

PRODUCT INFORMATION

C18 Ganglioside G_{D2}-d₇ (d18:1/18:0-d₇) (ammonium salt, synthetic) Item No. 43375

Formal Name: N-[(1S,2R,3E)-1-[[[O-(N-acetyl-α-neuraminosyl)-(2→8)-O-(N-acetyl-α-neuraminosyl)-(2→3)-O-[2-(acetylamino)-2-deoxy-β-D-galactopyranosyl-(1→4)]-O-β-D-galactopyranosyl-(1→4)-β-D-glucopyranosyl]oxy]methyl]-2-hydroxy-3-heptadecen-1-yl]-octadecanamide-16,16,17,17,18,18,18-d₇, diammonium salt

Synonyms: C18 Disialolactosylceramide-d₇, N-omega-CD₇-Octadecanoyl Disialoganglioside G_{D2}

MF: C₇₈H₁₂₉D₇N₄O₃₄ • 2NH₄

FW: 1,717.1

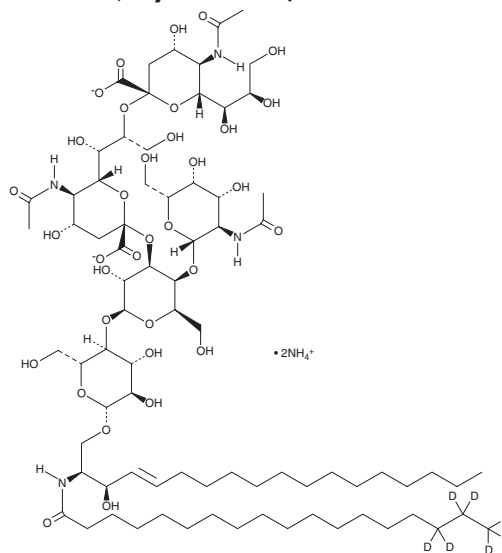
Chemical Purity: ≥98% (C18 ganglioside G_{D2})

Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀

Supplied as: A solution in methanol

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

C18 Ganglioside G_{D2}-d₇ (d18:1/18:0-d₇) (ammonium salt, synthetic) is intended for use as an internal standard for the quantification of C18 ganglioside G_{D2} 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

C18 Ganglioside G_{D2} is a ganglioside containing stearic acid (Item No. 10011298) as the acyl chain. Plasma levels of C18 ganglioside G_{D2} are increased in patients with neuroblastoma.¹ This product is fully synthetic and has no variations in the fatty acyl chain or sphingoid backbone.

Reference

1. Morini, M., Barco, S., Ardito, M., et al. Detection of plasma circulating GD2 ganglioside in patients with neuroblastoma and age-matched healthy children. Diagnostic and prognostic evaluation. *Oncologist* 30(2), oyaf008 (2025).

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, 05/06/2025

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