

## mFluor™ Violet 590-streptavidin conjugate

Catalog Number: 47260

Unit Size: 100 ug

### Product Details

---

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

### Chemical Properties

---

Appearance	Solid
Molecular Weight	~52 kDa
Soluble In	Water

### Spectral Properties

---

Excitation Wavelength	423 nm
Emission Wavelength	591 nm

### Applications

---

mFluor™ Violet 590-streptavidin conjugate consists of the tetrameric biotin-binding protein (streptavidin) covalently labeled with mFluor™ Violet 590 fluorescent dyes. Its high affinity for biotin conjugates is leveraged across a wide range of biochemical applications for the specific detection of various proteins, protein motifs, nucleic acids, and other molecules. This concept is central to the workflow of several detection protocols, including western blots, flow cytometry, imaging and microscopy, and microplate assays, and is crucial in purification workflows aimed at targeted fractionation. The conjugation of mFluor™ Violet 590 with streptavidin is carefully optimized to maximize fluorescence intensity while preserving the inherent binding properties of streptavidin. Upon excitation by a violet laser, mFluor™ Violet 590 exhibits a bright and photostable emission of orange fluorescence at 591 nm. mFluor™ Violet 590 streptavidin conjugates are highly effective as second-step reagents for indirect immunofluorescent staining when used in conjunction with biotinylated primary antibodies.