

Bodi Fluor™ 581/591 C11 *Lipid Peroxidation Probe*

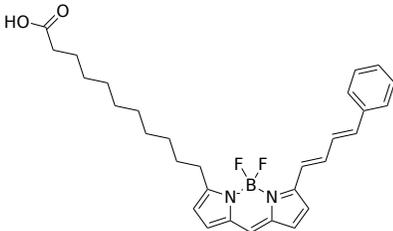
Catalog Number: 23290

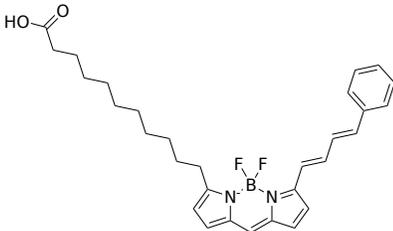
Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid deep blue
Molecular Weight	504.43
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	589 nm
Emission Wavelength	597 nm

Applications

Bodi Fluor™ 581/591 C11 is the same molecule as Bodipy 581/591 C11. It is a specific type of fluorescent dye used in various biological and biochemical applications. The "581/591" in the name refers to the wavelengths at which this particular Bodi Fluor dye exhibits maximum excitation (581 nm) and emission (591 nm). The "C11" part of the name typically indicates that the dye contains an 11-carbon aliphatic chain, which enhances its lipophilicity and makes it suitable for lipid-related studies. The lipophilic nature of the C11 moiety allows the dye to readily incorporate into cell membranes and other lipid-rich environments. One of the common applications of Bodi Fluor™ 581/591 C11 is to measure lipid peroxidation, a process involving the degradation of lipids through the action of reactive oxygen species (ROS). The dye's fluorescence changes with the oxidation state of lipids, and this unique property is exploited to assess lipid peroxidation levels in cells and tissues.