

# PRODUCT INFORMATION



## BODIPY 630/650X succinimide ester

Item No. 27758

**CAS Registry No.:** 380367-48-6  
**Formal Name:** (T-4)-[2,5-dioxo-1-pyrrolidinyl 6-[[2-[4-[2-[5-[[5-(2-thienyl)-2H-pyrrol-2-ylidene-κN]methyl]-1H-pyrrol-2-yl-κN]ethenyl]phenoxy]acetyl]amino] hexanoato]difluoro-boron

**Synonym:** BODIPY 630/650X SE

**MF:** C<sub>33</sub>H<sub>31</sub>BF<sub>2</sub>N<sub>4</sub>O<sub>6</sub>S

**FW:** 660.5

**Purity:** ≥90%

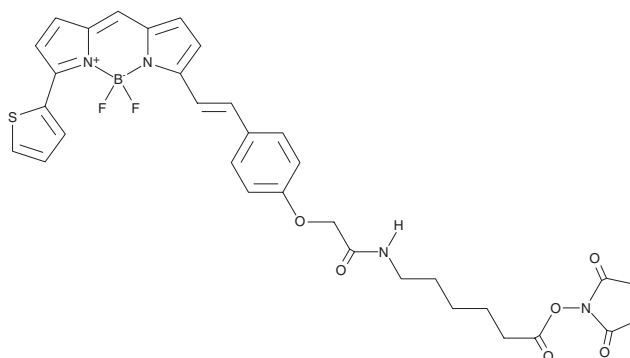
**UV/Vis.:** λ<sub>max</sub>: 628 nm

**Ex./Em. Max:** 630/650 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 630/650X succinimide ester (SE) is supplied as a solid. A stock solution may be made by dissolving the BODIPY 630/650X SE in the solvent of choice, which should be purged with an inert gas. BODIPY 630/650X SE is soluble in the organic solvent methanol.

### Description

BODIPY 630/650X SE is an amine-reactive fluorescent probe that contains an integrated aminohexanoyl linker (X).<sup>1,2</sup> It displays excitation/emission maxima of 630/650 nm, respectively. BODIPY 630/650X SE has been used, linked to adenosine A<sub>1</sub> receptor ligands, as a fluorescent probe to quantify ligand-receptor binding using fluorescence correlation spectroscopy.<sup>1,2</sup>

### References

1. Briddon, S.J., Middleton, R.J., Cordeaux, Y., *et al.* Quantitative analysis of the formation and diffusion of A<sub>1</sub>-adenosine receptor-antagonist complexes in single living cells. *Proc. Natl. Acad. Sci. U.S.A.* **101**(13), 4673-4678 (2004).
2. Dale, C.L., Hill, S.J., and Kellam, B. New potent, short-linker BODIPY-630/650™ labelled fluorescent adenosine receptor agonists. *Med. Chem. Commun.* **3**(3), 333-338 (2012).

#### 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, 12/05/2022

#### 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