



3050 Spruce Street  
Saint Louis, Missouri 63103 USA  
Telephone 800-325-5832 • (314) 771-5765  
Fax (314) 286-7828  
email: techserv@sial.com  
sigma-aldrich.com

## Product Information

### Deoxyribonuclease I

Catalog Number **D7291**

Storage Temperature  $-20^{\circ}\text{C}$

## TECHNICAL BULLETIN

CAS RN: 9003-98-9

EC 3.1.21.1

Synonym: DNase I

### Product Description

Deoxyribonuclease I is isolated from bovine pancreas and is processed to reduce RNase activity to below detectable levels.

DNase I digests single- and double-stranded DNA to a mixture of mono- and oligonucleotides carrying 5'-phosphates and 3'-OH termini. This catalytic activity is divalent ion-dependent. In the presence of  $\text{Mg}^{2+}$ , DNase I hydrolyzes each strand of double-stranded DNA randomly and independently. In the presence of  $\text{Mn}^{2+}$ , both strands can be cleaved. It is useful for nick translation, DNase footprinting, bisulfite-mediated mutagenesis, and RNA purification<sup>1,2</sup>, and is suitable for all but the most stringent RNA purification procedures.

However, RNA treated with this DNase I should not be used to generate a cDNA library or used for RT-PCR. For these sensitive applications, it is recommended to use DNase I, Amplification Grade, Catalog Number AMP-D1.

Deoxyribonuclease I is supplied as a solution in 50% (v/v) glycerol, 20 mM sodium acetate, pH 6.5, 5 mM  $\text{CaCl}_2$ , 0.1 mM PMSF.

Specific activity:  $\geq 10,000$  units/mg protein

Unit Definition: One unit will cause a  $\Delta A_{260}$  of 0.001 per minute per ml reaction mixture using calf thymus DNA as substrate.

RNase: none detected

### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

### Storage/Stability

Store at  $-20^{\circ}\text{C}$

### Procedure

#### Typical protocol to remove trace DNA

Add to an RNase-free PCR tube:

- 1  $\mu\text{g}$  of RNA sample in 8  $\mu\text{l}$  water
- 1  $\mu\text{l}$  of 10X reaction buffer (200 mM Tris-HCl, pH 8.3, with 20 mM  $\text{MgCl}_2$ )
- 1  $\mu\text{l}$  of DNase I, 1 unit/ $\mu\text{l}$ \*

\*Refer to the Certificate of Analysis for the lot specific specific activity and the number of mg protein per ml.

Incubate for 15 minutes at room temperature

To stop add 1  $\mu\text{l}$  of Stop Solution to bind calcium and magnesium ions and to inactivate the DNase I.

The Stop Solution (50 mM EDTA) must be added before heating to prevent metal (Mg/Ca) ion catalyzed hydrolysis of the RNA. Heat at  $70^{\circ}\text{C}$  for 10 minutes to denature both the DNase I and the RNA.

**Note:** This product should not be used for digestions longer than 15 minutes or for digestions at temperatures higher than room temperature, or the residual contaminating RNase activity will begin to degrade the RNA.

## Product Profile

### Activity Assay

Assay Buffer: 100 mM sodium acetate, pH 5.0 at 25 °C,  
containing 5 mM MgCl<sub>2</sub>  
50 µg/ml calf thymus DNA

5-20 µg of DNase I was added to 3 ml of reaction mixture at 25 °C and the  $\Delta A_{260}$  is monitored for 10 minutes. The maximum linear rate was used to calculate the activity.

### RNase Assay

Two µg transfer RNA were incubated with 2 µg DNase I in a 50 µl reaction mixture containing 30 mM Tris-HCl, pH 7.8, 50 mM NaCl and 10 mM MgCl<sub>2</sub> for 16 hours at 37 °C. No degradation of the tRNA was detected by polyacrylamide gel electrophoresis

## References

1. Galas, D.J., and Schmitz, A., DNase footprinting: a simple method for the detection of protein-DNA binding specificity. *Nucleic Acids Res.*, **9**, 3157-3170 (1978).
2. Greenfield, L., et al., Conversion of closed circular DNA molecules to single-nicked molecules by digestion with DNAase I in the presence of ethidium bromide. *Biochim. Biophys. Acta*, **407**, 365-375 (1975)

ND,PHC 07/05-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.