

XFD350-streptavidin conjugate

Catalog Number: 16890

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	~52000
Soluble In	Water

Spectral Properties

Excitation Wavelength	343 nm
Emission Wavelength	441 nm

Applications

XFD350 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 350 (Alexa Fluor® is the trademark of Thermo Fisher). The conjugates of streptavidin are commonly used together with a biotin conjugate for specific detection of a variety of proteins, protein motifs, nucleic acids, and other biomolecules in western blots, flow cytometry, imaging and microscopy, and microplate assays. XFD350-streptavidin conjugate is equivalent to Alexa Fluor® 350 streptavidin, comprised of a biotin-binding streptavidin covalently attached to XFD350.