

Recombinant c-Myc protein

Catalog No: 31117

Expressed In: *E. coli*

Quantity: 5 µg

Concentration: 0.12 µg/µl

Source: Human

Buffer Contents: 5 µg of Recombinant c-Myc protein in Dilution Buffer AM1 (20 mM Tris-Cl (pH 7.5), 20% glycerol, 100 mM KCl, 1 mM DTT and 0.2 mM EDTA).

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.

Protein Details: Recombinant c-Myc was expressed in a bacterial system (accession number NM 002467) with an N-terminal His-Tag and purified by an affinity column in combination with FPLC chromatography. The purified recombinant protein is greater than 75% homogeneous and contains no detectable protease, DNase and RNase activity.

Application Notes: Recombinant c-Myc protein is suitable for DNA and protein-protein interaction assays. 100 ng are sufficient for DNA-protein and protein-protein interaction studies. The standard curve for TransAM® c-Myc was generated using the range of 5 -0.625 ng of protein. The molecular weight of the protein is ~50 kDa.

NOTE: The presence of Poly [d(I-C)] in buffers may affect protein functionality and should be avoided.

References:

This product was used in the following publications:

Piccinni, E., *et. al.* (2011). "Direct interaction of Gas41 and Myc encoded by amplified genes in nervous system tumours." *Acta Biochim. Pol.* 58(4):529-34. PMID: 22068108.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

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