

Life Science Reagents for Western Blotting from Sigma-Aldrich®

Nucleic Acid
Blotting
DNA & RNA
(Southern) (Northern)

Nucleic Acid
Preparation
DNA & RNA

Nuclear Magnetic
Resonance (NMR)

Cloning

Nucleic Acid
Electrophoresis
DNA & RNA

Recombinant
Protein Expression
& Purification

**Protein
Detection**
Western Blots
& ELISAs

Proteomics

Transfection

Cell
Culture

Nucleic Acid
Amplification
PCR, QPCR, RT-PCR

Stem Cell
Culture

Cell
Signaling

Cell-Based
Assays

Protein
Electrophoresis
PAGE & 2D PAGE

Genotyping

We Ripple Through Everything You Do!

- Extensive reagent offering to support your Western Blot based research needs.
- Market leading manufacturer of biological buffers
- Quality and availability assured
- Market leading Technical Service
- Leading Web support tools:

Sigma Antibodies — sigma.com/antibodies

Molecular Weight Markers — sigma-aldrich.com/markers

Virtual Stockroom — sigma-aldrich.com/virtualstockroom

Life Science Reagents for Western Blotting from Sigma-Aldrich®

Sigma®'s New Prestige Antibodies™ are the most highly characterized antibodies in the industry.

- Designed and validated by the Human Proteome Resource (HDR) project
- Standardized in universal protocols
- Over 500 immunohistological images for every antibody
- Publicly available data on the Web
- Tested in multiple assurance steps to be the most specific products on the market

To learn more, please visit
sigma-aldrich.com/prestige.

Cat. No.	Description
Acids & Bases	
H1758	Hydrochloric acid Molecular Biology Grade, 36.5-38.0 %
S8045	Sodium hydroxide SigmaUltra, ≥98%, pellets (anhydrous)
T8657	Trichloroacetic acid for electrophoresis, suitable for fixing solution (for IEF and PAGE gels), ≥99%
Acrylamides	
A3553	Acrylamide for electrophoresis, ≥99%, powder
A9099	Acrylamide for molecular biology, ≥99%
A4058	Acrylamide solution 40%, for electrophoresis, sterile-filtered
A8887	Acrylamide suitable for electrophoresis, ≥99%
M7279	N,N'-Methylenebisacrylamide for electrophoresis, ≥98%, powder
M2022	N,N'-Methylenebisacrylamide for molecular biology, ≥98%
Amino Acids	
G8898	Glycine for electrophoresis, ≥99%
G6529	L-Glutathione reduced SigmaUltra, 98-100%
L8912	L-Leucine meets EP, JP, USP testing specifications, cell culture tested, ≥98.5%, from non-animal source
A8199	N-Acetyl-L-cysteine SigmaUltra, ≥99%
Buffers	
B8660	BICINE SigmaUltra, ≥99%
C6070	CAPS SigmaUltra, pH 3.0-7.0 (20 °C, 0.5 M in H ₂ O), ≥98.0% as NaOH
M8902	MES hemisodium salt ≥98%
M2933	MES hydrate Biotechnology Performance Certified, plant cell culture tested, ≥99.5%
M3058	MES sodium salt Biotechnology Performance Certified
M1317	MES solution Biotechnology Performance Certified, 1 M, cell culture tested
M3183	MOPS Biotechnology Performance Certified, cell culture tested, ≥99.5%
M9027	MOPS hemisodium salt ≥99%
M9024	MOPS sodium salt Biotechnology Performance Certified, cell culture tested, ≥99.5%
M1442	MOPS solution Biotechnology Performance Certified, cell culture tested, 1 M
P5493	Phosphate buffered saline 10× Concentrate, Biotechnology Performance Certified, cell culture tested
P3813	Phosphate buffered saline powder, pH 7.4
P3563	Phosphate buffered saline pH 7.4, contains Tween 20, dry powder
P4809	Phosphate-Citrate Buffer tablet
S3014	Sodium chloride for molecular biology, ≥98%
S6316	Sodium chloride solution 5 M, Biotechnology Performance Certified, cell culture tested
T1074	Tricine Biotechnology Performance Certified, 1 M in solution, cell culture tested, 0.2 µm filtered

