

# Assay Kits for Metabolite and Enzyme Analysis

Quality is our specialty,  
let us help you with **your** specialty



The Life Science  
business of Merck  
operates as  
MilliporeSigma in  
the U.S. and Canada.

**Sigma-Aldrich®**  
Lab & Production Materials

# The right assay kit for your specific needs

We provide a comprehensive portfolio of assay kits for a broad range of applications in life science research, drug discovery and development, and environmental monitoring and testing.

- easy-to-use, precise, and highly reliable assays
- spectrophotometric and/or fluorometric detection methods
- suitable for high-throughput analysis
- pre-tested for optimum quality to best serve your application needs

**Pulmonology**   
Page 12



**Rheumatology**  
Page 11



**Dermatology**  
Page 11



**Hematology**  
Page 10



**Obesity**  
Page 6



**Genitourinary**  
Page 10

**Orthopedics**  
Page 11





## Cardiology Cardiovascular

Page 6

## Neurology

Page 7

## Endocrinology- Diabetes

Page 9

## Gastroenterology

Page 7

### Additional Areas of Specialization

	Oncology	4
	Sample Preparation	4
	Geriatrics	11
	Pediatrics-Neonatology	11
	Nutrition	12
	Environmental Testing	12
	Food-Beverage-Cosmetic Testing	13
	Drug Discovery/Pharma	14



## Sample Preparation

Kit #	Product
MAK227	Chloramphenicol Extraction Kit
MAK175	Cholesterol Extraction Kit
MAK226	Enrofloxacin Extraction Kit
MAK174	Fatty Acid Extraction Kit

Kit #	Product
MAK338	Fatty Acid Extraction Kit, High Standard (450 ug/ml)
MAK225	Fatty Acid Methyl Ester Purification Kit Manufactured by Nacalai Tesque, Inc.
MAK224	Fatty Acid Methylation Kit Manufactured by Nacalai Tesque, Inc.



## Oncology

Kit #	Product
MAK474	Acetate Assay Kit
MAK566	Acetyl-Coenzyme A Assay Kit
MAK446	Acid Phosphatase Assay Kit
MAK525	Acid Phosphatase Fluorometric Assay Kit
MAK337	Aconitase Assay Kit
MAK433	Adenosine Quantification Assay Kit
MAK313	Adipolysis Assay Kit
MAK133	ADP Assay Kit
MAK518	ADP Assay Kit
MAK135	ADP/ATP Ratio Assay Kit
MAK436	AF Cholesterol Assay Kit
MAK500	Alanine Assay Kit
MAK327	Aldehyde Dehydrogenase Inhibitor Screening Kit
MAK447	Alkaline Phosphatase Assay Kit
MAK334	Antioxidant Assay Kit
MAK112	Arginase Activity Assay Kit
MAK473	ATP Assay Kit
MAK113	ATPase/GTPase Assay Kit
MAK138	Autophagy Assay Kit
MAK540	Beta-Hydroxybutyrate (Ketone Body) Assay Kit
MAK309	Bile Acid Assay Kit
MAK477	Calcium Assay Kit
MAK486	Carbonyl Assay Kit
MAK457	Caspase-3 Assay Kit
MAK325	Cholesterol Uptake Assay Kit
MAK333	Citrate Assay Kit
CS0720	Citrate Synthase Assay Kit
MAK504	Coenzyme A Assay Kit
MAK322	Collagen Assay Kit
MAK162	Colorimetric Maleimide Assay Kit
MAK152	Colorimetric Sphingomyelinase Assay Kit
MAK536	Cytotoxicity Assay Kit

Kit #	Product
MAK320	D-2-Hydroxyglutarate (D2HG) Assay Kit
MAK526	Diamine Oxidase Assay Kit
MAK174	Fatty Acid Extraction Kit
MAK338	Fatty Acid Extraction Kit (High Standard)
MAK156	Fatty Acid Uptake Kit
MAK509	Ferric Reducing Antioxidant Power (FRAP) Assay Kit
MAK552	Fluo-4 No Wash Calcium Assay Kit
MAK553	Fluorimetric cADP-Ribose Assay Kit
MAK150	Fluorometric Intracellular pH Assay Kit
MAK142	Fluorometric Intracellular Ros Kit (Deep Red)
MAK143	Fluorometric Intracellular Ros Kit (Green)
MAK144	Fluorometric Intracellular Ros Kit (Orange)
MAK145	Fluorometric Intracellular Ros Kit (Red)
MAK154	Fluorometric Sphingomyelin Assay Kit
MAK151	Fluorometric Thiol Assay Kit
MAK131	Formaldehyde Assay
MAK466	Free Fatty Acid Assay Kit
MAK492	Fumarate Assay Kit
MAK489	Glucose Uptake Assay Kit
MAK542	Glucose Uptake Assay Kit
MAK548	Glucose-6-Phosphate (G6P) Assay Kit
MAK451	Glucose-6-Phosphate Dehydrogenase Assay Kit
MAK452	Glucose-6-Phosphate Dehydrogenase Inhibitor Screening Kit
MAK330	Glutamate Assay Kit
MAK499	Glutamate Dehydrogenase (GDH) Assay Kit
MAK517	Glutathione (GSH) Assay Kit
MAK440	Glutathione GSH/GSSG Assay Kit
MAK437	Glutathione Peroxidase Assay Kit
MAK535	Glutathione Reductase Kit
MAK453	Glutathione S-transferase Assay Kit
MAK117	Glycerol Assay Kit
MAK439	Glycolysis Assay Kit

Kit #	Product
MAK114	Glyoxalase I Activity Assay Kit
MAK543	High Sensitivity Glucose Quantitation Kit
MAK458	Hyaluronidase Inhibitor Screening Assay Kit
MAK569	Hydroxyproline Assay Kit
MAK463	Hydroxyproline Assay Kit (Perchlorate-Free)
MAK556	Hypochlorite (Hypochlorous Acid) Assay Kit
MAK472	Iron Assay Kit
MAK524	Isocitrate Assay Kit
MAK493	Isocitrate Dehydrogenase (IDH) Assay Kit
MAK134	Ketone Body Assay
MAK570	Lactate Assay Kit
MAK464	Lactate Dehydrogenase Assay Kit
MAK487	Lactose Assay Kit
MAK490	L-Amino Acid Assay Kit
MAK529	LDH Cytotoxicity Assay Kit
MAK482	Lipase Assay Kit
MAK568	Lipid Peroxidation (MDA) Assay Kit
MAK329	L-Lactate Assay Kit (Colorimetric)
MAK443	L-Lactate Assay Kit (Fluorometric)
MAK511	Malate Assay Kit
MAK160	Mitochondrial Membrane Potential Kit (Flow Cytometry)
MAK159	Mitochondrial Membrane Potential Kit (Microplate)
MAK563	MPO Colorimetric Activity Assay Kit

Kit #	Product
MAK161	Multidrug Resistance Assay Kit
MAK516	Myeloperoxidase (MPO) Assay Kit
MAK468	NAD/NADH Assay Kit
MAK460	NAD <sup>+</sup> /NADH Assay Kit
MAK479	NADP/NADPH Assay Kit
MAK532	Nitric Oxide Synthase (NOS) Assay Kit
MAK515	Oxaloacetate Assay Kit
MAK506	Peroxidase Assay Kit
MAK137	Phospholipase D Activity Assay Kit
MAK461	pNPP Phosphatase Assay Kit
MAK172	Proteasome 20s Activity Assay Kit
MAK332	Pyruvate Assay Kit
MAK567	Pyruvate Dehydrogenase Activity Assay Kit
MAK485	Pyruvate Kinase Assay Kit
MAK314	Sialic Acid Assay Kit
MAK528	Superoxide Dismutase (SOD) Assay Kit
CS0170	Thioredoxin Reductase Assay Kit
MAK550	Tyrosinase Assay Kit
MAK173	Universal Fluorometric Kinase Assay Kit
MAK496	Zinc Assay Kit
MAK541	$\alpha$ -Ketoglutarate Assay Kit
MAK444	$\alpha$ -L-Fucosidase Assay Kit



