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 »

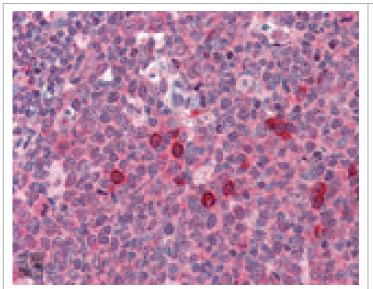
Ordering Information

Order Online »

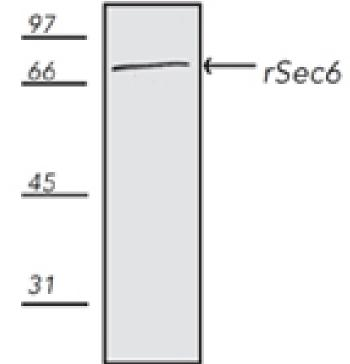
ADI-VAM-SV021-D	50µg
ADI-VAM-SV021-F	200µg

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.

lgG2a Isotype

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).

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com