


Recombinant DNase I (rDNase I)

Catalog Number AM2235

Pub. No. 4393899 Rev. B

Contents	Quantity	Storage conditions
rDNase I, 2 U/μL	1000 Units	Store at -20°C. <i>Do not store in a frost-free freezer.</i>
10X DNase I Buffer <ul style="list-style-type: none"> • 100 mM Tris-HCl, pH 7.5 • 25 mM MgCl₂ • 5 mM CaCl₂ 	1 mL	

 **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.

for RNA:DNA hybrids and single-stranded DNA is at least 2 orders of magnitude below that for double-stranded DNA. DNase I requires both Mg²⁺ and Ca²⁺ for maximal activity (Clark and Eichhorn, 1974; Junowicz and Spencer, 1973; Price, 1975).

Product description

Recombinant DNase I (rDNase I) is a highly purified DNase I derived from cloned bovine DNase I grown in a host with RNase levels that are 10⁷-fold lower than bovine pancreas.

Source: A non-animal host that overexpresses bovine DNase I.

Unit (U) definition: One unit is the amount of enzyme required to completely degrade 1 μg DNA in 10 min at 37°C, and it is equivalent to 0.04 Kunitz units.

Storage buffer (not included): 20 mM HEPES pH 7.5, 10 mM CaCl₂, 10 mM MgCl₂, 1 mM DTT and 50% (v/v) glycerol.

General information

Recombinant DNase I (rDNase I) is a recombinant form of bovine DNase I purified from an overexpressing host. rDNase I is bovine-free. In addition, the host expressing rDNase I contains 10⁷-fold less endogenous RNase activity than pancreatic tissues, the source for bovine DNase I. As a result, rDNase I can be more readily purified from contaminating RNases. The enzyme is provided free from any detectable RNase activity. rDNase I has a molecular weight of approximately 30 kDa.

DNase I is an endonuclease that hydrolyzes phosphodiester linkages yielding oligonucleotides with a 5'-phosphate and a free 3'-hydroxyl group (Kunitz, 1950). DNase I has been shown to act on single- and double-stranded DNA, chromatin, and RNA:DNA hybrids, although the specific activity of DNase I

Using rDNase I

The reaction kinetics of rDNase I are identical to that of native DNase I. Thus, rDNase I can be directly substituted for native DNase I with equivalent unit activity. Use rDNase I to degrade DNA in the presence of RNA when the absence of RNase is critical to maintain the integrity of the RNA. For example, DNase I is frequently used to remove template DNA after *in vitro* transcription reactions (Krieg et al, 1985), or to destroy genomic DNA in RNA preparations prior to reverse transcription-PCR (RT-PCR).

Removal of contaminating genomic DNA from RNA samples

1. If the nucleic acid solution concentration is >200 μg/mL, dilute it to 10 μg nucleic acid/50 μL.
2. Add 10X DNase I Buffer (supplied) to 1X concentration in the RNA sample.
3. Add 1 μL rDNase I (2 U) for up to 10 μg RNA in a 50 μL reaction, and incubate at 37°C for 30 minutes.
These reaction conditions will remove up to 2 μg of genomic DNA.
4. Extract the RNA sample with phenol/chloroform to inactivate the rDNase I.

Degradation of DNA template in a transcription reaction

1. After transcription, add 2 U of rDNase I to a 20 μ L transcription reaction. It is not necessary to add 10X DNase I Buffer to the transcription reaction.
2. Incubate at 37°C for 15 minutes.
 - If the transcript is to be gel purified, then gel loading buffer may be added directly to the rDNase I-treated transcription reaction.
 - If not, the rDNase I can be inactivated by phenol-chloroform extraction.

Conditions for complete DNA digestion

1. Add 10X DNase I Buffer to 1X concentration in the solution to be DNase-treated, and add approximately 1–2 U of rDNase I per 1 μ g DNA present.
2. Incubate at 37°C for 15–30 minutes.

Heat inactivation of rDNase I

Some protocols suggest heating at 75°C for 5 min to inactivate DNase I (Huang, Fasco, and Kaminsky, 1996). We recommend a 10-minute incubation at 75°C for complete inactivation of DNase I at a concentration of 0.1 U/ μ L. If this is the preferred method of inactivation, add EDTA to a final concentration of 5

mM before heating. If EDTA is not added, the RNA will undergo chemical scission when heated.

Note: For RNA samples that are to be used in reverse transcription reactions, the EDTA concentration in the RNA sample must be taken into account. Excess EDTA in an RNA sample may lower the free Mg^{2+} concentration and affect the efficiency of reverse transcription. After heat inactivation of DNase I, it may be necessary to add additional Mg^{2+} for maximum reverse transcriptase activity. Alternatively, DNase I can be inactivated and removed by phenol/chloroform extraction.

Gel analysis

Gel loading buffers should contain EDTA to eliminate DNase I activity; denaturing gel loading buffers containing mostly formamide or formaldehyde, such as Gel Loading Buffer II (Cat. no. AM8546G, AM8547) will also inactivate DNase I.

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.

References

- Clark, R. and Eichhorn, GL. (1974) *Biochem* 13, 5098.
- Huang Z, Fasco MJ, and Kaminsky LS (1996) *BioTechniques* 20,1012–1020.
- Junowicz, E. and Spencer, JH. (1973) *BBA* 312, 72.
- Krieg, PA, et al. (1985) *Genetic Engineering Principles and Methods* (Setlow JK, Hollaender A, ed.) Vol. 7, Plenum Press, New York, London.
- Kunitz, M. (1950) *J Gen Physiol* 33, 349.
- Price, P.A. (1975) *J Biol Chem* 250,1981–1986.

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

17 October 2012

