

TurboNuclease

Recombinant form of *Serratia marcescens* extracellular endonuclease

Conc.: 50,000 units @ 250 units/μl

Formulated in: Storage Buffer of 50 mM Tris-HCl, pH 8.0, 50 mM NaCl, 5 mM MgCl₂ and 50% glycerol.

Endotoxin Level: Total endotoxin level is <0.1EU/1,000 units of TurboNuclease as determined by the Endosafe® PTS™ LAL test system. No protease activity was detected.

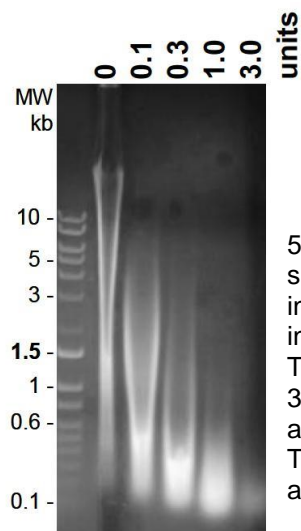
Stability: >6 months at -20°C

Description: Ultra-pure version of Benzonase. Nonspecific endonuclease, hydrolyzes both single and double-stranded nucleic acids (DNA and RNA) to 5'-phosphorylated oligonucleotides of 1-4 bases in length. MW = 27 kDa (homodimer). Expressed in *E. coli*.

Specific Activity: >1.3x10⁶ units/mg (>3x10⁶ Kunitz units/mg). One unit converts 1 OD260 of salmon sperm DNA into acid soluble nucleotides in 30 minutes at 37°C in a reaction buffer of 50 mM Tris-HCl, pH 8.0 and 1 mM MgCl₂. No detectable protease activity.

Application: Ideal for digesting nucleic acids during virus preparation and for reducing cell lysate viscosity during protein purification.

Quality Assurance



50 μg of salmon sperm DNA was incubated with the indicated units of TurboNuclease at 37°C for 30 min in a buffer of 50 mM Tris-HCl, pH 8.0 and 1 mM MgCl₂.

10% SDS-PAGE Coomassie staining

MW: 27 kDa
Purity: ≥99%

