

ZHX2 Antibody [#C11999]

Catalog Number:	C11999
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y6X8
Other Names:	AFP regulator 1, AFR1; RAF;, alpha-fetoprotein regulator 1, regulator of AFP, zinc finger and homeodomain protein 2
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 ZHX2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZHX2 antibody detects endogenous levels of total ZHX2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Kawata H., Biochem. J. 373:747-757(2003). The MGC Project Team; Genome Res. 14:2121-2127(2004). Nagase T., DNA Res. 5:355-364(1998).