

Recombinant AP-2α protein

Catalog No: 31331

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 AP-2α protein.

Protein Details:

AP-2α transcription factors activate transcription and regulate gene expression during embryogenesis. There are four different isoforms of AP-2 (AP-2α, AP-2β, AP-2γ and AP-2δ). The family members share a common structure in which the N-terminal region is involved with transcriptional activation and the C-terminal region is involved in DNA binding.

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 AP-2α 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.