

mFluor™ UV 460 Biotin ConjugateCatalog number: 3112
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	538.57
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	358 nm
Emission Wavelength	456 nm

Applications

mFluor™ UV 460 Biotin Conjugate contains mFluor™ UV 460 fluorophore that is well excited by the common 350 nm laser. It has a similar spectrum to BD's BV480, and demonstrates excellent avidin binding and strong blue/green fluorescence. It has avidin-binding properties similar to native biotin in terms of high affinity, fast association, and non-cooperative binding to avidin and streptavidin tetramers. mFluor™ UV 460 Biotin Conjugate can be used for detecting avidin, streptavidin or neutravidin conjugates. It is a water-soluble reagent and has pH independent fluorescence from pH 4 to pH 10. The flexible spacer between biotin moiety and fluorescent tag minimizes the steric hindrance involved in binding to avidin, streptavidin or neutravidin.