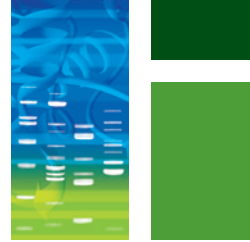


ELECTROPHORESIS

# Thermo Fisher Scientific (Invitrogen) Midi Precast Gels to Bio-Rad Criterion TGX and Criterion TGX Stain-Free Precast Gels

## Catalog Number Conversion Chart



Thermo Fisher Scientific (Invitrogen) Bis-Tris and Tris-Glycine Precast Gels, Midi Size (Boxes of 10)		
	Description	Catalog #
8% Bis-Tris (MOPS buffer)	NuPAGE 8% Bis-Tris Midi Gel, 12+2-well	WG1001BOX
	NuPAGE 8% Bis-Tris Midi Gel, 20-well	WG1002BOX
	NuPAGE 8% Bis-Tris Midi Gel, 26-well	WG1003BOX
10% Bis-Tris (MOPS buffer)	NuPAGE 10% Bis-Tris Midi Gel, 12+2-well	WG1201BOX
	NuPAGE 10% Bis-Tris Midi Gel, 20-well	WG1202BOX
	NuPAGE 10% Bis-Tris Midi Gel, 26-well	WG1203BOX
4-12% Bis-Tris (MOPS buffer)	NuPAGE 4-12% Bis-Tris Midi Gel, 12+2-well	WG1401BOX
	NuPAGE 4-12% Bis-Tris Midi Gel, 20-well	WG1402BOX
	NuPAGE 4-12% Bis-Tris Midi Gel, 26-well	WG1403BOX
10% Tris- Glycine	Novex 10% Tris-Glycine Midi Gel, 12+2-well	WXP01012BOX
	Novex 10% Tris-Glycine Midi Gel, 20-well	WXP01020BOX
	Novex 10% Tris-Glycine Midi Gel, 26-well	WXP01026BOX
12% Tris- Glycine	Novex 12% Tris-Glycine Midi Gel, 12+2-well	WXP01212BOX
	Novex 12% Tris-Glycine Midi Gel, 20-well	WXP01220BOX
	Novex 12% Tris-Glycine Midi Gel, 26-well	WXP01226BOX
4-12% Tris- Glycine	Novex 4-12% Tris-Glycine Midi Gel, 12+2-well	WXP41212BOX
	Novex 4-12% Tris-Glycine Midi Gel, 20-well	WXP41220BOX
	Novex 4-12% Tris-Glycine Midi Gel, 26-well	WXP41226BOX
4-20% Tris- Glycine	Novex 4-20% Tris-Glycine Midi Gel, 12+2-well	WXP42012BOX
	Novex 4-20% Tris-Glycine Midi Gel, 20-well	WXP42020BOX
	Novex 4-20% Tris-Glycine Midi Gel, 26-well	WXP42026BOX
	Novex 4-20% Tris-Glycine Midi Gel, 12+2-well	WXP42012BOX
	Novex 4-20% Tris-Glycine Midi Gel, 26-well	WXP42026BOX
8-16% Tris- Glycine	Novex 8-16% Tris-Glycine Midi Gel, 12+2-well	WXP81612BOX
	Novex 8-16% Tris-Glycine Midi Gel, 20-well	WXP81620BOX
	Novex 8-16% Tris-Glycine Midi Gel, 26-well	WXP81626BOX

