

MDA5 (human) polyclonal antibody (AT113)

MDA5 is a highly conserved helicase whose 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. The helicase domain, once freed from the CARD domains, then translocates from the cytoplasm to the nucleus where it acts on chromatin architecture. This could allow a more easy access of CAD, thereby accelerating the degradation of genomic DNA.

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

Citations: 25

[View Online »](#)

Ordering Information

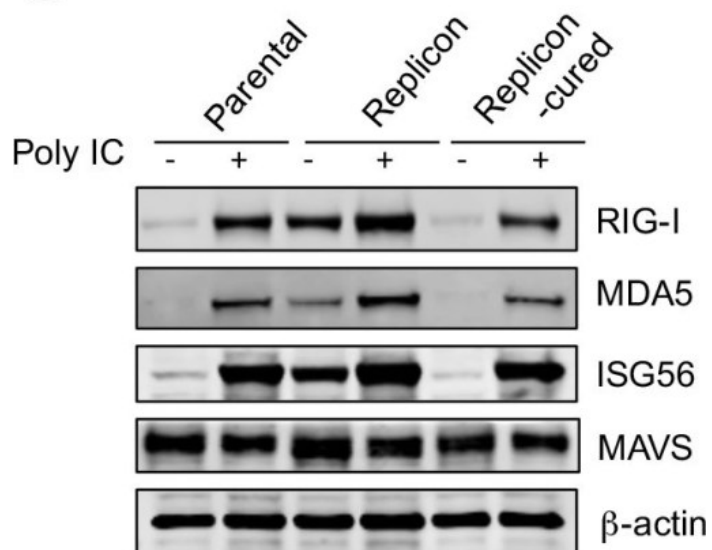
[Order Online »](#)

ALX-210-935-C100	100µg
------------------	-------

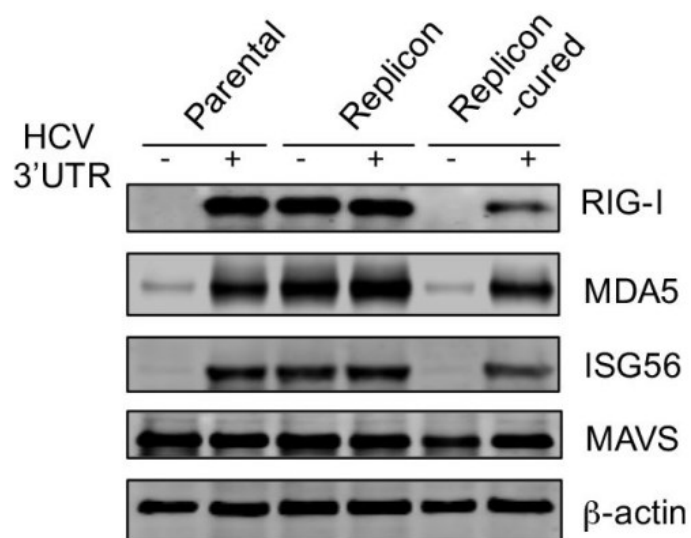
Manuals, SDS & CofA

[View Online »](#)

G



J



HEV does not target MAVS. (A) Confocal images showing MAVS and viral antigens in HepG2 cells infected with either HEV (top) or HAV (bottom). Cells were stained 5 days after infection with a rabbit anti-MAVS, chimpanzee 1313 serum (HEV), or a murine monoclonal antibody K24F2 (HAV). DAPI was used to stain the nuclei. Scale bar: 10 μ m. (B) Confocal images showing the mitochondrial localization of MAVS in HepG2 cells with or without HEV replicon. MAVS was stained with a rabbit antibody against MAVS (green). Mitochondria were visualized with MitoTracker (red). Nuclei were stained with DAPI. Scale bar: 10 μ m. (C) HepG2 cells with or without the HEV replicon were transfected with a MAVS-expressing plasmid along with a HAV 3ABC-expressing plasmid or an empty vector. The endogenous (closed arrowheads) and overexpressed MAVS (open arrowheads) were detected with a rabbit anti-MAVS antibody. Note that co-

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored as recommended.
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	IP, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Recombinant human MDA5 (melanoma differentiation-associated gene 5) (aa 78-555).
Purity Detail	Epitope-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	Q9BYX4

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)