

Synaptotagmin monoclonal antibody (ASV48)

Synaptotagmins (Syt) are membrane-trafficking proteins characterized by an N-terminal transmembrane region, a variable linker, and two C-terminal C2-domains, C2A and C2B. At least twelve Synaptotagmins are expressed in vertebrates and can be found in both vesicular and plasma membranes.

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

Citations: 5

[View Online »](#)

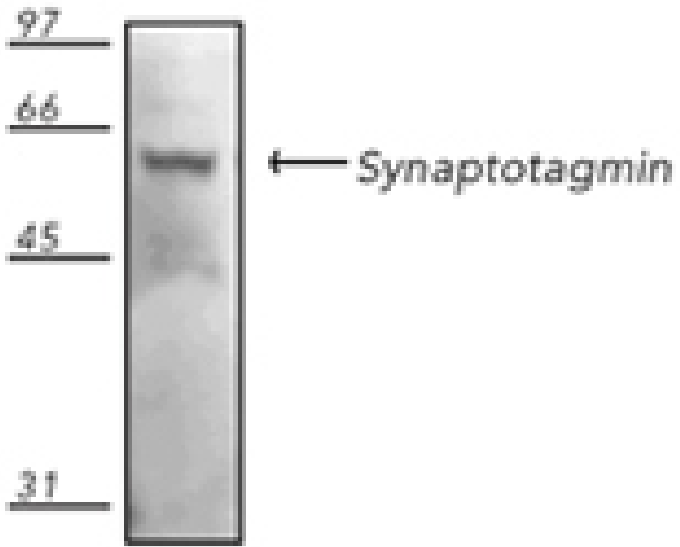
Ordering Information

[Order Online »](#)

ADI-SYA-148-D	50µg
---------------	------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of rat brain tissue extract, probed with Synaptotagmin, mAb (ASV48).

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	SYT
Application	ICC, IHC, IP, WB
Application Notes	Detects a band of ~65kDa by Western blot.
Clone	ASV48
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
GenBank ID	X52772
Host	Mouse
Immunogen	Rat brain synaptic junction protein complexes.
Isotype	IgG2b
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:500)Western Blot (0.1µg/ml, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from mouse ascites.
Species Reactivity	Monkey, Mouse, Rat, Shark
UniProt ID	P21707

Last modified: May 29, 2024



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)