

# PRODUCT INFORMATION



## C6 Biotin Ganglioside G<sub>D1b</sub> (d18:1/6:0) (ammonium salt)

Item No. 9004919

**Formal Name:** (2S,4S,5R,6R)-5-acetamido-6-((1S,2R)-2-(((2S,4S,5R,6R)-5-acetamido-2-carboxylato-4-hydroxy-6-((1R,2R)-1,2,3-trihydroxypropyl)tetrahydro-2H-pyran-2-yl)oxy)-1,3-dihydroxypropyl)-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-(6-(5-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanamido)hexanamido)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-hydroxytetrahydro-2H-pyran-2-carboxylate, diammonium salt

**MF:** C<sub>82</sub>H<sub>137</sub>N<sub>7</sub>O<sub>41</sub>S • 2NH<sub>4</sub>

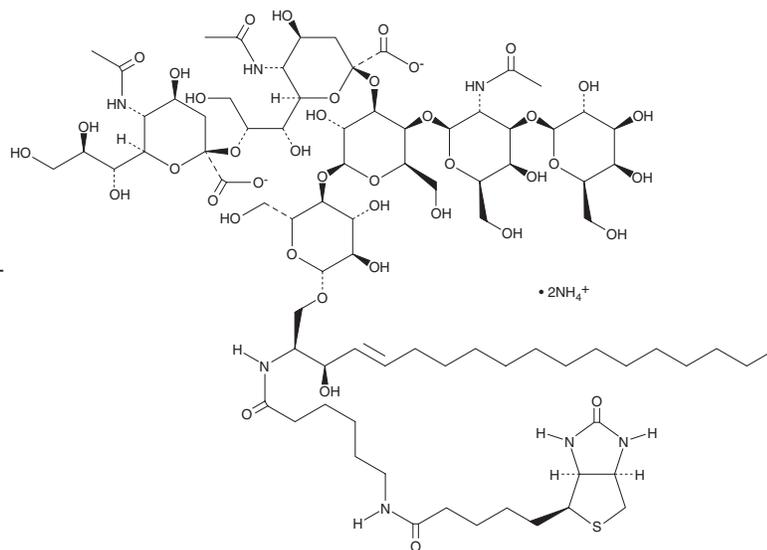
**FW:** 1,915.1

**Purity:** ≥90%

**Supplied as:** A neat 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

C6 Biotin ganglioside G<sub>D1b</sub> (d18:1/6:0) (ammonium salt) is supplied as a neat solid. A stock solution may be made by dissolving the C6 biotin ganglioside G<sub>D1b</sub> (d18:1/6:0) (ammonium salt) in the solvent of choice, which should be purged with an inert gas. C6 Biotin ganglioside G<sub>D1b</sub> (d18:1/6:0) (ammonium salt) is slightly soluble (0.1-1 mg/ml) in methanol.

### Description

C6 Biotin ganglioside G<sub>D1b</sub> (d18:1/6:0) is an affinity probe that allows ganglioside G<sub>D1b</sub> (Item Nos. 31708 | 31593 | 34778) to be detected or immobilized through interaction with the biotin ligand. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

#### 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, 01/12/2026

#### 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