

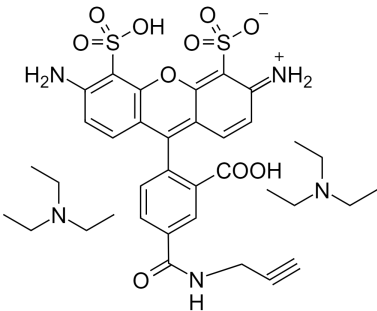
## Alpha Fluor™ 488 alkyne

Catalog number: 1700  
Unit size: 1 mg

### Product Details

Storage Conditions	Freeze (<math>< -15\text{ }^{\circ}\text{C}</math>), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

Appearance	Orange solid
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. Alpha Fluor™ 488 alkyne has the same chemical structure to Alexa Fluor® 488 alkyne (Alexa Fluor® is the Trademark of ThermoFisher Scientific). Alpha Fluor™ 488 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.