

New Precision Plus Protein™ standard plugs for all 2-D gels.

# A higher standard

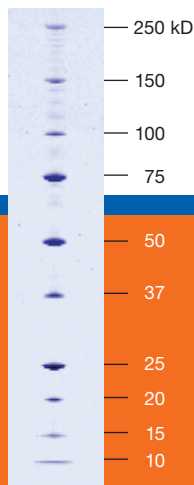
*Actual size*



Precision Plus Protein\* unstained standard plugs allow you to easily, quickly, and cleanly load molecular weight standards on a 2-D gel.

**BIO-RAD**

### Precision Plus Protein standard plug



## The Best Molecular Weight Standards for 2-D Gels

Precision Plus Protein unstained standard plugs allow you to easily, quickly, and cleanly load molecular weight standards on both vertical and horizontal 2-D gels without a reference well.

You no longer have to be creative to load standards when running IPG strips on 2-D gels. Techniques currently used, such as making your own plugs or wicking the standard into the gel, can be things of the past. You can now simply lay the plug next to the IPG strip and begin your run. No fuss, no mess.

### Innovative Precision Plus Protein Standard Plugs Features

- Absolute molecular weight accuracy, confirmed by mass spectrometry
- Ideal 10-band range from 10 kD to 250 kD, with 3 reference bands
- *Strep*-tag amino acid sequence for detection and molecular weight determination on western blots
- Molecular weights that never change between lots
- $r^2 > 0.995$

### Ordering Information

Catalog #	Description
161-0378	Precision Plus Protein Standard Plugs, unstained, 2 x 12
161-1511	PROTEAN® Plus Tris-HCl Gel, 12%, 24 cm IPG well, 25 x 27 cm

*Strep*-tag is a trademark of Institut für Bioanalytik GmbH. *Strep*-tag technology for western blot detection is covered by US patent 5,506,121 and by UK patent 2,272,698. Bio-Rad Laboratories 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.



Break off one plastic plug from the casing.



Use spatula to push plug out of plastic casing onto your gel.



Push plug into IPG well next to IPG strip.

**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 31 3689 6600  
**Canada** 905 364 3435 **China** 86 20 8732 2339 **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** 30 210 777 4396 **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** 0508 805 500  
**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 **United Kingdom** 020 8328 2000