



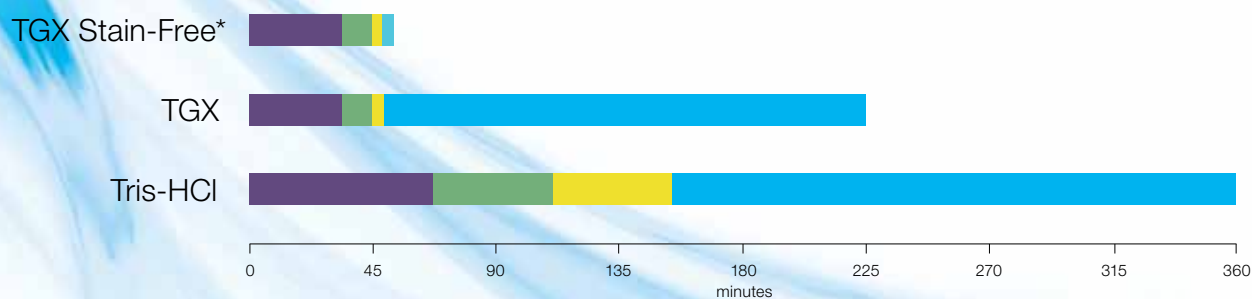
TGX™ and TGX Stain-Free™ FastCast™ Acrylamide Kits

The newest frontier for handcast gels



It's Time to Expect More from Handcast Gels!

FastCast kits provide unparalleled convenience and performance.



- Solution preparation and casting time (includes polymerization)
- Gel run time
- Transfer time (3-7 min transfers only with TGX or TGX Stain-Free gels)
- Visualizing/staining

*** Added benefits of TGX Stain-Free gels:**

- No stain needed to visualize protein bands
- ≤5 minute visualization and imaging

Advance Your Research with FastCast™ Acrylamide Solutions

FastCast acrylamide solutions bring the latest chemistry in gel electrophoresis to those who hand cast gels—providing increased convenience, speed, and performance over traditional handcast gels.

Long shelf life

Stain-free gel visualization

Confidence

Performance

Fast gel runs

Fast transfers

Easy casting

Efficient blotting

Convenience

Short casting time



Bring Convenience and Ease of Use to Your Lab



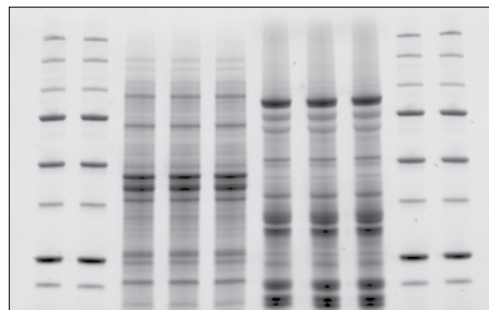
FastCast gels are ready to run in 30 minutes

- Cast stacker immediately after resolver with no overlay necessary
- Quickly mix 1:1, add TEMED and APS, and pour gels
- Use cost-effective Tris-glycine buffers
- Kits may be stored at room temperature for up to 1 year. Open and cast when needed

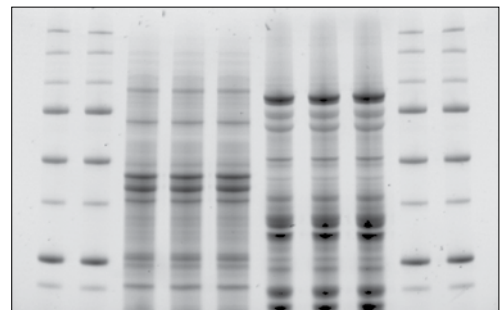


Cast gels once a month and use as experiments call for them

- Focus on your experimental plan, not your casting plan
- Cast less often with gels that maintain freshness and performance for 1 month
- Store at 4°C and use gels when needed



Freshly cast 10% TGX Stain-Free gel



After 1 month of storage at 4°C

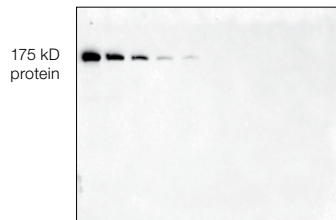
Exceed Your Performance Expectations



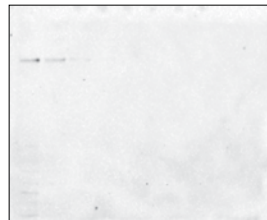
Get superior results faster

- Complete gel runs in 20 minutes
- Transfer proteins efficiently in 3 to 7 minutes
- Efficient transfer across high and low molecular weights

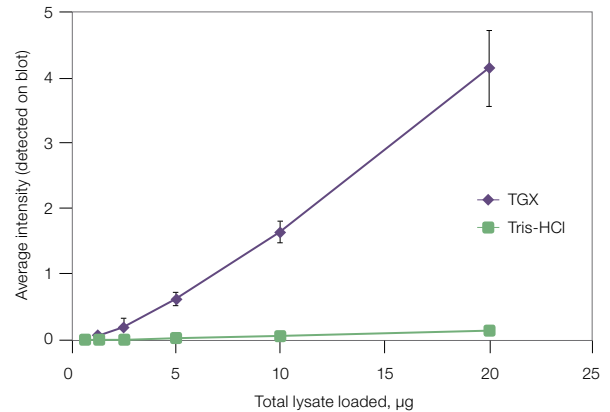
Blotting Efficiency Comparison: TGX gels enable more complete transfer



