

Product Information Sheet

Ordering Information

Product Number:	1077
Product Name:	iFluor™ 700 amine
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:	1049.02
Soluble in:	DMSO
Excitation Wavelength:	685 nm
Emission Wavelength:	710 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™ 700 dyes have fluorescence excitation and emission maxima of ~690 nm and ~710 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 700 labeling dyes (Alexa Fluor® is the trademark of Invitrogen). iFluor™ 700 amine is stable and used for modifying carbonyl groups (e.g., aldehyde and carboxy groups).