



## PROTEIN STANDARDS

# Precision Plus Protein™ WesternC™ Standards

- Colorimetric, chemiluminescent, and fluorescent properties in one standard
- Consistent, identical migration patterns within the Precision Plus Protein standards family
- Accurate molecular weight determination in gels and blots
- Visual gel monitoring and blot orientation

## Versatile, Multi-Application Protein Standards: Colorimetric, Chemiluminescent, and Fluorescent Properties All in One

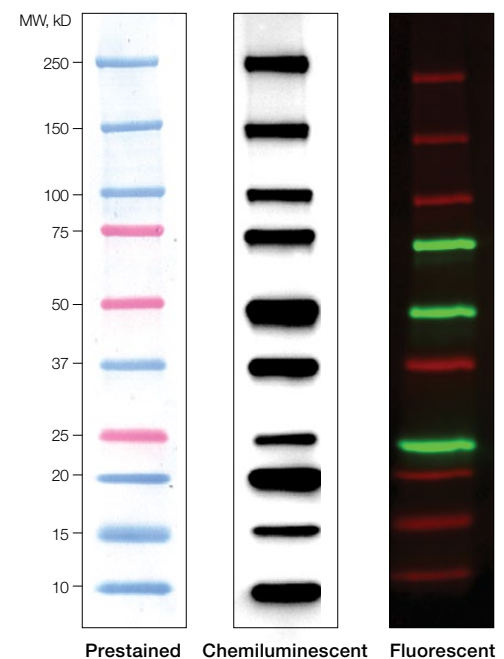
Standards are an integral part of every electrophoresis experiment because they help identify and characterize the molecules separated in a gel. Optimal protein standards will offer a combination of sharp and bright bands, accurate molecular weight estimation, reproducible migration pattern, and the ability to confirm electrophoretic separation. They can also verify transfer efficiency.

Traditional protein standards incorporate natural proteins of predetermined molecular weights for sizing. Bio-Rad's Precision Plus Protein standards incorporate recombinant proteins that exhibit reproducible and consistent migration regardless of staining, providing exceptional band sharpness and reproducibility not achievable with native proteins. The Precision Plus Protein standards family includes WesternC, All Blue, Dual Color, Dual Xtra, Kaleidoscope™, and Unstained standards for various experimental applications.

Precision Plus Protein WesternC standards are versatile, multi-application standards that contain ten prestained, recombinant bands from 10–250 kD. These bands contain an integrated *Strep*-tag sequence that enables parallel detection on western blots using a StrepTactin-horseradish peroxidase (HRP) conjugate. These features allow electrophoretic confirmation on a gel, verification of transfer efficiency on a blot, and detection and molecular weight estimation on chemiluminescent and fluorescent western blots.

### Features and benefits include:

- Reproducible and accurate migration pattern, allowing molecular weight estimation
- Exceptional band sharpness, providing clarity on electrophoretic separation



Precision Plus Protein WesternC standards.

- Integrated *Strep*-tag sequence, offering chemiluminescent detection and molecular weight determination on western blots
- Natural fluorescent properties with excitation spectra in the red (~635 nm) and green (~532 nm) channels, allowing single and multiplex fluorescent detection
- Prestained, dual-colored bands, providing easy gel monitoring and blot orientation
- Ready-to-use formulation



## Ordering Information

Catalog #	Description
161-0363	<b>Precision Plus Protein Unstained Standards,</b> 1,000 µl, 100 applications
161-0396	<b>Precision Plus Protein Unstained Standards Value Pack,</b> 5 x 1,000 µl, 500 applications
161-0373	<b>Precision Plus Protein All Blue Standards,</b> 500 µl, 50 applications
161-0393	<b>Precision Plus Protein All Blue Standards Value Pack,</b> 5 x 500 µl, 250 applications
161-0374	<b>Precision Plus Protein Dual Color Standards,</b> 500 µl, 50 applications
161-0394	<b>Precision Plus Protein Dual Color Standards Value Pack,</b> 5 x 500 µl, 250 applications
161-0377	<b>Precision Plus Protein Dual Xtra Standards,</b> 500 µl, 50 applications
161-0397	<b>Precision Plus Protein Dual Xtra Standards Value Pack,</b> 5 x 500 µl, 250 applications StrepTactin conjugates for Precision Plus Protein Standards
161-0375	<b>Precision Plus Protein Kaleidoscope Standards,</b> 500 µl, 50 applications
161-0395	<b>Precision Plus Protein Kaleidoscope Standards Value Pack,</b> 5 x 500 µl, 250 applications
161-0376*	<b>Precision Plus Protein WesternC Standards,</b> 250 µl, 50 applications
161-0399*	<b>Precision Plus Protein WesternC Standards Value Pack,</b> 5 x 250 µl, 250 applications
161-0385*	<b>Precision Plus Protein WesternC Standards + HRP Pack,</b> 250 µl WesternC standard, 125 µl HRP conjugate, 50 applications
161-0398*	<b>Precision Plus Protein WesternC Standards + HRP Value Pack,</b> 5 x 250 µl WesternC standard, 5 x 125 µl HRP conjugate, 250 applications
161-0380	<b>Precision Protein StrepTactin-HRP Conjugate,</b> 300 µl, 150 applications
161-0381	<b>Precision Protein StrepTactin-HRP Conjugate,</b> 125 µl, 50 applications
161-0382	<b>Precision Protein StrepTactin-AP Conjugate,</b> 300 µl, 150 applications
161-0378	<b>Precision Plus Protein Standard Plugs,</b> Unstained, 24 applications

\* Note that StrepTactin (-HRP or -AP) conjugate is needed for colorimetric or chemiluminescence blots.

StrepTactin and *Strep*-tag are trademarks of Institut für Bioanalytik GmbH.

*Strep*-tag technology for western blot detection is covered by U.S. patent 5,506,121 and by UK patent 2,272,698. StrepTactin is covered by German patent application P 19641876.3. Bio-Rad Laboratories, Inc. is licensed by Institut für Bioanalytik GmbH to sell these products for research use only.

### NOTICE TO PURCHASER: LIMITED LICENSE

\*This product is sold under license from Life Technologies Corporation, Carlsbad, CA for use only by the buyer of the product. The buyer is not authorized to sell or resell this product or its components.



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 11 5044 5699  
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00  
France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666  
New Zealand 64 9 415 2280 Norway 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 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55  
Taiwan 886 2 2578 7189 Thailand 800 88 22 88 United Kingdom 020 8328 2000