

Recombinant USF1 protein

Catalog No: 31333

Source: Human

Quantity: 10 µg

Expressed In: *E. coli*

Buffer: The protein is supplied at a concentration of 0.4 µg/µl in 20 mM Tris-HCl (pH7.3), 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Contents: 10 µg recombinant USF1 protein.

Protein Details:

USF1 (Upstream transcription factor 1) is a member of the basic helix-loop-helix leucine zipper family, and can function as a cellular transcription factor that binds to symmetrical DNA sequence (E-boxes) (5'-CACGTG-3') that are found in a variety of viral and cellular promoters. The USF1 gene has been linked to familial combined hyperlipidemia (FCHL). Two transcript variants encoding distinct isoforms have been identified for this gene.

Dilution of Protein:

Recombinant AP-2α is supplied in solution and should be kept as a solution in order to maintain full activity.

Application:

Recombinant USF1 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.