

PRODUCT INFORMATION



Ganglioside G_{D2} (bovine buttermilk) (ammonium salt)

Item No. 35774

Synonyms: Disialoganglioside G_{D2},
Ganglioside G_{D2}, Disialo

MF: C₈₃H₁₄₈N₄O₃₄ • 2NH₄ (for tricosanoyl)

FW: 1,746.1

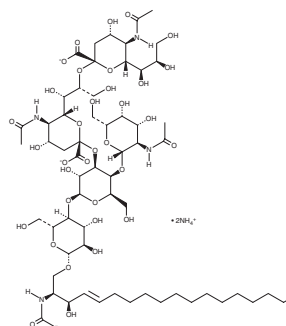
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years

Special Conditions: Forms micellar solution in water



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

Laboratory Procedures

Ganglioside G_{D2} (bovine buttermilk) (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the ganglioside G_{D2} (bovine buttermilk) (ammonium salt) in the solvent of choice. Ganglioside G_{D2} (bovine buttermilk) (ammonium salt) is soluble in a 2:1:0.1 solution of chloroform:methanol:DI water.

Description

Ganglioside G_{D2} is a glycosphingolipid that contains two sialic acid residues linked to an inner galactose unit. It is frequently expressed on tumor cells of neuroectodermal origin, including neuroblastomas.¹ This product is a semisynthetic mixture of ganglioside G_{D2} produced from buttermilk and is thus free from potential contamination with transmissible spongiform encephalopathies (TSE). It contains ganglioside G_{D2} molecular species with primarily C23:0, C22:0, and C24:0 fatty acyl chain lengths. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

Reference

1. Barker, E.L., Mueller, B.M., Handgretinger, R., *et al.* Effect of a chimeric anti-ganglioside G_{D2} antibody on cell-mediated lysis of human neuroblastoma cells. *Cancer Res.* **51(1)**, 144-149 (1991).

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, 01/16/2023

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