

Ready Gel System

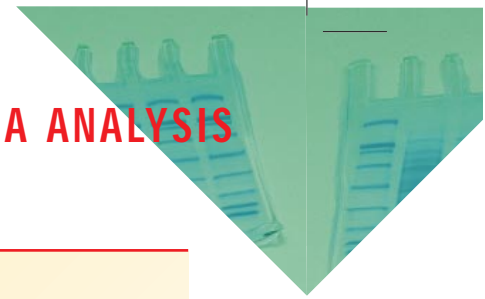
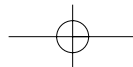
RESOURCE GUIDE ADDENDUM

This addendum includes new electrophoresis products introduced since the Ready Gel Resource Guide was published and updated information for a few products. Keep this addendum as a quick reference guide, and use it to find the electrophoresis application you're interested in and all related products available from Bio-Rad.






Applications:

- Protein Molecular Weight Analysis
- Native Protein Analysis
- Peptide Analysis
- Isoelectric Point Analysis/IEF
- Protease Analysis
- Oligosaccharide Analysis
- Glycoprotein Analysis
- dsDNA Analysis
- ssDNA, RNA Analysis
- Blot Processing Buffers

BIO-RAD

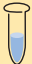



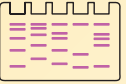


NUCLEIC ACIDS: FOR dsDNA ANALYSIS

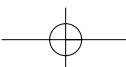
Step	Ready Gel System Product	Catalog #																																																	
 Sample Preparation	Nucleic Acid Sample Buffer, 5x, 10 ml	161-0733																																																	
	<table border="1"> <thead> <tr> <th rowspan="2">MW Range Resolved (bp)</th> <th rowspan="2">TBE Gel %T</th> <th colspan="5">Single gel catalog numbers - order as many as you need!</th> <th rowspan="2">Suggested Standard</th> <th rowspan="2">Standards Catalog #</th> </tr> <tr> <th>9 Well</th> <th>10 Well</th> <th>10 Well/50 µl</th> <th>12 Well</th> <th>15 Well</th> </tr> </thead> <tbody> <tr> <td>200-2,000</td> <td>5%</td> <td>161-1199</td> <td>161-1109</td> <td>161-1163</td> <td>161-1181</td> <td>161-1127</td> <td>NEW! EZ Load 20 bp Ruler,</td> <td>170-8351</td> </tr> <tr> <td>50-1,500</td> <td>10%</td> <td>161-1200</td> <td>161-1110</td> <td>161-1164</td> <td>161-1182</td> <td>161-1128</td> <td>NEW! EZ Load 100 bp Ruler,</td> <td>170-8352</td> </tr> <tr> <td>20-1,000</td> <td>15%</td> <td>161-1233</td> <td>161-1228</td> <td>161-1231</td> <td>161-1232</td> <td>161-1229</td> <td>AmpliSize DNA Size Standard, 50-2,000 bp</td> <td>170-8200</td> </tr> <tr> <td>10-2,000</td> <td>4-20%</td> <td>161-1239</td> <td>161-1234</td> <td>161-1237</td> <td>161-1238</td> <td>161-1235</td> <td>DNA Size Standard, 20 bp Ladder</td> <td>170-8201</td> </tr> </tbody> </table>	MW Range Resolved (bp)	TBE Gel %T	Single gel catalog numbers - order as many as you need!					Suggested Standard	Standards Catalog #	9 Well	10 Well	10 Well/50 µl	12 Well	15 Well	200-2,000	5%	161-1199	161-1109	161-1163	161-1181	161-1127	NEW! EZ Load 20 bp Ruler,	170-8351	50-1,500	10%	161-1200	161-1110	161-1164	161-1182	161-1128	NEW! EZ Load 100 bp Ruler,	170-8352	20-1,000	15%	161-1233	161-1228	161-1231	161-1232	161-1229	AmpliSize DNA Size Standard, 50-2,000 bp	170-8200	10-2,000	4-20%	161-1239	161-1234	161-1237	161-1238	161-1235	DNA Size Standard, 20 bp Ladder	170-8201
MW Range Resolved (bp)	TBE Gel %T			Single gel catalog numbers - order as many as you need!							Suggested Standard	Standards Catalog #																																							
		9 Well	10 Well	10 Well/50 µl	12 Well	15 Well																																													
200-2,000	5%	161-1199	161-1109	161-1163	161-1181	161-1127	NEW! EZ Load 20 bp Ruler,	170-8351																																											
50-1,500	10%	161-1200	161-1110	161-1164	161-1182	161-1128	NEW! EZ Load 100 bp Ruler,	170-8352																																											
20-1,000	15%	161-1233	161-1228	161-1231	161-1232	161-1229	AmpliSize DNA Size Standard, 50-2,000 bp	170-8200																																											
10-2,000	4-20%	161-1239	161-1234	161-1237	161-1238	161-1235	DNA Size Standard, 20 bp Ladder	170-8201																																											
 Precast Gel Selection																																																			
	<table border="1"> <thead> <tr> <th>10x TBE Buffer</th> <th></th> <th>Catalog #</th> </tr> </thead> <tbody> <tr> <td></td> <td>1 L</td> <td>161-0733</td> </tr> <tr> <td></td> <td>5 L NEW!</td> <td>161-0770</td> </tr> <tr> <td></td> <td>6 x 1 L</td> <td>161-0756</td> </tr> </tbody> </table>	10x TBE Buffer		Catalog #		1 L	161-0733		5 L NEW!	161-0770		6 x 1 L	161-0756																																						
10x TBE Buffer		Catalog #																																																	
	1 L	161-0733																																																	
	5 L NEW!	161-0770																																																	
	6 x 1 L	161-0756																																																	
 Running Buffer																																																			
 Electrophoresis Cell and Power Supply	Mini-PROTEAN 3 Electrophoresis Module	165-3302																																																	
	PowerPac 300 (100/120 V)	165-5050																																																	
 Gel Staining	Ethidium Bromide Solution, 10 ml, 10 mg/ml	161-0433																																																	
	Ethidium Bromide Tablets, 10 x 11 mg	161-0430																																																	
	Silver Stain Plus Kit	161-0449																																																	
	Silver Stain Kit	161-0443																																																	

