

XFD555 acid

Catalog Number: 1796

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	946.99
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	553 nm
Emission Wavelength	568 nm

Applications

XFD555 acid is the same molecule to the Alexa Fluor® 555 acid (Alexa Fluor® is the trademark of ThermoFisher). It is a cyanine dye with bright yellow-fluorescent dye optimal for labeling antibodies. XFD555 dye is water soluble and pH-insensitive from pH 4 to 10. XFD555 acid is the most economic amine-reactive form for conjugating this dye to a peptide.