

Recombinant RAP74 protein

Catalog No: 31315

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 0.3 µg/µl

Source: Human

Buffer Contents: 10 µg recombinant RAP74 protein supplied at a concentration of 0.3 µg/µl in 20 mM Tris-HCl, pH 7.3, 500 mM KCl, 0.2 mM EDTA and 20% glycerol.

Background: **RAP74** (General Transcription Factor IIF, Polypeptide 1, 74kDa) TFIIF is a general transcription initiation factor that binds to RNA polymerase II and helps to recruit it to the initiation complex in collaboration with TFIIB. It promotes transcription elongation.

Protein Details: Full length recombinant RAP74 was expressed from (accession number NP_002087.2) human protein with a C-terminal polyhistidine tag.

Application Notes: Recombinant RAP74 is suitable for *in vitro* transcription, protein-protein interactions and protein-DNA interaction studies. A recommended starting point for *in vitro* transcription assays is 20 ng of recombinant protein per reaction. The molecular weight of the protein is ~74 kDa.

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 guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.