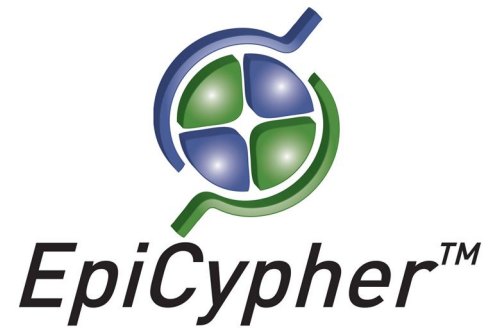


HP1 β Chromodomain, Recombinant Human

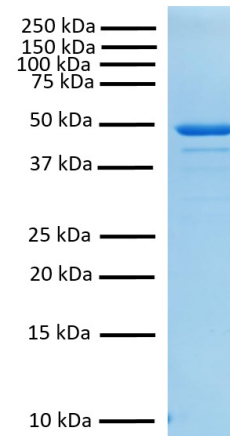
Catalog No. 15-0058
Lot No. 17006001
Pack Size 100 μ g



Type Chromodomain **Expressed In** *E. coli*
Mol. Wgt. 49 kDa **Epitope Tag** GST

Product Description:

Recombinant human HP1 β (CBX1, M31, Heterochromatin protein 1 beta, accession P83916, amino acids 1-185) containing an N-terminal GST 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 lysing 9 (H3K9me).



Protein Gel Data: HP1 β Chromodomain (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.

Formulation:

Recombinant GST-tagged protein at 1 mg/ml in 50 mM Tris-HCl pH 7.0, 300 mM NaCl, 1 mM DTT, 10% 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:

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