

iFluor™ 546 maleimideCatalog number: 1059
Unit size: 1 mg**Product Details**

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

Chemical Properties

Appearance	Solid
Molecular Weight	1082.10
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	541 nm
Emission Wavelength	557 nm

Applications

AAT Bioquest's iFluor™ dyes are optimized for labeling proteins, in particular, antibodies. These dyes are bright, photostable and have minimal quenching on proteins. iFluor™ 546 is optimized to be well excited by the 546 nm laser or LED light sources, enabling new biological applications for fluorescence instruments that is equipped with the 546 nm 546 nm laser or LED light sources. iFluor™ 546 dye might be the brightest fluorophore available that is well excited by the 546 nm laser or LED light sources. iFluor™ 546 maleimide is stable and shows good reactivity and selectivity with thiol group.