## Cardiology - Cardiovascular

Kit #	Product
MAK435	Acetylcholine Assay Kit
MAK337	Aconitase Assay Kit
MAK433	Adenosine Quantification Assay Kit
MAK313	Adipolysis Assay Kit
MAK436	AF Cholesterol Assay Kit
MAK331	AF HDL and LDL/VLDL Assay Kit
MAK500	Alanine Assay Kit
CS0002	Angiotensin I Converting Enzyme (ACE) Activity Assay Kit (Fluorometric)
MAK334	Antioxidant Assay Kit
MAK467	Aspartate Transaminase (AST) Assay Kit
MAK309	Bile Acid Assay Kit
MAK551	Butyrylcholinesterase (BChE) Activity Assay Kit
MAK486	Carbonyl Assay Kit
MAK531	Catalase Assay Kit
MAK106	CETP Activity Assay Kit
MAK305	CETP RP Activity Assay Kit
CS0005	Cholesterol Quantification Assay kit
MAK325	Cholesterol Uptake Assay Kit
MAK116	Creatine Kinase Assay Kit
MAK509	Ferric Reducing Antioxidant Power (FRAP) Assay Kit
MAK553	Fluorimetric cADP-Ribose Assay Kit
MAK165	Fluorimetric Hydrogen Peroxide Assay Kit
MAK166	Fluorimetric Hydrogen Peroxide Assay Kit
MAK535	Glutathione Reductase Kit
MAK546	Glycerol 3-Phosphate (G3P) Assay Kit
MAK565	High Sensitivity Triglyceride Fluorometric Assay Kit
MAK569	Hydroxyproline Assay Kit
MAK463	Hydroxyproline Assay Kit (Perchlorate-Free)
MAK164	Intracellular Hydrogen Peroxide Assay

Kit #	Product
MAK306	LCFC-LCAT Acyltransferase Activity Assay
MAK560	Lipid (Oil Red O) Staining Kit
MAK568	Lipid Peroxidation (MDA) Assay Kit
MAK329	L-Lactate Assay Kit (Colorimetric)
MAK443	L-Lactate Assay Kit (Fluorometric)
MAK109	Lpl Activity Assay Kit
MAK507	Magnesium Assay Kit
MAK307	Malachite Green Phosphate Assay Kit
MAK563	Mpo Colorimetric Activity Assay Kit
MAK110	MTP Activity Assay Kit
MAK516	Myeloperoxidase (MPO) Assay Kit
MAK454	Nitric Oxide Assay Kit
MAK532	Nitric Oxide Synthase (NOS) Assay Kit
MAK323	Nitric Oxide Synthase Inhibitor Screening Kit
MAK506	Peroxidase Assay Kit
MAK311	Peroxide Assay Kit
MAK308	Phosphate Assay Kit
MAK108	Pltp Activity Assay Kit
MAK461	pNPP Phosphatase Assay Kit
MAK445	Protein Creatinine Ratio Assay Kit
MAK157	Renin Assay Kit
MAK314	Sialic Acid Assay Kit
MAK335	Succinate Assay Kit
MAK528	Superoxide Dismutase (SOD) Assay Kit
CS0170	Thioredoxin Reductase Assay Kit
MAK564	Triglyceride Quantification Kit
MAK173	Universal Fluorometric Kinase Assay Kit
MAK483	Uric Acid Assay Kit
MAK497	Xanthine Oxidase Assay Kit



## Obesity

Kit #	Product
MAK313	Adipolysis Assay Kit
MAK500	Alanine Assay Kit
MAK551	Butyrylcholinesterase (BChE) Activity Assay Kit


Kit #	Product
MAK325	Cholesterol Uptake Assay Kit
MAK546	Glycerol 3-Phosphate (G3P) Assay Kit
MAK560	Lipid (Oil Red O) Staining Kit



## Gastroenterology

Kit #	Product
MAK321	Acetaldehyde Assay Kit
MAK446	Acid Phosphatase Assay Kit
MAK525	Acid Phosphatase Fluorometric Assay Kit
MAK498	Alcohol Dehydrogenase (ADH) Assay Kit
MAK447	Alkaline Phosphatase Assay Kit
MAK571	ALT Activity Assay Kit
MAK310	Ammonia Assay Kit
MAK538	Ammonia Assay Kit
MAK112	Arginase Assay Kit
MAK467	Aspartate Transaminase (AST) Assay Kit
MAK124	BCG Albumin Assay Kit
MAK125	BCP Albumin Assay Kit
MAK540	Beta-Hydroxybutyrate (Ketone Body) Assay Kit
MAK309	Bile Acid Assay Kit
MAK126	Bilirubin Assay Kit
MAK562	Branched Chain Amino Acid Kit
MAK486	Carbonyl Assay Kit
MAK322	Collagen Assay Kit
MAK127	Copper Assay Kit
MAK526	Diamine Oxidase Assay Kit