Run conditions per 2 gels:
 Begin - 100 V (C), 26 mA, 45-105 minutes
 End - 100 V (C), 22 mA

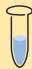
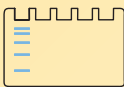


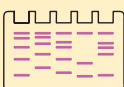
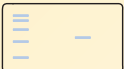
NUCLEIC ACIDS: ssDNA, RNA ANALYSIS

Step	Ready Gel System Product	Catalog #																																								
 Sample Preparation	TBE-Urea Sample Buffer, 30 ml	161-0768																																								
	<table border="1"> <thead> <tr> <th rowspan="2">MW Range Resolved (bases)</th> <th rowspan="2">TBE-Urea Gel %T</th> <th colspan="5">Single gel catalog numbers - order as many as you need!</th> <th rowspan="2">Suggested Standard</th> <th rowspan="2">Standards Catalog #</th> </tr> <tr> <th>9 Well</th> <th>10 Well</th> <th>12 Well</th> <th>15 Well</th> <th>2-D/Prep</th> </tr> </thead> <tbody> <tr> <td>50-1,000</td> <td>5%</td> <td>161-1205</td> <td>161-1115</td> <td>161-1187</td> <td>161-1133</td> <td>161-1151</td> <td>AmpliSize DNA Size Standard, 50-2,000 bp</td> <td>170-8200</td> </tr> <tr> <td>25-200</td> <td>10%</td> <td>161-1206</td> <td>161-1116</td> <td>161-1188</td> <td>161-1134</td> <td>161-1152</td> <td>DNA Size Standard, 20 bp Ladder</td> <td>170-8201</td> </tr> <tr> <td>10-50</td> <td>15%</td> <td>161-1207</td> <td>161-1117</td> <td>161-1189</td> <td>161-1135</td> <td>161-1153</td> <td>Oligo Standard, 8-32 mer</td> <td>161-0430</td> </tr> </tbody> </table>	MW Range Resolved (bases)	TBE-Urea Gel %T	Single gel catalog numbers - order as many as you need!					Suggested Standard	Standards Catalog #	9 Well	10 Well	12 Well	15 Well	2-D/Prep	50-1,000	5%	161-1205	161-1115	161-1187	161-1133	161-1151	AmpliSize DNA Size Standard, 50-2,000 bp	170-8200	25-200	10%	161-1206	161-1116	161-1188	161-1134	161-1152	DNA Size Standard, 20 bp Ladder	170-8201	10-50	15%	161-1207	161-1117	161-1189	161-1135	161-1153	Oligo Standard, 8-32 mer	161-0430
MW Range Resolved (bases)	TBE-Urea Gel %T			Single gel catalog numbers - order as many as you need!							Suggested Standard	Standards Catalog #																														
		9 Well	10 Well	12 Well	15 Well	2-D/Prep																																				
50-1,000	5%	161-1205	161-1115	161-1187	161-1133	161-1151	AmpliSize DNA Size Standard, 50-2,000 bp	170-8200																																		
25-200	10%	161-1206	161-1116	161-1188	161-1134	161-1152	DNA Size Standard, 20 bp Ladder	170-8201																																		
10-50	15%	161-1207	161-1117	161-1189	161-1135	161-1153	Oligo Standard, 8-32 mer	161-0430																																		
 Precast Gel and Standard Selection																																										
	<table border="1"> <thead> <tr> <th>10x TBE Buffer</th> <th></th> <th>Catalog #</th> </tr> </thead> <tbody> <tr> <td></td> <td>1 L</td> <td>161-0733</td> </tr> <tr> <td></td> <td>5 L NEW!</td> <td>161-0770</td> </tr> <tr> <td></td> <td>6 x 1 L</td> <td>161-0756</td> </tr> </tbody> </table>	10x TBE Buffer		Catalog #		1 L	161-0733		5 L NEW!	161-0770		6 x 1 L	161-0756																													
10x TBE Buffer		Catalog #																																								
	1 L	161-0733																																								
	5 L NEW!	161-0770																																								
	6 x 1 L	161-0756																																								
 Running Buffer																																										
 Electrophoresis Cell and Power Supply	Mini-PROTEAN 3 Electrophoresis Module	165-3302																																								
	PowerPac 300 (100/120 V)	165-5050																																								
 Gel Staining	Ethidium Bromide Solution, 10 ml, 10 mg/ml	161-0433																																								
	Ethidium Bromide Tablets, 10 x 11 mg	161-0430																																								
	Radiant Red® RNA Stain Solution	170-3122																																								

