## Recombinant MAX protein



## Catalog No: 31244 Expressed In: *E. coli*

Quantity: 10 µg Concentration: 0.15 µg/µl Source: Human

**Buffer Contents:** 10 µg of recombinant protein supplied at a concentration of 0.15 µg/µl in 20 mM Tris-HCI, pH 7.3, 300 mM KCI, 0.2 mM EDTA and 20% glycerol.

**Background:** Max protein, isoform b (also known as Myc associated factor X or MGC10775) belongs to a basic helix-loop-helix leucine zipper (bHLHZ) family of transcription factors. Other family members include Mad, Mxi1, and Myc, which bind the E box DNA site 5'-CAC(GA)TG-3'. c-Myc and Mad form heterodimers with Max to regulate transcription. Mad-Max heterodimers are associated with transcriptional activation, while Myc-Max heterodimers are associated with transcriptional activation. This variant lacks an in-frame segment, resulting in a shorter protein known as isoform b.

**Protein Details:** Recombinant MAX protein was expressed in *E. coli* as the full-length human protein (accession number NP\_660087) with an N-terminal FLAG tag. The molecular weight of the protein is ~18 kDa.

**Application Notes:** Recombinant MAX protein is suitable for use in *in vitro* transcription assays. We recommended to start with a range of protein concentrations (*e.g.* 5 ng, 50 ng, and 500 ng) to determine the optimal amount needed for each *in vitro* transcription assay.

**Storage and Guarantee:** 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.