# 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 »

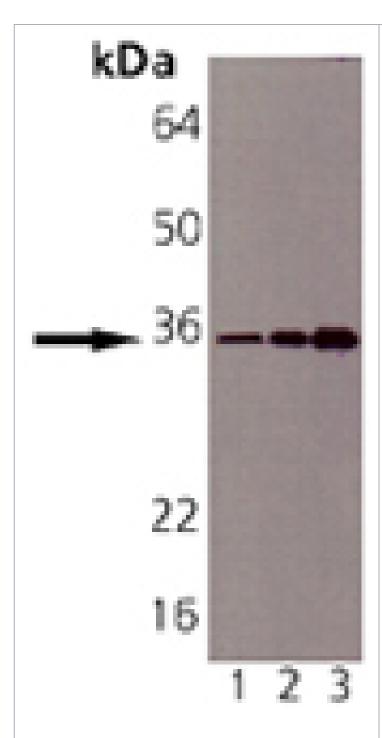
## **Ordering Information**

**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.

lgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Immunoprecipitation (12µg/ml)Western Blot (1:1,000,

ECL)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 Hamster, Human, Mouse, Rat

UniProt ID Q63635

Last modified: May 29, 2024



info-

eu@enzolifesciences.com