

PRODUCT INFORMATION



BODIPY 480/508-Cholesteryl Oleate

Item No. 39685

Formal Name: (3S,8S,9S,10R,13R,14S,17R)-17-((R)-4-(5,5-difluoro-1,3,7,9-tetramethyl-5H-4λ⁴,5λ⁴-dipyrrolo[1,2-c:2',1'-f][1,3,2]diazaborinin-10-yl)butan-2-yl)-10,13-dimethyl-2,3,4,7,8,9,10,11,12,13,14,15,16,17-tetradecahydro-1H-cyclopenta[a]phenanthren-3-yl oleate

Synonym: BODIPY 480/508-Cholesterol Oleate

MF: C₅₄H₈₃BF₂N₂O₂

FW: 841.1

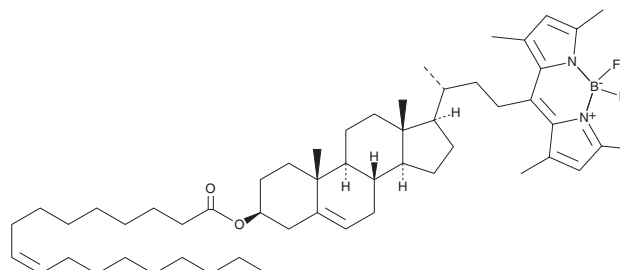
Purity: ≥95%

Ex./Em. Max: 480/508 nm

Supplied as: A solution in methyl acetate

Storage: -80°C

Stability: ≥1 year



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

Laboratory Procedures

BODIPY 480/508-cholesteryl oleate is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as chloroform purged with an inert gas can be used. The solubility of BODIPY 480/508-cholesteryl oleate in these solvents is approximately 1 mg/ml.

Description

BODIPY 480/508-cholesteryl oleate is a cell-permeable and fluorescently tagged derivative of cholesteryl oleate (Item No. 26472). BODIPY 480/508-cholesteryl oleate displays excitation/emission maxima of 480/508 nm, respectively, and has been used to monitor cholesterol trafficking in HeLa cells.¹

Reference

1. Ikonen, E. and Blom, T. Lipoprotein-mediated delivery of BODIPY-labeled sterol and sphingolipid analogs reveals lipid transport mechanisms in mammalian cells. *Chem. Phys. Lipids* **194**, 29-36 (2016).

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, 01/29/2024

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