
Product Information Sheet

Ordering Information

Product Number:	1895
Product Name:	AF647 maleimide [equivalent to Alexa Fluor™ C2 647 maleimide]
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:	1095.15
Soluble In:	DMSO
Excitation Wavelength:	N/A
Emission Wavelength:	N/A

Application Notes

AF647 C2 Maleimide is the same molecule to the Alexa Fluor® 647 C2 Maleimide (Alexa Fluor® is the trademark of ThermoFisher). It is a bright red-fluorescent dye optimal for use with the frequency-doubled He-Ne laser line. Alpha Fluor™ 594 dye is water soluble and pH-insensitive from pH 4 to pH 10. The maleimide of AF647 is the most convenient thiol-reactive form for conjugating this dye to a protein or antibody.