.
Poor transfer of positively charged proteins	Protein net charge in the transfer buffer is positive; proteins move toward cathode	Reverse the transfer stack such that the Immobilon® transfer membrane is on the cathode side of the gel.
Poor semi-dry transfer	Current bypasses the gel stack	Make sure the membrane and blotting paper are cut exactly to the gel size and there are no overhangs.
Poor transfer of a wide range of protein sizes	Different conditions required to transfer large and small proteins	<p>Refer to A two-step procedure for efficient electrotransfer of both high molecular weight (> 400,000) and low molecular weight (<20,000) proteins²².</p> <p>Use three-buffer system for semi-dry transfer (Protocol 1.2)</p>

3. Protein Visualization

Symptom	Possible Cause	Remedy
Poor detection by transillumination	Inappropriate membrane	Transillumination works best with Immobilon®-P transfer membrane. It is not recommended for nitrocellulose or Immobilon®-P ^{SQ} transfer membrane.
	Membrane wasn't completely dried prior to wetting with methanol	Be sure that the membrane was dried completely after the transfer prior to immersing it in the 20% methanol solution. Make sure to use a 20% methanol solution.
	Blot saturated with water only	Saturate the blot with 20% methanol.
	Insufficiently saturated blot	Increase concentration of methanol up to 40%.
Weak or uneven stain	Membrane wasn't wetted in methanol prior to staining	The membrane must be pre-wet with methanol; the entire membrane should change uniformly from opaque to semi-transparent.
Uneven/splotchy results	Use sufficient volume of incubation solutions and ensure that the membrane is completely covered with these solutions during incubation.	The container used should be large enough to allow solution to move freely across the blot. Do not incubate more than one blot at a time in that same container. In addition, the protein side of the blot should be facing up so as not to be interacting with the bottom surface of the container.
	Air bubbles	The blot should not have any air bubbles on the surface. Gently pull the membrane across the edge of the container to remove bubbles.
	Poor reagent quality	All of the buffers and reagents should be fresh and free of particulates and contaminants. Filtration of buffers with Millex® syringe filters or Stericup®, Steritop® or Steriflip® filter units and centrifugation of antibody stocks may be required.
High background staining	Nonspecific protein binding to the membrane	Make sure to use clean electrotransfer equipment and components and high-quality reagents and Milli-Q® water.

4. Immunodetection

Symptom	Possible Cause	Remedy
Weak signal	Improper blocking reagent	The blocking agent may have an affinity for the protein of interest and thus obscure the protein from detection. Try a different blocking agent and/or reduce both the amount or exposure time of the blocking agent.
	Insufficient antibody reaction time	Increase the incubation time.
	Antibody concentration is too low or antibody is inactive	Multiple freeze-thaw or bacterial contamination of antibody solution can change antibody titer or activity. Increase antibody concentration or prepare it fresh.
	Outdated detection reagents	Use fresh substrate and store properly. Outdated substrate can reduce sensitivity.
	Protein transfer problems	Optimize protein transfer (see above).
	Dried blot in chromogenic detection	If there is poor contrast using a chromogenic detection system, the blot may have dried. Try rewetting the blot in water to maximize the contrast.
	Tap water inactivates chromogenic detection reagents	Use Milli-Q® water for reagent preparation.
	Azide inhibits HRP	Do not use azide in the blotting solutions.
	Antigen concentration is too low	Load more antigen on the gel prior to the blotting.
No signal	Antibody concentration too low	Increase concentration of primary and secondary antibodies.
	HRP inhibition	HRP-labeled antibodies should not be used in solutions containing sodium azide.
	Primary antibody was raised against native protein	Separate proteins in non-denaturing gel or use antibody raised against denatured antigen.
Uneven blot	Fingerprints, fold marks or forceps imprints on the blot	Avoid touching or folding membrane; use gloves and blunt end forceps.
Speckled background	Aggregates in the blocking reagent	Filter blocking reagent solution through 0.2 µm or 0.45 µm Millex® syringe filter unit.
	Aggregates in HRP-conjugated secondary antibody	Filter secondary antibody solution through 0.2 µm or 0.45 µm Millex® syringe filter unit.

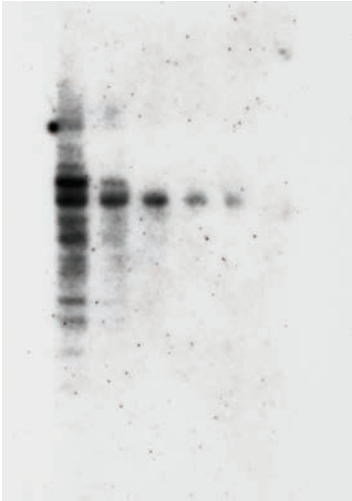
4. Immunodetection (continued)

Symptom	Possible Cause	Remedy
High background	Insufficient washes	Increase washing volumes and times. Pre-filter all of your solutions, including the transfer buffer, using Millex [®] syringe filter units , Stericup [®] , or Steriflip [®] filter units.
	Secondary (enzyme-conjugated) antibody concentration is too high	Increase antibody dilution.
	Protein-protein interactions	Use Tween [®] -20 detergent (0.05%) in the wash and detection solutions to minimize protein-protein interactions and increase the signal to noise ratio.
	Immunodetection on Immobilon [®] -P ^{5Q} transfer membrane	Increase the concentration or volume of the blocking agent used to compensate for the greater surface area of the membrane. Persistent background can be reduced by adding up to 0.5M NaCl and up to 0.2% SDS to the wash buffer and extending the wash time to 2 hours.
	Poor-quality reagents	Use high-quality reagents and Milli-Q [®] water.
	Cross-reactivity between blocking reagent and antibody	Use different blocking agent or use Tween [®] -20 detergent in the washing buffer.
	Film overexposure	Shorten exposure time.
	Membrane drying during incubation process	Use volumes sufficient to cover the membrane during incubation.
	Poor-quality antibodies	Use high-quality affinity purified antibodies.
Excess detection reagents	Drain blots completely before exposure.	
Persistent background	Nonspecific binding	Use High Salt Wash. (Protocol 3.5)
High background (rapid immunodetection)	Membrane wets out during rapid immunodetection	Reduce the Tween [®] -20 (< 0.04%) detergent in the antibody diluent. Use gentler agitation during incubations. Rinse the blot in Milli-Q [®] water after electrotransfer to remove any residual SDS carried over from the gel. Be sure to dry the blot completely prior to starting any detection protocol.
	Membrane was wet in methanol prior to the immunodetection	Do not pre-wet the membrane.
	Membrane wasn't completely dry prior to the immunodetection	Make sure the membrane is completely dry prior to starting the procedure.
Nonspecific binding	Primary antibody concentration too high	Increase primary antibody dilution.
	Secondary antibody concentration too high	Increase secondary antibody dilution.
	Antigen concentration too high	Decrease amount of protein loaded on the gel.
Reverse images on film (white bands on dark background)	Too much HRP-conjugated secondary antibody	Reduce concentration of secondary HRP-conjugated antibody.
Poor detection of small proteins	Small proteins are masked by large blocking molecules such as BSA	Consider casein or a low molecular weight polyvinylpyrrolidone (PVP). Surfactants such as Tween [®] detergent and Triton [®] X-100 detergent may have to be minimized.
		Avoid excessive incubation times with antibody and wash solution.

