

Recombinant TFCP2 protein

Catalog No: 31250

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 0.1 µg/µl

Source: Human

Buffer Contents: 10 µg of recombinant protein supplied at a concentration of 0.1 µg/µl in 20 mM Tris-HCl, pH 7.3, 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Background: TFCP2 (also known as transcription factor CP2, alpha globin, transcription factor LSF, TFCP2C, SAA3 enhancer factor or LBP-1c, isoform 1) is a component of the stage selector protein (SSP) complex, which consists of TFCP2 and two copies of NFE4. It aids in the interaction of gamma-globin genes with enhancer elements in fetal erythroid cells. TFCP2 has also been shown to interact with APBB1 and RNF2.

Protein Details: Recombinant TFCP2 protein was expressed in *E. coli* as the full-length human protein (accession number NP_005644) with an N-terminal polyhistidine and FLAG tag. The molecular weight of the protein is ~57 kDa.

Application Notes: Recombinant TFCP2 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: Lyophilized proteins can be stored at -20°C or -80°C, preferably desiccated. 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.