



Product Information Sheet

| | |
|----------------------------------|--|
| Product Name: | Streptavidin, B-Phycoerythrin Conjugated |
| Catalog Number: | 60662 |
| Amount: | 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. DO NOT FREEZE. |
| Instructions: | The recommended concentration for most immunofluorescent staining is 1-10 µg/mL. |

For *in vitro* research use only