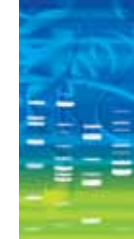


ELECTROPHORESIS

Ready Gel® to Mini-PROTEAN® TBE, TBE-Urea, and Tris-Tricine Precast Gels

Catalog Number Conversion Chart



Ready Gel Precast Gels (1 gel/pack)		
	Description	Catalog #
16.5% Tris-Tricine	Ready Gel Tris-Tricine, 16.5%, 10-well, 30 µl	161-1107
	Ready Gel Tris-Tricine, 16.5%, 10-well, 50 µl	161-1161
	Ready Gel Tris-Tricine, 16.5%, 12-well, 20 µl	161-1179
	Ready Gel Tris-Tricine, 16.5%, 15-well, 15 µl	161-1125
10-20% Tris-Tricine	Ready Gel Tris-Tricine, 10-20%, 10-well, 30 µl	161-1108
	Ready Gel Tris-Tricine, 10-20%, 10-well, 50 µl	161-1162
	Ready Gel Tris-Tricine, 10-20%, 12-well, 20 µl	161-1180
	Ready Gel Tris-Tricine, 10-20%, 15-well, 15 µl	161-1126
5% TBE	Ready Gel TBE, 5%, 10-well, 30 µl	161-1109
	Ready Gel TBE, 5%, 10-well, 50 µl	161-1163
	Ready Gel TBE, 5%, 12-well, 20 µl	161-1181
	Ready Gel TBE, 5%, 15-well, 15 µl	161-1127
10% TBE	Ready Gel TBE, 10%, 10-well, 30 µl	161-1110
	Ready Gel TBE, 10%, 10-well, 50 µl	161-1164
	Ready Gel TBE, 10%, 12-well, 20 µl	161-1182
	Ready Gel TBE, 10%, 15-well, 15 µl	161-1128
15% TBE	Ready Gel TBE, 15%, 10-well, 30 µl	161-1228
	-	-
	Ready Gel TBE, 15%, 12-well, 20 µl	161-1232
	Ready Gel TBE, 15%, 15-well, 15 µl	161-1229
4-20% TBE	Ready Gel TBE, 4-20%, 10-well, 30 µl	161-1234
	Ready Gel TBE, 4-20%, 10-well, 50 µl	161-1237
	-	-
	Ready Gel TBE, 4-20%, 15-well, 15 µl	161-1235
10% TBE-Urea	Ready Gel TBE-Urea, 10%, 10-well, 30 µl	161-1116
	Ready Gel TBE-Urea, 10%, 15-well, 15 µl	161-1134
15% TBE-Urea	Ready Gel TBE-Urea, 15% 10-well, 30 µl	161-1117
	Ready Gel TBE-Urea, 15% 12-well, 20 µl	161-1189
	Ready Gel TBE-Urea, 15%, 15-well, 15 µl	161-1135

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

→

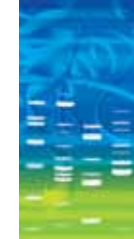
→

Mini-PROTEAN Precast Gels (1 box of 2 gels/pack)		
Catalog #		Description
456-3063		Mini-PROTEAN Tris-Tricine, 16.5%, 10-well, 30 µl
456-3064		Mini-PROTEAN Tris-Tricine, 16.5%, 10-well, 50 µl
456-3065		Mini-PROTEAN Tris-Tricine, 16.5%, 12-well, 20 µl
456-3066		Mini-PROTEAN Tris-Tricine, 16.5%, 15-well, 15 µl
456-3113		Mini-PROTEAN Tris-Tricine, 10-20%, 10-well, 30 µl
456-3114		Mini-PROTEAN Tris-Tricine, 10-20%, 10-well, 50 µl
456-3115		Mini-PROTEAN Tris-Tricine, 10-20%, 12-well, 20 µl
456-3116		Mini-PROTEAN Tris-Tricine, 10-20%, 15-well, 15 µl
456-5013		Mini-PROTEAN TBE, 5%, 10-well, 30 µl
456-5014		Mini-PROTEAN TBE, 5%, 10-well, 50 µl
456-5015		Mini-PROTEAN TBE, 5%, 12-well, 20 µl
456-5016		Mini-PROTEAN TBE, 5%, 15-well, 15 µl
456-5033		Mini-PROTEAN TBE, 10%, 10-well, 30 µl
456-5034		Mini-PROTEAN TBE, 10%, 10-well, 50 µl
456-5035		Mini-PROTEAN TBE, 10%, 12-well, 20 µl
456-5036		Mini-PROTEAN TBE, 10%, 15-well, 15 µl
456-5053		Mini-PROTEAN TBE, 15%, 10-well, 30 µl
456-5054		Mini-PROTEAN TBE, 15%, 10-well, 50 µl
456-5055		Mini-PROTEAN TBE, 15%, 12-well, 20 µl
456-5056		Mini-PROTEAN TBE, 15%, 15-well, 15 µl
456-5093		Mini-PROTEAN TBE, 4-20%, 10-well, 30 µl
456-5094		Mini-PROTEAN TBE, 4-20%, 10-well, 50 µl
456-5095		Mini-PROTEAN TBE, 4-20%, 12-well, 20 µl
456-5096		Mini-PROTEAN TBE, 4-20%, 15-well, 15 µl
456-6033		Mini-PROTEAN TBE-Urea, 10%, 10-well, 30 µl
456-6036		Mini-PROTEAN TBE-Urea, 10%, 15-well, 15 µl
456-6053		Mini-PROTEAN TBE-Urea, 15%, 10-well, 30 µl
456-6055		Mini-PROTEAN TBE-Urea, 15%, 12-well, 20 µl
456-6056		Mini-PROTEAN TBE-Urea, 15%, 15-well, 15 µl



