
Product Information Sheet

Ordering Information

Product Number:	1790
Product Name:	Alpha Fluor™ 488 acid [equivalent to Alexa Fluor™ 488 acid]
Unit Size:	5 mg
Storage Conditions:	Freeze (<-15 °C), Desiccated, Avoid Light
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Appearance:	Solid
Molecular Weight:	626.70
Soluble In:	DMSO
Excitation Wavelength:	530
Emission Wavelength:	555

Application Notes

Alpha Fluor™ 488 acid is the same molecule to the Alexa Fluor® 4882 acid (Alexa Fluor® is the trademark of ThermoFisher). It is a bright yellow-fluorescent dye optimal for use with the 488 nm Argon laser line. Alpha Fluor™ 488 dye is water soluble and pH-insensitive from pH 4 to pH 10. Alpha Fluor™ 488 is the most economic amine-reactive form for conjugating this dye to a peptide.