

## TAF1(C-term) Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # T28-35BG**

Lot # F648-1

### Product Description

Recombinant human TAF1 (1032-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The TAF1 gene accession number is [NM\\_138923](#).

### Gene Aliases

BA2R; CCG1; CCGS; DYT3; DYT3/TAF1; KAT4; N-TAF1; NSCL2; OF; P250; TAF2A

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

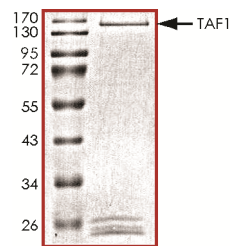
### Scientific Background

TAF1 is a DNA-binding protein complex required for RNA polymerase II-mediated transcription of many, if not all, protein-encoding genes in eukaryotic cells. TAF1 plays a key role in the initiation process, since it binds to the TATA element to form a complex that nucleates the assembly of the other components into a preinitiation complex and that may be stable through multiple rounds of transcription. TAF1 may be targeted to specific chromatin-bound promoters and may play a key role in chromatin recognition (1). TAF1 also serves to link the control of transcription to the cell cycle (2).

### References

- Jacobson, R. H. et.al: Structure and function of a human TAFII250 double bromodomain module. *Science* 288: 1422-1425, 2000.
- Starr, D. B. et.al: TFIID binds in the minor groove of the TATA box. *Cell* 67: 1231-1240, 1991.

### Purity



The purity of TAF1 was determined to be **>70%** by densitometry. Approx. MW **155 kDa**.

## TAF1(C-term) Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number T28-35BG

Specific Lot Number F648-1

Purity	>70%
Concentration	0.1 µg/µl
Stability	1yr At $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**