

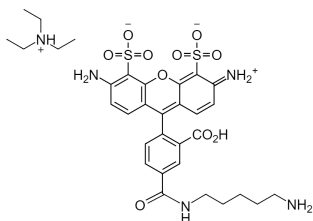
## Alpha Fluor™ 488 amine

Catalog number: 1705  
Unit size: 1 mg

### Product Details

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

### Chemical Properties

Appearance	Orange solid
Molecular Weight	719.83
Soluble In	DMSO
Chemical Structure	 <p>5- and 6-isomers</p>

### Spectral Properties

Excitation Wavelength	499 nm
Emission Wavelength	520 nm

### Applications

Alpha Fluor™ 488 amine is the same molecule to Alexa Fluor® 488 C5 amine. The fluorescent dye amines are reactive molecules that can be used to add a fluorescent label to biomolecules containing a carbonyl group. Alpha Fluor™ 488 amine reacts with carboxylic acids, aldehydes and ketones. It is a very useful building block for developing Alpha Fluor™ 488 probes.