

RNase Inhibitor – recombinant, L pack

RNase inhibitor is a recombinant protein which completely inhibits a broad spectrum of eukaryotic RNases, including RNase A, B and C. It inhibits RNases by binding noncovalently in a 1:1 ratio with high affinity ($4 \times 10^{-14}M$). It does not inhibit the RNases I, T1, T2, H, U1, U2 and C13. In addition, RNase inhibitor shows no inhibition of polymerase or reverse transcriptase activity and so can be used for cDNA synthesis and in one-step RT-PCR reactions. The murine version of RNase inhibitor lacks the pair of cysteines identified in human version, therefore it has significantly improved resistance to oxidation. Nuclease, DNase and RNase free

Ordering Information

[Order Online »](#)

JBS-PCR-392L	5x2000U
--------------	---------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes RNase inhibitor should be applied in a working concentration of 40 units (1 µl) per 50 µl reaction volume. The application of RNase inhibitor is recommended for:

- cDNA synthesis / RT-PCR
- In vitro transcription/translation
- RNA purification
- RNA protection assays
- Separation and identification of specific ribonuclease activities
- Other applications where the integrity of RNA is essential

Concentration 40 units/µl

Formulation liquid (Supplied in 20 mM HEPES-KOH pH 7.5, 50 mM KCl, 8 mM DTT and 50 % [v/v] glycerol)

Source Mus musculus, recombinant

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)