

Bodi Fluor™ 488 acid [equivalent to Bodipy® FL] *CAS 165599-63-3*

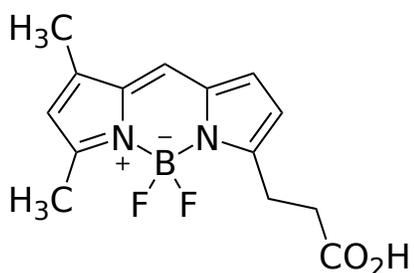
Catalog Number: 700

Unit Size: 10 mg

Product Details

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

Chemical Properties

Appearance	Solid red
Molecular Weight	292.09
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	504 nm
Emission Wavelength	510 nm

Applications

Bodi Fluor™ 488 is the same molecule to Bodipy® FL (Bodipy® is the trademark of Invitrogen). Bodi Fluor™ 488 has the excitation and emission spectra similar to those of fluorescein. Compared to fluorescein, Bodi Fluor™ 488 is a neutral molecule, and its fluorescence is not pH-dependent. Bodi Fluor™ 488 gives bright green fluorescence with a sharp emission peak. Bodi Fluor™ 488 has been used for developing fluorescent intracellular tracers. Its conjugates are cell-permeable as long as its carriers are cell-permeable.