

Recombinant BRPF3 (576-701) protein

Catalog No: 31487, 31787**Expressed In:** *E. coli***Quantity:** 100, 1000 µg**Concentration:** 3.5 µg/µl**Source:** Human

Buffer Contents: Recombinant BRPF3 (576-701) protein expressed in *E. coli* cells at a concentration of 3.5 µg/µl in 25 mM Tris pH 7.4, 150 mM NaCl, 5% glycerol.

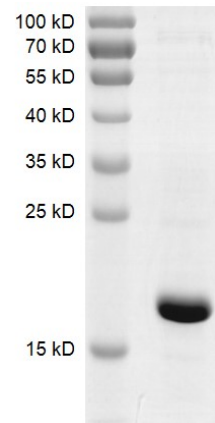
Background: BRPF3 (Bromodomain and PHD finger containing 3) is a component of the MOZ/MORF histone acetyltransferase (HAT) complex. The acetylation of histone lysine residues plays a crucial role in the epigenetic regulation of gene transcription. Acetylated lysine residues can be recognized by bromodomain. The bromodomains serve as “readers” of histone acetylation marks regulating the transcription of target promoters. There is less study on its function now.

Protein Details: The peptide corresponding to amino acids 576-701 that contains the bromodomain sequences of BRPF3 (accession number NM_015695.2) was expressed in *E. coli* and contains an N-terminal His tag and C-terminal FLAG tag with an observed molecular weight of 19.7 kDa. The recombinant protein is >90% pure by SDS-PAGE.

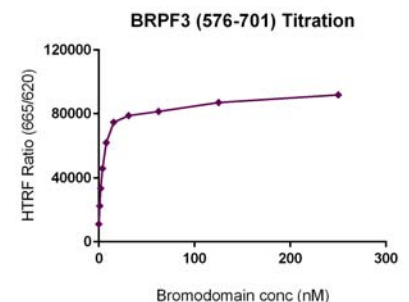
Application Notes: Recombinant BRPF3 (576-701) is suitable for use in binding assays, inhibitor screening, and selectivity profiling.

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.

**Recombinant BRPF3 (576-701) protein gel.**

BRPF3 (576-701) protein was run on an SDS-PAGE gel and stained with Coomassie blue.

**Recombinant BRPF1 (576-701) HTRF activity assay.**

3.3 µM histone peptide H4K5/8/12/16 (4Ac) was incubated with BRPF1 (576-701) protein in reaction buffer including 50mM HEPES-NaOH pH 7.0, 0.1% BSA overnight at room temperature. Anti-Flag antibody was used to detect reaction products.