

Bodi Fluor™ R6G NHS Ester [equivalent to Bodipy® R6G, NHS Ester]

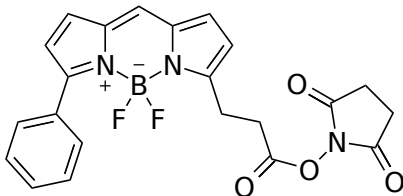
Catalog Number: 703

Unit Size: 5 mg

Product Details

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

Chemical Properties

Appearance	Solid dark red
Molecular Weight	437.21
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	528 nm
Emission Wavelength	547 nm

Applications

Bodi Fluor™R6G, NHS ester is the same molecule to Bodipy® R6G, NHS ester (Bodipy® is the trademark of Invitrogen). Bodi Fluor™R6G has the excitation and emission spectra similar to rhodamine 6G. Compared to rhodamine 6G, Bodi Fluor™R6G is a neutral molecule, and its fluorescence is not pH-dependent. Bodi Fluor™R6G gives bright green fluorescence with a sharp emission peak. Bodi Fluor™R6G has been used for developing fluorescent intracellular tracers. Its conjugates are cell-permeable as long as its carriers are cell-permeable.