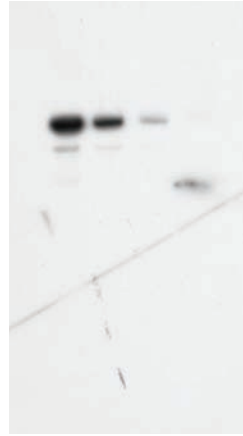
5. Fluorescent Detection

Symptom	Possible Cause	Remedy
High overall background	High background fluorescence from the blotting membrane	Use Immobilon®-FL PVDF blotting membrane.
	Blocking reagents not optimized for fluorescence detection	Use Block-FL noise-cancelling reagents, optimized for fluorescent Western blots.
Multiplexing problems	Experimental design	The two antibodies must be derived from different host species so that they can be differentiated by secondary antibodies of different specificities. Before combining the two primary antibodies, test the banding pattern on separate blots to determine where bands will appear. Use cross-adsorbed secondary antibodies in two-color detection.
Speckled background	Dust/powder particles on the Handle blots	Handle blots with powder-free gloves and clean surface of the scanner.
Low signal	Wet blot	Drying the blot may enhance signal strength. The blot can be scanned after re-wetting. Do not wrap the blot in plastic/Saran™ wrap while scanning.
	Blot photo-bleached	While fluorescent dyes usually provide long-lasting stable signal, some fluorescent dyes can be easily photo-bleached. To prevent photo-bleaching, protect the membrane from light during secondary antibody incubations and washes, and until the membrane is ready to be scanned. Store developed blots in the dark for subsequent imaging.
	Wrong excitation wavelength	Follow dye manufacturers' instructions for blot imaging or use an emission filter.

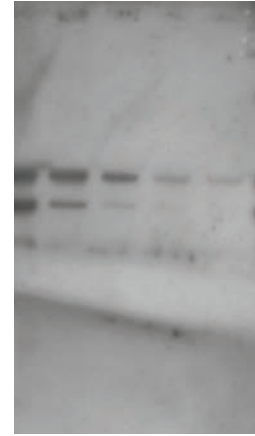
Examples and Causes of Blot Failure



Fingerprints

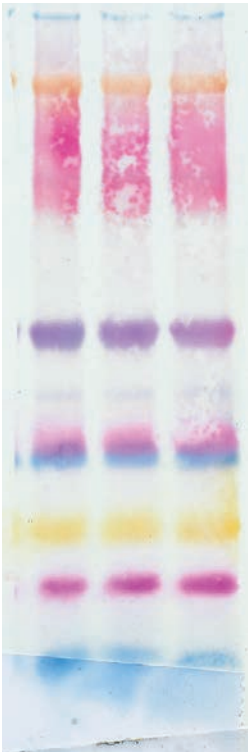


Scratches and forceps imprints

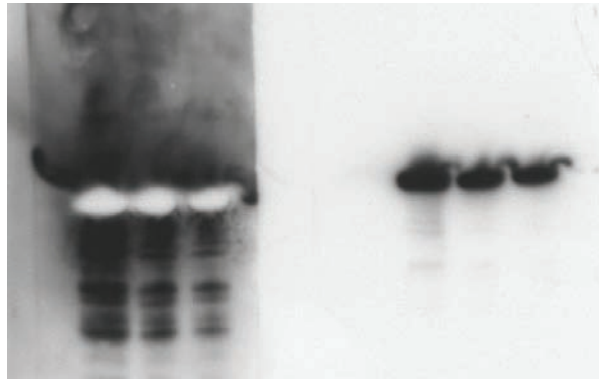


Folded membrane

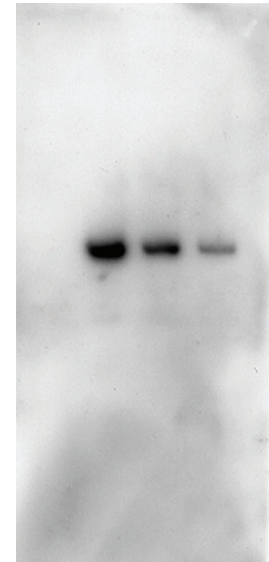
Spotchy background caused by insufficient washes and/or unfiltered blocking solution. Can be improved by filtering the blocking solution through 0.2 μm Millex®-GP syringe filter and adding extra washes.



Bubbles between the membrane and SDS-PAGE gel introduced during the transfer.



Negative (reverse) image on film caused by excess secondary HRP-conjugated antibody. The detection was improved by increasing secondary antibody dilution 10-fold. Blot on left was incubated with a 1:5,000 dilution of secondary antibody, and blot on right was incubated with a 1:50,000 dilution.



High background and dark edges caused by insufficient washing of the blot. Can be alleviated by additional or/and longer washes or a high salt wash (Protocol 3.5)

Glossary

1-D

One-dimensional

2-D

Two-dimensional

Adsorption

The process whereby molecules or particles (e.g., protein) bind to a surface (e.g., membrane).

Anode

Positively charged electrode in an electrophoresis system.

Blocking

Technique used to reduce nonspecific binding of antibodies during immunodetection; unoccupied membrane sites are blocked with an inert protein, nonionic detergent, or protein-free blocking reagent.

Blot

A microporous membrane with biomolecules adsorbed onto the polymer.

Blotting

Process of transferring proteins, nucleic acids or biomolecules from a gel to a membrane. A membrane with proteins immobilized on it is called a Western blot.

Cathode

Negatively charged electrode in an electrophoresis system.

Chemiluminescent detection

Immunodetection technique that results in the production of visible light at the site of the target protein.

Chromogenic detection

Immunodetection technique that results in the deposition of a colored precipitate at the site of the target protein.

Dot blot

A blot prepared by filtration of liquid samples through a membrane using a dot blot manifold (see "Filtration").

Edman degradation

A process that uses the Edman reagent, phenyl isothiocyanate (PITC), to sequentially remove one amino acid from the N-terminus of a protein. The chemically derivatized amino acid is then analyzed after it is cleaved from the protein. Sequential processing of the protein provides the amino acid sequence of the protein.

Electrotransfer

Common method for the transfer of proteins from a gel to a membrane. Proteins elute from the gel and get absorbed onto the membrane when an electrical field is applied perpendicular to the plane of the gel.

ELISA

Enzyme-Linked Immunosorbent Assay; a rapid method to determine the presence and quantity of a specific substance. It is based on an antigen-antibody interaction where the antibody or antigen is linked to an enzyme as a means of detecting its presence.

Filtration

Direct application of sample onto a membrane.

A dissolved sample is pulled through the membrane by applying a vacuum; proteins bind to the membrane and the other sample components pass through.

Fluorescent detection

Immunodetection of a protein using a primary or secondary antibody conjugated to a fluorescent dye.

Gel

The substrate, usually polyacrylamide, on which sample proteins have been separated.

Immunoblot

A Western blot that has been analyzed for a target protein using a specific antibody.

Immunodetection

Method of protein detection using a specific antibody to identify the location of a target membrane-bound protein. The specificity of antigen-antibody binding permits the identification of a specific protein in a complex sample.

Isoelectric focusing

Method of protein separation on the basis of isoelectric points. Usually achieved by electrophoresis of proteins in a stabilized pH gradient where proteins migrate to the pH corresponding to their isoelectric points.

Isoelectric point (pI)

The pH value at which the net electric charge of a molecule, such as a protein or amino acid, is zero. It is usually the first dimension of a 2-D gel.

Polyacrylamide

A crosslinked polymer of acrylamide used in gel electrophoresis.

Primary antibody

The first antibody used in an immunodetection protocol. The primary antibody is specific for the target antigen.

