

# Cancer Research Primary Antibodies You Can Rely On



Generate reproducible western blots with PrecisionAb™ Antibodies

## High Specificity and Sensitivity

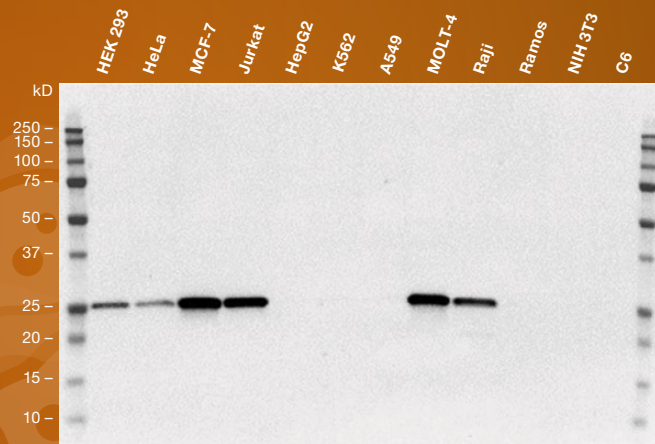
- Screened using 12 cell lysates expressing endogenous protein levels

## 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 mouse anti-human Bcl-2 antibody (Bio-Rad, catalog #VMA00017) followed by detection with HRP-conjugated goat anti-mouse IgG (1:10,000, Bio-Rad, catalog #1705047) and visualized on the ChemiDoc™ MP Imaging System with 67 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/CancerPrecisionAbs](http://bio-rad.com/web/CancerPrecisionAbs)



## Ordering Information

Catalog #	Description	Catalog #	Description
<b>Polyclonal Anti-Human Antibody</b>		<b>Monoclonal Anti-Human Antibody</b>	
VPA00068	Goat Anti-ABCC4	VMA00012	Human Anti-Cytokeratin 14
VPA00027	Goat Anti-ATF2	VMA00010	Human Anti-MEK 1
VPA00116	Goat Anti-Catenin Alpha 1	VMA00008	Human Anti-Vimentin
VPA00132	Goat Anti-Cytochrome B245	VMA00253	Mouse Anti-AKT1
VPA00061	Goat Anti-DDX6	VMA00066	Mouse Anti-ANAPC7
VPA00003	Goat Anti-ErbB3	VMA00017	Mouse Anti-Bcl-2
VPA00059	Goat Anti-GRP94	VMA00164	Mouse Anti-Calpain-2
VPA00057	Goat Anti-MIF	VMA00129	Mouse Anti-Catalase
VPA00110	Goat Anti-NADPH Oxidase 1	VMA00098	Mouse Anti-Cdc2
VPA00067	Goat Anti-PEBP1	VMA00022	Mouse Anti-Cdk6
VPA00056	Goat Anti-Peroxiredoxin 1	VMA00023	Mouse Anti-Cdk7
VPA00194	Goat Anti-PPP2R1A	VMA00156	Mouse Anti-Cytochrome P450 2E1
VPA00028	Goat Anti-PPP2R5D	VMA00117	Mouse Anti-Cytochrome P450 3A4
VPA00155	Goat Anti-PSMB8	VMA00002	Mouse Anti-Cytokeratin 19
VPA00069	Goat Anti-Superoxide Dismutase (Mn)	VMA00061	Mouse Anti-EGF Receptor
VPA00146	Goat Anti-TPI1	VMA00134	Mouse Anti-EIF2S1
VPA00111	Goat Anti-UBE2 L3	VMA00120	Mouse Anti-EIF4B
VPA00017	Rabbit Anti-AIF	VMA00147	Mouse Anti-ErbB2
VPA00192	Rabbit Anti-AKT2	VMA00157	Mouse Anti-Fibronectin
VPA00077	Rabbit Anti-Annexin V	VMA00131	Mouse Anti-Heat Shock Protein 40
VPA00086	Rabbit Anti-ATG16	VMA00041	Mouse Anti-Heat Shock Protein 60
VPA00230	Rabbit Anti-ATG7	VMA00042	Mouse Anti Heat Shock Protein 70
VPA00164	Rabbit Anti-Catenin Beta	VMA00081	Mouse Anti-Heat Shock Protein 90 Alpha
VPA00254	Rabbit Anti-COP9 Signalosome Complex Subunit 1	VMA00082	Mouse Anti-Heat Shock Protein 90 Beta
VPA00026	Rabbit Anti-DNA Methyltransferase 1	VMA00102	Mouse Anti-H-FABP
VPA00167	Rabbit Anti-ErbB3	VMA00113	Mouse Anti-hnRNP F
VPA00076	Rabbit Anti-Galectin-3	VMA00261	Mouse Anti-hnRNP L
VPA00205	Rabbit Anti-GSK3 Alpha	VMA00260	Mouse Anti-hnRNP M
VPA00175	Rabbit Anti-Histone Deacetylase 1	VMA00040	Mouse Anti-H-Ras
VPA00171	Rabbit Anti-Histone Deacetylase 6	VMA00078	Mouse Anti-IGFBP-7
VPA00133	Rabbit Anti-IGFBP-7	VMA00140	Mouse Anti-KEAP1
VPA00084	Rabbit Anti-IKK Gamma	VMA00266	Mouse Anti-Laminin Beta 3
VPA00252	Rabbit Anti-MMP-14 Hemopexin	VMA00268	Mouse Anti-Myocyte-Specific Enhancer Factor 2D
VPA00174	Rabbit Anti-mTOR	VMA00150	Mouse Anti-NF Kappa B Inhibitor Alpha
VPA00193	Rabbit Anti-Myeloperoxidase	VMA00262	Mouse Anti-NUMB
VPA00190	Rabbit Anti-POLB	VMA00019	Mouse Anti-p53
VPA00227	Rabbit Anti-Proto-Oncogene c-Rel	VMA00018	Mouse Anti-PCNA
VPA00210	Rabbit Anti-PSMB5	VMA00263	Mouse Anti-PI-3 Kinase C2 Subunit Alpha
VPA00247	Rabbit Anti-RPL7	VMA00016	Mouse Anti-Poly (ADP-Ribose) Polymerase 1
VPA00248	Rabbit Anti-RPL9	VMA00111	Mouse Anti-Proliferation Associated Protein 2G4
VPA00246	Rabbit Anti-RPS14	VMA00076	Mouse Anti-PSMD1
VPA00087	Rabbit Anti-Smac/DIABLO	VMA00124	Mouse Anti-RAF-1
VPA00182	Rabbit Anti-SMAD2	VMA00168	Mouse Anti-RALB
VPA00166	Rabbit Anti-SMAD3	VMA00158	Mouse Anti-RPA1
VPA00234	Rabbit Anti-SMURF1	VMA00027	Mouse Anti-SOCS-3
VPA00250	Rabbit Anti-Splicing Factor 3b Subunit 2	VMA00242	Mouse Anti-STAT3
VPA00008	Rabbit Anti-STAT1 Alpha	VMA00306	Mouse Anti-SUMO-1

Visit [bio-rad.com/web/CancerPrecisionAbs](http://bio-rad.com/web/CancerPrecisionAbs) to view our complete listing.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [bio-rad.com](http://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 41 026 674 55 05  
Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

