

# 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](#).

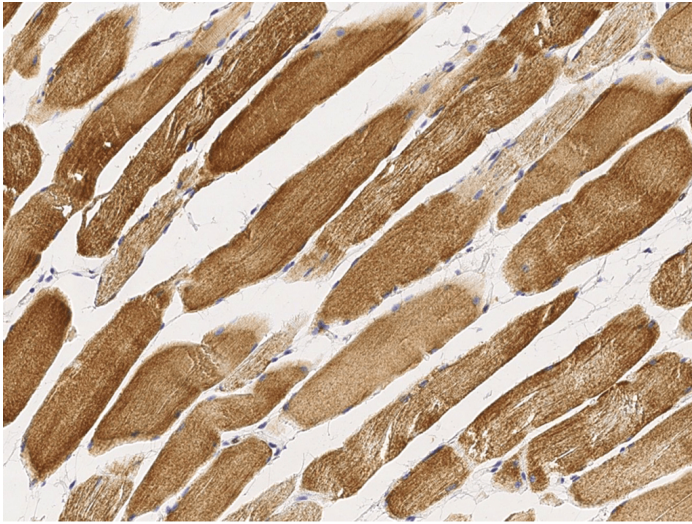
## Ordering Information

[Order Online »](#)

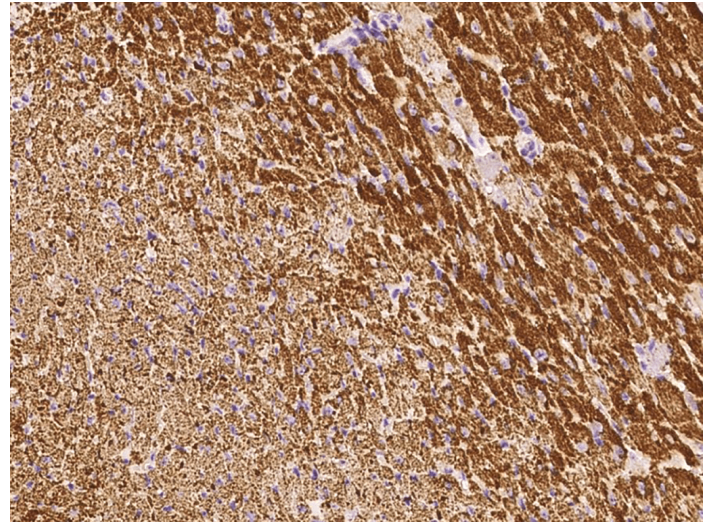
ENZ-ABS1065-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



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



Immunochemical 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 Dilutions/Conditions	Western Blot (1:500-1:2'000) 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 <a href="#">Worry-Free Guarantee</a>



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)