

## Recombinant BRPF1 (629-742), GST-tag

---

**Catalog No:** 31547, 31947

**Lot No:** 22716001

**Expressed In:** *E. coli*

**Quantity:** 100 µg

**Concentration:** 3 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant BRPF1 (629-742), GST-tag is supplied at a concentration of 10 µg/µl in 25 mM Tris pH 8.0, 300 mM NaCl, 5% glycerol.

**Background:** Bromodomain and PHD finger containing (BRPF1) protein is a component of the MOZ/MORF complex which has histone H3 acetyltransferase activity. The bromodomain of BRPF1 recognizes acetylated histone lysine residues and function as a 'reader' of these epigenetic histone marks to regulate chromatin structure and gene expression by linking associated proteins to the recognized acetylated nucleosomal targets. BRPF1 is known to positively regulate the transcription of RUNX1 and RUNX2.

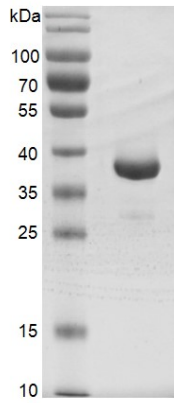
**Protein Details:** The peptide corresponding to amino acids 629 - 742 that contains the bromodomain sequences of BRPF1 (accession number NP\_004625.2) was expressed in *E. coli* cells and contains an N-terminal GST tag with a molecular weight of 39.4 kDa. BRPF1 (629-742), GST-tag is >92% pure by SDS-PAGE.

**Application Notes:** Recombinant BRPF1 (629-742), GST-tag 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.

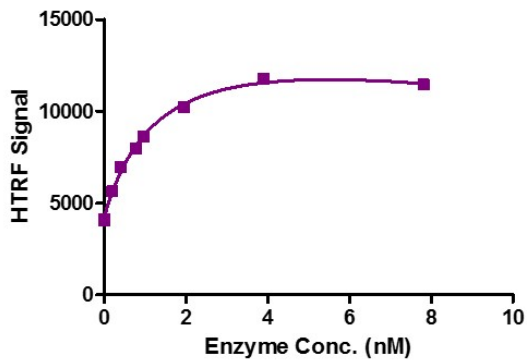
### BRPF1 (629-742)



### Recombinant BRPF1 (629-742), GST-tag gel.

BRPF1 (629-742), GST-tag protein was run on a 12% SDS-PAGE gel and stained with Coomassie Blue.

### GST-BRPF1 (629-742) Titration

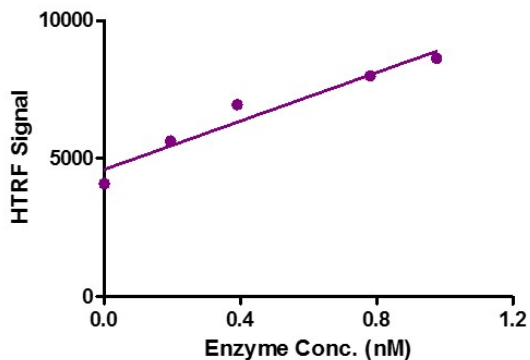


### HTRF assay for BRPF1 (629-742), GST-tag activity

3  $\mu$ M histone peptide H4K5/8/12/16(4Ac) was incubated with BRPF1 (629-742), GST-tag in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature.

Anti-GST antibody was used to detect reaction products.

### GST-BRPF1 (629-742) Titration



### HTRF assay for BRPF1 (629-742), GST-tag activity

3  $\mu$ M histone peptide H4K5/8/12/16(4Ac) was incubated with BRPF1 (629-742), GST-tag in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature.

Anti-GST antibody was used to detect reaction products.