PVDF

Polyvinylidene fluoride; the polymer used to make Immobilon®-P, Immobilon®-P^{5Q} and Immobilon®-FL transfer membranes. (Occasionally this acronym is erroneously defined in the blotting literature as polyvinylidene difluoride.)

Rapid immunodetection

Faster method of immunodetection which eliminates the need for (and the risks of) blocking.

Reprobing

The process of sequentially cycling a blot through multiple rounds of immunodetection.

Retention

In the context of stripping and reprobing immunoblots, retention refers to the ability of a protein to remain adsorbed to a membrane surface under conditions that disrupt immunocomplexes.

SDS

Sodium dodecyl sulfate. SDS is a detergent that binds to proteins, giving them a net negative charge. It is used in denaturing protein gel electrophoresis. It is also widely used to disrupt cell walls and dissociate protein complexes. Lithium dodecyl sulfate (LDS) can be substituted for SDS.

SDS-PAGE

Sodium dodecyl sulfate-polyacrylamide gel electrophoresis

Secondary antibody

The second antibody used in an immunodetection protocol. The secondary antibody is specific for the primary antibody and is typically conjugated to an enzyme used for signal amplification.

Semi-dry transfer

Electrotransfer technique where the traditional buffer reservoir is replaced by layers of filter paper soaked in buffer; an equally effective, but faster technique than tank transfer. Multiple buffer systems can be used during the transfer.

Slot blot

A blot prepared by the filtration of liquid samples through a membrane using a slot blot manifold.

Dot blotting

A blot prepared by manually applying sample to the membrane.

Staining

Technique used to make protein bands visible on a gel or blot. The colored stain may be reversible or non-reversible.

Stripping

Process of removing an antibody from a membrane prior to a subsequent round of immunodetection.

Substrate

The compound that interacts with the enzyme conjugated to the secondary antibody to yield a detectable signal in immunodetection. The signal is typically in the form of a colored precipitate (chromogenic substrates) or electromagnetic radiation (light from chemiluminescent or fluorescent substrates).

Tank transfer

Traditional electrotransfer technique where the gel and membrane are completely immersed in a reservoir of buffer; an effective but slow technique.

Transfer buffer(s)

The buffer(s) used as the chemical environment to facilitate the transfer of biomolecules from a gel onto a membrane.

Transillumination

Non-destructive, reversible visualization technique used to make protein bands visible on a blot. The protein bands appear as clear areas in dark field when placed on a light box.

Vacuum blotting

The process of transferring biomolecules from a gel to a membrane using vacuum as the driving force; not typically used for protein gels.

Western blot

A blot prepared by transferring protein from a polyacrylamide gel to a membrane.

References

1. Southern, E.M. Detection of specific sequences among DNA fragments separated by gel electrophoresis. *J Mol Biol.* 1975, 5;98(3):503-17.
2. Pluskal, M.; Przekop, M.; Kavonian, M.; Hicks, D.; Vecoli, C. Immobilon® PVDF transfer membrane: a new membrane substrate for Western blotting. *Biotechniques.* 1986;4(3):272-82.
3. Kurien BT, Scofield RH. Protein blotting: a review. *J Immunol Methods.* 2003; 274(1-2):1-15.
4. Lauritzen, E. and Pluskal, M. Improved HIV antiglycoprotein antibody detection by immunoblotting on a hydrophobic membrane. *J Acquir Immune Defic Syndr.* 1988;1(4):333-9.
5. Mansfield, M. Protein blotting using polyvinylidene fluoride membranes. In: Dunbar B, editor. *Protein blotting: a practical approach.* Oxford: IRL Press; 1994. p 33-52.
6. Matsudaira, P. Sequence from picomole quantities of proteins electroblotted onto polyvinylidene difluoride membranes. *Biol Chem.* 1987;262(21):10035-8.
7. Iwamatsu, A. S-carboxymethylation of proteins transferred onto polyvinylidene difluoride membranes followed by in situ protease digestion and amino acid sequencing. *Electrophoresis.* 1992;13:142-7.
8. Fernandez, J.; DeMott, M.; Atherton, D.; Mische, S.M. Internal protein sequence analysis: Enzymatic digestion for less than 10 µg of protein bound to polyvinylidene difluoride or nitrocellulose membranes. *Anal Biochem.* 1992;201(2):255-64.
9. McKeon, T.A. and Lyman, M.L. Calcium improves electrophoretic transfer of calmodulin and other small proteins. *Anal Biochem.* 1991; 193:125-30.
10. Reese, G.; Schmechel, D.; Ayuso, R.; Lehrer, S.B. Grid-immunoblotting: a fast and simple technique to test multiple allergens with small amounts of antibody for cross-reactivity. *J Chromatogr B Biomed Sci Appl.* 2001;756(1-2):151-6.
11. Celis, J.E. and Gromov, P. High-resolution two-dimensional gel electrophoresis and protein identification using Western blotting and ECL detection. *EXS.* 2000;88:55-67.
12. Poland, J.; Bohme, A.; Schubert, K.; Sinha, P. Revisiting electroblotting of immobilized pH gradient gels: a new protocol for studying post-translational modification of proteins. *Electrophoresis.* 2002;23(24):4067-71.
13. Eto, M.; Watanabe, K.; Moriyama, T.; Makino, I. Apolipoprotein E phenotyping from plasma by isoelectric focusing and immunoblotting. *Tohoku J Exp Med.* 1990;160(4):301-9.
14. Kurien, B.T. and Scofield, R.H. Multiple immunoblots after non-electrophoretic bidirectional transfer of a single SDS-PAGE gel with multiple antigens. *J Immunol Methods.* 1997;205(1):91-4.
15. Chen, H. and Chang G.D. Simultaneous immunoblotting analysis with activity gel electrophoresis in a single polyacrylamide gel. *Electrophoresis.* 2001;22(10):1894-9.
16. Peferoen, M.; Huybrechts, R.; De Loof, A. Vacuum-blotting: a new simple and efficient transfer of proteins from sodium dodecyl sulfate-polyacrylamide gels to nitrocellulose. *FEBS Letters.* 1982;145(2):369-72.
17. Towbin, H.; Staehelin, T.; Gordon, J. Electrophoretic transfer of proteins from polyacrylamide gels to nitrocellulose sheets: procedure and some applications. *Proc Natl Acad Sci.* 1979; 76(9):4350-4.
18. Kyhse-Anderson, J. Electroblotting of multiple gels: a simple apparatus without buffer tank for rapid transfer of proteins from polyacrylamide to nitrocellulose. *J Biophys Biochem Methods.* 1984;10:203-209.
19. Schafer-Nielsen, C.; Svendsen, P.J.; Rose, C. Separation of macromolecules in isotachopheresis involving single or multiple counterions. *J Biochem Biophys Methods.* 1980; 3(2):97-128.
20. Mozdzanowski, J.; Speicher, D.W. Microsequence analysis of electroblotted proteins. *Anal Biochem.* 1992;207(1):11-8.
21. Towbin, H. and Gordon, J. Immunoblotting and dot immunoblotting – current status and outlook. *J Immunol Methods.* 1984;72(2):313-40.
22. Otter, T.; King, S.M.; Witman, G.B. A two-step procedure for efficient electrotransfer of both high molecular weight (> 400,000) and low molecular weight (<20,000) proteins. *Anal Biochem.* 1987;162:370-7.
23. Bolt, M.W. and Mahoney, P.A. High-efficiency blotting of proteins of diverse sizes following sodium dodecyl sulfate-polyacrylamide gel electrophoresis. *Anal Biochem.* 1997;247(2):185-92.
24. Erickson P.F.; Minier, L.N.; Lasher, R.S. Quantitative electrophoretic transfer of polypeptides from SDS polyacrylamide gels to nitrocellulose sheets: a method for their re-use in immunoradiographic detection of antigens. *J Immunol Methods.* 1982;51(2):241-9.
25. Elkon, K.B.; Jankowski, P.W.; Chu, J.L. Blotting intact immunoglobulins and other high molecular weight proteins after composite agarose-polyacrylamide gel electrophoresis. *Anal Biochem.* 1984;140(1):208-13.
26. Thompson, D. and Larson, G. Western blots using stained protein gels. *BioTechniques.* 1992; 12(5):656-8.
27. Reig, J.A. and Klein, D.C. Submicron quantities of unstained proteins are visualized on polyvinylidene difluoride membranes by transillumination. *Applied and Theoretical Electrophoresis.* 1988;1:59-60.
28. Berggren, K.; Steinberg, T.H.; Lauber, W.M.; Carroll, J.A.; Lopez, M.F.; Chernokalskaya, E.; Zieske, L.; Diwu, Z.; Haugland, R.P.; Patton, W.F. A luminescent ruthenium complex for ultrasensitive detection of proteins immobilized on membrane supports. *Anal Biochem.* 1999;276(2):129-143.
29. Haugland, R.P. *Handbook of fluorescent probes and research products.* 9th edition. Eugene (OR): Molecular Probes, Inc.; 2002.
30. Dunn, M.J. Detection of total proteins on Western blots of 2-D polyacrylamide gels. *Methods Mol Biol.* 1999;112:319-29.
31. Bickar, D. and Reid, P.D. A high-affinity protein stain for Western blots, tissue prints, and electrophoretic gels. *Anal Biochem.* 1992; 203(1):109-15.
32. Van Dam, A.P.; Van den Brink, H.G.; Smeenk, R.J.T. Technical problems concerning the use of immunoblots for the detection of antinuclear antibodies. *J Immunol Methods.* 1990;129(1):63-70.
33. Zampieri, S.; Ghirardello, A.; Doria, A.; Tonello, M.; Bendo, R.; Rossini, K.; Gambari, P.F. The use of Tween®-20 in immunoblotting assays for the detection of autoantibodies in connective tissue diseases. *J Immunol Methods.* 2000; 239(1-2):1-11.
34. Sharma, S.K. and Carew, T.J. Inclusion of phosphatase inhibitors during Western blotting enhances signal detection with phosphospecific antibodies. *Anal Biochem.* 2002; 307(1):187-9.
35. Lasne, F. Double blotting: a solution to the problem of nonspecific binding of secondary antibodies in immunoblotting procedures. *J Immunol Methods.* 2001;253:125-131.
36. Lasne, F. Double blotting: a solution to the problem of nonspecific binding of secondary antibodies in immunoblotting procedures. *J Immunol Methods.* 2003;276:223-6.
37. Leary, J.J.; Brigati, D.J.; Ward, D.C. Rapid and sensitive colorimetric method for visualizing biotin-labeled DNA probes hybridized to DNA or RNA immobilized on nitrocellulose: bio-blots. *Proc Natl Acad Sci.* 1983;80(13):4045-9.
38. Cortese, J.D. Beyond film: laboratory imagers. *The Scientist.* 2002;16(7):41.
39. Hengen, P.N. Chemiluminescent detection methods. *TIBS.* 1997; 22:313-14.
40. Van Dyke, K.K.; Van Dyke, R., editors. *Luminescence immunoassay and molecular applications,* Boca Raton (FL): CRC Press; 1990. p 61-5.
41. Gharahdaghi, F.; Kirchner, M.; Fernandez, J.; Mische, S.M. Peptide-mass profiles of polyvinylidene difluoride-bound proteins by matrix-

