

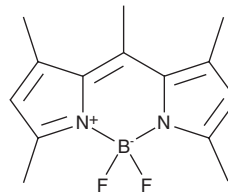
PRODUCT INFORMATION



BODIPY 493/503

Item No. 25892

CAS Registry No.: 121207-31-6
Formal Name: (T-4)-[2-[1-(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)ethyl]-3,5-dimethyl-1H-pyrrolato-κN]difluoro-boron
MF: C₁₄H₁₇BF₂N₂
FW: 262.1
Purity: ≥98%
Ex./Em. Max: 493/503 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 493/503 is supplied as a solid. A stock solution may be made by dissolving the BODIPY 493/503 in the solvent of choice, which should be purged with an inert gas. BODIPY 493/503 is soluble in organic solvents such as methanol, chloroform, DMSO, and dimethyl formamide.

Description

BODIPY 493/503 is a lipophilic fluorescent probe that localizes to polar lipids and can be used to label cellular neutral lipid contents, particularly those localized to lipid droplets, in live and fixed cells.^{1,2} BODIPY 493/503 is compatible with epifluorescent, confocal, and two-photon microscopy, as well as flow cytometry. It displays excitation/emission maxima of 493/503 nm, respectively, and can be used for live and fixed cell applications.

References

1. Listenberger, L.L. and Brown, D.G. Fluorescent detection of lipid droplets and associated proteins. *Curr. Protoc. Cell. Biol. Supplement* **35**, 24.2.1-24.2.11 (2007).
2. Qiu, B. and Simon, M.C. BODIPY 493/503 staining of neutral lipid droplets for microscopy and quantification by flow cytometry. *Bio. Protoc.* **6**(17), e1912 (2016).

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, 10/13/2022

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