

5M Ammonium Acetate

Store at room temperature or -20°C .

Catalog # (P/N):	AM9070G	AM9071	AM9077G
Amount:	100 mL	500 mL	1 mL
Product Description:	A nuclease-free solution for routine precipitation of nucleic acids.		
Appearance:	Colorless liquid		
Storage Conditions:	Store at room temperature or -20°C .		
Safety Information:	Warning! Causes eye, skin, and respiratory tract irritation. May be harmful if swallowed, inhaled, or absorbed through skin. Avoid breathing vapor. Use with adequate ventilation. Avoid contact with eyes and skin. Wear appropriate protective eyewear, clothing, and gloves. Read Material Safety Data Sheet.		

USER INFORMATION

General Information:	Ammonium acetate is used for routine precipitation of nucleic acids, and is especially useful for reducing the coprecipitation of unwanted dNTPs and contaminating oligosaccharides. This solution is ready for use and requires no extra preparation. Note: Do not autoclave Ammonium Acetate Buffer. If precipitates form, warm solution to 37°C to resuspend.
Applications:	Routine Precipitation of DNA or RNA at Concentrations ≥ 20 ng/mL <ol style="list-style-type: none"> Use 5 M Ammonium Acetate to bring the final concentration of ammonium acetate in the solution to 2.5 M. Add 2 volumes of 100% ethanol for DNA or 2.5 volumes of 100% ethanol for RNA. Chill the solution at or below -20°C for ≥ 30 min. Centrifuge at top speed in a microcentrifuge for 15 min. Discard the supernatant and wash the pellet with ice cold 70–75% ethanol to remove residual salt. Resuspend in nuclease-free water (P/N AM9937) or 10 mM Tris, pH 7–8, 1 mM EDTA. <p>Note: Ammonium acetate should not be used when the nucleic acid will be phosphorylated using T4 polynucleotide kinase, because this enzyme is inhibited by ammonium ions.</p>

QUALITY CONTROL

	5 M Ammonium Acetate is tested in the following quality control assays before and after final packaging. The product is filtered through a $0.2\ \mu\text{m}$ filter into the final package.
Nonspecific Endonuclease Activity:	A sample is incubated for 14–16 hr with supercoiled plasmid DNA and analyzed by agarose gel electrophoresis.
Exonuclease Activity:	A sample is incubated for 14–16 hr with labeled double-stranded DNA, followed by PAGE analysis.
RNase Activity:	A sample is incubated for 14–16 hr with labeled RNA, followed by PAGE analysis.

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. 4386629B