

Products for Electrophoresis

Catalog #	Description
161-0373	Precision Plus Protein Standards, all blue prestained, 500 µl
161-0374	Precision Plus Protein Standards, dual color prestained, 500 µl
161-0737	Laemmli Sample Buffer, 30 ml
161-0786	Bio-Safe™ Coomassie Stain, 1 L

Products for Blot Detection

Catalog #	Description
161-0380	StrepTactin-HRP Conjugate, 0.3 ml, 150 applications
161-0382	StrepTactin-AP Conjugate, 0.3 ml, 150 applications
170-5040	Immun-Star™ HRP Chemiluminescent Substrate
170-5012	Immun-Star AP Substrate Pack

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.

2000 Alfred Nobel Dr., Hercules, CA 94547
(510) 741-1000

4110023 Rev B

Instruction Manual

Precision Plus Protein™ Standards

Unstained

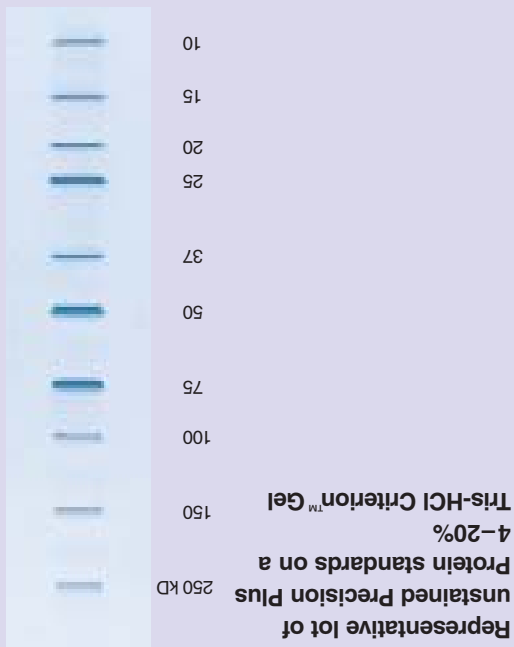
Catalog #161-0363

Product shipped at room temperature.

Store at -20°C upon arrival.

BIO-RAD

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 (catalog #161-0374 only). All Precision Plus Protein standards contain three reference bands (the 25 kD, 50 kD, and 75 kD bands are three times as intense as the other bands). These standards are available in unstained, prestained all blue, and prestained dual color formulations.



Product Description

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

The 25 kD, 50 kD, and 75 kD bands are three times the intensity of the other bands and serve as reference markers.

The molecular weights of the Precision Plus Protein standards, unstained, are confirmed by sequencing, mass spectrometry, and migration in a Laemmli SDS-PAGE system.

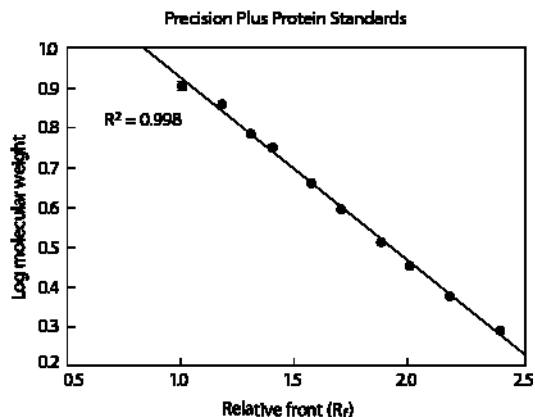
Precision Plus Protein standards, unstained, contain an integral *Strep*-tag sequence, enabling parallel detection on western blots using StrepTactin-HRP (catalog #161-0380) or StrepTactin-AP (catalog #161-0382).

Instructions for Use

Precision Plus Protein standards, unstained, are provided in loading buffer, ready to load, with no dilution required for most applications. Allow the tube to reach room temperature and thoroughly mix before use. This will ensure that any solids that may have precipitated at -20°C have returned to solution. Do not heat the product above room temperature.

Specifications

Catalog #	161-0363
Volume	1 ml
Number of applications	100
Protein quantity	50 kD band, 750 ng; 20 kD and 100 kD bands, 150 ng each*
	* These values should be used only to determine a rough approximation of quantity of a protein of interest.
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	1 year at -20°C



The curve generated by plotting the relative mobility (R_f) vs. the log of the molecular weight can be used to determine the molecular weight of unknown proteins.

Recommended Loading Volumes for Mini Gel Systems

Load Volume	PAGE Application
10 μ l	Coomassie Blue G-250 or R-250 staining
1–3 μ l	Silver staining; SYPRO Ruby staining
1–4 μ l	Blotting: Colorimetric AP development
1–6 μ l	Blotting: Colorimetric HRP development
5–10 μ l	Blotting: Chemiluminescent AP development*
5–10 μ l	Blotting: Chemiluminescent HRP development**

* Dilute Precision Plus Protein standards 1:15–1:30 in Laemmli buffer prior to use.

**Dilute Precision Plus Protein standards 1:30–1:60 in Laemmli buffer prior to use.

Coomassie is a trademark of Imperial Chemical Industries. *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 UK patent 22 72 698. StrepTactin is covered by German patent application P 19641876.3. Bio-Rad is licensed to sell these products for research use only. SYPRO is a trademark of Molecular Probes, Inc.