

PRODUCT INFORMATION



Ganglioside G_{D1b} Polyclonal Antibody Item No. 28642

Overview and Properties

Contents:	This vial contains 100 µl polyclonal antibody to ganglioside G _{M1} .
Immunogen:	Purified ganglioside G _{D1b} 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 optimal working concentration/dilution should be determined empirically.

Description

Ganglioside G_{D1b} (Item No. 19569) is an acidic glycosphingolipid that contains two sialic acid residues linked to an inner galactose unit. It is a component of plasma membranes where it packs densely with cholesterol to form lipid microdomains that modulate both intra- and intercellular signaling events.¹ The concentration of ganglioside G_{D1b} in human brain increases with age, constituting 7.85% of total sialic acid in the brain of 0- to 10-year-old subjects and 20.29% in 11- to 30-year-old subjects.² Ganglioside G_{D1b} levels are positively correlated with pilocytic astrocytoma tumor grade, and G_{D1b} has been detected in various other gliomas, including primitive neuroectodermal tumors, glioblastomas, and anaplastic astrocytomas.³ Ganglioside G_{D1b} Polyclonal Antibody can be used for ELISA and TLC immunoblotting applications. [Matreya, LLC. Catalog No. 1964]

References

1. Kolter, T. Ganglioside biochemistry. *ISRN Biochem.* 506160 (2012).
2. Riboni, L., Sonnino, S., Acquotti, D., *et al.* Natural occurrence of ganglioside lactones. Isolation and characterization of G_{D1b} inner ester from adult human brain. *J. Biol. Chem.* **261(18)**, 8514-8519 (1986).
3. Comas, T.C., Tai, T., Kimmel, D., *et al.* Immunohistochemical staining for ganglioside GD1b as a diagnostic and prognostic marker for primary human brain tumors. *Neuro Oncol.* **1(4)**, 261-267 (1999).

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