

rSec6 monoclonal antibody (9H5)

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: 5

[View Online »](#)

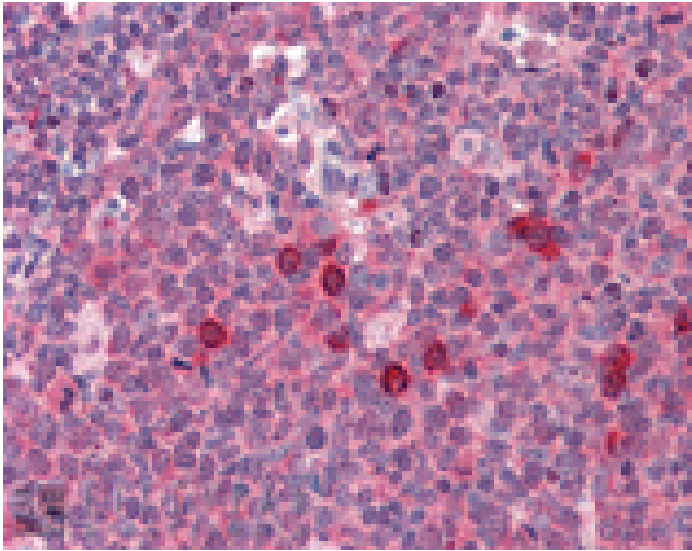
Ordering Information

[Order Online »](#)

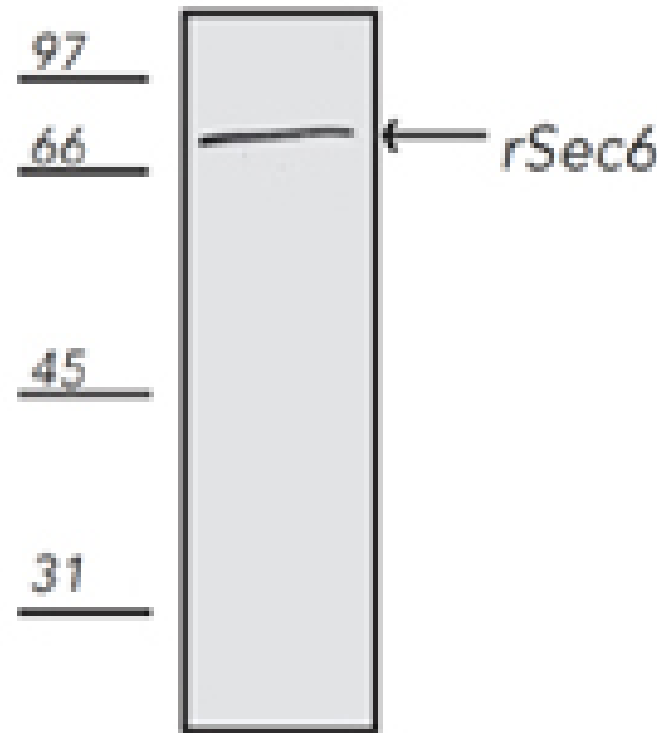
ADI-VAM-SV021-D	50µg
ADI-VAM-SV021-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human tonsil tissue stained with rSec6, mAb (9H5) at 10µg/ml.



Western blot analysis of pig ESK-4 cell lysate, probed with Prod. No. ADI-VAM-SV021.

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 3
Application	ICC, IHC (PS), WB
Application Notes	Detects a band of ~87kDa by Western blot.
Clone	9H5
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Host	Mouse
Immunogen	Recombinant rat Sec6.
Isotype	IgG2a
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:250, ECL)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, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep
Technical Info / Product Notes	Recommended by the Human Protein Atlas Organization for IHC (Ensembl No. ENSG00000180104).

UniProt ID

Q62825

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)