- assisted laser desorption/ionization time-of-flight mass spectrometry in the presence of nonionic detergents. *Anal Biochem.* 1996;233(1):94–9.
42. Bienvenut, W.V.; Sanchez, J.C.; Karmime, A.; Rouge, V.; Rose, K.; Binz, P.A.; Hochstrasser, D.F. Toward a clinical molecular scanner for proteome research: parallel protein chemical processing before and during Western blot. *Anal Chem.* 1999;71(21):4800–7.
43. Bunai, K.; Nozaki, M.; Hamano, M.; Ogane, S.; Inoue, T.; Nemoto, T.; Nakanishi, H.; Yamane, K. Proteomic analysis of acrylamide gel separated proteins immobilized on polyvinylidene difluoride membranes following proteolytic digestion in the presence of 80% acetonitrile. *Proteomics.* 003;3(9):1738–49.
44. Binz, P.A.; Muller, M.; Walther, D.; Bienvenut, W.V.; Gras, R.; Hoogland, C.; Bouchet, G.; Gasteiger, E.; Fabbretti, R.; Gay, S.; et al. A molecular scanner to automate proteomic research and to display proteome images. *Anal Chem.* 1999;71(21):4981–8.
45. Bienvenut, W.V.; Déon, C.; Pasquarello, C.; Campbell, J.M.; Sanchez, J.C.; Vestal, M.L.; Hochstrasser, D.F. Matrix-assisted laser desorption/ionization-tandem mass spectrometry with high resolution and sensitivity for identification and characterization of proteins. *Proteomics.* 2002; 2(7), 868–76
46. Sloane, A.J.; Duff, J.L.; Wilson, N.L.; Gandhi, P.S.; Hill, C.J.; Hopwood, F.G.; Smith, P.E.; Thomas, M.L.; Cole, R.A.; Packer, N.H.; et al. High throughput peptide mass fingerprinting and protein macroarray analysis using chemical printing strategies. *Mol Cell Proteomics.* 2002;1(7):490–9.
47. Mansfield, M. Rapid immunodetection of polyvinylidene fluoride membrane blots without blocking. *Anal Biochem.* 1995;229(1):140–3.
48. Towbin, H.; Staehelin, T.; Gordon, J. Immunoblotting in the clinical laboratory. *J Clin Chem Clin Biochem.* 1989;27(8):495–501.
49. Stahl, D.; Lacroix-Desmazes, S.; Mouthon, L.; Kaveri, S.V.; Kazatchkine, M.D. Analysis of human self-reactive antibody repertoires by quantitative immunoblotting. *J Immunol Methods.* 2000; 240(1–2):1–14.
50. Bauer, G. Simplicity through complexity: immunoblot with recombinant antigens as the new gold standard in Epstein-Barr virus serology. *Clin Lab.* 2001;47(5–6):223–30.
51. Mylonakis, E.; Paliou, M.; Lally, M.; Flanigan, T.P.; Rich, J.D. Laboratory testing for infection with the human immunodeficiency virus: established and novel approaches. *Am J Med.* 2000; 109(7):568–76.
52. Heermann, K.H.; Gultekin, H.; Gerlich, W.H. Protein blotting: techniques and application in virus hepatitis research. *Ric Clin Lab.* 1988;18(2–3):193–221.
53. Ghaemmaghami, S.; Huh, W.K.; Bower, K.; Howson, R.W.; Belle, A.; Dephoure, N.; O'Shea, E.K.; Weissman, J.S. Global analysis of protein expression in yeast. *Nature.* 2003;425(6959):737–41
54. Rudolph, C.; Adam, G.; Simm, A. Determination of copy number of c-Myc protein per cell by quantitative Western blotting. *Anal Biochem.* 1999;269(1):66–71.
55. Wang, R. and Thompson, J.E. Detection of ATP competitive protein kinase inhibition by Western blotting. *Anal Biochem.* 2001;299:110–2.
56. Ahmed, F.E. Detection of genetically modified organisms in foods. *Trends Biotechnol.* 2002;20(5):215–23.
57. Grässer, F.A.; Sauder, C.; Haiss, P.; Hille, A.; König, S.; Göttel, S.; Kremmer, E.; Leinenbach, H.P.; Zeppezauer, M.; Mueller-Lantzsch, N. Immunological detection of proteins associated with the Epstein-Barr virus nuclear antigen 2A. *Virology.* 1993;195(2):550–60.
58. Silva, C.M.; Tully, D.B.; Petch, L.A.; Jewell, C.M.; Cidlowski, J.A. Application of a protein-blotting procedure to the study of human glucocorticoid receptor interactions with DNA. *Proc Natl Acad Sci.* 1987;84(7):1744–8.
59. Aksamitiene, E.; Hoek, J.B.; Kholodenko, B.; Kiyatkin, A. Multistrip Western blotting to increase quantitative data output. *Electrophoresis.* 2007;28(18):3163–3173.
60. Swank, M.W.; Kumar, V.; Zhao, J.; Wu, G.Y. A novel method of loading samples onto mini-gels for SDS-PAGE: Increased sensitivity and Western blots using sub-microgram quantities of protein. *J Neurosci Methods.* 2006;158:224–233.
61. Schilling, M.; Maiwald, T.; Bohl, S.; et al. Quantitative data generation for systems biology: the impact of randomisation, calibrators and normalisers. *Syst Biol (Stevenage).* 2005;152(4):193–200
62. Schilling, M.; Maiwald, T.; Bohl, S.; et al. Computational processing and error reduction strategies for standardized quantitative data in biological networks. *Febs J.* 2005;272(24):6400–6411.

