

Turbo Nuclease, S pack

Ordering Information

[Order Online »](#)

| | |
|-------------|--------|
| JBS-EN-180S | 10000U |
|-------------|--------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Benzonase®

Application Notes Turbo Nuclease is very effective in degrading nucleic acid from protein samples and it is used in a variety of application where complete hydrolysis of DNA/RNA is required:
Reduction of viscosity from cell lysates
Prevention of cell clumping
Elimination of unspecific protein-DNA complexes prior 2D- or native gel electrophoresis
Removal of nucleic acids from large-scale protein preparations

Formulation liquid (Supplied in 20 mM Tris-HCl pH 8.0, 20 mM NaCl, 2 mM MgCl₂, 1 mM DTT and 50 % [v/v] glycerol)

Source Serratia marcescens, recombinant, E. coli

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com