AM7940 Catalog #:



Store at or below -70°C. Do not store in a frost-free freezer.

Catalog #: AM7940 Amount: 200 µg Concentration: 1 mg/mL 1.7 - 2.1Absorbance 260/280:

Storage Conditions: Store at or below -70°C. Avoid multiple freeze-thaw cycles. Aliquots of the product may be stored short-term at

–20°C. Do not store in a frost-free freezer.

Storage Buffer: THE RNA Storage Solution (1 mM Sodium Citrate, pH 6.4; Cat #AM7000)

USER INFORMATION

Product Description: E. coli Total RNA is prepared from DH5α cells. Total RNA is prepared using Ambion RNA isolation reagents. The

purified RNA undergoes a rigorous DNase treatment. It has been precisely quantified and is provided at 1 mg/mL in THE RNA Storage Solution. The integrity of the RNA is verified by capillary electrophoresis using the Agilent® 2100

bioanalyzer.

E. coli Total RNA provides the researcher with access to RNA isolated from a source that might otherwise be inconvenient or expensive to isolate. In addition, the RNA may serve as a positive control when a particular mRNA is known to be expressed in this strain. E. coli Total RNA is ready for use in RT-PCR, ribonuclease protection assays,

S1 nuclease assays, cDNA synthesis, Northern analysis, and in vitro translation.

Handling Instructions: RNA is very sensitive to degradation by exogenous ribonucleases introduced during handling. Wear gloves when

handling this product. Use RNase-free reagents, tubes, and barrier pipette tips.

Thawing Instructions

Thaw just to completion at 37°C, vortex for a few seconds when fully thawed, and place on ice. Aliquot the RNA, if

necessary, to minimize freeze-thaw cycles (≤5).

QUALITY CONTROL

Nonspecific Endonuclease

Activity:

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.

Meets or exceeds specification when a sample is incubated for 14–16 hr with 40 ng 32P-labeled Sau3A fragments of **Exonuclease Activity:**

pUC19 and analyzed by PAGE.

All Total RNA products undergo accelerated stability testing. RNA integrity is checked on the Agilent bioanalyzer **Functional Testing:**

both before and after a 16–18 hr incubation at 37°C. DNA contamination is checked by real-time PCR using 16S

primers and probe.

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. Alternatively, e-mail your request to 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 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, Patents, and Licensing:

Applied Biosystems, AB (Design), and Ambion are registered trademarks of Applera Corporation or its subsidiaries in the US and/or

certain other countries. All other trademarks are the sole property of their respective owners.

© 2007 Ambion, Inc. All rights reserved. 4386600A

Applied Biosystems Ambion