Kit #	Product
MAK336	D-Lactate Assay Kit
MAK156	Fatty Acid Uptake Kit
MAK499	Glutamate Dehydrogenase (GDH) Assay Kit
MAK546	Glycerol 3-Phosphate (G3P) Assay Kit
MAK547	High Sensitivity Beta-Hydroxybutyrate (Ketone Body) Assay Kit
MAK128	Indican Assay Kit
MAK326	Indole Assay Kit
MAK482	Lipase Assay Kit
MAK109	LPL Activity Assay Kit
MAK307	Malachite Green Phosphate Assay Kit
MAK308	Phosphate Assay Kit
MAK461	pNPP Phosphatase Assay Kit
MAK332	Pyruvate Assay Kit
MAK317	Sorbitol Dehydrogenase Assay Kit
MAK558	Trypsin Activity Assay Kit
MAK120	Urease Assay Kit
MAK497	Xanthine Oxidase Assay Kit
MAK478	$\alpha$ -Amylase Activity Assay Kit
MAK444	$\alpha$ -L-Fucosidase Assay Kit
MAK129	$\beta$ -Glucosidase Assay Kit



## Neurology

Kit #	Product
MAK435	Acetylcholine Assay Kit
MAK119	Acetylcholinesterase Assay Kit
MAK324	Acetylcholinesterase Inhibitor Screening Kit
MAK566	Acetyl-Coenzyme A Assay Kit
MAK337	Aconitase Assay Kit
MAK433	Adenosine Quantification Assay Kit
MAK436	AF Cholesterol Assay Kit
CS0002	Angiotensin I Converting Enzyme (ACE) Activity Assay Kit (Fluorometric)
MAK334	Antioxidant Assay Kit
MAK533	Arginase Assay Kit
MAK495	Aspartate Assay Kit
MAK473	ATP Assay Kit
MAK171	Biotin Assay Kit
MAK562	Branched Chain Amino Acid Kit
MAK477	Calcium Assay Kit
MAK486	Carbonyl Assay Kit
MAK457	Caspase-3 Assay Kit
MAK531	Catalase Assay Kit

Kit #	Product
MAK508	Choline Assay Kit
MAK139	Colorimetric Aldehyde Assay Kit
MAK140	Colorimetric Aldehyde Assay Kit, Blue
MAK127	Copper Assay Kit
MAK494	Creatine Assay Kit
MAK475	Creatinine Assay Kit
MAK320	D-2-Hydroxyglutarate (D2HG) Assay Kit
MAK336	D-Lactate Assay Kit
MAK509	Ferric Reducing Antioxidant Power (FRAP) Assay Kit
MAK552	Fluo-4 No Wash Calcium Assay Kit
MAK554	Fluorimetric Acetylcholinesterase Assay Kit
MAK553	Fluorimetric cADP-Ribose Assay Kit
MAK165	Fluorimetric Hydrogen Peroxide Assay Kit
MAK166	Fluorimetric Hydrogen Peroxide Assay Kit
MAK150	Fluorometric Intracellular pH Assay Kit
MAK142	Fluorometric Intracellular Ros Kit (Deep Red)
MAK143	Fluorometric Intracellular Ros Kit (Green)
MAK144	Fluorometric Intracellular Ros Kit (Orange)
MAK145	Fluorometric Intracellular Ros Kit (Red)



## Neurology (continued)

Kit #	Product
MAK521	Fumarase Assay Kit
MAK330	Glutamate Assay Kit
MAK440	Glutathione GSH/GSSG Assay Kit
MAK535	Glutathione Reductase Kit
MAK453	Glutathione S-transferase Assay Kit
MAK546	Glycerol 3-Phosphate (G3P) Assay Kit
MAK565	High Sensitivity Triglyceride Fluorometric Assay Kit
MAK569	Hydroxyproline Assay Kit
MAK463	Hydroxyproline Assay Kit (Perchlorate-Free)
MAK164	Intracellular Hydrogen Peroxide Assay
MAK568	Lipid Peroxidation (MDA) Assay Kit
MAK512	Malate Dehydrogenase Assay Kit
MAK557	Melanin Assay Kit
MAK520	Monoamine Oxidase (MAO) Inhibitor Screening Kit
MAK136	Monoamine Oxidase Assay Kit
MAK460	NAD <sup>+</sup> /NADH Assay Kit

Kit #	Product
MAK454	Nitric Oxide Assay Kit
MAK506	Peroxidase Assay Kit
MAK311	Peroxide Assay Kit
MAK484	Phenylalanine Assay Kit
MAK461	pNPP Phosphatase Assay Kit
MAK172	Proteasome 20S Activity Assay Kit
MAK567	Pyruvate Dehydrogenase Activity Assay Kit
MAK561	Succinate Dehydrogenase Assay Kit
MAK528	Superoxide Dismutase (SOD) Assay Kit
MAK564	Triglyceride Quantification Kit
MAK523	Tryptophan Assay Kit
MAK173	Universal Fluorometric Kinase Assay Kit
MAK496	Zinc Assay Kit
MAK123	$\alpha$ -Glucosidase Activity Assay Kit
MAK318	$\alpha$ -Mannosidase Assay Kit
MAK129	$\beta$ -Glucosidase Assay Kit







