

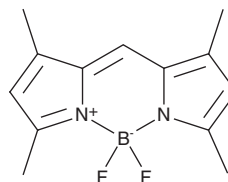
# PRODUCT INFORMATION



## BODIPY 505/515

Item No. 25893

**CAS Registry No.:** 21658-70-8  
**Formal Name:** (T-4)-[2-[(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)methyl]-3,5-dimethyl-1H-pyrrolato-κN]difluoro-boron  
**MF:** C<sub>13</sub>H<sub>15</sub>BF<sub>2</sub>N<sub>2</sub>  
**FW:** 248.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 228, 363, 504 nm  
**Ex./Em. Max:** 505/515 nm  
**Supplied as:** A crystalline solid  
**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 505/515 is supplied as a crystalline solid. A stock solution may be made by dissolving the BODIPY 505/515 in the solvent of choice, which should be purged with an inert gas. BODIPY 505/515 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of BODIPY 505/515 in these solvents is approximately 1.7 and 25 mg/ml, respectively.

BODIPY 505/515 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BODIPY 505/515 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BODIPY 505/515 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

BODIPY 505/515 is a lipophilic fluorescent probe that localizes to intracellular lipid bodies and has been used to label lipid droplets.<sup>1</sup> BODIPY 505/515 has been used for electron, epifluorescent, and confocal microscopy, as well as flow cytometry applications in various algae species.<sup>1,2</sup> It displays excitation/emission maxima of 505/515 nm, respectively, and has been used for live and fixed cell applications.

### References

1. Velmurugan, N., Sathishkumar, Y., Yim, S.S., *et al.* Study of cellular development and intracellular lipid bodies accumulation in the thraustochytrid *Aurantiochytrium* sp. KRS101. *Bioresour. Technol.* **161**, 149-154 (2014).
2. Rumin, J., Bonnefond, H., Saint-Jean, B., *et al.* The use of fluorescent Nile red and BODIPY for lipid measurement in microalgae. *Biotechnol. Biofuels* **8:42** (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, 01/11/2021

#### 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