

Recombinant MAX protein, His-Tag

Catalog No: 81026, 81726

Lot No: 01318001

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1.5 µg/µl

Source: Human

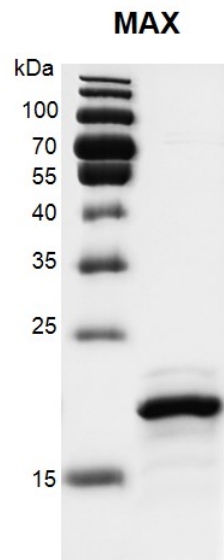
Buffer Contents: Recombinant MAX protein, His-tag is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol and 0.5 mM TCEP.

Background: MAX (MYC Associated Factor X, also known as BHLHd4) is a member of the basic helix-loop-helix leucine zipper (bHLHZ) family of transcription factors. It is able to form homodimers and heterodimers with other family members, which include Mad, Mxi1 and Myc. MAX forms a sequence-specific DNA-binding protein complex with MYC or MAD which recognizes the core sequence 5-CAC[GA]TG-3. The MYC:MAX complex is a transcriptional activator, whereas the MAD:MAX complex is a repressor. May repress transcription via the recruitment of a chromatin remodeling complex containing H3 Lys-9 histone methyltransferase activity.

Protein Details: Recombinant MAX protein, His-tag was expressed in *E. coli* cells as the full length protein (accession number NP_660087.1) with an N-terminal 6×His tag. The molecular weight of this protein is 20.7 kDa.

Application Notes: Recombinant MAX protein, His-tag is suitable for use in *in vitro* transcription assays and protein interaction.

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.



Recombinant MAX protein, His-tag, protein gel

12.5% SDS-PAGE Coomassie staining

MW: 20.7 kDa

Purity: > 85%