

RNase T1, *Aspergillus oryzae*, Recombinant, *E. coli*

Cat.-No.	Amount
EN-155S	1 mg
EN-155L	5 mg

Supplied as lyophilized powder.

Ribonuclease T1 (RNase T1) is an endoribonuclease that specifically cuts RNA at the 3'-end of guanosine residues through a 2', 3'-cyclic phosphate intermediate mechanism. RNase T1 is cloned from *Aspergillus oryzae* and over expressed in *E. coli* to produce a highly pure enzyme without contaminating DNase or non-specific RNase activity.

RNase T1 is used for RNA structure and mapping studies and for RNA protection assays, as well as for removal of RNA from DNA samples.

AVOID FREEZE/THAW CYCLES.

For in vitro use only!

Purity: > 95% by SDS-PAGE.

Absence of contaminants: Tested for the absence of endo- and exodeoxyribonucleases.

Selected references:

Walz et al. (1979) Base-group specificity at the primary recognition site of ribonuclease T for minimal RNA substrates. *Arch. Biochem. Biophys.* **195**:95.

McGregor et al. (1999) Secondary structure mapping of an RNA ligand that has high affinity for the MetJ repressor protein and interference modification analysis of the protein-RNA complex. *J. Biol. Chem.* **274**:2255.

Krebs et al. (2003) RNaseCut: a MALDI mass spectrometry-based method for SNP discovery. *Nucleic Acids Res.* **31**:e37.

Unit definition: One unit converts 100 ng of *E. coli* ribosomal RNA into acid-soluble nucleotides in 5 seconds at 37°C in 50 mM Tris-HCl (pH 7.0), 2 mM EDTA.

1 mg corresponds to 380.000 Units.

Store: 4 °C