

## Data Sheet

### TAF1 (1519-1651)

Human, recombinant, N-terminal HIS-tag  
**Catalog #:** 31110  
Lot #: 120323-GF      Conc.: 0.79 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04 % Tween-20, 20% glycerol.

**Stability:** >6 months at -80°C

#### References:

1. Sako W, *et al.*, *Neuroscience*. 2011 Aug 25;189:100-7.
2. Zaborowska J, *et al.*, *Transcription*. 2012 Mar 1;3(2).

#### Description:

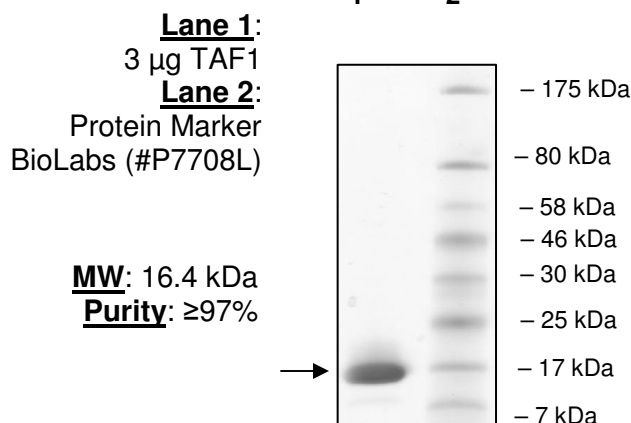
Human Bromodomain TAF1, amino-acids 1519 – 1651 (GenBank Accession No. NM\_004606) with N-terminal HIS-tag, MW = 16.4 kDa, expressed in an *E. coli* expression system.

#### Application:

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

## Quality Assurance

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



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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Protein Sequence:**

MHHHHHLLDDDDQVAFSFDNIIVTQKMMAVPDSWPFHH  
PVNKKFVPDYYKVIVNPM DLETIRKNISKHKYQSRESFLD  
DVNLILANSVKYNGPESQYTKTAQEIVNVCYQTLTEYDEH  
LTQLEKDICTAKEAALEEEAE

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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)