

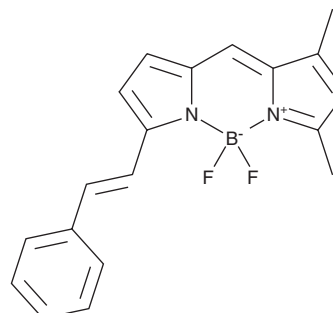
PRODUCT INFORMATION



STY-BODIPY

Item No. 27089

CAS Registry No.: 2383063-37-2
Formal Name: (T-4)-[2-[(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)methyl]-5-[(1E)-2-phenylethenyl]-1H-pyrrolato-κN]difluoro-boron
Synonyms: Styrene-BODIPY, Styrene-Conjugated BODIPY
MF: C₁₉H₁₇BF₂N₂
FW: 322.2
Purity: ≥98%
UV/Vis.: λ_{max}: 240, 313, 562 nm
Abs. Max: 571 nm
Supplied as: A solution in benzene
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

STY-BODIPY is a styrene-conjugated fluorogenic probe for radical-trapping antioxidant (RTA) activity.¹ Co-autoxidation of the STY-BODIPY signal carrier and a hydrocarbon co-substrate can be quantified by monitoring the loss of absorbance at 571 nm. STY-BODIPY has been used to measure the activity of RTAs, as well as the kinetics and stoichiometry of RTA reactions in cell-free assays.¹⁻³

References

1. Haidasz, E.A., Van Kessel, A.T.M., and Pratt, D.A. A continuous visible light spectrophotometric approach to accurately determine the reactivity of radical-trapping antioxidants. *J. Org. Chem.* **81**(3), 737-744 (2016).
2. Shah, R., Shchepinov, M.S., and Pratt, D.A. Resolving the role of lipoxygenases in the initiation and execution of ferroptosis. *ACS Cent. Sci.* **4**(3), 387-396 (2018).
3. Chauvin, J.-P.R., Haidasz, E.A., Griesser, M., *et al.* Polysulfide-1-oxides react with peroxy radicals as quickly as hindered phenolic antioxidants and do so by a surprising concerted homolytic substitution. *Chem. Sci.* **7**(10), 6347-6356 (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, 11/18/2020

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