

XFD750 acid

Catalog Number: 1798

Unit Size: 10 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	999.07
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	752 nm
Emission Wavelength	776 nm

Applications

XFD750 acid is the same molecule to the Alexa Fluor® 750 acid (Alexa Fluor® is the trademark of ThermoFisher). It is a cyanine dye with bright NIR-fluorescence optimal for labeling antibodies. It has the identical spectra to Cy7 dye. XFD750 dye is water soluble and pH-insensitive from pH 4 to 10. XFD750 acid is the most economic amine-reactive form for conjugating this dye to a peptide.