

# 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 »](#)

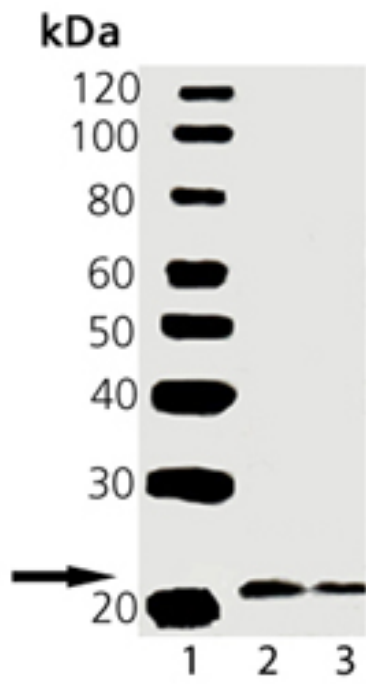
Ordering Information

[Order Online »](#)

ADI-VAS-SV006-D	50µg
-----------------	------

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

<b>Alternative Name</b>	Synaptobrevin, Vesicle-associated membrane protein 2
<b>Application</b>	ICC, IP, WB
<b>Application Notes</b>	Detects a band of ~18kDa by Western blot.
<b>Formulation</b>	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
<b>GenBank ID</b>	M24105
<b>Host</b>	Rabbit
<b>Immunogen</b>	Synthetic peptide corresponding to a portion of rat VAMP2. The sequence is completely conserved in mouse, human, bovine, and Xenopus.
<b>Purity Detail</b>	Protein A affinity purified.
<b>Recommendation Dilutions/Conditions</b>	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.
<b>Source</b>	Purified from rabbit serum.
<b>Species Reactivity</b>	Bovine, Guinea pig, Human, Mouse, Rabbit, Rat
<b>UniProt ID</b>	P63045
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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)