

XFD488-streptavidin conjugate

Catalog Number: 16891

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--------------------------------------------

Expiration Date	12 months upon receiving
-----------------	--------------------------

Unit Details

Units	16891 (1 mg)
-------	--------------

Reconstitution Volume	1 mL ddH ₂ O
-----------------------	-------------------------

Chemical Properties

Appearance	Solid orange
------------	--------------

Molecular Weight	N/A
------------------	-----

Soluble In	Water
------------	-------

Spectral Properties

Excitation Wavelength	499 nm
-----------------------	--------

Emission Wavelength	520 nm
---------------------	--------

Applications

XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (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. XFD488-streptavidin conjugate is equivalent to Alexa Fluor® 488 streptavidin. The conjugate is comprised of a biotin-binding streptavidin covalently attached to XFD488.