

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.

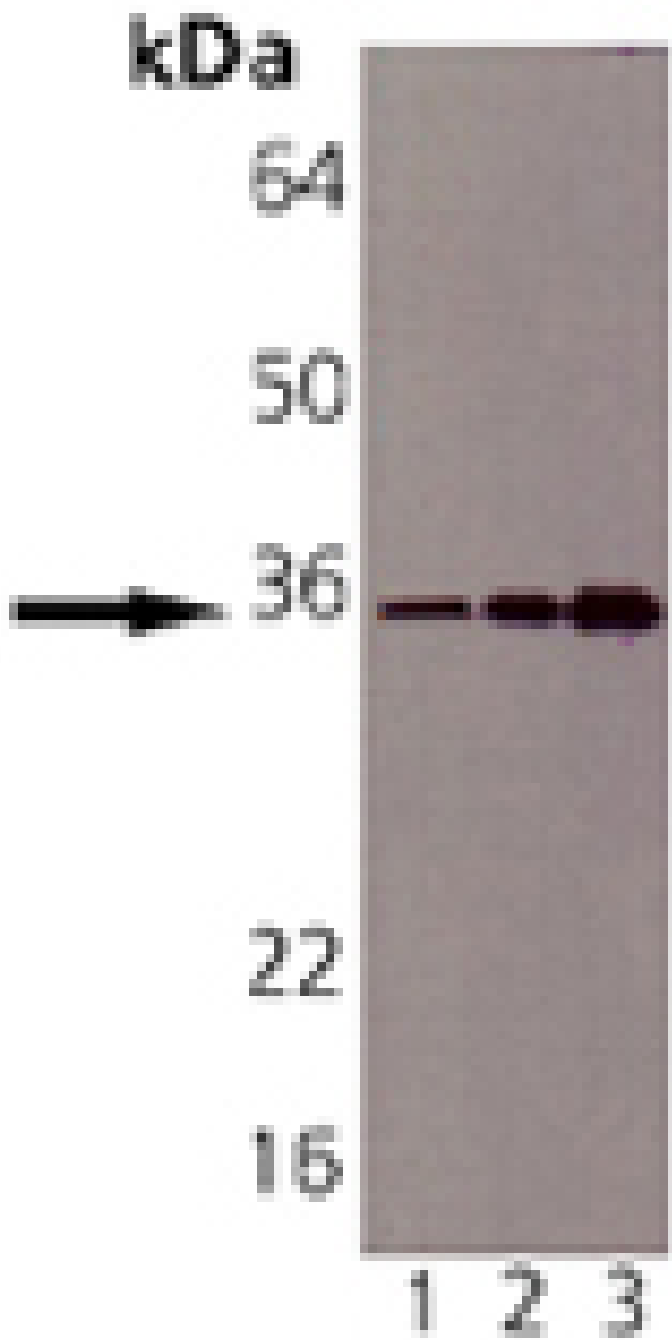
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 4

[View Online »](#)

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 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



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)