



Product Information Sheet

Product Name:	Streptavidin, 5-TAMRA Conjugated
Catalog Number:	60670
Amount:	1 mg
Concentration:	1mg/mL
Degree of Substitution: (DOS)	See label on vial
Fluorescence:	Red fluorescence. Excitation/Emission wavelength= 555 nm/580 nm
Buffer:	10 mM phosphate, 150 mM NaCl, pH 7.2, with BSA and 2 mM NaN ₃
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 endogenous biotin before staining samples with the fluorophore-streptavidin conjugate. If non-specific binding background is high, the concentration may be decreased to minimize background. The conjugate may also be centrifuged briefly using the supernatant only for staining.

For *in vitro* research use only