

Recombinant TCEA1 protein

Catalog No: 31242

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 0.25 µg/µl

Source: Human

Buffer Contents: 10 µg of recombinant protein supplied at a concentration of 0.25 µg/µl in 20 mM Tris-HCl, pH 7.3, 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Background: TCEA1 is transcription elongation factor A protein 1 (also known as transcription elongation factor S-II, TFIIS, TF2S, GTF2S or RNA polymerase II elongation factor). RNA polymerase II is an enzyme found in eukaryotic cells composed of 12 subunits that catalyze transcription of DNA. Transcription consists of three phases: initiation, elongation and termination. TCEA1 (TFIIS) is a transcription elongation factor that stimulates RNA transcript cleavage and polymerase read-through past template-encoded arresting sites.

Protein Details: Recombinant TCEA1 protein was expressed in *E. coli* as the full-length human protein (accession number AAA61138) with an N-terminal FLAG tag. The molecular weight of the protein is ~36 kDa.

Application Notes: Recombinant TCEA1 protein is suitable for use in *in vitro* transcription assays. We recommended 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 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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.