

Histone H2B

CATALOG NO.: HMT-11-147

LOT NO.:

DESCRIPTION: Full-length, untagged human recombinant histone H2B (residues 1-126; Genbank Accession # NM_080593; MW = 13.9 kDa) expressed in *E. coli*. One of the four, core nucleosomal histones, H2B forms dimers with histone H2A, two of which associate with the H3-H4 tetramer (H3-H4)₂, forming the core octamer of the nucleosome (see review¹). Histone H2B is subject to several regulatory post-translational modifications including monoubiquitination², phosphorylation³, acetylation⁴, lysine methylation⁴ and crotonylation⁵. Monoubiquitination of H2B lysine-120 (H2BK120ub) stimulates H3K4 methylation in chromatin⁶, an effect which recent evidence suggests is mediated in trans via ubiquitin binding by Ash2L, a component of the human MLL/SET1 family methyltransferase complexes⁷. Monoubiquitination at both H2BK34⁸ and H2BK120⁹ has been shown to enhance H3K79 methylation by Dot1L. RBC's Histone useful for the assay of histone acetyltransferases (HATs; e.g. p300, CBP; see Figure, below) either by radiolabeling with [³H]-S-adenosylmethionine (e.g. gel electrophoresis/autoradiography or filterplate/scintillation counting) or by non-radiolabeled methods (e.g. detection with site-specific anti-acetyllysine antibodies).

PURITY: >90% by SDS-PAGE.

ASSAY CONDITIONS: Histone acetyltransferase assays (see figure, below) were performed with histone H2B and [³H]-Acetyl-CoA as substrates. Activity was determined as TCA-precipitated counts in a scintillation/filter plate assay (Multiscreen FB, Topcount). Reaction conditions: 50 mM Tris-HCl, pH 8.0, 50 mM NaCl, 1 mM DTT, 0.1 mM EDTA, 0.1 mM PMSF, 30°C, 60 min.

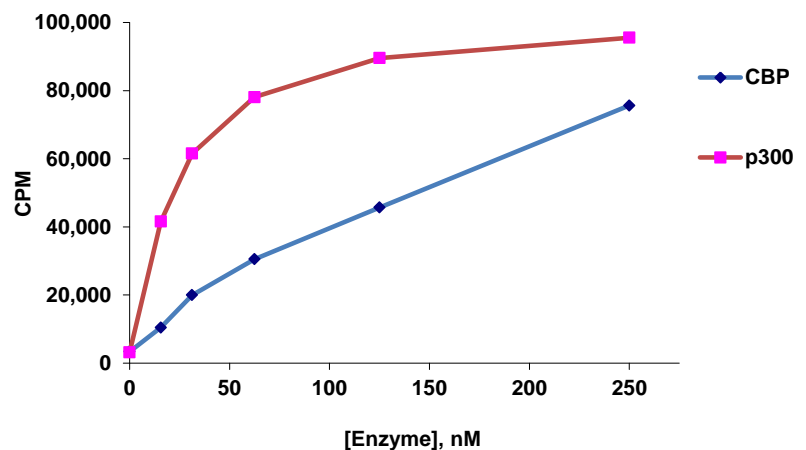
SUPPLIED AS: ___ µg/µl total protein in in 20 mM sodium phosphate pH 7.5, 300 mM NaCl, 1 mM DTT, 1 mM EDTA, 10% glycerol (v/v) as determined by OD₂₈₀

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.

REFERENCES: 1) J.J. Wyrick *et al. Biochim. Biophys. Acta* 2012 **1819** 892; 2) B. Zhu *et al. Mol. Cell.* 2005 **20** 601; 3) W.L. Cheung *et al. Cell* 2003 **113** 507; 4) H.C. Beck *et al. Mol. Cell. Proteomics* 2006 **5** 1314; 5) M. Tan *et al. Cell* 2011 **146** 1016; 6) J. Kim *et al. Cell* 2009 **137** 459; 7) L. Wu *et al. Mol. Cell* 2013 doi: 10.1016/j.molcel.2013.01.033. [Epub ahead of print]; 8) L. Wu *et al. Mol. Cell* 2011 **43** 132; 9) S.J. Whitcomb *et al. J. Biol. Chem.* 2012 **287** 23718



Coomassie blue stained SDS-PAGE (16% acrylamide) of 2 µg of purified Histone H2B. MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



Assay of Histone Acetyltransferases with Histone H2B. The indicated concentrations of either p300 or CBP were assayed as described above in 25 µL reactions, 60 min., 30°C, with 3 µM [³H]-Acetyl-CoA and 5 µM histone H2B as substrates.