

Data Sheet

TAF1L (1398-1516)

Human, recombinant, N-terminal GST tag

Catalog #: 31105

Lot #: 120221 Conc.: 0.41 mg/ml

Formulated in: 50 mM Tris-HCl, pH 8.0,
138 mM NaCl, 2.7 mM KCl, 20 mM
Glutathione, 3 mM EDTA, 10% glycerol
and 2 mM DTT.

Stability: >6 months at -80°C

Reference:

Wang, P.J., *et. al.*, *Hum. Mol. Genet.* 2002
Sep 15; **11(19)**: 2341-2346.

Description:

Human TAF1 RNA polymerase II, TAT box binding protein (TBP)-associated factor, also known as TAF1L, amino-acids 1398 – 1516 (GenBank Accession No. NM_153809) with N-terminal GST tag, MW = 40.6 kDa, expressed in an *E. coli* expression system.

Application:

Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

Quality Assurance

**4-20% SDS-PAGE
Coomassie staining**

Lane 1:

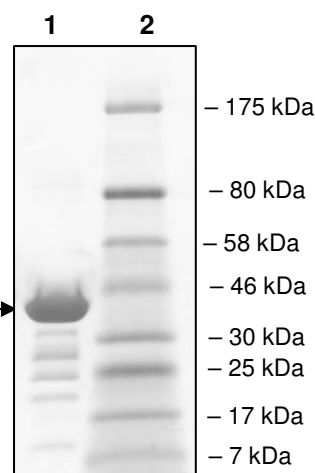
1 µg TAF1L

Lane 2:

Protein Marker
BioLabs (#P7708L)

MW: 40.6 kDa

Purity: ≥82%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com