General references

- Beisiegel, U. Protein blotting. *Electrophoresis.* 1986;7:1–18.
- Gershoni, J.M. Protein blotting: a manual. *Meth Biochem Analysis.* 1987;33:1–58.
- Stott, D.I. Immunoblotting and dot blotting. *J Immunol Methods.* 1989;119(2):153–87.

Patents

- U.S. Pat. No. 6,632,339, issued to W.V. Bienvenut et al. on October 14, 2003.
- U.S. Pat. App. Pub. No. 2002/0136668 (published September 26, 2002), filed by D. Wallace et al. on December 18, 2001.
- International Patent Application No. PCT/AU98/00265 of A. Gooley et al. (published October 22, 1998 as International Publication No. WO 98/47006).
- International Patent Application No. PCT/FR01/01331 of Laboratoire National de Dépistage du Dopage with Hospices Civils de Lyon which also holds the patent (2 786 273).

7. Ordering Information

Electrophoresis

Description	Cat. No.
mPAGE™ Precast Gels	
mPAGE™ 4-12% Bis-Tris, 10x8, 10-well 10 gels	MP41G10
mPAGE™ 4-12% Bis-Tris, 10x8, 12-well 10 gels	MP41G12
mPAGE™ 4-12% Bis-Tris, 10x8, 15-well 10 gels	MP41G15
mPAGE™ 4-20% Bis-Tris, 10x8, 10-well	MP42G10
mPAGE™ 4-20% Bis-Tris, 10x8, 12-well	MP42G12
mPAGE™ 4-20% Bis-Tris, 10x8, 15-well	MP42G15
mPAGE™ 8-16% Bis-Tris, 10x8, 10-well	MP81G10
mPAGE™ 8-16% Bis-Tris, 10x8, 12-well	MP81G12
mPAGE™ 8-16% Bis-Tris, 10x8, 15-well	MP81G15
mPAGE™ 8% Bis-Tris, 10x8, 10-well	MP8W10
mPAGE™ 8% Bis-Tris, 10x8, 12-well	MP8W12
mPAGE™ 8% Bis-Tris, 10x8, 15-well	MP8W15
mPAGE™ 10% Bis-Tris, 10x8, 10-well	MP10W10
mPAGE™ 10% Bis-Tris, 10x8, 12-well	MP10W12
mPAGE™ 10% Bis-Tris, 10x8, 15-well	MP10W15
mPAGE™ 12% Bis-Tris, 10x8, 10-well	MP12W10
mPAGE™ 12% Bis-Tris, 10x8, 12-well	MP12W12
mPAGE™ 12% Bis-Tris, 10x8, 15-well	MP12W15
mPAGE™ Buffers	
mPAGE™ 4X LDS Sample Buffer	MPSB-10ML
mPAGE™ 4X LDS Sample Buffer	MPSB-250ML
mPAGE™ MES SDS Running Buffer Powder	MPMES
mPAGE™ MOPS SDS Running Buffer Powder	MPM0PS
mPAGE™ Transfer Buffer Powder	MPTRB
mPAGE™ Trial Packs	
mPAGE™ Trial Kit, 4-12%, 12-well, MOPS	MP41G12TR1
mPAGE™ Trial Kit, 4-20%, 12-well, MOPS	MP42G12TR1
mPAGE™ Trial Kit, 10%, 12-well, MOPS	MP10W12TR1
mPAGE™ Trial Kit, 4-12%, 12-well, MES	MP41G12TR2
mPAGE™ Trial Kit, 10%, 12-well, MES	MP10W12TR2
mPAGE™ Trial Kit, 12%, 12-well, MES	MP12W12TR2
mPAGE™ Accessories	
mPAGE™ Adapter Plates	MPTA
mPAGE™ Gel Cassette Opener	MPC0
mPAGE™ Buffer Dam	MPBD
MW Markers	
mPAGE™ Unstained Protein Standard	MPSTD3
mPAGE™ Color Protein Standard	MPSTD4
mPAGE™ Western Protein Standard	MPSTD2
Biotinylated Molecular Weight Marker, mol wt 6,500–180,000 Da	B2787
Bovine Serum Albumin, dimer mol wt ~132 kDa	A8654
Bovine Serum Albumin, For use as a marker in SDS-PAGE	A7517
Bromophenol Blue sodium salt, for molecular biology, for electrophoresis	B8026
Color Marker Ultra-low Range (M.W. 1,060–26,600)	C6210
ColorBurst™ Electrophoresis Marker, mol wt 8,000–220,000 Da	C1992
ECL™ DualVue Western Markers, GE Healthcare, RPN810, sufficient for 25 blots	GERPN810
ECL Plex™ Fluorescent Rainbow™ Markers, GE Healthcare, RPN850E, pack of 120 µL	GERPN850E
Full Range Molecular Weight Marker Rainbow™, GE Healthcare, RPN800E	GERPN800E
High Range Molecular Weight Marker Rainbow™, GE Healthcare, RPN756E	GERPN756E
Low Range Molecular Weight Marker Rainbow™, GE Healthcare, RPN755E	GERPN755E
Prestained Molecular Weight Marker, mol wt 26,600–180,000 Da	SDS7B2
Protein Gel Stains	
ReadyBlue™ Protein Gel Stain, 1L	RSB-1L
Brilliant Blue G solution, Concentrate	B8522
Brilliant Blue R, pure	B7920
EZBlue™ Gel Staining Reagent	G1041
Fast Green FCF, Dye content ≥85 %	F7252
InstantBlue™, Ultrafast Protein Stain	ISB1L
SYPRO® Orange Protein Gel Stain	S5692-500UL
ProteoSilver™ Silver Stain Kit	PROTSIL1-1KT
ProteoSilver™ Plus Silver Stain Kit	PROTSIL2-1KT

