

## Recombinant RAP30 protein

---

**Catalog No:** 31316

**Source:** Human

**Quantity:** 10 µg

**Expressed In:** E. coli

**Buffer:** The recombinant protein is supplied at a concentration of 0.4 µg/µl in 20 mM Tris-HCl, pH 7.3, 100 mM KCl, 0.2 mM EDTA and 20% glycerol.

**Contents:** 10 µg recombinant RAP30 protein

### Protein Details:

RAP30 is a component of human TFIIF, which is involved in initiation and elongation of RNA polymerase II-dependent transcription. Recombinant RAP30 was expressed from full-length (accession number NP\_004119.1) human protein with an N-terminal polyhistadine tag.

### Dilution of Protein:

Recombinant RAP30 is supplied in solution and should be kept as a solution in order to maintain full activity.

### Application:

Recombinant RAP30 is suitable for *in vitro* transcription, protein-protein interactions and protein-DNA interaction studies. A recommended starting point for *in vitro* transcription assays is 8 ng of recombinant protein per reaction. The molecular weight of the protein is ~30 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.