



## Mouse IgG<sub>1</sub> conjugated to SureLight® Allophycocyanin (APC) (Isotype control)

Product Number D3-1601  
Amount 100 µg  
Store at 2-8°C

### Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 2-8°C in the dark, do not freeze. Phycobiliproteins are sensitive to freeze-thaw cycles.

### Handling

Reconstitute to 1.0 ml with distilled deionized water vortex gently and allow vial to sit on ice for 20 minutes. We recommend that the investigator determine the appropriate working concentration for their specific application.

Avoid exposure to heat and light.

### Buffer

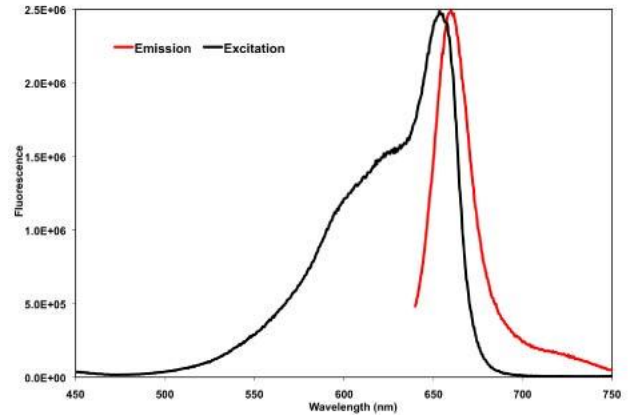
Upon rehydration with 1.0 ml distilled deionized water; the product is in 10 mM Tris (pH 8.2), 50 mM sucrose, 150 mM sodium chloride, 0.1% BSA as a stabilizer, 1µg/ml pentachlorophenol and 50mm sucrose. The concentration of the conjugate is 100 µg/ml

### Stability

Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within one year. If further dilution of the conjugate is required, use diluted material within one month.

### Note

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



Fluorescence excitation and emission spectra of APC in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. The emission scan was taken with excitation at 630 nm. The excitation scan was taken with an emission at 660 nm. The curves were normalized to equalize peak heights.

### Spectral Characteristics

Visible absorption maxima	650 nm
Emission maximum	660 nm

Fluor:Protein	~1:1
---------------	------



**COLUMBIA BIOSCIENCES**