

# Ribonuclease H

Catalog Number 18021-071

Doc. Part No. 18021071.pps Pub. No. MAN0001330 Rev. A.0



**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](http://thermofisher.com/support).

## Product description

The Invitrogen™ Ribonuclease H (RNase H) is an endoribonuclease that specifically degrades the RNA strand of an RNA-DNA hybrid to produce 5' phosphate-terminated oligoribonucleotides and single-stranded DNA. It is suitable for removing mRNA during second-strand cDNA synthesis, for removing the poly(A) sequences from mRNA in the presence of oligo(dT), and for oligodeoxyribonucleotide-directed cleavage of RNA. The RNase H is purified from *E. coli* expressing the *E. coli* RNase H gene from a plasmid.

## Contents and storage

Contents	Amount	Storage
Ribonuclease H (2 U/μL)	4 × 30 U	Store at -20°C in a non-frost-free freezer.

**Note: Unit (U) definition:** One unit is defined as the amount of RNase H that degrades 1 nmol of poly(A) in 20 minutes at 37°C

### Storage buffer

20 mM Tris-HCl (pH 7.5), 100 mM KCl, 10 mM MgCl<sub>2</sub>, 0.1 mM EDTA, 0.1 mM DTT, 50 μg/mL BSA, 50% (v/v) glycerol

### 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.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

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

---

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

## DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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.

**Revision history:** Pub. No. MAN0001330

Revision	Date	Description
A.0	15 June 2016	Format, style, and legal updates
—	25 May 2007	Baseline for this revision history

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

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA  
1 800 955 6288

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit [thermofisher.com/support](http://thermofisher.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[thermofisher.com](http://thermofisher.com)

15 June 2016

**ThermoFisher**  
SCIENTIFIC