

# GS28 monoclonal antibody (HFD9)

Golgi SNARE of 28 kDa (GS28), also known as p28 or GOS28, is a 28 kDa integral membrane protein on the surface of the Golgi apparatus that serves as a t-SNARE in ER to Golgi transport. The amino-terminal of GS28 is exposed to the cytosol and is anchored to the cis Golgi via a 20 amino acid carboxyl-terminal hydrophobic tail. GS28 co-immunoprecipitates complexes consisting of Syntaxin 5, Rbet1, Membrin, Rsec22, and Rsly1, and is therefore implicated in ER to-Golgi or intra-Golgi vesicle transport.

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

Citations: 4

[View Online »](#)

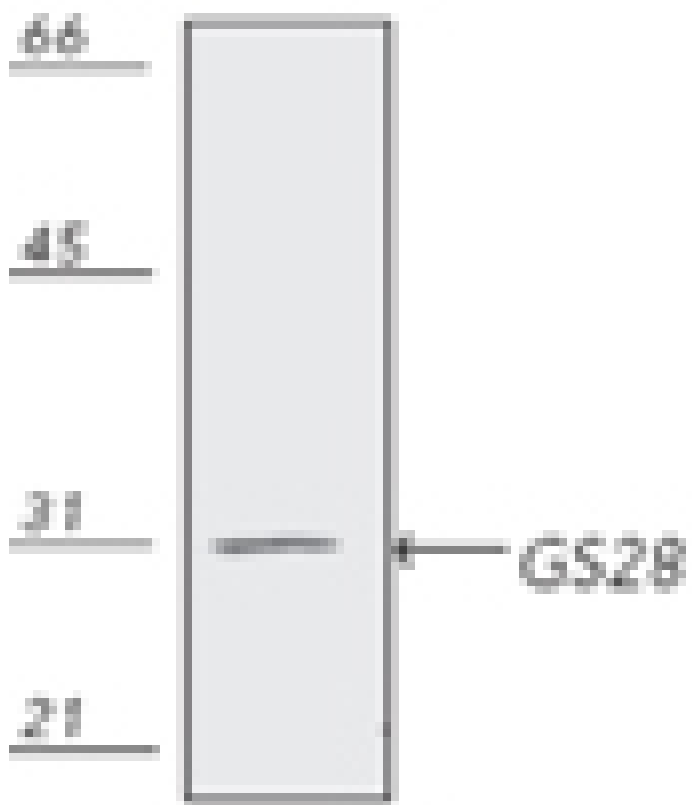
## Ordering Information

[Order Online »](#)

ADI-VAM-PT047-D	50µg
ADI-VAM-PT047-F	200µg

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of pig kidney ESK-4 cell lysate, probed with GS28 mAb (HFD9).

# 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                      GOSR1, Golgi SNAP receptor complex member 1

Application                      Electron microscopy, WB

Application Notes                      Detects a band of ~28kDa by Western blot.

Clone                      HFD9

Formulation                      Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

Host                      Mouse

Immunogen                      Rat liver Golgi-enriched membranes.

Isotype                      IgG1

Purity Detail                      Protein G affinity purified.

Recommendation Dilutions/Conditions                      Western Blot (1:1,000, 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, Mouse, Porcine, Rabbit, Rat, Sheep

UniProt ID                      Q62931

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)