

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 »](#)

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

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)