

c-Myc antibody (mAb)

Catalog Nos: 61075, 61076

RRID: AB_2793497

Clone: 6A10

Isotype: IgG2a

Application(s): EMSA, IP, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 68 kDa

Background: **c-Myc** (Myc) is a sequence specific DNA binding transcription factor that regulates gene expression, especially genes involved in cell division. **Myc** family members contain a basic Helix-Loop-Helix leucine zipper (bHLHZ) domain. Myc family proteins associated with thousands of genes, especially those involved in cell growth and proliferation. **c-Myc** serves as an integrator of signal transduction, linking extracellular signals to an acute cellular growth response. Mutations and overexpression of c-Myc are associated with cell proliferation and cancer.

Immunogen: This c-Myc antibody was raised against a recombinant protein corresponding to amino acids 1-262 of human c-Myc protein.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

This antibody is also available as an AbFlex[®] engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91213.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.