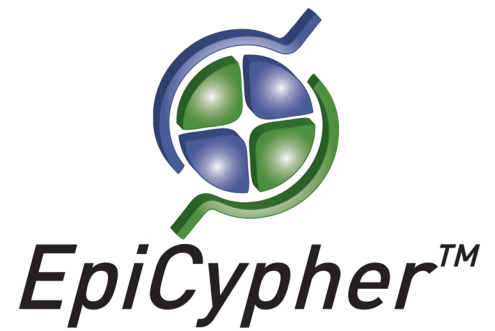


CBP Bromodomain, His-Tagged

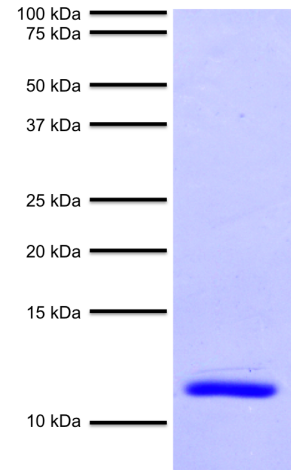
Catalog No. 15-0048
Lot No. 15119001
Pack Size 100 µg



Type Bromodomain **Expressed In** *E. coli*
Mol. Wgt. 12 kDa **Epitope Tag** His

Product Description:

Recombinant human CBP Bromodomain (CREB Binding protein, CREBBP, accession number Q92793, amino acids 1091 to 1190) containing an N-terminal His-tag, expressed in *E. coli*. CBP is an important regulator of cell function and serves to integrate a number of key cell signaling pathways. CBP contains intrinsic histone acetyltransferase (HAT) activity and serves as a co-activator for a large number of transcription factors. The bromodomain of CBP specifically recognizes a number of acetylated proteins, including histones and transcription factors. The CBP gene is mutated in Rubinstein-Taybi Syndrome, a rare developmental disorder.



Protein Gel Data: CBP 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 CBP 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.