

Product Information Sheet

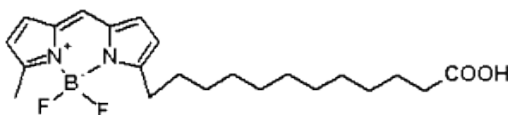
Ordering Information

Product Number:	22120
Product Name:	C1-Bodi Fluor 500/510-C12 [equivalent to BODIPY [®] 500/510 C1, C12]
Unit Size:	5 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 418.33

Chemical Structure:



Soluble in: DMSO

Excitation Wavelength: 503

Emission Wavelength: 512

Application Notes

The green fluorescent BODIPY FL C12 fatty acid can be used as a synthetic precursor to a wide variety of fluorescent phospholipids. BODIPY dyes are relatively environment insensitive and is fluorescent in both aqueous and lipid environments with sharp emission peak.