

PRODUCT INFORMATION

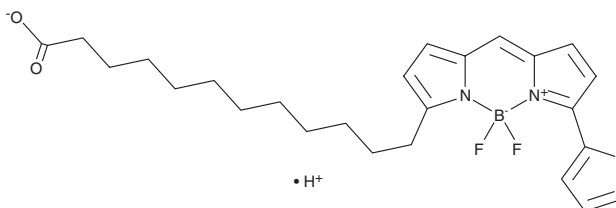


BODIPY 558/568 C12

Item No. 27014

CAS Registry No.: 158757-84-7
Formal Name: (T-4)-difluoro[5-[[5-(2-thienyl)-2H-pyrrol-2-ylidene-κN]methyl]-1H-pyrrole-2-dodecanoato(2-)-κN¹]-borate(1-), monohydrogen

Synonym: Red C12
MF: C₂₅H₃₀BF₂N₂O₂S • H
FW: 472.4
Purity: ≥95%
Ex./Em. Max: 558/568 nm
Supplied as: An oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

BODIPY 558/568 C12 is supplied as an oil. A stock solution may be made by dissolving the BODIPY 558/568 C12 in the solvent of choice, which should be purged with an inert gas. BODIPY 558/568 C12 is soluble in DMSO and dimethyl formamide.

Description

BODIPY 558/568 C12 is a fatty acid-conjugated fluorescent probe for lipid droplets.^{1,2} It displays excitation/emission maxima of 558/568 nm, respectively, and has been used to monitor the localization and dynamics of lipid droplets in live cells.

References

1. Targett-Adams, P., Chambers, D., Gledhill, S., *et al.* Live cell analysis and targeting of the lipid droplet-binding adipocyte differentiation-related protein. *J. Biol. Chem.* **278(18)**, 15998-6007 (2003).
2. Rambold, A.S., Cohen, S., and Lippincott-Schwartz, J. Fatty acid trafficking in starved cells: Regulation by lipid droplet lipolysis, autophagy, and mitochondrial fusion dynamics. *Dev. Cell.* **32(6)**, 678-692 (2015).

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, 12/19/2022

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