Run conditions per 2 gels:
 Begin - 200 V (C), 30 mA
 End - 200 V (C), 20 mA, 40-90 minutes

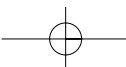


SDS-PAGE: PROTEIN MOLECULAR WEIGHT ANALYSIS


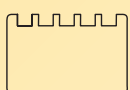


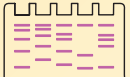
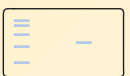
Step	Ready Gel System Product	Catalog #							
 Sample Preparation	Laemmli Sample Buffer, 30 ml	161-0737							
	2-mercaptoethanol, 25 ml	161-0710							
 Ready Gel and Standards Selection	Single gel catalog numbers - order as many as you need!						Suggested SDS-PAGE Standard	Standard Catalog #	
	MW Range Resolved	Tris-HCl Gel %T	9 Well	10 Well	10 Well/50 µl	12 Well	15 Well	2-D/Prep	
	100-250 kDa	5%	161-1215	161-1210	161-1213	161-1214	161-1211	161-1212	High range 161-0303
	40-200 kDa	7.5%	161-1190	161-1100	161-1154	161-1172	161-1118	161-1136	High range 161-0303
	30-150 kDa	10%	161-1191	161-1101	161-1155	161-1173	161-1119	161-1137	Broad range 161-0317
	20-120 kDa	12%	161-1192	161-1102	161-1156	161-1174	161-1120	161-1138	Low range 161-0304
	10-100 kDa	15%	161-1193	161-1103	161-1157	161-1175	161-1121	161-1139	Low range 161-0304
	6-50 kDa	18%	161-1221	161-1216	161-1219	161-1220	161-1217	161-1218	Low range 161-0304
	20-250 kDa	4-15%	161-1194	161-1104	161-1158	161-1176	161-1122	161-1140	Broad range 161-0317
	10-200 kDa	4-20%	161-1195	161-1105	161-1159	161-1177	161-1123	161-1141	Broad range 161-0317
	6-70 kDa	8-16%	161-1227	161-1222	161-1225	161-1226	161-1223	161-1224	Broad range 161-0317
10-100 kDa	10-20%	161-1196	161-1106	161-1160	161-1178	161-1124	161-1142	Broad range 161-0317	
 Running Buffer			Catalog #						
	10x Tris/Glycine/SDS	1 L	161-0732						
		5 L NEW!	161-0772						
		6 x 1 L	161-0755						
 Electrophoresis Cell and Power Supply	Mini-PROTEAN 3 Electrophoresis Module	165-3302							
	PowerPac 300 (100/120 V)	165-5050							
	PowerPac 300 (220/240 V)	165-5051							
Run conditions per 2 gels: Begin - 200 V (C), 100 mA End - 200 V (C), 60 mA, 35 minutes									
 Protein Stain	Bio-Safe Coomassie	1 L NEW!	161-0786						
		5 L NEW!	161-0787						
	Premixed Coomassie R-250 Kit		161-0435						
	Silver Stain Plus Kit		161-0449						
	Silver Stain Kit		161-0443						
	SYPRO Orange Fluorescent Stain		170-3120						
Zinc Stain Kit		161-0440							
Transfer Equipment, Membranes, and Detection Reagents (see Ready Gel Resource Guide page 54)									
 Alternative Detection: Western Blotting	Trans-Blot/Mini Trans-Blot, Transfer Buffer:								
	10x Tris/Glycine	1 L	161-0734						
		5 L NEW!	161-0771						
		6 x 1 L	161-0757						
	Semi-Dry Transfer, Transfer Buffer:								
10x Tris/CAPS	1 L NEW!	161-0778							



