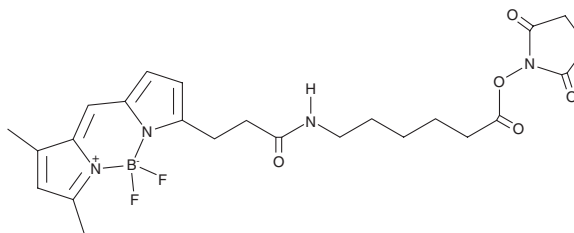


PRODUCT INFORMATION



3-BODIPY-propanoylaminocaproic Acid N-hydroxysuccinimide ester Item No. 20672

CAS Registry No.: 217190-09-5
Formal Name: (T-4)-[5-[(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)methyl]-N-[6-[(2,5-dioxo-1-pyrrolidinyl)oxy]-6-oxohexyl]-1H-pyrrole-2-propanamidato-κN¹]difluoro-boron
Synonyms: BODIPY FL-X SE, BODIPY FL-X N-succinimidyl ester
MF: C₂₄H₂₉BF₂N₄O₅
FW: 502.3
Purity: ≥98%
UV/Vis.: λ_{max}: 504 nm
Supplied as: A crystalline 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

3-BODIPY-propanoylaminocaproic acid N-hydroxysuccinimide ester (BODIPY FL-X SE) is supplied as a crystalline solid. A stock solution may be made by dissolving the BODIPY FL-X SE in the solvent of choice, which should be purged with an inert gas. BODIPY FL-X SE is soluble in organic solvents such as methanol and chloroform. The solubility of BODIPY FL-X SE in these solvents is approximately 1 mg/ml.

Description

BODIPY FL-X SE is an amine-reactive fluorescent probe.¹ BODIPY FL-X SE has been used to visualize proteins immobilized on PVDF membranes with minor excitation peaks at 300-360 nm, a primary excitation peak at 504 nm, and an emission maximum at 511 nm. It has also been used to increase the hydrophobicity and detectability of amino acids in transient trapping-micellar electrokinetic chromatography.²

References

1. Top, K.P., Hatleberg, G., Berggren, K.N., *et al.* Green/red dual fluorescence detection of total protein and alkaline phosphate-conjugated probes on blotting membranes. *Electrophoresis* **22(5)**, 896-905 (2001).
2. Sueyoshi, K., Hashiba, K., Kawai, T., *et al.* Hydrophobic labeling of amino acids: Transient trapping-capillary/microchip electrophoresis. *Electrophoresis* **32(10)**, 1233-1240 (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, 11/16/2022

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