

# Syntaxin 6

## monoclonal antibody

### (3D10)

Syntaxins are type II integral membrane proteins localized to membranes that regulate vesicle transport, protein maturation, and exocytosis. Syntaxin 1A resides in the nerve terminals of sensory neurons and nerve fibers reaching small blood vessels. Syntaxin 2 localizes to the apical plasma membrane and regulates epithelial-mesenchymal interactions. Syntaxin 6 is primarily localized to the trans-Golgi network, and partially co-localizes with AP-1 on clathrin-coated membranes. Syntaxin 13 is ubiquitously expressed and localizes to the tubular early and recycling endosomes, as well as endosomal vacuoles.

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

Citations: 4

[View Online »](#)

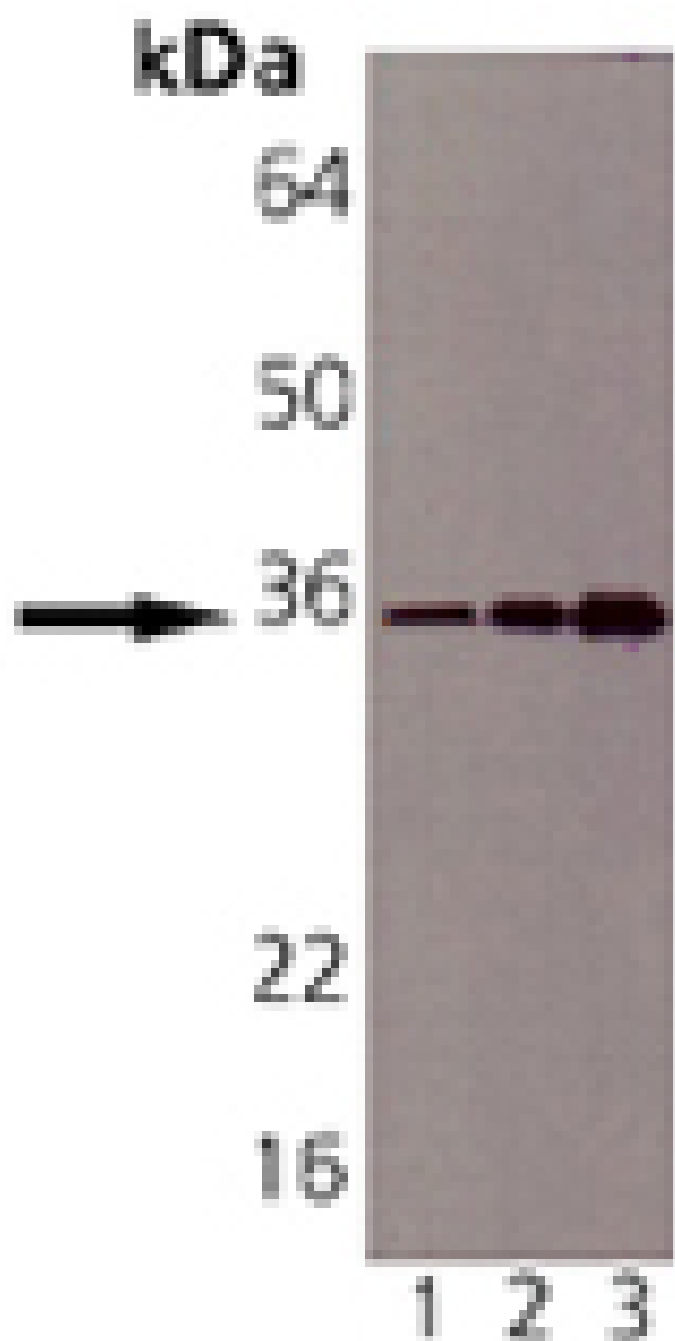
### Ordering Information

[Order Online »](#)

ADI-VAM-SV025-D	50µg
ADI-VAM-SV025-F	200µg

### Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Syntaxin 6: Lane 1: CHO-K1,  
Lane 2: Mouse Brain, Lane 3: Rat Brain.

## 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** STX6

**Application** IP, WB

**Application Notes** Detects a band of ~35kDa by Western blot.

**Clone** 3D10

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

**Host** Mouse

**Immunogen** Recombinant rat Syntaxin 6.

**Isotype** IgG1

**Purity Detail** Protein G affinity purified.

**Recommendation** Immunoprecipitation (12µg/ml) Western Blot (1:1,000, ECL) Suggested  
**Dilutions/Conditions** 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** Hamster, Human, Mouse, Rat

**UniProt ID** Q63635



ENZO LIFE SCIENCES,  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 1 42 44 55 55  
[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)

This antibody is covered by our **Worry Free Guarantee**