

Recombinant TFIIIB protein

Catalog No: 31240

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 0.4 µg/µl

Source: Human

Buffer Contents: 10 µg of recombinant protein supplied at a concentration of 0.4 µg/µl in 20 mM Tris-HCl, pH 7.3, 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Background: TFIIIB, transcription factor II B (also known as general transcription factor IIB, GTF2B) is required by RNA polymerase II for initiation of transcription. TFIIIB forms a complex with transcription factors IID and IIA (DAB complex) within the nucleus.

Protein Details: Recombinant TFIIIB protein was expressed in *E. coli* as the full-length human protein (accession number AAA61149) with an N-terminal FLAG tag. The molecular weight of the protein is ~36.8 kDa.

Application Notes: Recombinant TFIIIB protein is suitable for use in *in vitro* transcription assays. A recommended starting point for the *in vitro* transcription assay is 10-50 ng recombinant protein per reaction.

Storage and Guarantee: 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.