

# RIG-I (human) polyclonal antibody (AT111)

RIG-I (retinoic acid-inducible gene I; Ddx58), a member of the DEAD-box RNA helicases, is activated by cytoplasmic dsRNA and 5'-ppp RNA produced during the viral replication. The protein is characterized by a N-terminal region with two caspase recruitment domains (CARD) and a C-terminal region harboring potential ATP-dependent RNA helicase activity. RIG-1 recruits the CARD adaptor inducing IFN- $\beta$  (Cardif) in a CARD-CARD-dependent manner resulting in NF- $\kappa$ B and IRF3 activation.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 10

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-210-932-C100	100 $\mu$ g
------------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	RIG-1, Retinoic acid-inducible gene 1 protein, DEAD-box protein 58, DEAD (Asp-Glu-Ala-Asp) box polypeptide 58 M, Probable ATP-dependent RNA helicase DDX58
Application	IP, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Recombinant human RIG-I (aa 201-713).
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (1:100)Western blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Human
UniProt ID	O95786
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



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)