



## Goat anti-Human IgG conjugated to PerCP

Product Number D11-110-1MG  
Amount 1mg  
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. PerCP is 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.

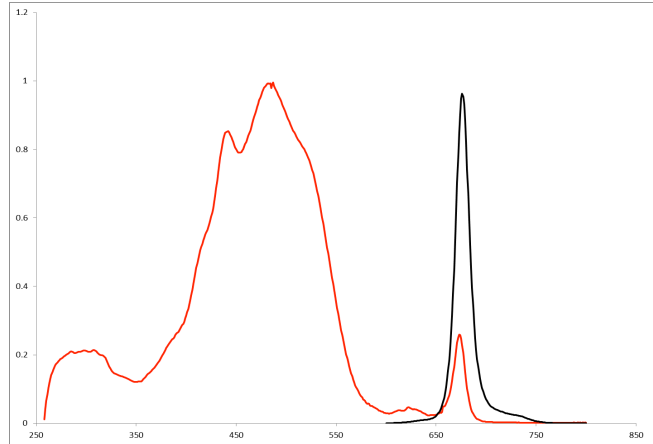
### Buffer

Upon rehydration with 1.0 ml distilled deionized water; the product is in 100 mM Sodium Phosphate (pH 7.4), 100 mM NaCl, 50 mM Sucrose, 0.1% BSA and 2 mM sodium azide. The concentration of the conjugate is 1mg/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 6 months. If further dilution of the

conjugate is required, use diluted material within



one month.

Absorbance and emission spectra of Peridinin- Chlorophyll Protein Complex (PerCP) in 10 mM Sodium Phosphate (pH 7.4). Emission scan was taken with an excitation at 488nm.

### Note

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

### Spectral Characteristics

Visible absorption maxima	483 nm
Emission maximum	673 nm



**COLUMBIA BIOSCIENCES**