

Poly (A) carrier RNA, L pack

Poly (A) carrier RNA is used for quantitative precipitation/purification of RNA and DNA. It improves recovery of short fragments (< 200 bp) or low amounts of nucleic acids. Chain length: 2100 to 10000 nucleotides
Recommended working concentration: 0.5 mg/ml

Ordering Information

[Order Online »](#)

| | |
|---------------|----|
| JBS-RNT-017-L | 1g |
|---------------|----|

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 Polyadenylic acid, potassium salt

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



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)