See page 44 of the Ready Gel Resource Guide for more on stains and stain kits

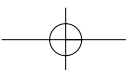


NATIVE PAGE: FOR NATIVE PROTEIN ANALYSIS

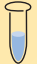



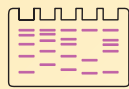
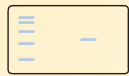
Step	Ready Gel System Product	Catalog #							
 Sample Preparation	Native Sample Buffer, 30 ml	161-0738							
	Single gel catalog numbers - order as many as you need!								
	Tris-HCl Gel %T	9 Well	10 Well	10 Well/50 µl	12 Well	15 Well	2-D/Prep	Suggested SDS-PAGE Standard	Standard Catalog #
	5%	161-1215	161-1210	161-1213	161-1214	161-1211	161-1212	Prestained Blue, Broad	161-0318
	7.5%	161-1190	161-1100	161-1154	161-1172	161-1118	161-1136	Kaleidoscope, Broad	161-0324
	10%	161-1191	161-1101	161-1155	161-1173	161-1119	161-1137		
	12%	161-1192	161-1102	161-1156	161-1174	161-1120	161-1138		
	15%	161-1193	161-1103	161-1157	161-1175	161-1121	161-1139		
	18%	161-1221	161-1216	161-1219	161-1220	161-1217	161-1218		
	4-15%	161-1194	161-1104	161-1158	161-1176	161-1122	161-1140		
4-20%	161-1195	161-1105	161-1159	161-1177	161-1123	161-1141			
8-16%	161-1227	161-1222	161-1225	161-1226	161-1223	161-1224			
10-20%	161-1196	161-1106	161-1160	161-1178	161-1124	161-1142			
 Ready Gel Selection							Use Prestained standards for reference only, not for molecular weight determination.		
 Running Buffer	10x Tris/Glycine Buffer	1 L	161-0734						
		5 L NEW!	161-0771						
		6 x 1 L	161-0757						
 Electrophoresis Cell and Power Supply	Mini-PROTEAN 3 Electrophoresis Module		165-3302						
	PowerPac 300 (100/120 V)		165-5050						
	PowerPac 300 (220/240 V)		165-5051						
		Run conditions per 2 gels:							
		Begin - 200 V (C), 100 mA							
		End - 200 V (C), 60 mA, 35 minutes							
 Protein Stain	Bio-Safe Coomassie	1 L NEW!	161-0786	Tip: Use any stain except Cu and Zn stains for native separations. Cu and Zn stains require SDS in the gel. SYPRO also requires SDS, however, the SYPRO instruction manual includes a protocol for pre-treating native gels with an SDS containing solution prior to staining.					
		5 L NEW!	161-0787						
	Premixed Coomassie R-250 Kit		161-0435						
	Silver Stain Plus Kit		161-0449						
	Bio-Rad Silver Stain Kit		161-0443						
Transfer Equipment, Membranes, and Detection Reagents (see Ready Gel Resource Guide page 55)									
 Alternative Detection: Western Blotting	Trans-Blot/Mini Trans-Blot, Transfer Buffer:								
	10x Tris/Glycine	1 L	161-0734						
		5 L NEW!	161-0771						
		6 x 1 L	161-0757						
Semi-Dry Transfer, Transfer Buffer:									
10 x Tris/CAPS	1 L NEW!	161-0778							



See page 55 of the Ready Gel Resource Guide for more on transfer equipment

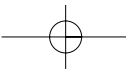


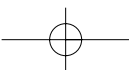
SDS-PAGE: FOR PEPTIDE AND SMALL PROTEIN ANALYSIS

Step	Ready Gel System Product	Catalog #							
 Sample Preparation	Tricine Sample Buffer, 30 ml	161-0739							
	2-mercaptoethanol, 25 ml	161-0710							
 Ready Gel and Standard Selection	Single gel catalog numbers - order as many as you need!						Suggested Standard	Standard Catalog #	
	MW Range Resolved	Tris-Tricine Gel %T	9 Well	10 Well	10 Well/50 µl	12 Well	15 Well	2-D/Prep	
	4-30 kDa	16.5%	161-1197	161-1107	161-1161	161-1179	161-1125	161-1143	Polypeptide SDS-PAGE
5-40 kDa	10-20%	161-1198	161-1108	161-1162	161-1180	161-1126	161-1144	Kaleidoscope Polypeptide Standard	
 Running Buffer			Catalog #						
	10x Tris/Tricine/SDS Buffer	1 L	161-0744						
 Electrophoresis Cell and Power Supply	6 x 1 L	161-0760							
	Mini-PROTEAN 3 Electrophoresis Module	165-3302							
	PowerPac 300 (100/120 V)	165-5050							
	PowerPac 300 (220/240 V)	165-5051							
 Protein Stain	Run conditions per 2 gels: Begin - 100 V (C), 65 mA End - 100 V (C), 35 mA, 100 minutes								
	Bio-Safe Coomassie	1 L NEW!	161-0786						
		5 L NEW!	161-0787						
	Coomassie G-250, 10 g		161-0406	Recommended					
	Silver Stain Plus Kit		161-0449	Silver Stains recommended for peptides larger than 3.5 Kd					
 Alternative Detection: Western Blotting	Silver Stain Kit		161-0443						
	Transfer Equipment, Membranes, and Detection Reagents (see Ready Gel Resource Guide page 54)								
	Trans-Blot/Mini Trans-Blot, Transfer Buffer:								
	10x Tris/Glycine	1 L	161-0734						
		5 L NEW!	161-0771						
	6 x 1 L	161-0757							
	Semi-Dry Transfer, Transfer Buffer:								
10x Tris/CAPS	1 L NEW!	161-0778							


