

Acrylamide/Bis 19:1 40% (w/v) Solution

Store at 4°C.

Catalog #:	AM9022	AM9024
Amount:	500 mL	2 x 500 mL
Product Description:	A 40% (w/v) solution of Ultrapure Acrylamide (38%) and bis-Acrylamide (2%)	
Appearance:	Clear liquid	
Molecular Weight:	Acrylamide component – 71.08 N,N' Methylene-bis-Acrylamide component – 154.17	
Caution:	Poison. Acrylamide is a neurotoxin and a suspected carcinogen. Avoid contact with eyes, skin, and clothing. Wash thoroughly with soap and water after handling. See MSDS.	
Storage Conditions:	Store at 4°C. Keep tightly sealed. If high resolution is required (such as sequencing), fresh solution (less than 2 months) should be used.	

USER INFORMATION

General Information: Ambion® Acrylamide/Bis Solution is provided in a ready-to-use form, eliminating the dust and inhalation hazard associated with weighing and preparing acrylamide and bis-acrylamide powders and solutions. Acrylamide/Bis 19:1 is a 40% (w/v) solution of acrylamide (38%) and bis-acrylamide (2%). Ideal for use in ribonuclease protection assays, general small nucleic acid electrophoresis (such as gel purification of RNA probes, PCR product analysis), or sequencing applications.

QUALITY CONTROL

Quality Verification: Specifications listed in this document have been verified to pass all criteria.

OTHER INFORMATION

Material Safety Data Sheets: Material Safety Data Sheets (MSDSs) for any chemical product supplied by Applied Biosystems or Ambion are available 24 hours a day. At www.appliedbiosystems.com, select Support, then MSDS. Search by chemical name, product name, product part number, or MSDS part number. Right-click to print or download the MSDS of interest. At www.ambion.com, go to the web catalog page for the product of interest. Select MSDS, then right-click to print or download. Or, e-mail (MSDS_Inquiry_CCRM@appliedbiosystems.com), telephone (650-554-2756; USA), or fax (650-554-2252; USA) your request, specifying the catalog or part number(s) and the name of the product(s). We will e-mail the associated MSDSs unless you request fax or postal delivery. Requests for postal delivery require 1-2 weeks for processing.

Warranty and Liability: *For research use only. Not for use in diagnostic procedures.*

Ambion, a part of Life Technologies Corporation, is committed to delivering superior product quality and performance, supported by industry-leading global service and technical support teams. Warranty information for the accompanying consumable product is available at www.ambion.com/info/warranty in "Limited Warranty for Consumables," which is subject to the exclusions, conditions, exceptions, and limitations set forth under the caption "EXCLUSIONS, CONDITIONS, EXCEPTIONS, AND LIMITATIONS" in the full warranty statement. Please contact Ambion if you have any questions about our warranties or would like information about post-warranty support.

Information in this document is subject to change without notice. Ambion, a part of Life Technologies Corporation, assumes no responsibility for any errors that may appear in this document.

Ambion disclaims all warranties with respect to this document, expressed or implied, including but not limited to those of merchantability or fitness for a particular purpose. In no event shall Ambion be liable, whether in contract, tort, warranty, or under any statute or on any other basis for special, incidental, indirect, punitive, multiple or consequential damages in connection with or arising from this document, including but not limited to the use thereof.

Trademarks, Patents, and Licensing:

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

© 2009 Life Technologies Corporation. All rights reserved. 4386625B