

XFD488 TCO

Catalog Number: 4601

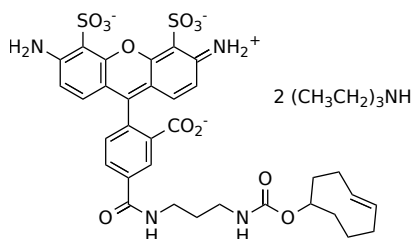
Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid orange-red
Molecular Weight	945.16
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	499 nm
Emission Wavelength	520 nm

Applications

XFD488 contains the same fluorophore of Alexa Fluor® 488. It is a bright, green-fluorescent dye with excitation ideally suited for the 488 nm laser line. It is used for stable signal generation in imaging and flow cytometry. AF488 TCO demonstrates exceptionally fast reaction with tetrazines, the fastest kinetics ever reported for any biorthogonal reaction. It provides a convenient tool for making biological conjugates with tetrazine-modified substrates. The resulted conjugates provide the bright, green fluorescence of Alexa Fluor® 488 (Alexa Fluor® is the trademark of ThermoFisher Scientific).