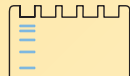



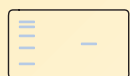


See page 38 of the Ready Gel Resource Guide for more on power supplies





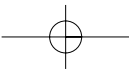
IEF: ISOELECTRIC POINT ANALYSIS

Step	Ready Gel System Product	Catalog #																																								
 Sample Preparation	IEF Sample Buffer, 30 ml	161-0763																																								
	 Ready Gel and Standard Selection																																									
	<table border="1"> <thead> <tr> <th>IEF Ready Gels %T</th> <th>pH Range Resolved</th> <th colspan="6">Single gel catalog numbers - order as many as you need!</th> <th>Suggested SDS-PAGE Standard</th> <th>Standard Catalog #</th> </tr> </thead> <tbody> <tr> <td>5%</td> <td>pH 3-10</td> <td>9 Well</td> <td>10 Well</td> <td>10 Well/50 µl</td> <td>12 Well</td> <td>15 Well</td> <td>2-D/Prep</td> <td>IEF Standard, pI 4.6-9.6</td> <td>161-0310</td> </tr> <tr> <td>5%</td> <td>pH 5-8</td> <td>161-1201</td> <td>161-1111</td> <td>161-1165</td> <td>161-1183</td> <td>161-1129</td> <td>161-1147</td> <td>IEF Standard, pI 4.6-9.6</td> <td>161-0310</td> </tr> <tr> <td></td> <td></td> <td>161-1202</td> <td>161-1112</td> <td>161-1166</td> <td>161-1184</td> <td>161-1130</td> <td>161-1148</td> <td></td> <td></td> </tr> </tbody> </table>	IEF Ready Gels %T	pH Range Resolved	Single gel catalog numbers - order as many as you need!						Suggested SDS-PAGE Standard	Standard Catalog #	5%	pH 3-10	9 Well	10 Well	10 Well/50 µl	12 Well	15 Well	2-D/Prep	IEF Standard, pI 4.6-9.6	161-0310	5%	pH 5-8	161-1201	161-1111	161-1165	161-1183	161-1129	161-1147	IEF Standard, pI 4.6-9.6	161-0310			161-1202	161-1112	161-1166	161-1184	161-1130	161-1148			
IEF Ready Gels %T	pH Range Resolved	Single gel catalog numbers - order as many as you need!						Suggested SDS-PAGE Standard	Standard Catalog #																																	
5%	pH 3-10	9 Well	10 Well	10 Well/50 µl	12 Well	15 Well	2-D/Prep	IEF Standard, pI 4.6-9.6	161-0310																																	
5%	pH 5-8	161-1201	161-1111	161-1165	161-1183	161-1129	161-1147	IEF Standard, pI 4.6-9.6	161-0310																																	
		161-1202	161-1112	161-1166	161-1184	161-1130	161-1148																																			
 Running Buffer	10x Anode Buffer, 250 ml	161-0761																																								
	10x Cathode Buffer, 250 ml	161-0762																																								
 Electrophoresis Cell and Power Supply	Mini-PROTEAN 3 Electrophoresis Module	165-3302																																								
	PowerPac 1000 (100/120 V)	165-5054																																								
	PowerPac 1000 (220/240 V)	165-5055																																								
	Run conditions per 2 gels stepwise: 100 V (C), 5-25 mA 1 hour, 250 V (C), 5-25 mA 1 hour, 500 V (C), 5-25 mA, 30 minutes																																									
 Gel Stain	Premixed IEF Stain, 1 L	161-0434																																								
	Silver Stain Plus Kit	161-0449																																								
	Silver Stain Kit	161-0443																																								
 Alternative Detection: Western Blotting	Transfer Equipment, Membranes, and Detection Reagents (see Ready Gel Resource Guide page 54)																																									
	Trans-Blot/Mini Trans-Blot, Transfer Buffer: 0.7% Acetic Acid, see Tip #25, page 55 in the Ready Gel Resource Guide																																									

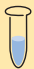





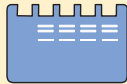
1x Formulations:
Anode Buffer: 7 mM Phosphoric Acid
Cathode Buffer: 20 mM Lysine/20 mM Arginine

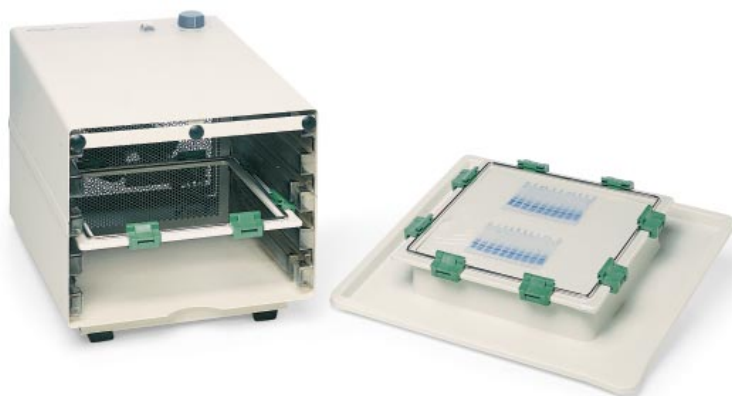


See page 42 of the Ready Gel Resource Guide for more on premixed sample and running buffers



