

## Micrococcal Nuclease

**Cat. No.** C06070001

**Source:** E. coli

**Lot #:** 001

**Size:** 250, 500, 1000 µl

**Concentration:** 500 U/ml

**Source:** Staphylococcus aureus

**Purity:** Purified using glycerol gradient and ion exchange chromatography, >98% purity as determined by SDS-PAGE.

**Storage buffer:** 20mM Tris-Cl pH 8.0, 50mM NaCl, 1mM DTT and 50% glycerol.

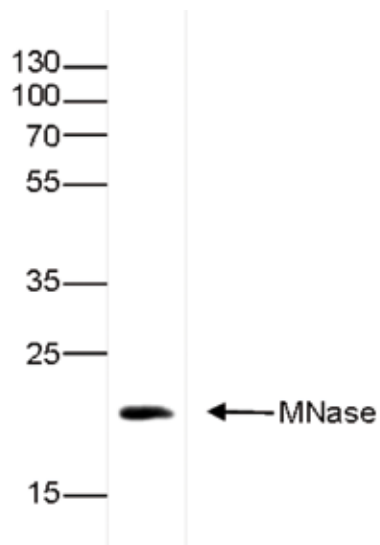
**Unit definition:** One unit will produce 1.0 µmole of acid soluble polynucleotides from native DNA per min at pH 8 at 37 °C, based on EM/260 = 10,000 for the mixed nucleotides.

**Storage:** Store at -20°C; guaranteed stable for 2 years from date of receipt when stored properly.

**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

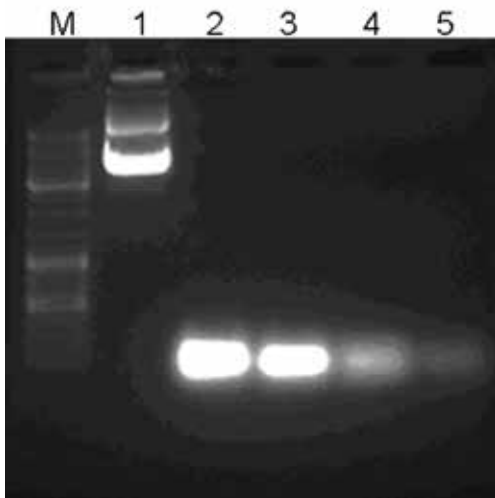
**Description:** Micrococcal nuclease or MNase is a 16.9 kDa endonuclease derived from Staphylococcus aureus. It is purified from an E. coli strain expressing an N-terminal 6XHIS tagged micrococcal nuclease. Purified protein exhibit a strong endonuclease activity against single-stranded, double-stranded, circular and linear nucleic acids. The enzyme is active in the pH range of 7.0 - 10.0, with optimal activity at pH 9.2 for both RNA and DNA substrates. The rate of cleavage is 30 times greater at the 5' side of A or T than at G or C and results in the production of mononucleotides and oligonucleotides with terminal 3'-phosphates. MNase is suitable for removing nucleic acids from cell lysates, releasing chromatin-bound proteins and shearing chromatin for use in chromatin immunoprecipitation (ChIP) experiments.

### Quality control



**Figure 1.**

SDS page of the micrococcal MNase. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.



**Figure 2.**

MNase activity assay. 0.5  $\mu$ g plasmid DNA was digested with the Diagenode MNase (concentration 5U/ml) in a buffer containing 20 mM Tris-HCl Ph 7.6, 3 mM CaCl<sub>2</sub> and 0.01% BSA and analyzed by agarose gel electrophoresis. The digestion was carried out for 0.5, 1, 2 and 4 minutes (lane 2, 3, 4 and 5, respectively). An undigested control is shown in lane 1.

**Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

**Diagenode Inc. USA | NORTH AMERICA**

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**Last update:** September 10, 2015