

RNAlater® Tissue Collection: RNA Stabilization Solution

IMPORTANT! Before using this product, read and understand the “Safety Information” appendix in this document.

- Product Description 1
- Product Information 2
- Guidelines for Use of RNAlater® Solution 4
- Storage in RNAlater® Solution 5
- RNA Isolation from Samples in RNAlater® Solution 7
- Quality Control 8
- Appendix A Safety Information 9
- Documentation and Support 11

Product Description

RNAlater® Tissue Collection: RNA Stabilization Solution is an aqueous tissue storage reagent that rapidly permeates most tissues to stabilize and protect RNA in fresh specimens. It eliminates the need to immediately process or freeze samples; the specimen can simply be submerged in RNAlater® Solution and stored for analysis at a later date.

Samples in RNAlater® Solution can be stored for extended periods under conditions where RNA degradation would normally take place rapidly (Figure 1). Tissues can be stored indefinitely in RNAlater® Solution at -20°C or below.

Figure 1 RNA from Tissue Stored in RNAlater® Solution. RNA was extracted from mouse tissues stored in RNAlater® Solution as shown. The top panel is an ethidium bromide-stained denaturing agarose gel; the bottom panel shows a Northern blot of the same gel.

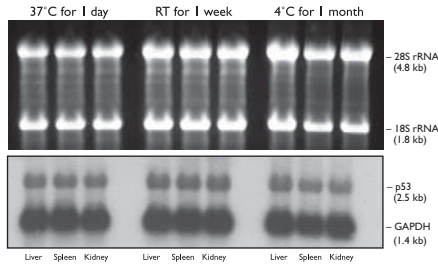
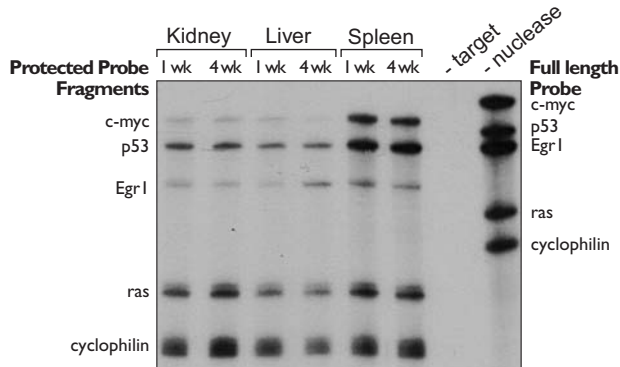


Figure 2 mRNA Profiles of Mouse Tissues Stored in RNAlater® Solution. Mouse tissues were stored in RNAlater® Solution for 1 or 4 weeks at 4°C. RNA was isolated from each tissue and analyzed using the Ambion® RPA III™ Kit. The data demonstrate the stability of expression profiles in tissue stored in RNAlater® Solution.



Product Information

Storage and stability

- Store RNAlater® Solution at room temperature.
- If any precipitation of RNAlater® Solution is seen, heat it to 37°C and agitate to redissolve it.

Sample types compatible with RNAlater® Solution

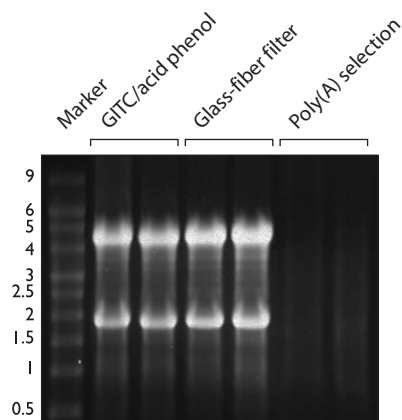
RNAlater® Solution can be used for RNA preservation with most tissues, cultured cells, bacteria, and yeast. It may not be effective in tissues that are poorly penetrated by the solution, such as waxy plant tissue and bone.

RNAlater® Solution has been extensively tested with animal tissues, including brain, heart, kidney, spleen, liver, testis, skeletal muscle, fat, lung, and thymus. It has also been proven effective for RNA preservation in *E. coli*, *Drosophila*, tissue culture cells, white blood cells, and some plant tissues. For more information, go to www.invitrogen.com/site/us/en/home/support/technical-support.html.

