

## Mouse Anti-c-myc IgG conjugated to PerCP

Product Number D11-1714  
Amount 0.1 mg total protein  
Clone 9E10 (sequence EQKLISEEDL)  
Store at 2-8°C

### Form/Shipping & Storage

Supplied lyophilized. Upon receipt, store at -20 °C. Reconstitute with 1 ml of ddH<sub>2</sub>O and store at 2-8 °C. Phycobiliproteins are sensitive to freeze-thaw cycles. Concentration after resuspension is 01.mg/mL

### Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

### Buffer

Upon reconstitution, the product is in 10 mM tris (pH 8.2) + 150 mM NaCl + 50 mM Sucrose, 0.01% NaN<sub>3</sub> + 0.1% BSA.

### Stability

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.

### Note

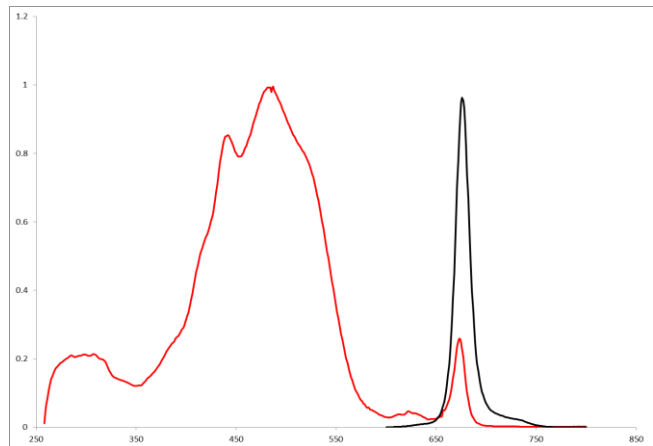
For research use only, not for diagnostic or therapeutic use.

### 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.



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.

### Spectral Characteristics

Visible absorption maxima	483
Emission maximum	673



**COLUMBIA BIOSCIENCES**