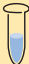


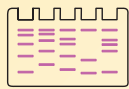
ZYMOGRAM: FOR PROTEASE ANALYSIS

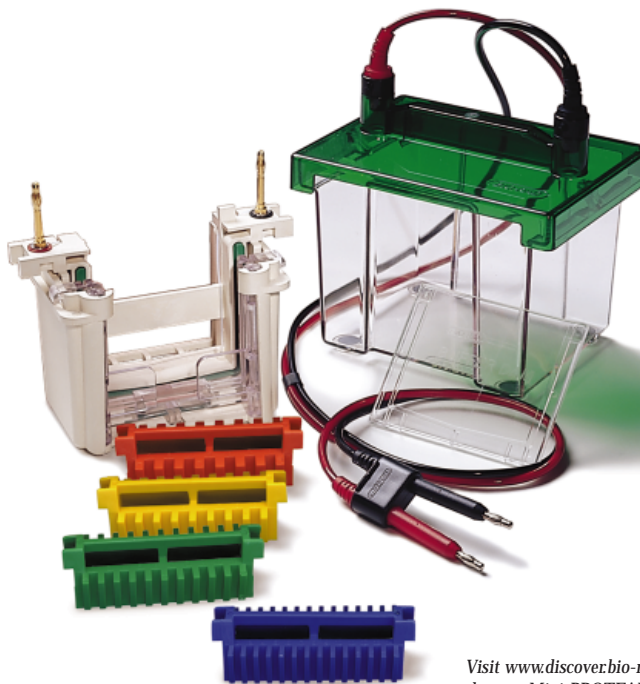
Step	Ready Gel System Product	Catalog #																																									
 Sample Preparation	Zymogram Sample Buffer, 30 ml	161-0764																																									
 Ready Gel and Standard Selection	<table border="1"> <thead> <tr> <th rowspan="2">MW Range Resolved</th> <th rowspan="2">Zymogram Gel %T With Contents</th> <th colspan="5">Single gel catalog numbers - order as many as you need!</th> <th rowspan="2">Suggested SDS-PAGE Standard</th> <th rowspan="2">Standard Catalog #</th> </tr> <tr> <th>9 Well</th> <th>10 Well</th> <th>10 Well/50 µl</th> <th>12 Well</th> <th>15 Well</th> </tr> </thead> <tbody> <tr> <td>30-150 kDa</td> <td>10% with Gelatin</td> <td>161-1203</td> <td>161-1113</td> <td>161-1167</td> <td>161-1185</td> <td>161-1131</td> <td>Prestained, low range</td> <td>161-0315</td> </tr> <tr> <td>20-120 kDa</td> <td>12% with β-Casein</td> <td>161-1204</td> <td>161-1114</td> <td>161-1168</td> <td>161-1186</td> <td>161-1132</td> <td>Prestained, broad range</td> <td>161-0318</td> </tr> <tr> <td></td> <td></td> <td></td> <td></td> <td></td> <td></td> <td></td> <td>Prestained, high range</td> <td>161-0309</td> </tr> </tbody> </table>	MW Range Resolved	Zymogram Gel %T With Contents	Single gel catalog numbers - order as many as you need!					Suggested SDS-PAGE Standard	Standard Catalog #	9 Well	10 Well	10 Well/50 µl	12 Well	15 Well	30-150 kDa	10% with Gelatin	161-1203	161-1113	161-1167	161-1185	161-1131	Prestained, low range	161-0315	20-120 kDa	12% with β-Casein	161-1204	161-1114	161-1168	161-1186	161-1132	Prestained, broad range	161-0318								Prestained, high range	161-0309	
MW Range Resolved	Zymogram Gel %T With Contents			Single gel catalog numbers - order as many as you need!							Suggested SDS-PAGE Standard	Standard Catalog #																															
		9 Well	10 Well	10 Well/50 µl	12 Well	15 Well																																					
30-150 kDa	10% with Gelatin	161-1203	161-1113	161-1167	161-1185	161-1131	Prestained, low range	161-0315																																			
20-120 kDa	12% with β-Casein	161-1204	161-1114	161-1168	161-1186	161-1132	Prestained, broad range	161-0318																																			
							Prestained, high range	161-0309																																			
 Running Buffer	<table border="1"> <thead> <tr> <th></th> <th></th> <th>Catalog #</th> </tr> </thead> <tbody> <tr> <td rowspan="3">10x Tris/Glycine/SDS Buffer</td> <td>1 L</td> <td>161-0733</td> </tr> <tr> <td>5 L NEW!</td> <td>161-0770</td> </tr> <tr> <td>6 x 1 L</td> <td>161-0756</td> </tr> </tbody> </table>			Catalog #	10x Tris/Glycine/SDS Buffer	1 L	161-0733	5 L NEW!	161-0770	6 x 1 L	161-0756																																
		Catalog #																																									
10x Tris/Glycine/SDS Buffer	1 L	161-0733																																									
	5 L NEW!	161-0770																																									
	6 x 1 L	161-0756																																									
 Electrophoresis Cell and Power Supply	<table border="1"> <tbody> <tr> <td>Mini-PROTEAN 3 Electrophoresis Module</td> <td>165-3302</td> </tr> <tr> <td>PowerPac 300 (100/120 V)</td> <td>165-5050</td> </tr> <tr> <td>PowerPac 300 (220/240 V)</td> <td>165-5051</td> </tr> </tbody> </table> <p>Run conditions per 2 gels: Begin - 100 V (C), 25 mA, 80-100 minutes End - 100 V (C), 12 mA</p>	Mini-PROTEAN 3 Electrophoresis Module	165-3302	PowerPac 300 (100/120 V)	165-5050	PowerPac 300 (220/240 V)	165-5051																																				
Mini-PROTEAN 3 Electrophoresis Module	165-3302																																										
PowerPac 300 (100/120 V)	165-5050																																										
PowerPac 300 (220/240 V)	165-5051																																										
 Renaturation	Zymogram Renaturation Buffer, 10x, 125 ml	161-0765																																									
 Development	Zymogram Development Buffer, 10x, 125 ml	161-0766																																									
 Gel Staining	<p>Note: Document gels prior to staining since stain will obscure standards.</p> <table border="1"> <tbody> <tr> <td>Coomassie R-250, 10 g</td> <td>161-0400</td> </tr> </tbody> </table> <p>(For Zymogram Staining, use as a solution of: 0.5% Coomassie R250, 10% Acetic Acid, 40% methanol)</p>	Coomassie R-250, 10 g	161-0400																																								
Coomassie R-250, 10 g	161-0400																																										



