

## User Guide

# Immobilon® ECL UltraPlus

Western HRP Substrate

**WBULP-10ML, WBULP-20ML, WBULP-100ML**

## Introduction

Immobilon® ECL UltraPlus detection reagent is a non-isotopic, luminol-based chemiluminescence substrate, designed for the chemiluminescent detection of immobilized proteins conjugated with horseradish peroxidase (HRP). It is compatible with both PVDF and nitrocellulose blotting membranes, as well as commonly used buffers and blocking reagents.

Immobilon® ECL UltraPlus substrate is for research use only. It is not for use in diagnostic procedures.

## Package Contents

Solution A: Luminol/enhancer solution (amber bottle)  
Solution B: Peroxide solution (white bottle)

Cat. No.	Solution Volumes	Membrane Area
WBULP-10ML	5 mL Solution A 5 mL Solution B	Sufficient for 200 cm <sup>2</sup>
WBULP-20ML	10 mL Solution A 10 mL Solution B	Sufficient for 400 cm <sup>2</sup>
WBULP-100ML	50 mL Solution A 50 mL Solution B	Sufficient for 2000 cm <sup>2</sup>

## Storage

Store at room temperature for up to one year from date of receipt.

## Usage Guidelines

- Immobilon® ECL UltraPlus substrate is a two-part highly stable chemiluminescent substrate.
- Immobilon® ECL UltraPlus substrate is extremely sensitive and the concentration of the primary and secondary antibody may need to be reduced for optimal signal-noise ratio.
- Suggested antibody dilutions to start is 1:10,000 - 1:200,000 for primary antibody concentration and 1:300,000 - 1:1,000,000 for secondary antibody.

- Use of blocking buffer to dilute antibodies may reduce background and increase sensitivity.
- To avoid high background, always wear gloves when handling membrane.
- Handle the membrane with blunt tip forceps (Cat. No. XX6200006P) to avoid tearing it.
- Do not use sodium azide in any blocking buffers or wash solutions, since it inhibits HRP activity.

## Chemiluminescent Detection Protocol

Approximately 0.1 mL of Immobilon® ECL UltraPlus substrate is required per cm<sup>2</sup> of membrane area. The volumes of substrate required for some common membrane sizes are indicated below:

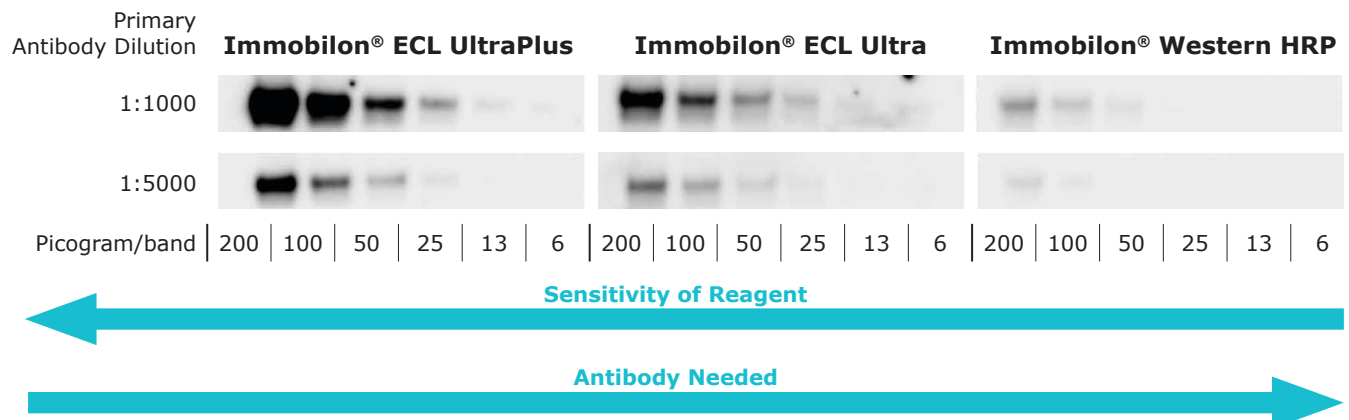
Blot Size	Immobilon® ECL UltraPlus Substrate Volume
7 × 8.4 cm	6 mL
10 × 10 cm	10 mL
8.5 × 13.5 cm	12 mL

Prepare working solution by gently mixing solutions A and B in a 1:1 ratio. Avoid vigorous agitation. Store protected from light until ready to use.

