

# PRODUCT INFORMATION



## C11 BODIPY 581/591

Item No. 27086

**CAS Registry No.:** 217075-36-0  
**Formal Name:** (T-4)-difluoro[5-[[5-[(1E,3E)-4-phenyl-1,3-butadien-1-yl]-2H-pyrrol-2-ylidene-κN]methyl]-1H-pyrrole-2-undecanoato(2-)-κN<sup>1</sup>]-borate(1-), monohydrogen

**MF:** C<sub>30</sub>H<sub>34</sub>BF<sub>2</sub>N<sub>2</sub>O<sub>2</sub> • H

**FW:** 504.4

**Purity:** ≥95%

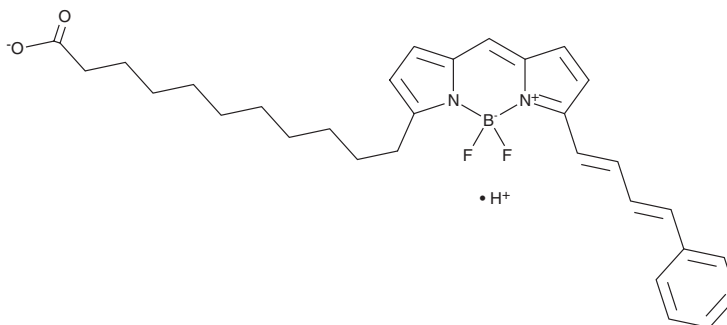
**UV/Vis.:** λ<sub>max</sub>: 581 nm

**Ex./Em. Max:** 581/591 nm, respectively, when reduced;  
500/510 nm, respectively, when oxidized

**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

C11 BODIPY 581/591 is supplied as a solid. A stock solution may be made by dissolving the C11 BODIPY 581/591 in the solvent of choice, which should be purged with an inert gas. C11 BODIPY 581/591 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml. C11 BODIPY 581/591 is slightly soluble in methanol.

### Description

C11 BODIPY 581/591 is a lipid-soluble ratiometric fluorescent indicator of lipid oxidation.<sup>1</sup> Upon oxidation, its excitation maximum shifts from 581 to 500 nm and the emission maximum shifts from 591 to 510 nm. C11 BODIPY 581/591 has been used in the quantitation of ferroptosis.<sup>1,2</sup>

### References

1. Pap, E.H.W., Drummen, G.P.C., Winter, V.J., *et al.* Ratio-fluorescence microscopy of lipid oxidation in living cells using C11-BODIPY<sup>581/591</sup>. *FEBS Lett.* **453(3)**, 278-282 (1999).
2. Naguib, Y.M.A. A fluorometric method for measurement of peroxy radical scavenging activities of lipophilic antioxidants. *Anal. Biochem.* **265(2)**, 290-298 (1998).

#### 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, 10/07/2022

#### 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