Cat. No.	Description
T5816	Tricine Biotechnology Performance Certified, cell culture tested, ≥99%
T8793	Tris Buffered Saline, pH 8.0, with 3% nonfat Milk powder
T6789	Tris Buffered Saline, with BSA, pH 8.0
T6664	Tris Buffered Saline pH 8.0, powder
T9039	Tris Buffered Saline, with TWEEN® 20, pH 8.0 powder
T4904	Tris-Glycine Buffer 10× Concentrate
T6066	Trizma® base Biotechnology Performance Certified, meets EP, USP testing specifications, cell culture tested, ≥99.9%
T5941	Trizma hydrochloride for molecular biology, Biotechnology Performance Certified, cell culture tested, ≥99.0%

Detergents

L3771	Sodium dodecyl sulfate for electrophoresis, ≥98.5%
L4390	Sodium dodecyl sulfate for molecular biology, ≥98.5%
T8787	Triton® X-100 for molecular biology
P5927	TWEEN 20 for electrophoresis, suitable for solubilizing agent of membrane proteins and as a blocking reagent in immunoblotting
P9416	TWEEN 20 for molecular biology, viscous liquid

General Reagents

M7154	2-Mercaptoethanol for electrophoresis
A2504	6-Aminocaproic acid ≥99% (titration), powder
B4287	Albumin from bovine serum for molecular biology, powder
B2518	Albumin, Acetylated from bovine serum lyophilized powder, for molecular biology
A9647	Albumin from bovine serum Heat-shocked fractionate, pH ~7 (1 % (w/v) in H ₂ O), ≥96% (agarose gel electrophoresis), lyophilized powder
A3678	Ammonium persulfate for electrophoresis, ≥98%
B4639	Biotin cell culture tested, insect cell culture tested, ≥99%, powder
D9163	DL-Dithiothreitol for electrophoresis, ≥99%
D9779	DL-Dithiothreitol for molecular biology, ≥98% (TLC), ≥99% (titration)
E5134	EDTA disodium salt dihydrate for molecular biology, ≥99%
E1510	Ethidium bromide solution 10 mg/mL, for molecular biology, aqueous solution
F3526	Fluorescent Molecular Weight Marker mol wt. 20,100-205,000 Da
G7663	Gelatin blocking buffer for molecular biology, powder blend
G4004	Glutaraldehyde solution 25 wt. % in H ₂ O
G8773	Glycerol for electrophoresis, ≥99%
G3272	Guanidine hydrochloride for molecular biology, ≥99%
216763	Hydrogen peroxide solution ACS reagent, 30 wt. % in H ₂ O
494437	Methanol biotech. grade, 99.93%
M1770	Methanol (for solid phase protein sequence analysis)
S7920	Sodium fluoride SigmaUltra, ≥99%
T9281	TEMED for electrophoresis, ~99%

Cat. No.	Description
U6504	Urea for electrophoresis
W3513	Water Biotechnology Performance Certified
W4502	Water Molecular Biology Reagent
W0138	Western Blocker™ Solution for HRP detection systems

Markers

SDS7B2	Prestained Molecular Weight Marker mol wt. 26,600-180,000 Da
C2242	Chemichrome™ Western Control
B2787	Biotinylated Molecular Weight Marker mol wt. 6,500-180,000 Da
C4861	ColorBurst™ Electrophoresis Marker High Range, mol wt. 30,000-220,000 Da
C1992	ColorBurst Electrophoresis Marker mol wt. 8,000-220,000 Da
C3187	Color Marker Low Range, mol wt. 6,500-45,000 Da
C6210	Color Marker Ultra-low Range (M.W. 1,060-26,600)

Membranes

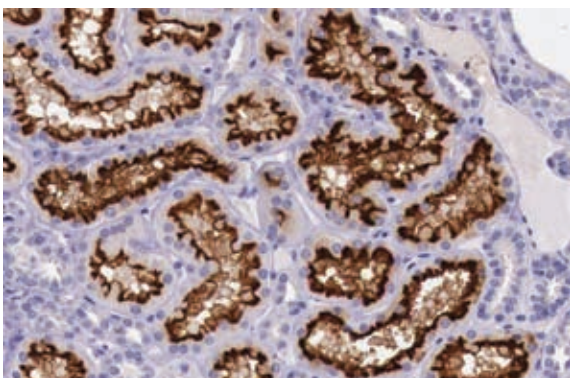
P7796	Blotting paper, pure cellulose Extra Thick, size 7 cm × 10 cm
Z671088	Whatman® Westran® PVDF membranes Westran Clear Signal, Rolls, pore size 0.45 µm, size 30 cm × 3 m
Z671045	Whatman Westran PVDF membranes Westran Clear Signal, Sheets, pore size 0.45 µm, size 10 cm × 10 cm
Z671053	Whatman Westran PVDF membranes Westran Clear Signal, Sheets, pore size 0.45 µm, size 15 cm × 15 cm
Z671061	Whatman Westran PVDF membranes Westran Clear Signal, Sheets, pore size 0.45 µm, size 20 cm × 20 cm