## Endocrinology - Diabetes

Kit #	Product
MAK337	Aconitase Assay Kit
MAK313	Adipolysis Assay Kit
CS0002	Angiotensin I Converting Enzyme (ACE) Activity Assay Kit (Fluorometric)
MAK334	Antioxidant Assay Kit
MAK495	Aspartate Assay Kit
MAK540	Beta-Hydroxybutyrate (Ketone Body) Assay Kit
MAK562	Branched Chain Amino Acid Kit
MAK551	Butyrylcholinesterase (BChE) Activity Assay Kit
MAK477	Calcium Assay Kit
MAK531	Catalase Assay Kit
CS0720	Citrate Synthase Assay Kit
MAK504	Coenzyme A Assay Kit
MAK156	Fatty Acid Uptake Kit
MAK552	Fluo-4 No Wash Calcium Assay Kit
MAK167	Fluorometric Maleimide Assay Kit
MAK151	Fluorometric Thiol Assay Kit
MAK466	Free Fatty Acid Assay Kit
MAK476	Glucose Assay Kit
MAK450	Glucose Dehydrogenase Assay Kit
MAK501	Glucose Oxidase Activity Assay Kit
MAK542	Glucose Uptake Assay Kit (Colorimetric)
MAK489	Glucose Uptake Assay Kit (Fluorometric)
MAK437	Glutathione Peroxidase Assay Kit
MAK453	Glutathione S-transferase Assay Kit

Kit #	Product
MAK546	Glycerol 3-Phosphate (G3P) Assay Kit
MAK465	Glycogen Assay Kit
MAK439	Glycolysis Assay Kit
MAK547	High Sensitivity Beta-Hydroxybutyrate (Ketone Body) Assay Kit
MAK543	High Sensitivity Glucose Quantitation Kit
MAK134	Ketone Body Assay Kit
MAK570	Lactate Assay Kit
MAK560	Lipid (Oil Red O) Staining Kit
MAK568	Lipid Peroxidation (MDA) Assay Kit
MAK329	L-Lactate Assay Kit (Colorimetric)
MAK443	L-Lactate Assay Kit (Fluorometric)
MAK507	Magnesium Assay Kit
MAK513	Maltose Assay Kit
MAK172	Proteasome 20S Activity Assay Kit
MAK445	Protein Creatinine Ratio Assay Kit
MAK442	Sorbitol Assay Kit
MAK317	Sorbitol Dehydrogenase Assay Kit
MAK173	Universal Fluorometric Kinase Assay Kit
MAK471	Urea Assay Kit III
MAK483	Uric Acid Assay Kit
MAK478	$\alpha$ -Amylase Activity Assay Kit
MAK123	$\alpha$ -Glucosidase Assay Kit
MAK444	$\alpha$ -L-Fucosidase Assay Kit
MAK456	$\beta$ -N-Acetylglucosaminidase Assay Kit



## Immunology

Kit #	Product
MAK112	Arginase Assay Kit
MAK138	Autophagy Assay Kit
MAK545	Beta-Lactamase Activity Assay Kit
MAK127	Copper Assay Kit
MAK536	Cytotoxicity Assay Kit
MAK526	Diamine Oxidase Assay Kit

Kit #	Product
MAK438	Glutamine Assay Kit
MAK557	Melanin Assay Kit
MAK454	Nitric Oxide Assay Kit
MAK532	Nitric Oxide Synthase (NOS) Assay Kit
MAK335	Succinate Assay Kit
MAK496	Zinc Assay Kit



## Hematology

Kit #	Product
MAK133	ADP Assay Kit
MAK518	ADP Assay Kit
MAK126	Bilirubin Assay Kit
MAK477	Calcium Assay Kit
MAK127	Copper Assay Kit
MAK536	Cytotoxicity Assay Kit
MAK554	Fluorimetric Acetylcholinesterase Assay Kit
MAK450	Glucose Dehydrogenase Assay Kit
MAK548	Glucose-6-Phosphate (G6P) Assay Kit
MAK503	Glucose-6-Phosphate Assay Kit
MAK451	Glucose-6-Phosphate Dehydrogenase Assay Kit
MAK452	Glucose-6-Phosphate Dehydrogenase Inhibitor Screening Kit
MAK535	Glutathione Reductase Kit

