

Histone H4 (Ab-8) Antibody [#D0032]

Catalog Number:	D0032
Concentration:	1mg/ml
Swiss-Prot No.:	P62805
Other Names:	H4F2, HIST, HIST1H4A, Histone H4
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 non-acetylated peptide derived from human Histone H4 around the acetylation site of lysine 8 (G-G-K ^P -G-L).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Histone H4 (Ab-8) antibody detects endogenous levels of total Histone H4 protein.
Reactivity:	Human, Mouse, Rat
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	KW Marvin, J. Biol. Chem., Nov 1990; 265: 19839 - 19847. TA Chen, J. Biol. Chem., Apr 1991; 266: 6489 - 6498. Michael J. Hendzel, Mol. Biol. Cell, Sep 1998; 9: 2491.