



## Product Information Sheet

---

Product Name:	Streptavidin, B-Phycoerythrin Conjugated
Catalog Number:	AS-60662
Size:	0.5 mg
Concentration:	1 mg/mL
Degree of Substitution: (DOS)	See label on vial.
Fluorescence:	Excitation = 545, 563 nm. Emission = 578 nm.
Buffer:	10 mM phosphate, 150 mM NaCl, pH 7.2, with 2 mM sodium azide.
Storage:	Store the conjugate at 4°C and keep away from light. The product is stable for 6 months. DO NOT FREEZE.
Instructions:	The recommended concentration for most immunofluorescent staining is 1-10 µg/mL.

For *in vitro* research use only