

## Product Information Sheet

### Ordering Information

Product Number:	1079
Product Name:	iFluor™ 750 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:	1047.07
Soluble in:	DMSO
Excitation Wavelength:	749 nm
Emission Wavelength:	775 nm

### Application Notes

AAT Bioquest's iFluor™ dyes are developed for labeling proteins, in particular, antibodies. These dyes are optimized to have minimal fluorescence quenching effect on proteins and nucleic acids. Our iFluor™ 750 dyes have fluorescence excitation and emission maxima close to 750 nm and 780 nm respectively with good photostability. Our in-house comparable studies indicated that our iFluor™ 750 dyes are significantly brighter than the corresponding Cy7® and Alexa Fluor® 750. These spectral characteristics make them a superior alternative to Cy7® and Alexa Fluor® 750 (Cy7® and Alexa Fluor® are the trademarks of GE Healthcare and Invitrogen respectively). iFluor™ 750 conjugates have been widely used in fluorescence animal imaging applications. The iFluor™ 750 amine is stable and used for modifying carbonyl groups (e.g., aldehyde and carboxy groups).