

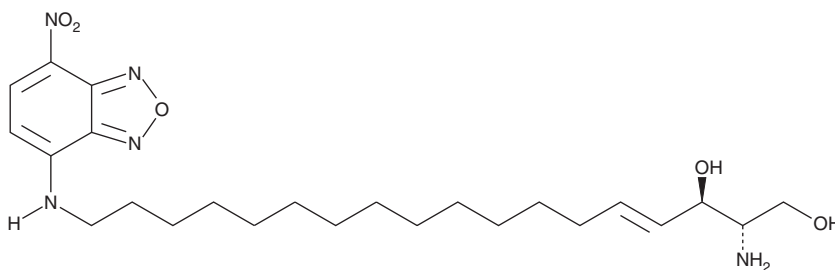
PRODUCT INFORMATION



NBD Sphingosine

Item No. 25348

CAS Registry No.: 1449370-25-5
Formal Name: (2S,3R,4E)-2-amino-18-[(7-nitro-2,1,3-benzoxadiazol-4-yl)amino]-4-octadecene-1,3-diol
MF: C₂₄H₃₉N₅O₅
FW: 477.6
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

NBD Sphingosine is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice.

Description

NBD sphingosine is a biologically active derivative of sphingosine (Item No. 10007907) that is tagged with a fluorescent nitrobenzoxadiazole (NBD) group. It has been used to study the cellular localization and metabolism of sphingosine in HeLa cells.¹ NBD sphingosine inhibits proliferation of HCT116, PC3, HEK293T, and HFE-145 cells (IC₅₀s = 25.6, 40.5, 10.4, and 12.5 μM, respectively).

References

1. Lee, Y.M., Lim, C., Lee, H.S., *et al.* Synthesis and biological evaluation of a polyene-containing sphingoid base probe as a chemical tool. *Bioconjug. Chem.* **24(8)**, 1324-1331 (2013).

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, 11/16/2018

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