



# Antibodies You Can Rely On



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

## Expect High Specificity and Sensitivity

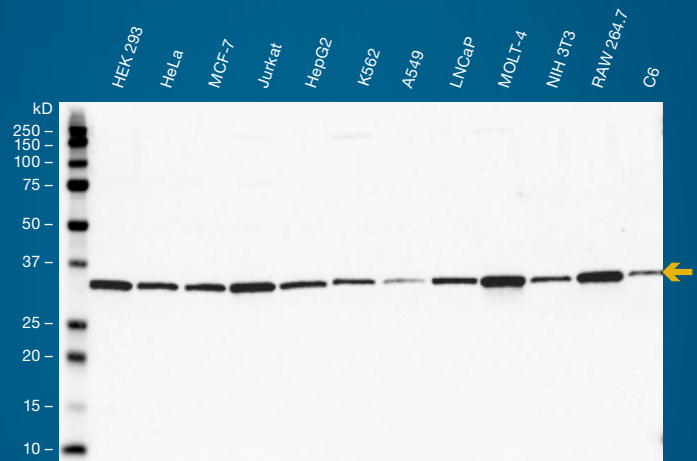
- Screened using 12 cell lysates expressing endogenous protein levels

## 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/PrecisionAbs](https://www.bio-rad.com/PrecisionAbs)



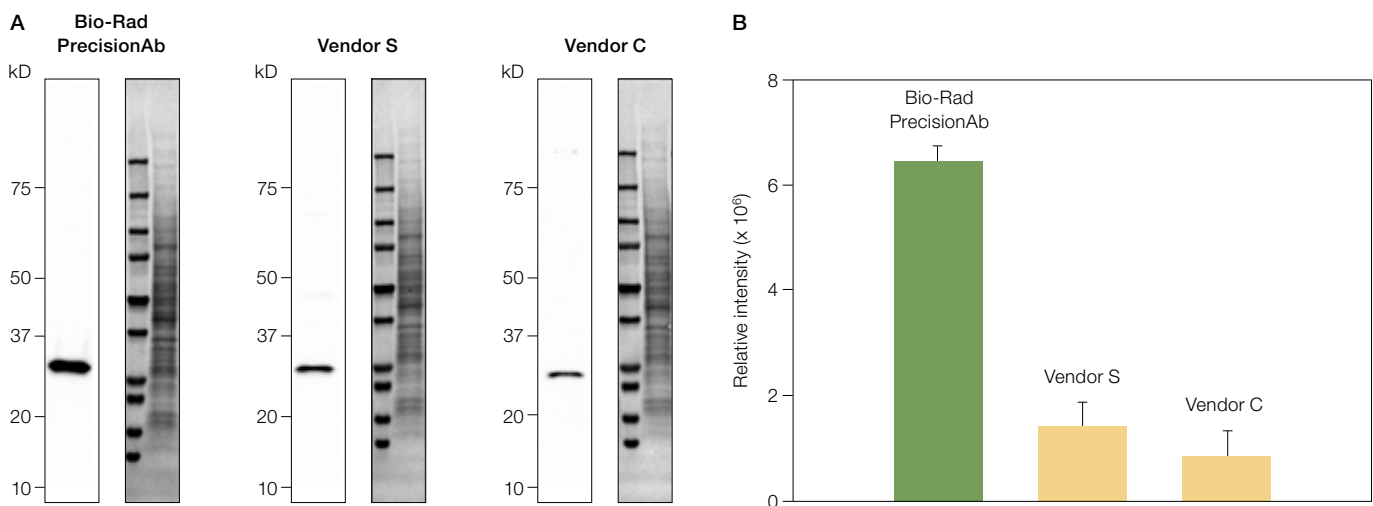
## Customer Testimonial



“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/PrecisionAbs](http://bio-rad.com/PrecisionAbs)



**Trial size**  
Minimum two western blots (20 µl)



**Antibody with control**  
Minimum ten western blots (100 µl)



**Antibody only**  
(100 µl)

**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 66 662 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

