

# mFluor™ Violet 450-PEG4-Biotin Conjugate

Catalog Number: 3116

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	942.26
Soluble In	DMSO

## Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

## Applications

---

mFluor™ Violet 450-PEG4-Biotin Conjugate contains mFluor™ Violet 450 fluorophore that is well excited by the common 405 nm violet laser. It demonstrates excellent avidin binding and strong blue fluorescence. Its avidin-binding properties is similar to the native biotin in terms of high affinity, fast association, and non-cooperative binding to avidin and streptavidin tetramers. mFluor™ Violet 450 biotin conjugate can be used for detecting and quantifying biotin binding sites of avidin, streptavidin or neutravidin. It is a water-soluble reagent and has pH independent fluorescence from pH 4 to pH 10. The flexible PEG4 spacer between biotin moiety and fluorescent tag minimizes the steric hindrance involved in binding to avidin, streptavidin or neutravidin.