

# 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 »](#)

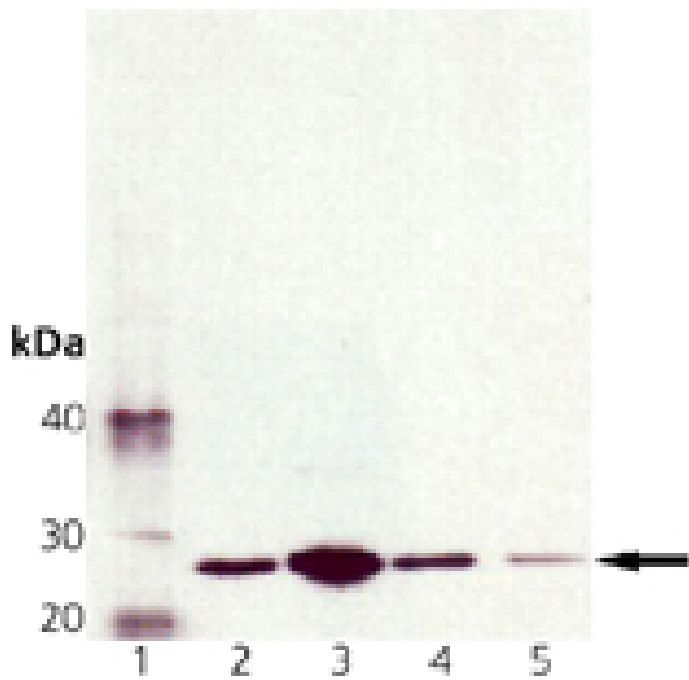
## Ordering Information

[Order Online »](#)

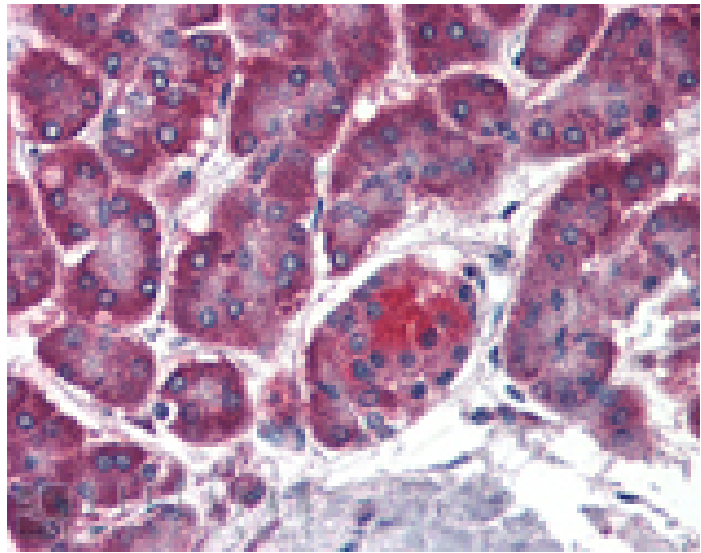
|                 |       |
|-----------------|-------|
| ADI-VAM-PT046-D | 50µg  |
| ADI-VAM-PT046-F | 200µg |

## 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 µg/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 <a href="#">Worry-Free Guarantee</a> |