

## Recombinant TAF1 (1398-1524) protein

---

**Catalog No:** 31403, 31803

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

**Quantity:** 100, 1000 µg

**Concentration:** 1 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant TAF1 (1398-1524) protein is supplied at a concentration of 1 µg/µl in 25 mM Tris pH 8.0, 500 mM NaCl, 20% glycerol.

**Background:** TAF1 RNA polymerase II, TATA box binding protein (TBP)-associated factor (TAF1) is the largest component and core scaffold of the TFIID basal transcription factor complex. TAF1 contains novel N- and C-terminal Ser/Thr kinase domains which can autophosphorylate or transphosphorylate other transcription factors. TAF1 also contains bromodomains for interaction with other proteins. The bromodomain functions as a 'reader' of epigenetic histone marks and regulates chromatin structure and gene expression by linking associated proteins to the recognized acetylated nucleosomal targets. TAF1 has been shown to phosphorylate threonine 55 of p53 which leads to MDM2-mediated degradation of p53. TAF1 also phosphorylates GTF2A1 and GTF2F1 on serine residues. In addition, TAF1 possesses DNA-binding activity and is essential for progression of the G1 phase of the cell cycle.

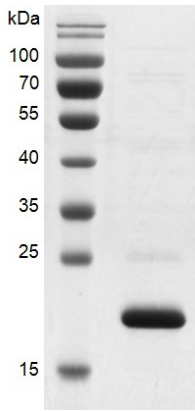
**Protein Details:** Recombinant TAF1 (1398-1524) protein that includes amino acids 1398-1524 of human TAF1 protein (accession number NP\_004597.2) was expressed in *E. coli* and contains an N-terminal His tag and C-terminal FLAG tag with a molecular weight of 21 kDa. The purity of the protein is >90% by SDS-PAGE.

**Application Notes:** Recombinant TAF1 (1398-1524) is suitable for use in binding assays, inhibitor screening, and selectivity profiling.

**Storage and Guarantee:** 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 guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

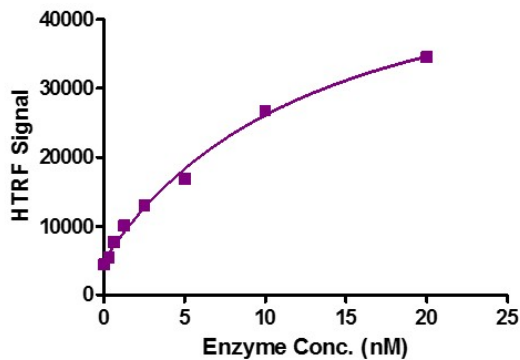
### TAF1 (1398-1524)



### Recombinant TAF1 (1398-1524) protein gel.

TAF1 (1398-1524) protein was run on a 12% SDS-PAGE gel and stained with Coomassie Blue.

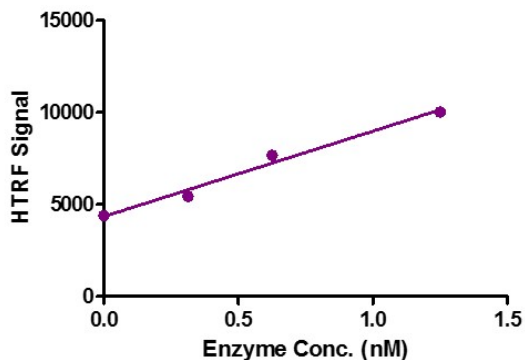
### TAF1 (1398-1524) Titration



### HTRF Assay for Recombinant TAF1 (1398-1524) protein activity.

3  $\mu$ M histone peptide H4K5/K8/K12/K16 (4ac) was incubated with TAF1 (1398-1524) protein in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-Flag antibody was used to detect reaction products.

### TAF1(1398-1524) Titration



### HTRF Assay for Recombinant TAF1 (1398-1524) protein activity.

3  $\mu$ M histone peptide H4K5/K8/K12/K16 (4ac) was incubated with TAF1 (1398-1524) protein in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-Flag antibody was used to detect reaction products.