

PRODUCT INFORMATION



BODIPY 645/660-Cholesterol

Item No. 38027

Formal Name: (3S,8S,9S,10R,13R,14S,17R)-17-((R)-4-(5,5-difluoro-3-((E)-4-methoxystyryl)-7-(1H-pyrrol-2-yl)-5H-4H,5H-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-ol

Synonyms: BODIPY 645/660-Provitamin D, Red BODIPY-Cholesterol

MF: $C_{45}H_{54}BF_2N_3O_2$

FW: 717.7

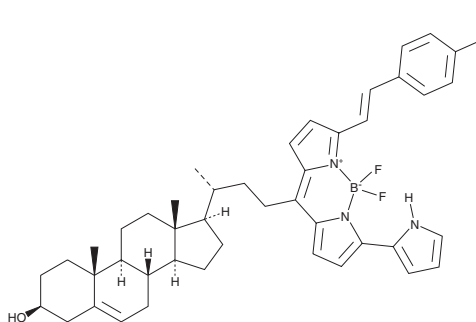
Purity: $\geq 90\%$

Ex./Em. Max: 645/660 nm

Supplied as: A 1 mg/ml solution in methyl acetate

Storage: -20°C

Stability: ≥ 1 year



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

Description

BODIPY 645/660-cholesterol is a cell-permeable and fluorescently tagged derivative of cholesterol (Item Nos. 9003100 | 39088 | 39448). BODIPY 645/660-cholesterol displays excitation/emission maxima of 645/660 nm, respectively, and has been used to monitor cholesterol trafficking in A549 cells.

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, 01/26/2026

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