



ELECTROPHORESIS

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

Tips & Tricks

Tip #1

Pour your desired number of gels by taking out the appropriate volumes (in a 1:1 ratio) from each bottle.

Tip #2

Use Tris-glycine buffers when running the FastCast gels. The buffers are optimized to complement the gel chemistry.

Tip #3

It helps to insert the comb at a slight diagonal before pressing down when casting the gel. This helps to prevent air bubbles in the gel.

Tip #4

Use water instead of ethanol to test glass plates for leaks. Ethanol may cause erosion and cracking of the glass plates over time, which would result in leaks.

Tip #5

Kits can be stored between 4°C and 28°C, but be sure to bring kits back to room temperature before use.

Tip #6

Carefully position the pipet to be angled 20 degrees to the lab bench as you distribute the solution into the glass plates or cassette. Hold the pipet steady and aim for the center of the glass plates or cassette to ensure even distribution of the solution.

Tip #7

Pour the stacker immediately after the resolver with TGX FastCast acrylamide kits. Using an overlay to make sure the top of the resolver is straight is not necessary.

Tip #8

If using glass plates, be sure to check that they are not cracked or chipped. This will help avoid leaks.

Tip #9

Do not run different gel types (chemistry) or percentages in the same tank at the same time.

Tip #10

Using too much APS or TEMED may cause inconsistent gel polymerization and could compromise your results.

Tip #11

If using glass plates, be sure to replace the casting stand gaskets regularly. This will help to prevent leaks.

Tip #12

Mark the cassette (or glass plates) to ensure the resolving solution is poured to the desired distance below the comb.

Tip #13

Make fresh 10% APS solution on the same day that gels are cast for optimal polymerization.

Tip #14

Cover gels with a paper towel wetted with deionized water and store horizontally in a zipper sealed plastic bag at 4°C for up to 1 month.

Tip #15

Check polymerization by pouring extra resolving and/or stacking solution into a Falcon tube.

For more product information, go to www.bio-rad.com/go/fastcast.

BIO-RAD

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.

Falcon is a trademark of Becton, Dickinson and Company.



BIO-RAD

**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 026674 55 05
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000