

PRODUCT INFORMATION



BODIPY-Palmitate

Item No. 26749

CAS Registry No.: 1246809-50-6
Formal Name: (T-4)-[o-(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)-3,5-dimethyl-1H-pyrrole-2-hexadecanoato(2-)-κN¹] difluoro-borate(1-)

Synonym: C16 BODIPY, FA16:0-BODIPY

MF: C₂₈H₄₃BF₂N₂O₂

FW: 488.5

Purity: ≥98%

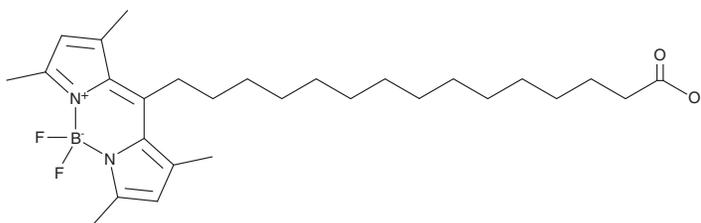
UV/Vis.: λ_{max}: 497 nm

Ex./Em. Max: 488/508 nm

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

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

Description

BODIPY-Palmitate is a fluorescently tagged form of palmitic acid (Item No. 10006627) that displays excitation/emission maxima of 488/508 nm, respectively.^{1,2} It has been used to monitor fatty acid uptake and metabolism in cultured cells.

References

1. Palanivel, R. and Sweeney, G. Regulation of fatty acid uptake and metabolism in L6 skeletal muscle cells by resistin. *FEBS Lett.* **579(22)**, 5049-5054 (2005).
2. Nealon, J.R., Blanksby, S.J., Donaldson, P.J., *et al.* Fatty acid uptake and incorporation into phospholipids in the rat lens. *Invest. Ophthalmol. Vis. Sci.* **52(2)**, 804-809 (2011).

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, 03/07/2024

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