

ELECTROPHORESIS

Life Technologies Midi Precast Gels to Bio-Rad Criterion™ TGX™ and Criterion™ TGX Stain-Free™ Precast Gels  
Catalog Number Conversion Chart



Life Technologies Bis-Tris and Tris-Glycine Precast Gels Midi Size (Box of 10)		
	Description	Catalog #
8% Bis-Tris (MOPS buffer)	NuPAGE Novex 8% Bis-Tris Midi Gel 12+2-well	WG1001BOX →
	NuPAGE Novex 8% Bis-Tris Midi Gel 20-well	WG1002BOX →
	NuPAGE Novex 8% Bis-Tris Midi Gel 26-well	WG1003BOX →
10% Bis-Tris (MOPS buffer)	NuPAGE Novex 10% Bis-Tris Midi Gel 12+2-well	WG1201BOX →
	NuPAGE Novex 10% Bis-Tris Midi Gel 20-well	WG1202BOX →
	NuPAGE Novex 10% Bis-Tris Midi Gel 26-well	WG1203BOX →
4-12% Bis-Tris (MOPS buffer)	NuPAGE Novex 4-12% Bis-Tris Midi Gel, 12+2-well	WG1401BOX →
	NuPAGE Novex 4-12% Bis-Tris Midi Gel, 20-well	WG1402BOX →
	NuPAGE Novex 4-12% Bis-Tris Midi Gel, 26-well	WG1403BOX →
8% Tris-Glycine	Novex 8% Tris-Glycine Midi Gel 12+2-well	WT0081BOX →
	Novex 8% Tris-Glycine Midi Gel 20-well	WT0082BOX →
	Novex 8% Tris-Glycine Midi Gel 26-well	WT0083BOX →
10% Tris-Glycine	Novex 10% Tris-Glycine Midi Gel 12+2-well	WT0101BOX →
	Novex 10% Tris-Glycine Midi Gel 20-well	WT0102BOX →
	Novex 10% Tris-Glycine Midi Gel 26-well	WT0103BOX →
12% Tris-Glycine	Novex 12% Tris-Glycine Midi Gel 12+2-well	WT0121BOX →
	Novex 12% Tris-Glycine Midi Gel 20-well	WT0122BOX →
	Novex 12% Tris-Glycine Midi Gel 26-well	WT0123BOX →
4-12% Tris-Glycine	Novex 4-12% Tris-Glycine Midi Gel, 12+2-well	WT4121BOX →
	Novex 4-12% Tris-Glycine Midi Gel, 20-well	WT4122BOX →
	Novex 4-12% Tris-Glycine Midi Gel, 26-well	WT4123BOX →
4-20% Tris-Glycine	Novex 4-20% Tris-Glycine Midi Gel, 12+2-well	WT4201BOX →
	Novex 4-20% Tris-Glycine Midi Gel, 20-well	WT4202BOX →
	Novex 4-20% Tris-Glycine Midi Gel, 26-well	WT4203BOX →
	Novex 4-20% Tris-Glycine Midi Gel, 12+2-well	WT4201BOX →
	Novex 4-20% Tris-Glycine Midi Gel, 20-well	WT4202BOX →
8-16% Tris-Glycine	Novex 8-16% Tris-Glycine Midi Gel 12+2-well	WT8161BOX →
	Novex 8-16% Tris-Glycine Midi Gel 20-well	WT8162BOX →
	Novex 8-16% Tris-Glycine Midi Gel 26-well	WT8163BOX →

