

## Products for Electrophoresis

Catalog #	Description
161-0363	Precision Plus Protein Unstained Standards, 1,000 µl
161-0373	Precision Plus Protein All Blue Standards, 500 µl
161-0375	Precision Plus Protein Kaleidoscope Standards, 500 µl
161-0737	Laemmli Sample Buffer, 30 ml
161-0786	Bio-Safe™ Coomassie Stain, 1 L

Coomassie is a trademark of BASF Aktiengesellschaft.

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

4110025 Rev C

**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 but the Precision Plus Protein Kaleidoscope standards contain three reference bands (the 25 kD, 50 kD, and 75 kD bands are three times as intense as the other bands).

# Instruction Manual

## Precision Plus Protein™ Standards

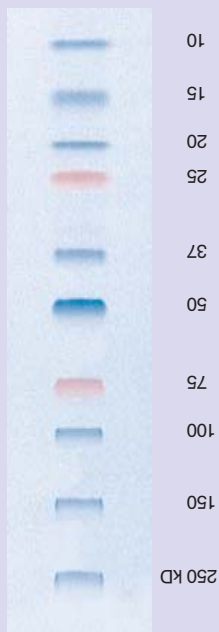
### Dual Color

Catalog #161-0374

Product shipped at room temperature.

Store at –20°C upon arrival.

**BIO-RAD**



**Representative lot of Precision Plus Protein Dual Color standards on a 4–20% Tris-HCl Criterion™ Gel**

## Product Description

Precision Plus Protein Dual Color standards contain ten 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 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 kD and 75 kD) and a more intense blue reference band (50 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 only be used for molecular weight estimation. For molecular weight determination on XT gels, use the Precision Plus Protein Unstained or All Blue standards.

## Instructions for Use

Precision Plus Protein Dual Color 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 at  $-20^{\circ}\text{C}$  have returned to solution. Do not heat the product above room temperature.

## Recommended Loading Volumes

Load Volume	PAGE Application
10 $\mu\text{l}$	Mini gels
5 $\mu\text{l}$	Blotting, to monitor transfer (mini gels)
20 $\mu\text{l}$	Large gels

## Specifications

Catalog #	161-0374
Volume	500 $\mu\text{l}$
Number of applications	50
Loading buffer	30% (w/v) glycerol, 2% SDS, 62.5 mM Tris, pH 6.8, 50 mM DTT, 5 mM EDTA, 0.02% $\text{NaN}_3$
Shelf life	1 year at $-20^{\circ}\text{C}$ *

\*If the product has been stored as directed for one year, its shelf life may be extended by adding dithiothreitol (DTT) to approximately 50 mM.