

RNase A (DNase free), L pack

Ordering Information

[Order Online »](#)

JBS-EN-179L	2x250mg
-------------	---------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Endoribonuclease

Appearance dry powder

Application Notes Isolation of DNA Plasmid Preparation Analytical purposes

CAS 9001-99-4

MW 13.7 kDa (monomer)

Purity ≥ 90 % (ion exchange chromatography), salt free, chromatographically homogeneous lyophilisate

Source Bovine pancreas

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