

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](#).

Manuals, SDS & CofA

[View Online »](#)

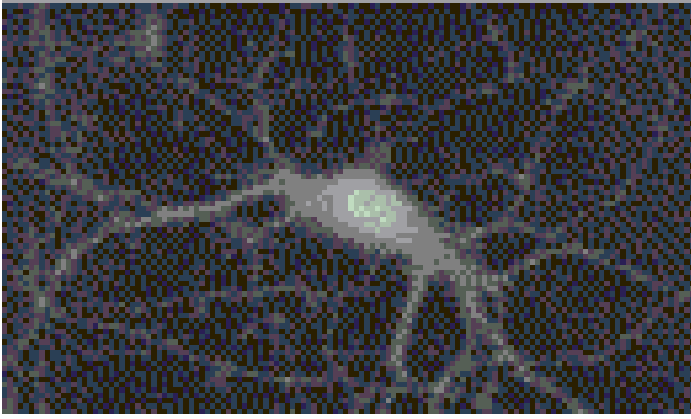


Figure: Immunocytochemistry of SNAP-25 in cultured rat hippocampal cells using PAb to SNAP-25 (Prod. No. [ALX-210-279](#)).¹, SNAP-25 polyclonal antibody
Immunocytochemistry, SNAP-25 polyclonal antibody
Immunocytochemistry, yes
ALX-210-302, ALX-210-302_WB.gif, http://static.enzolifesciences.com/fileadmin/files/image/ALX-210-302_WB.gif, **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