Final blot of 10% TGX Stain-Free gel demonstrating excellent transfer efficiency of a high molecular weight protein

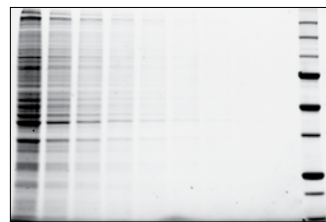


Final blot of 10% Tris-HCl gel demonstrating reduced transfer efficiency versus the TGX Stain-Free gel



Increase confidence in your results

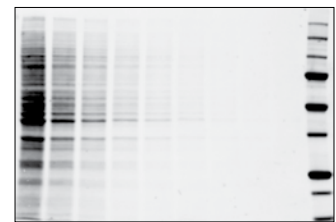
- With TGX Stain-Free FastCast kits, see your results in less than 5 minutes after gel run
- No more fixing, staining, and destaining necessary
- Monitor success of electrophoresis and transfers through stain-free visualization of gels and membranes



Stain-free image upon completion of gel electrophoresis confirming separation completed successfully



Stain-free image of gel post-transfer showing that the majority of proteins transferred from the gel to the membrane



Stain-free image of membrane post-transfer demonstrating successful transfer of proteins to the membrane

Ordering Information

Catalog # Description

Gel Casting Solutions and Reagents

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**

Electrophoresis Cells

165-8000 **Mini-PROTEAN® Tetra Cell, 10-well, 0.75 mm thickness**
 165-8001 **Mini-PROTEAN Tetra Cell, 10-well, 1.0 mm thickness**
 165-8006 **Mini-PROTEAN Tetra Cell, 10-well, 1.5 mm thickness**
 165-4100 **Mini-PROTEAN 3 Dodeca™ cell**
 165-4130 **Criterion™ Dodeca™ cell**
 165-4139 **Criterion Dodeca cell and PowerPac™ Universal Power Supply**

Mini-PROTEAN and Criterion Empty Cassettes and Combs

456-0003 **Mini-PROTEAN Empty Cassettes, 1.0 mm thick, 10-well, pkg of 50**
 456-0005 **Mini-PROTEAN Empty Cassettes, 1.0 mm thick, 12-well, pkg of 50**
 456-0006 **Mini-PROTEAN Empty Cassettes, 1.0 mm thick, 15-well, pkg of 50**
 456-0001 **Mini-PROTEAN Empty Cassettes, 1.0 mm thick, IPG well, pkg of 50**
 456-0013 **Mini-PROTEAN Combs, 10-well, pkg of 50, for use with 456-0003**
 456-0015 **Mini-PROTEAN Combs, 12-well, pkg of 50, for use with 456-0005**
 456-0016 **Mini-PROTEAN Combs, 15-well, pkg of 50, for use with 456-0006**
 456-0011 **Mini-PROTEAN Combs, IPG well, pkg of 50, for use with 456-0001**
 345-9901 **Criterion Empty Cassettes, 1.0 mm, 12+2-well, pkg of 10**
 345-9902 **Criterion Empty Cassettes, 1.0 mm, 18-well, pkg of 10**
 345-9903 **Criterion Empty Cassettes, 1.0 mm, 26-well, pkg of 10**
 345-9904 **Criterion Empty Cassettes, 1.0 mm, prep+2-well, pkg of 10**
 345-9906 **Criterion Empty Cassettes, 1.0 mm, IPG+1-well, pkg of 10**

Catalog # Description

AnyGel™ Stands

165-4131 **AnyGel Stand, Pkg of 1, single-row gel stand, holds 1 PROTEAN® gel, 2 Criterion gels, or 3 Mini-PROTEAN or Ready Gel® precast mini gels**
 165-5131 **AnyGel Stand, Pkg of 1, 6-row gel stand, holds 1 PROTEAN gel, 12 Criterion gels, or 18 Mini-PROTEAN or Ready Gel precast mini gels**

Detection Reagent

170-5060 **Clarity™ Western ECL Substrate, 200 ml**

Blotting Systems

170-3935 **Mini Trans-Blot® Module, without lower buffer tank and lid, for use with the Mini-PROTEAN Tetra cell**
 170-4155 **Trans-Blot® Turbo™ Transfer Starter System, blotting instrument, includes base, 2 cassettes to hold 1–2 midi or up to 4 mini blotting sandwiches, blot roller, and starter consumable kit**
 170-4270 **Trans-Blot Turbo RTA Transfer Kit, ready-to-assemble, mini, nitrocellulose, pkg of 40**
 170-4271 **Trans-Blot Turbo RTA Transfer Kit, ready-to-assemble, midi, nitrocellulose, pkg of 40**
 170-4272 **Trans-Blot Turbo RTA Transfer Kit, ready-to-assemble, mini, PVDF, pkg of 40**
 170-4273 **Trans-Blot Turbo RTA Transfer Kit, ready-to-assemble, midi, PVDF, pkg of 40**
 170-4274 **Trans-Blot Turbo RTA Transfer Kit, ready-to-assemble, mini, LF PVDF, pkg of 40**
 170-4275 **Trans-Blot Turbo RTA Transfer Kit, ready-to-assemble, midi, LF PVDF, pkg of 40**

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**

Buffers

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**

Power Supplies

164-5050 **PowerPac Basic Power Supply**
 164-5052 **PowerPac HC Power Supply**
 164-5056 **PowerPac HV Power Supply**
 164-5070 **PowerPac Universal Power Supply**

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

**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