Substrates

T0565	3,3',5,5'-Tetramethylbenzidine (TMB) Liquid Substrate System for Membranes
D7304	3,3'-Diaminobenzidine (DAB) Liquid Substrate System tetrahydrochloride
C8302	4-Chloro-1-naphthol solution
AEC101	AEC Staining Kit liquid
90101	Ampliflu Red BioChemika, for fluorescence, ≥98.0%
B1911	BCIP®/NBT Liquid Substrate System
B6404	BCIP/NBT solution, premixed
C0712	CDP-Star® Chemiluminescent Substrate 0.25 mM in H ₂ O
CPS1120	Chemiluminescent Peroxidase Substrate-1 sufficient for 120 mL substrate
CPS1300	Chemiluminescent Peroxidase Substrate-1 sufficient for 300 mL substrate
CPS160	Chemiluminescent Peroxidase Substrate-1 sufficient for 60 mL substrate
D4418	SIGMAFAST™ 3,3'-Diaminobenzidine tablets tablet, To prepare 15 mL
B5655	SIGMAFAST BCIP/NBT tablet
D0426	SIGMAFAST DAB with Metal Enhancer tablet
F4523	SIGMAFAST Fast Red TR/Naphthol AS-MX Tablets tablet, To prepare 10 mL

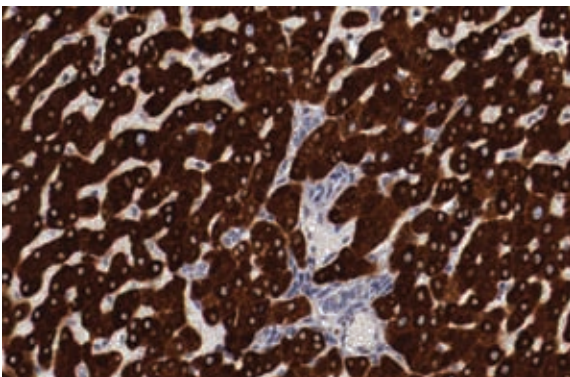
Prestige Antibodies™ from Sigma-Aldrich®

Antibodies that work, right out of the box!



Anti-ACE2 shows strong cytoplasmic and/or membranous staining of cells in tubuli in kidney.

Cat. No. HPA000288



Anti-MBL2 shows strong cytoplasmic and/or membranous staining of hepatocytes in liver.

Cat. No. HPA000207

The Most highly characterized antibodies in the industry:

- Designed & validated by the Human Proteome Resource (HPR)
- Standardized in universal protocols
- Over 500 immunohistology images for every antibody
- Publicly available database on the web



The Human Protein Atlas (*proteinatlas.org*) displays a collection of over 2.9 million IHC images of normal and cancer tissues, cell lines and primary cells. Prestige Antibodies, targeting approximately 1,600 human proteins, have been screened for protein profiling.

To search or browse by Gene name or ID, please visit **Prestige Antibodies at sigma.com/prestige.**

World Headquarters

3050 Spruce St., St. Louis, MO 63103
(314) 771-5765
sigma-aldrich.com

Order/Customer Service (800) 325-3010 • Fax (800) 325-5052

Technical Service (800) 325-5832 • sigma-aldrich.com/techservice

Development/Bulk Manufacturing Inquiries SAFC™ (800) 244-1173

Accelerating Customers' Success through Leadership in Life Science, High Technology and Service

©2008 Sigma-Aldrich Co. All rights reserved. SIGMA, SAFC, SAFC, SIGMA-ALDRICH, ALDRICH, FLUKA, and SUPELCO are trademarks belonging to Sigma-Aldrich Co. and its affiliate Sigma-Aldrich Biotechnology, L.P. Sigma brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.

Trizma, BCIP, REDTaq, Sigma and Sigma-Aldrich are registered trademarks of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co. JumpStart, AccuTaq, ColorBurst, Chemichrome, SIGMAFAST and Western Blocker are trademarks of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co. Triton is a registered trademark of Union Carbide Corp. TWEEN is a registered trademark of Uniqema, a business unit of ICI Americas, Inc. CDP-Star is a registered trademark of Tropix, Inc. Whatman and Westran are registered trademarks of Whatman International Ltd. Prestige Antibodies is a trademark of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co. Sigma and Sigma-Aldrich are registered trademarks of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co.

KSS
04646
0078