

## Product Information Sheet

### Ordering Information

Product Number: 1078  
Product Name: iFluor<sup>™</sup> 710 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: 1001.97  
Soluble in: DMSO  
Excitation Wavelength: 710 nm  
Emission Wavelength: 736 nm

### Application Notes

AAT Bioquest's iFluor<sup>™</sup> 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<sup>™</sup> 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<sup>®</sup> 700 labeling dyes (Alexa Fluor<sup>®</sup> is the trademark of Invitrogen). iFluor<sup>™</sup> 710 amine is stable and used for modifying carbonyl groups (e.g., aldehyde and carboxy groups).