

SNAP-25 polyclonal antibody

SNAP-25 (synaptosomal associated protein of 25kDa) is a soluble protein of 25kDa which plays a key role in vesicle membrane fusion events with the plasma membrane. This protein was originally discovered to be enriched in subsets of neurons in the brain and displays a presynaptic pattern of expression. The membrane proteins SNAP-25, synaptobrevin, and syntaxin form the core of a ubiquitous membrane fusion machine that interacts with the soluble proteins NSF and α -SNAP. Functional interactions have been demonstrated to exist between SNAP-25, syntaxin, and the synaptic protein interaction site on voltage-sensitive L- and N-type Ca^{2+} channels. SNAP-25 has been shown to be required for synaptic vesicle fusion with the plasma membrane during release of NGF. Regulated exocytosis of cortical granule secretion in fertilized eggs and membrane fusion events in neurons and endocrine cells is mediated by SNAP-25 in a Ca^{2+} -dependent mechanism. SNAP-25 protein levels have been shown to be elevated in prolactinoma and GH/PRL tumor cells while reduced SNAP-25 protein expression has been observed between schizophrenic and normal hippocampal cells. Altered patterns of SNAP-25 expression have also been observed in the inferior temporal cortex and prefrontal association cortex between normal brains and brains from individuals affected with schizophrenia. SNAP-25 accumulation due to cytoskeletal dysfunction is also observed in the swollen axons of the white matter of individuals with severe Alzheimer's dementia.

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

Ordering Information

[Order Online »](#)

ALX-210-279-R100	100 μ l
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

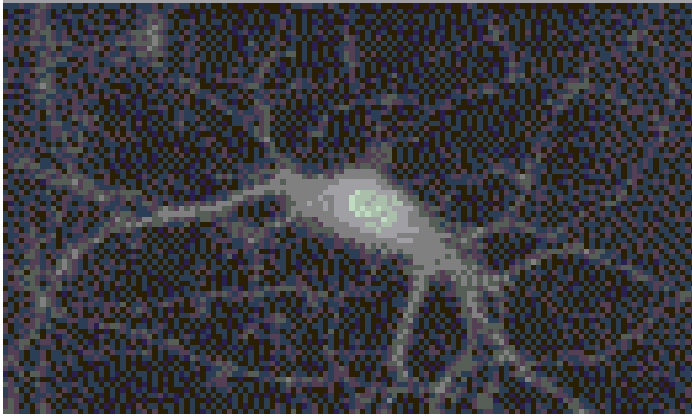


Figure: Immunocytochemistry of SNAP-25 in cultured rat hippocampal cells using PAb to SNAP-25 (Prod. No. [ALX-210-279](#)).^{""}, SNAP-25 polyclonal antibody

Immunocytochemistry, SNAP-25 polyclonal antibody

Immunocytochemistry, yes

ALX-210-302, ALX-210-302_WB.gif, http://static.enzolifesciences.com/fileadmin/files/image/ALX-210-302_WB.gif, **Figure:** Western blot using rabbit polyclonal anti PARP-1.

Lane 1: recombinant human PARP-1 (ALX-201-053

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 Synaptosomal-associated protein 25

Application ICC, WB

Application Notes Detects a band of ~25kDa by Western blot.

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

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 195-206 (A195NQRATKMLGSG206) of rat SNAP-25 (synaptosomal-associated protein 25).

Purity Detail Epitope-affinity purified IgG.

Recommendation Dilutions/Conditions Immunocytochemistry (1:100)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Bovine, Mouse, Rat, Zebrafish

UniProt ID P60881

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-](mailto:info-usa@enzolifesciences.com)
[usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-](mailto:info-eu@enzolifesciences.com)
[eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-](mailto:info-fr@enzolifesciences.com)
[fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-](mailto:info-fr@enzolifesciences.com)
[fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-](mailto:info-de@enzolifesciences.com)
[de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-](mailto:info-uk@enzolifesciences.com)
[uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

Last modified: May 29, 2024