



Product Information Sheet

Product Name:	Streptavidin, Cross-linked Allophycocyanin Conjugated
Catalog Number:	60663
Amount:	0.2 mg
Concentration:	1 mg/mL
Degree of Substitution: (DOS)	See label on vial.
Fluorescence:	Excitation = 595, 633, 650 nm. Emission = 660 nm. Maximum absorbance is 650 nm.
Buffer:	10 mM phosphate, 150 mM NaCl, pH 7.2, with 2 mM sodium azide.
Storage:	Store conjugate at 4°C and keep away from light. DO NOT FREEZE.
Instructions:	The recommended concentration for most immunofluorescent staining is 1-10 µg/mL.

For *in vitro* research use only