Electrophoresis

Description	Cat. No.
Acrylamides	
Acrylamide, for electrophoresis, ≥99% (HPLC), powder	A3553
Acrylamide, suitable for electrophoresis, ≥99%	A8887
Acrylamide solution, 40%, for electrophoresis, sterile-filtered	A4058
N,N'Methylenebisacrylamide solution, 2%, for electrophoresis	M1533
Acrylamide/Bis-acrylamide, 30% solution, BioReagent, suitable for electrophoresis, 29:1	A3574
Acrylamide/Bis-acrylamide, 30% solution, BioReagent, suitable for electrophoresis, 37.5:1	A3699
Acrylamide/bis-acrylamide, 40% solution, BioReagent, suitable for electrophoresis, 19:1	A9926
Acrylamide/bis-acrylamide, 40% solution, BioReagent, suitable for electrophoresis, 29:1	A7802
Acrylamide/bis-acrylamide, 40% solution, BioReagent, suitable for electrophoresis, 37.5:1	A7168
Buffers	
BICINE, BioXtra, ≥99% (titration)	B8660
Ficoll® 400, BioXtra, for molecular biology, lyophilized powder	F2637
Glycine, for electrophoresis, ≥99%	G8898
Laemmli 2x Concentrate Sample Buffer	S3401
MES solution, BioPerformance Certified, 1 M, suitable for cell culture	M1317
Phenol Red, powder, BioReagent, suitable for cell culture	P3532
Sodium bisulfite, ACS reagent, mixture of NaHSO ₃ and Na ₂ S ₂ O ₅	243973
Trichloroacetic acid, for electrophoresis, suitable for fixing solution (for IEF and PAGE gels), ≥99%	T8657
Tricine, BioPerformance Certified, cell culture tested, ≥99% (titration)	T5816
Triethanolamine, ≥99.0% (GC)	90279
Tris-Glycine Buffer, 10x Concentrate	T4904
Tris-Glycine-SDS Buffer, 10x Concentrate	T7777
Tris-Tricine-SDS Buffer, 10x Concentrate	T1165
Trizma® hydrochloride solution, BioPerformance Certified, pH 7.5, 1 M, suitable for cell culture	T2319
Trizma® base, anhydrous, free-flowing, Redi-Dri™, ≥99.9%	RDD008
Detergents	
3-(Benzyltrimethylammonio)propanesulfonate, BioXtra, ≥99.0% (HPCE)	17236
IGEPAL® CA-630, CMC 0.083 mM	56741
Octyl β-D-glucopyranoside solution, ≥98% (HPLC), 50 % (w/v) in H ₂ O	O3757
Tween®-20 detergent	P1379
Gel Casting Reagents	
Ammonium persulfate, for molecular biology, for electrophoresis, ≥98%	A3678
1,4-Bis(acryloyl)piperazine, BioXtra, suitable for electrophoresis, ≥99.0% (TLC)	14470
(-)-Riboflavin, BioReagent, suitable for cell culture, suitable for insect cell culture, ≥98%	R9504
N,N,N',N'-Tetramethylethylenediamine, BioReagent, suitable for electrophoresis, ~99%	T9281
SDS Gel Preparation Kit	8091
Gel Cross-linkers	
N,N'-Bis(acryloyl)cystamine, BioReagent, suitable for electrophoresis	A4929
N,N'-Methylenebisacrylamide, powder, for molecular biology, for electrophoresis, ≥99.5%	M7279
N,N'-Methylenebisacrylamide solution, for electrophoresis, 2% in H ₂ O	M1533
Piperazine diacrylamide, For acrylamide gel electrophoresis	D1538
Reducing Agents	
DL-Dithiothreitol, for electrophoresis, ≥99%	D9163
Lithium dodecyl sulfate, BioReagent, for molecular biology, suitable for electrophoresis	L9781
Sodium dodecyl sulfate, BioReagent, suitable for electrophoresis, for molecular biology, ≥98.5% (GC)	L3771
5-Sulfosalicylic acid dihydrate, for electrophoresis, ≥99%	S3147
Tris(2-carboxyethyl)phosphine hydrochloride, BioUltra, for electrophoresis, SDS-PAGE tested	68957

Transfer

Description	Cat. No.
Blotting Papers	
Immobilon® Blotting Filter Paper, 7 x 8.4 cm sheet	IBFP0785C
Immobilon® Blotting Filter Paper, 8.5 x 13.5 cm sheet	IBFP0813C
Hybond blotting paper, L x W 20 cm x 20 cm, sheets, pack of 100 ea	GERPN6101M
Whatman® gel blotting paper, Grade GB003, GB003 blotting sheets, 15 x 20 cm, 100/pk	WHA10427812
Whatman® gel blotting paper, Grade GB003, GB003 blotting sheets, 20 x 20 cm, 100/pk	WHA10427818
Whatman® membrane marking pen	WHA10499001

Transfer

Description	Format	Cat. No.
PVDF Transfer Membrane		
Immobilon®-E PVDF Transfer Membrane, 0.45 µm (no alcohol pre-wet step required)	8.5 cm x 10 m roll	IEVH85R
	26.5 cm x 1.875 m roll	IEVH00005
	7 cm x 8.4 cm sheets, 50/pk	IEVH07850
	8 cm x 10 cm sheets, 10/pk	IEVH08100
	Blotting Sandwich*, 7 cm x 8.4 cm, 20/pk	IESN07852
	Blotting Sandwich*, 8.5 cm x 13.5 cm, 20/pk	IESN08132
	7 cm x 8.4 cm sheets, 4/pk (trial size)	IEVH07804
Immobilon®-P PVDF Transfer Membrane, 0.45 µm	8.5 cm x 10 m roll	IPVH85R
	26.5 cm x 3.75 m roll	IPVH00010
	26.5 cm x 1.875 m roll	IPVH00005
	7 cm x 8.4 cm sheets, 50/pk	IPVH07850
	8.5 cm x 13.5 cm, 10/pk	IPVH08130
	Blotting Sandwich*, 7 cm x 8.4 cm, 20/pk	IPSN07852
	Blotting Sandwich*, 8.5 cm x 13.5 cm, 20/pk	IPSN08132
Immobilon®-FL PVDF Transfer Membrane, 0.45 µm	8.5 cm x 10 m roll	IPFL85R
	26.5 cm x 3.75 m roll	IPFL00010
	26.5 cm x 1.875 m roll	IPFL00005
	7 cm x 8.4 cm sheets, 10/pk	IPFL07810
	10 cm x 10 cm sheets, 10/pk	IPFL10100
Immobilon®-PSQ PVDF Transfer Membrane, 0.2 µm	8.5 cm x 10 m roll	ISEQ85R
	26.5 cm x 3.75 m roll	ISEQ00010
	26.5 cm x 1.875 m roll	ISEQ00005
	7 cm x 8.4 cm sheets, 50/pk	ISEQ07850
	8.5 cm x 13.5 cm, 10/pk	ISEQ08130
	20 cm x 20 cm, 10/pk	ISEQ20200
Immobilon® NOW Dispenser	For 8.5 cm x 10 cm rolls	IMDISP

