



## **Mouse anti-c-Myc IgG conjugated to Biotin**

Product Number    B1-1714  
Amount             100 µg  
Store at             4°C

### Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at -20° Celsius.

### Handling

Prior to use, reconstitute to 1 ml with distilled water, vortex gently and allow material to reconstitute for 20 minutes. After reconstitution store the product at 4°C.

### Buffer

Upon rehydration with 1 ml distilled water the product is in PBS (pH 7.4) and 50 mM sucrose.

### Stability

Lyophilized material is stable for up to one year.

### Western blot detection

Recommended dilution of 1:500 - 1:10000 depending on target protein concentration.

### Antibody/Antigen Information

Mouse IgG<sub>1</sub> clone 9E10 against sequence EQKLISEEDL.

### 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.



**COLUMBIA BIOSCIENCES**