

Recombinant TAF1L (1392-1654) protein

Catalog No: 31436, 31806

Lot No: 12318003

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1.2 µg/µl

Source: Human

Buffer Contents: Recombinant TAF1L (1392-1654) protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, and 10% glycerol.

Background: TAF1 RNA polymerase II, TATA box binding protein (TBP)-associated factor, 210 kDa-like (TAF1L) is homologous to TAF1 and contains bromodomains for interaction with other proteins. The bromodomain functions as a 'reader' of epigenetic histone marks and regulates chromatin structure and gene expression by linking associated proteins to the recognized acetylated nucleosomal targets. TAF1L is specifically expressed in the testis in male germ cells. It may act as a functional substitute for TAF1 during male meiosis when sex chromosomes are transcriptionally silenced.

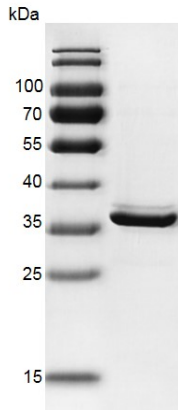
Protein Details: The peptide corresponding to amino acids 1392-1654 that contains the bromodomain sequences of TAF1L (accession number NP_722516.1) was expressed in *E. coli* and contains an N-terminal His-Tag and C-terminal FLAG-Tag with an observed molecular weight of 33.9 kDa.

Application Notes: Recombinant TAF1L (1392-1654) is suitable for use in binding assays, inhibitor screening, and selectivity profiling.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

TAF1L (1392-1654)



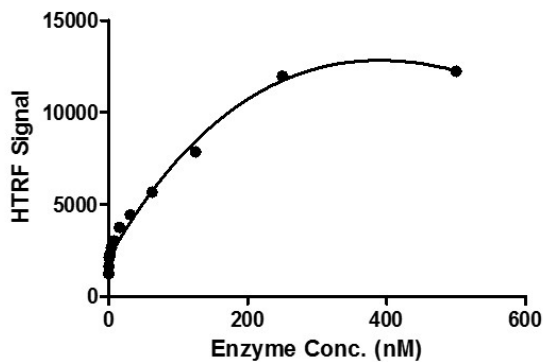
Recombinant TAF1L (1392-1654) protein gel.

TAF1L (1392-1654) protein was run on a 12.5% SDS-PAGE gel and stained with Coomassie Blue.

MW: 33.9 kDa

Purity: > 90%

TAF1L (1392-1654) Titration



Recombinant TAF1L (1392-1654) activity assay.

3 μ M histone peptide H4K5/8/12/16 (4ac) was incubated with TAF1L (1392-1654) in reaction buffer including 50 mM HEPES-NaOH pH 7.4, 0.1% BSA for 1 hour at room temperature. FLAG antibody and SA-XL665 mixture (each 1:100 dilution in Binding Buffer) was used to detect reaction products. All operations were performed at RT. HTRF assay was used for detection.

The file cannot be found:
31436-12318003-3.jpg

Recombinant TAF1L (1392-1654) activity assay.

3 μ M histone peptide H4K5/8/12/16 (4ac) was incubated with TAF1L (1392-1654) in reaction buffer including 50 mM HEPES-NaOH pH 7.4, 0.1% BSA for 1 hour at room temperature. FLAG antibody and SA-XL665 mixture (each 1:100 dilution in Binding Buffer) was used to detect reaction products. All operations were performed at RT. HTRF assay was used for detection.