



Mouse anti-c-myc IgG conjugated to R-Phycocerythrin

Product Number D5-1714
Lot Number XXXX
Amount 100 µg
Store at 2-8°C
Clone 9E10 (sequence EQKLISEEDL)
IgG Class mouse IgG_{1,κ}

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

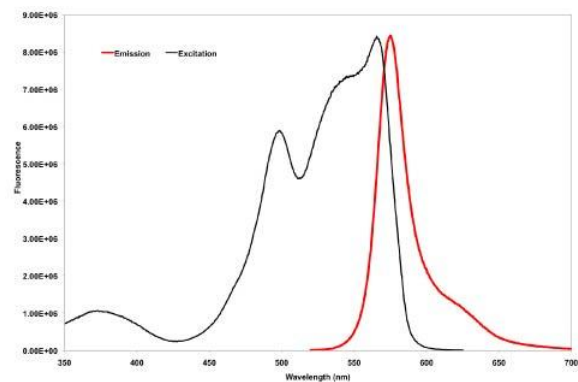
Upon rehydration with 1.0 ml distilled deionized water; the product is in 100 mM sodium phosphate (pH 7.4), 50 mM sucrose, 100 mM sodium chloride, 0.1% BSA, and 2 mM sodium azide. The concentration of the conjugate is 100 µg/ml

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: 100 µg/ml

References:

Gazitt Y, He YJ, Erdos GW, Chang L, Ashktorab H, Cohen RJ. Development of a two color immunofluorescence stain and immunolocalization method for N-myc and c-myc oncoproteins with a newly generated mouse IgM anti N-myc antibody. J Immunol Methods. 1992 Apr 8;148(1-2):159-69.

Kieke MC, Cho BK, Boder ET, Kranz DM, Wittrup KD. Isolation of anti-T cell receptor scFv mutants by yeast surface display. Protein Eng. 1997 Nov;10(11):1303-10.

Lincoln ST, Bauer KD. Limitations in the measurement of c-myc oncoprotein and other nuclear antigens by flow cytometry. Cytometry. 1989 Jul;10(4):456-62.

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com