

DNase I

Catalog Number 18047-019

Doc. Part No. 18047019.pps Pub. No. MAN0001337 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.

Product description

The Invitrogen™ DNase I digests single- and double-stranded DNA to oligodeoxyribonucleotides that contain a 5' phosphate. Ribonuclease has been reduced to non-detectable levels. DNase I is suitable for removing DNA from protein preparations, nick translating DNA, and generating random fragments for dideoxy sequencing. DNase I is purified from bovine pancreas. The specific activity of DNase I is typically in the range of 10,000–25,000 units/mg. To remove DNA from RNA preparations, use DNase I, Amplification Grade (Cat. No. 18068-015).

Contents and storage

Contents	Amount	Storage
DNase I (deoxyribonuclease; 50–375 U/μL)	20,000 U	Store at –20°C in a non-frost-free freezer.

Note: Unit (U) definition: One unit increases the absorbance of a high molecular weight DNA solution at a rate of 0.001 A_{260} units/minute/mL of reaction mixture at 25°C.

Storage buffer

- 20 mM sodium acetate (pH 6.5)
- 5 mM CaCl₂
- 0.1 mM PMSF
- 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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

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. MAN0001337

Revision	Date	Description
A.0	15 June 2016	Format, style, and legal updates
—	1 October 2008	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 or email techsupport@lifetech.com

thermofisher.com

15 June 2016

ThermoFisher
SCIENTIFIC