



Product Information Sheet

Product Name: Streptavidin, R-phycoerythrin conjugated

Catalog Number: 60669

Amount: 0.5 mg

Concentration: 1 mg/mL

Degree of Substitution: See label on vial (DOS)

Fluorescence: Excitation = 498, 539, 565 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