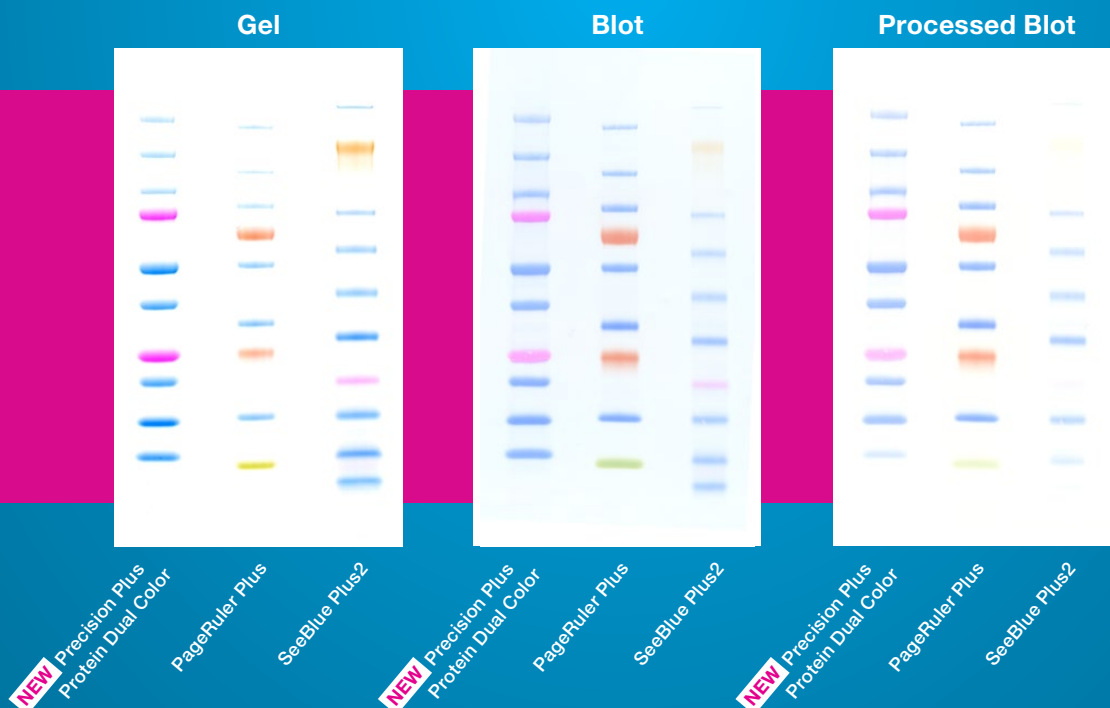




# Brighter, Sharper, Stronger Bands

## Bio-Rad's New and Improved Precision Plus Protein Dual Color Standard

- Brighter for easy identification of target protein
- Sharper for accurate molecular weight estimation
- Stronger band intensity after blot development



# Brighter, Sharper, Stronger Bands



## Ordering Information

Catalog #	Description
<b>Protein Standards</b>	
161-0374	<b>Precision Plus Protein Dual Color Standards</b> , 500 µl, 50 applications
161-0394	<b>Precision Plus Protein Dual Color Standards Value Pack</b> , 2.5 ml (5 x 500 µl), 250 applications

Buffers	
161-0732	<b>10x Tris/Glycine/SDS</b> , 1 L
161-0747	<b>4x Laemmli Sample Buffer</b> , 10 ml

Electrophoresis Cells	
165-6001	<b>Criterion™ Cell</b> , includes electrophoresis buffer tank, lid with power cables, 3 sample loading guides (12+2-well, 18-well, 26-well)
165-8004	<b>Mini-PROTEAN® Tetra Cell for Mini Precast Gels</b> , 4-gel vertical electrophoresis system, includes electrode assembly, companion running module, tank, lid with power cables, mini cell buffer dam

Power Supplies	
164-5050	<b>PowerPac™ Basic Power Supply</b> , 100–120/220–240 V
164-5052	<b>PowerPac HC Power Supply</b> , 100–120/220–240 V
164-5056	<b>PowerPac HV Power Supply</b> , 100–120/220–240 V
164-5070	<b>PowerPac Universal Power Supply</b> , 100–120/220–240 V

Blotting System	
170-4155	<b>Trans-Blot® Turbo™ Transfer Starter System</b> , blotting instrument, includes base, 2 cassettes to hold 1–2 midi or up to 4 mini blotting sandwiches, blot roller, and starter consumable kit




