

ELECTROPHORESIS

NuPAGE Bis-Tris Precast Gels to Mini-PROTEAN® TGX™
and Mini-PROTEAN® TGX Stain-Free™ Precast Gels

Catalog Number Conversion Chart



NuPAGE Bis-Tris Precast Gels (Box of 10)			
	Description	Well Volume	Catalog #
10% (MOPS buffer)	NuPAGE Novex 10% Bis-Tris, 1.0 mm, 9-well	28 µl	NP0307BOX →
	NuPAGE Novex 10% Bis-Tris, 1.0 mm, 10-well	25 µl	NP0301BOX →
	NuPAGE Novex 10% Bis-Tris, 1.5 mm, 10-well	37 µl	NP0315BOX →
	NuPAGE Novex 10% Bis-Tris, 1.0 mm, 12-well	20 µl	NP0302BOX →
	NuPAGE Novex 10% Bis-Tris, 1.0 mm, 15-well	15 µl	NP0303BOX →
	NuPAGE Novex 10% Bis-Tris, 1.5 mm, 15-well	25 µl	NP0316BOX →
	NuPAGE Novex 10% Bis-Tris, 1.0 mm, 1-well	700 µl	NP0304BOX →
12% (MOPS buffer)	NuPAGE Novex 12% Bis-Tris, 1.0 mm, 10-well	25 µl	NP0341BOX →
	NuPAGE Novex 12% Bis-Tris, 1.0 mm, 10-well	25 µl	NP0341BOX →
	NuPAGE Novex 12% Bis-Tris, 1.0 mm, 12-well	20 µl	NP0342BOX →
	NuPAGE Novex 12% Bis-Tris, 1.0 mm, 15-well	15 µl	NP0343BOX →
	NuPAGE Novex 12% Bis-Tris, 1.0 mm, 17-well	15 µl	NP0349BOX →
	NuPAGE Novex 12% Bis-Tris, 1.0 mm, 1-well	700 µl	NP0344BOX →
4–12% (MOPS buffer)	NuPAGE Novex 4–12% Bis-Tris, 1.0 mm, 9-well	28 µl	NP0327BOX →
	NuPAGE Novex 4–12% Bis-Tris, 1.0 mm, 10-well	25 µl	NP0321BOX →
	NuPAGE Novex 4–12% Bis-Tris, 1.5 mm, 10-well	37 µl	NP0335BOX →
	NuPAGE Novex 4–12% Bis-Tris, 1.0 mm, 12-well	20 µl	NP0322BOX →
	NuPAGE Novex 4–12% Bis-Tris, 1.0 mm, 15-well	15 µl	NP0323BOX →
	NuPAGE Novex 4–12% Bis-Tris, 1.5 mm, 15-well	25 µl	NP0336BOX →
	NuPAGE Novex 4–12% Bis-Tris, 1.0 mm, 17-well	15 µl	NP0329BOX →
	NuPAGE Novex 4–12% Bis-Tris, 1.0 mm, 1-well	700 µl	NP0324BOX →

Mini-PROTEAN TGX Precast Gels (Box of 10)			
	Catalog #	Well Volume	Description
12%	456-1049	30 µl	Mini-PROTEAN TGX, 12%, 8+1-well*
	456-1043	30 µl	Mini-PROTEAN TGX, 12%, 10-well
	456-1044	50 µl	Mini-PROTEAN TGX, 12%, 10-well
	456-1045	20 µl	Mini-PROTEAN TGX, 12%, 12-well
	456-1046	15 µl	Mini-PROTEAN TGX, 12%, 15-well
	456-1046	15 µl	Mini-PROTEAN TGX, 12%, 15-well
	456-1041	7 cm strip/450 µl	Mini-PROTEAN TGX, 12%, IPG/prep
Any kD™	456-9039	30 µl	Mini-PROTEAN® TGX Any kD™, 8+1-well
	456-9033	30 µl	Mini-PROTEAN TGX Any kD, 10-well
	456-9035	20 µl	Mini-PROTEAN TGX Any kD, 12-well
	456-9036	15 µl	Mini-PROTEAN TGX Any kD, 15-well
	456-9036	15 µl	Mini-PROTEAN TGX Any kD, 15-well
	456-9031	7 cm strip/450 µl	Mini-PROTEAN TGX Any kD, IPG/prep
4–20%	456-1099	30 µl	Mini-PROTEAN TGX, 4–20%, 8+1-well
	456-1093	30 µl	Mini-PROTEAN TGX, 4–20%, 10-well
	456-1094	50 µl	Mini-PROTEAN TGX, 4–20%, 10-well
	456-1095	20 µl	Mini-PROTEAN TGX, 4–20%, 12-well
	456-1096	15 µl	Mini-PROTEAN TGX, 4–20%, 15-well
	456-1096	15 µl	Mini-PROTEAN TGX, 4–20%, 15-well
	456-1096	15 µl	Mini-PROTEAN TGX, 4–20%, 15-well
	456-1091	7 cm strip/450 µl	Mini-PROTEAN TGX, 4–20%, IPG/prep

