



Pure Convenience

Acrylamide
Premixed Acrylamide/Bis Powders
Premixed Acrylamide/Bis Solutions
Premixed Buffers

BIO-RAD

Convenience Reagents

Bio-Rad's full line of convenience reagents allows you to obtain reproducible electrophoresis results faster while maintaining the highest standards of purity.

Choose from a variety of reagents tailored to any application or budget. Contact your local sales office for pricing information.

Table 1. Polyacrylamide Gel Composition and Separation Ranges

Resolving Gel or Gradient %T	Approximate Separation Range
SDS-PAGE	
7.5%	45,000 to 200,000 daltons
10%	25,000 to 200,000 daltons
12%	14,400 to 120,000 daltons
15%	6,500 to 100,000 daltons
4–15%	20,000 to 200,000 daltons
4–20%	6,500 to 200,000 daltons
10–20%	6,500 to 120,000 daltons
TBE	
5%	200 to 2,000 base pairs
10%	50 to 1,750 base pairs
Tris-Tricine	
16.5%	2,000 to 70,000 daltons
10–20%	2,000 to 120,000 daltons

Table 2. How Many Gels Can I Run?

Approximate number of 12% mini format gels that can be run with acrylamide solutions, premixed powders, and premixed buffers:

Product Description	Number of 12% Mini Gels
500 ml 40% Acrylamide/Bis Solution	300
500 ml 30% Acrylamide/Bis Solution	225
150 g Premixed Acrylamide/Bis Powder	225
30 g Premixed Acrylamide/Bis Powder	45
100 g Acrylamide Powder	150
500 g Acrylamide Powder	750
1 liter of 10x Premixed Buffer	66

Table 3. Premixed Buffer Composition
(All buffers are made with 18 megaohm water)

Description	1x Buffer Concentration						
	Tris	Glycine	Tricine	SDS	Boric Acid	Acetic Acid	EDTA
10x Tris/Glycine/SDS	25 mM	192 mM		0.1%			
10x Tris/Glycine	25 mM	192 mM					
10x Tris/Tricine/SDS	100 mM		100 mM	0.1%			
10x Tris/Boric Acid/EDTA	89 mM				89 mM		2 mM
10x TBE Extended Range*	130 mM				45 mM		2.5 mM
50x Tris/Acetic Acid/EDTA	40 mM					20 mM	1 mM

Description	Tris-HCl, pH 6.8	SDS	Glycerol	Bromophenol Blue	Coomassie® G-250 Dye
Laemmli Sample Buffer	62.5 mM	2%	25%	0.01%	
Native Sample Buffer	62.5 mM		40%	0.01%	
Tricine Sample Buffer	200 mM	2%	40%		0.04%

* Extended buffer capacity for longer DNA sequencing runs.

Ordering Information

Catalog Number	Product Description	Quantity
Acrylamide Powder		
161-0100	Acrylamide, 99.9%	100 g
161-0101		500 g
161-0107		1 kg
161-0103		2 kg
161-0108		5 kg
Premixed Acrylamide/Bis Powders		
161-0120	Acrylamide/Bis Powder, 19:1 mixture, (5% C)	30 g
161-0123		150 g
161-0121	Acrylamide/Bis Powder, 29:1 mixture, (3.3% C)	30 g
161-0124		150 g
161-0122	Acrylamide/Bis Powder, 37.5:1 mixture, (2.67% C)	30 g
161-0125		150 g
Acrylamide/Bis Solutions		
161-0154	30% Acrylamide/Bis Solution, 19:1 mixture, (5% C)	500 ml
161-0156	30% Acrylamide/Bis Solution, 29:1 mixture, (3.3% C)	500 ml
161-0158	30% Acrylamide/Bis Solution, 37.5:1 mixture, (2.6% C)	500 ml
161-0144	40% Acrylamide/Bis Solution, 19:1 mixture, (5% C)	500 ml
161-0146	40% Acrylamide/Bis Solution, 29:1 mixture, (3.3% C)	500 ml
161-0148	40% Acrylamide/Bis Solution, 37.5:1 mixture, (2.6% C)	500 ml
161-0140	40% Acrylamide Solution	500 ml
161-0142	2% Bis Solution	500 ml

Ordering Information

Catalog Number	Product Description	Quantity
Catalysts		
161-0501	Riboflavin-5'-Phosphate	10 g
161-0700	Ammonium Persulfate	10 g
161-0800	TEMED	5 ml
Premixed Buffers - see Table 3 for exact formulations		
161-0732	10x Tris/Glycine/SDS	1 L
161-0755		6 x 1 L
161-0734	10x Tris/Glycine	1 L
161-0757		6 x 1 L
161-0744	10x Tris/Tricine/SDS	1 L
161-0760		6 x 1 L
161-0733	10x Tris/Boric Acid/EDTA	1 L
161-0756		6 x 1 L
161-0741	10x TBE Extended Range	1 L
161-0758		6 x 1 L
161-0743	50x Tris/Acetic Acid/EDTA	1 L
161-0759		6 x 1 L
Premixed Sample Buffers		
161-0737	Laemmli Sample Buffer	30 ml
161-0738	Native Sample Buffer	30 ml
161-0739	Tricine Sample Buffer	30 ml
Detergents		
161-0300	SDS (Sodium dodecyl sulfate)	25 g
161-0301		100 g
161-0302		1 kg
161-0416	SDS Solution 10% (w/v)	250 ml
161-0418	SDS Solution 20% (w/v)	1,000 ml
161-0407	Triton X-100 Detergent	500 ml
170-6531	Tween 20 EIA Grade	100 ml
161-0460	CHAPS	1 g
161-0465	CHAPSO	1 g

Coomassie is a trademark of I C I Organics, Inc.



**Bio-Rad
Laboratories**

1-800-541-9913
www.bio-rad.com

**Life Science
Group**

U.S. (800) 424-0700 Australia (02-94542300) Austria (51-3778903)
 Belgium (04-38855555) Canada (905) 752-2775 China (86-30) 2048622
 Denmark (89) 370047 Finland (00) 3042200 France (6) 48 00 48 00
 Germany (089) 8884-0 Hong Kong (852) 2308300 India (01) 55-065-0508
 Israel (03) 055 4327 Italy (02) 25800.3 Japan (03-28) 55-6270
 The Netherlands (0338) 38-290888 New Zealand (09-448) 8000
 Singapore (65) 222-4877 Spain (91) 685 70 85
 Sweden (08) 038 62 750 00 Switzerland (01) 800 55 55
 United Kingdom (0800) 385384