Kit #	Product
MAK316	Heme Assay Kit
MAK115	Hemoglobin Assay Kit
MAK472	Iron Assay Kit
MAK464	Lactate Dehydrogenase Assay Kit
MAK529	LDH Cytotoxicity Assay Kit
MAK563	MPO Colorimetric Activity Assay Kit
MAK516	Myeloperoxidase (MPO) Assay Kit
MAK461	pNPP Phosphatase Assay Kit
MAK485	Pyruvate Kinase Assay Kit
MAK314	Sialic Acid Assay Kit
MAK455	Whole Blood Hemoglobin (Hb) Kit
MAK496	Zinc Assay Kit
MAK129	$\beta$ -Glucosidase Assay Kit



## Genitourinary

Kit #	Product
MAK525	Acid Phosphatase Fluorometric Assay Kit
MAK530	Alkaline Phosphatase (ALP) Assay Kit
MAK310	Ammonia Assay Kit
MAK124	BCG Albumin Assay Kit
MAK125	BCP Albumin Assay Kit
MAK477	Calcium Assay Kit
MAK494	Creatine Assay Kit
MAK116	Creatine Kinase Assay Kit
MAK475	Creatinine Assay Kit
MAK537	D-Amino Acid Assay Kit
MAK552	Fluo-4 No Wash Calcium Assay Kit

Kit #	Product
MAK553	Fluorimetric cADP-Ribose Assay Kit
MAK150	Fluorometric Intracellular Ph Assay Kit
MAK169	High Sensitivity Pyrophosphate Assay Kit
MAK556	Hypochlorite (Hypochlorous Acid) Assay Kit
MAK490	L-Amino Acid Assay Kit
MAK315	Oxalate Assay Kit
MAK132	Sulfate Assay Kit
MAK471	Urea Assay Kit III
MAK483	Uric Acid Assay Kit
MAK496	Zinc Assay Kit
MAK456	$\beta$ -N-Acetylglucosaminidase Assay Kit



## Rheumatology

Kit #	Product
MAK337	Aconitase Assay Kit
MAK334	Antioxidant Assay Kit
MAK322	Collagen Assay Kit
MAK152	Colorimetric Sphingomyelinase Assay Kit
MAK169	High Sensitivity Pyrophosphate Assay Kit

Kit #	Product
MAK569	Hydroxyproline Assay Kit
MAK463	Hydroxyproline Assay Kit (Perchlorate-Free)
MAK534	Salicylate Assay Kit
MAK483	Uric Acid Assay Kit
MAK497	Xanthine Oxidase Assay Kit



## Orthopedics

Kit #	Product
MAK530	Alkaline Phosphatase (ALP) Assay Kit
MAK127	Copper Assay Kit
MAK116	Creatine Kinase Assay Kit
MAK166	Fluorimetric Hydrogen Peroxide Assay Kit

Kit #	Product
MAK165	Fluorimetric Hydrogen Peroxide Assay Kit
MAK164	Intracellular Hydrogen Peroxide Assay
MAK555	Lysyl Oxidase (LOX) Assay Kit
MAK311	Peroxide Assay Kit



## Geriatrics

Kit #	Product
MAK127	Copper Assay Kit
MAK116	Creatine Kinase Assay Kit
MAK165	Fluorimetric Hydrogen Peroxide Assay Kit
MAK166	Fluorimetric Hydrogen Peroxide Assay Kit
MAK440	Glutathione GSH/GSSG Assay Kit

Kit #	Product
MAK164	Intracellular Hydrogen Peroxide Assay
MAK460	NAD+/NADH Assay Kit
MAK532	Nitric Oxide Synthase (NOS) Assay Kit
MAK311	Peroxide Assay Kit



## Ped-Neonatology

Kit #	Product
MAK502	Galactose Assay Kit
MAK465	Glycogen Assay Kit
MAK134	Ketone Body Assay Kit

Kit #	Product
MAK487	Lactose Assay Kit
MAK311	Peroxide Assay Kit
MAK484	Phenylalanine Assay Kit



## Dermatology

Kit #	Product
MAK171	Biotin Assay Kit
MAK531	Catalase Assay Kit
MAK322	Collagen Assay Kit
MAK127	Copper Assay Kit

