

RNase-Free Microcentrifuge Tubes

Catalog Numbers AM12300, AM12350, AM12400, AM12425, AM12450, AM12475, AM12480

Pub. No. MAN1001375 Rev. A



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 [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

Ambion™ RNase-free microcentrifuge tubes are marked with graduations for convenience, have flat lids for easy labeling, and have a positive seal design which prevents caps from coming open in boiling water baths. These polypropylene tubes are autoclave-safe.

The 2.0 mL Elution Tubes can accommodate the RNAqueous™ Kit (Cat. no. [AM1912](#)).

Contents

Catalog numbers that appear as links open the web pages for those products.

Product	Cat. No.	Amount
Nonstick, RNase-Free Microfuge Tubes, 0.5 mL	AM12350	500
RNase-Free Microcentrifuge Tubes, 0.5 mL	AM12300	1000
Nonstick, RNase-free Microfuge Tubes, 1.5 mL	AM12450	250
RNase-Free Microcentrifuge Tubes, 1.5 mL	AM12400	500
Nonstick, RNase-Free Microfuge Tubes, 2.0 mL	AM12475	250
RNase-Free Microcentrifuge Tubes, 2.0 mL	AM12425	500
Elution Tubes, RNase-Free Microfuge Tubes, 2.0 mL	AM12480	100

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.

[thermofisher.com/support](https://www.thermofisher.com/support) | [thermofisher.com/askaquestion](https://www.thermofisher.com/askaquestion)



Life Technologies Corporation | 2130 Woodward Street | Austin, Texas 78744 USA

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

Revision history: Pub. No. MAN1001375 A

Revision	Date	Description
A	11 March 2025	<ul style="list-style-type: none"> First release with new publication number format. Supersedes publication number 4386463 Rev. C. The product names for Cat. No. AM12300, AM12400, and AM12425 were updated. The quality control section was removed.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2025 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

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