


RNase A (Biochemistry Grade)

Catalog Number AM2274

Pub. No. 4393904 Rev. B

Contents	Quantity	Storage conditions
RNase A (Biochemistry Grade), 1 µg/mL	200 µL	Store at -20°C. <i>Do not store in a frost-free freezer.</i>
10X RNA Structure Buffer	1 mL	
1X Alkaline Hydrolysis Buffer	1 mL	
1X RNA Sequencing Buffer	1 mL	
Precipitation/Inactivation Buffer (add 3.2 mL 100% ethanol before use)	4.8 mL	
Gel Loading Buffer II	1.4 mL	
Yeast RNA (10 mg/mL)	100 µL	

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

Product description

RNase A (Biochemistry Grade) is isolated by an extensive series of purification steps including affinity chromatography. RNase A is an endonuclease that specifically cleaves 3' of U and C residues. RNase A is used in rapid analysis of the physical structure of RNA, in combination with single-strand specific RNase T1 and double-strand specific RNase V1.

Source: Manufactured from Bovine pancreas by a USDA-approved source.

Unit (U) definition: One unit is the amount of RNase A required to give an increase of 0.0146 A_{286} absorbance units per minute in a 1-mL volume at room temperature and is equivalent to 0.1177 Kunitz Units. Assay Conditions: 100 mM Tris-acetate (pH 6.5), 1 mM EDTA and 1 mM cyclic 2', 3'-CMP.

Storage buffer (not included): 10 mM HEPES (pH 7.2), 20 mM NaCl, 0.1% Triton X-100, and 50% glycerol (v/v).

General information about biochemistry-grade ribonucleases

Biochemistry-grade ribonucleases are optimized for the study of RNA structure, RNA sequencing, protein foot-printing, and

boundary experiments. They are also tested for purity to ensure the absence of nonspecific nuclease or other contaminating ribonuclease activities that could introduce unexpected cleavage sites and interfere with RNA structure studies.

In addition to RNA structural analysis, biochemistry-grade ribonucleases can be used to map protein binding sites on RNAs by comparing cleavage patterns in the presence and absence of an RNA binding protein.

Using RNase A (Biochemistry Grade)

To visualize the RNase A digestion products, the RNA should be end-labeled either at its 5' end using a kinase reaction with [γ - 32 P]ATP (KinaseMax™ Kit, Cat. no. AM1520), or at its 3' end using an RNA ligation reaction with [32 P]pCp (T4 RNA Ligase, Cat. no. AM2140).

RNA structure analysis

In the following procedure, each RNA is digested with three sequential 10-fold dilutions of ribonuclease, to quickly optimize conditions for distinguishing nucleotides that are structurally constrained from those that are not. Further dilution may be necessary to achieve the optimal digestion ladder.

Note: Before you begin, complete the Precipitation/Inactivation Buffer by adding 3.2 mL of 100% ethanol to the bottle supplied.

1. Thaw the 10X RNA Structure Buffer and end-labeled RNA at room temperature.

2. Mix 4 μL of 10X RNA Structure Buffer, 0.2–4 μg of end-labeled RNA, 4 μg of yeast RNA, and Nuclease-free Water to bring the final volume to 36 μL .
3. Distribute 9 μL aliquots of the RNA/buffer/water mixture into 4 microcentrifuge tubes numbered 1–4. To Sample 2 add 1 μL of ribonuclease. Mix thoroughly by pipetting.
4. Transfer 1 μL from Sample 2 to Sample 3. Mix thoroughly by pipetting.
5. Transfer 1 μL from Sample 3 to Sample 4. Mix thoroughly by pipetting.
6. Incubate all samples at room temperature for 15 minutes.
7. Add 20 μL of completed Inactivation/Precipitation Buffer, vortex and incubate at -20°C for 15 minutes.
8. Microcentrifuge at maximum speed for 15 minutes at 4°C , aspirate the supernatant, and wash the pellet with 70% ethanol.
9. Air dry the pellet and dissolve it in 7 μL of Gel Loading Buffer II.
10. Heat the samples at 95°C for 5 minutes, then fractionate 3 μL of each sample using a 6–20% acrylamide/7 M Urea sequencing gel.
11. Use autoradiography to assess the digestion products.

Expected results: Sample 1 is the end-labeled RNA before nuclease treatment, and any non-full length bands represent cleavage products already present within the RNA sample. These bands will also be present in the nuclease-treated samples and should be disregarded in your analysis. Samples 2, 3, and 4 were digested with decreasing amounts of ribonuclease. Bands evident in the samples with lower amounts of ribonuclease typically represent nucleotides that are most accessible to the ribonuclease.

Alkaline hydrolysis

This procedure provides a "ladder" of hydrolyzed RNA fragments for gel electrophoresis. In the procedure, three

hydrolysis times are used to provide RNA samples that are increasingly hydrolyzed. Select the ladder that provides the best distribution of nucleic acids over the range of lengths needed for your experiments.

1. Add 0.1–3 μg end-labeled RNA and 3 μg yeast tRNA in a volume not to exceed 5 μL .
2. Add sufficient 1X Alkaline Hydrolysis Buffer to bring the final reaction volume to 15 μL .
3. Distribute 5 μL aliquots of the RNA/buffer mixture into 3 tubes labeled 1–3.
4. Heat each sample to 95°C .
5. After 2 minutes, remove Sample 1 and place it in an ice bucket.
6. After 5 minutes, remove Sample 2 and place it in an ice bucket.
7. After 15 minutes, remove Sample 3 and place it in an ice bucket.
8. Add 10 μL of Gel Loading Buffer II to each of the 3 samples. Prepare an untreated sample by mixing 1 μL of 5' end-labeled RNA with 8 μL of Gel Loading Buffer II.
9. Fractionate 3 μL of each sample (4 samples total) using a 6–20% acrylamide/7 M urea sequencing gel.
10. Use autoradiography to visualize the fractionated RNA products.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

For Research Use Only. Not for use in diagnostic procedures.

The information in this guide is subject to change without notice.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) 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.

NOTICE TO PURCHASER: 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.

© 2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com

9 October 2012