Kit #	Product
MAK569	Hydroxyproline Assay Kit
MAK555	Lysyl Oxidase (LOX) Assay Kit
MAK557	Melanin Assay Kit



## Pulmonology

Kit #	Product
MAK510	Chloride Assay Kit
MAK322	Collagen Assay Kit
MAK165	Fluorimetric Hydrogen Peroxide Assay Kit
MAK166	Fluorimetric Hydrogen Peroxide Assay Kit
MAK453	Glutathione S-transferase Assay Kit

Kit #	Product
MAK164	Intracellular Hydrogen Peroxide Assay
MAK482	Lipase Assay Kit
MAK555	Lysyl Oxidase (LOX) Assay Kit
MAK311	Peroxide Assay Kit
MAK528	Superoxide Dismutase (SOD) Assay Kit



## Nutrition

Kit #	Product
MAK313	Adipolysis Assay Kit
MAK505	Ascorbic Acid Assay Kit
MAK138	Autophagy Assay Kit
MAK124	BCG Albumin Assay Kit
MAK125	BCP Albumin Assay Kit
MAK171	Biotin Assay Kit
MAK504	Coenzyme A Assay Kit
MAK494	Creatine Assay Kit
MAK475	Creatinine Assay Kit
MAK537	D-Amino Acid Assay Kit

Kit #	Product
MAK526	Diamine Oxidase Assay Kit
MAK544	Enterokinase (Enteropeptidase) Activity Assay Kit
MAK509	Ferric Reducing Antioxidant Power (FRAP) Assay Kit
MAK128	Indican Assay Kit
MAK490	L-Amino Acid Assay Kit
MAK513	Maltose Assay Kit
MAK522	Starch Assay Kit
MAK559	Total Carbohydrate Assay Kit
MAK471	Urea Assay Kit III



## Environmental Testing

Kit #	Product
MAK434	Acetaldehyde Assay Kit
MAK334	Antioxidant Assay Kit
MAK448	Boron Assay Kit
MAK139	Colorimetric Aldehyde Assay Kit
MAK140	Colorimetric Aldehyde Assay Kit, Blue
MAK459	Detergent Assay Kit
MAK491	Formate Assay Kit

Kit #	Product
MAK117	Glycerol Assay Kit
MAK556	Hypochlorite (Hypochlorous Acid) Assay Kit
MAK118	Invertase Activity Assay Kit
MAK488	Phosphate Assay Kit
MAK442	Sorbitol Assay Kit
MAK120	Urease Assay Kit



## Food-Beverage - Cosmetic Testing

Kit #	Product
MAK434	Acetaldehyde Assay Kit (Colorimetric)
MAK321	Acetaldehyde Assay Kit (Fluorometric)
MAK474	Acetate Assay Kit
MAK331	AF HDL and LDL/VLDL Assay Kit
MAK500	Alanine Assay Kit
MAK334	Antioxidant Assay Kit
MAK505	Ascorbic Acid Assay Kit
MAK477	Calcium Assay Kit
MAK227	Chloramphenicol Extraction Kit
MAK510	Chloride Assay Kit
MAK508	Choline Assay Kit
MAK333	Citrate Assay Kit
MAK322	Collagen Assay Kit
MAK494	Creatine Assay Kit
MAK537	D-Amino Acid Assay Kit
MAK336	D-Lactate Assay Kit
MAK226	Enrofloxacin Extraction Kit
MAK480	Ethanol Assay Kit
MAK481	Ethanol Assay Kit
MAK338	Fatty Acid Extraction Kit (High Standard)
MAK174	Fatty Acid Extraction Kit, Low Standard (150 ug/ml)
MAK225	Fatty Acid Methyl Ester Purification Kit
MAK224	Fatty Acid Methylation Kit
MAK509	Ferric Reducing Antioxidant Power (FRAP) Assay Kit
MAK142	Fluorometric Intracellular Ros Kit (Deep Red)
MAK144	Fluorometric Intracellular Ros Kit (Orange)
MAK491	Formate Assay Kit
MAK449	Free Amino Nitrogen Assay Kit
MAK466	Free Fatty Acid Assay Kit
MAK502	Galactose Assay Kit
MAK476	Glucose Assay Kit

