

# Recombinant RNasin® Ribonuclease Inhibitor

Don't see what you are looking for?  
[Click here for custom options.](#)



RNases are ubiquitous, cause RNA degradation and can compromise RNA integrity. Recombinant RNasin® Inhibitor is a 50kDa protein that inhibits RNase A family and human placental RNases by noncovalently binding to RNases in a 1:1 ratio. Recombinant RNasin® Inhibitor does not inhibit RNase T1, S1 nuclease, RNase from *Aspergillus*, RNase H, RNase ONE™ Ribonuclease and enzymes for downstream applications such as GoScript™ Reverse Transcriptase, AMV/M-MLV reverse transcriptases, SP6, T7/T3 RNA polymerase, and *Taq* DNA polymerases. Learn more about our custom options for this product at: [www...](#)

## Expand to Read More »



[Learn more about the Helix program](#)



Prices valid for US customers only

PRODUCT	SIZE	CONC.	CATALOG #	*LIST PRICE	QTY	
Recombinant RNasin® Ribonuclease Inhibitor	2,500u	20–40u/ μl	N2511	\$ 114.00	<input type="text"/>	<a href="#">Add to cart</a>
Protocol <a href="#">MSDS</a> <a href="#">Components</a> <a href="#">Add to Helix</a>						
Recombinant RNasin® Ribonuclease Inhibitor	10,000u	20–40u/ μl	N2515	\$ 277.00	<input type="text"/>	<a href="#">Add to cart</a>
Protocol <a href="#">MSDS</a> <a href="#">Components</a> <a href="#">Add to Helix</a>						

### QC Tests

SDS-PAGE/purity, DNase, RNase, latent RNase, endonuclease.

### Source

Recombinant strain of *E. coli* (Cat.# N2511, N2515).

### Storage Buffer

20mM HEPES-KOH (pH 7.6 at 4°C), 50mM KCl, 8mM DTT and 50% (v/v) glycerol.

### Storage Conditions

Store at -20°C.

### Unit Definition

One unit is defined as the amount of RNasin® Ribonuclease Inhibitor required to inhibit the activity of 5ng of ribonuclease A by 50%. Activity is measured by the inhibition of hydrolysis of cytidine 2,3'-cyclic monophosphate by ribonuclease A.

For product intended use please see Patents & Disclaimers tab.

### You Also May Be Interested In



[GoTaq® Real-Time qPCR and RT-qPCR Systems for Dye-Based Detection](#)

[GoScript™ Reverse Transcription System](#)

[QuantiFluor® RNA System](#)

[M-MLV Reverse Transcriptase, RNase H Minus, Point Mutant](#)