

Data Sheet

TAF1L (1517-1649)

Human, recombinant, N-terminal HIS tag

Catalog #: 31104

Lot #: 120216G Conc.: 0.30 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.05% Tween-
20, and 20% glycerol.

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 1517 – 1649 (GenBank Accession No. NM_153809) with N-terminal HIS tag, MW = 16.4 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:

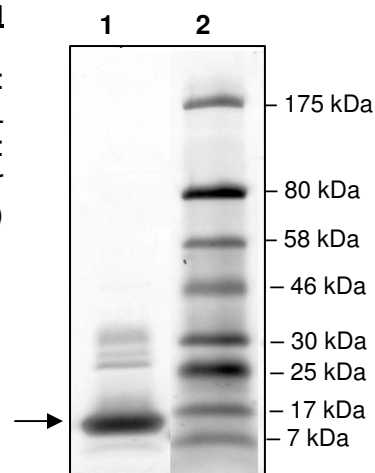
3 µg TAF1L

Lane 2:

Protein Marker
BioLabs (#P7708L)

MW: 16.4 kDa

Purity: ≥79%



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