

P300 Bromodomain, His-Tagged

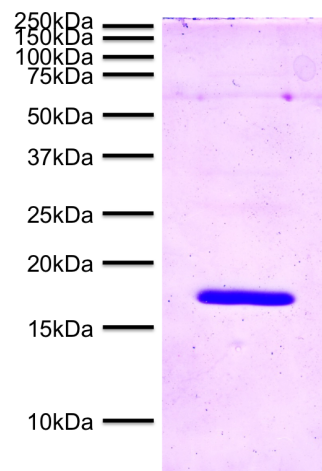
Catalog No. 15-0046
Lot No. 15018001
Pack Size 100 µg



Type Bromodomain **Expressed In** *E. coli*
Mol. Wgt. 18 kDa **Epitope Tag** His

Product Description:

Recombinant 6xHis-tagged P300 Bromodomain (E1A binding protein p300, EP300, accession number Q09472, amino acids 1031 to 1170) containing an N-terminal His-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.



Protein Gel Data: P300 Bromodomain, His-tagged (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

Formulation:

Recombinant 6xHis-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, His-tagged is useful for protein binding and screening experiments examining acetylated protein substrates.

References:

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