Description	Cat. No.
Protran® Nitrocellulose Membranes	
Amersham™ Protran® Western blotting membranes, nitrocellulose pore size 0.2 µm, roll W x L 300 mm x 4 m, pkg of 1 ea	GE10600001
Amersham™ Protran® Western blotting membranes, nitrocellulose pore size 0.2 µm, sheet W x L 200 mm x 200 mm, pkg of 25 ea	GE10600043
Amersham™ Protran® Western blotting membranes, nitrocellulose pore size 0.2 µm, sheet W x L 200 mm x 200 mm, pkg of 10 ea	GE10600044
Amersham™ Protran® Western blotting membranes, nitrocellulose pore size 0.2 µm, roll W x L 200 mm x 4 m, pkg of 1 ea	GE10600006
Amersham™ Protran® Premium Western blotting membranes, nitrocellulose pore size 0.2 µm, roll W x L 300 mm x 4 m, pkg of 1 ea	GE10600004
Amersham™ Protran® Supported Western blotting membranes, nitrocellulose pore size 0.2 µm, roll W x L 300 mm x 4 m, pkg of 1 ea	GE10600015
Amersham™ Protran® Western blotting membranes, nitrocellulose pore size 0.45 µm, sheet W x L 200 mm x 200 mm, pkg of 25 ea	GE10600041
Amersham™ Protran® Western blotting membranes, nitrocellulose pore size 0.45 µm, roll W x L 300 mm x 4 m, pkg of 1 ea	GE10600002
Amersham™ Protran® Western blotting membranes, nitrocellulose pore size 0.45 µm, roll W x L 200 mm x 4 m, pkg of 1 ea	GE10600007
Amersham™ Protran® Western blotting membranes, nitrocellulose pore size 0.45 µm, roll W x L 150 mm x 4 m, pkg of 1 ea	GE10600012
Amersham™ Protran® Premium Western blotting membranes, nitrocellulose pore size 0.45 µm, sheet W x L 80 mm x 90 mm, pkg of 25 ea	GE10600096
Amersham™ Protran® Premium Western blotting membranes, nitrocellulose pore size 0.45 µm, sheet W x L 200 mm x 200 mm, pkg of 10 ea	GE10600048
Amersham™ Protran® Premium Western blotting membranes, nitrocellulose pore size 0.45 µm, roll W x L 300 mm x 4 m, pkg of 1 ea	GE10600003
Amersham™ Protran® Premium Western blotting membranes, nitrocellulose pore size 0.45 µm, roll W x L 200 mm x 4 m, pkg of 1 ea	GE10600008
Amersham™ Protran® Supported Western blotting membranes, nitrocellulose pore size 0.45 µm, roll W x L 300 mm x 4 m, pkg of 1 ea	GE10600016
Amersham™ Protran® Supported Western blotting membranes, nitrocellulose pore size 0.45 µm, roll W x L 200 mm x 4 m, pkg of 1 ea	GE10600018
Protein Blot Stains	
Alcian Blue 8GX, certified by the Biological Stain Commission	A3157
Amido Black Staining Solution 2x, electrophoresis reagent	A8181
Glycoprotein Detection Kit	GLYCOPRO
Gold solution, colloidal, 0.0065% gold basis (AAS)	50755
Ponceau S, BioReagent, suitable for electrophoresis	P3504
Ponceau S solution, BioReagent, suitable for electrophoresis, 0.1 % (w/v) in 5% acetic acid	P7170
Reversible Protein Detection Kit for Membranes and Polyacrylamide Gels	RPROB

Blocking & Antibody Addition

Description	Cat. No.
Immobilon® GO immunodetection device	
Trial Pack: Immobilon® GO Device 2/pk	IMGDV002
Immobilon® GO Device 10/pk	IMGDV010
SNAP i.d.® 2.0 systems	
SNAP i.d.® 2.0 System - Mini (7.5 x 8.4 cm)	SNAP2MINI
SNAP i.d.® 2.0 System - Midi (8.5 x 13.5 cm)	SNAP2MIDI
SNAP i.d.® 2.0 System - MultiBlot (4.5 x 8.4 cm)	SNAP2MB3
SNAP i.d.® 2.0 System - Mini and Midi (7.5 x 8.4 cm and 8.5 x 13.5 cm)	SNAP2MM
SNAP i.d.® 2.0 System - Mini and MultiBlot (7.5 x 8.4 cm and 4.5 x 8.4 cm)	SNAP2MB1
SNAP i.d.® 2.0 System - Midi and MultiBlot (8.5 x 13.5 cm and 4.5 x 8.4 cm)	SNAP2MB2
SNAP i.d.® 2.0 consumables	
SNAP i.d.® 2.0 Mini Blot Holders (7.5 x 8.4 cm) 100/pk	SNAP2BHMN0100
SNAP i.d.® 2.0 Midi Blot Holders (8.5 x 13.5 cm) 100/pk	SNAP2BHMD0100
SNAP i.d.® 2.0 MultiBlot Holders (4.5 x 8.4 cm) 50/pk	SNAP2BHMB050
SNAP i.d.® 2.0 accessories	
SNAP i.d.® 2.0 Antibody Collection Tray 20/pk	SNAPABTR
SNAP i.d.® 2.0 Blot Roller 1/pk	SNAP2RL
SNAP i.d.® 2.0 Mini Blot Holding Frames (double pack) 2/pk	SNAP2FRMN02
SNAP i.d.® 2.0 Midi Blot Holding Frames (double pack) 2/pk	SNAP2FRMD02
SNAP i.d.® 2.0 Mini Blot Holding Frame (single pack) 1/pk	SNAP2FRMN01
SNAP i.d.® 2.0 Midi Blot Holding Frame (single pack) 1/pk	SNAP2FRMD01
SNAP i.d.® 2.0 MultiBlot Frame (single pack) 1/pk	SNAP2FRMB01
Signal enhancer reagents	
Immobilon® Signal Enhancer for Immunodetection	WBSH0500
Blocking buffers and reagents	
BløK®-CH Reagent Chemiluminescence detection 500 mL/bottle	WBAVDCH01
BløK®-FL Reagent Fluorescence detection 500 mL/bottle	WBAVDL01
BløK®-PO Reagent Phosphoprotein detection 500 mL/bottle	WBAVDP001
Western Blocker™ Solution, for HRP detection systems	W0138
Bovine Serum Albumin, heat shock fraction, protease free, fatty acid free, essentially globulin free, pH 7, ≥98%	A7030
Bovine Serum Albumin, heat shock fraction, pH 7, ≥98%	A7906
Gelatin blocking buffer, powder blend	G7663
Phosphate buffered saline, 10x concentrate, BioPerformance Certified, suitable for cell culture	P5493
Phosphate buffered saline, dry powder, pH 7.4, contains 3% nonfat milk	P2194
Phosphate buffered saline, pH 7.4, contains BSA, powder	P3688
Phosphate buffered saline, powder, pH 7.4	P3813
Tris Buffered Saline, with BSA, pH 8.0, powder	T6789
Wash buffers	
Phosphate buffered saline, BioUltra, pH 7.4 (in solution), contains TWEEN 20, tablets	08057
Phosphate buffered saline, pH 7.4, contains TWEEN® 20, dry powder	P3563
Tris Buffered Saline, pH 8.0, powder	T6664
Tris buffered saline with TWEEN® 20, BioUltra, tablets (for 500 mL), pH 7.6	91414
Tris Buffered Saline, with TWEEN® 20, pH 8.0, powder	T9039
Peroxidase conjugates	
Anti-Mouse IgG (Fab specific)-Peroxidase antibody produced in goat, affinity isolated antibody buffered aqueous solution	A3682
Anti-Mouse IgG (Fab specific)-Peroxidase antibody produced in goat, affinity isolated antibody buffered aqueous solution	A9917
Anti-Mouse IgG (Fc specific)-Peroxidase antibody produced in goat, affinity isolated antibody	A0168
Anti-Mouse IgG (Fc specific)-Peroxidase antibody produced in goat, affinity isolated antibody buffered aqueous solution	A2554
Anti-Mouse IgG (whole molecule)-Peroxidase antibody produced in rabbit, IgG fraction of antiserum buffered aqueous solution	A9044
Anti-Rabbit IgG (whole molecule)-Peroxidase antibody produced in goat, affinity isolated antibody	A0545
Avidin-Peroxidase, lyophilized powder	A3151
Anti-Goat IgG (whole molecule)-Peroxidase antibody produced in rabbit, affinity isolated antibody buffered aqueous solution	A5420
Anti-Mouse IgG (whole molecule)-Peroxidase antibody produced in goat, affinity isolated antibody buffered aqueous solution	A4416
Anti-Rabbit IgG (whole molecule)-Peroxidase antibody produced in goat, affinity isolated antibody buffered aqueous solution	A6154
Anti-Goat IgG Antibody, HRP conjugate; produced in rabbit; Purified rabbit IgG conjugated to horseradish peroxidase in buffer containing 0.01 M Sodium Phosphate, 0.15 M NaCl, pH 7.1 with 15 mg/mL BSA and 0.01% thimerosal. Lyophilized.	AP106P

