

PRODUCT INFORMATION



FBBBE

Item No. 14606

CAS Registry No.: 1522117-83-4
Formal Name: 3',6'-bis((4-(4,4,5,5-tetramethyl-1,3,2-dioxaborolan-2-yl)benzyl)oxy)-3H-spiro[isobenzofuran-1,9'-xanthen]-3-one

Synonym: Fluorescein bis (benzyl boronic ester)

MF: C₄₆H₄₆B₂O₉

FW: 764.5

Purity: ≥95%

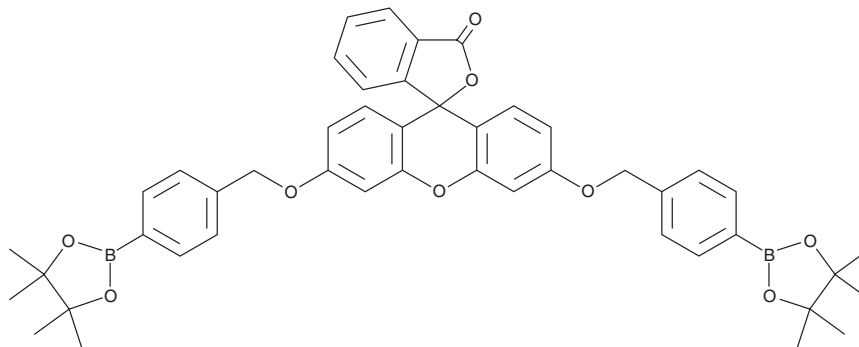
UV/Vis.: λ_{max}: 230, 275 nm

Ex./Em. Max: 480/512 nm

Supplied as: A crystalline 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

FBBBE is supplied as a crystalline solid. A stock solution may be made by dissolving the FBBBE in the solvent of choice. FBBBE is soluble in organic solvents such as DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of FBBBE in these solvents is approximately 25 and 30 mg/ml, respectively.

Description

FBBBE is a derivatized fluorescein compound that is used to detect the production hydrogen peroxide (H₂O₂) by cells. Intracellular H₂O₂, but not extracellular H₂O₂, initiates the conversion of FBBBE to fluorescein, resulting in an increase in intracellular fluorescence (excitation/emission maxima: 480/512 nm).¹

Reference

1. Daniel, K.B., Agrawal, A., Manchester, M., *et al.* Readily accessible fluorescent probes for sensitive biological imaging of hydrogen peroxide. *ChemBioChem* **14**(5), 593-598 (2013).

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, 01/16/2024

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