

c-myc

Proto-oncogene

human, recombinant, Sf9 insect cells

Cat. No.	Amount
PR-773	5 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 20 mM Tris-HCl pH 8.0, 100 mM KCl, 0.2 mM EDTA, 1 mM DTT, 20% glycerol.

Activity

100 ng are sufficient for a protein-protein interaction assay.

Application

c-myc has been applied in DNA and protein-protein interaction assays.

Molecular Weight

64 kDa

Purity

>95% by SDS-PAGE

Description

The universal deregulation of c-myc gene expression in tumor cells suggests that this oncogene represents an attractive target for cancer therapeutic purposes. The c-myc promoter integrates diverse mitogenic signalling cascades, which are constitutively activated in tumor cells, and translates them into expression of the c-MYC transcription factor, which promotes cell proliferation, growth, differentiation, and apoptosis by regulating the expression of numerous target genes. The structural and biochemical features of the MYC family (MYC, N-MYC, and L-MYC) mark them as direct regulators of gene expression. As basic helix-loop-helix leucine zipper proteins (bHLH-ZIP), the MYCs acquire the capacity to bind the DNA sequence CACGTG (E-box) when dimerized with MAX (another bHLH-ZIP). A head-to-tail pair of MYC-MAX dimers may, in turn, form a heterotetramer capable of bridging distant E-boxes. Among the broadly distributed positive enforcers of MYC action that are often recruited to target genes are chromatin remodeling (SWI/SNF relatives) and modifying complexes (TRAPP/GCN5 and relatives); these complexes mobilize nucleosomes and acetylate histones and/or other targets to activate gene expression. MYC binds TBP along an auxiliary pathway to control gene expression. MAD and MNT generally oppose MYC action by enlisting histone deacetylase complexes. Besides acting at the level of chromatin, MYC may also operate at later stages of the transcription cycle, after pre-initiation complex formation. In addition to using generic chromatin complexes to up- or down-regulate transcription, the MYC network also conscripts individual factors to modify expression locally on an ad hoc basis. For example, YY1, AP2, MIZ1, SP1, BRCA1, and other proteins interact directly with MYC, and so may directly modify the output of the MYC network.

Recombinant c-myc was expressed in a baculovirus system and purified by an affinity column in combination with FPLC chromatography.

The purified recombinant protein is greater than 90% homogeneous and contains no detectable protease, DNase, and RNase activity.

Selected References:

O'Donnell *et al.* (2005) c-Myc-regulated microRNAs modulate E2F1 expression. *Nature*. **435**:839.

Dang *et al.* (1999) Function of the c-Myc oncogenic transcription factor.

c-myc

Proto-oncogene

human, recombinant, Sf9 insect cells

Exp. Cell Res. **253**:63.

Grandori *et al.* (2000) The Myc/Max/Mad network and the transcriptional control of cell behavior. *Annu. Rev. Cell Dev. Biol.* **16**:653.

Eisenman (2001) Deconstructing myc. *Genes & Dev.* **15**:2023.

McMahon *et al.* (2000) The essential cofactor TRRAP recruits the histone acetyltransferase hGCN5 to c-Myc. *Mol. Cell. Biol.* **20**:556.

Bouchard *et al.* (2001) Regulation of cyclin D2 gene expression by the Myc/Max/Mad network: Myc-dependent TRRAP recruitment and histone acetylation at the cyclin D2 promoter. *Genes & Dev.* **15**:2042.

Nair *et al.* (2003) X-ray structures of Myc-Max and Mad-Max recognizing DNA. Molecular bases of regulation by proto-oncogenic transcription factors. *Cell* **112**:193.

Cheng *et al.* (1999) c-MYC interacts with INI1/hSNF5 and requires the SWI/SNF complex for transactivation function. *Nat. Genet.* **22**:102.

Luscher (2001) Function and regulation of the transcription factors of the Myc/Max/Mad network. *Gene* **277**:1.

Eberhardy *et al.* (2001) c-Myc mediates activation of the cad promoter via a post-RNA polymerase II recruitment mechanism. *J. Biol. Chem.* **276**:48562.

Shrivastava *et al.* (1996) YY1 and c-Myc associate *in vivo* in a manner that depends on c-Myc levels. *Proc. Natl. Acad. Sci.* **93**:10638.

Sakamuro *et al.* (1999) New Myc-interacting proteins: a second Myc network emerges. *Oncogene* **18**:2942.