



Anti-c-myc IgG conjugated to Dylight® 650

Product Number	D10-1714
Amount	0.1 mg total protein
Clone	9E10 (sequence EQKLISEEDL)
Lot Number	XXXX
Approximately	5-7
Conc	1.0 mg/ml
F:P Ratio	2.08
Store at	2-8°C

Form/Shipping & Storage

Supplied liquid. Upon receipt, store at 2-8 °C. Fluorescent dyes are sensitive to light. Please store in the dark and avoid exposure to sunlight.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application.

Buffer

0.1M sodium phosphate, 0.1M NaCl, pH7.4, 2 mM NaN₃

Stability

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.

Note

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

Spectral Characteristic

Visible absorption maxima	655 nm
Emission maximum	672 nm

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