

# electrophoresis

## Mini-PROTEAN® Empty Cassette Gel Casting Instructions

### Instructions

1. Prepare the resolving gel acrylamide solution by combining all reagents except ammonium persulfate (APS) and TEMED. Degas the solution under a vacuum for at least 15 min.
2. The Mini-PROTEAN cassette should be held vertically for gel casting. The AnyGel™ stand (catalog #165-4131) or a test tube rack works well to stabilize the cassette during casting.
3. Add APS and TEMED to the degassed acrylamide solution and pour the solution into the cassette. Pour the solution smoothly to prevent it from mixing with air. Fill the cassette until the solution is 1 cm below the teeth of the comb.
4. Immediately overlay the acrylamide solution with distilled water. Apply the overlay slowly, with a steady, even rate of delivery to prevent mixing. DO NOT use alcohol overlay solutions, as these may discolor the cassette.
5. Allow the gel to polymerize for 45 min to 1 hr. Rinse off the overlay solution with distilled water.
6. Prepare the stacking gel acrylamide solution. Combine all reagents except APS and TEMED and degas under vacuum for at least 15 min.
7. Dry the area above the separating gel with filter paper before pouring the stacking gel.
8. Place a comb in the cassette and tilt it so that the teeth are at a slight (~10°) angle. This will prevent air from being trapped under the comb teeth while the acrylamide solution is being poured.
9. Add APS and TEMED to the degassed acrylamide solution and pour the solution down the edge of the cassette nearest to the upturned side of the comb. Pour until all the teeth have been covered by solution. Then properly align the comb in the cassette and add solution to fill completely.
10. Allow the gel to polymerize 30–45 min.

### Stock Solutions

#### 30% Acrylamide/bis (37.5:1)

- 29.2 g acrylamide
- 0.8 g N,N'-bis-methylene-acrylamide

Make to 100 ml with deionized water. Filter and store at 4°C in the dark.

Or, use

30% Acrylamide/Bis Solution, 37.5:1, 500 ml (catalog #161-0158)  
Preweighed Acrylamide/Bis Powder, 37.5:1, 30 g (catalog #161-0122)

#### 1.5 M Tris-HCl, pH 8.8

- 18.15 g Tris base
- ~60 ml deionized water

Adjust to pH 8.8 with 6 N HCl.

Make to 100 ml with deionized water and store at 4°C in the dark.

Or, use

1.5 M Tris-HCl, pH 8.8, 1 L (catalog #161-0798)

#### 0.5 M Tris-HCl, pH 6.8

- 6 g Tris base
- ~60 ml deionized water

Adjust to pH 6.8 with 6 N HCl.

Make to 100 ml with deionized water and store at 4°C in the dark.

Or, use

0.5 M Tris-HCl, pH 6.8, 1 L (catalog #161-0799)

#### 10% (w/v) SDS

Dissolve 10 g SDS in 90 ml water and bring to 100 ml with deionized water.

Or, use

SDS Solution, 10% (w/v), 250 ml (catalog #161-0416)

	Stacking Gel		Resolving Gel	
	4%	7.5%	12%	X%
30% Acrylamide/bis	1.32 ml	2.5 ml	4.0 ml	0.33•X ml
0.5 M Tris-HCl, pH 6.8	2.52 ml	—	—	—
1.5 M Tris-HCl, pH 8.8	—	2.5 ml	2.5 ml	2.5 ml
10% SDS	100 µl	100 µl	100 µl	100 µl
Distilled deionized water	6 ml	4.85 ml	3.35 ml	7.35 – (0.33•X) ml
TEMED	10 µl	5 µl	5 µl	5 µl
10% APS	50 µl	50 µl	50 µl	50 µl
<b>Total Volume</b>	<b>10 ml</b>	<b>10 ml</b>	<b>10 ml</b>	<b>10 ml</b>



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

*Life Science  
Group*

**Web site** [www.bio-rad.com](http://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