

ZNF24 Antibody [#C10397]

Catalog Number:	C10397
Concentration:	1mg/ml
Swiss-Prot No.:	P17028
Other Names:	Zinc finger protein 18, Zinc finger protein KOX11, Heart development-specific gene 1 protein, Zinc finger protein with KRAB and SCAN domains 6, ZNF18, HDSG1, KOX11, ZKSCAN6
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 N-terminal of human ZNF24.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZNF24 antibody detects endogenous levels of total ZNF24 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Guo L.L., Yi Chuan 27:523-530(2005). Ota T., Nat. Genet. 36:40-45(2004). Thiesen H.-J., New Biol. 2:363-374(1990).