

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 1000 μl
161-0374	Precision Plus Protein Dual Color Standards, 500 μl
161-0394	Precision Plus Protein Dual Color 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

Coomassie is a trademark of BASF Aktiengesellschaft.

Tween is a trademark of ICI Americas Inc.

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.

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

Precision Plus Protein™ All Blue Standards Instruction Manual

Single unit
Catalog #161-0373

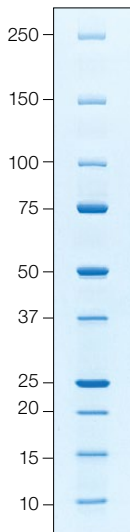
5-Pack value pack
Catalog #161-0393

Product shipped at
room temperature.

Store at 4°C for 6 months
or -20°C for 12 months.

BIO-RAD

Representative lot of
Precision Plus Protein
All Blue prestained
standards on a 4–20%
Tris-HCl Criterion™ 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 (Dual Color and Dual Xtra only).

Note: Prestained standards may fade during blocking, antibody incubation, or washing steps due to the use of Tween 20 in buffers. For other options, contact Bio-Rad Technical Support at 1-800-424-6723.

Product Description

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

Precision Plus Protein All Blue standards represent a breakthrough in protein staining, with molecular weight shifts of <5% and minimal band broadening. This feature, along with three reference bands of 25, 50, and 75 kD, and the broad range of molecular weights, makes location and size approximation of proteins of interest quick and easy.

The molecular weights of the Precision Plus Protein All Blue 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.

Instructions for Use

Precision Plus Protein All Blue standards are provided in loading 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 have returned to solution. Do not heat the product above room temperature.

Recommended Loading Volumes

Load Volume	Application
1–3 µl	Blotting mini (8.3 W x 6.4 L cm) and midi (13.3 W x 8.7 L cm) gels, monitoring transfer
10 µl	Mini and midi gel electrophoresis
20 µl	Large (18.3 W x 19.3 L cm) gel electrophoresis

Specifications

Catalog #	161-0373, single unit 161-0393, 5-pack value pack
Volume	500 µl
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 , 0.01% bromophenol blue
Shelf life	6 months at 4°C or 12 months at –20°C; if the product has been stored as directed for longer than 12 months, the shelf life may be extended by adding dithiothreitol (DTT) to 50 mM