

XFD555 amine

Catalog Number: 1707

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	1103.10
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	553 nm
Emission Wavelength	568 nm

Applications

XFD555 amine is the same molecule to the Alexa Fluor® 555 amine (Alexa Fluor® is the trademark of ThermoFisher). It is a cyanine dye with bright yellow-fluorescence optimal for labeling antibodies. XFD555 dye is water soluble and pH-insensitive from pH 4 to 10. The amine of XFD555 is a versatile carbonyl-reactive form for conjugating this fluorophore to peptides and other biological molecules.