

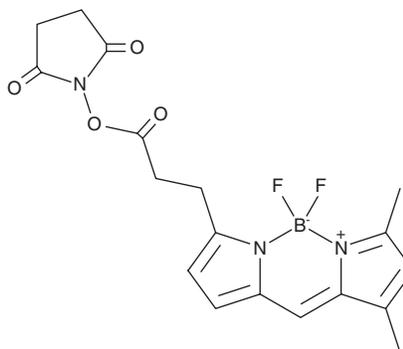
PRODUCT INFORMATION



BODIPY FL succinimide ester

Item No. 29508

CAS Registry No.: 146616-66-2
Formal Name: (T-4)-[1-[3-[5-[(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)methyl]-1H-pyrrol-2-yl-κN]-1-oxopropoxy]-2,5-pyrrolidinedionato]difluoro-boron
Synonyms: BODIPY FL NHS ester, BODIPY FL SE, EverFluor FL, SE
MF: C₁₈H₁₈BF₂N₃O₄
FW: 389.2
Purity: ≥95%
Em./Ex. Max: 502/511 nm
Supplied as: A solid
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 FL succinimide ester is supplied as a solid. A stock solution may be made by dissolving the BODIPY FL succinimide ester in the solvent of choice, which should be purged with an inert gas. BODIPY FL succinimide ester is soluble in methanol.

Description

BODIPY FL succinimide ester (BODIPY FL SE) is an amine-reactive fluorescent probe.¹ It displays excitation/emission maxima of 502/511 nm, respectively. BODIPY FL SE has been used in the synthesis of protease substrates that are non-fluorescent until unquenched following proteolytic cleavage.² It has also been used in the synthesis of a fluorescent verapamil derivative and of fluorescent inhibitors of cholesterol absorption.^{3,4}

References

1. D'Amore, C., Orso, G., Forgiarini, A., *et al.* Synthesis and biological characterization of a new norbormide derived BODIPY FL-conjugated fluorescent probe for cell imaging. *Front. Pharmacol.* **9**, 1055 (2018).
2. Jones, L.J., Upson, R.H., Haugland, R.P., *et al.* Quenched BODIPY dye-labeled casein substrates for the assay of protease activity by direct fluorescence measurement. *Anal. Biochem.* **251(2)**, 144-152 (1997).
3. Kubo, Y., Nakazawa, A., Akanuma, S.I., *et al.* Blood-to-retina transport of fluorescence-labeled verapamil at the blood-retinal barrier. *Pharm. Res.* **35(5)**, 93 (2018).
4. Burnett, D.A., Caplen, M.A., Browne, M.E., *et al.* Synthesis of fluorescent biochemical tools related to the 2-azetidinone class of cholesterol absorption inhibitors. *Bioorg. Med. Chem. Lett.* **12(3)**, 315-318 (2002).

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, 12/13/2022

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