

## Recombinant BRPF1 (1077-1205), GST-tag

**Catalog No:** 31484, 31884

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

**Quantity:** 100, 1000 µg

**Concentration:** 2 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant BRPF1 (1077-1205), GST-tag protein expressed in *E. coli* at a concentration of 2 mg/ml in 25 mM Tris pH 7.4, 150 mM NaCl and 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.

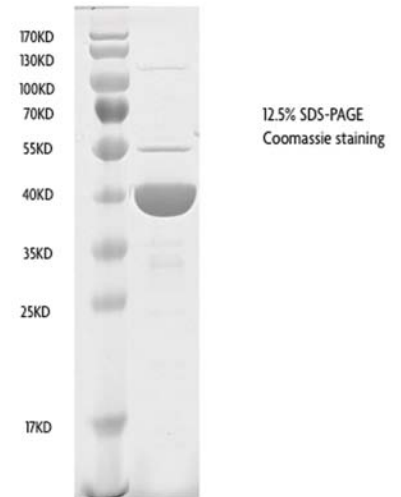
BRPF1 possesses a **PWWP (Pro-Trp-Trp-Pro) domain** that functions in the association of PWWP domain-containing proteins with chromatin. PWWP domains have been shown to bind DNA and also specifically recognized methylated lysine residues of histones. In doing so they are suspected to be involved in the mediation of the cross-talk between DNA and histone epigenetic marks. Mutations in PWWP domains have been shown to be associated with various human diseases, including cancer.

**Protein Details:** The peptide corresponding to amino acids 1077 - 1205 that contains the PWWP domain sequences of BRPF1 (accession number NP\_001003694.1) was expressed in *E. coli* and contains an N-terminal GST tag with an observed molecular weight of 42.8 kDa. The recombinant protein is >92% pure by SDS-PAGE.

**Application Notes:** Recombinant BRPF1 (1077-1205), 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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

BRPF1 (1077-1205)



**Recombinant BRPF1 (1077-1205),  
GST-tag protein gel.**

BRPF1 (1077-1205), GST-tag protein was run on an SDS-PAGE gel and stained with Coomassie blue.