

PRODUCT INFORMATION



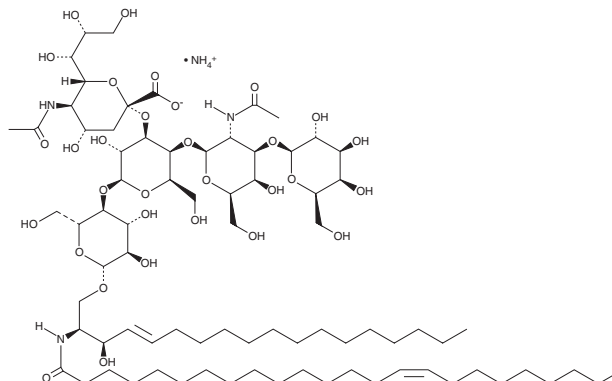
C24:1 Ganglioside G_{M1} (d18:1/24:1) (ammonium salt)

Item No. 42981

Formal Name: (2S,4S,5R,6R)-5-acetamido-2-(((2R,3S,4R,5R,6S)-3-(((2S,3R,4R,5R,6R)-3-acetamido-5-hydroxy-6-(hydroxymethyl)-4-(((2R,3R,4S,5R,6R)-3,4,5-trihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)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_{M1}, N-Nervonoyl Monosialoganglioside G_{M1}, N-Selacholeoyl Monosialoganglioside G_{M1}, N-Tetracosenoyl-(cis-15)-monosialoganglioside G_{M1}

MF: C₇₉H₁₄₀N₃O₃₁ • NH₄⁺
FW: 1,646.0
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

C24:1 Ganglioside G_{M1} (d18:1/24:1) (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the C24:1 ganglioside G_{M1} (d18:1/24:1) (ammonium salt) in the solvent of choice, which should be purged with an inert gas. C24:1 Ganglioside G_{M1} (d18:1/24:1) (ammonium salt) is soluble in a 60:40:9 solution of chloroform:methanol:2.5N ammonium hydroxide.

Description

Ganglioside G_{M1} is a monosialylated ganglioside and the prototypic ganglioside for those containing one sialic acid residue.^{1,2} C24:1 Ganglioside G_{M1} is a monosialylated ganglioside that contains nervonic acid (Item No. 13940). As this product is derived from a natural source, there may be variations in the sphingoid backbone.

References

1. Kolter, T. Ganglioside biochemistry. *ISRN Biochem.* 506160 (2012).
2. Mocchetti, I. Exogenous gangliosides, neuronal plasticity and repair, and the neurotrophins. *Cell Mol. Life Sci.* **62**(19-20), 2283-2294 (2005).

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, 01/27/2025

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