

Recombinant MAX protein

Catalog No: 81017, 81717

Lot No: 20817001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.16 µg/µl

Source: Human

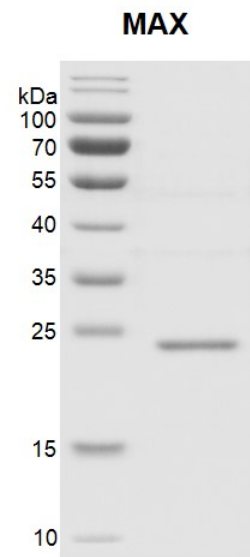
Buffer Contents: Recombinant MAX protein is supplied at a concentration of 0.16 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

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 was expressed in a baculovirus expression system as the full length protein (accession number NP_660087.1) with an N-terminal FLAG tag. The molecular weight of MAX protein is 18.5 kDa.

Application Notes: Recombinant MAX protein 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 gel
12.5% SDS-PAGE gel stained with Coomassie blue.
MW: 18.5 kDa
Purity: > 90%