PRODUCT INFORMATION



PBD-BODIPY

Item No. 27945

CAS Registry No.:	148185-52-8	/
Formal Name:	(T-4)-[3,5-dimethyl-2-[[5-(4-phenyl-1,3-	
	butadienyl)-2H-pyrrol-2-ylidene]methyl]-1H- pyrrolato-N ¹ ,N ²]difluoro-boron	
MF:	$C_{21}H_{19}BF_2N_2$	
FW:	348.2	
Purity:	≥95%	
UV/Vis.:	λ _{max} : 255, 329, 543, 580 nm	/-
Abs. Max:	591 nm	
Supplied as:	A solution in methyl acetate	L 11
Storage:	-20°C	
Stability:	≥2 years	

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

Description

PBD-BODIPY is a probe for the spectrophotometric measurement of autoxidation reactions.¹ Co-autoxidation of the PBD-BODIPY signal carrier and a hydrocarbon co-substrate can be quantified by monitoring loss of absorbance at 591 nm.^{1,2} PBD-BODIPY has been used to measure the activity of radicaltrapping antioxidants in cell-free assays. It has also been used as a fluorescent probe for the detection of epoxidation activity.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. Zilka, O., Shah, R., Li, B., et al. On the mechanism of cytoprotection by ferrostatin-1 and liproxstatin-1 and the role of lipid peroxidation in ferroptotic cell death. ACS Cent. Sci. 3(3), 232-243 (2017).
- 3. De Cremer, G., Bartholomeeusen, E., Pescarmona, P.P., et al. The influence of diffusion phenomena on catalysis: A study at the single particle level using fluorescence microscopy. Catal. Today 157(1-4), 236-242 (2010).

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.

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, 04/25/2024

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