

# Cardif/MAVS (human) monoclonal antibody (58N3B6)

RIG-I (retinoic acid-inducible gene I; Ddx58) and MDA5 (melanoma differentiation-associated gene 5, also known as Ifih1 or Helicard) are proteins that sense viral replication intermediates, such as double-stranded RNA and triggers the host antiviral programs. These molecules signal the downstream activation of NF- $\kappa$ B and IFN regulatory factor (IRF) -3, which coordinately regulate the expression of type-I interferons. Cardif (also called VISA/IPS-1/MAVS) is a CARD (caspase activation and recruitment domain)-containing adaptor protein that interacts with the CARD domain of RIG-I and MDA5, leading to the activation of NF- $\kappa$ B and IRF3. Cardif is located to the mitochondrial outer membrane. Removal of the mitochondrial-targeting domain of cardif abolishes its ability to induce IFNs. Cardif is cleaved and inactivated by NS3-4A, a serine protease from hepatitis C virus known to block interferon- $\beta$  production.

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

Citations: 3

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-ABS259-0100	100 $\mu$ g
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

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	CARD adapter inducing interferon-β, IPS-1, Interferon-β promoter stimulator protein 1, Mitochondrial antiviral signalling protein, VISA, Virus-induced signalling adapter
Application	IHC (PS), WB
Clone	58N3B6
Formulation	Liquid. In 0.2ml PBS containing 0.05% sodium azide.
Gene/Protein Identifier	57506 (Entrez GeneID)
Host	Mouse
Immunogen	Human MAVS (IPS-1)
Isotype	IgG2bk
Purity Detail	Protein G purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:10-1:2,000) (paraffin sections 3-5µg/ml)Western Blot (2-5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q7Z434

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-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)