



Product Information Sheet

Ordering Information

Product Number: 1071

Product Name: iFluor™ 405 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: 748.68

Soluble In: DMSO

Excitation Wavelength: 402

Emission Wavelength: 425

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™ 405 dyes have fluorescence excitation and emission maxima of ~400 nm and ~420 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 405 labeling dye (Alexa Fluor® is the trademark of Invitrogen). iFluor™ 405 amine is stable and used for modifying carbonyl groups (e.g., aldehyde and carboxy groups).