

ELECTROPHORESIS

Criterion Tris-HCl to Criterion TGX and Criterion TGX Stain-Free Precast Gels

Catalog Number Conversion Chart (contd)

Criterion Tris-HCl Precast Gels			Criterion TGX Precast Gels		Criterion TGX Stain-Free Precast Gels		
	Description	Catalog #	Catalog #	Description	Catalog #	Description	
4-20%	Criterion, 4-20% Tris-HCl, IPG+1-well, 11 cm IPG Strip	345-0104	→	567-1091	Criterion TGX, 4-20%, IPG+1-well, 11 cm IPG Strip	567-8091	Criterion TGX Stain-Free, 4-20%, IPG+1-well, 11 cm IPG Strip
	Criterion, 4-20% Tris-HCl, prep+2-well, 800 µl	345-0035	→	567-1092	Criterion TGX, 4-20%, prep+2-well, 800 µl	567-8092	Criterion TGX Stain-Free, 4-20%, prep+2-well, 800 µl
	Criterion, 4-20% Tris-HCl, 12+2-well, 45 µl	345-0032	→	567-1093	Criterion TGX, 4-20%, 12+2-well, 45 µl	567-8093	Criterion TGX Stain-Free, 4-20%, 12+2-well, 45 µl
	Criterion, 4-20% Tris-HCl, 18-well, 30 µl	345-0033	→	567-1094	Criterion TGX, 4-20%, 18-well, 30 µl	567-8094	Criterion TGX Stain-Free, 4-20%, 18-well, 30 µl
	Criterion, 4-20% Tris-HCl, 26-well, 15 µl	345-0034	→	567-1095	Criterion TGX, 4-20%, 26-well, 15 µl	567-8095	Criterion TGX Stain-Free, 4-20%, 26-well, 15 µl
8-16%	Criterion, 8-16% Tris-HCl, IPG+1-well, 11 cm IPG Strip	345-0105	→	567-1101	Criterion TGX, 8-16%, IPG+1-well, 11 cm IPG Strip	567-8101	Criterion TGX Stain-Free, 8-16%, IPG+1-well, 11 cm IPG Strip
	Criterion, 8-16% Tris-HCl, prep+2-well, 800 µl	345-0040	→	567-1102	Criterion TGX, 8-16%, prep+2-well, 800 µl	567-8102	Criterion TGX Stain-Free, 8-16%, prep+2-well, 800 µl
	Criterion, 8-16% Tris-HCl, 12+2-well, 45 µl	345-0037	→	567-1103	Criterion TGX, 8-16%, 12+2-well, 45 µl	567-8103	Criterion TGX Stain-Free, 8-16%, 12+2-well, 45 µl
	Criterion, 8-16% Tris-HCl, 18-well, 30 µl	345-0038	→	567-1104	Criterion TGX, 8-16%, 18-well, 30 µl	567-8104	Criterion TGX Stain-Free, 8-16%, 18-well, 30 µl
	Criterion, 8-16% Tris-HCl, 26-well, 15 µl	345-0039	→	567-1105	Criterion TGX, 8-16%, 26-well, 15 µl	567-8105	Criterion TGX Stain-Free, 8-16%, 26-well, 15 µl
10-20%	Criterion, 10-20% Tris-HCl, IPG+1-well, 11 cm IPG Strip	345-0107	→	567-1111	Criterion TGX, 10-20%, IPG+1-well, 11 cm IPG Strip	567-8111	Criterion TGX Stain-Free, 10-20%, IPG+1-well, 11 cm IPG Strip
	Criterion, 10-20% Tris-HCl, prep+2-well, 800 µl	345-0045	→	567-1112	Criterion TGX, 10-20%, prep+2-well, 800 µl	567-8112	Criterion TGX Stain-Free, 10-20%, prep+2-well, 800 µl
	Criterion, 10-20% Tris-HCl, 12+2-well, 45 µl	345-0042	→	567-1113	Criterion TGX, 10-20%, 12+2-well, 45 µl	567-8113	Criterion TGX Stain-Free, 10-20%, 12+2-well, 45 µl
	Criterion, 10-20% Tris-HCl, 18-well, 30 µl	345-0043	→	567-1114	Criterion TGX, 10-20%, 18-well, 30 µl	567-8114	Criterion TGX Stain-Free, 10-20%, 18-well, 30 µl
	Criterion, 10-20% Tris-HCl, 26-well, 15 µl	345-0044	→	567-1115	Criterion TGX, 10-20%, 26-well, 15 µl	567-8115	Criterion TGX Stain-Free, 10-20%, 26-well, 15 µl