Blocking & Antibody Addition

Description	Cat. No.
Peroxidase conjugates (continued)	
Anti-Rabbit IgG Antibody, HRP-conjugate; produced in rabbit; Goat polyclonal conjugated with horseradish peroxidase IgG in buffer containing 50% storage buffer 0.02 M potassium phosphate, 0.15 M NaCl, pH 7.2, 10 mg/mL BSA, 0.01% (w/v) gentamicin sulfate and 50% glycerol. Lyophilized.	12-348
Anti-Mouse IgG Antibody, HRP conjugate; produced in goat; Purified goat polyclonal IgG in buffer containing conjugated to horseradish peroxidase lyophilized from 0.02 M Potassium Phosphate, 0.15 M NaCl, pH 7.2, 10 mg/mL BSA, and 0.01% gentamicin sulfate.	12-349
Anti-Mouse IgG Antibody, (H+L) HRP conjugate; produced in goat; Purified Goat IgG in buffer containing liquid in PBS, pH 7.1 with 15 mg/mL BSA and 0.01% thimerosal.	AP308P
Anti-Rabbit light chain, HRP conjugate Antibody; produced in mouse; Purified from antiserum by chromatographic procedures. Lyophilized in 0.01M Sodium Phosphate, 0.25M NaCl, pH 7.6. 15 mg/mL BSA as stabilizer.	MAB201P
Anti-Mouse light chain Antibody, HRP conjugate; produced in goat; Lyophilized in 0.01M Sodium Phosphate, 0.25M NaCl, pH 7.6. 15 mg/mL BSA as stabilizer. Purified by immunoaffinity chromatography.	AP200P
Fluorophore conjugates	
ECL Plex™ G-A-M IgG, Cy [®] 3, GE Healthcare, PA43009, pack of 150 µg	GEPA43009
ECL Plex™ G-A-M IgG, Cy [®] 5, GE Healthcare, PA45010, pack of 600 µg	GEPA45010
ECL Plex™ Gar IgG Cy [®] 3, GE Healthcare, 28-9011-06, pack of 150 µg	GE28-9011-06
ECL Plex™ G-A-R IgG, Cy [®] 5, GE Healthcare, PA45011, pack of 150 µg	GEPA45011

Blot Stripping Reagents

Description	Cat. No.
ReBlot™ Plus Mild Antibody Stripping Solution, 10 x 50 mL	2502
ReBlot™ Plus Strong Antibody Stripping Solution, 10 x 50 mL	2504

Detection

Description	Cat. No.
Immobilon® Western Chemiluminescent HRP Substrate	
2 x 25 mL	WBKLS0050
2 x 50 mL	WBKLS0100
2 x 250 mL	WBKLS0500
Immobilon® ECL Ultra Western HRP Substrate	
2 x 50 mL	WBULS0100
2 x 250 mL	WBULS0500
Immobilon® Classico Western HRP Substrate	
100 mL	WBLUC0100
500 mL	WBLUC0500
Immobilon® Crescendo Western HRP Substrate	
100 mL	WBLUR0100
500 mL	WBLUR0500
Immobilon® Forte Western HRP Substrate	
100 mL	WBLUF0100
500 mL	WBLUF0500
Amersham ECL Western Blotting Detection Reagent	
Amersham™ ECL™ Western Blotting Reagent Pack GE Healthcare, RPN2124, pack of 1 kit	GERPN2124
ECL™ Western Blotting Analysis System GE Healthcare, RPN2108	GERPN2108
ECL™ Western Blotting Detection Reagents GE Healthcare, RPN2209	GERPN2209
ECL™ Western Blotting Reagents GE Healthcare, RPN2106	GERPN2106
ECL™ Western Blotting Reagents GE Healthcare, RPN2134	GERPN2134
ECL™ Western Blotting Reagents GE Healthcare, RPN2109	GERPN2109
ECL™ Start Western Blotting Reagent GE Healthcare, RPN2109	GERPN3243
ECL™ Start Western Blotting Reagent GE Healthcare, RPN2109	GERPN3244

Detection

Description	Cat. No.
Amersham ECL Prime Western Blotting Detection Reagent	
ECL™ Prime Western Blotting System GE Healthcare, RPN2232	GERPN2232
ECL™ Prime Western Blotting Detection Reagent GE Healthcare, RPN2236	GERPN2236
Amersham ECL Select Western Blotting Detection Reagent	
ECL™ Select Western Blotting Detection Reagent GE Healthcare, RPN2235	GERPN2235
Detection Films	
Hyperfilm™ ECL™, W x L 18 cm x 24 cm, pack of 100 ea	GE28-9068-37
Hyperfilm™ MP, W x L 8 in. x 10 in., pack of 50 ea	GE28-9068-45
Chromogenic Substrates	
3,3',5,5'-Tetramethylbenzidine (TMB) Liquid Substrate System for Membranes, ready-to-use solution	T0565
3,3'-Diaminobenzidine (DAB) Enhanced Liquid Substrate System tetrahydrochloride, for Membrane Applications	D6815

Accessories

Description	Cat. No.
Scienceware® gel incubating tray W x L 1.6 cm (0.63 in.) x 11.5 cm (4.5 in.)	BAF451000000-5EA
Scienceware® gel incubating tray W x L 9.0 cm (3.5 in.) x 9.0 cm (3.5 in.)	BAF451000001-5EA
Scienceware® gel incubating tray W x L 16.5 cm (6.5 in.) x 16.5 cm (6.5 in.)	BAF451000002-5EA
Scienceware® electrophoresis plate rack, polypropylene	E1028-1EA
Gel Handler™ gel support	Z376957-1PAK

To learn more, please visit: SigmaAldrich.com/westernblot

To place an order or receive technical assistance in Europe, please call Customer Service:

France: 0825 045 645

Spain: 901 516 645 Option 1

Germany: 069 86798021

Switzerland: 0848 645 645

Italy: 848 845 645

United Kingdom: 0870 900 4645

For other countries across Europe, please call: +44 (0) 115 943 0840

Or visit: MerckMillipore.com/offices

For Technical Service visit: MerckMillipore.com/techservice

© 2020 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. Merck and the vibrant are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.