

Membrin monoclonal antibody (4HAD6)

Membrin, also known as GS27 (Golgi SNARE of 27K), is an integral membrane protein present on the surface of the Golgi apparatus. GS27 is a member of a protein complex that consists of cis-Golgi vesicle receptor syntaxin 5 and GS28, v-SNARE rbet1 and Rsec22, and Rsly1. These complexes are postulated to direct the fusion of ER-derived vesicles with vesicular tubular cluster (VTC), and the fusion of VTCs to form cis-Golgi compartment.

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

Citations: 3

[View Online »](#)

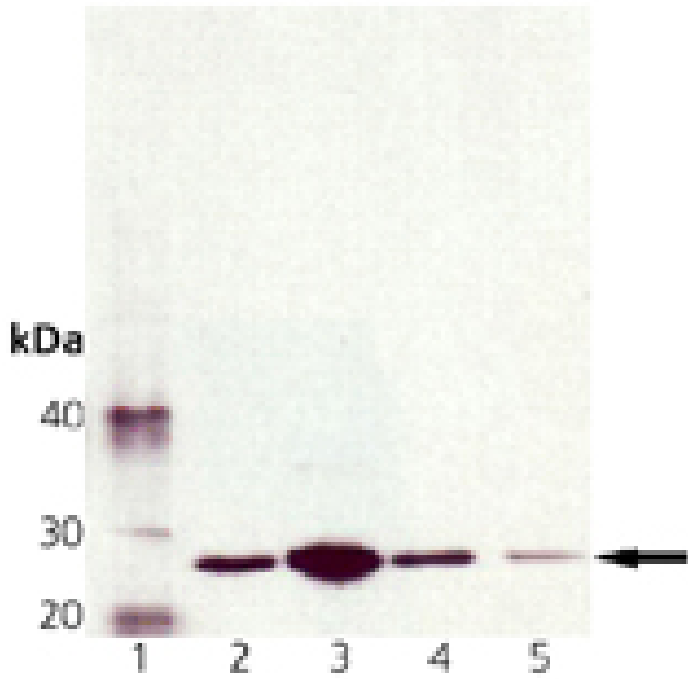
Ordering Information

[Order Online »](#)

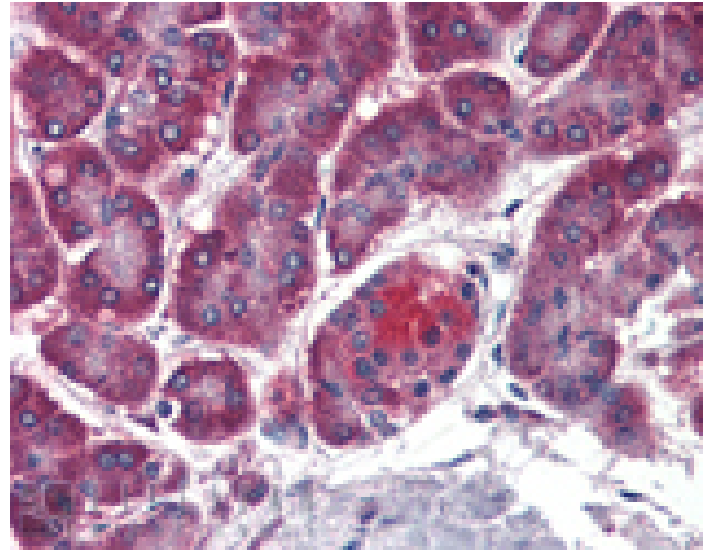
ADI-VAM-PT046-D	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of VAM-PT046: Lane 1: MW marker, Lane 2: HeLa, Lane 3: PC-12, Lane 4: RK13, Lane 5: Vero.



Immunohistochemistry analysis of human pancreas tissue stained with Membrin, mAb (4HAD6) at 20 μ g/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	GOSR2, golgi SNAP receptor complex member
Application	ICC, IHC (PS), WB
Application Notes	Detects a band of ~27kDa by Western blot.
Clone	4HAD6
Formulation	Liquid. In PBS, pH 7.2, containing 0.1MM PMSF and 50% glycerol.
Host	Mouse
Immunogen	Recombinant rat Membrin protein with a His tag.
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Source	Purified from ascites.
Species Reactivity	Chicken, Dog, Hamster, Human, Monkey, Porcine, Rabbit, Rat
UniProt ID	O35165
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee