

VDAC2 polyclonal antibody

This antibody detects Voltage-Dependent Anion Channel 2 (VDAC2), which is involved in mitochondrial permeability and apoptosis regulation.

VDAC2 is a mitochondrial outer membrane protein involved in metabolic regulation and apoptosis. It interacts with apoptotic proteins like Bcl-2 and Bax, making it an important target in cancer and neurodegenerative diseases.

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

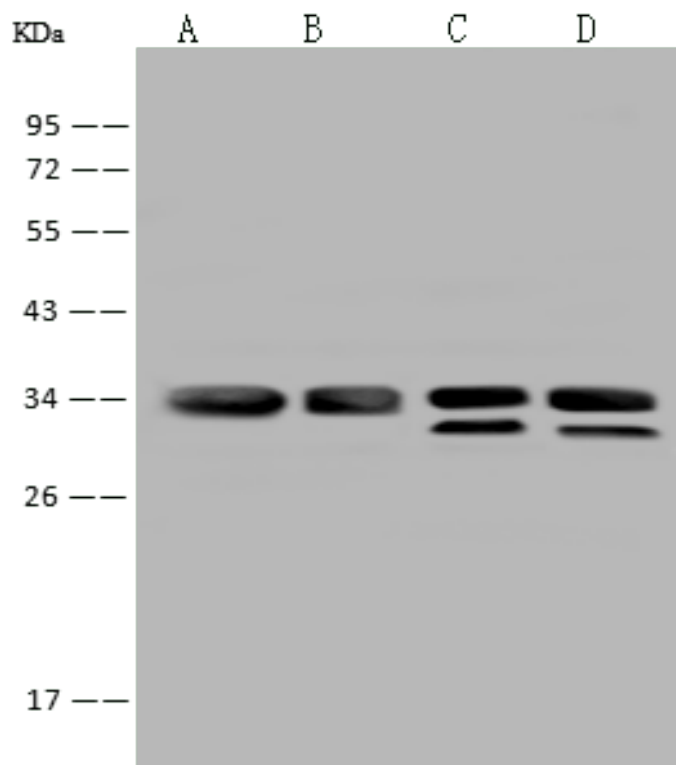
Ordering Information

[Order Online »](#)

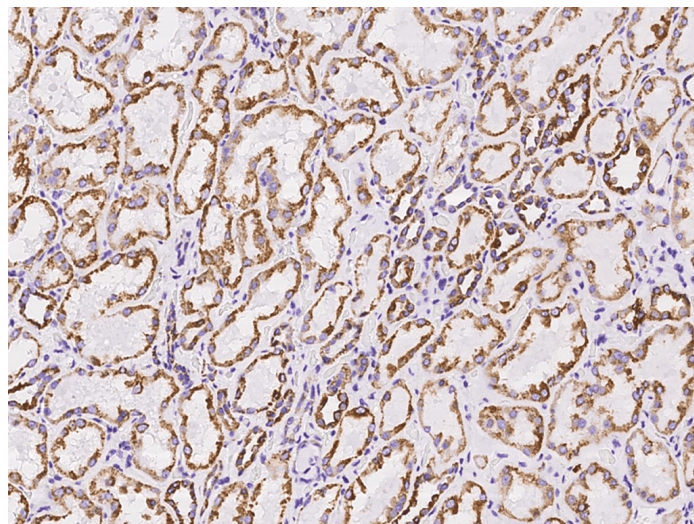
ENZ-ABS1072-0100	100µl
------------------	-------

