

Syntaxin 13

monoclonal antibody

(15G2)

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: 1

[View Online »](#)

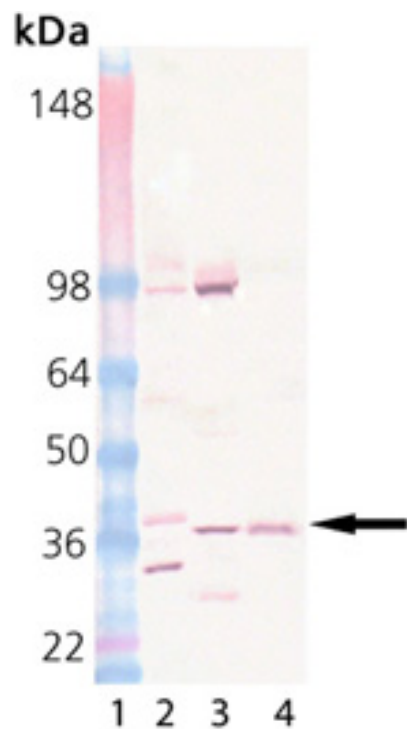
Ordering Information

[Order Online »](#)

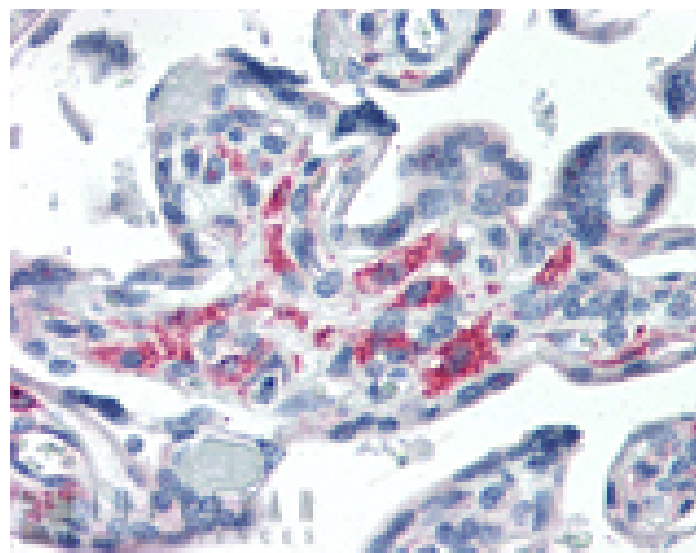
ADI-VAM-SV026-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Syntaxin 13 (15G2), mAb (Prod. No. ADI-VAM-SV026): Lane 1: MW Marker, Lane 2: Rat Brain Lysate (Prod. No. ADI-LYT-RB100), Lane 3: Mouse Brain Lysate (Prod. No. ADI-LYT-MB100), Lane 4: CHO-K1 Cell Lysate (Heat Shocked)



Immunohistochemistry analysis of human placenta tissue stained with Syntaxin 13, mAb (15G2) at 20 µg/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	STX13, Syntaxin 12
Application	IHC (PS), WB
Application Notes	Detects a band of ~38kDa by Western blot.
Clone	15G2
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
GenBank ID	AF044581
Host	Mouse
Immunogen	Recombinant cytoplasmic domain of rat syntaxin 13.
Isotype	IgG
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 ascites.
Species Reactivity	Dog, Hamster, Human, Mouse, Rat
UniProt ID	G3V7P1

Worry-free Guarantee

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com