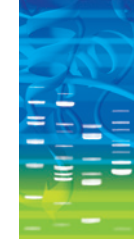
Bio-Rad Laboratories, Inc.				
Criterion TGX Precast Gels*		Criterion TGX Stain-Free Precast Gels*		
	Catalog #	Description	Catalog #	Description
7.5%	5671023	Criterion TGX, 7.5%, 12+2-well	5678023	Criterion TGX Stain-Free, 7.5%, 12+2-well
	5671024	Criterion TGX, 7.5%, 18-well	5678024	Criterion TGX Stain-Free, 7.5%, 18-well
	5671025	Criterion TGX, 7.5%, 26-well	5678025	Criterion TGX Stain-Free, 7.5%, 26-well
10%	5671033	Criterion TGX, 10%, 12+2-well	5678033	Criterion TGX Stain-Free, 10%, 12+2-well
	5671034	Criterion TGX, 10%, 18-well	5678034	Criterion TGX Stain-Free, 10%, 18-well
	5671035	Criterion TGX, 10%, 26-well	5678035	Criterion TGX Stain-Free, 10%, 26-well
4-15%	5671083	Criterion TGX, 4-15%, 12+2-well	5678083	Criterion TGX Stain-Free, 4-15%, 12+2-well
	5671084	Criterion TGX, 4-15%, 18-well	5678084	Criterion TGX Stain-Free, 4-15%, 18-well
	5671085	Criterion TGX, 4-15%, 26-well	5678085	Criterion TGX Stain-Free, 4-15%, 26-well
10%	5671033	Criterion TGX, 10%, 12+2-well	5678033	Criterion TGX Stain-Free, 10%, 12+2-well
	5671034	Criterion TGX, 10%, 18-well	5678034	Criterion TGX Stain-Free, 10%, 18-well
	5671035	Criterion TGX, 10%, 26-well	5678035	Criterion TGX Stain-Free, 10%, 26-well
12%	5671043	Criterion TGX, 12%, 12+2-well	5678043	Criterion TGX Stain-Free, 12%, 12+2-well
	5671044	Criterion TGX, 12%, 18-well	5678044	Criterion TGX Stain-Free, 12%, 18-well
	5671045	Criterion TGX, 12%, 26-well	5678045	Criterion TGX Stain-Free, 12%, 26-well
4-15%	5671083	Criterion TGX, 4-15%, 12+2-well	5678083	Criterion TGX Stain-Free, 4-15%, 12+2-well
	5671084	Criterion TGX, 4-15%, 18-well	5678084	Criterion TGX Stain-Free, 4-15%, 18-well
	5671085	Criterion TGX, 4-15%, 26-well	5678085	Criterion TGX Stain-Free, 4-15%, 26-well
4-20%	5671093	Criterion TGX, 4-20%, 12+2-well	5678093	Criterion TGX Stain-Free, 4-20%, 12+2-well
	5671094	Criterion TGX, 4-20%, 18-well	5678094	Criterion TGX Stain-Free, 4-20%, 18-well
	5671095	Criterion TGX, 4-20%, 26-well	5678095	Criterion TGX Stain-Free, 4-20%, 26-well
	5671123	Criterion TGX, Any kD, 12+2-well	5678123	Criterion TGX Any kD Stain-Free, 12+2-well
	5671124	Criterion TGX, Any kD, 18-well	5678124	Criterion TGX Any kD Stain-Free, 18-well
8-16%	5671125	Criterion TGX, Any kD, 26-well	5678125	Criterion TGX Any kD Stain-Free, 26-well
	5671103	Criterion TGX, 8-16%, 12+2-well	5678103	Criterion TGX Stain-Free, 8-16%, 12+2-well
	5671104	Criterion TGX, 8-16%, 18-well	5678104	Criterion TGX Stain-Free, 8-16%, 18-well
	5671105	Criterion TGX, 8-16%, 26-well	5678105	Criterion TGX Stain-Free, 8-16%, 26-well

\* Criterion TGX and TGX Stain-Free Gels are also available in IPG/prep+1-well and Prep+2-well sizes but are not listed here as there are no equivalent Thermo Fisher Scientific midi precast gels available.

