

# Bodi Fluor™ 488 SE [equivalent to Bodipy® FL, SE] \*CAS 146616-66-2\*

Catalog Number: 701

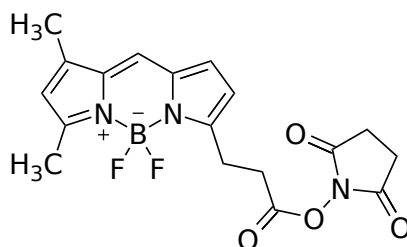
Unit Size: 5 mg

## Product Details

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

## Chemical Properties

Appearance	Solid red
Molecular Weight	389.16
Soluble In	DMSO
Chemical Structure	



## Spectral Properties

Excitation Wavelength	504 nm
Emission Wavelength	510 nm

## Applications

Bodi Fluor™ 488, SE is the same molecule to Bodipy® FL, SE (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.