

MDA5 polyclonal antibody

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: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-ABS299-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

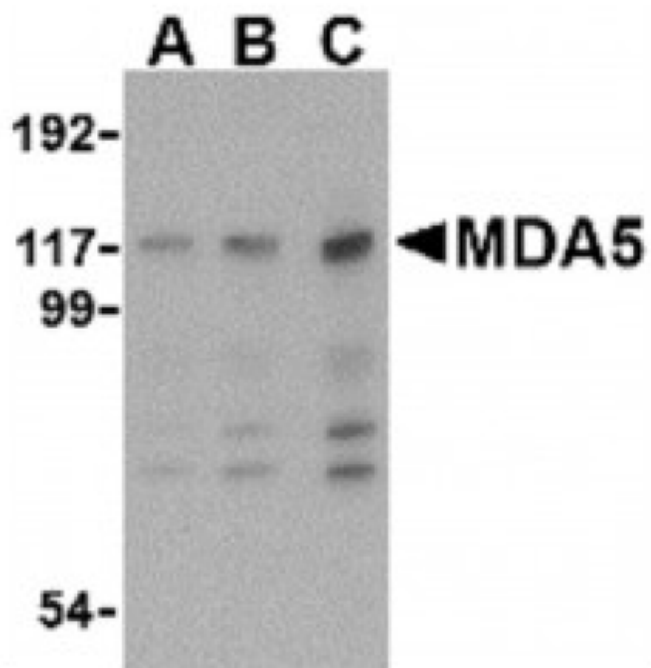


Figure 1: Western blot analysis of MDA5 in Daudi cell lysate with MDA5 antibody at (A) 1, (B) 2 and (C) 4 $\mu\text{g/mL}$.

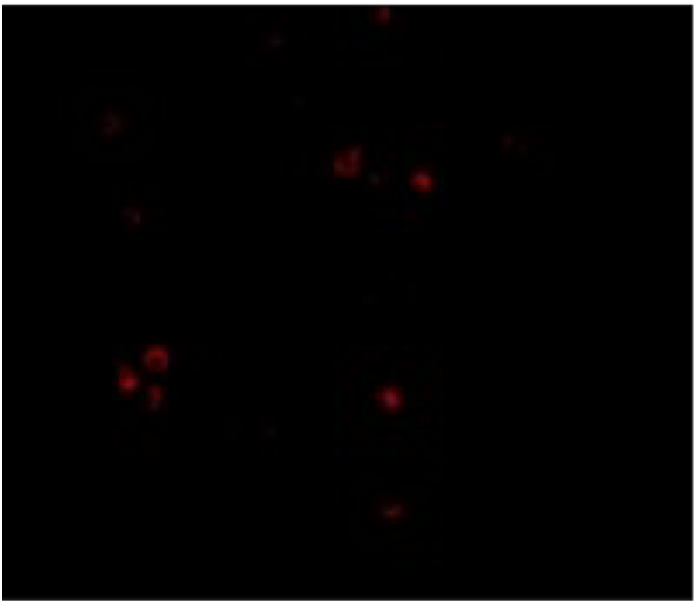


Figure 2: Immunofluorescence of MDA5 in Human Lymph Node cells with MDA5 antibody at 20 $\mu\text{g/mL}$.

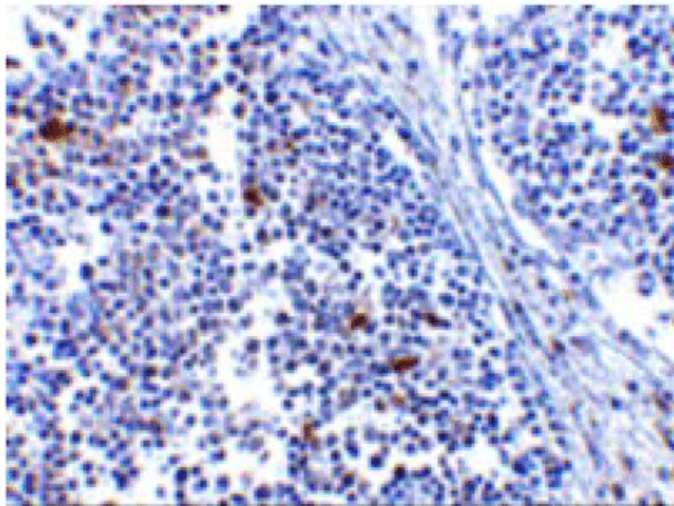


Figure 3: Immunohistochemistry of MDA5 in human lymph node tissue with MDA5 antibody at 5 $\mu\text{g/mL}$.

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	ELISA, IF, IHC (PS), WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to 16 amino acids from near the center of human MDA5.
Isotype	IgG
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Immunofluorescence (10µg/ml)Immunohistochemistry (5µg/ml)Western Blot (1-4µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	Q9BYX4 (human), Q8R5F7 (mouse)

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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