

## Histone H2B (Ab-12) Antibody [#D0024]

<b>Catalog Number:</b>	D0024
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P57053
<b>Other Names:</b>	Histone cluster 1, H2bb,H2B.1; H2B/f; H2BFF, H2B histone family, member F,histone 1, H2bb
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-acetylated peptide derived from human Histone H2B around the acetylation site of lysine 12 (P-K-K <sup>A</sup> -G-S).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Histone H2B (Ab-12) antibody detects endogenous levels of total Histone H2B protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Michael A. Parra, Mol. Cell. Biol., May 2006; 26: 3842 - 3852. Fiona A. Myers, J. Biol. Chem., Sep 2003; 278: 36315 - 36322. Débora Bonenfant, Mol. Cell. Proteomics, Mar 2006; 5: 541 - 552.