

## 5M Ammonium Acetate

Store at room temperature.

<b>Catalog #:</b>	AM9070G	AM9071	AM9077G
<b>Amount:</b>	100 mL	500 mL	1 mL
<b>Product Description:</b>	A nuclease-free solution for routine precipitation of nucleic acids.		
<b>Appearance:</b>	Colorless liquid		
<b>Storage Conditions:</b>	Store at room temperature.		

---

### 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**

1. Use 5 M Ammonium Acetate to bring the final concentration of ammonium acetate in the solution to 2.5 M.
2. Add 2 volumes of 100% ethanol for DNA or 2.5 volumes of 100% ethanol for RNA.
3. Chill the solution at or below -20°C for ≥30 min.
4. Centrifuge at top speed in a microcentrifuge for 15 min.
5. Discard the supernatant and wash the pellet with ice cold 70–75% ethanol to remove residual salt.
6. Resuspend in nuclease-free water (Cat #AM9937) or 10 mM Tris, pH 7–8, 1 mM EDTA.

**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.

---

### 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 µm filter into the final package.
<b>Nonspecific Endonuclease Activity:</b>	Meets or exceeds specification when a sample is incubated for 14–16 hr with 300 ng supercoiled plasmid DNA and analyzed by agarose gel electrophoresis.
<b>Exonuclease Activity:</b>	Meets or exceeds specification when a sample is incubated for 14–16 hr with 40 ng labeled <i>Sau3A</i> fragments of pUC19 and analyzed by PAGE.
<b>RNase Activity:</b>	Meets or exceeds specification when a sample is incubated for 14–16 hr with 25 ng labeled RNA and analyzed by PAGE.

---

### OTHER INFORMATION

---

**Material Safety Data Sheets:** Material Safety Data Sheets (MSDSs) can be printed or downloaded from product-specific links on our website at the following address: [www.ambion.com/techlib/msds](http://www.ambion.com/techlib/msds). Alternatively, e-mail your request to [MSDS\\_Inquiry\\_CCRM@appliedbiosystems.com](mailto:MSDS_Inquiry_CCRM@appliedbiosystems.com). Specify the catalog or part number(s) of the product(s), and we will e-mail the associated MSDSs unless you specify a preference for fax delivery. For customers without access to the internet or fax, our technical service department can fulfill MSDS requests placed by telephone or postal mail. (Requests for postal delivery require 1–2 weeks for processing.)

**Warranty and Liability:**

**For research use only. Not for use in diagnostic procedures.**

Applied Biosystems 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](http://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 Applied Biosystems 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. Applied Biosystems assumes no responsibility for any errors that may appear in this document.

Applied Biosystems 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 Applied Biosystems 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:**

Applied Biosystems, AB (Design), and Ambion are registered trademarks of Applied Biosystems Corporation or its subsidiaries in the US and/or certain other countries. All other trademarks are the sole property of their respective owners.

© 2008 Ambion, Inc. All rights reserved. 4386629A