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

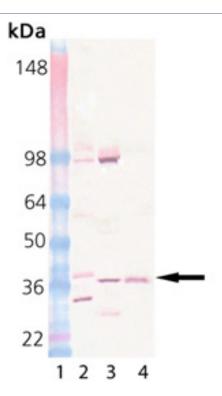
Order Online »

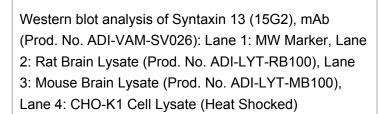
ADI-VAM-SV026-E

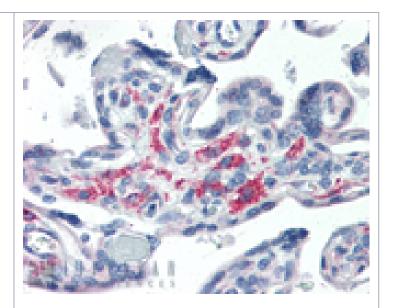
100µg

Manuals, SDS & CofA

View Online »







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

Handling & Storage

-20°C **Long Term Storage**

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

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

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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com