

iFluor® 670 maleimide

Catalog Number: 72502

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	971.03
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

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. They can be well excited by the major laser lines of fluorescence instruments (e.g., 350, 405, 488, 514, 555, 633, 647 and 808 nm). iFluor® 670 dyes have fluorescence excitation and emission maxima of ~670 nm and ~690 nm respectively. These spectral characteristics make them a unique acceptor for preparing PE, APC and PerCP tandems that are widely used in flow cytometry applications. iFluor® 670 dyes have excellent photostability. iFluor® 670 Maleimide is stable and shows good reactivity and selectivity with protein thiol group.