1. Place the blot, protein-side up, in a clean container or on a clear plastic sheet protector, and add Immobilon® ECL UltraPlus substrate onto the blot.
2. Incubate the blot for 1 to 5 minutes at room temperature.
3. Drain the excess substrate, transfer to a clean sheet protector, and cover the blot with plastic wrap or equivalent.
4. Visualize blots using a CCD-based imaging system or place blots into film cassette and expose to suitable x-ray film for required amount of time.

## Performance

Detection of Recombinant Human IKB alpha Protein (38 kDa)



Two-fold dilution series of recombinant human IKB alpha (Abcam ab113133) was resolved by SDS-PAGE and transferred to Immobilon®-P membrane using a fast semi-dry blotting system. The blots were blocked with 0.5% non-fat dry milk, probed with primary antibody rabbit anti-IKB alpha (Abcam ab32518) diluted 1:1,000 and 1:5,000 as well as secondary antibody goat anti-rabbit IgG HRP conjugated (Cat. No. AP187P) diluted 1:300,000. The blot was then incubated with Immobilon® ECL UltraPlus substrate for 2 minutes and exposed to a digital imager for 60 seconds.

## Product Ordering

Description	Membrane Area	Cat. No.
Immobilon® ECL UltraPlus Western HRP Substrate		
5 mL Solution A		
5 mL Solution B	200 cm <sup>2</sup>	WBULP-10ML
10 mL Solution A		
10 mL Solution B	400 cm <sup>2</sup>	WBULP-20ML
50 mL Solution A		
50 mL Solution B	2000 cm <sup>2</sup>	WBULP-100ML
Blotting Membranes	Qty/Pk	
Immobilon®-P PVDF, 0.45 µm, 26.5 x 375 cm roll	1	IPVH00010
Immobilon®-P PVDF, 0.45 µm, 7 x 8.4 cm sheet	50	IPVH07850
Immobilon®-P PVDF, 0.45 µm, 8.5 x 13.5 cm sheet	10	IPVH08130
Immobilon®-P <sup>SQ</sup> PVDF, 0.2 µm, 26.5 x 375 cm roll	1	ISEQ00010
Immobilon®-P <sup>SQ</sup> PVDF, 0.2 µm, 7 x 8.4 cm sheet	50	ISEQ07850
Accessories	Qty/Pk	
Filter forceps, blunt end	3	XX6200006P

For primary and secondary antibodies go to [SigmaAldrich.com/antibodies](https://www.sigmaaldrich.com/antibodies).

## Disposal

Collect and dispose of used material according to all applicable international, federal, state, and local regulations.

## Safety Data Sheet

Safety Data Sheets (SDS) are available on the product page at [SigmaAldrich.com](https://www.sigmaaldrich.com).

## Technical Assistance

Visit the tech service page on our web site at [SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice).

## Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://www.sigmaaldrich.com/terms).

## Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://www.sigmaaldrich.com/offices).

## Troubleshooting

Symptom	Possible Cause	Solution
High background	Concentration of HRP-conjugated antibody too high	Increase dilution of secondary antibody.
	Inefficient blocking	Optimize blocking conditions.
	Insufficient washing	Increase wash buffer volume and/or increase number of washes.
Negative staining (white bands on black background)	Substrate depleted due to high antibody concentration	Increase antibody dilution and/or decrease antigen concentration.
Highly speckled background	High concentration of antibody and/or protein	Increase dilution of primary and secondary antibodies.
	Formation of aggregates in the HRP conjugate	Reduce protein load in the gel. Filter conjugate through 0.2 µm filter.
Nonspecific bands	High concentration of primary antibody	Increase dilution of primary antibody.
Signal disappears quickly in a blot that initially had a very high signal	High HRP-antibody concentration exhausted the substrate prematurely	Increase dilution of antibody significantly and/or decrease antigen concentration.
Weak or no signal	Antibody concentration too low	Increase exposure time. Decrease antibody dilution and/or increase antigen concentration.
	Protein did not transfer to membrane	Verify protein transfer.
	Antibody did not bind to target	Verify that antibody is specific for intended target.
	Expired substrate	Replace substrate with new lot.

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

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

Merck, Millipore, Immobilon, and Sigma-Aldrich 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.

© 2021-2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

Document Template 00035533 Ver 1.0

20595716 Ver 3.0, Rev 05JUL2024, AU

