

RIG-I (mouse) monoclonal antibody (SS1A)

RIG-1/DDX58 (Retinoic acid-inducible gene 1 protein/ DEAD-box protein 58) is an RNA helicase that belongs to the DEAD/H box. Contains 2 CARD domains, a helicase ATP-binding domain and a helicase C-terminal domain. Predicted molecular weight approximately 107 kDa. RIG-1 has an essential function in double-stranded RNA-induced innate antiviral responses. RIG-1 and MDA5 (melanoma differentiation-associated gene 5) are key intracellular receptors that sense intracellular viral infection and trigger a signal for innate antiviral responses including the production of type I IFN. Clone SS1A has been shown to be useful for Western blotting of the mouse RIG-1 protein. Involved in innate immune defense against viruses. RIG-1 has an essential function in double-stranded RNA-induced innate antiviral responses.

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

Citations: 3

[View Online »](#)

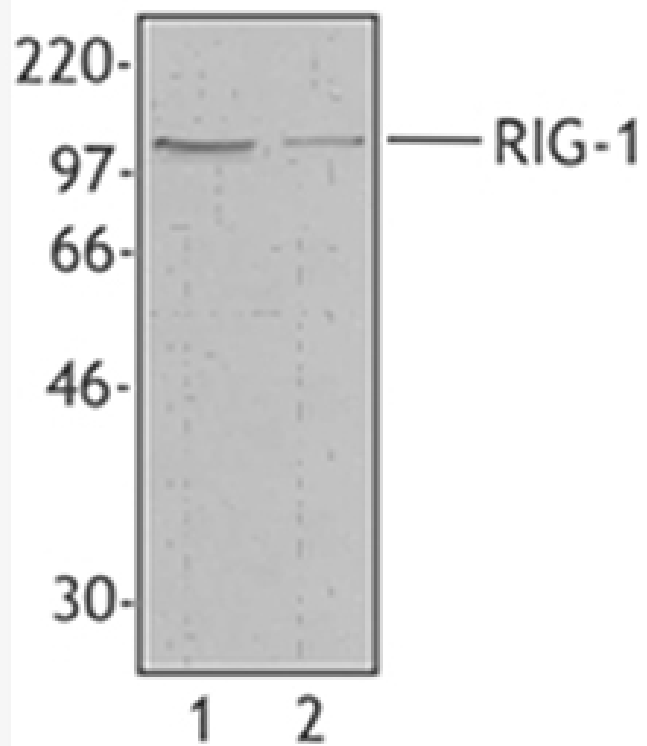
Ordering Information

[Order Online »](#)

ALX-804-960-0100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of RIG-I (mouse), mAb (SS1A) (Prod. No. ALX-804-960) using NIH-3T3 cells extracts (lane 1) and mouse splenocytes (lane 2).

Handling & Storage

Long Term Storage +4°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	WB
Clone	SS1A
Formulation	Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.
Gene/Protein Identifier	109906 (Entrez GeneID)
Host	Rat
Immunogen	Synthetic peptide corresponding to C-terminal mouse RIG-I conjugated to KLH.
Isotype	IgG2a
Purity Detail	Affinity purified.
Species Reactivity	Mouse
UniProt ID	Q6Q899
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)