

# XFD488 alkyne

Catalog Number: 1700

Unit Size: 1 mg

## Product Details

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

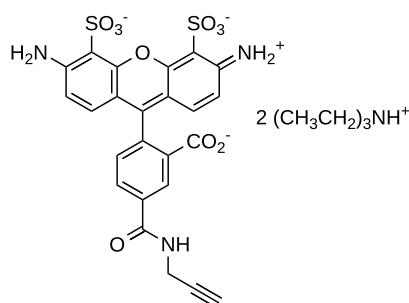
## Chemical Properties

Appearance Solid orange

Molecular Weight 773.92

Soluble In DMSO

Chemical Structure



## Spectral Properties

Excitation Wavelength 499 nm

Emission Wavelength 520 nm

## Applications

The fluorescent dye alkynes react with azides via a copper-catalyzed click reaction (CuAAC). This click chemistry is increasingly being used in a variety of biological applications. XFD488 alkyne has a chemical structure similar to Alexa Fluor® 488 alkyne (Alexa Fluor® is the Trademark of Thermo Fisher Scientific). XFD488 alkyne 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.