

**mFluor™ Violet 500 Biotin Conjugate**Catalog number: 3113  
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	742.82
Soluble In	DMSO

**Spectral Properties**

---

Excitation Wavelength	410 nm
Emission Wavelength	501 nm

**Applications**

---

mFluor™ Violet 500 Biotin Conjugate contains mFluor™ Violet 500 fluorophore that is well excited by the common 405 nm violet laser. It has a similar spectrum to BD Horizon™ V500 and demonstrates excellent avidin binding and strong 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™ Violet 500 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.