



## Mouse anti-6 Histidine IgG conjugated to PerCP

Product Number D11-1711  
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. 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), 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

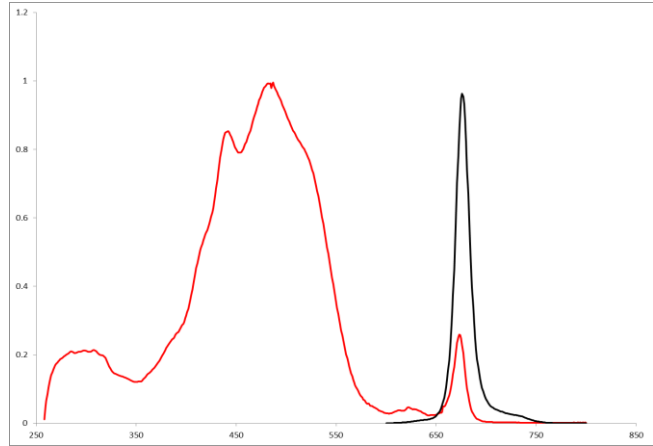
Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 6 months. If further dilution of the conjugate is required, use diluted material within one month.

### References

Tang YQ, Han SY, Zheng H, Wu L, Ueda M, Wang XN, Lin Y. Construction of cell surface-engineered yeasts displaying antigen to detect antibodies by immunofluorescence and yeast-ELISA. *Appl Microbiol Biotechnol*. 2008 Jul;79(6):1019-26.

Kojima T, Takei Y, Ohtsuka M, Kawarasaki Y, Yamane T, Nakano H. PCR amplification from single DNA molecules on magnetic beads in emulsion: application for high-throughput screening of transcription factor targets. *Nucleic Acids Res*. 2005 Oct 6;33(17):e150.

Tran TM, Moreno A, Yazdani SS, Chitnis CE, Barnwell JW, Galinski MR. Detection of a *Plasmodium vivax* erythrocyte binding protein by flow cytometry. *Cytometry A*. 2005;63(1):59-66.



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**