

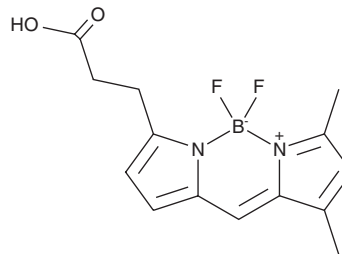
PRODUCT INFORMATION



BODIPY 503/512

Item No. 28800

CAS Registry No.: 165599-63-3
Formal Name: (T-4)-[5-[(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)methyl]-1H-pyrrole-2-propanoato(2-)-κN¹]difluoroborate(1-), monohydrogen
Synonyms: BDP FL, 3-BODIPY-Propanoic Acid
MF: C₁₄H₁₅BF₂N₂O₂
FW: 292.1
Purity: ≥95%
Ex./Em. Max: 503/512 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 503/512 is supplied as a solid. A stock solution may be made by dissolving the BODIPY 503/512 in the solvent of choice, which should be purged with an inert gas. BODIPY 503/512 is soluble in the organic solvent methanol.

Description

BODIPY 503/512 is a lipophilic, amine-reactive fluorescent probe.¹ It has been used to label oligonucleotide probes and primers for the quantitation of DNA and RNA by PCR and to monitor the uptake and trafficking of BODIPY-labeled proteins and other compounds within cells by fluorescence microscopy.¹⁻⁴ BODIPY 503/512 has also been used to identify and sort adipocytes from mouse white and brown adipose tissue by flow cytometry.⁵ It displays excitation/emission maxima of 503/512 nm, respectively.¹

References

1. Murakami, M., Cabral, H., Matsumoto, Y., *et al.* Improving drug potency and efficacy by nanocarrier-mediated subcellular targeting. *Sci. Transl. Med.* **3(64)**, 64ra2 (2011).
2. Kurata, S., Kanagawa, T., Yamada, K., *et al.* Fluorescent quenching-based quantitative detection of specific DNA/RNA using a BODIPY® FL-labeled probe or primer. *Nucleic Acids Res.* **29(6)**, E34 (2001).
3. Bittel, A.M., Saldívar, I.S., Dolman, N.J., *et al.* Superresolution microscopy with novel BODIPY-based fluorophores. *PLoS One* **13(10)**, e0206104 (2018).
4. Surewaard, B.G.J., Deniset, J.F., Zemp, F.J., *et al.* Identification and treatment of the *Staphylococcus aureus* reservoir in vivo. *J. Exp. Med.* **213(7)**, 1141-1151 (2016).
5. Hagberg, C.E., Li, Q., Kutschke, M., *et al.* Flow cytometry of mouse and human adipocytes for the analysis of browning and cellular heterogeneity. *Cell Rep.* **24(10)**, 2746-2756 (2018).

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, 11/30/2022

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