

Product Information Sheet

Ordering Information

Product Number: 1833

Product Name: AF647 NHS Ester [equivalent to Alexa Fluor™ 647 NHS Ester]

Unit Size: 1 mg

Storage Conditions: Freeze (<-15 °C), Minimize light exposure

Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance: Solid

Molecular Weight: ~1250

Soluble In: DMSO

Excitation Wavelength: 651

Emission Wavelength: 672

Application Notes

AF647 NHS Ester (Succinimidyl Ester) is the same molecule to the Alexa Fluor® 647 NHS Ester (Alexa Fluor® is the trademark of ThermoFisher). It is a bright red fluorescent dye. AF647 dye is water soluble and pH-insensitive from pH 4 to pH 10. The NHS ester (or succinimidyl ester) of AF647 is the most convenient amine-reactive form for conjugating this dye to a protein or antibody.