

VDAC1 polyclonal antibody

This antibody detects Voltage-Dependent Anion Channel 1 (VDAC1), a key protein involved in mitochondrial function and apoptosis.

VDAC1 is a mitochondrial outer membrane protein that regulates metabolite exchange between the mitochondria and cytoplasm. It plays a crucial role in cell survival, apoptosis, and metabolic regulation. VDAC1 is associated with cancer, neurodegeneration, and metabolic disorders.

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

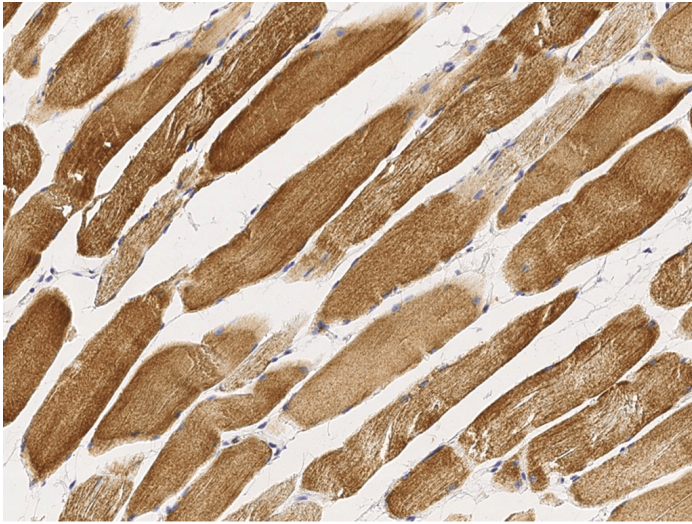
Ordering Information

[Order Online »](#)

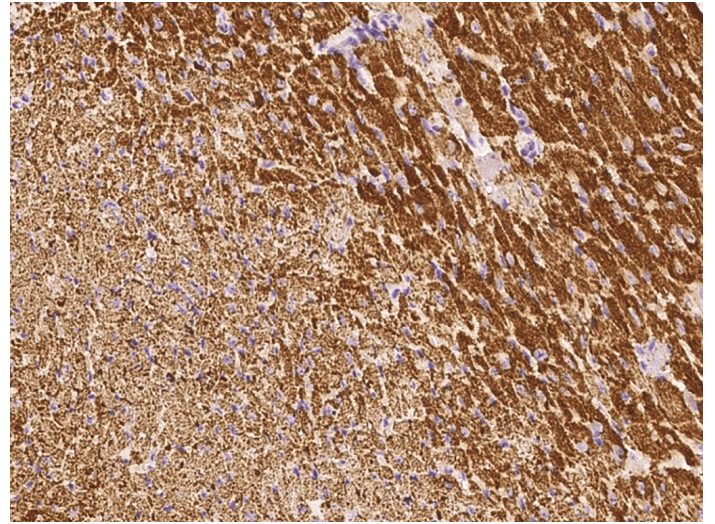
ENZ-ABS1065-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining VDAC1 in human skeletal muscle at 1:2'000 dilution, formalin-fixed paraffin embedded sections.



Immunohistochemical staining of VDAC1 in mouse heart at 1:1'000 dilution, formalin-fixed paraffin embedded sections.

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 VDAC1

Application IHC (PS), IP, WB

Crossreactivity Cross-reacts with human VDAC2 and VDAC3, mouse VDAC1 and VDAC3 and rat VDAC1.

Formulation Liquid. 0.2 µm filtered solution in PBS. Carrier-free. Preservative-free.

Host Rabbit

Immunogen Synthetic peptide corresponding to the center region of human VDAC1.

Isotype IgG

Purity Detail Protein A & antigen affinity purified.

Recommendation Western Blot (1:500-1:2'000)

Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:500-1:3'000)

Immunoprecipitation (1-4 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, Mouse

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



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)