

## ZNF7 Antibody [#C10399]

<b>Catalog Number:</b>	C10399
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P17097
<b>Other Names:</b>	Zinc finger protein 7, Zinc finger protein KOX4, Zinc finger protein HF.16, ZNF7, KOX4
<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 ZNF7.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ZNF7 antibody detects endogenous levels of total ZNF7 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 78%, Positives = 85%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Lania L., Genomics 6:333-340(1990). The MGC Project Team, Genome Res. 14:2121-2127(2004). Thiesen H.-J., New Biol. 2:363-374(1990).