

Recombinant TAF1L (1402-1521) protein

Catalog No: 31451, 31851

Expressed In: *E. coli*

Quantity: 100 µg

Concentration: 1.5 µg/µl

Source: Human

Buffer Contents: Recombinant TAF1L (1402-1521) protein is supplied at a concentration of 1.5 µg/µl in 25 mM Tris pH 8.0, 500 mM NaCl, 20% 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.

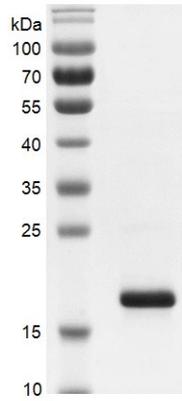
Protein Details: Recombinant TAF1L (1402-1521) protein that includes amino acids 1402-1521 of human TAF1L protein (accession number NP_722516.1) was expressed in *E. coli* and contains an N-terminal His tag and C-terminal FLAG tag with a molecular weight of 20.08 kDa. The purity of the protein is > 92% by SDS-PAGE.

Application Notes: Recombinant TAF1L (1402-1521) 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 (1402-1521)



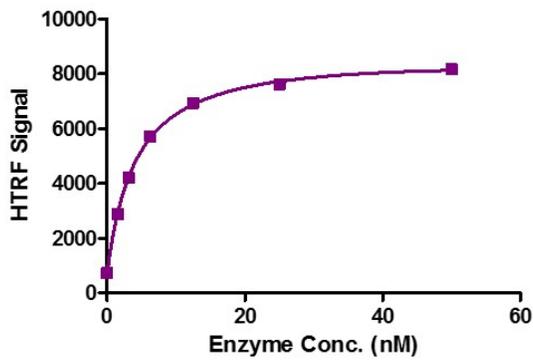
Recombinant TAF1L (1402-1521) protein gel.

TAF1L (1402-1521) protein was run on a 12% SDS-PAGE gel and stained with Coomassie Blue.

MW: 20.1 kDa

Purity: > 92%

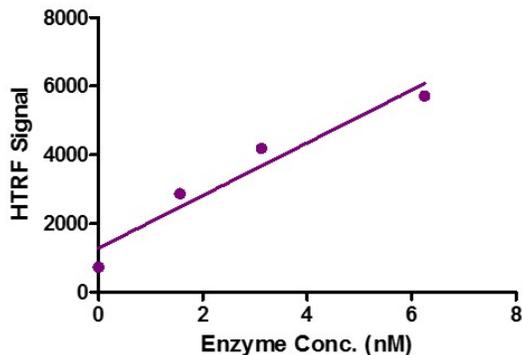
TAF1L (1402-1521) Titration



HTRF Assay for Recombinant TAF1L (1402-1521) protein activity.

3 μ M histone peptide H4K5/8/12/16(4ac) was incubated with TAF1L (1402-1521) in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-FLAG antibody was used to detect reaction products.

TAF1L (1402-1521) Titration



HTRF Assay for Recombinant TAF1L (1402-1521) protein activity.

3 μ M histone peptide H4K5/8/12/16(4ac) was incubated with TAF1L (1402-1521) in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-FLAG antibody was used to detect reaction products.