Mini-PROTEAN TGX Stain-Free Precast Gels (Box of 10)			
	Catalog #	Well Volume	Description
12%	456-8049	30 µl	Mini-PROTEAN TGX Stain-Free, 12%, 8+1-well
	456-8043	30 µl	Mini-PROTEAN TGX Stain-Free, 12%, 10-well
	456-8044	50 µl	Mini-PROTEAN TGX Stain-Free, 12%, 10-well
	456-8045	20 µl	Mini-PROTEAN TGX Stain-Free, 12%, 12-well
	456-8046	15 µl	Mini-PROTEAN TGX Stain-Free, 12%, 15-well
	456-8046	15 µl	Mini-PROTEAN TGX Stain-Free, 12%, 15-well
	456-8041	7 cm strip/450 µl	Mini-PROTEAN TGX Stain-Free, 12%, IPG/prep
Any kD™	456-8129	30 µl	Mini-PROTEAN® TGX Any kD Stain-Free™, 8+1-well
	456-8123	30 µl	Mini-PROTEAN TGX Any kD Stain-Free, 10-well
	456-8125	20 µl	Mini-PROTEAN TGX Any kD Stain-Free, 12-well
	456-8126	15 µl	Mini-PROTEAN TGX Any kD Stain-Free, 15-well
	456-8126	15 µl	Mini-PROTEAN TGX Any kD Stain-Free, 15-well
	456-8121	7 cm strip/450 µl	Mini-PROTEAN TGX Any kD Stain-Free, IPG/prep
4–20%	456-8099	30 µl	Mini-PROTEAN TGX Stain-Free, 4–20%, 8+1-well
	456-8093	30 µl	Mini-PROTEAN TGX Stain-Free, 4–20%, 10-well
	456-8094	50 µl	Mini-PROTEAN TGX Stain-Free, 4–20%, 10-well
	456-8095	20 µl	Mini-PROTEAN TGX Stain-Free, 4–20%, 12-well
	456-8096	15 µl	Mini-PROTEAN TGX Stain-Free, 4–20%, 15-well
	456-8096	15 µl	Mini-PROTEAN TGX Stain-Free, 4–20%, 15-well
	456-8096	15 µl	Mini-PROTEAN TGX Stain-Free, 4–20%, 15-well
	456-8091	7 cm strip/450 µl	Mini-PROTEAN TGX Stain-Free, 4–20%, IPG/prep

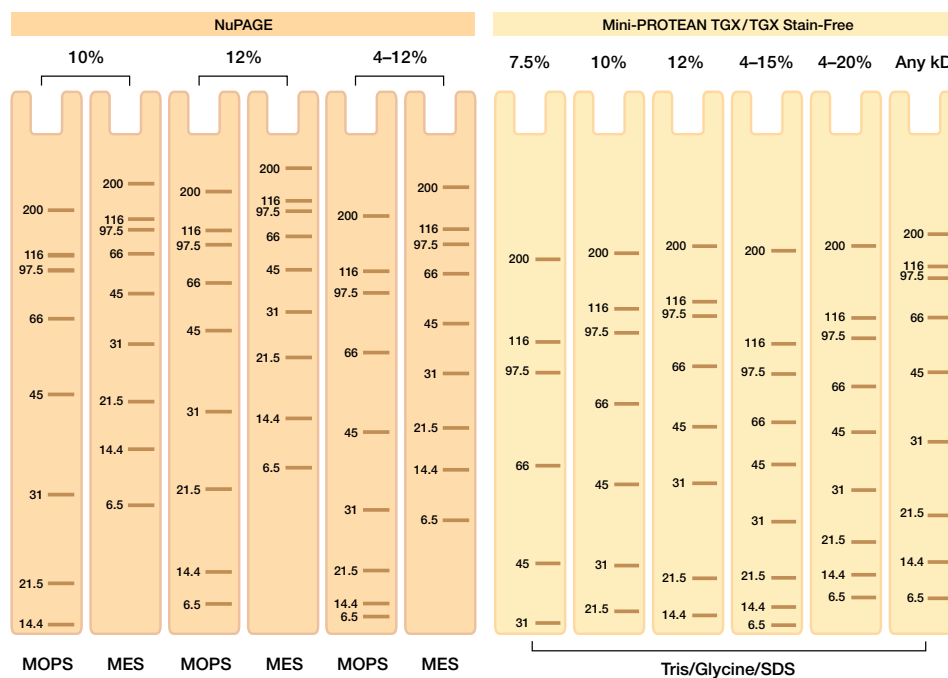
* 8+1-well gels are compatible with multichannel pipets.



Mini-PROTEAN TGX and TGX Stain-Free Precast Gels

The Mini-PROTEAN TGX and TGX Stain-Free (Tris-Glycine eXtended) gels deliver accelerated run times with an extended shelf life of 12 months

- Run times as short as 15 min
- Transfer in as little as 15 min
- Formulated for use with Laemmli system
- Inexpensive buffer system, low running costs
- Robust system for difficult samples
- Bottom-open cassette design for ease of gel handling and blotting
- Compatible with Mini-PROTEAN Tetra, Mini-PROTEAN 3, and Mini-PROTEAN® 3 Dodeca™ systems



Side-by-side comparison of migration patterns of broad range SDS-PAGE unstained standards (catalog #161-0317) run on NuPAGE () and Mini-PROTEAN TGX and TGX Stain-Free precast gels ().

Related Literature

- 1658100 Mini-PROTEAN Precast Gels Instruction Manual and Application Guide
- 5535 Mini-PROTEAN Tetra Cell Brochure
- 5871 Mini-PROTEAN TGX Precast Gels Product Information Sheet
- 5910 Mini-PROTEAN TGX Precast Gel: A Gel for SDS-PAGE with Improved Stability — Comparison with Standard Laemmli Gels Tech Note
- 5911 Mini-PROTEAN TGX Gel: A Versatile and Robust Laemmli-Like Precast Gel for SDS-PAGE Tech Note
- 6133 ChemiDoc™ MP Imaging System Brochure
- 5976 Gel Doc™ EZ Imaging System Brochure
- 6088 V3 Western Workflow™ Brochure

For further information visit www.miniprotean.com.

NuPAGE and Novex are registered trademarks of Life Technologies.