



Mouse anti-c-Myc IgG conjugated to SureLight™ P-3

Product Number D7-1714
Amount 100 µg
Lot# G022AC01
Clone 9E10 (sequence EQKLISEEDL)
Store at 4°C

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 4°C in the dark. Phycobiliproteins are sensitive to freeze-thaw cycles; after reconstitution, store at 2-8°C in the dark, do not freeze. The concentration after resuspension is 50 µg/mL

Handling

Prior to use, reconstitute to 200 µl with distilled deionized water, vortex gently and allow vial to sit on ice for 20 minutes. Avoid exposure to heat and light.

Buffer

Upon rehydration with 200 µl distilled deionized water, the product is in 100 mM sodium phosphate (pH 7.4), 200 mM sucrose, 150 mM sodium chloride, 0.1% BSA as a stabilizer, and 2 mM sodium azide as a preservative.

Stability

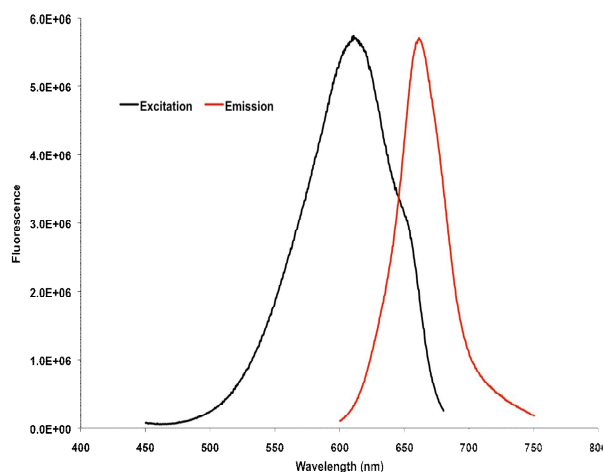
Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

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.

Kieckhefer 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.



Note

For research use only, not for diagnostic or therapeutic use.

Spectral Characteristics

Visible absorption maxima 614>650 (sh) nm
Emission maximum 662 nm

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com