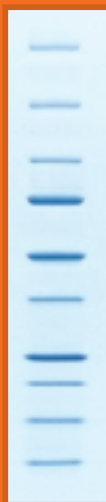


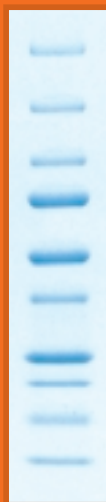
Precision Plus Protein™ Standards

The Sharpest Standards in the Box

Now with dual color and more reference bands.



Precision Plus Protein
Unstained
Standards



Precision Plus Protein
All Blue
Standards



Precision Plus Protein
Dual Color
Standards

BIO-RAD

Protein standards have never looked better.

New Precision Plus Protein standards offer unsurpassed band sharpness, accurate molecular weight estimations, lot-to-lot consistency, and instant orientation for gel electrophoresis and western blotting. Precision Plus Protein standards contain ten highly purified recombinant proteins, including three reference bands, in an ideal range of bands from 10 kD to 250 kD. Precision Plus Protein standards are available in unstained, all blue, and dual color options, providing three new ways to get sharper looking results.



New! Precision Plus Protein dual color standards offer:

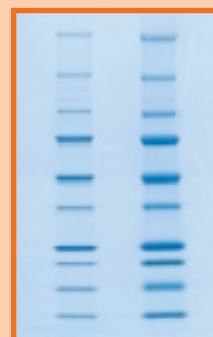
- > Prestained pink and blue bands
- > Unsurpassed band sharpness
- > Superior lot-to-lot consistency
- > Ideal ten-band range from 10 kD to 250 kD
- > Three high-intensity reference bands: 25 kD, 50 kD, and 75 kD
- > Molecular weights that never change, so band sizes are easy to remember
- > Migration pattern that matches unstained standards
- > A ready-to-use, load-and-go format



Dual color standards are distinct and easy to recognize

New! Precision Plus Protein all blue standards offer:

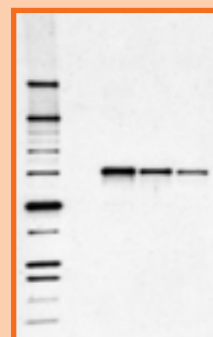
- > Prestained blue bands
- > Unsurpassed band sharpness
- > Superior lot-to-lot consistency
- > Ideal ten-band range from 10 kD to 250 kD
- > Three high-intensity reference bands: 25 kD, 50 kD, and 75 kD
- > Molecular weights that never change, so band sizes are easy to remember
- > Migration pattern that matches unstained standards
- > A ready-to-use, load-and-go format



Perfectly comigrating unstained (left) and all blue (right) standards

New! Precision Plus Protein unstained standards offer:

- > Absolute molecular weight accuracy, confirmed by mass spectrometry
- > Ideal ten-band range from 10 kD to 250 kD
- > Three high-intensity reference bands: 25 kD, 50 kD, and 75 kD
- > Quantitation capability
- > *Strep*-tag amino acid sequence for detection and molecular weight determinations on western blots
- > Superior lot-to-lot consistency
- > Molecular weights that never change, so band sizes are easy to remember
- > A ready-to-use, load-and-go format



Sharp, clear bands on gels and western blots

Precision Plus Protein unstained standards (left lane) incubated with StrepTactin-HRP and detected with the Immun-Star™ HRP chemiluminescent kit.

Ordering Information

Catalog #	Description
161-0374	Precision Plus Protein Dual Color Standards, 50 applications
161-0373	Precision Plus Protein All Blue Standards, 50 applications
161-0363	Precision Plus Protein Unstained Standards, 100 applications
161-0380	Precision Protein™ StrepTactin-HRP conjugate, 0.3 ml
161-0382	Precision Protein StrepTactin-AP conjugate, 0.3 ml

Strep-tag and StrepTactin are trademarks of Institut für Bioanalytik GmbH. *Strep*-tag technology for western blot detection is covered by US patent 5,506,121 and by UK patent 2,272,698.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191 Canada (905) 712-2771 China (86-21) 63052255 Denmark 45 44 52-1000 Finland 358 (0)9 804 2200 France 01 47 95 69 65 Germany 089 318 84-177 Hong Kong 852-2789-3300 India (91-124) 6398112/113/114, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 47-23-38-41-30 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-2729877 South Africa 00 27 11 4428508 Spain 590 5200 Sweden 46 (0)8-55 51 27 00 Switzerland 061 717-9555 United Kingdom 0800-181134