



Primary Antibodies You Can Rely On



Experience superior and reproducible western blot performance with PrecisionAb™ Antibodies.

Expect High Specificity and Sensitivity

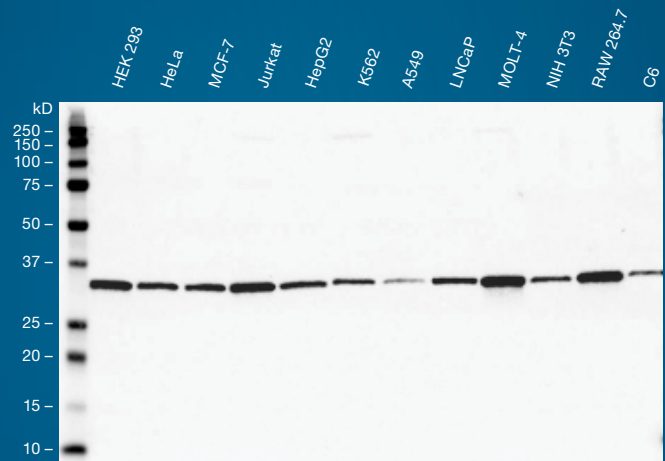
- Screened using 12 cell lysates expressing endogenous protein levels

Reliably Replicate Data


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

Save Time and Money Optimizing Experiments

- Trial size for each antibody
- Positive controls and protocols available to help set up experiments



Western blot analysis of 12 whole cell lysates probed with mouse anti-human PCNA primary antibody followed by detection with goat anti-mouse secondary antibody and visualized on the ChemiDoc™ MP Imaging System with Clarity™ Western ECL Substrate.

 Order your trial size antibody at [bio-rad.com/web/PrecisionAbs](https://www.bio-rad.com/web/PrecisionAbs)



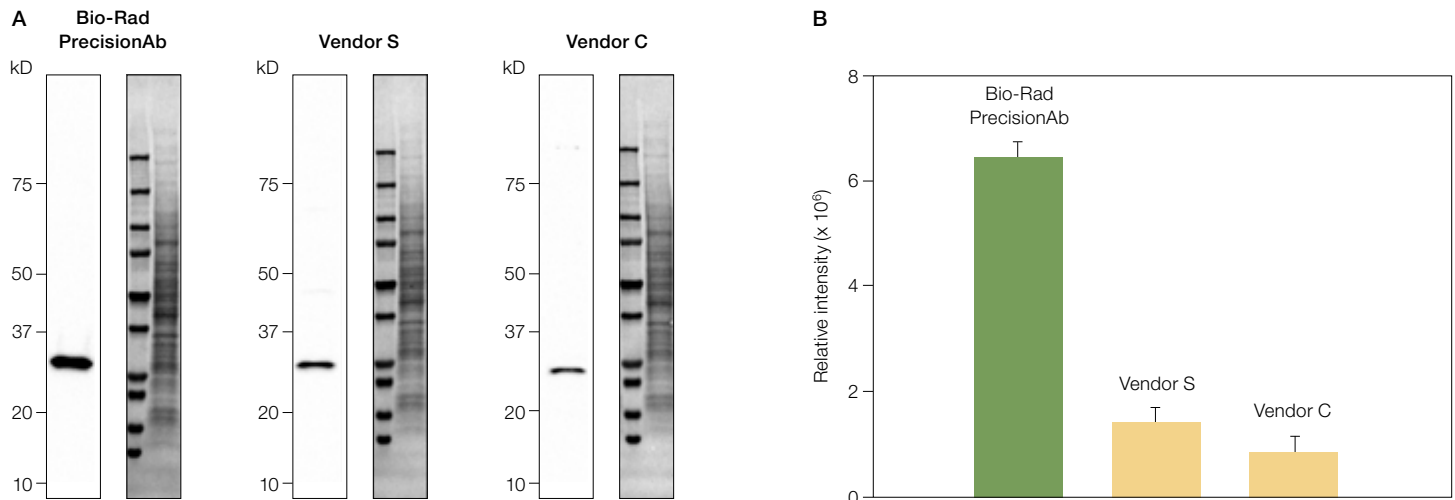
Customer Testimonial



“Antibodies provided from Bio-Rad were of very high quality for 80% of the antibodies tested, suggesting most antibodies will be great for western blotting.”

—Aldrin Gomes, PhD
Associate Professor, UC Davis

Comparative Data



Side-by-side testing of anti-human BCL-2 antibodies on Jurkat cell lysate. **A**, western blot and stain-free total protein imaging data for each antibody; **B**, relative band intensities for each of the western blots. The PrecisionAb Antibody shows ~5x greater signal intensity even at a lower dilution. Primary antibody dilutions: PrecisionAb, 1:1,000; Vendor S, 1:1,000; Vendor C, 1:500. Secondary antibody dilution was 1:10,000 for all blots.

Product Formats

bio-rad.com/web/PrecisionAbs



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



Antibody only
Minimum 10 western blots (100 μ l)



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 41 026 674 55 05 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311
United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

