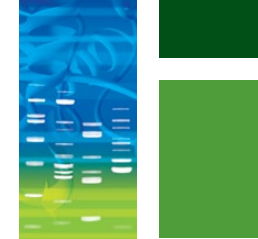


ELECTROPHORESIS

Ready Gel® to Mini-PROTEAN® TGX™ and Mini-PROTEAN® TGX Stain-Free™ Precast Gels
Catalog Number Conversion Chart



Ready Gel Precast Gels (1 gel/pack)		
	Description	Catalog #
7.5% Tris-HCl	Ready Gel, 7.5% Tris-HCl, 10-well, 30 µl	161-1100
	Ready Gel, 7.5% Tris-HCl, 10-well, 30 µl	161-1100
	Ready Gel, 7.5% Tris-HCl, 10-well, 50 µl	161-1154
	Ready Gel, 7.5% Tris-HCl, 12-well, 20 µl	161-1172
	Ready Gel, 7.5% Tris-HCl, 15-well, 15 µl	161-1118
	Ready Gel, 7.5% Tris-HCl, 2-D prep/450 µl	161-1136
10% Tris-HCl	Ready Gel, 10% Tris-HCl, 9-well, 30 µl	161-1191
	Ready Gel, 10% Tris-HCl, 10-well, 30 µl	161-1101
	Ready Gel, 10% Tris-HCl, 10-well, 50 µl	161-1155
	Ready Gel, 10% Tris-HCl, 12-well, 20 µl	161-1173
	Ready Gel, 10% Tris-HCl, 15-well, 15 µl	161-1119
	Ready Gel, 10% Tris-HCl, 2-D prep/450 µl	161-1137
12% Tris-HCl	Ready Gel, 10% Tris-HCl, IPG, 7 cm strip	161-1390
	Ready Gel, 12% Tris-HCl, 10-well, 30 µl	161-1102
	Ready Gel, 12% Tris-HCl, 10-well, 30 µl	161-1102
	Ready Gel, 12% Tris-HCl, 10-well, 50 µl	161-1156
	Ready Gel, 12% Tris-HCl, 12-well, 20 µl	161-1174
	Ready Gel, 12% Tris-HCl, 15-well, 15 µl	161-1120
8-16% Tris-HCl	Ready Gel, 12% Tris-HCl, 2-D prep/450 µl	161-1138
	Ready Gel, 12% Tris-HCl, IPG, 7 cm strip	161-1391
	Ready Gel, 8-16% Tris-HCl, 10-well, 30 µl	161-1222
	Ready Gel, 8-16% Tris-HCl, 10-well, 30 µl	161-1222
	Ready Gel, 8-16% Tris-HCl, 10-well, 50 µl	161-1225
	Ready Gel, 8-16% Tris-HCl, 12-well, 20 µl	161-1226
4-15% Tris-HCl	Ready Gel, 8-16% Tris-HCl, 15-well, 15 µl	161-1223
	Ready Gel, 8-16% Tris-HCl, IPG, 7 cm strip	161-1394
	Ready Gel, 4-15% Tris-HCl, 9-well, 30 µl	161-1194
	Ready Gel, 4-15% Tris-HCl, 10-well, 30 µl	161-1104
	Ready Gel, 4-15% Tris-HCl, 10-well, 50 µl	161-1158
	Ready Gel, 4-15% Tris-HCl, 12-well, 20 µl	161-1176
4-20% Tris-HCl	Ready Gel, 4-15% Tris-HCl, 15-well, 15 µl	161-1122
	Ready Gel, 4-15% Tris-HCl, 2-D prep/450 µl	161-1140
	Ready Gel, 4-15% Tris-HCl, IPG, 7 cm strip	161-1392
	Ready Gel, 4-20% Tris-HCl, 10-well, 30 µl	161-1105
	Ready Gel, 4-20% Tris-HCl, 10-well, 30 µl	161-1105
	Ready Gel, 4-20% Tris-HCl, 10-well, 50 µl	161-1159
	Ready Gel, 4-20% Tris-HCl, 12-well, 20 µl	161-1177
	Ready Gel, 4-20% Tris-HCl, 15-well, 15 µl	161-1123
	Ready Gel, 4-20% Tris-HCl, 2-D prep/450 µl	161-1141
	Ready Gel, 4-20% Tris-HCl, IPG, 7 cm strip	161-1393

