



Anti-c-myc conjugated to SureLight™ APC

Product Number D3-1714
Amount 100 µg 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 ddH₂O and store at 2-8 °C. Phycobiliproteins are sensitive to freeze-thaw cycles.

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, 1 µg/mL pentachlorophenol + 0.1% BSA.

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

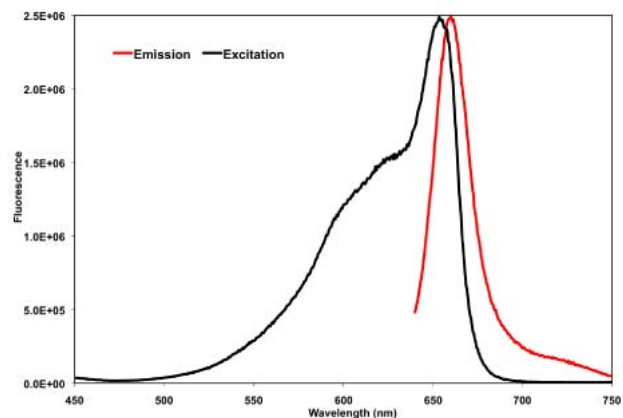
For research use only, not for diagnostic or therapeutic use.

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.



Fluorescence excitation and emission spectra of APC in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. The emission scan was taken with excitation at 630 nm. The excitation scan was taken with an emission at 660 nm. The curves were normalized to equalize peak heights.

Spectral Characteristics

Visible absorption maxima 652
Emission maximum 657

Surelight APC 0.441 µM
Anti-c-myc 0.407 µM

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com



Anti-c-myc conjugated to SureLight™ APC

Product Number D3-1714
Amount 1.0 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 ddH₂O and store at 2-8°C. Phycobiliproteins are sensitive to freeze-thaw cycles.

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, 1 µg/mL pentachlorophenol + 0.1% BSA.

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

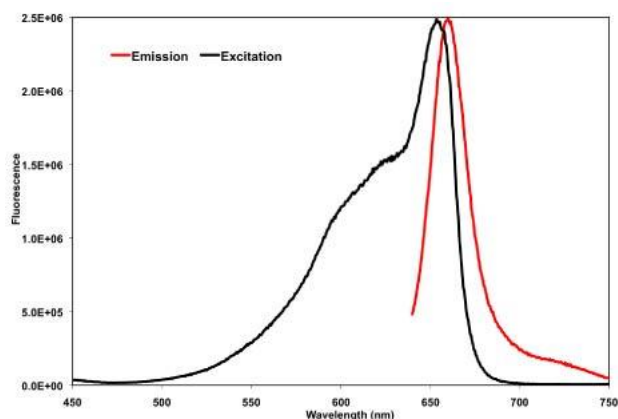
For research use only, not for diagnostic or therapeutic use.

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.



Fluorescence excitation and emission spectra of APC in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. The emission scan was taken with excitation at 630 nm. The excitation scan was taken with an emission at 660 nm. The curves were normalized to equalize peak heights.

Spectral Characteristics

Visible absorption maxima 652
Emission maximum 657

Surelight APC 4.81 µM
Anti-c-myc 4.21 µM
Fluor Protein -1.14:1

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com