

HOXB9 Antibody [#C10404]

Catalog Number:	C10404
Concentration:	1mg/ml
Swiss-Prot No.:	P17482
Other Names:	Homeobox protein Hox-B9, Hox-2E, Hox-2.5, HOXB9, HOX2E
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 HOXB9.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HOXB9 antibody detects endogenous levels of total HOXB9 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Kosaki K., Teratology 65:50-62(2002). Giampaolo A., Differentiation 40:191-197(1989). Boncinelli E., Genome 31:745-756(1989).