

## Recombinant c-Myc / MAX complex

---

**Catalog No:** 81087, 81787

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.4 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant c-Myc / MAX Complex is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 20% glycerol, 0.04% Triton X-100 and 0.5 mM TCEP.

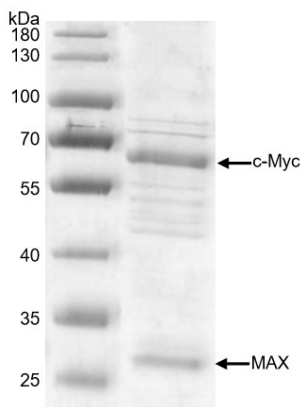
**Background:** c-Myc (MYC Proto-Oncogene, BHLH Transcription Factor), also known as MYCC or MRTL, is a transcription factor that binds DNA in a non-specific manner, yet also specifically recognizes the core sequence 5-CAC[GA]TG-3. Diseases associated with MYC include Burkitt Lymphoma and Precursor T-Cell Acute Lymphoblastic Leukemia. Among its related pathways are Common Cytokine Receptor Gamma-Chain Family Signaling Pathways and IL-2 Pathway. c-Myc can activate the transcription of growth-related genes. MAX (MYC Associated Factor X), also known as BHLHd4, is a member of transcription regulators. MAX forms a sequence-specific DNA-binding protein complex with MYC or MAD, and the MYC / MAX complex is a transcriptional activator, whereas the MAD / MAX complex is a repressor. It may repress transcription via the recruitment of a chromatin remodeling complex containing H3 Lys-9 histone methyltransferase activity. c-Myc / MAX complex can specifically recognize the core sequence 5-CAC[GA]TG-3 and bind to the E box DNA consensus sequence and regulates the transcription of specific target genes.

**Protein Details:** Recombinant c-Myc / MAX Complex that includes human c-Myc amino acids 2 - 454 (accession number NP\_002458.2) with an N-terminal FLAG tag and full length human MAX protein (accession number NP\_660087.1) with an N-terminal 6xHis Tag was expressed in Sf9 cells. The molecular weights of c-Myc (2 - 454) and MAX are 51.8 kDa and 20.6 kDa, respectively.

**Application Notes:** This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

**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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

### c-Myc / MAX Complex



### Recombinant c-Myc / MAX Complex

10% SDS-PAGE Coomassie staining

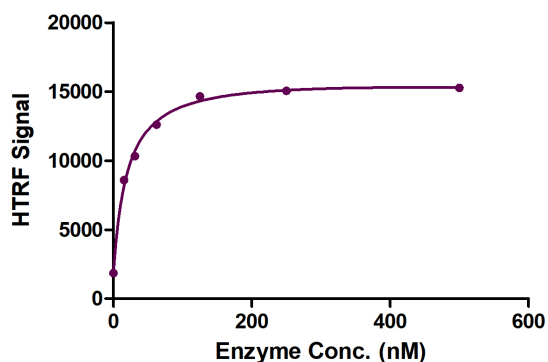
Molecular Weights:

c-Myc: 51.8 kDa

MAX: 20.6 kDa

Purity: >85%

### c-Myc / MAX Complex Titration



### Recombinant c-Myc / MAX Complex

1  $\mu$ M substrate DNA (5'-Biotin GGGACCACGTGGTCCC-3') were incubated with different concentrations of c-Myc / MAX Complex protein in HTRF binding buffer contain 50mM HEPES-NaOH pH 7.4, 0.1% BSA for 1 hour, then 10  $\mu$ l detection reagents contain anti-His antibody (1:100) and SA-XL665 (1:100) diluted with 1 $\times$  Detection Buffer were added and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.