

Product Information Sheet

Ordering Information

Product Number: 1823

Product Name: Alpha Fluor™ 568 NHS Ester [equivalent to Alexa Fluor™ 568 NHS Ester]

Unit Size: 1 mg

Storage Conditions:

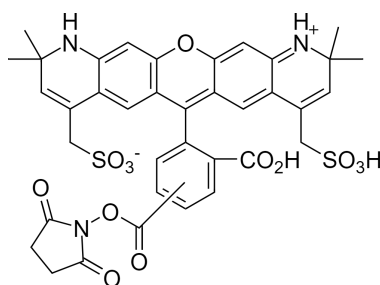
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance: Solid

Molecular Weight: 791.80

Chemical Structure:



Soluble In: DMSO

Excitation Wavelength: 579

Emission Wavelength: 603

Application Notes

Alpha Fluor™ 568 NHS Ester (Succinimidyl Ester) is the same molecule to the Alexa Fluor® 568 NHS Ester (Alexa Fluor® is the trademark of ThermoFisher). It is a bright orange-fluorescent dye optimal for use with the frequency-doubled Nd:YAG laser line. Alpha Fluor™ 568 dye is water soluble and pH-insensitive from pH 4 to pH 10. The NHS ester (or succinimidyl ester) of Alpha Fluor™ 568 is the most convenient amine-reactive form for conjugating this dye to a protein or antibody.