Voltage-dependent anion channel polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

Ordering Information

Order Online »

ALX-210-785-C100

100µg

Manuals, SDS & CofA

View Online »

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 VDAC, Mitochondrial porin

Application IP, WB

Application NotesDetects a band of ~31kDa by Western blot.

Formulation Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium

azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to a part

(CNDGTEFGGSIYQK) of C-terminal human VDAC (voltage-dependent anion channel) with an added N-terminal cysteine. This sequence is completely conserved between isoforms 1, 2 and 3 of human, mouse and rat

VDAC.

Purity Detail Epitope-affinity purified IgG.

Recommendation Dilutions/Conditions Western Blot (2µg/ml)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Monkey, Mouse, Rat

UniProt ID P21796

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com