

PRODUCT INFORMATION



Ganglioside G_{M2} Asialo Polyclonal Antibody Item No. 28639

Overview and Properties

Contents:	This vial contains 50 µl of polyclonal antibody ganglioside GM2 asialo.
Immunogen:	Purified ganglioside G _{M2} asialo and complete Freund's adjuvant
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥2 years
Host:	Rabbit
Isotype:	IgG/IgM
Applications:	ELISA and TLC immunoblotting; the recommended starting dilution is 1:200 for ELISA and 1:100 for TLC immunoblotting. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

Ganglioside G_{M2} asialo (asialo-G_{M2}; Item No. 24629) is a glycosphingolipid containing three monosaccharide residues and a fatty acid of variable chain length but lacking the sialic acid residue present on ganglioside M₂. Asialo-G_{M2} levels are low-to-undetectable in normal human brain, but it accumulates in the brain of patients with Tay-Sachs and Sandhoff disease, which are neurodegenerative disorders characterized by deficiency of lysosomal β-hexosaminidase A and B, respectively.¹ It also binds to various bacteria, including *Pseudomonas* isolates derived from cystic fibrosis patients.² Ganglioside G_{M2} Asialo Polyclonal Antibody can be used for ELISA and TLC immunoblotting.

References

1. Baek, R.C., Martin, D.R., Cox, N.R., *et al.* Comparative analysis of brain lipids in mice, cats, and humans with Sandhoff disease. *Lipids* **44**(3), 197-205 (2009).
2. Krivan, H.C., Ginsburg, V., and Roberts, D.D. *Pseudomonas aeruginosa* and *Pseudomonas cepacia* isolated from cystic fibrosis patients bind specifically to gangliotetraosylceramide (asialo GM1) and gangliotriaosylceramide (asialo GM2). *Arch. Biochem. Biophys.* **260**(1), 493-496 (1988).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/02/2021

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM