

Histone 1F0 Antibody [#C10293]

Catalog Number:	C10293
Concentration:	1mg/ml
Swiss-Prot No.:	P07305
Other Names:	H1, H1 histone family, member 0, H10, H1FV, histone H1', histone H1.0
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 peptide derived from internal of human Histone 1F0.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Histone 1F0 antibody detects endogenous levels of total Histone 1F0 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Doenecke D., J. Mol. Biol. 187:461-464(1986). Collins J.E., Genome Biol. 5:RESEARCH84.1-RESEARCH84.11(2004). Dunham I., Nature 402:489-495(1999).