

XFD488 azide

Catalog Number: 1701

Unit Size: 1 mg

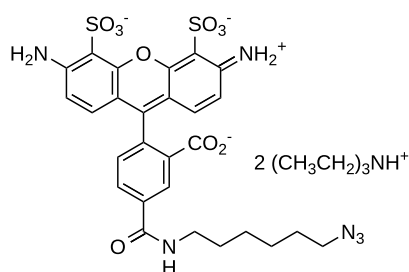
Product Details

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

Chemical Properties

Appearance	Solid orange
Molecular Weight	861.04
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	499 nm
Emission Wavelength	520 nm

Applications

The fluorescent dye azides react with alkynes via a copper-catalyzed click reaction (CuAAC). This click chemistry is increasingly being used in a variety of biological applications. XFD488 azide has a chemical structure similar to Alexa Fluor® 488 azide (Alexa Fluor® is the Trademark of Thermo Fisher Scientific). XFD488 azide is reactive with terminal alkynes via a copper-catalyzed click reaction as a bio-orthogonal or biologically unique hapten for use in applications requiring signal amplification.