Mini-PROTEAN TGX Gels (1 box of 10 gels/pack)		
	Catalog #	Description (Option 1)
7.5%	456-1029	Mini-PROTEAN TGX, 7.5%, 8+1-well, 30 µl*
	456-1023	Mini-PROTEAN TGX, 7.5%, 10-well, 30 µl
	456-1024	Mini-PROTEAN TGX, 7.5%, 10-well, 50 µl
	456-1025	Mini-PROTEAN TGX, 7.5%, 12-well, 20 µl
	456-1026	Mini-PROTEAN TGX, 7.5%, 15-well, 15 µl
	456-1021	Mini-PROTEAN TGX, 7.5%, IPG/prep, 7 cm strip/450 µl
10%	456-1039	Mini-PROTEAN TGX, 10%, 8+1-well, 30 µl
	456-1033	Mini-PROTEAN TGX, 10%, 10-well, 30 µl
	456-1034	Mini-PROTEAN TGX, 10%, 10-well, 50 µl
	456-1035	Mini-PROTEAN TGX, 10%, 12-well, 20 µl
	456-1036	Mini-PROTEAN TGX, 10%, 15-well, 15 µl
	456-1031	Mini-PROTEAN TGX, 10%, IPG/prep, 7 cm strip/450 µl
12%	456-1031	Mini-PROTEAN TGX, 10%, IPG/prep, 7 cm strip/450 µl
	456-1049	Mini-PROTEAN TGX, 12%, 8+1-well, 30 µl
	456-1043	Mini-PROTEAN TGX, 12%, 10-well, 30 µl
	456-1044	Mini-PROTEAN TGX, 12%, 10-well, 50 µl
	456-1045	Mini-PROTEAN TGX, 12%, 12-well, 20 µl
	456-1046	Mini-PROTEAN TGX, 12%, 15-well, 15 µl
8-16%	456-1041	Mini-PROTEAN TGX, 12%, IPG/prep, 7 cm strip/450 µl
	456-1041	Mini-PROTEAN TGX, 12%, IPG/prep, 7 cm strip/450 µl
	456-1109	Mini-PROTEAN TGX, 8-16%, 8+1-well, 30 µl
	456-1103	Mini-PROTEAN TGX, 8-16%, 10-well, 30 µl
	456-1104	Mini-PROTEAN TGX, 8-16%, 10-well, 50 µl
	456-1105	Mini-PROTEAN TGX, 8-16%, 12-well, 20 µl
4-15%	456-1106	Mini-PROTEAN TGX, 8-16%, 15-well, 15 µl
	456-1101	Mini-PROTEAN TGX, 8-16%, IPG/prep, 7 cm strip/450 µl
	456-1089	Mini-PROTEAN TGX, 4-15%, 8+1-well, 30 µl
	456-1083	Mini-PROTEAN TGX, 4-15%, 10-well, 30 µl
	456-1084	Mini-PROTEAN TGX, 4-15%, 10-well, 50 µl
	456-1085	Mini-PROTEAN TGX, 4-15%, 12-well, 20 µl
4-20%	456-1086	Mini-PROTEAN TGX, 4-15%, 15-well, 15 µl
	456-1081	Mini-PROTEAN TGX, 4-15%, IPG/prep, 7 cm strip/450 µl
	456-1081	Mini-PROTEAN TGX, 4-15%, IPG/prep, 7 cm strip/450 µl
	456-1099	Mini-PROTEAN TGX, 4-20%, 8+1-well, 30 µl
	456-1093	Mini-PROTEAN TGX, 4-20%, 10-well, 30 µl
	456-1094	Mini-PROTEAN TGX, 4-20%, 10-well, 50 µl
	456-1095	Mini-PROTEAN TGX, 4-20%, 12-well, 20 µl
	456-1096	Mini-PROTEAN TGX, 4-20%, 15-well, 15 µl
	456-1091	Mini-PROTEAN TGX, 4-20%, IPG/prep, 7 cm strip/450 µl
	456-1091	Mini-PROTEAN TGX, 4-20%, IPG/prep, 7 cm strip/450 µl

