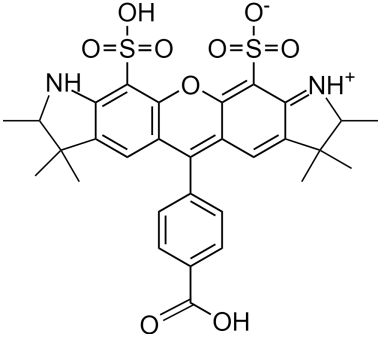
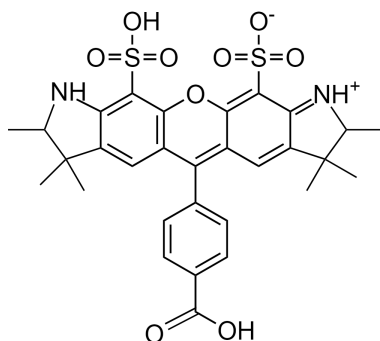


**Alpha Fluor™ 532 acid [equivalent to Alexa Fluor™ 532 acid]**Catalog number: 1795  
Unit size: 10 mg**Product Details**

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

**Chemical Properties**

Appearance	Red solid
Molecular Weight	626.70
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	534 nm
Emission Wavelength	553 nm

**Applications**

Alpha Fluor™ 532 acid is the same molecule to the Alexa Fluor® 532 acid (Alexa Fluor® is the trademark of ThermoFisher). It is a bright yellow-fluorescent dye optimal for use with the frequency-doubled Nd:YAG laser line. Alpha Fluor™ 532 dye is water soluble and pH-insensitive from pH 4 to pH 10. Alpha Fluor™ 532 is the most economic amine-reactive form for conjugating this dye to a peptide.