



## Goat anti-Llama IgG conjugated to R-Phycoerythrin

Product Number D5-111  
Amount 100 ug  
Lot # RPE019-12-002  
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

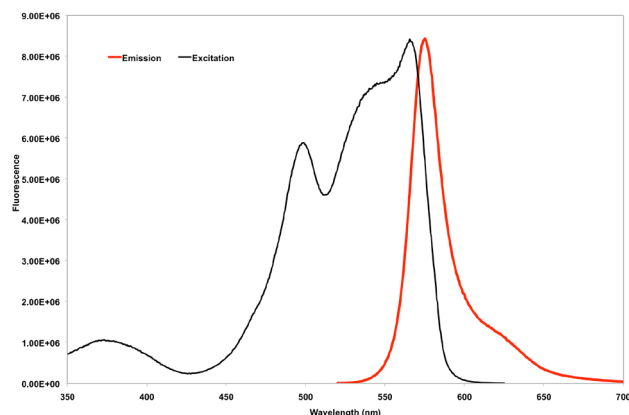
Upon rehydration with 1.0 ml distilled deionized water; the product is in 100 mM sodium phosphate (pH 7.4), 50 mM sucrose, 150 mM sodium chloride, 0.1% BSA as a stabilizer, and 0.05% sodium azide as a preservative. The concentration of the conjugate is 100 µg/ml.

### 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 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.



### Spectral Characteristics

Visible absorption maxima 565>540>498  
Emission maximum 578

### Concentration

After reconstitution to 1.0 ml

Fluor:Protein = ~1.2:1

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com