

Data Sheet

TAF1 (1519-1651)

Human, recombinant, N-terminal GST-tag

Catalog #: 31126

Lot #: 121127 **Conc.:** 3.6 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, and 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 GST-tag, MW = 42.2 kDa, expressed in an *E. coli* expression system.

Application:

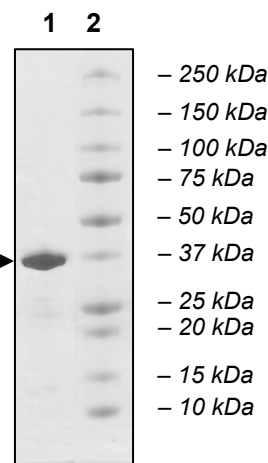
Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2.7 µg TAF1
Lane 2:
Protein Marker

MW: 42.2 kDa →
Purity: ≥90%



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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

*MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKL TQSM IIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSR IAYS KDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKR IEAIPQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDLEVL FQGPLG SLLDDD
DQVAFS FILDNIVTQKMMAVPDSWPFHHPVNKKFVPDYYKVIVNPMDLETIRKNISKHK
YQSRESFLDDVNLILANSVKYNGPESQYTKTAQEIVNVCYQTLTEYDEHLTQLEKDICT
AKEAALEAE*

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