

XFD488 tetrazine

Catalog Number: 4600

Unit Size: 1 mg

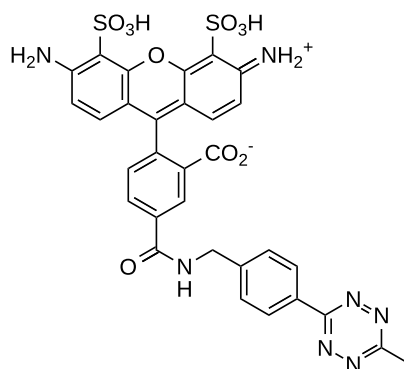
Product Details

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

Chemical Properties

Appearance	Solid dark red
Molecular Weight	717.68
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	499 nm
Emission Wavelength	520 nm

Applications

XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). XFD488 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. XFD488 tetrazine demonstrates exceptionally fast reaction with trans-cyclooctenes (TCO) as the dienophile, the fastest kinetics ever reported for any biorthogonal reaction. It provides a convenient tool for making biological conjugates with TCO-modified substrates. The resulted conjugates provide the bright, green fluorescence of XFD488.