

# PRODUCT INFORMATION

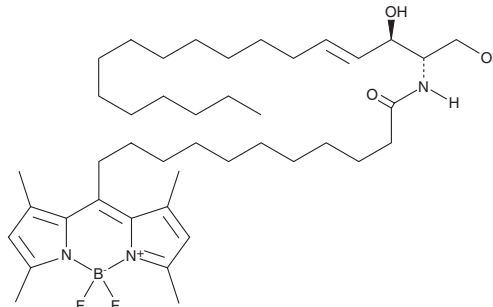


## BODIPY-C12 Ceramide (d18:1/12:0)

Item No. 25997

**CAS Registry No.:** 1246355-58-7  
**Formal Name:** (T-4)-[λ-(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)-N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-3,5-dimethyl-1H-pyrrole-2-dodecanamidato-κN<sup>1</sup>]difluoro-boron  
**Synonyms:** BODIPY-C12 Ceramide, BODIPY-Ceramide (d18:1/12:0), BODIPY-Cer(d18:1/12:0), BODIPY-N-Lauroyl-D-erythro-Sphingosine, B12Cer

**MF:** C<sub>42</sub>H<sub>70</sub>BF<sub>2</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 713.8  
**Purity:** ≥98%  
**Ex./Em. Max:** 505/540 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

BODIPY-C12 Ceramide (B12Cer) (d18:1/12:0) is supplied as a solid. A stock solution may be made by dissolving the B12Cer (d18:1/12:0) in the solvent of choice, which should be purged with an inert gas. B12Cer (d18:1/12:0) is slightly soluble in chloroform and methanol.

### Description

B12Cer is a fluorescently tagged form of C12 ceramide (Item No. 22530) that displays excitation/emission maxima of 505/540 nm, respectively.<sup>1</sup> It is formed when acid sphingomyelinase hydrolyzes BODIPY-C12 sphingomyelin *in vitro* and has been used to quantify acid sphingomyelinase activity in plasma from patients with Niemann-Pick disease.

### Reference

1. He, X., Chen, F., Dagan, A., *et al.* A fluorescence-based, high-performance liquid chromatographic assay to determine acid sphingomyelinase activity and diagnose types A and B Niemann-Pick disease. *Anal. Biochem.* **314**(1), 116-120 (2003).

#### 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, 12/19/2022

#### 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