

Recombinant BPTF / FALZ (2791-2911) protein

Catalog No: 31447, 31847

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1 µg/µl

Source: Human

Buffer Contents: Recombinant BPTF / FALZ (2791-2911) protein is supplied at a concentration of 1 µg/µl in 25 mM Tris pH 8.0, 500 mM NaCl, 20% glycerol.

Background: BPTF (Bromodomain PHD finger transcription factor) is a histone-binding component of NuRF (nucleosome-remodeling factor), a complex which catalyzes ATP-dependent nucleosome sliding and facilitates transcription of chromatin. BPTF protein contains a DNA-binding domain, a zinc finger motif and a bromodomain. Bromodomains recognize acetylated histone lysine residues and function as 'readers' of these epigenetic histone marks to regulate chromatin structure and gene expression by linking associated proteins to the acetylated nucleosomal targets. BPTF also specifically recognizes H3 tails trimethylated on lysine 4 (H3K4me3), which mark transcription start sites of virtually all active genes. It may also regulate transcription through direct binding to DNA or transcription factors.

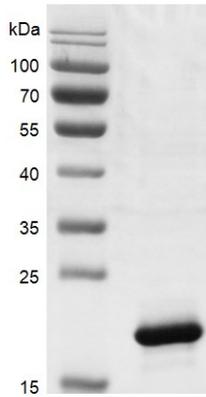
Protein Details: Recombinant BPTF / FALZ (2791-2911) protein that includes amino acids 2791-2911 of human FALZ protein (accession number NP_872579.2) was expressed in *E. coli* and contains an N-terminal His tag and C-terminal FLAG tag with a molecular weight of 20.3 kDa. The purity of the protein is >90% by SDS-PAGE.

Application Notes: Recombinant BPTF / FALZ (2791-2911) 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.

BPTF / FALZ (2791-2911)



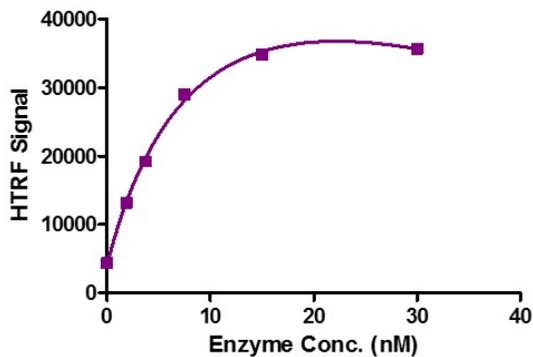
Recombinant BPTF / FALZ (2791-2911) protein gel.

BPTF / FALZ (2791-2911) protein was run on a 12% SDS-PAGE gel and stained with Coomassie Blue.

MW: 20.3 kDa

Purity: > 90%

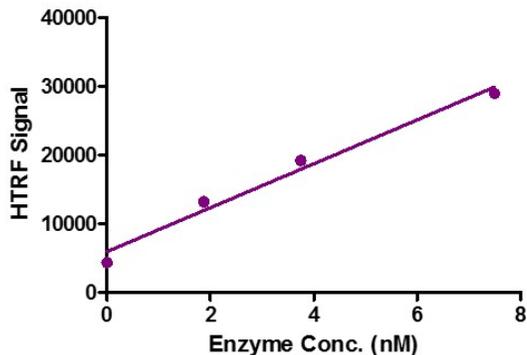
FALZ (2791-2911) Titration



HTRF Assay for Recombinant BPTF / FALZ (2791-2911) activity.

3 μ M histone peptide H4K5/8/12/16(4Ac) was incubated with BPTF / FALZ (2791-2911) protein 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.

FALZ (2791-2911) Titration



HTRF Assay for Recombinant BPTF / FALZ (2791-2911) activity.

3 μ M histone peptide H4K5/8/12/16(4Ac) was incubated with BPTF / FALZ (2791-2911) protein 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.