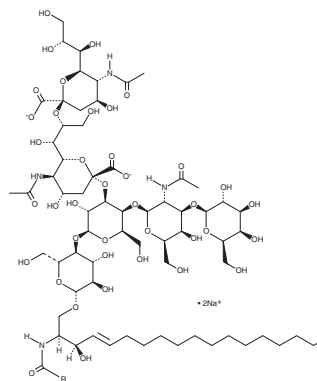


PRODUCT INFORMATION



Ganglioside G_{D1b} (porcine) (sodium salt) Item No. 19569

CAS Registry No.: 19553-76-5
Synonyms: Disialoganglioside G_{D1b},
Ganglioside C₁, Ganglioside G₂
MF: C₈₄H₁₄₆N₄O₃₉ • 2Na (for stearoyl)
FW: 1,882.1
Purity: ≥98%
Supplied as: A lyophilized powder
Storage: -20°C
Stability: ≥4 years



Special Conditions: Hygroscopic. Protect from moisture.

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ganglioside G_{D1b} (porcine) (sodium salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the ganglioside G_{D1b} (porcine) (sodium salt) in the solvent of choice, which should be purged with an inert gas. Ganglioside G_{D1b} (porcine) (sodium salt) is soluble in a 2:1 solution of chloroform:methanol. Ganglioside G_{D1b} (porcine) (sodium salt) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Ganglioside G_{D1b} 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 This product contains ganglioside G_{D1b} molecular species with primarily C18:0 fatty acyl chain lengths.

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, 11/29/2022

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