RNA isolation from RNAlater® Solution

RNAlater® Solution is compatible with most RNA isolation methods. Samples stored in RNAlater® Solution have been used successfully with TRI Reagent® Solution (P/N AM9738), and all of Ambion® RNA isolation kits and reagents, including: the TōTALLY RNA™ Kit, the PARIS™ Kit, the *mirVana*™ miRNA Isolation Kit, and the RNAqueous® and Poly(A)Purist™ product families.

Figure 3 RNA isolated from tissue stored in RNAlater® Solution using different methods. Whole mouse hearts (left lane of each set) and livers (right lane of each set) were dissected, and stored in RNAlater® Solution for 3 days at 4°C. RNA was isolated from equal mass amounts of each tissue using the indicated methods. RNA (5 µg) was run on denaturing agarose and stained with ethidium bromide.



Isolating genomic DNA from RNAlater® Solution-stored samples

DNA can be isolated from RNAlater® Solution-stored samples. For more information, go to www.invitrogen.com/site/us/en/home/support/technical-support.html.

Isolating protein from RNAlater® Solution-stored samples

Proteins are also preserved in RNAlater® Solution. RNAlater® Solution will denature proteins; therefore, protein obtained from samples stored in it will be suitable for applications such as Western blotting or 2D gel electrophoresis, but not for applications that require native protein.

Guidelines for Use of RNAlater® Solution

- Use RNAlater® Solution with *fresh tissue only*; do not freeze tissues before immersion in RNAlater® Solution.
- Before immersion in RNAlater® Solution, cut large tissue samples to ≤ 0.5 cm in any single dimension.
- Place the fresh tissue in 5–10 volumes of RNAlater® Solution.
- Most samples in RNAlater® Solution can be stored at room temperature for 1 week without compromising RNA quality, or at -20°C or -80°C indefinitely.
- Do not freeze samples in RNAlater® Solution immediately; store at 4°C overnight (to allow the solution to thoroughly penetrate the tissue), remove supernatant, then move to -20°C or -80°C for long-term storage.

Note: We offer RNAlater®-ICE (P/N AM7030) to recover tissues that have already been frozen. RNAlater®-ICE renders frozen tissues pliant enough for homogenization while maintaining the low temperatures needed to protect the RNA from degradation.

Animal Tissue

RNAlater® Solution does not disrupt the structure of tissues; thus, tissue that has been equilibrated in RNAlater® Solution can be removed from the solution, sectioned into smaller pieces, and returned to RNAlater® Solution, if desired.

Small organs such as mouse liver, kidney and spleen can be stored whole in RNAlater® Solution.

Plant Tissue	Plant tissues that have natural barriers to diffusion, such as waxy coatings on leaves, will often require disruption to allow RNAlater® Solution access to the tissue. However, many plant tissues can simply be submerged in RNAlater® Solution whole; we have successfully isolated intact RNA from tobacco leaf explants, entire <i>Arabidopsis</i> and alfalfa seedlings, and from potato shoot tips.
Tissue Culture Cells	Pellet cells according to the protocols followed by your laboratory. Remove supernatant and then add 5–10 volumes RNAlater® Solution. The cells can be washed in PBS before resuspending in RNAlater® Solution, if desired.
Blood and Plasma	White blood cells can be effectively preserved in RNAlater® Solution when separated from the red blood cells and sera and treated as tissue culture cells. RNAlater® Solution can also be added to small volumes of anticoagulated whole blood, sera, and plasma; however, the procedure is not presented here—see the Ambion RiboPure™-Blood Kit (P/N AM1928) protocol for detailed instructions.
Yeast	<p>Pellet up to 3×10^8 cells (centrifuge at $12,000 \times g$ for 2 min). Remove supernatant and immediately resuspend the pellet in 0.5–1 mL of RNAlater® Solution. Yeast cells can be stored in RNAlater® Solution for up to 8 hr at 25°C, or up to a week at 4°C.</p> <p>For long-term storage, incubate the cells in RNAlater® Solution for 1 hr. Repellet the cells (centrifuge at $>12,000 \times g$ for 5 min), remove supernatant, flash freeze, and store at –80°C.</p>
Bacteria	RNAlater® Solution is bacteriostatic; although bacteria do not grow in it, the cells remain intact. <i>E. coli</i> stored in RNAlater® Solution for 1 month at 4°C are intact and yield undegraded RNA.

