

Neuroscience Research Primary Antibodies You Can Rely On



Generate reproducible western blots with PrecisionAb™ Antibodies

High Specificity and Sensitivity

- Screened using 12 cell lysates expressing endogenous protein levels

Superior Data Reproducibility

- Stringent lot-to-lot QC requirements ensure reproducible performance

Less Time and Money for Optimization

- Trial size available
- Positive controls and protocols available



Western blot analysis of 12 whole cell lysates probed with goat anti-human superoxide dismutase (Cu-Zn) antibody (1:1,000, Bio-Rad, catalog #VPA00070) followed by detection with HRP-conjugated donkey anti-sheep/goat IgG (1:10,000, bio-rad-antibodies.com, #STAR88P) and visualized on the ChemiDoc™ MP Imaging System with 12 second exposure.

Product Formats



Trial size with positive control lysate
Minimum 2 western blots (20 µl)



Antibody only
Minimum 10 western blots (100 µl)



Order your trial size antibody at bio-rad.com/web/NeuroPrecisionAbs



Ordering Information

Catalog #	Description	Catalog #	Description
Polyclonal Anti-Human Antibody		Monoclonal Anti-Human Antibody	
VPA00009	Rabbit Anti-Lysine-Specific Histone Demethylase	VMA00003	Mouse Anti-Beta 2 Microglobulin
VPA00017	Rabbit Anti-AIF	VMA00014	Human Anti-Actin Beta
VPA00019	Goat Anti-Fyn	VMA00037	Mouse Anti-CD71 (Soluble)
VPA00028	Goat Anti-PPP2R5D	VMA00046	Mouse Anti-GAPDH
VPA00065	Goat Anti-DCTN1	VMA00048	Mouse Anti-Actin Beta
VPA00096	Goat Anti-Calnexin	VMA00049	Mouse Anti-Actin Gamma
VPA00103	Goat Anti-VPS35	VMA00054	Mouse Anti-EBP50
VPA00118	Goat Anti-LRRK2	VMA00065	Mouse Anti-C9orf72
VPA00119	Goat Anti-Lamin B1	VMA00069	Mouse Anti-Neuron Specific Enolase
VPA00138	Goat Anti-Clathrin Light Chain A/B	VMA00073	Mouse Anti-Synaptophysin
VPA00143	Goat Anti-SDCBP	VMA00074	Mouse Anti-Lysosome Membrane Protein 2
VPA00159	Goat Anti-MTHFD1L	VMA00076	Mouse Anti-PSMD1
VPA00173	Rabbit Anti-Histone Deacetylase 5	VMA00079	Mouse Anti-Spectrin Alpha Chain
VPA00189	Rabbit Anti-GGA3	VMA00080	Mouse Anti-Adenosine A2A Receptor
VPA00216	Rabbit Anti-FUSE-Binding Protein 2	VMA00089	Mouse Anti-Haptoglobulin
VPA00243	Rabbit Anti-FUSE-Binding Protein 1	VMA00109	Mouse Anti-AKR1A1
VPA00014	Rabbit Anti-CD54	VMA00129	Mouse Anti-Catalase
VPA00030	Goat Anti-TDP-43	VMA00131	Mouse Anti-Heat Shock Protein 40
VPA00044	Goat Anti-CD195	VMA00135	Mouse Anti-USP5
VPA00067	Goat Anti-PEBP1	VMA00141	Mouse Anti-DLD
VPA00068	Goat Anti-ABCC4	VMA00142	Mouse Anti-PSMB4
VPA00069	Goat Anti-Superoxide Dismutase (Mn)	VMA00149	Mouse Anti-PSMF1
VPA00070	Goat Anti-Superoxide Dismutase (Cu-Zn)	VMA00152	Mouse Anti-Neuron Specific Enolase
VPA00097	Goat Anti-Apolipoprotein A1	VMA00154	Mouse Anti-NTRK3
VPA00098	Goat Anti-FADS1	VMA00161	Mouse Anti-EPN2
VPA00099	Goat Anti-CD20	VMA00184	Mouse Anti-TCF12
VPA00152	Goat Anti-Plasminogen	VMA00189	Mouse Anti-Neurofilament
VPA00154	Goat Anti-SNAP-25	VMA00196	Mouse Anti-UQCRC1
VPA00168	Rabbit Anti-ROBO1	VMA00203	Mouse Anti-ALAS1
VPA00187	Rabbit Anti-GAPDH	VMA00290	Mouse Anti-Lysine-Specific Demethylase 3A
VPA00193	Rabbit Anti-Myeloperoxidase	VMA00303	Mouse Anti-PIN1
VPA00200	Sheep Anti-Alpha 2 Macroglobulin	VMA00316	Mouse Anti-Superoxide Dismutase (Cu-Zn)
VPA00240	Rabbit Anti-CD271	VMA00346	Mouse Anti-AKR1C2
VPA00244	Rabbit Anti-Histone Acetyltransferase 1	VMA00226	Mouse Anti-C14orf166
VPA00090	Rabbit Anti-UBC13	VMA00042	Mouse Anti-Heat Shock Protein 70
VPA00234	Rabbit Anti-SMURF1	VMA00155	Mouse Anti-SATB1
VPA00203	Rabbit Anti-TRIM28	VMA00177	Mouse Anti-SUPT3H
VPA00206	Rabbit Anti-UBE2K	VMA00249	Mouse Anti-Survival Motor Neuron Protein
VPA00051	Goat Anti-TRAF2	VMA00268	Mouse Anti-Myocyte-Specific Enhancer Factor 2D
VPA00026	Rabbit Anti-DNA Methyltransferase 1	VMA00300	Mouse Anti-USP14
VPA00190	Rabbit Anti-POLB	VMA00333	Mouse Anti-14-3-3 Beta/Alpha
VPA00023	Rabbit Anti-MCM6	VMA00335	Mouse Anti-Creatine Phosphokinase
VPA00060	Goat Anti-APEX1	VMA00342	Mouse Anti-GSK3 Alpha/Beta
VPA00195	Rabbit Anti-UBA1	VMA00349	Mouse Anti-Cystatin B
VPA00151	Goat Anti-HR23B	VMA00350	Mouse Anti-Carbonic Anhydrase VIII
VPA00130	Goat Anti-POU2F1	VMA00247	Mouse Anti-E3 Ubiquitin-Protein Ligase CBL

Visit bio-rad.com/web/NeuroPrecisionAbs to view our complete listing.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 177 Belgium 32 (0)3 710 53 00 Brazil 55 11 3065 7550
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 45 44 52 10 00 Finland 358 09 804 22 00
France 33 01 47 95 69 65 Germany 49 89 31 884 0 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 972 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 31 (0)318 540 666
New Zealand 64 9 415 2280 Norway 47 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04
Singapore 65 6415 3188 South Africa 27 (0) 861 246 723 Spain 34 91 590 5200 Sweden 46 08 555 12700 Switzerland 41 026674 55 05
Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

