

Cardif (human) polyclonal antibody (AT107)

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- κ 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- κ 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- β production.

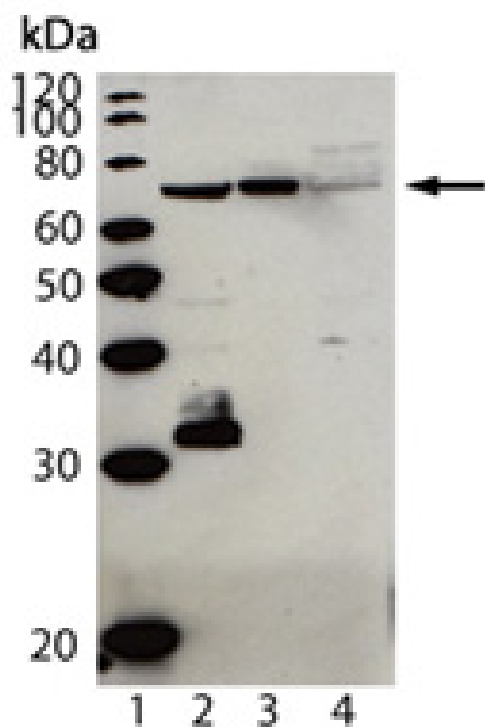
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 40 [View Online »](#)

Ordering Information [Order Online »](#)

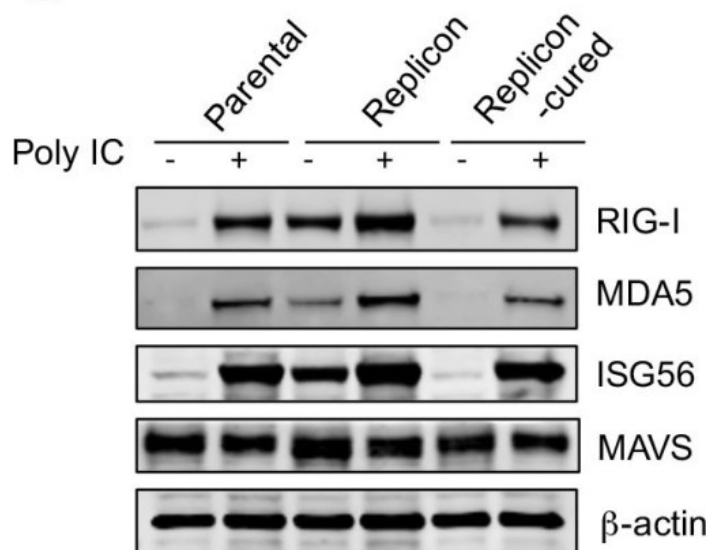
ALX-210-929-C100	100 μ g
------------------	-------------

Manuals, SDS & CofA [View Online »](#)

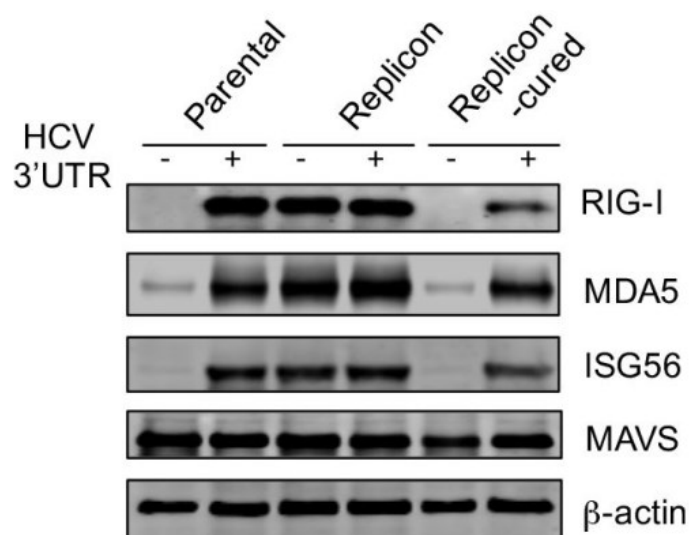


Western blot analysis of Cardif (human), pAb (AT107) (Prod. No. ALX-210-929): Lane 1: MW marker; Lane 2: HepG2; Lane 3: PALA; and Lane 4: HeLa. Additional bands observed probably represent isoforms or cleaved products of Cardif.

G

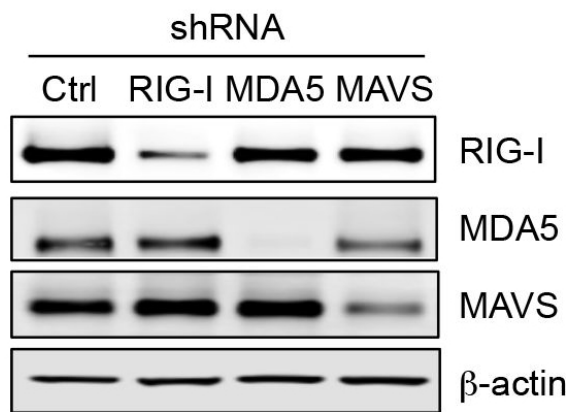


J



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 was 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-

A



Signaling pathways involved in HEV-induced IFN response. Immunoblots of RIG-I, MDA5, MAVS and β -actin in HepG2 cells expressing gene-specific shRNA, or GFP (Ctrl). (B-C) IFN- β promoter activity in HepG2 cells with different gene knockdown following Sendai virus (SeV) infection (B) or poly IC transfection (C). Cells were transfected with IFN- β -Luc and TK-RLuc (for normalization of transfection efficiency) 20 h prior to SeV infection or poly IC transfection. Cells were lysed and luciferase activity was determined 20 h after SeV infection or 12 h after poly IC transfection. Data are presented as fold changes relative to non-treated cells. Shown are representative results from two independent experiments each performed in triplicate. (D-F) Effect of RIG-I, MDA5 or MAVS knockdown on HEV replication and host IFN responses. Control and shRNA-expressing HepG2 cells were inoculated with HEV (1000 GE/cell). IFN- λ mRNA expression (D), HEV RNA abundance (E), and HEV-positive foci (F) were determined at 5 days after infection. The results show the mean \pm SEM of 2 independent experiments performed in duplicate each. * $P < 0.05$; ** $P < 0.01$. (G) Immunoblots of IRF-3, IRF-7 and β -actin in HepG2 cells transduced with lentiviruses expressing GFP (Ctrl) or gene-specific shRNA. (H-J) Effect of IRF-3 or IRF-7 knockdown on HEV replication and IFN responses. Control and shRNA-expressing cells were inoculated with HEV (1,000 GE/cell). IFN- λ mRNA expression (H), HEV RNA abundance (I), and HEV-positive foci (J) in different cells were determined after 5 days. The results show the mean \pm SEM of 2 independent experiments each performed in duplicate wells. * $P < 0.05$; ** $P < 0.01$. Scale bar (upper panel in F), 100 μ m.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Hepatitis E virus persists in the presence of a type III interferon response. *PLoS Pathog* (2017)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
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	CARD adapter inducing interferon- β , IPS-1, Interferon- β promoter stimulator protein 1, MAVS, Mitochondrial antiviral signalling protein, VISA, Virus-induced signalling adapter
Application	ICC, IP, WB
Application Notes	Detects a band of ~65kDa by Western blot.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Recombinant human Cardif (CARD adapter inducing interferon- β) (aa 160-450).
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:500)Immunoprecipitation (1:100)Western Blot (1:2,000)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

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)