

Sample Preparation Buffers

1 M Tris-HCl, pH 7.6 (100 ml)

Tris base	12.11 g
Deionized H ₂ O (diH ₂ O)	80 ml
Adjust pH to 7.6 with HCl	
diH ₂ O	to 100 ml

0.5 M Tris-HCl, pH 6.8 (100 ml)

(catalog #161-0799)

Tris base	6.06 g
diH ₂ O	~60 ml
Adjust to pH 6.8 with HCl	
diH ₂ O	to 100 ml
Store at 4°C	

10% SDS (10 ml)

(catalog #161-0416)

SDS	1.00 g
diH ₂ O	to 10 ml

1.0% Bromophenol Blue (10 ml)

(catalog #161-0404)

Bromophenol blue	100 mg
diH ₂ O	to 10 ml

RIPA Solubilization Buffer (100 ml)

25 mM Tris-HCl pH 7.6, 150 mM NaCl, 5 mM EDTA, 1% NP-40 or 1% Triton X-100, 1% sodium deoxycholate, 0.1% SDS

NaCl	0.88 g
EDTA	0.15 g
NP-40 or Triton X-100	1 g
Sodium deoxycholate	1 g
SDS	0.10 g
diH ₂ O	80 ml
1 M Tris-HCl, pH 7.6	2.5 ml
diH ₂ O	to 100 ml

Phosphate Buffered Saline (PBS, 1 L)

0.9% (w/v) sodium chloride in 10 mM phosphate buffer, pH 7.4

NaCl	8.00 g
KCl	0.20 g
Na ₂ HPO ₄	1.44 g
KH ₂ PO ₄	0.24 g
diH ₂ O	800 ml
Adjust pH to 7.4 with HCl or NaOH	
diH ₂ O	to 1 L

SDS-PAGE Sample Buffer (2x, 8 ml)

(catalog #161-0737, 30 ml)

62.5 mM Tris-HCl, pH 6.8, 25% glycerol, 2% SDS, 0.01% bromophenol blue

0.5 mM Tris-HCl, pH 6.8	1.0 ml
25% Glycerol	2.0 ml
1.0% Bromophenol blue	0.08 ml
10% SDS	1.6 ml
diH ₂ O	2.92 ml

Store as 1–2 ml aliquots at –70°C and add β-mercaptoethanol (0.4 ml) or 3% DTT immediately before use.

Protein Precipitation Solution (100 ml)

20% Trichloroacetic acid (TCA), 0.2% DTT in ice-cold acetone (–20°C)

TCA	20.00 g
DTT	0.20 g
Acetone	80 ml
Dissolve	
Acetone	to 100 ml

Store at –20°C

Wash Solution (100 ml)

0.2% DTT in ice-cold acetone (–20°C)

DTT	0.20 g
Acetone	80 ml
Dissolve	
Acetone	to 100 ml

Store at –20°C

Lysis Buffer (50 ml)

2 M thiourea, 7 M urea, 4% (w/v) CHAPS, 1% (w/v) DTT, 2% (v/v) carrier ampholytes (pH 3–10)

Urea	21.00 g
Thiourea	7.60 g
diH ₂ O	to 45 ml
CHAPS	2.00 g
Bio-Lyte® ampholytes	1.0 ml
DTT	0.50 g
diH ₂ O	to 50 ml

SDS Sample Solubilization Buffer (50 ml)

1% (w/v) SDS, 100 mM Tris-HCl (pH 9.5)

SDS	0.50 g
Tris base	0.60 g
diH ₂ O	40 ml
Titrate to pH 9.5 with diluted HCl	
diH ₂ O	to 50 ml

Gel Casting Reagents

Acrylamide/Bis (30% T, 2.67% C)

Acrylamide (29.2 g/100 ml)	87.60 g
N'N'-bis-methylene-acrylamide	2.40 g
diH ₂ O	to 300 ml

Filter and store at 4°C in the dark (30 days).

Premade alternatives:

Catalog #161-0125, 37.5:1 acrylamide/bis powder

Catalog #161-0158, 30% acrylamide/bis solution

1.5 M Tris-HCl, pH 8.8 (150 ml)

(catalog #161-0798, 1 L)

Tris base (18.15 g/100 ml)	27.23 g
diH ₂ O	80 ml
Adjust to pH 8.8 with 6 N HCl	
diH ₂ O	to 150 ml

Store at 4°C

0.5 M Tris-HCl, pH 6.8

(catalog #161-0799, 1 L)

Tris base	6.00 g
diH ₂ O	60 ml
Adjust to pH 6.87 with 6 N HCl	
diH ₂ O	to 100 ml

Store at 4°C

10% (w/v) SDS (100 ml)

(catalog #161-0416)

SDS	10.00 g
diH ₂ O	90 ml
Dissolve with gentle stirring	
diH ₂ O	to 100 ml

10% (w/v) APS (fresh daily)

(catalog #161-0416)

Ammonium persulfate	0.10 g
diH ₂ O	1 ml

Water-Saturated *n*-Butanol

<i>n</i> -Butanol	50 ml
diH ₂ O	5 ml

Combine in a bottle and shake. Use the top phase only. Store at room temperature.

Sample Buffers

2× SDS-PAGE (Laemmli, 30 ml)

(catalog #161-0737)

62.5 mM Tris-HCl, pH 6.8, 2% SDS, 25% glycerol, 0.01% bromophenol blue, 5% β-mercaptoethanol (added fresh)

0.5 M Tris-HCl, pH 6.8	3.75 ml
50% Glycerol	15.0 ml
1.0% Bromophenol blue	0.3 ml
10% SDS	6.0 ml
diH ₂ O	to 30 ml

β-mercaptoethanol (50 μl to 950 μl sample buffer) before use.

2× Native PAGE (30 ml)

(catalog #161-0738)

62.5 mM Tris-HCl, pH 6.8, 40% glycerol, 0.01% bromophenol blue

0.5 M Tris-HCl, pH 6.8	3.75 ml
50% Glycerol	24.0 ml
1.0% Bromophenol blue	0.3 ml
diH ₂ O	to 30 ml

2× Tricine (30 ml)

(catalog #161-0739)

200 mM Tris-HCl, pH 6.8, 2% SDS, 40% glycerol, 0.04% Coomassie Brilliant Blue G-250, 2% β-mercaptoethanol (added fresh)

0.5 M Tris-HCl, pH 6.8	12.0 ml
50% Glycerol	24.0 ml
10% SDS	6.0 ml
Coomassie Blue G-250	12.00 mg
diH ₂ O	to 30 ml

β-mercaptoethanol (20 μl to 980 μl sample buffer) before use.

Running Buffers

10× SDS-PAGE (1 L)

(catalog #161-0732)

250 mM Tris, 1.92 M glycine, 1% SDS, pH 8.3	
Tris base	30.30 g
Glycine	144.10 g
SDS	10.00 g
diH ₂ O	to 1 L

Do not adjust the pH (~pH 8.3).

10× Native PAGE (1 L)

(catalog #161-0734)

250 mM Tris, 1.92 M glycine, pH 8.3	
Tris base	30.30 g
Glycine	144.10 g
diH ₂ O	to 1 L

Do not adjust the pH (~pH 8.3).

10× Tris-Tricine (1 L)

(catalog #161-0744)

1 M Tris, 1 M Tricine, 1% SDS, pH 8.3	
Tris base	121.10 g
Tricine	179.20 g
SDS	10.00 g
diH ₂ O	to 1 L

Do not adjust the pH (~pH 8.3).

Buffer Components

0.5 M Tris-HCl, pH 6.8 (1 L)

(catalog #161-0799)

Tris base	60.60 g
diH ₂ O	~900 ml
Adjust to pH 6.8 with HCl	
diH ₂ O	to 1 L

Store at 4°C

10% SDS (250 ml)

(catalog #161-0416)

SDS	25.00 g
diH ₂ O	to 250 ml

1.0% Bromophenol Blue (10 ml)

(catalog #161-0404)

Bromophenol blue	100 mg
diH ₂ O	to 10 ml

This is an excerpt from Bio-Rad's comprehensive Electrophoresis Guide (Bulletin 6640).



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 11 5044 5699
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00
France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666
New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04
Singapore 65 6415 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55
Taiwan 886 2 2578 7189 Thailand 800 88 22 88 United Kingdom 020 8328 2000