

## Product Information Sheet

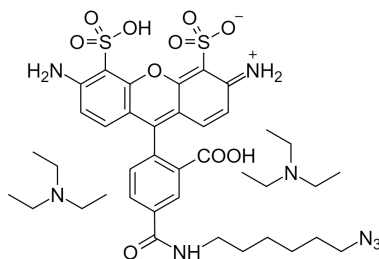
### Ordering Information

Product Number: 1701  
Product Name: Alpha Fluor™ 488 azide  
Unit Size: 1 mg  
Storage Conditions: Freeze (<-15 °C), Desiccated, Avoid Light  
Expiration Date: 12 months upon receiving

### Chemical and Spectral Properties

Appearance: Solid  
Molecular Weight: 861.04

Chemical Structure:



Soluble In: DMSO  
Excitation Wavelength: 494  
Emission Wavelength: 517

### Application Notes

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. 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.