

VAMP8 polyclonal antibody

This polyclonal antibody detects Vesicle-Associated Membrane Protein 8 (VAMP8), a key regulator of vesicle trafficking.

VAMP8 is a SNARE protein that mediates intracellular vesicle fusion and trafficking. It plays a role in exocytosis, autophagy, and immune cell activation, making it an important target in studies of inflammation and neurodegeneration.

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

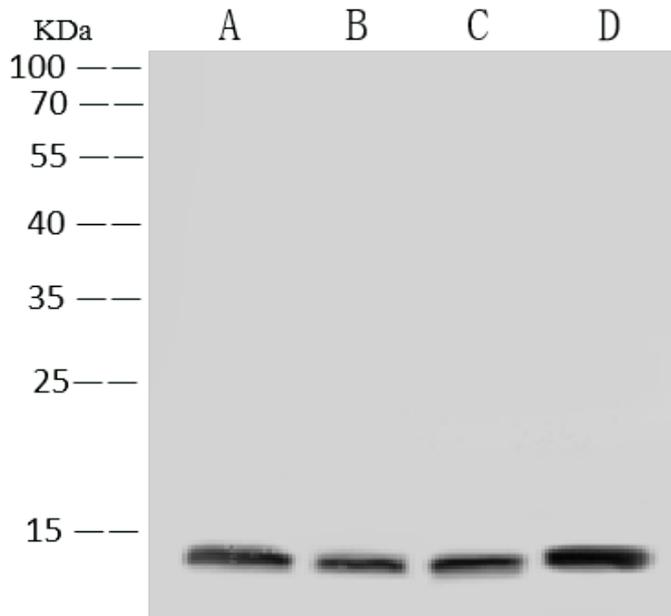
Ordering Information

[Order Online »](#)

ENZ-ABS1073-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of various lysates using VAMP8 rabbit polyclonal antibody (ENZ-ABS1073) at 1:500 dilution. Secondary Antibody: Goat Anti-Rabbit IgG (H+L)/HRP at 1:10'000 dilution. Lane A: A431 Whole Cell Lysate, Lane B: HEK293 Whole Cell Lysate, Lane C: Hela Whole Cell Lysate, Lane D: THP-1 Whole Cell Lysate. Lysates/proteins at 30 μ g per lane.

Handling & Storage

Handling Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Vesicle-associated membrane protein 8, Endobrevin

Application IP, WB

Formulation Liquid. In PBS, pH 7.0, with 0.03% Proclin300. Carrier-free.

Host Rabbit

Immunogen E. coli-derived human VAMP8 fragment.

Isotype IgG

Purity Detail Protein A & antigen affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:500-1:2'000)
Immunoprecipitation (1-5 µL/mg of lysate)

Suggested dilutions/conditions may not be available for all applications.

Optimal conditions must be determined individually for each application.

Species Reactivity Human

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



ENZO LIFE SCIENCES,
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