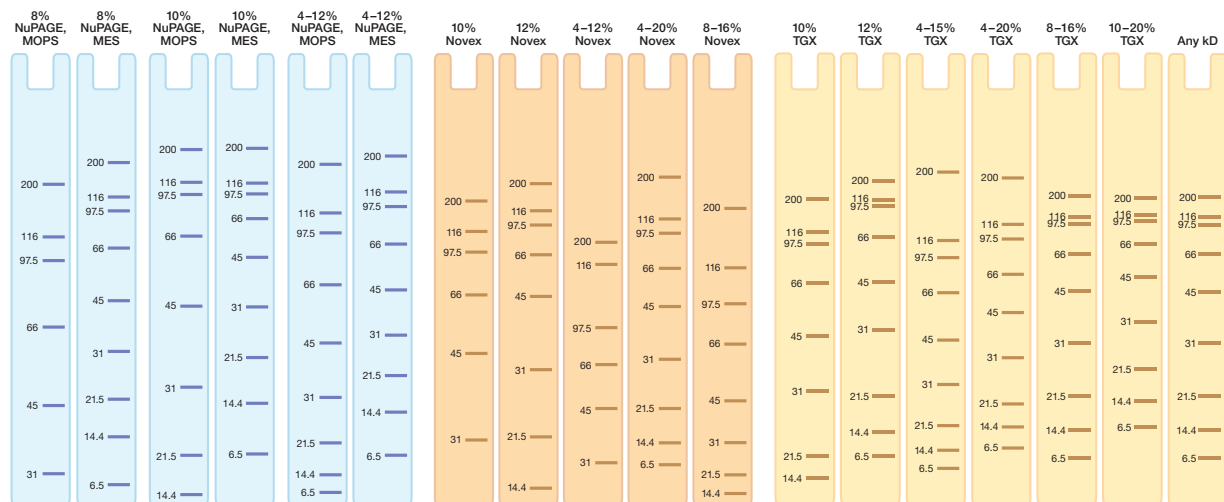


# Criterion TGX and TGX Stain-Free Precast Gels

## Accelerated Electrophoresis with Faster Run Times



### Migration Charts Comparing Criterion TGX and TGX Stain-Free with NuPAGE and Novex Gels



Side-by-side comparison of migration patterns of broad range SDS-PAGE unstained standards run on Thermo Fisher Scientific (Invitrogen) midi precast gels: NuPAGE Bis-Tris Gels (■), Novex Tris-Glycine Gels (■), Criterion TGX and TGX Stain-Free Precast Gels (■). The 6.5 kD protein standard band is visible only on the Criterion TGX and not on the TGX Stain-Free Precast Gels because it lacks tryptophan residues that are required for Stain-Free detection. Migration patterns are shown for NuPAGE Bis-Tris Gels run using both MOPS and MES buffers.



### Criterion TGX and TGX Stain-Free Precast Gels deliver the fastest separation and extended shelf life using the Laemmli PAGE system

- Run times as short as 15 min
- Transfer in as little as 8 min
- Inexpensive Laemmli buffer system, low running costs
- One-year shelf life

### Criterion TGX Stain-Free Precast Gels deliver

- Built-in checkpoints throughout electrophoresis and western blotting workflow
- Complete protein separation, gel imaging, and data analysis in 25 min
- Comparable sensitivity to Coomassie stain
- Better reproducibility and quantitation compared to staining procedures
- Use of the same gel for downstream western blotting, standard staining methods, and mass spectrometry analysis

### Related Literature

- 3668 ChemiDoc™ Go Imaging System
- 4110001 Criterion Precast Gels Instruction Manual and Application Guide
- 5934 NuPAGE Bis-Tris Precast Gels to Mini-PROTEAN™ TGX and Mini-PROTEAN TGX Stain-Free Precast Gels Catalog Number Conversion Chart
- 5974 Mini-PROTEAN and Criterion TGX Stain-Free Precast Gels
- 5981 Criterion Tris-HCl to Criterion TGX and Criterion TGX Stain-Free Precast Gels Conversion Chart
- 6007 Criterion TGX Precast Gels Product Information Sheet
- 6040 A Guide to Polyacrylamide Gel Electrophoresis and Detection
- 6873 ChemiDoc and ChemiDoc MP Imaging System Brochure
- 7284 GelDoc™ Go Imaging System Brochure

Visit [bio-rad.com/CriterionGel](http://bio-rad.com/CriterionGel) for more information.

Still unsure what gels would work best for your experiment? Try the [Free Western Blotting Selector Tool](#).

BIO-RAD, CHEMIDOC, GELDOC, and MINI-PROTEAN are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2024 Bio-Rad Laboratories, Inc.

TGX Stain-Free Precast Gels are covered by U.S. Patent Numbers 7,569,130 and 8,007,646.