Criterion Tris-HCl Stain Free to Criterion TGX Stain-Free Precast Gels

Catalog Number Conversion Chart

Criterion Tris-HCl Stain Free Precast Gels			Criterion TGX Stain-Free Precast Gels		
	Description	Catalog #	Catalog #	Description	
10%	Criterion Stain Free, 10% Tris-HCl, 12+2-well, 45 µl	345-1012	→	567-8033	Criterion TGX Stain-Free, 10%, 12+2-well, 45 µl
	Criterion Stain Free, 10% Tris-HCl, 18-well, 30 µl	345-1018	→	567-8034	Criterion TGX Stain-Free, 10%, 18-well, 30 µl
4-20%	Criterion Stain Free, 4-20% Tris-HCl, 12+2-well, 45 µl	345-0412	→	567-8093	Criterion TGX Stain-Free 4-20%, 12+2-well, 45 µl
	Criterion Stain Free, 4-20% Tris-HCl, 18-well, 30 µl	345-0418	→	567-8094	Criterion TGX Stain-Free 4-20%, 18-well, 30 µl
	Criterion Stain Free, 4-20% Tris-HCl, 26-well, 15 µl	345-0426	→	567-8095	Criterion TGX Stain-Free 4-20%, 26-well, 15 µl
8-16%	Criterion Stain Free, 8-16% Tris-HCl, IPG+1, 11 cm IPG Strip	345-8161	→	567-8101	Criterion TGX Stain-Free, 8-16%, IPG+1-well, 11 cm IPG Strip
	Criterion Stain Free, 8-16% Tris-HCl, 12+2-well, 45 µl	345-8162	→	567-8103	Criterion TGX Stain-Free, 8-16%, 12+2-well, 45 µl
	Criterion Stain Free, 8-16% Tris-HCl, 26-well, 15 µl	345-8166	→	567-8105	Criterion TGX Stain-Free, 8-16%, 26-well, 15 µl

ELECTROPHORESIS

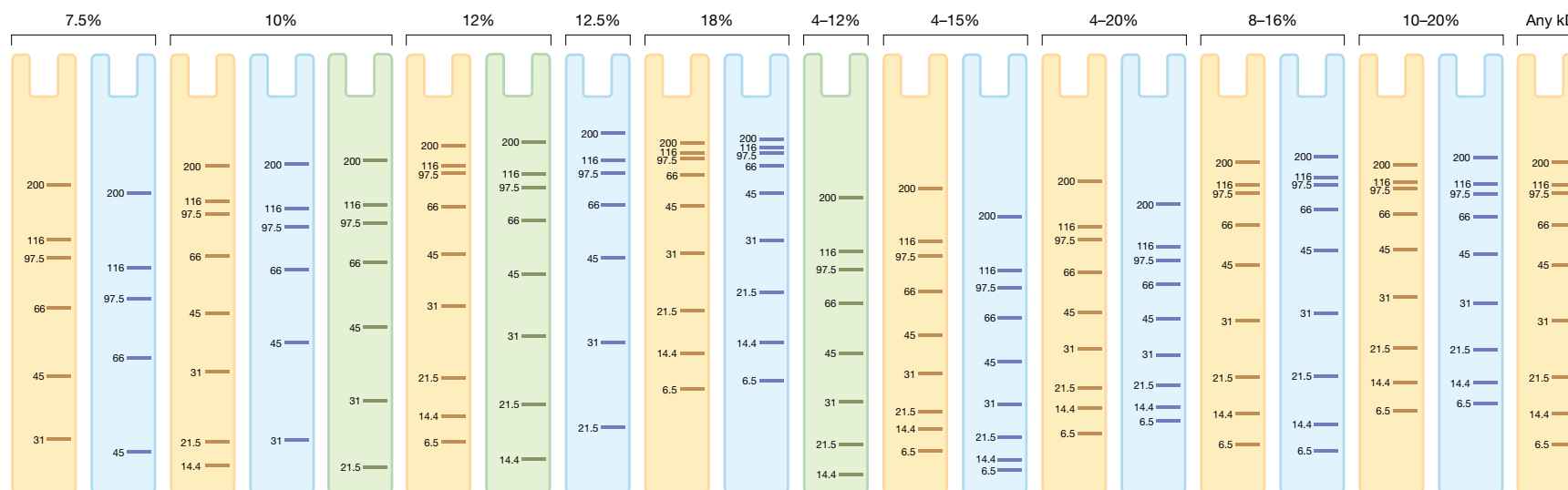
Criterion XT to Criterion TGX and Criterion TGX Stain-Free Precast Gels

