

RNaseZap[®] Solution

RNase Decontamination Solution

Part Number AM9780-AM9784



A. Background

RNaseZap[®] RNase Decontamination Solution completely removes RNase contamination from glass and plastic surfaces. It contains three different ingredients known to be active against RNase. It effectively removes high levels of RNase contamination that similar products cannot. RNaseZap Solution has been used to remove RNase contamination from reaction vessels. With thorough rinsing, RNaseZap Solution leaves no residues that are inhibitory to enzymatic reactions.

B. Instructions for use

RNaseZap Solution is ready to use

Do not dilute because dilution will reduce its effectiveness. If there is a precipitate (as may happen at low temperatures), shake and/or heat at 37°C to bring the precipitate back into solution.



CAUTION

Always wear gloves while using RNaseZap Solution because prolonged contact with skin may cause irritation. When using the spray bottle attachment, always work in a fume hood as inhalation of aerosol or vapor may cause irritation to lungs and mucous membranes.



CAUTION

Please note that RNaseZap Solution should not be used on corrodible metal surfaces.

For cleaning work surfaces:

Apply RNaseZap Solution directly to surface to be cleaned, wipe thoroughly with paper towel, rinse with water and then dry with clean paper towel.

For cleaning lab apparatus:

Apply RNaseZap Solution liberally to a paper towel and wipe all exposed surfaces of the apparatus thoroughly. Rinse with water and then wipe dry. Some small parts may be cleaned by briefly soaking them in RNaseZap Solution, rinsing them with water and then drying.

For cleaning plastic and glass vessels:

Add or spray enough RNaseZap Solution so that the entire surface of the vessel can be coated with the solution upon swirling, or vortexing in the case of centrifuge and microfuge tubes. After discarding the solution, rinse vessels thoroughly two times with distilled water.

For cleaning pipettors:

Remove shaft from pipettor according to manufacturers instructions. Remove seals and gaskets from the shaft and then soak shaft for one minute in RNaseZap Solution. Rinse the shaft thoroughly with water and then reassemble pipettor.

C. Safety Information

Chemical safety guidelines

To minimize the hazards of chemicals:

- Read and understand the Material Safety Data Sheets (MSDS) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the MSDS.

- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the MSDS.
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended on the MSDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.

About MSDSs

Chemical manufacturers supply current Material Safety Data Sheets (MSDSs) with shipments of hazardous chemicals to new customers. They also provide MSDSs with the first shipment of a hazardous chemical to a customer after an MSDS has been updated. MSDSs provide the safety information you need to store, handle, transport, and dispose of the chemicals safely.

Each time you receive a new MSDS packaged with a hazardous chemical, be sure to replace the appropriate MSDS in your files.

Obtaining the MSDS

To obtain Material Safety Data Sheets (MSDSs) for any chemical product supplied by Applied Biosystems or Ambion:

- 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. Click **MSDS**, then right-click to print or download.
- E-mail (MSDS_Inquiry_CCRM@appliedbiosystems.com) or telephone (650-554-2756; USA) your request, specifying the catalog or part number(s) and the name of the prod-

uct(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.

For the MSDSs of chemicals not distributed by Applied Biosystems or Ambion, contact the chemical manufacturer.



Manual 9780M Revision C

Revision Date: September 30, 2008

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

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.

Literature Citation: When you are describing a procedure utilizing this product in a Materials and Methods Section for publication, we would appreciate that you refer to it as RNaseZap[®] Solution.

Warranty and Liability: 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 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.

Trademarks: Applied Biosystems, AB (Design), Ambion, and RNaseZap are registered trademarks of Applied Biosystems, Inc. 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.

AB Applied Biosystems

Ambion[®]

U.S./Canada
(English) 800-327-3002
(French) 800-668-6913
Fax +1-512-651-0201

Europe
tel +44 (0)1480-373-020
fax +44 (0)1480-373-010

Japan
tel 81 (0) 3 5566 6230
fax 81 (0) 3 5566 6507

Direct free phone numbers, distributors:
www.appliedbiosystems.com