

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

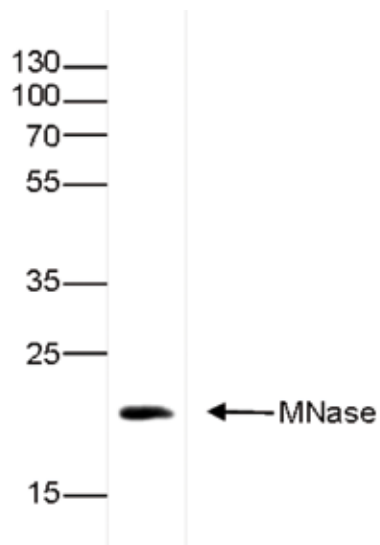


Figure 1.

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.

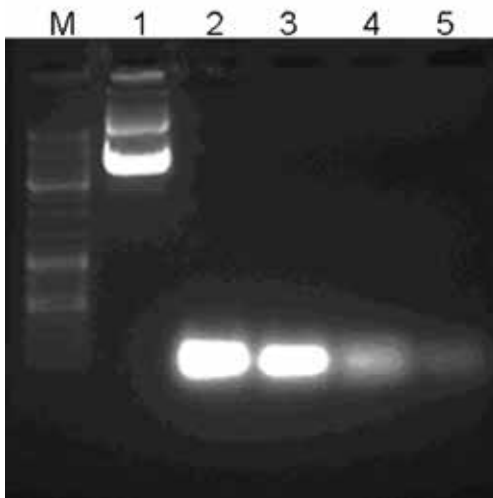


Figure 2.

MNase activity assay. 0.5 µg plasmid DNA was digested with the Diagenode MNase (concentration 5U/ml) in a buffer containing 20 mM Tris-HCl Ph 7.6, 3 mM CaCl₂ 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