



Mouse anti-HA (haemagglutinin) tag (YPYDVPDYA) IgG conjugated to R-Phycoerythrin

Product Number D5-1722
Amount 100 µg
Clone 16B12
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

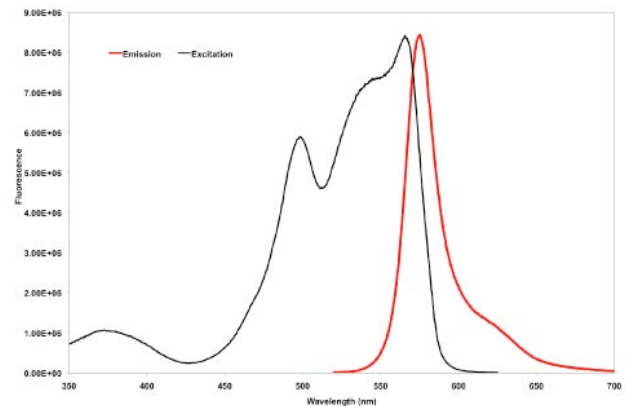
Upon rehydration with 1.0 ml distilled deionized water; the product is in 100 mM sodium phosphate (pH 7.4), 50 mM sucrose, 150 mM sodium chloride, 0.1% BSA as a stabilizer, and 0.05% sodium azide as a preservative. The concentration of the conjugate is 100 µg/ml

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.



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

Fluor:Protein = ~1.3:1



COLUMBIA BIOSCIENCES