Catalog Number Conversion Chart



Criterion XT Precast Gels		
	Description	Catalog #
10%	Criterion XT Bis-Tris, 10%, 12+2-well, 45 µl	345-0111
	Criterion XT Bis-Tris, 10%, 18-well, 30 µl	345-0112
	Criterion XT Bis-Tris, 10%, 26-well, 15 µl	345-0113
12%	Criterion XT Bis-Tris, 12%, 12+2-well, 45 µl	345-0117
	Criterion XT Bis-Tris, 12%, 18-well, 30 µl	345-0118
	Criterion XT Bis-Tris, 12%, 26-well, 15 µl	345-0119
4-12%	Criterion XT Bis-Tris, 4-12%, IPG+1-well, 11 cm IPG Strip	345-0127
	Criterion XT Bis-Tris, 4-12%, prep+2-well, 800 µl	345-0126
	Criterion XT Bis-Tris, 4-12%, 12+2-well, 45 µl	345-0123
	Criterion XT Bis-Tris, 4-12%, 18-well, 30 µl	345-0124
	Criterion XT Bis-Tris, 4-12%, 26-well, 15 µl	345-0125

Criterion TGX Precast Gels			Criterion TGX Stain-Free Precast Gels	
	Catalog #	Description	Catalog #	Description
10%	567-1033	Criterion TGX, 10%, 12+2-well, 45 µl	567-8033	Criterion TGX Stain-Free, 10%, 12+2-well, 45 µl
	567-1034	Criterion TGX, 10%, 18-well, 30 µl	567-8034	Criterion TGX Stain-Free, 10%, 18-well, 30 µl
	567-1035	Criterion TGX, 10%, 26-well, 15 µl	567-8035	Criterion TGX Stain-Free, 10%, 26-well, 15 µl
12%	567-1043	Criterion TGX, 12%, 12+2-well, 45 µl	567-8043	Criterion TGX Stain-Free, 12%, 12+2-well, 45 µl
	567-1044	Criterion TGX, 12%, 18-well, 30 µl	567-8044	Criterion TGX Stain-Free, 12%, 18-well, 30 µl
	567-1045	Criterion TGX, 12%, 26-well, 15 µl	567-8045	Criterion TGX Stain-Free, 12%, 26-well, 15 µl
Any kD	567-1121	Criterion TGX Any kD, IPG+1-well, 11 cm IPG Strip	567-8121	Criterion TGX Any kD Stain-Free, IPG+1-well, 11 cm IPG Strip
	567-1122	Criterion TGX Any kD, prep+2-well, 800 µl	567-8122	Criterion TGX Any kD Stain-Free, prep+2-well, 800 µl
	567-1123	Criterion TGX Any kD, 12+2-well, 45 µl	567-8123	Criterion TGX Any kD Stain-Free, 12+2-well, 45 µl
	567-1124	Criterion TGX Any kD, 18-well, 30 µl	567-8124	Criterion TGX Any kD Stain-Free, 18-well, 30 µl
	567-1125	Criterion TGX Any kD, 26-well, 15 µl	567-8125	Criterion TGX Any kD Stain-Free, 26-well, 15 µl



Side-by-side comparison of migration patterns of broad range SDS-PAGE unstained standards run on Criterion TGX/TGX Stain-Free* gels (■), Criterion Tris-HCl gels (■), and Criterion XT gels (MOPS) (■).

* 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, which are required for stain-free detection.

Criterion TGX and Criterion TGX Stain-Free Precast Gels

Criterion TGX and TGX Stain-Free precast gels deliver the fastest separations and an 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

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

Related Literature

6007	Criterion TGX Precast Gels Product Information Sheet
5974	Criterion TGX Stain-Free Precast Gels Product Information Sheet
6008	Comparison of the Criterion TGX Stain-Free Precast Gel System and Standard Coomassie Staining Procedures for Running and Imaging Protein Gels
4110001	Criterion Gel Application Guide
5976	Gel Doc™ EZ Imaging System Brochure

Coomassie is a trademark of BASF Aktiengesellschaft.

Purchase of Criterion XT Bis-Tris gels, XT MOPS running buffer, XT MES running buffer, XT MOPS buffer kit, and XT MES buffer kit is accompanied by a limited license under U.S. Patent Numbers 6,143,154; 6,096,182; 6,059,948; 5,578,180; 5,922,185; 6,162,338; and 6,783,651, and corresponding foreign patents.

