

# Synaptotagmin monoclonal antibody (ASV30)

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: 3

[View Online »](#)

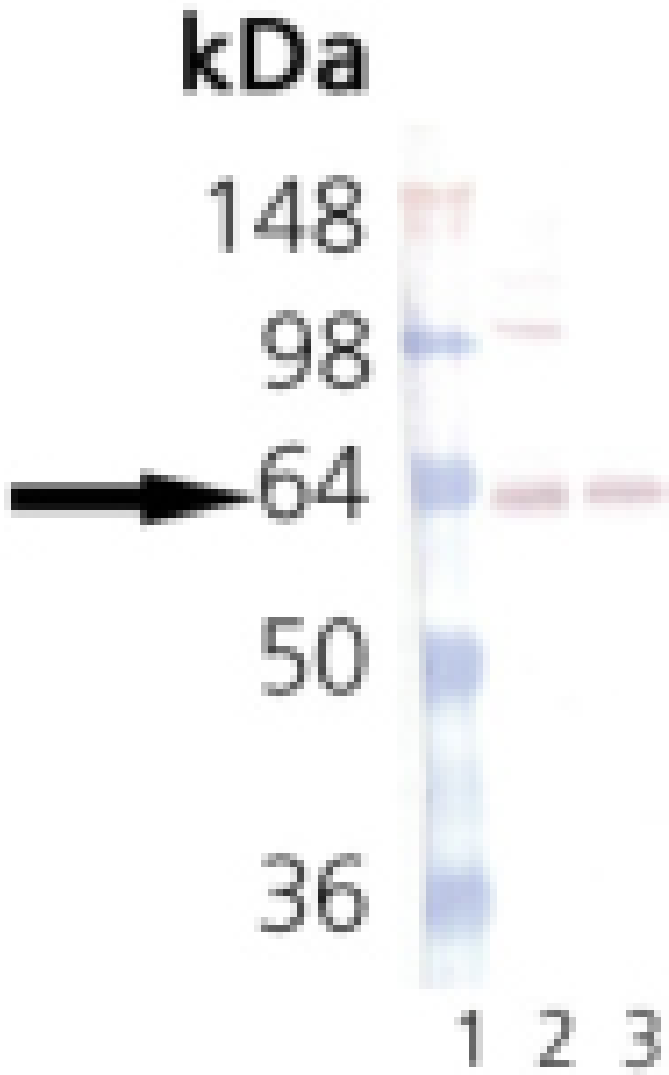
Ordering Information

[Order Online »](#)

ADI-SYA-130-F	200µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Synaptotagmin: Lane 1: MW marker, Lane 2: Mouse Brain Tissue Lysate (Prod No. ADI-LYT-MB100), Lane 3: Rat Brain Tissue Lysate (Prod No. ADI-LYT-RB100).

## 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** IHC, WB

**Application Notes** Detects a band of ~65kDa by Western blot.

**Clone** ASV30

**Formulation** Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

**GenBank ID** X52772

**Host** Mouse

**Immunogen** Rat brain synaptic junction protein complexes .

**Isotype** IgG2a

**Purity Detail** Protein G affinity purified.

**Recommendation Dilutions/Conditions** Western Blot (1:1,000, 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** Human, Mouse, Rat

**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)