

HBP1 Antibody [#C10135]

Catalog Number:	C10135
Concentration:	1mg/ml
Swiss-Prot No.:	O60381
Other Names:	HMG box-containing protein 1, HMG box-containing protein 1, High mobility group box transcription factor 1, HBP1
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 HBP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HBP1 antibody detects endogenous levels of total HBP1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Lemerrier C., Mol. Cell. Biol. 20:6627-6637(2000). Hillier L.W., Nature 424:157-164(2003). Ota T., Nat. Genet. 36:40-45(2004).