Kit #	Product
MAK501	Glucose Oxidase Activity Assay Kit
MAK330	Glutamate Assay Kit
MAK117	Glycerol Assay Kit
MAK547	High Sensitivity Beta-Hydroxybutyrate (Ketone Body) Assay Kit
MAK169	High Sensitivity Pyrophosphate Assay Kit
MAK565	High Sensitivity Triglyceride Fluorometric Assay Kit
MAK432	Histamine Quantification Assay Kit
MAK326	Indole Assay Kit
MAK118	Invertase Activity Assay Kit
MAK472	Iron Assay Kit
MAK524	Isocitrate Assay Kit
MAK570	Lactate Assay Kit
MAK487	Lactose Assay Kit
MAK490	L-Amino Acid Assay Kit
MAK507	Magnesium Assay Kit
MAK511	Malate Assay Kit
MAK513	Maltose Assay Kit
MAK514	Mannitol Assay Kit
MAK454	Nitric Oxide Assay Kit
MAK488	Phosphate Assay Kit
MAK122	Phospholipid Assay Kit
MAK168	Pyrophosphate Assay Kit
MAK442	Sorbitol Assay Kit
MAK522	Starch Assay Kit
MAK175	Sterol Extraction Kit
MAK528	Superoxide Dismutase (SOD) Assay Kit
MAK559	Total Carbohydrate Assay Kit
MAK550	Tyrosinase Assay Kit
MAK471	Urea Assay Kit III
MAK496	Zinc Assay Kit

# Drug Discovery/Pharma

Kit #	Product
MAK435	Acetylcholine Assay Kit
MAK313	Adipolysis Assay Kit
MAK135	ADP/ATP Ratio Assay Kit
MAK331	AF HDL and LDL/VLDL Assay Kit
MAK500	Alanine Assay Kit
MAK327	Aldehyde Dehydrogenase Inhibitor Screening Kit
MAK530	Alkaline Phosphatase (ALP) Assay Kit
MAK334	Antioxidant Assay Kit
MAK545	Beta-Lactamase Activity Assay Kit
MAK457	Caspase-3 Assay Kit
MAK106	CETP Activity Assay Kit
MAK305	CETP RP Activity Assay Kit
MAK325	Cholesterol Uptake Assay Kit
MAK536	Cytotoxicity Assay Kit
MAK450	Glucose Dehydrogenase Assay Kit
MAK451	Glucose-6-Phosphate Dehydrogenase Assay Kit

Kit #	Product
MAK452	Glucose-6-Phosphate Dehydrogenase Inhibitor Screening Kit
MAK330	Glutamate Assay Kit
MAK441	Kinase Assay Kit
MAK306	LCFC-LCAT Activity Assay Kit
MAK329	L-Lactate Assay Kit (Colorimetric)
MAK443	L-Lactate Assay Kit (Fluorometric)
MAK109	LPL Activity Assay Kit
MAK110	MTP Activity Assay Kit
MAK161	Multidrug Resistance Assay Kit
MAK121	Neuraminidase Assay Kit
MAK323	Nitric Oxide Synthase Inhibitor Screening Kit
MAK122	Phospholipid Assay Kit
MAK108	PLTP Activity Assay Kit
MAK332	Pyruvate Assay Kit

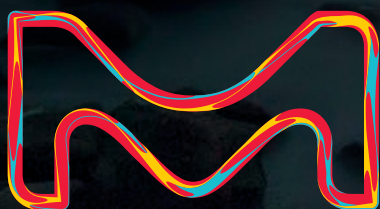


Navigate toward  
**Metabolomic  
 Discovery**

**Explore our portfolio:**

- Comprehensive metabolomic resources, reagents, libraries and kits
- Large collection of amino acid metabolites and carbohydrate metabolites
- Enzymatic-based kits for the quantitation of nutrients and metabolites

[SigmaAldrich.com/metabolomics](https://SigmaAldrich.com/metabolomics)



© 2019 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. Merck and the Merck logo 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.

**Sigma-Aldrich®**  
 Lab & Production Materials

The Life Science  
 business of Merck  
 operates as  
 MilliporeSigma in  
 the U.S. and Canada.

# Sigma-Aldrich®

Lab & Production Materials

Merck KGaA  
Frankfurter Strasse 250  
64293 Darmstadt, Germany

**SigmaAldrich.com**

**To place an order or receive technical assistance**

Order/Customer Service: **[SigmaAldrich.com/order](https://SigmaAldrich.com/order)**

Technical Service: **[SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice)**

Safety-related Information: **[SigmaAldrich.com/safetycenter](https://SigmaAldrich.com/safetycenter)**

© 2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. Merck and the vibrant M 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.

LIL No. MK\_CA3044EN Ver. 4.0 50422 05/2024