

RNase I (DNase free), S pack

Endonuclease

Ordering Information

[Order Online »](#)

JBS-EN-176S	2000U
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes Degrades RNA to cyclic nucleotide monophosphates leaving a 5'-OH and 2', 3'-cyclic monophosphate
Ribonuclease Protection Assays
Mapping or Quantitation of RNA by selective cleavage of single-strand regions
Eliminates RNA from DNA- and protein purification

Formulation liquid (Supplied in 10 mM Tris-HCl pH 8.0, 200 mM NaCl and 50 % [v/v] glycerol)

MW 27.0 kDa

Purity > 95 % (SDS-PAGE)

Source 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](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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