

# Membrin polyclonal antibody

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: 4

[View Online »](#)

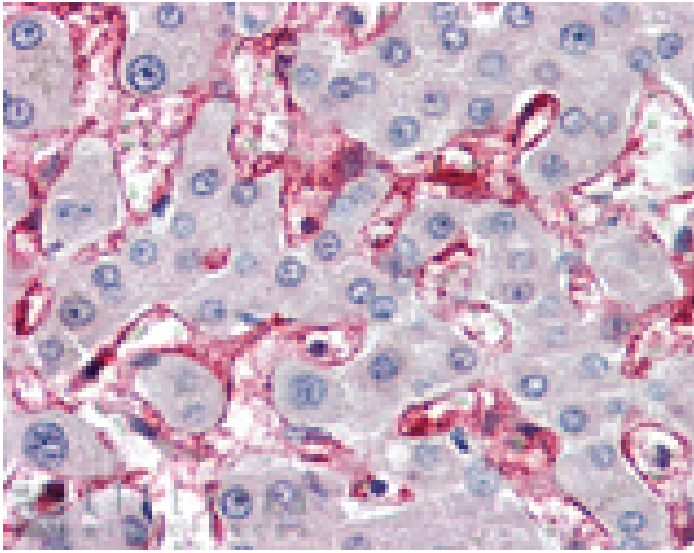
## Ordering Information

[Order Online »](#)

ADI-VAP-PT049-C	25µg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human liver tissue stained with Membrin, pAb at 10µg/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	GOSR2, golgi SNAP receptor complex member
Application	IHC (PS), WB
Application Notes	Detects a band of ~27kDa by Western blot. Cited applications include Immunofluorescence.
Formulation	Liquid. In borate buffered saline, pH 8.4, containing 25mM sodium borate, 100mM boric acid, 75mM sodium chloride, 5mM EDTA (disodium salt).
GenBank ID	U91539
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of rat membrin.
Purity Detail	Peptide immunoaffinity purified.
Recommendation Dilutions/Conditions	Western Blot (2µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Bovine, Chicken, Dog, Drosophila, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep
Technical Info / Product Notes	Recommended by the Human Protein Atlas Organization for IHC (Ensembl No. ENSG00000108433).

UniProt ID

O35165

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)