



## Product Information

### Acrylamide/Bis-acrylamide, 30% solution (Mix Ratio 29:1)

Catalog Number **A3574**  
Storage Temperature 2–8 °C

#### Product Description

The Acrylamide/Bis-acrylamide solution is used in protein and nucleic acid electrophoresis. The solution concentration (30%) is based on the total weight of both the acrylamide and bis-acrylamide. The mix or feed ratio (w/w) of acrylamide:bis-acrylamide is 29:1. The solution is prepared from electrophoresis grade acrylamide and bis-acrylamide in ultrapure water. The product is passed through a 0.2 µm filter.

The product is suitable for electrophoresis.

#### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

#### Storage/Stability

Store the product at 2–8 °C.

#### Procedure

The following formula may be used to calculate the volume of stock solution required for a gel concentration of <30%:

$$V_r = \frac{c \times V_t}{30\%}$$

Where:

**V<sub>r</sub>** is the required volume of Acrylamide/Bis-acrylamide, 30% Solution

**c** is the final desired acrylamide concentration (%)

**V<sub>t</sub>** is the total volume of final solution to be prepared

To prepare gels for use in the Laemmli system, the following solutions may be combined as indicated in Tables 1 and 2.

Solution A - Acrylamide/Bis-acrylamide 30% Solution, Catalog Number A3574

Solution B - Combine:

	Catalog No.	Amount
Trizma <sup>®</sup> Base	T6066	7.28 g
TEMED	T9281	92 µl

Adjust the pH of the solution to 8.9 with 1 M HCl and add ultrapure water to 40 ml final volume.

Solution C - Combine:

Trizma Base	T6066	1.21 g
TEMED	T9281	92 µl

Adjust the pH of the solution to 6.9 with 1 M HCl and add ultrapure water to 20 ml final volume.

Solution D - Dissolve 1 g of Sodium Dodecyl Sulfate (Catalog Number L3771) in 10 ml of ultrapure water and filter.

Solution E - Dissolve 60 mg of Ammonium Persulfate (Catalog Number A3678) in 5 ml of ultrapure water. Prepare fresh each day.

See Tables 1 and 2 for mixing instructions.

**Table 1.**

Separating Gel Solution (Final volume = 20 ml)

Component	Final Acrylamide Concentration									
	5%	6%	7%	8%	9%	10%	11%	12%	15%	20%
Ultrapure Water (ml)	10.47	9.80	9.13	8.47	7.80	7.13	6.47	5.80	3.80	0.47
Solution A (ml)	3.33	4.00	4.67	5.33	6.00	6.67	7.33	8.00	10.00	13.33
Solution B (ml)	5.0	5.0	5.0	5.0	5.0	5.0	5.0	5.0	5.0	5.0
Solution D (ml)	0.2	0.2	0.2	0.2	0.2	0.2	0.2	0.2	0.2	0.2
Solution E (ml)	1.0	1.0	1.0	1.0	1.0	1.0	1.0	1.0	1.0	1.0

**Table 2.**

Stacking Gel Solution (Final volume = 8 ml)

Component	Final Acrylamide Concentration
	3%
Ultrapure Water (ml)	5.52
Solution A (ml)	0.80
Solution C (ml)	1.00
Solution D (ml)	0.08
Solution E (ml)	0.60

**References**

1. Laemmli, U.K., Nature, **227**, 680 (1970).

Trizma is a registered trademark of Sigma-Aldrich™  
Biotechnology LP and Sigma-Aldrich Co.

KR,MAM 09/07-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.