

Streptavidin-iFluor 750®-extra conjugate

Catalog Number: 46008, 46009

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	Liquid
Molecular Weight	N/A
Soluble In	PBS

Spectral Properties

Excitation Wavelength	757 nm
Emission Wavelength	779 nm

Applications

Streptavidin-iFluor 750®-extra Conjugate is a highly sensitive fluorescent reagent designed for detecting biotinylated biomolecules with near-infrared fluorescence. It features a streptavidin core for strong, specific binding to biotin, and is conjugated to iFluor® 750, a photostable fluorophore with excitation/emission maxima around 755/778 nm.

The near-infrared emission of iFluor® 750-extra minimizes background autofluorescence interference and avoids spectral overlap with visible-range dyes, making it ideal for multicolor fluorescence detection in microscopy, flow cytometry, Western blotting, and in vivo imaging.