

Histone H4 (Acetyl-Lys5) Antibody [#D0011]

Catalog Number:	D0011
Concentration:	1mg/ml
Swiss-Prot No.:	P62805
Other Names:	H4F2, HIST, HIST1H4A, Histone H4
All Sites:	Human: Lys5; Mouse: Lys5; Rat: Lys5
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized acetylated peptide derived from human Histone H4 around the acetylation site of lysine 5 (R-G-K ^A -G-G).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific acetylated peptide. The antibody against non-acetylated peptide was removed by chromatography using non-acetylated peptide corresponding to the acetylation site.
Specificity:	Histone H4 (Acetyl-Lys5) antibody detects endogenous levels of Histone H4 only when acetylated at lysine 5.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	M Couppez, J. Biol. Chem., Feb 1987; 262: 2854 - 2860. BM Turner, J. Cell Sci., Jun 1990; 96: 335. KW Marvin, J. Biol. Chem., Nov 1990; 265: 19839 - 19847.