

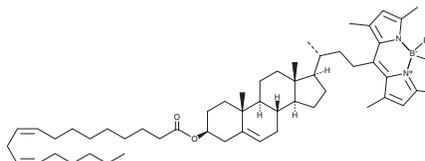
# PRODUCT INFORMATION



## BODIPY 480/508-Cholesteryl Linoleate

Item No. 39684

**CAS Registry No.:** 2260795-60-4  
**Formal Name:** (T-4)-[(3β)-24-(3,5-dimethyl-1H-pyrrol-2-yl-κN)-24-(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)chol-5-en-3-yl, 9Z,12Z-octadecadienoato]difluoro-boron  
**Synonym:** BODIPY 480/508-Cholesterol Linoleate  
**MF:** C<sub>54</sub>H<sub>81</sub>BF<sub>2</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 839.0  
**Purity:** ≥95%  
**Ex./Em. Max:** 480/508 nm  
**UV/Vis.:** λ<sub>max</sub>: 496 nm  
**Supplied as:** A 1 mg/ml solution in methyl acetate  
**Storage:** -80°C  
**Stability:** ≥2 years



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

### Description

BODIPY 480/508-cholesteryl linoleate is a cell-permeable and fluorescently tagged derivative of cholesteryl linoleate (Item No. 22597). BODIPY 480/508-cholesteryl linoleate displays excitation/emission maxima of 480/508 nm, respectively, and has been used to monitor cholesterol trafficking in isolated human monocyte-derived macrophages.<sup>1</sup>

### Reference

1. Jin, X., Dimitriadis, E.K., Liu, Y., *et al.* Macrophages shed excess cholesterol in unique extracellular structures containing cholesterol microdomains. *Arterioscler. Thromb. Vasc. Biol.* **38(7)**, 1504-1518 (2018).

#### 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, 01/27/2026

#### 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