

TAF1(N-term) Protein

Recombinant human protein expressed in Sf9 cells

Catalog # T28-35G

Lot # X731-2

Product Description

Recombinant human TAF1 (1-561) 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°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.

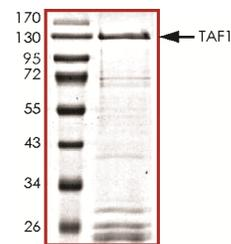
Scientific Background

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 **130 kDa**.

TAF1(N-term) Protein

Recombinant human protein expressed in Sf9 cells

| | |
|---------------------|---|
| Catalog Number | T28-35G |
| Specific Lot Number | X731-2 |
| Purity | >70% |
| Concentration | 0.1 µg/µl |
| Stability | 1yr At -70°C from date of shipment |
| Storage & Shipping | Store product at -70°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
www.signalchem.com

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