

SAP97 monoclonal antibody (RPI 197.4)

Synapse-associated proteins SAP97 and SAP102 display extensive sequence homology to the PSD 95 family of proteins that facilitate ion channel clustering at synaptic terminals. In addition to the CNS, SAP97 is also detected at the basal lateral membrane between a variety of epithelial cells. SAP102 associates with NMDA receptors through the C-terminal NR2B subunit of the receptor. The N-terminal SAP97 PDZ domain mediates its interaction with the synaptic ras-GTPase activating protein SynGAP.

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

Citations: 9

[View Online »](#)

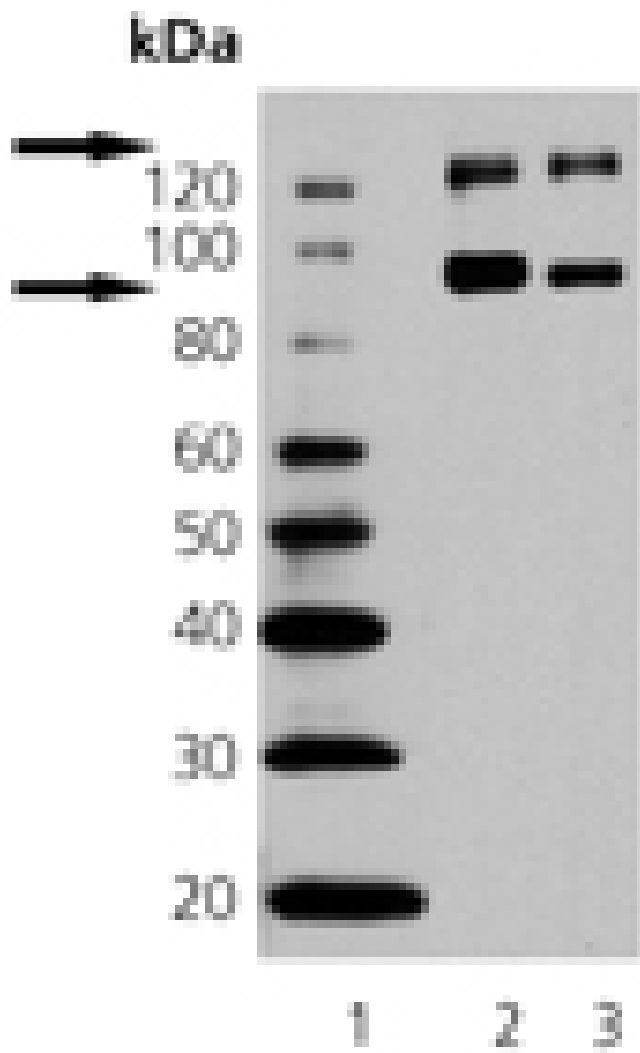
Ordering Information

[Order Online »](#)

ADI-VAM-PS005-D	50µg
ADI-VAM-PS005-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: Lane 1: MW marker, Lane 2: Rat Brain Tissue Extract, Lane 3: Mouse Brain Tissue Extract.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Synapse-associated protein 97, Disks large homolog 1
Application	Electron microscopy, ICC, IHC, IP, WB
Application Notes	Detects bands of ~97kDa and ~140kDa by Western blot.
Clone	RPI 197.4
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Host	Mouse
Immunogen	Recombinant rat SAP97 (N-terminal fragment).
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	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 ascites.
Species Reactivity	Dog, Human, Mouse, Rat
UniProt ID	Q62696
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)