

mFluor™ UV455-streptavidin conjugate

Catalog Number: 47200, 47201

Unit Size: 100 ug, 1 mg

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	357 nm
Emission Wavelength	461 nm

Applications

mFluor™ UV455 streptavidin conjugate consists of the tetrameric biotin-binding protein (streptavidin) covalently labeled with mFluor™ UV 455 fluorescent dyes. Its high affinity for biotinylated 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™ UV455 with streptavidin is carefully optimized to maximize fluorescence intensity while preserving the inherent binding properties of streptavidin. Upon excitation by an ultraviolet laser, mFluor™ UV455 exhibits a bright and photostable emission of blue fluorescence at 461 nm. mFluor™ UV455 streptavidin conjugates are highly effective as second-step reagents for indirect immunofluorescent staining when used in conjunction with biotinylated primary antibodies.