See page 40 of the Ready Gel Resource Guide for more on gel drying equipment





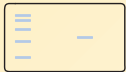
OLIGOSACCHARIDE ANALYSIS

Step	Ready Gel System Product / Catalog #		
 Sample Preparation Kit Selection	N-Linked Oligosaccharide Analysis Glycoprotein in non-Tris buffer Profiling Kit / 170-6501 Sequencing Kit / 170-6510	O-Linked Oligosaccharide Analysis Glycoprotein in non-Tris buffer Profiling Kit / 170-6815	Monosaccharide Analysis Glycoprotein in non-Tris buffer Composition Kit / 170-6811
 Running Buffer	Oligosaccharide Buffer Pack/ 170-6503	Oligosaccharide Buffer Pack/ 170-6503	Monosaccharide Buffer Pack/ 170-6814
 Electrophoresis Cell and Power Supply	Mini-PROTEAN 3 Electrophoresis Module / 165-3302 Carbohydrate Gaskets 165-3157 PowerPac 300 (100/120 V) 165-5050 PowerPac 300 (220/240 V) 165-5051 Run conditions per 2 gels: Begin: 100-200 V, 30 mA (C), 60-90 minutes End: 200-300 V, 30 mA (C)	Mini-PROTEAN 3 Electrophoresis Module / 165-3302 Carbohydrate Gaskets 165-3157 PowerPac 300 (100/120 V) 165-5050 PowerPac 300 (220/240 V) 165-5051 Run conditions per 2 gels: Begin: 200-300 V, 40 mA (C), 50-60 minutes End: 600-700 V, 40 mA (C)	Mini-PROTEAN 3 Electrophoresis Module / 165-3302 Carbohydrate Gaskets 165-3157 PowerPac 300 (100/120 V) 165-5050 PowerPac 300 (220/240 V) 165-5051 Run conditions per 2 gels: Begin: 300-400 V, 60 mA (C), 50-60 minutes End: 500-600 V, 60 mA (C)
 Detection Method Imaging Options	ANTS Fluorophore Gel Doc 2000 or mini transilluminator	ANTS Fluorophore Gel Doc 2000 or mini transilluminator	AMAC Fluorophore Gel Doc 2000 or mini transilluminator



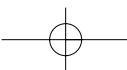
Visit www.discover.bio-rad.com for information on the new Mini-PROTEAN 3 Electrophoresis System

