

Recombinant MAD1 protein

Catalog No: 31329

Source: Human

Quantity: 10 µg

Expressed In: *E. coli*

Buffer: The protein is supplied at a concentration of 0.1 µg/µl in 20 mM Tris-HCl (pH7.3), 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Contents: 10 µg recombinant MAD1 protein.

Protein Details:

MAD1 (also known as MXD1, MAX dimerization protein 1) is a member of the MAD family of proteins. MAD and MAX proteins form heterodimers that bind to the DNA sequence 5'-CAC(GA)TG-3' to repress transcription, while MAX and MYC proteins form heterodimers that activate transcription.

Recombinant MAD1 protein was expressed in *E. coli* as the full-length human protein (accession number AAA36194) with an N-terminal FLAG tag. The molecular weight of the protein is ~27 kDa.

Dilution of Protein:

Recombinant MAD1 is supplied in solution and should be kept as a solution in order to maintain full activity.

Application:

Recombinant MAD1 protein is suitable for use in *in vitro* transcription assays. We recommend 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:

Recombinant proteins in solution are temperature sensitive and should be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage.

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.