

# PB1 Bromodomain 1

**Catalog No.** 15-0027  
**Lot No.** 13228001  
**Pack Size** 100 µg



**Type** Bromodom **Expressed In** *E. coli*  
**Mol. Wgt.** 40 kDa **Epitope Tag** GST

## Product Description:

Recombinant human PB1 Bromodomain 1 (BAF180, PBRM1, accession number Q86U86, amino acids 32 to 170) containing an N-terminal GST tag, expressed in *E. coli*. PB1 is a member of the BAF transcriptional activation complex and contains six discrete bromodomains. PB1 appears to play a central role, regulating many functions of the BAF complex including chromatin localization, recruitment of protein subunits and alteration of chromatin architecture.

## Formulation:

Recombinant GST-tagged protein at 1 mg/ml in 20 mM Tris-HCl pH 8.0, 200 mM NaCl, 1 mM DTT, 20% glycerol.

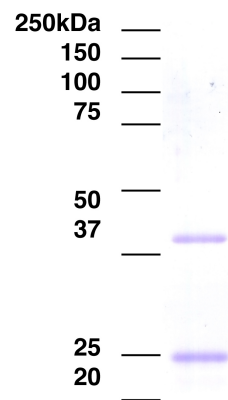
## Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

## Application Notes:

Recombinant PB1 Bromodomain 1 is useful for protein binding and screening experiments examining acetylated protein substrates.

## References:



**Protein Gel Data:** PB1 Bromodomain 1 (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

This product is for *in vitro* research use only and is not intended for use in humans or animals.