

AMPIGENE[®] RNase Inhibitor

Recombinant protein that blocks ribonuclease activity to protect samples from RNase digestion

AMPIGENE[®] RNase Inhibitor works to inhibit RNase activity for sensitive applications such as RT-qPCR, cDNA synthesis, and RNA sequencing.

Ordering Information

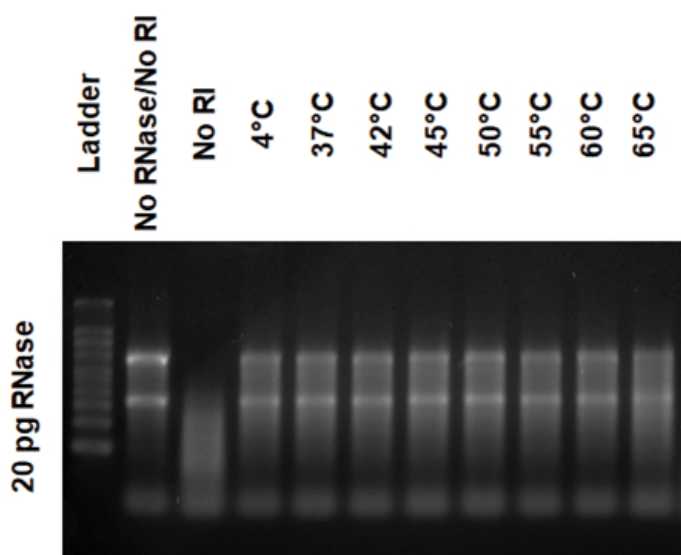
[Order Online »](#)

ENZ-NUC136-2500	2500U
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Inhibits eukaryotic RNases, including RNase A, B, and C
- Compatible with reverse transcriptase, Taq DNA polymerase, and RNA polymerase
- Ribonuclease and phosphatase free
 - Stable up to 65°C for up to 30 minutes



AMPIGENE® RNase Inhibitor is thermostable up to 65°C for 30 minutes

AMPIGENE® RNase Inhibitor (40U) was incubated at the indicated temperatures for 30 minutes. After the incubation, the enzyme was added to the mix containing 1 ug RNA, 20 pg RNase, and reaction buffer (1x final concentration). Samples were then incubated at 37°C for 30 min (to allow RNase to degrade RNA, if active), then loaded on a 1% agarose gel. AMPIGENE® RNase Inhibitor was stable up to 65°C ensuring RNA protection from RNase degradation.

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application PCR

Contents AMPIGENE[®] RNase Inhibitor (40U/μL)



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)