

PRODUCT INFORMATION



C24:1 Ganglioside G_{M2} (d18:1/24:1) (ammonium salt)

Item No. 42983

Formal Name: (2S,4S,5R,6R)-5-acetamido-2-(((2R,3S,4R,5R,6S)-3-(((2S,3R,4R,5R,6R)-3-acetamido-4,5-dihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)-6-(((2R,3S,4R,5R,6R)-4,5-dihydroxy-6-(((2S,3R,E)-3-hydroxy-2-((Z)-tetracos-15-enamido)octadec-4-en-1-yl)oxy)-2-(hydroxymethyl)tetrahydro-2H-pyran-3-yl)oxy)-5-hydroxy-2-(hydroxymethyl)tetrahydro-2H-pyran-4-yl)oxy)-4-hydroxy-6-((1R,2R)-1,2,3-trihydroxypropyl)tetrahydro-2H-pyran-2-carboxylate, monoammonium salt

Synonyms: C24:1 G_{M2},
N-Nervonoyl-(cis-15)-monosialoganglioside G_{M2},
N-Selacholeoyl-(cis-15)-monosialoganglioside G_{M2},
N-Tetracosenoyl-(cis-15)-monosialoganglioside G_{M2}

MF: C₇₃H₁₃₀N₃O₂₆ • NH₄

FW: 1,483.9

Purity: ≥98%

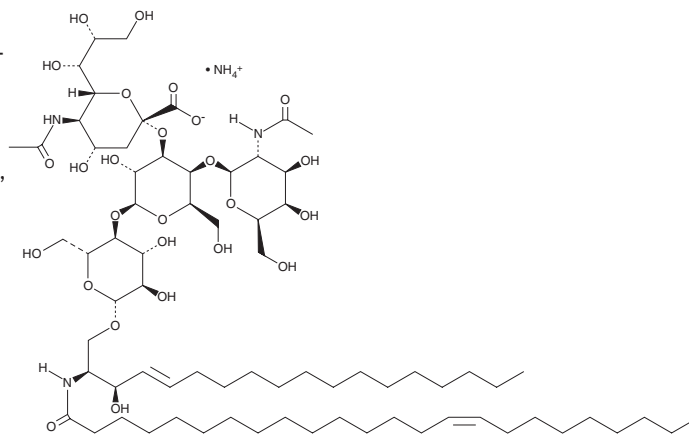
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special conditions Forms a micellar solution in water

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



Laboratory Procedures

C24:1 Ganglioside G_{M2} (d18:1/24:1) (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the C24:1 Ganglioside G_{M2} (d18:1/24:1) (ammonium salt) in the solvent of choice, which should be purged with an inert gas. C24:1 Ganglioside G_{M2} (d18:1/24:1) (ammonium salt) is soluble in a 2:1:0.1 solution of chloroform:methanol:water. We do not recommend storing the aqueous solution for more than one day.

Description

C24:1 Ganglioside G_{M2} is an endogenous monosialylated ganglioside.¹ It is a self-lipid that binds to CD1d in HEK293T cells.² C24:1 Ganglioside G_{M2} has been found as a minor ganglioside G_{M2} species in human brain from a patient with variant B Tay-Sachs disease.¹ As this product is derived from a natural source, there may be variations in the sphingoid backbone.

References

1. Mauri, L., Valsecchi, M., Casellato, R., et al. Procedure for separation of GM2 ganglioside species with different ceramide structures by a flash reversed-phase silica gel liquid chromatography. *J. Chromatogr. B. Analyt. Technol. Biomed. Life Sci.* **796**(1), 1-10 (2003).
2. Haig, N.A., Guan, Z., Li, D., et al. Identification of self-lipids presented by CD1c and CD1d proteins. *J. Biol. Chem.* **286**(43), 37692-37701 (2011).

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, 01/21/2025

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