Storage in RNAlater® Solution

If refrigeration is available:	<p>Storage at –80°C</p> <p>Storage at –80°C is recommended for archival samples and will provide optimal preservation. Samples can be stored at –80°C indefinitely. RNAlater® Solution will freeze at –80°C.</p>
---------------------------------------	---

To prepare samples for storage at -80°C , first incubate the samples in RNAlater® Solution overnight at 4°C to allow thorough penetration of the tissue, then transfer to -80°C . To expedite thawing of the samples, we recommend removing the tissue, or pelleting cells, from the RNAlater® Solution before freezing at -80°C .

Samples can subsequently be thawed at room temperature and refrozen without significantly affecting the amount or the integrity of the recoverable RNA.

Storage at -20°C

Storage at -20°C can also be used for archival samples. Samples will not freeze at -20°C , but crystals may form; this will not affect subsequent RNA isolation. Samples can be stored at -20°C indefinitely.

To prepare samples for storage at -20°C , first incubate the samples in RNAlater® Solution overnight at 4°C to allow thorough penetration of the tissue, then transfer to -20°C .

Samples can subsequently be thawed at room temperature and refrozen without affecting the amount or the integrity of the recoverable RNA.

Storage at 4°C

Most samples can be stored in RNAlater® Solution at 4°C for up to 1 month without significant RNA degradation.

If refrigeration is not available:

Place samples in the coolest environment available. If ambient temperature is above 25°C , incubate the samples in RNAlater® Solution on ice for a few hours, if possible, before storing at ambient temperature.

Storage at 25°C (room temperature)

Most samples can be stored at 25°C in RNAlater® Solution for up to 1 week without significant loss of RNA quality. After 2 weeks at 25°C , RNA generally appears slightly degraded (marginally acceptable for Northern analysis, but still of sufficient quality for nuclease protection assays or RT-PCR analysis).

Storage at 37°C

RNA isolated from samples stored at 37°C is intact after a 24 hour incubation, but is partially degraded after 3 days.

RNA Isolation from Samples in RNAlater® Solution

Remove RNAlater® Solution from samples

RNase inactivation is reversible; do not rinse RNAlater® Solution from samples before using. Blot tissues with a wipe, or pellet cells to remove excess RNAlater® Solution.

Tissue

Retrieve tissue from RNAlater® Solution with sterile forceps, quickly blot away excess RNAlater® Solution with an absorbent lab wipe or paper towel, and then submerge the sample in RNA isolation lysis solution. Homogenize tissue promptly after placing it in lysis/denaturation solution.

Cells

There are two options for isolating RNA from cells stored in RNAlater® Solution. The preferred method is to remove the solution from the cells prior to extraction. Alternatively, cells in RNAlater® Solution can be used directly for RNA extraction. Because of the greater volume that the cells are in, this method generally requires additional lysis solution.

- **Removal of RNAlater® Solution prior to extraction**

Because of the density of RNAlater® Solution, greater centrifugal forces are required to pellet cells from RNAlater® Solution than from normal media. Generally, cells become much less fragile when stored in RNAlater® Solution and can be centrifuged at high speed without lysis. Most cell types can be centrifuged at 5000 x g without damage to the cells. Since different cell types vary in their ability to withstand centrifugal forces, we recommend testing the centrifugal speed with an expendable sample. Alternatively, dilute the RNAlater® Solution by adding an equal volume of ice cold PBS (or other buffered solution) immediately before centrifugation to reduce the density of the solution, then centrifuge at normal speeds.

