

PSD-95 monoclonal antibody (7E3-1B8)

Postsynaptic density protein 95 kDa (PSD-95), also known as Synapse associated protein 90 kDa (SAP90), is a brain specific protein that is highly similar to the Drosophila dlg tumor suppressor protein. PSD-95 is a member of membrane associated proteins that are localized beneath the postsynaptic membrane of synapses in the CNS. PSD-95 interacts with NMDA receptor and Shaker-type K⁺ channels and contributes to their clustering and localization at the synaptic spines in hippocampal neurons and the pinceau terminal of cerebellar basket cells, respectively.

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

Citations: 5

[View Online »](#)

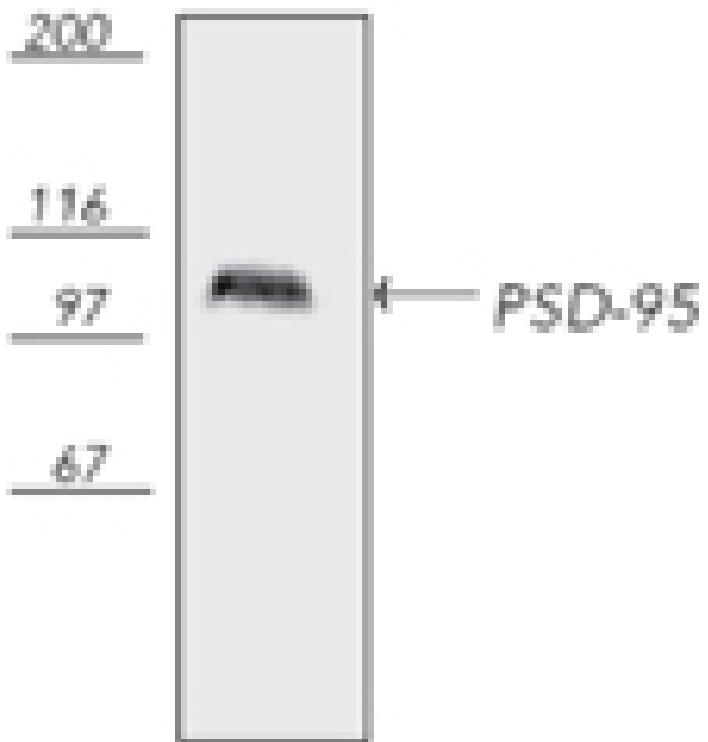
Ordering Information

[Order Online »](#)

ADI-VAM-PS001-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of bovine brain tissue extract, probed with PSD-95, mAb (7E3-1B8).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Postsynaptic density protein 95, Disks large homolog 4, Synapse-associated protein 90, SAP90
Application	WB
Application Notes	Detects a band of ~95kDa by Western blot.
Clone	7E3-1B8
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Host	Mouse
Immunogen	Recombinant rat PSD-95.
Isotype	IgG2a
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	Bovine, Mouse, Rat
UniProt ID	Q62108
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)