

PRODUCT INFORMATION



BODIPY-Linoleate

Item No. 29795

Formal Name: (9Z,12Z)-17-(5,5-difluoro-1,3,7,9-tetramethyl-5H-4(1,3,2-diazaborinin-10-yl)heptadeca-9,12-dienoic acid

MF: C₃₀H₄₃BF₂N₂O₂

FW: 512.5

Purity: ≥95%

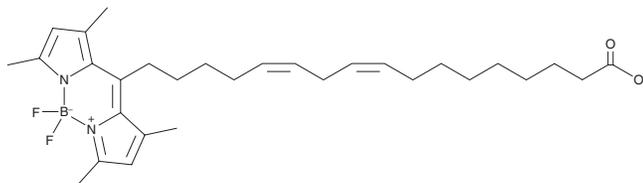
Supplied as: A 1 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥4 years

Special Conditions: Light and temperature sensitive

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



Laboratory Procedures

BODIPY-Linoleate is supplied as a solution in ethanol. A stock solution may be made by dissolving the BODIPY-linoleate in the solvent of choice, which should be purged with an inert gas. BODIPY-Linoleate is slightly soluble (0.1-1 mg/ml) in methanol and sparingly soluble (1-10 mg/ml) in chloroform.

Description

BODIPY-Linoleate is a fluorescently tagged derivative of linoleic acid (Item Nos. 90150 | 90150.1 | 21909).¹ BODIPY displays excitation/emission maxima of 470/503 nm, respectively.

Reference

- Schmitt, A., Hinkeldey, B., Wild, M., *et al.* Synthesis of the core compound of the BODIPY dye class: 4,4'-Difluoro-4-bora-(3a,4a)-diazas-indacene. *J. Fluoresc.* **19**(4), 755-758 (2009).

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, 01/05/2026

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