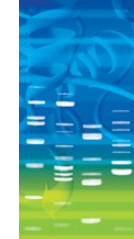
Mini-PROTEAN TGX Stain-Free Gels (1 box of 10 gels/pack)		
	Catalog #	Description (Option 2)
7.5%	456-8029	Mini-PROTEAN TGX Stain-Free, 7.5%, 8+1-well, 30 µl
	456-8023	Mini-PROTEAN TGX Stain-Free, 7.5%, 10-well, 30 µl
	456-8024	Mini-PROTEAN TGX Stain-Free, 7.5%, 10-well, 50 µl
	456-8025	Mini-PROTEAN TGX Stain-Free, 7.5%, 12-well, 20 µl
	456-8026	Mini-PROTEAN TGX Stain-Free, 7.5%, 15-well, 15 µl
	456-8021	Mini-PROTEAN TGX Stain-Free, 7.5%, IPG/prep, 7 cm strip/450 µl
10%	456-8039	Mini-PROTEAN TGX Stain-Free, 10%, 8+1-well, 30 µl
	456-8033	Mini-PROTEAN TGX Stain-Free, 10%, 10-well, 30 µl
	456-8034	Mini-PROTEAN TGX Stain-Free, 10%, 10-well, 50 µl
	456-8035	Mini-PROTEAN TGX Stain-Free, 10%, 12-well, 20 µl
	456-8036	Mini-PROTEAN TGX Stain-Free, 10%, 15-well, 15 µl
	456-8031	Mini-PROTEAN TGX Stain-Free, 10%, IPG/prep, 7 cm strip/450 µl
12%	456-8031	Mini-PROTEAN TGX Stain-Free, 10%, IPG/prep, 7 cm strip/450 µl
	456-8049	Mini-PROTEAN TGX Stain-Free, 12%, 8+1-well, 30 µl
	456-8043	Mini-PROTEAN TGX Stain-Free, 12%, 10-well, 30 µl
	456-8044	Mini-PROTEAN TGX Stain-Free, 12%, 10-well, 50 µl
	456-8045	Mini-PROTEAN TGX Stain-Free, 12%, 12-well, 20 µl
	456-8046	Mini-PROTEAN TGX Stain-Free, 12%, 15-well, 15 µl
8-16%	456-8041	Mini-PROTEAN TGX Stain-Free, 12%, IPG/prep, 7 cm strip/450 µl
	456-8041	Mini-PROTEAN TGX Stain-Free, 12%, IPG/prep, 7 cm strip/450 µl
	456-8109	Mini-PROTEAN TGX Stain-Free, 8-16%, 8+1-well, 30 µl
	456-8103	Mini-PROTEAN TGX Stain-Free, 8-16%, 10-well, 30 µl
	456-8104	Mini-PROTEAN TGX Stain-Free, 8-16%, 10-well, 50 µl
	456-8105	Mini-PROTEAN TGX Stain-Free, 8-16%, 12-well, 20 µl
4-15%	456-8106	Mini-PROTEAN TGX Stain-Free, 8-16%, 15-well, 15 µl
	456-8101	Mini-PROTEAN TGX Stain-Free, 8-16%, IPG/prep, 7 cm strip/450 µl
	456-8089	Mini-PROTEAN TGX Stain-Free, 4-15%, 8+1-well, 30 µl
	456-8083	Mini-PROTEAN TGX Stain-Free, 4-15%, 10-well, 30 µl
	456-8084	Mini-PROTEAN TGX Stain-Free, 4-15%, 10-well, 50 µl
	456-8085	Mini-PROTEAN TGX Stain-Free, 4-15%, 12-well, 20 µl
4-20%	456-8086	Mini-PROTEAN TGX Stain-Free, 4-15%, 15-well, 15 µl
	456-8081	Mini-PROTEAN TGX Stain-Free, 4-15%, IPG/prep, 7 cm strip/450 µl
	456-8081	Mini-PROTEAN TGX Stain-Free, 4-15%, IPG/prep, 7 cm strip/450 µl
	456-8099	Mini-PROTEAN TGX Stain-Free, 4-20%, 8+1-well, 30 µl
	456-8093	Mini-PROTEAN TGX Stain-Free, 4-20%, 10-well, 30 µl
	456-8094	Mini-PROTEAN TGX Stain-Free, 4-20%, 10-well, 50 µl
	456-8095	Mini-PROTEAN TGX Stain-Free, 4-20%, 12-well, 20 µl
	456-8096	Mini-PROTEAN TGX Stain-Free, 4-20%, 15-well, 15 µl
	456-8091	Mini-PROTEAN TGX Stain-Free, 4-20%, IPG/prep, 7 cm strip/450 µl
	456-8091	Mini-PROTEAN TGX Stain-Free, 4-20%, IPG/prep, 7 cm strip/450 µl

