

mFluor™ Violet 545-streptavidin conjugate

Catalog Number: 47250

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	393 nm
Emission Wavelength	543 nm

Applications

The mFluor™ Violet 545-streptavidin conjugate is a tetrameric biotin-binding protein (streptavidin) covalently labeled with mFluor™ Violet 545 fluorescent dyes. It has a strong binding affinity for biotin conjugates, making it useful in various biochemical applications for detecting specific proteins, protein motifs, nucleic acids, and other molecules. This concept is essential for workflows such as western blots, flow cytometry, imaging and microscopy, and microplate assays, as well as for purification workflows aimed at targeted fractionation. The conjugation of mFluor™ Violet 545 with streptavidin is carefully optimized to maximize fluorescence intensity while preserving the inherent binding properties of streptavidin. When excited by a violet laser, mFluor™ Violet 545 emits bright and photostable green fluorescence at 543 nm. mFluor™ Violet 545 streptavidin conjugates are particularly useful as second-step reagents for indirect immunofluorescent staining when used with biotinylated primary antibodies.