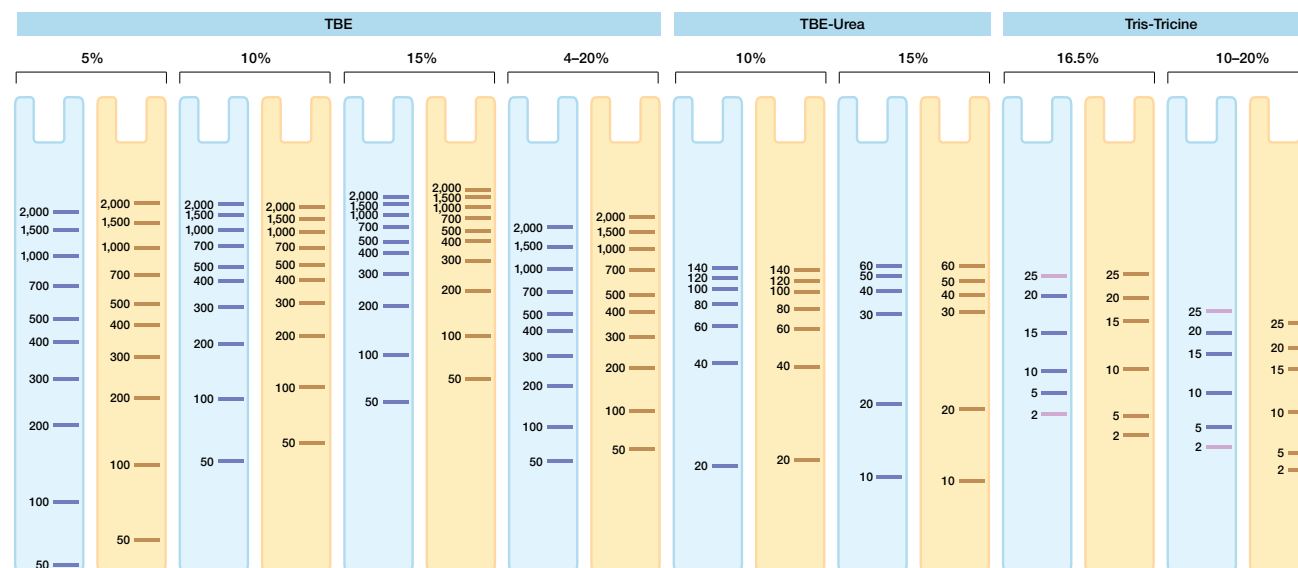
# Mini-PROTEAN Precast Gels

## Beyond Ready Gels



### Mini-PROTEAN precast gels advance the convenience of Ready Gel precast gels to the next level.

- Same gel chemistry as the Ready Gel precast gels
- Easy-to-open cassette for faster downstream processing
- Compatibility with Mini-PROTEAN Tetra, Mini-PROTEAN 3, and Mini-PROTEAN® Dodeca™ systems



Side-by-side comparison of migration patterns on Mini-PROTEAN (■) and Ready Gel (■) precast gels. Numbers indicate base pairs (TBE and TBE-Urea) or molecular weights in kD (Tris-Tricine).



### Related Literature

- Bulletin 1658100 Mini-PROTEAN Precast Gels Instruction Manual and Application Guide
- Bulletin 5535 Mini-PROTEAN Tetra Cell Brochure
- Bulletin 6062 Mini-PROTEAN TBE, TBE-Urea, and Tris-Tricine Precast Gel Quick Start Guide