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



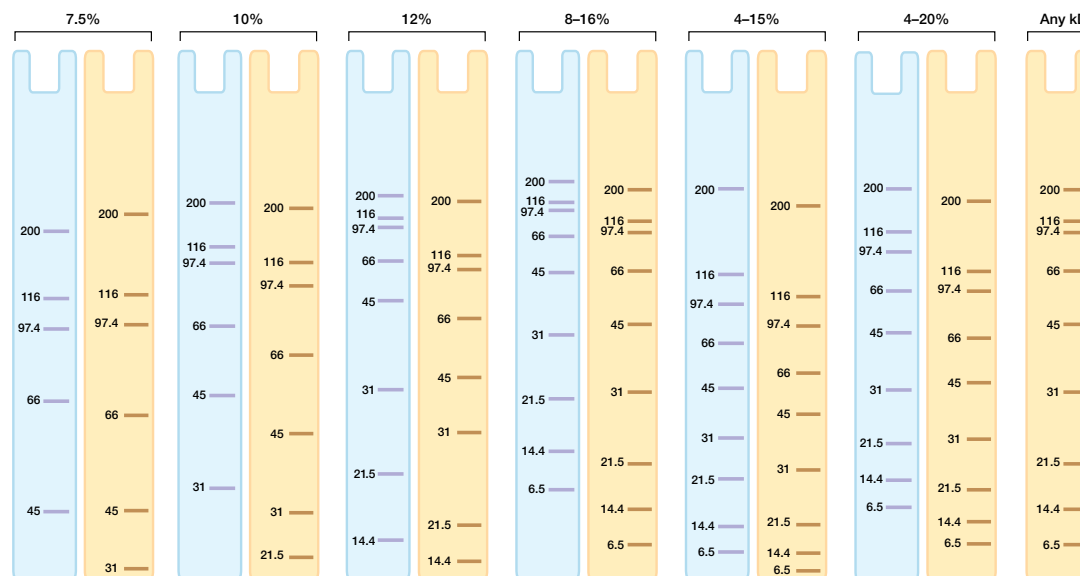
Mini-PROTEAN TGX and TGX Stain-Free Precast Gels

Beyond Ready 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

- Direct drop-in for Laemmli system
- Inexpensive buffer system, low running costs
- Run times as short as 15 min
- Transfer in as little as 15 min
- Accurate molecular weight estimation
- Bottom-open cassette design for ease of gel handling and blotting
- Easy-to-open cassette for faster downstream processing
- 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 Ready Gel precast gels (■) and Mini-PROTEAN TGX and TGX Stain-Free gels (■).

Related Literature

1658100	Mini-PROTEAN Precast Gels Instruction Manual and Application Guide
5535	Mini-PROTEAN Tetra Cell Brochure
5871	Mini-PROTEAN TGX 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
5976	Gel Doc™ EZ Imaging System Brochure
6133	ChemiDoc™ MP Imaging System Brochure
6088	V3 Western Workflow™ Brochure

For further information visit www.miniprotean.com.