GLYCOPROTEIN ANALYSIS

Step	Ready Gel System Product	Catalog #								
 Sample Preparation	Laemmli Sample Buffer, 30 ml	161-0737								
	2-mercaptoethanol, 25 ml	161-0710								
 Ready Gel and Standard Selection	Single gel catalog numbers - order as many as you need!						Suggested Standard	Standard Catalog #		
	MW Range Resolved	Tris-HCl Gel %T	9 Well	10 Well	10 Well/50 µl	12 Well	15 Well	2-D/Prep		
	100-250 kDa	5%	161-1215	161-1210	161-1213	161-1214	161-1211	161-1212	High range	161-0303
	40-200 kDa	7.5%	161-1190	161-1100	161-1154	161-1172	161-1118	161-1136	High range	161-0303
	30-150 kDa	10%	161-1191	161-1101	161-1155	161-1173	161-1119	161-1137	Broad range	161-0317
	20-120 kDa	12%	161-1192	161-1102	161-1156	161-1174	161-1120	161-1138	Low range	161-0304
	10-100 kDa	15%	161-1193	161-1103	161-1157	161-1175	161-1121	161-1139	Low range	161-0304
	6-50 kDa	18%	161-1221	161-1216	161-1219	161-1220	161-1217	161-1218	Low range	161-0304
	20-250 kDa	4-15%	161-1194	161-1104	161-1158	161-1176	161-1122	161-1140	Broad range	161-0317
	10-200 kDa	4-20%	161-1195	161-1105	161-1159	161-1177	161-1123	161-1141	Broad range	161-0317
	6-70 kDa	8-16%	161-1227	161-1222	161-1225	161-1226	161-1223	161-1224	Broad range	161-0317
	10-100 kDa	10-20%	161-1196	161-1106	161-1160	161-1178	161-1124	161-1142	Broad range	161-0317
	 Running Buffer	10x Tris-Glycine/SDS Buffer	1 L	161-0732						
10x Tris-Glycine/SDS Buffer		5 L NEW!	161-0772							
Mini-PROTEAN 3 Electrophoresis Module			165-3302							
PowerPac 300 (100/120 V)			165-5050							
 Electrophoresis Cell and Power Supply	PowerPac 300 (220/240 V)		165-5051							
	Run conditions per 2 gels: Begin - 200 V (C), 100 mA End - 200 V (C), 60 mA, 35 minutes									
 Detection Method	Glycoprotein Detection Colorimetric Western Blot (BCIP/NBT)		170-6490							
	Deglycosylation Detection Coomassie R-250 Staining Solutions Kit		170-6500 170-0435							



See page 38 of the Ready Gel Resource Guide for more on power supplies



BLOT PROCESSING BUFFERS

Sample	Method of Transfer	Buffer Stock	Volume	Catalog #
Proteins	Trans-Blot/Mini Trans-Blot	10x Tris/Glycine	1 L	161-0734
			5 L NEW!	161-0771
			6 x 1 L	161-0757
Proteins	Trans-Blot Semi-Dry	10x Tris/CAPS (requires SDS)	1 L	161-0778
			250 ml	161-0416
DNA/RNA	Trans-Blot/Mini Trans-Blot	10x TBE	1 L	161-0733
			5 L NEW!	161-0770
			6 x 1 L	161-0756
		50x TAE	1 L	161-0743
			5 L NEW!	161-0773
6 x 1 L	161-0759			
DNA/RNA	Trans-Blot Semi-Dry	10x TBE	1 L	161-0733
			5 L NEW!	161-0770
			6 x 1 L	161-0756
DNA/RNA	Capillary transfer	20x SSC	1 L NEW!	161-0774
			5 L NEW!	161-0775

Item	Product	Package Size	Catalog #
Wash buffers	10x TBS 10x PBS 20x SSC	1 L NEW!	170-6435
		1 L NEW!	161-0780
		1 L NEW!	161-0774
		5 L NEW!	161-0775
Blocking Buffers	1x TBS/1% Casein 1x PBS/1% Casein 10% Tween	1 L NEW!	161-0782
		1 L NEW!	161-0783
		1 L NEW!	161-0781

Ready Gels Accelerate Your Research!



Bio-Rad
Laboratories

**Life Science
Group**

Website www.bio-rad.com Bio-Rad Laboratories Main Office 2000 Alfred Nobel Drive, Hercules, CA 94547, Ph. (510) 741-1000, Fx. (510)741-5800
 Also in: Australia Ph. 02 9914 2800, Fx. 02 9914 2889 Austria Ph. (01) 877 89 01, Fx. (01) 876 56 29 Belgium Ph. 09-385 55 11, Fx. 09-385 65 54
 Canada Ph. (905) 712-2771, Fx. (905) 712-2990 China Ph. 86-10-62051850/51, Fx. 86-10-62051876 Denmark Ph. 45 39 17 99 47, Fx. 45 39 27 16 98
 Finland Ph. 358 (0)9 804 2200, Fx. 358 (0)9 804 1100 France Ph. 01 43 90 46 90, Fx. 01 46 71 24 67 Germany Ph. 089 318 84-0, Fx. 089 318 84-100
 Hong Kong Ph. 852-2789-3300, Fx. 852-2789-1257 India Ph. (91-11) 461-0103, Fx. (91-11) 461-0765 Israel Ph. 03 951 4127, Fx. 03 951 4129
 Italy Ph. 39-02-216091, Fx.39-02-21609-399 Japan Ph. 03-5811-6270, Fx. 03-5811-6272 Korea Ph. 82-2-3473-4460, Fx. 82-2-3472-7003
 Latin America Ph. 305-894-5950, Fx. 305-894-5960 Mexico Ph. 514-2210, Fx. 514-2209 The Netherlands Ph. 0318-540666, Fx. 0318-542216
 New Zealand Ph. 64-9-4152280, Fx. 64-9-4152284 Russia Ph. 7-095-4585822, Fx. 7-095-4585782 Singapore Ph. 65-2729877, Fx. 65-2734835
 Spain Ph. 34-91-661-7085, Fx. 34-91-661-9698 Sweden Ph. 46 (0)8 627 50 00, Fx. 46 (0)8 627 54 00 Switzerland Ph. 01-809 55 55, Fx. 01-809 55 00
 United Kingdom Ph. 0800-181134, Fx. 01442-259118