Criterion TGX Precast Gels*		Criterion TGX Stain-Free Precast Gels*		
	Catalog #	Description	Catalog #	Description
7.5%	567-1023	Criterion TGX 7.5%, 12+2-well	567-8023	Criterion TGX Stain-Free, 7.5%, 12+2-well
	567-1024	Criterion TGX 7.5%, 18-well	567-8024	Criterion TGX Stain-Free, 7.5%, 18-well
	567-1025	Criterion TGX 7.5%, 26-well	567-8025	Criterion TGX Stain-Free, 7.5%, 26-well
10%	567-1033	Criterion TGX 10%, 12+2-well	567-8033	Criterion TGX Stain-Free, 10%, 12+2-well
	567-1034	Criterion TGX 10%, 18-well	567-8034	Criterion TGX Stain-Free, 10%, 18-well
	567-1035	Criterion TGX 10%, 26-well	567-8035	Criterion TGX Stain-Free, 10%, 26-well
4-15%	567-1083	Criterion TGX, 4-15%, 12+2-well	567-8083	Criterion TGX Stain-Free, 4-15%, 12+2-well
	567-1084	Criterion TGX, 4-15%, 18-well	567-8084	Criterion TGX Stain-Free, 4-15%, 18-well
	567-1085	Criterion TGX, 4-15%, 26-well	567-8085	Criterion TGX Stain-Free, 4-15%, 26-well
7.5%	567-1023	Criterion TGX, 7.5%, 12+2-well	567-8023	Criterion TGX Stain-Free, 7.5%, 12+2-well
	567-1024	Criterion TGX, 7.5%, 18-well	567-8024	Criterion TGX Stain-Free, 7.5%, 18-well
	567-1025	Criterion TGX, 7.5%, 26-well	567-8025	Criterion TGX Stain-Free, 7.5%, 26-well
10%	567-1033	Criterion TGX, 10%, 12+2-well	567-8033	Criterion TGX Stain-Free, 10%, 12+2-well
	567-1034	Criterion TGX, 10%, 18-well	567-8034	Criterion TGX Stain-Free, 10%, 18-well
	567-1035	Criterion TGX, 10%, 26-well	567-8035	Criterion TGX Stain-Free, 10%, 26-well
12%	567-1043	Criterion TGX, 12%, 12+2-well	567-8043	Criterion TGX Stain-Free, 12%, 12+2-well
	567-1044	Criterion TGX, 12%, 18-well	567-8044	Criterion TGX Stain-Free, 12%, 18-well
	567-1045	Criterion TGX, 12%, 26-well	567-8045	Criterion TGX Stain-Free, 12%, 26-well
4-15%	567-1083	Criterion TGX, 4-15%, 12+2-well	567-8083	Criterion TGX Stain-Free, 4-15%, 12+2-well
	567-1084	Criterion TGX, 4-15%, 18-well	567-8084	Criterion TGX Stain-Free, 4-15%, 18-well
	567-1085	Criterion TGX, 4-15%, 26-well	567-8085	Criterion TGX Stain-Free, 4-15%, 26-well
4-20%	567-1093	Criterion TGX, 4-20%, 12+2-well	567-8093	Criterion TGX Stain-Free, 4-20%, 12+2-well
	567-1094	Criterion TGX, 4-20%, 18-well	567-8094	Criterion TGX Stain-Free, 4-20%, 18-well
	567-1095	Criterion TGX, 4-20%, 26-well	567-8095	Criterion TGX Stain-Free, 4-20%, 26-well
Any kD™	567-1123	Criterion TGX Any kD™, 12+2-well	567-8123	Criterion™ TGX Any kD Stain-Free™, 12+2-well
	567-1124	Criterion TGX Any kD, 18-well	567-8124	Criterion TGX Any kD Stain-Free, 18-well
	567-1125	Criterion TGX Any kD, 26-well	567-8125	Criterion TGX Any kD Stain-Free, 26-well
8-16%	567-1103	Criterion TGX, 8-16% 12+2-well	567-8103	Criterion TGX Stain-Free, 8-16% 12+2-well
	567-1104	Criterion TGX, 8-16% 18-well	567-8104	Criterion TGX Stain-Free, 8-16% 18-well
	567-1105	Criterion TGX, 8-16% 26-well	567-8105	Criterion TGX Stain-Free, 8-16% 26-well

