

Recombinant PC4 protein

Catalog No: 31317

Source: Human

Quantity: 10 µg

Expressed In: *E. coli*

Buffer: The recombinant protein is supplied at a concentration of 1 µg/µl in 20 mM Tris-HCl, pH 7.3, 100 mM KCl, 0.2 mM EDTA and 20% glycerol.

Contents: 10 µg of recombinant PC4 protein

Protein Details:

PC4, activated RNA polymerase II transcription cofactor 4, is a non-specific DNA/chromatin binding factor implicated in transcription and DNA replication. Recombinant PC4 was expressed from full-length (accession number NP_006704.3) human protein.

Dilution of Protein:

Recombinant PC4 is supplied in solution and should be kept as a solution in order to maintain full activity.

Application:

Recombinant PC4 is suitable for *in vitro* transcription, protein-protein interactions and protein-DNA interaction studies. A recommended starting point for *in vitro* transcription assays is 100-200 ng recombinant protein per reaction. The molecular weight of the protein is ~15 kDa.

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.