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 Synaptotagamins 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 »

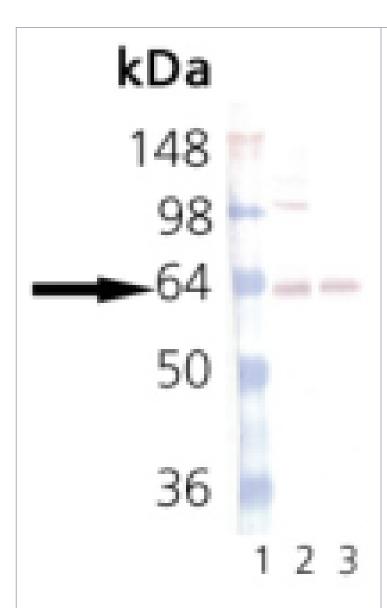
Ordering Information

Order Online »

ADI-SYA-130-D	50µg
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 NotesDetects 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 .

lgG2a

Purity Detail Protein G affinity purified.

Recommendation Dilutions/ConditionsWestern 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



info-

eu@enzolifesciences.com