

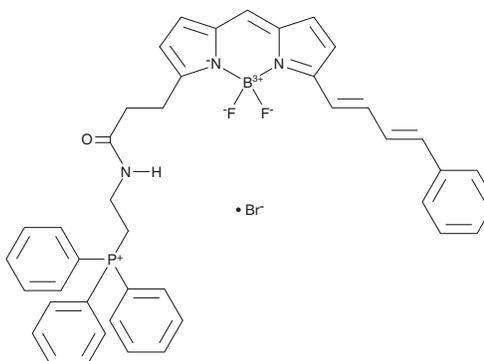
# PRODUCT INFORMATION



## MitoPerOx

Item No. 18798

**CAS Registry No.:** 1392820-50-6  
**Formal Name:** (2-(3-(5,5-difluoro-7-((1E,3E)-4-phenylbuta-1,3-dien-1-yl)-5H-5λ<sup>4</sup>,6λ<sup>4</sup>-dipyrrolo[1,2-c:2',1'-f][1,3,2]diazaborinin-3-yl)propanamido)ethyl)triphenylphosphonium, monobromide  
**MF:** C<sub>42</sub>H<sub>38</sub>BF<sub>2</sub>N<sub>3</sub>OP • Br  
**FW:** 760.5  
**Purity:** ≥90%  
**Ex./Em. Max:** 495/520-590 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

MitoPerOx is supplied as a solid. A stock solution may be made by dissolving the MitoPerOx in the solvent of choice, which should be purged with an inert gas. MitoPerOx is soluble in the organic solvent DMSO at a concentration of approximately 25 mM.

### Description

MitoPerOx is a ratiometric fluorescent probe that can be used to assess changes in lipid peroxidation within mitochondria.<sup>1</sup> It is composed of a BODIPY fluorophore conjugated *via* a dienyl link to a triphenylphosphonium cation component that drives its accumulation in mitochondria. MitoPerOx displays an excitation maximum of 495 nm and exhibits a shift in emission maxima from 590 to 520 nm upon mitochondrial lipid peroxidation, enabling determination of ratiometric measurements of lipid peroxidation in live cells.

### Reference

1. Prime, T.A., Forkink, M., Logan, A., *et al.* A ratiometric fluorescent probe for assessing mitochondrial phospholipid peroxidation within living cells. *Free Radic. Biol. Med.* **53(3)**, 544-553 (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, 07/11/2023

#### 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