

## HOXA7 Antibody [#C10524]

<b>Catalog Number:</b>	C10524
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P31268
<b>Other Names:</b>	Homeobox protein Hox-A7, Hox-1A, Hox 1.1, HOXA7, HOX1A
<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 HOXA7.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	HOXA7 antibody detects endogenous levels of total HOXA7 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	La Celle P.T., J. Biol. Chem. 276:32844-32853(2001). McIlhatton M.A., Biochim. Biophys. Acta 1442:329-333(1998). Kim M.H., Mol. Biotechnol. 14:19-24(2000).