

# 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 »](#)

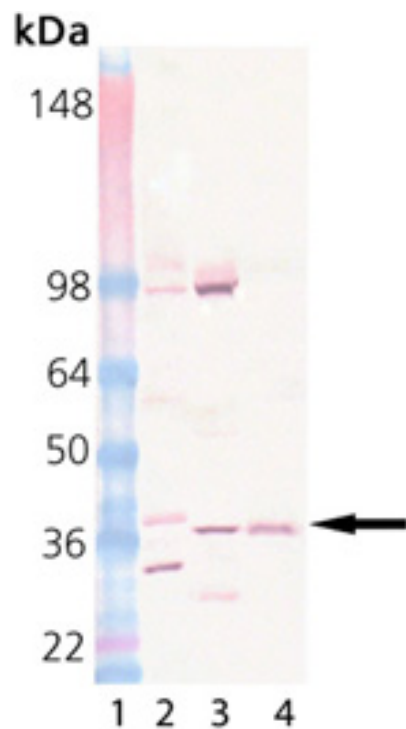
### Ordering Information

[Order Online »](#)

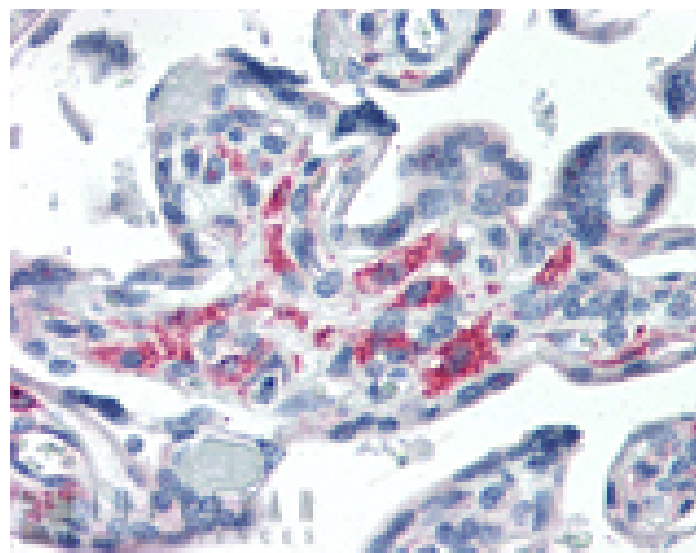
|                 |       |
|-----------------|-------|
| ADI-VAM-SV026-E | 100µg |
|-----------------|-------|

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](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)