



## Streptavidin conjugated to B-Phycoerythrin

Product Number D4-2212  
Amount 1 mg  
Store at 2-8°C

### Form/Shipping & Storage

Supplied lyophilized. Upon receipt, store at -20°C. Reconstitute with 1 ml of ddH<sub>2</sub>O and store at 2-8°C. Phycobiliproteins are sensitive to freeze-thaw cycles.

### Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

### Buffer

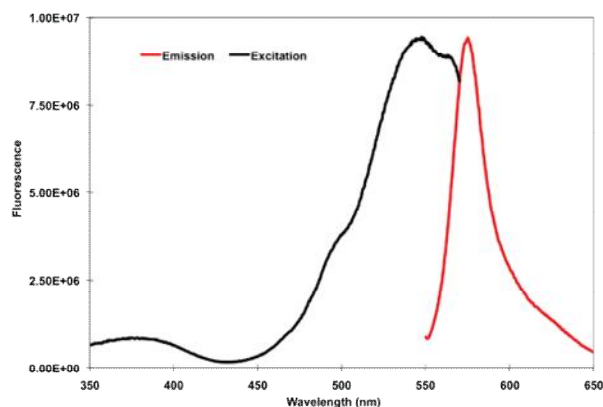
The product is in 100 mM sodium phosphate (pH 7.4) + 100 mM NaCl + 2 mM NaN<sub>3</sub> + 0.1% BSA + 50 mM Sucrose.

### Stability

Product should be stored at 2-8°C in the dark and be used within 3 months. If further dilution of the conjugate is required, use diluted material within one week.

### Note

For research use only, not for diagnostic or therapeutic use.



Fluorescence excitation and emission spectra of B-phycoerythrin in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. Emission scan was made with excitation at 498nm. Excitation scan was taken with emission at 573 nm. Scans were normalized to equalize peak heights.

### Spectral Characteristics

Visible absorption maxima	545
Emission maximum	572

### Concentrations

Streptavidin	10.07 µM
B-Phycoerythrin	2.52 µM