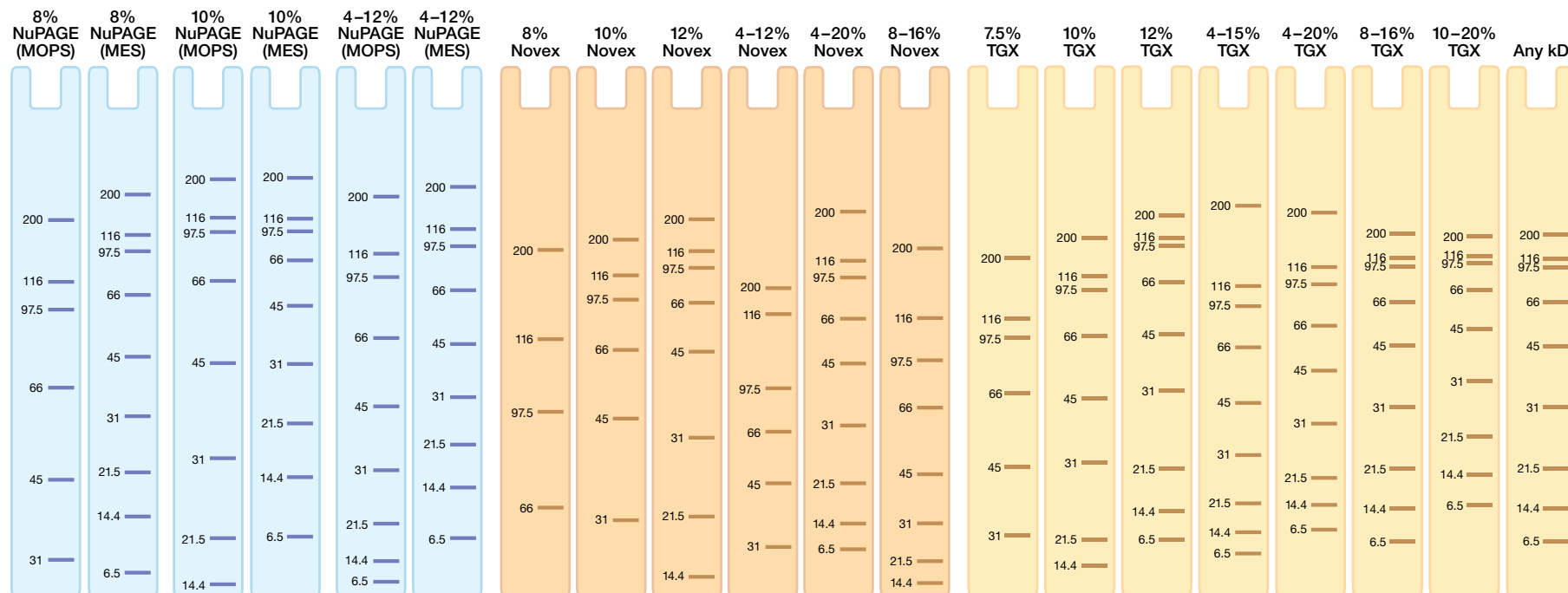
\* Criterion TGX and TGX Stain-Free gels are also available in the 10-20% format but are not listed here as there are no equivalent Life Technologies midi precast gels available.



# Criterion TGX and TGX Stain-Free Precast Gels

Accelerated Electrophoresis with Faster Run Times

## Criterion TGX/TGX Stain-Free and Novex Gel Migration Charts



Side-by-side comparison of migration patterns of broad range SDS-PAGE unstained standards run on Life Technologies midi precast gels: NuPAGE Novex Bis-Tris gels (■), Novex Tris-Glycine gels (■), Criterion TGX/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 Novex Bis-Tris gels run using both MOPS and MES buffers.



### Criterion TGX and TGX Stain-Free precast gels deliver the fastest separations and extended shelf life using the Laemmli PAGE system

- Run times as short as 20 min
- Transfer in as little as 15 min
- Inexpensive Laemmli buffer system, low running costs
- 1 year shelf life

### Related Literature

6007 Criterion TGX Precast Gels Product Information Sheet  
 5974 TGX Stain-Free Precast Gels Product Information Sheet  
 5976 Gel Doc™ EZ Imaging System Brochure

### Criterion TGX Stain-Free precast gels deliver

- Complete protein separations, 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

6133 ChemiDoc™ MP Imaging System Brochure  
 6088 V3 Western Workflow™ Brochure  
 4110001 Criterion Precast Gels Instruction Manual and Applications Guide

Coomassie is a trademark of BASF Aktiengesellschaft.  
 NuPAGE and Novex are trademarks of Life Technologies.