

# MDA5 (human) monoclonal antibody (Hely-1)

MDA5 and RIG-I are highly conserved helicases involved in the innate immune response to virus. MDA-5 ATPase activity can be stimulated by RNA and less potently by DNA. It contains two N-terminal CARD domains and a C-terminal helicase domain. After binding nonself-ligand, RIG-I and MDA5 interact via their CARD domains with the signaling-adaptor molecule, IPS-1 (also known as MAVS, VISA and Cardif), which leads to the induced expression of IFN- $\beta$ , IRF3-target genes and NF- $\kappa$ B target genes to drive antiviral and inflammatory responses.

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

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-863-C100

100 $\mu$ g

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Helicard, Melanoma differentiation-associated gene 5, Interferon-induced helicase C domain-containing protein 1, RH116, CADM-140 autoantigen, RIG-I-like receptor 2, RLR-2
Application	ELISA, IP, WB
Application Notes	ELISA Immunoprecipitation (1:100) Western Blot (1:1'000)
Clone	Hely-1
Formulation	Liquid. In PBS containing 10% glycerol and 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human MDA5 (melanoma differentiation-associated gene 5) (aa 78-555).
Isotype	IgG1
Purity Detail	Protein G-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.

<b>Source</b>	Purified from concentrated hybridoma tissue cell culture supernatant.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	Q9BYX4
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

**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)