

HP1 β , Recombinant Human, His-Tagged



EpiCypher[®]

Catalog No 15-0074
Lot No 19198001
Pack Size 100 μ g

Type Chromodomain **Expressed In** *E. coli*
Mol. Wgt. 24.4 kDa **Epitope Tag** 6xHis

Product Description:

Recombinant human HP1 β (CBX1, M31, Heterochromatin protein 1 beta, accession P83916, amino acids 1-184) containing an N-terminal 6xHis tag, expressed in *E. coli*. HP1 β is a component of constitutive (transcriptionally silent) heterochromatin, recruited through the interaction of the Chromodomain with histone H3 methylated at lysine 9 (H3K9me).

Formulation:

Recombinant 6xHis-tagged protein at 0.94 mg/mL in 20 mM Tris HCl pH 7.5, 100 mM NaCl, 1 mM DTT, 5% glycerol.

Storage and Stability:

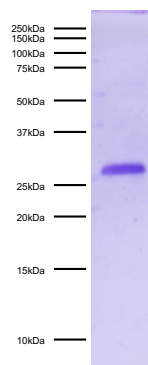
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

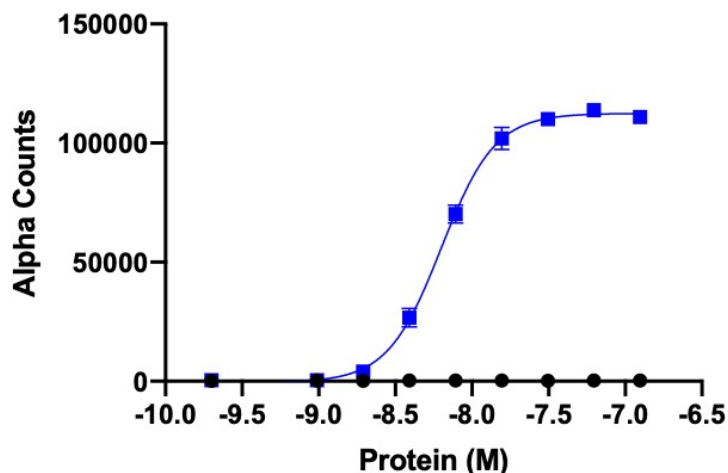
Recombinant HP1 β Chromodomain is useful for protein binding and screening experiments examining methylated protein and peptide substrates.

References:

Li Y et al (2014) Cell 159: 558-571.



Protein Gel Data: HP1 β , Recombinant Human, His-Tagged (1 μ g) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.



AlphaScreen Data: HP1 β , Recombinant Human, His-Tagged used in nucleosome binding studies to detect interactions with specific post-translational modifications. The identities of the nucleosomes tested are as follows: rNuc unmodified (black), H3K9me3 (blue).

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

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

EpiCypher, Inc • Phone: 855-374-2461 • Fax: 855-420-6111 • www.epicypher.com