



ELECTROPHORESIS

It's Time to Expect More from Handcast Gels!

Introducing Bio-Rad's TGX™ and TGX Stain-Free™ FastCast™ Acrylamide Kits



- Cast gels and use as needed for up to a month
- Gels are ready to run in 30 minutes
- Enables more efficient transfer than Tris-HCl gels

With the TGX Stain-Free kits, you can also:

- Monitor success of your electrophoresis and transfer
- Visualize your proteins in less than 5 minutes after your gel run
- Validate your data by total protein normalization and analysis

How it works:

1. Quickly mix solutions 1:1
2. Add APS and TEMED
3. Pour stacker immediately after pouring resolver
4. 30 minute polymerization time
5. 20 minute run time
6. More efficient transfers in just 3–7 minutes

Starter and full kit sizes now available in 7.5%, 10%, and 12% formulations.

Visit www.bio-rad.com or contact your sales representative for more information.

Having a four-week shelf life is extremely useful.

— FastCast Customer
Chris

Using the stain-free kit, we can activate gels and look at the loading efficiency while the gel is being transferred. Once the gel is transferred to the membrane, we can look at the transfer efficiency and conduct total protein normalization using the same gel.

— FastCast Customer
Isabelle

With the new FastCast system I can breeze through what typically is the tedious task of running a gel and move on to the next step much more quickly.

— FastCast Customer
Aaron

BIO-RAD

Approximate number of gels per kit.

	Mini Gels (~9 x 7 cm)				Midi Gels (~13 x 9 cm)
	Mini-PROTEAN® Cassettes	0.75 mm Glass Plates	1.0 mm Glass Plates	1.5 mm Glass Plates	Criterion™ Cassettes
Starter Kit	11 gels	19 gels	13 gels	9 gels	6 gels
Kit	55 gels	95 gels	65 gels	45 gels	30 gels

Ordering Information

Catalog #	Description
161-0170	TGX FastCast Acrylamide Starter Kit, 7.5%
161-0171	TGX FastCast Acrylamide Kit, 7.5%
161-0172	TGX FastCast Acrylamide Starter Kit, 10%
161-0173	TGX FastCast Acrylamide Kit, 10%
161-0174	TGX FastCast Acrylamide Starter Kit, 12%
161-0175	TGX FastCast Acrylamide Kit, 12%
161-0180	TGX Stain-Free FastCast Acrylamide Starter Kit, 7.5%
161-0181	TGX Stain-Free FastCast Acrylamide Kit, 7.5%
161-0182	TGX Stain-Free FastCast Acrylamide Starter Kit, 10%
161-0183	TGX Stain-Free FastCast Acrylamide Kit, 10%
161-0184	TGX Stain-Free FastCast Acrylamide Starter Kit, 12%
161-0185	TGX Stain-Free FastCast Acrylamide Kit, 12%
161-0800	TEMED, 5 ml
161-0700	Ammonium Persulfate (APS), 10 g

Kits Including TEMED and APS:

161-0171TA	TGX FastCast Acrylamide Kit, 7.5% with TEMED, APS
161-0173TA	TGX FastCast Acrylamide Kit, 10% with TEMED, APS
161-0175TA	TGX FastCast Acrylamide Kit, 12% with TEMED, APS
161-0181TA	TGX Stain-Free FastCast Acrylamide Kit, 7.5% with TEMED
161-0183TA	TGX Stain-Free FastCast Acrylamide Kit, 10% with TEMED
161-0185TA	TGX Stain-Free FastCast Acrylamide Kit, 12% with TEMED

Protein Standards

161-0396	Precision Plus Protein™ Unstained Standards Value Pack, 10–250 kD, 5 x 1,000 µl, 500 applications
161-0393	Precision Plus Protein All Blue Standards Value Pack, 10–250 kD, 5 x 500 µl, 250 applications
161-0394	Precision Plus Protein Dual Color Standards Value Pack, 10–250 kD, 5 x 500 µl, 250 applications
161-0395	Precision Plus Protein™ Kaleidoscope™ Standards Value Pack, 10–250 kD, 5 x 500 µl, 250 applications
161-0397	Precision Plus Protein Dual Xtra Value Pack, 2–250 kD, 5 x 500 µl, 250 applications
161-0398	Precision Plus Protein™ WesternC™ Pack, 10–250 kD, 5 x 250 µl WesternC standard, 5 x 250 µl HRP conjugate, 250 applications
161-0378	Precision Plus Protein Standard Plugs, unstained, 10–250 kD, 24 applications

Buffer

161-0747	4x Laemmli Sample Buffer, 10 ml
161-0737	2x Laemmli Sample Buffer, 30 ml
161-0732	10x Tris/Glycine/SDS, 1 L
161-0734	10x Tris/Glycine, 1 L

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 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