

XFD488 acid

Catalog Number: 1790

Unit Size: 5 mg

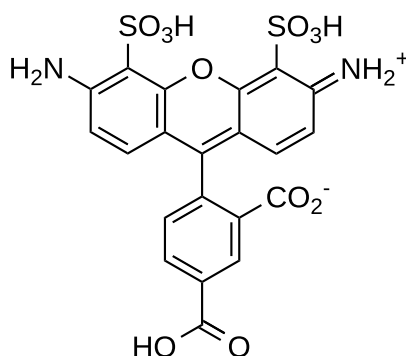
Product Details

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

Chemical Properties

Appearance	Solid orange
Molecular Weight	534.47
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 acid has a similar molecule as Alexa Fluor® 488 acid. It is a bright yellow-fluorescent dye optimal for use with the 488 nm Argon laser line. XFD488 dye is water soluble and pH-insensitive from pH 4 to pH 10. XFD488 acid is the most economic amine-reactive form for conjugating this dye to a peptide.