

## Product Information Sheet

### **Ordering Information**

Product Number: 1077

Product Name: iFluor™ 700 amine

Unit Size: 1 mg

Storage Conditions: Freeze (<-15 °C), Minimize light exposure

Expiration Date: 12 months upon receiving

### **Chemical and Spectral Properties**

Appearance: Solid

Molecular Weight: 1049.02

Soluble In: DMSO

Excitation Wavelength: 685

Emission Wavelength: 710

### **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).