

## c-Myc antibody (pAb) (65 kDa form)

**Catalog No:** 39012

**RRID:** AB\_2793149

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human, Monkey, Mouse

**Volume:** 100  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Concentration:** 0.337  $\mu$ g/ $\mu$ l

**Molecular Weight:** 65 kDa

**Background:** c-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. 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 synthetic peptide corresponding to amino acid residues 9-22 of human c-Myc; it recognizes a 65 kDa form of c-Myc.

**Buffer:** PBS containing 0.2% gelatin, 0.01% thimersol and 0.1% sodium azide. Thimersol and sodium azide are highly toxic.

### Application Notes:

Validated Applications:

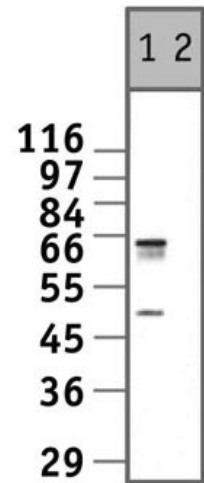
WB: 1:500 - 1:1,000 dilution

Published Applications: see references

For optimal results, primary antibody incubations should be performed overnight at 37°C. Individual optimization may be required.

**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.



### c-Myc antibody (65 kDa form) tested by Western blot.

Detection of c-Myc by Western blot analysis. A 65 kDa form of the transcription factor c-Myc is detected in nuclear extracts derived from 3T6 Swiss albino cells using c-Myc rabbit polyclonal antibody at a 1:1,000 dilution (lane 1). The signal is eliminated when the c-Myc immunizing peptide is added to the reaction mixture (lane 2).