

## P300 Bromodomain

**Catalog No.** 15-0037  
**Lot No.** 14231001  
**Pack Size** 100 µg

**Type** Bromodom **Expressed In** *E. coli*  
**Mol. Wgt.** 46 kDa **Epitope Tag** GST

### Product Description:

Recombinant GST-tagged P300 Bromodomain (E1A binding protein p300, EP300, accession number Q09472, amino acids 1031 to 1170) containing an N-terminal GST tag, expressed in *E. coli*. P300 is potent transcriptional co-activator and binds directly to a number of transcription factors including CREB, HIF1A and TFAP2A. P300 contains histone acetyltransferase activity, as well as a Bromodomain, through which it interacts with acetylated histones.

### 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.

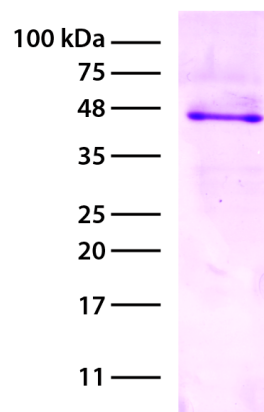
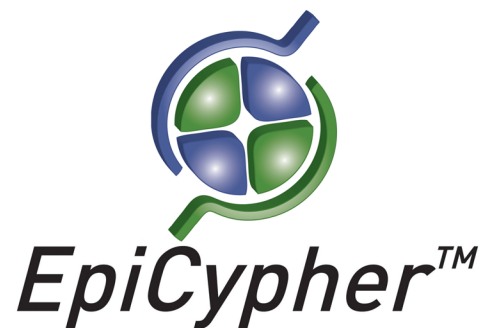
### Storage and Stability:

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

### Application Notes:

Recombinant P300 Bromodomain is useful for protein binding and screening experiments examining acetylated protein substrates.

### References:



**Protein Gel Data:** P300 Bromodomain (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.