

VAMP2 polyclonal antibody

Vesicle-associated membrane protein (VAMP), also known as synaptobrevin, is an 18 kDa integral membrane protein localized to the cytoplasmic surface of synaptic vesicles. Two VAMP homologs, VAMP1 and VAMP2, are differentially expressed in the nervous system. VAMP2 is found in neurons associated with autonomic, sensory and integrative functions, and in non-neuronal Langerhans islets and glomerular cells. VAMP is a component of a 20S complex implicated in vesicle docking and fusion, the formation of which involves the interaction of VAMP with syntaxin through an ATP-dependent reaction modulated by synaptotagmin, SNAP-25, alpha-SNAP, and NSF.

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

Citations: 8

[View Online »](#)

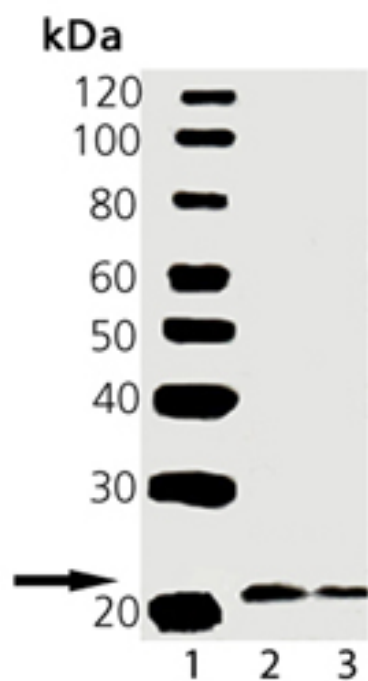
Ordering Information

[Order Online »](#)

ADI-VAS-SV006-D	50µg
ADI-VAS-SV006-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of VAMP2, pAb: Lane 1: Molecular Weight Marker, Lane 2: Mouse Brain Tissue Extract (Prod. No.: ADI-LYT-MB100), Lane 3: Rat Brain Tissue Extract (Prod. No.: ADI-LYT-RB100)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Synaptobrevin, Vesicle-associated membrane protein 2

Application ICC, IP, WB

Application Notes Detects a band of ~18kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

GenBank ID M24105

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of rat VAMP2. The sequence is completely conserved in mouse, human, bovine, and Xenopus.

Purity Detail Protein A affinity purified.

Recommendation Western Blot (1:1,000, ECL) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Guinea pig, Human, Mouse, Rabbit, Rat

UniProt ID P63045

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



ENZO LIFE SCIENCES,
INC.
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