

## HOXA1/B1/D1 Antibody [#C10372]

<b>Catalog Number:</b>	C10372
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Homeobox protein Hox-B1, Hox-2l, HOXB1, HOX2l
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human HOXA1/B1/D1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	HOXA1/B1/D1 antibody detects endogenous levels of total HOXA1/B1/D1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:20000
<b>References:</b>	Acampora D., Nucleic Acids Res. 17:10385-10402(1989). Boncinelli E., Genome 31:745-756(1989). Piper D.E., Cell 96:587-597(1999).