

# PRODUCT INFORMATION



## C3 BODIPY Ganglioside G<sub>M1</sub> (d18:1/3:0) (ammonium salt)

Item No. 33799

**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)-6-(((2S,3R,E)-2-(3-(5,5-difluoro-7,9-dimethyl-5H-5λ<sup>4</sup>,6λ<sup>4</sup>-dipyrrolo[1,2-c:2',1'-f][1,3,2]diazaborinin-2-yl)propanamido)-3-hydroxyoctadec-4-en-1-yl)oxy)-4,5-dihydroxy-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:**

C3-BODIPY G<sub>M1</sub>,  
N-Propanoyl-BODIPY-monosialoganglioside G<sub>M1</sub>

**MF:** C<sub>69</sub>H<sub>109</sub>BF<sub>2</sub>N<sub>5</sub>O<sub>31</sub> • NH<sub>4</sub>

**FW:** 1,571.5

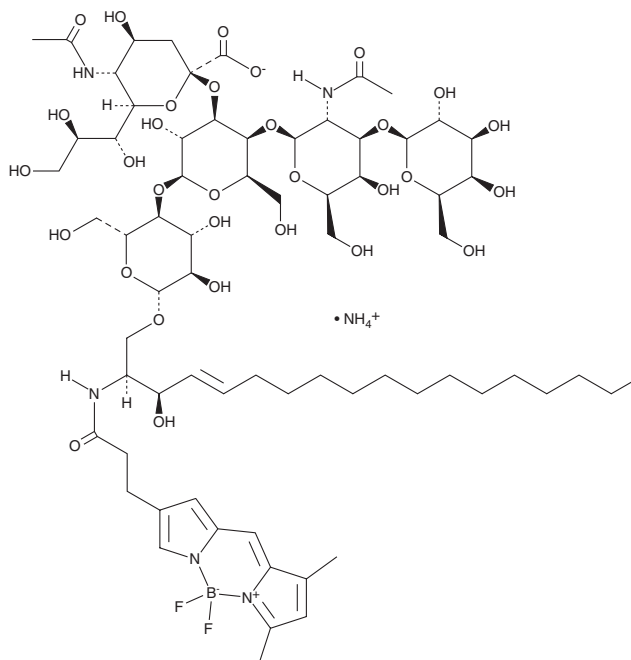
**Purity:** ≥98%

**Ex./Em. Max:** 503/512 nm

**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

C3 BODIPY Ganglioside G<sub>M1</sub> (d18:1/3:0) (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the C3 BODIPY ganglioside G<sub>M1</sub> (d18:1/3:0) (ammonium salt) in the solvent of choice, which should be purged with an inert gas. C3 BODIPY Ganglioside G<sub>M1</sub> (d18:1/3:0) (ammonium salt) is soluble in a 2:1:0.2 solution of chloroform:methanol:DI water. We do not recommend storing the aqueous solution for more than one day.

### Description

C3 BODIPY Ganglioside G<sub>M1</sub> (d18:1/3:0) is a fluorescently labeled form of ganglioside G<sub>M1</sub>. It is composed of ganglioside G<sub>M1</sub> with BODIPY 503/512 (Item No. 28800), which displays excitation/emission maxima of 503/512 nm, respectively, incorporated at the N-acyl chain position.<sup>1</sup> As this product is derived from a natural source, there may also be variations in the sphingoid backbone.

### Reference

1. Rasmussen, J.-A.M. and Hermetter, A. Chemical synthesis of fluorescent glycerol- and sphingolipids. *Prog. Lipid Res.* **47**(6), 436-460 (2008).

**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, 12/13/2022

**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