



Streptavidin conjugated to R-Phycoerythrin

Product Number D5-2212
Amount 1 mg
Store at 2-8°C

Form/Shipping & Storage

Upon receipt, store at 2-8 °C. DO NOT FREEZE.
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

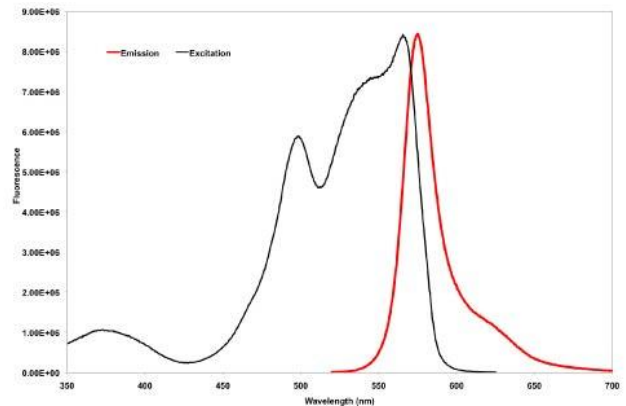
The product is in 100 mM sodium phosphate (pH 7.4) + 100 mM NaCl + 2mM NaN₃.

Stability

Product should be stored at 2-8°C in the dark and be used within 6 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 R-phycoerythrin in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. Emission scan was taken with excitation at 498 nm. Excitation scan was taken with emission at 575 nm. Scans were normalized to equalize peak heights.

Spectral Characteristics

Visible absorption maxima 565>540>498
Emission maximum 572

Fluor:Protein = ~1:1



COLUMBIA BIOSCIENCES