



Mouse anti-HRP IgG2b (clone 3P/2H11) conjugated to R-Phycoerythrin

Product Number D5-1901
Lot number RPE058-16-001
Amount 100 µg total protein
Store at 4°C

Form/ Storage

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

Handling

Avoid exposure to heat and light. Prior to use reconstitute to 1 ml with distilled deionized water, vortex and allow it to sit on ice for 20 minutes.

Buffer

Upon reconstitution, the product is in 100 mM sodium phosphate (pH 7.4), 50 mM sucrose, 100 mM sodium chloride, 0.1% BSA as a stabilizer, and 2 mM sodium azide as a preservative.

Stability

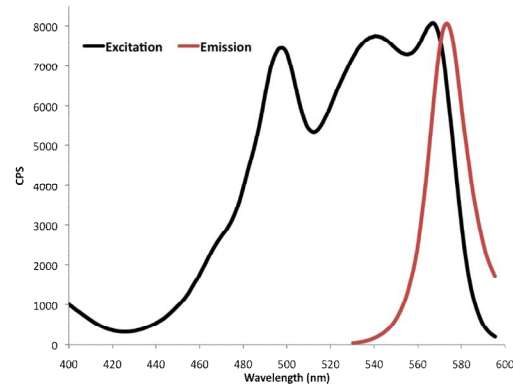
Lyophilized material is stable for one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 3 months.

Antigen Info

Peroxidase (Horseradish) (HRP)

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.

Spectral Characteristics

| | |
|---------------------------|-------------|
| Visible absorption maxima | 565>540>498 |
| Emission maximum | 578 nm |

Concentration

After reconstitution to 1.0 ml
0.1 mg/mL

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com
For sales inquiries: sales@columbiabiosciences.com