

# AMPIGENE<sup>®</sup> RNase Inhibitor

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

AMPIGENE<sup>®</sup> RNase Inhibitor works to inhibit RNase activity for sensitive applications such as RT-qPCR, cDNA synthesis, and RNA sequencing.

## Ordering Information

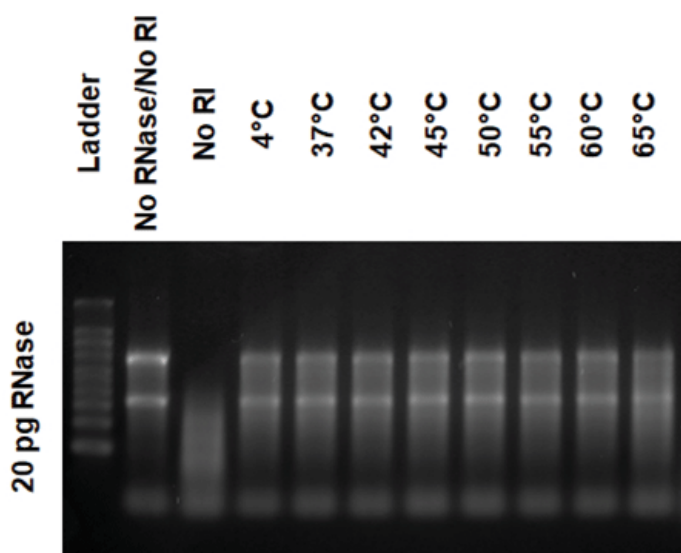
[Order Online »](#)

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<sup>®</sup> RNase Inhibitor (40U/μL)

Last modified: May 29, 2024



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)