



Product Information Sheet

Product Name:	Streptavidin, HiLyte Fluor™ 555-labeled
Catalog Number:	AS-60666
Size:	1 mg
Concentration	1mg/mL
Degree of Substitution: (DOS)	See label on vial
Fluorescence:	Orange fluorescence. Excitation/Emission wavelength= 554 nm/570 nm.
Buffer:	10 mM phosphate, 150 mM NaCl, pH 7.2, with BSA, and 2 mM Sodium azide
Storage:	The dye-protein conjugate is stable for 2~3 months at 4°C. For long-term storage, divide the solution into aliquots and store at -20°C or add an equal volume of glycerol (ACS grade or higher) and store the solution at -20°C without aliquoting. Avoid multiple thaw-freeze cycles. The product is stable for 1 year at -20°C.
Instructions:	The recommended concentration for most immunofluorescent staining is 1-10 µg/mL. It is necessary to block the endogenous biotin before staining your sample with fluorophore-streptavidin conjugate. If the non-specific binding background is high, you may decrease the concentration to minimize the background. You may also centrifuge the conjugate briefly and use the supernatant only for staining.

For *in vitro* research use only