- ***RNA extraction from cells in RNAlater[®] Solution***

One-step phenol-based disruption/extraction solutions, such as Ambion TRI Reagent[®] Solution or RNAwiz[™] Reagent (available only in Japan), can be used to purify RNA from cells suspended in RNAlater[®] Solution. This can be done by adding ten volumes of the one-step solution to the cell mixture, and proceeding normally. When RNAwiz Reagent is used in this way, it may be necessary to dilute the aqueous phase before the RNA precipitation step. See below for more information.

Tips for RNA isolation

Glass fiber-based extraction

Lysates from RNAlater[®] Solution-treated samples often require more force to pass through glass-fiber filters than lysates from untreated samples. Therefore, it may be necessary to use centrifugation instead of vacuum pressure to pass lysates through glass-fiber filters.

One-step disruption/extraction solutions

When using one-step RNA isolation products such as TRI Reagent Solution or RNAWIZ Reagent (available only in Japan), on RNAlater[®] Solution-preserved samples, the aqueous phase will occasionally appear cloudy; this will not adversely affect RNA recovery or quality.

With RNAWIZ Reagent, there may be a problem getting the aqueous phase to mix with isopropanol at the precipitation step because of RNAlater[®] Solution carryover. If this occurs, add a mixture of 50% water, 50% isopropanol until the solution becomes clear and the two phases mix. The amount of water/isopropanol required will depend on how much RNAlater[®] Solution was carried over; if the sample was mostly RNAlater[®] Solution, as much as an equal volume may be needed.

Quality Control

RNAlater[®] Solution undergoes quality assurance testing to verify that its composition is invariant from lot to lot.

Appendix A Safety Information



WARNING! GENERAL SAFETY. Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
 - Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs), and use appropriate personal protective equipment (gloves, gowns, eye protection, etc.). To obtain SDSs, see the “Documentation and Support” section in this document.
-

Chemical safety



WARNING! GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below, and consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the “Documentation and Support” section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer’s cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary

container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)

- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.

IMPORTANT! Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

Biological hazard safety



WARNING! Depending on the samples used on the instrument, the surface may be considered a biohazard. Use appropriate decontamination methods when working with biohazards



WARNING! BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Follow all applicable local, state/provincial, and/or national regulations. Wear appropriate protective equipment, which includes but is not limited to: protective eyewear, face shield, clothing/lab coat, and gloves. All work should be conducted in properly equipped facilities using the appropriate safety equipment (for example, physical containment devices). Individuals should be trained according to applicable regulatory and company/institution requirements before working with potentially infectious materials. Read and follow the applicable guidelines and/or regulatory requirements in the following:

- U.S. Department of Health and Human Services guidelines published in Biosafety in Microbiological and Biomedical Laboratories found at: www.cdc.gov/biosafety
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030; www.access.gpo.gov/nara/cfr/waisidx_01/29cfr1910a_01.html).

- Your company's/institution's Biosafety Program protocols for working with/handling potentially infectious materials.
 - Additional information about biohazard guidelines is available at: www.cdc.gov
-

In the EU:

Check local guidelines and legislation on biohazard and biosafety precaution and refer to the best practices published in the World Health Organization (WHO) Laboratory Biosafety Manual, third edition, found at: www.who.int/csr/resources/publications/biosafety/WHO_CDS_CSR_LYO_2004_11/en/

Documentation and Support

Obtaining SDSs

Safety Data Sheets (SDSs) are available from:

- www.invitrogen.com/sds

or

- www.appliedbiosystems.com/sds

Note: For the SDSs of chemicals not distributed by Life Technologies, contact the chemical manufacturer.

Obtaining support

For the latest services and support information for all locations, go to:

- www.invitrogen.com

or

- www.appliedbiosystems.com

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches



7020MF

For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.

Information in this document is subject to change without notice.

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. TO THE FULLEST EXTENT ALLOWED BY LAW, 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, WHETHER OR NOT FORESEEABLE AND WHETHER OR NOT APPLIED BIOSYSTEMS IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

LIMITED USE LABEL LICENSE: RESEARCH USE ONLY

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

© 2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. TRI Reagent is a registered trademark of Molecular Research Center, Inc.

Part Number 7020M Rev. F 03/2011

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA
Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit www.appliedbiosystems.com/support

www.lifetechnologies.com

