

# rSec8 monoclonal antibody (14G1)

rSec6 and rSec8 are rat homologs of the yeast Sec6 and Sec8 proteins that are essential for protein transport between the Golgi apparatus and the plasma membrane. In yeast, Sec8 is part of a 19.5S complex that includes Sec6 and Sec15, that also associate with the ras-like GTP binding protein Sec4. An analogous 17S complex in rat brain was found to contain rSec6 and rSec8, along with eight proteins with a combined molecular weight of 743 kDa. The membrane-associated rSec6/8 complex co-immunoprecipitates with syntaxin, providing evidence that this complex is involved in synaptic vesicle docking and fusion at the presynaptic terminal.

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

Citations: 13

[View Online »](#)

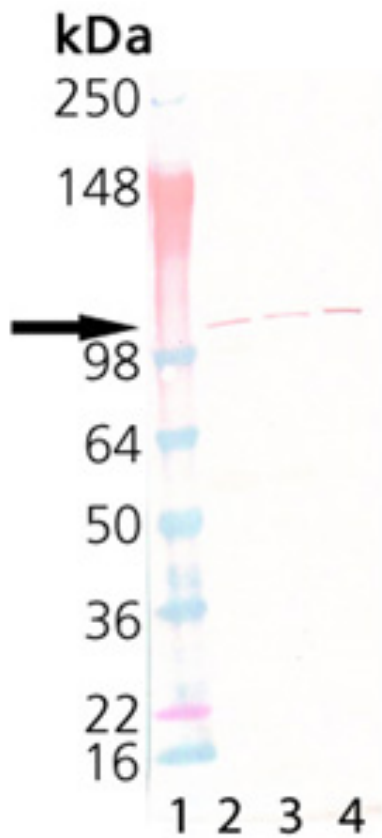
## Ordering Information

[Order Online »](#)

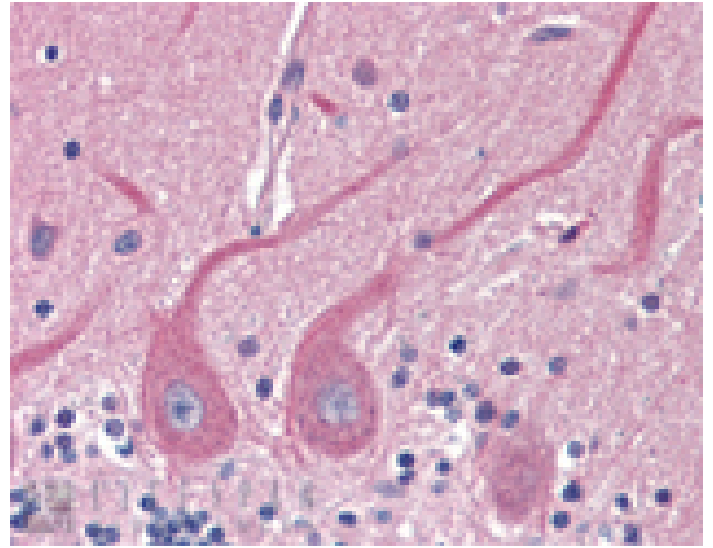
ADI-VAM-SV016-F	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of rSec8: Lane 1: MWM, Lane 2: Brain (rat), (tissue extract) (Prod No. ADI-LYT-RB100), Lane 3: Brain (mouse), (tissue extract) (Prod No. ADI-LYT-MB100), Lane 4: MDBK Cell Lysate.



Immunohistochemistry analysis of human brain, cerebellum tissue stained with rSec8, mAb (14G1) at 5µg/ml.

## 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** Exocyst complex component 4

**Application** ICC, IHC (PS), IP, WB

**Application Notes** Detects a band of ~110kDa by Western blot.

**Clone** 14G1

**Formulation** Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

**GenBank ID** U32498

**Host** Mouse

**Immunogen** Recombinant rat rSec8.

**Isotype** IgG1

**Purity Detail** Protein G affinity purified.

**Recommendation Dilutions/Conditions** Immunoprecipitation (12.5µg/ml)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** Bovine, Chicken, Dog, Fish, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep

UniProt ID

Q62824

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)