## 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 typel 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** 

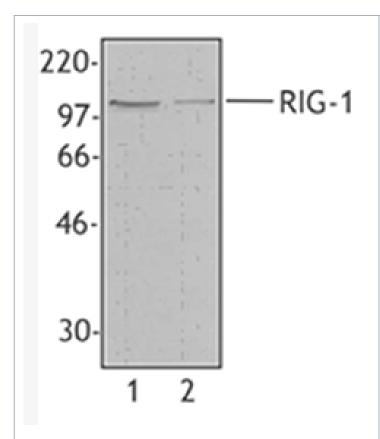
**Order Online** »

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.

lgG2a

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com