

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 »](#)

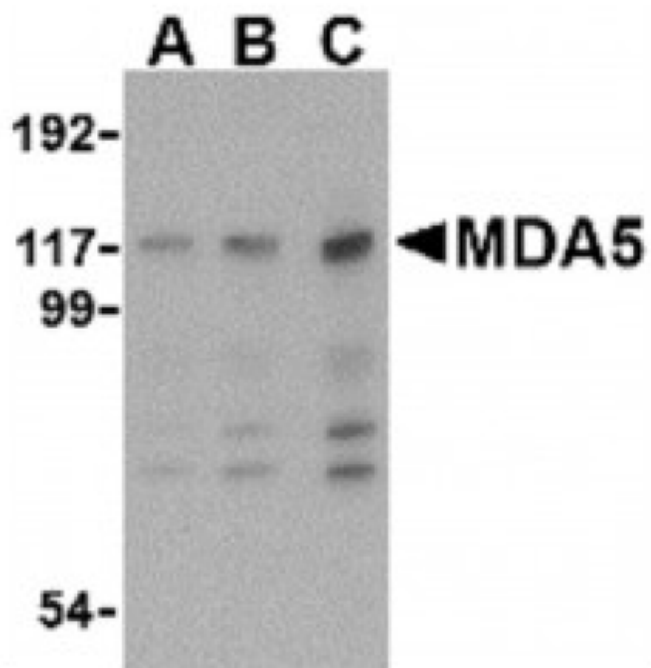


Figure 1: Western blot analysis of MDA5 in Daudi cell lysate with MDA5 antibody at (A) 1, (B) 2 and (C) 4 $\mu\text{g/mL}$.

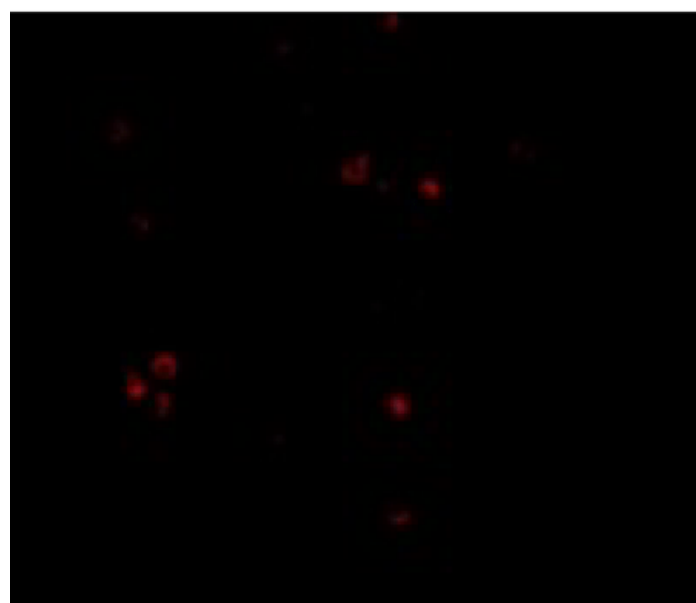


Figure 2: Immunofluorescence of MDA5 in Human Lymph Node cells with MDA5 antibody at 20 $\mu\text{g/mL}$.

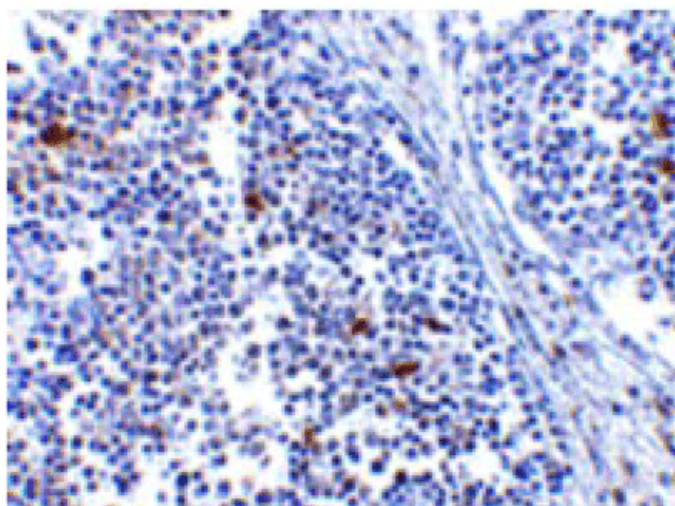


Figure 3: Immunohistochemistry of MDA5 in human lymph node tissue with MDA5 antibody at 5 $\mu\text{g/mL}$.

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)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
Info-
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELSI) 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

This antibody is covered by our **Worry-Free Guarantee**.