

**mFluor™ Violet 610-streptavidin conjugate**

Catalog Number: 47272

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 | 421 nm |
| Emission Wavelength   | 612 nm |

**Applications**

---

mFluor™ Violet 610-streptavidin conjugate consists of the tetrameric biotin-binding protein (streptavidin) covalently labeled with mFluor™ Violet 610 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 610 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 610 exhibits a bright and photostable emission of red fluorescence at 612 nm. mFluor™ Violet 610 streptavidin conjugates are highly effective as second-step reagents for indirect immunofluorescent staining when used in conjunction with biotinylated primary antibodies.