

XFD488 amine

Catalog Number: 1705

Unit Size: 1 mg

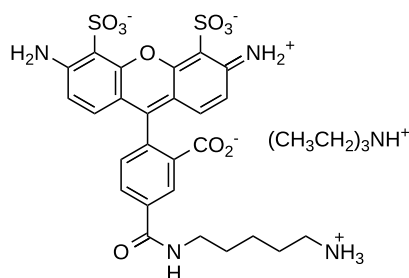
Product Details

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

Chemical Properties

Appearance	Solid orange
Molecular Weight	719.83
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	499 nm
Emission Wavelength	520 nm

Applications

XFD488 amine is the same molecule to Alexa Fluor® 488 C5 amine. The fluorescent dye amines are reactive molecules that can be used to add a fluorescent label to biomolecules containing a carbonyl group. XFD488 amine reacts with carboxylic acids, aldehydes and ketones. It is a very useful building block for developing XFD488 probes.