

## Recombinant TFIIA2 protein

---

**Catalog No:** 31249

**Expressed In:** *E. coli*

**Quantity:** 10 µg

**Concentration:** 0.2 µg/µl

**Source:** Human

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

**Background:** TFIIA2 (also known as TF2A2, GTF2A2, general transcription factor IIA subunit 2, TFIIS, T2AG or TFIIA-gamma) is a component of RNA polymerase II transcription complex. TFIIA is a heterodimeric protein consisting of the large unprocessed subunit 1, and the small subunit gamma (TFIIA2). TFIIA associates with TATA-binding subunit (TBP) to regulate transcriptional activity.

**Protein Details:** Recombinant TFIIA2 protein was expressed in *E. coli* as the full-length human protein (accession number AAB58247) with an N-terminal FLAG tag. The molecular weight of the protein is ~12 kDa.

**Application Notes:** Recombinant TFIIA2 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:** 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.