



## Rabbit anti-T7 tag IgG conjugated to R-Phycoerythrin

Product Number D5-1832  
Lot# XXXX  
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 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

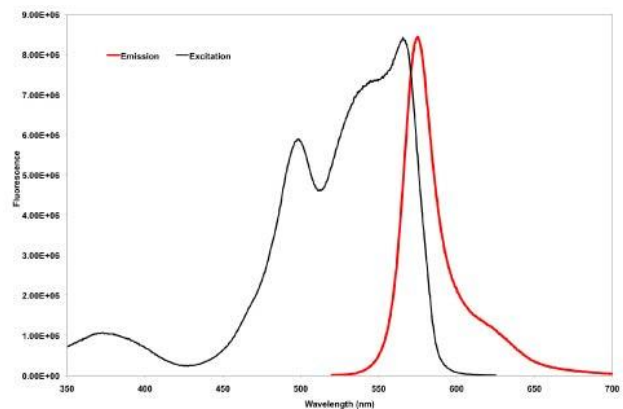
Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 12 months. If further dilution of the conjugate is required, use diluted material within one week.

### Antigen Info

Antibody was raised against T7-tag with amino acid sequence MASMTGGQMG.

### 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 578

### Concentration

After reconstitution to 1.0 ml is 0.1mg/mL

### References

Itoh T, Fujita N, Kanno E, Yamamoto A, Yoshimori T, Fukuda M. Golgi-resident Small GTPase Rab33B Interacts with Atg16L and Modulates Autophagosome Formation. *Mol Biol Cell*. 2008 Jul; 19(7): 2916-2925.

Day B, Dahlbeck D, Staskawicz BJ. NDR1 Interaction with RIN4 Mediates the Differential Activation of Multiple Disease Resistance Pathways in Arabidopsis. *Plant Cell*. 2006 Oct; 18(10): 2782-2791.

Rayala SK, Martin E, Sharina IG, Molli PR, Wang X, Jacobson R, Murad F, Kumar R. Dynamic interplay between nitration and phosphorylation of tubulin cofactor B in the control of microtubule dynamics. *PNAS*. 2007 Dec 4; 104(49): 19470-19475

4985 Winchester Blvd. Frederick, MD 21703  
301.732.5415  
www.columbiabiosciences.com  
For technical inquiries: info@columbiabiosciences.com  
For sales inquiries: sales@columbiabiosciences.com