

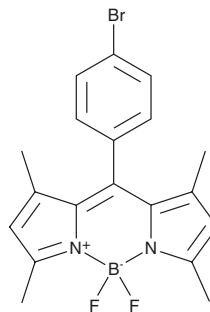
PRODUCT INFORMATION



8(4'-bromophenyl)-1,3,5,7-tetramethyl-BODIPY

Item No. 40123

CAS Registry No.: 850534-66-6
Formal Name: (T-4)-[2-[(4-bromophenyl)(3,5-dimethyl-2H-pyrrol-2-ylidene-κN)methyl]-3,5-dimethyl-1H-pyrrolato-κN]difluoro-boron
MF: C₁₉H₁₈BBrF₂N₂
FW: 403.1
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years
Special Conditions: Light and oxygen sensitive



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

Laboratory Procedures

8(4'-bromophenyl)-1,3,5,7-tetramethyl-BODIPY is supplied as a solid. A stock solution may be made by dissolving the 8(4'-bromophenyl)-1,3,5,7-tetramethyl-BODIPY in the solvent of choice, which should be purged with an inert gas. 8(4'-bromophenyl)-1,3,5,7-tetramethyl-BODIPY is soluble in acetonitrile, chloroform, and methanol.

Description

8(4'-bromophenyl)-1,3,5,7-tetramethyl-BODIPY is an intermediate in the synthesis of BODIPY-containing fluorescent dyes.¹ It has also been used as a precursor in the synthesis of BODIPY-containing photosensitizer dyes with antibacterial activity.²

References

1. Guo, Y., Liang, Y., Wang, Y., *et al.* Orthogonal conformation regulation enables highly efficient phosphorescence emission in BODIPY-fluorene derivatives without heavy atoms. *Dyes and Pigments* **208**, 110791 (2023).
2. Sen, P., Sindelo, A., Nnaji, N., *et al.* Diiodinated mono- and dipyritylvinyl BODIPY dyes: Photophysical properties, *in vitro* antibacterial studies, molecular docking and theoretical calculations. *Photochem. Photobiol.* **99(3)**, 947-956 (2023).

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, 11/16/2023

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