

# PRODUCT INFORMATION



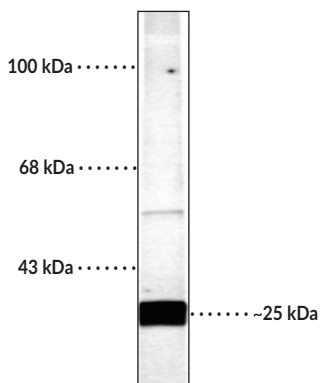
## SNAP-25 Polyclonal Antibody (Clone D005)

Item No. 29293

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µl of affinity-purified rabbit polyclonal antibody.
<b>Synonyms:</b>	SUP, Super Protein, Synaptosomal-associated 25 kDa Protein
<b>Immunogen:</b>	Peptide corresponding to amino acid residues from the C-terminal region of rat SNAP-25
<b>Species Reactivity:</b>	(+) Rat; other species not tested
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 µg/ml BSA, and 50% glycerol
<b>Host:</b>	Rabbit
<b>Applications:</b>	Western blot (WB); the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



WB of rat cortical lysate showing specific labeling of the ~25 kDa SNAP-25 protein.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/30/2023

### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

Synaptosomal-associated protein 25 (SNAP-25) is a member of the SNARE complex, which also includes syntaxin and VAMP, that is responsible for fusing synaptic vesicles with the presynaptic plasma membrane to facilitate neurotransmitter release.<sup>1</sup> Two SNAP-25 isoforms, SNAP-25a and SNAP-25b, are generated through alternative splicing, with SNAP-25b expressed only during the postnatal period and as the predominant isoform in the brain.<sup>2</sup> SNAP-25 contains two  $\alpha$  helices, as well as one large and several smaller intrinsically disordered domains.<sup>1</sup> It also contains a cluster of closely spaced cysteine residues that are subject to palmitoylation, a modification that regulates the intracellular distribution of SNAP-25, as well as residues subject to phosphorylation, modifications that modulate the rate of vesicle recruitment and its interactions with syntaxin.<sup>1,3,4</sup> SNAP-25 is located primarily on the intracellular side of the presynaptic plasma membrane in neurons and interacts with a variety of proteins to orchestrate vesicle fusion in a calcium-triggered manner and to mediate spine development.<sup>1</sup> It is also found in the pancreas, enteroendocrine cells, and the chromaffin cells of the adrenal medulla where it is involved in hormone secretion.<sup>2</sup> Decreased expression of SNAP-25 in a heterozygous mouse model leads to moderate hyperactivity, impaired learning and memory, and increased susceptibility to seizures induced by kainate. SNPs in SNAP-25 are associated with attention deficient/hyperactivity disorder, as well as bipolar disorder, schizophrenia, and autism spectrum disorder.<sup>5</sup> Protein levels of SNAP-25 are decreased in postmortem brain from patients with Down syndrome and Alzheimer's disease and in postmortem hippocampus from patients with schizophrenia.<sup>6,7</sup> Cayman's SNAP-25 Polyclonal Antibody (Clone D005) can be used for Western blot (WB) applications. The antibody recognizes SNAP-25 at 25 kDa from rat samples.

## References

1. Karmakar, S., Sharma, L.G., Roy, A., *et al.* Neuronal SNARE complex: A protein folding system with intricate protein-protein interactions, and its common neuropathological hallmark, SNAP25. *Neurochem. Int.* **122**, 196-207 (2019).
2. Kádiová, A., Radecke, J., and Sørensen, J.B. The SNAP-25 protein family. *Neuroscience* **420**, 50-71 (2019).
3. Nagy, G., Matti, U., Nehring, R.B., *et al.* Protein kinase C-dependent phosphorylation of synaptosome-associated protein of 25 kDa at Ser<sup>187</sup> potentiates vesicle recruitment. *J. Neurosci.* **22(21)**, 9278-9286 (2002).
4. Greaves, J. and Chamberlain, L.H. Differential palmitoylation regulates intracellular patterning of SNAP25. *J. Cell. Sci.* **124(Pt 8)**, 1351-1360 (2010).
5. Safari, M.R., Omrani, M.D., Noroozi, R., *et al.* Synaptosome-associated protein 25 (SNAP25) gene association analysis revealed risk variants for ASD, in Iranian population. *J. Mol. Neurosci.* **61(3)**, 305-311 (2017).
6. Greber, S., Lubec, G., Cairns, N., *et al.* Decreased levels of synaptosomal associated protein 25 in the brain of patients with Down Syndrome and Alzheimer's disease. *Electrophoresis* **20(4-5)**, 928-934 (1999).
7. Fatemi, S.H., Earle, J.A., Stary, J.M., *et al.* Altered levels of the synaptosomal associated protein SNAP-25 in hippocampus of subjects with mood disorders and schizophrenia. *Neuroreport* **12(15)**, 3257-3262 (2001).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM