

Product Information Sheet

Ordering Information

Product Number:	1092
Product Name:	iFluor™ 555 alkyne
Unit Size:	1 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	24 months upon receiving

Chemical and Spectral Properties

Molecular Weight:	781.86
Soluble in:	DMSO
Excitation Wavelength:	555 nm
Emission Wavelength:	565 nm

Application Notes

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, 555 and 633 nm). iFluor™ 555 dyes have fluorescence excitation and emission maxima of ~550 nm and ~570 nm respectively. iFluor™ 555 family has the spectral properties essentially identical to those of Cy3® (Cy3® is the trademark of GE Healthcare). Compared to Cy3 probes iFluor™ 555 family has much stronger fluorescence and higher photostability. Their fluorescence is pH-independent from pH 3 to 11. These spectral characteristics make this new dye family a superior alternative to Cy3®. iFluor™ 555 family has become an excellent replacement for Cy3® and Alexa Fluor® 555 labeling dye (Cy3® and Alexa Fluor® are the trademark of Invitrogen and GE Health Care). iFluor™ 555 alkyne is reasonably stable and shows good reactivity and selectivity with azido group in click chemistry.