Manuals, SDS & CofA

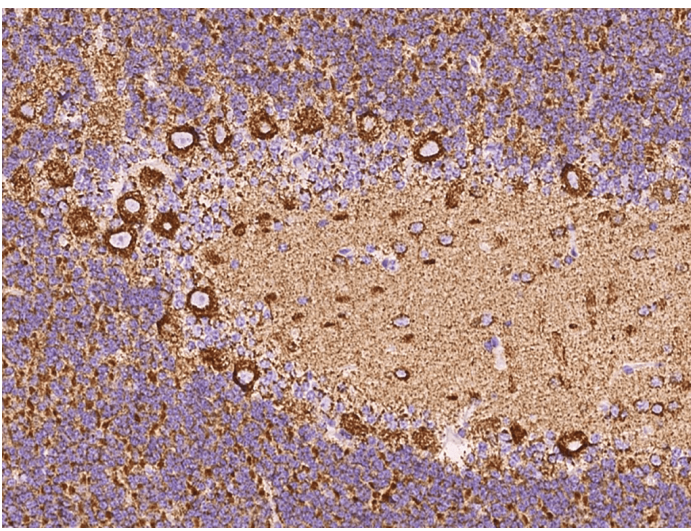
[View Online »](#)



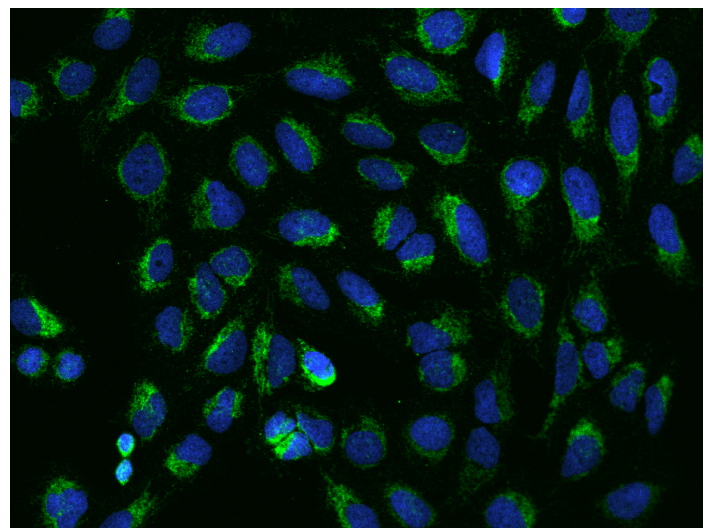
Western blot analysis of various lysates using VDAC2 rabbit polyclonal antibody (ENZ-ABS1072) at 1:500 dilution. Secondary Antibody: Goat Anti-Rabbit IgG (H+L)/HRP at 1:10'000 dilution. Lane A: HeLa Whole Cell Lysate, Lane B: HepG2 Whole Cell Lysate, Lane C: HCT116 Whole Cell Lysate, Lane D: U-251 MG Whole Cell Lysate. Lysates/proteins at 30 µg per lane.



Immunohistochemical staining of human VDAC2 in human kidney with VDAC2 rabbit polyclonal antibody at 1:100 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of human VDAC2 in cynomolgus cerebellum with rabbit polyclonal antibody at 1:100 dilution, formalin-fixed paraffin embedded sections.



Immunofluorescence staining of VDAC2 in U2OS cells. Cells were fixed with 4% PFA, permeabilized with 0.1% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit VDAC2 polyclonal antibody (dilution ratio 1:200) at 4°C overnight. Then cells were stained with the Alexa Fluor®488-conjugated Goat Anti-rabbit IgG secondary antibody (green) and counterstained with DAPI (blue). Positive staining was localized to Cytoplasm.

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 Non-selective voltage-gated ion channel VDAC2

Application ICC, IF, IHC (PS), WB

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

Host Rabbit

Immunogen E. coli-derived human VDAC2 fragment.

Isotype IgG

Purity Detail Protein A & antigen affinity purified.

Recommendation Dilutions/Conditions

Western Blot (1:500-1:2'000)
Immunocytochemistry/Immuofluorescence (1:100-1:500)
Immunohistochemistry (paraffin sections) (1:50-1:200)

Suggested dilutions/conditions may not be available for all applications.
Optimal conditions must be determined individually for each application.

Species Reactivity Cynomolgus, Human

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

.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
Last modified: **March 28, 2025**
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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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