

iFluor™ 405 azideCatalog number: 1094
Unit size: 1 mg**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	660.64
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	403 nm
Emission Wavelength	427 nm

Applications

iFluor® 405 azide is reasonably stable and shows good reactivity and selectivity with alkyne group in click chemistry. 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 ~405 nm and ~420 nm respectively. iFluor® 405 family has the spectral properties essentially identical to those of Alexa Fluor® 405 (Alexa Fluor® 405 is the trademark of ThermoFisher). Its fluorescence is pH-independent from pH 3 to 11. These spectral characteristics make this new dye family a great alternative to Alexa Fluor® 405.