

MEOX2 Antibody [#C10662]

Catalog Number:	C10662
Concentration:	1mg/ml
Swiss-Prot No.:	P50222
Other Names:	Mesenchyme homeobox 2 Growth arrest-specific homeobox
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 MEOX2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MEOX2 antibody detects endogenous levels of total MEOX2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Grigoriou M., Genomics 26:550-555(1995). Lepage D.F., Genomics 24:535-540(1994). Hillier L.W., Nature 424:157-164(2003).