Trans-Blot Turbo Transfer Packs/Kits	
170-4156	<b>Trans-Blot Turbo Transfer Pack, Mini, PVDF</b> , pkg of 10
170-4157	<b>Trans-Blot Turbo Transfer Pack, Midi, PVDF</b> , pkg of 10
170-4158	<b>Trans-Blot Turbo Transfer Pack, Mini, Nitrocellulose</b> , pkg of 10
170-4159	<b>Trans-Blot Turbo Transfer Pack, Midi, Nitrocellulose</b> , pkg of 10
170-4270	<b>Trans-Blot Turbo RTA Transfer Kit, Mini, Nitrocellulose</b> , for 40 blots
170-4271	<b>Trans-Blot Turbo RTA Transfer Kit, Midi, Nitrocellulose</b> , for 40 blots
170-4272	<b>Trans-Blot Turbo RTA Transfer Kit, Mini, PVDF</b> , for 40 blots
170-4273	<b>Trans-Blot Turbo RTA Transfer Kit, Midi, PVDF</b> , for 40 blots
170-4274	<b>Trans-Blot Turbo RTA Transfer Kit, Mini, LF PVDF</b> , for 40 blots
170-4275	<b>Trans-Blot Turbo RTA Transfer Kit, Midi, LF PVDF</b> , for 40 blots

Substrates	
170-5060	<b>Clarity™ Western ECL Substrate</b> , 200 ml, contains 100 ml Clarity Western Peroxide Reagent and 100 ml Clarity Western Luminol/Enhancer Reagent
170-5061	<b>Clarity Western ECL Substrate</b> , 500 ml, contains 250 ml Clarity Western Peroxide Reagent and 250 ml Clarity Western Luminol/Enhancer Reagent

Description			
	10-Well 30 µl	10-Well 50 µl	12-Well 20 µl
<b>Mini-PROTEAN® TGX Stain-Free™ Precast Gels*</b>			
10%	456-8033	456-8034	456-8035
4–15%	456-8083	456-8084	456-8045
4–20%	456-8093	456-8094	456-8085
Any kD™	456-8123	456-8124	456-8125

All formats are available as both ten packs (catalog numbers listed) and two packs. To order as a two pack, add an "S" to the end of the catalog number for the corresponding ten pack.

\* Mini-PROTEAN TGX Stain-Free Gels are available in additional gel percentages.

Description			
	12+2-Well 45 µl	18-Well 30 µl	26-Well 15 µl
<b>Criterion™ TGX Stain-Free™ Precast Gels**</b>			
10%	567-8033	567-8034	567-8035
4–15%	567-8083	567-8084	567-8085
4–20%	567-8093	567-8094	567-8095
Any kD	567-8123	567-8124	567-8125

\* Reference well accommodates 15 µl of markers/standards.

\*\* Criterion TGX Stain-Free Gels are sold as a single gel and available in additional gel percentages.

Catalog #	Description
-----------	-------------

### Gel Casting Solutions

161-0170	<b>TGX™ FastCast™ Acrylamide Starter Kit</b> , 7.5%
161-0171	<b>TGX FastCast Acrylamide Kit</b> , 7.5%
161-0172	<b>TGX FastCast Acrylamide Starter Kit</b> , 10%
161-0173	<b>TGX FastCast Acrylamide Kit</b> , 10%
161-0174	<b>TGX FastCast Acrylamide Starter Kit</b> , 12%
161-0175	<b>TGX FastCast Acrylamide Kit</b> , 12%
161-0180	<b>TGX Stain-Free™ FastCast™ Acrylamide Starter Kit</b> , 7.5%
161-1081	<b>TGX Stain-Free FastCast Acrylamide Kit</b> , 7.5%
161-1082	<b>TGX Stain-Free FastCast Acrylamide Starter Kit</b> , 10%
161-1083	<b>TGX Stain-Free FastCast Acrylamide Kit</b> , 10%
161-1084	<b>TGX Stain-Free FastCast Acrylamide Starter Kit</b> , 12%
161-1085	<b>TGX Stain-Free FastCast Acrylamide Kit</b> , 12%

PageRuler is a trademark of Thermo Fisher Scientific Inc.

SeeBlue is a trademark of Life Technologies Corporation.

Precision Plus Protein Standards are 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  
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 11 3065 7550  
Canada 905 364 3435 China 86 21 6169 8500 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 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666  
New Zealand 64 9 415 2280 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 026 674 55 05  
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000