

PRODUCT INFORMATION



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

Item No. 42982

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-16,17,17,18,18,19,19,20,20,21,21,22,22,23,23,24,24,24-d₁₈)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}-d₁₈, N-Nervonoyl Monosialoganglioside G_{M1}-d₁₈, N-omega-CD₁₈-Tetracosenoyl-(cis-15)-monosialoganglioside G_{M1}, N-Selacholeoyl Monosialoganglioside G_{M1}-d₁₈, N-omega-CD₁₈-Tetracosenoyl-(cis-15)-monosialoganglioside G_{M1}

MF: C₇₉H₁₂₂D₁₈N₃O₃₁ • NH₄

FW: 1,664.1

Chemical Purity: ≥98% (C24:1 ganglioside G_{M1})

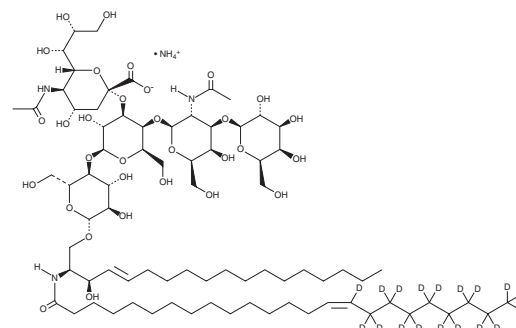
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₁₈); ≤1% d₀

Supplied as: A solid

Storage: -20°C

Stability: ≥4 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 is intended for use as an internal standard for the quantification of C24:1 ganglioside G_{M1} (Item No. 42981) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

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. Egerton, J.R., Yakstis, J.J., and Campbell, W.C. Efficacy of rafoxanide (3,5-diiodo-3'-chloro-4'-(p-chlorophenoxy) salicylanilide) against *Haemonchus contortus* in sheep. *Res. Vet. Sci.* **11**(4), 382-388 (1970).

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, 05/22/2025

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