

Recombinant TFIIE β protein

Catalog No: 31341

Source: Human

Quantity: 10 μ g

Expressed In: *E. coli*

Buffer: Recombinant TFIIE β protein supplied at a concentration of 0.1 μ g/ μ l in 20 mM Tris-HCl (pH7.3), 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Contents: 10 μ g recombinant TFIIE β protein.

Protein Details:

Recombinant TFIIE β protein was expressed in *E. coli* as the full-length human protein (accession number NP_002086) with an N-terminal FLAG tag. The molecular weight of the protein is 35 kDa.

Dilution of Protein:

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

Application:

Recombinant TFIIE β 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.