

Related Products for Electrophoresis

Catalog #	Description
161-0363	Precision Plus Protein Unstained Standards, 1,000 μl
161-0396	Precision Plus Protein Unstained Standards Value Pack, 5 x 1,000 μl
161-0373	Precision Plus Protein All Blue Standards, 500 μl
161-0393	Precision Plus Protein All Blue Standards Value Pack, 5 x 500 μl
161-0375	Precision Plus Protein™ Kaleidoscope™ Standards, 500 μl
161-0395	Precision Plus Protein Kaleidoscope Standards Value Pack, 5 x 500 μl
161-0377	Precision Plus Protein Dual Xtra Standards, 500 μl
161-0397	Precision Plus Protein Dual Xtra Standards Value Pack, 5 x 500 μl
161-0385	Precision Plus Protein™ WesternC™ Pack, Precision Plus Protein WesternC Standards and StrepTactin-HRP, 50 applications
161-0398	Precision Plus Protein WesternC Value Pack, 250 applications
161-0747	4x Laemmli Sample Buffer, 10 ml
161-0786	Bio-Safe™ Coomassie Stain, 1 L

For more information, go to www.bio-rad.com.

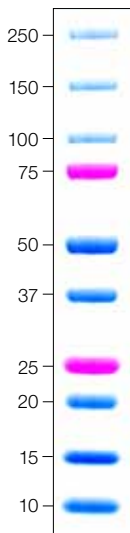
Precision Plus Protein™ Dual Color Standards Instruction Manual

Single unit
Catalog #161-0374
5-Pack value pack
Catalog #161-0394

Product shipped at
room temperature.

Store at 4°C for 6 months
or -20°C until expiration date.

Representative lot of
Precision Plus Protein
Dual Color standards
on a 4–20% Criterion™
TGX™ gel



Precision Plus Protein standards are the latest generation of recombinant proteins designed to provide a ladder of convenient, consistent sizes. These standards continue the tradition of accuracy and ease of use begun by Precision Protein™ standards, while also providing instant orientation, even greater accuracy, particularly in the 10–25 kD range, and two-color convenience.

Product Description

Precision Plus Protein Dual Color standards contain ten protein bands of 10, 15, 20, 25, 37, 50, 75, 100, 150, and 250 kD.

The Precision Plus Protein Dual Color standards represent a breakthrough in protein staining, with molecular weight shifts of <5% and minimal band broadening. Two pink reference bands (25 and 75 kD) and a more intense low molecular weight region (<25 kD) provide instant orientation of the gel pattern to quickly and easily identify proteins of interest.

The molecular weights of the Precision Plus Protein Dual Color standards are confirmed by migration in a Laemmli SDS-PAGE system. For electrophoretic determination of molecular weights at an accuracy of >95%, Precision Plus Protein Unstained standards (catalog #161-0363) should be used.

These standards are also compatible with the XT buffer system, but should be used only for molecular weight estimation. For molecular weight determination on XT gels, use the Precision Plus Protein Unstained (catalog #161-0363) or All Blue (catalog #161-0373) standards.

Instructions for Use

Precision Plus Protein Dual Color standards are provided in sample buffer, ready to load, with no dilution required. Allow the tube to reach room temperature and thoroughly mix before use. This will ensure that any solids that may have precipitated return to solution. Do not heat the product above room temperature.

Recommended Loading Volumes

Load Volume	Application
7 μ l	Blotting mini (8.3 W x 6.4 L cm) and midi (13.3 W x 8.7 L cm) gels, monitoring transfer
10 μ l	Mini and midi gel electrophoresis
20 μ l	Large (18.3 W x 19.3 L cm) gel electrophoresis

Specifications

Catalog #	161-0374, single unit 161-0394, 5-pack value pack
Volume	500 μ l
Number of applications	50, single unit 250, 5-pack value pack
Loading buffer	30% (w/v) glycerol, 2% SDS, 62.5 mM Tris, pH 6.8, 50 mM DTT, 5 mM EDTA, 0.02% NaN_3
Shelf life	6 months at 4°C or -20°C until expiration date; if the product has been stored as directed for longer than 12 months, the shelf life may be extended by adding additional dithiothreitol (DTT) to 50 mM

Coomassie is a trademark of BASF Aktiengesellschaft.

Tween is a trademark of ICI Americas Inc.

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.

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.