

ADH7 Antibody [#C14383]

Catalog Number:	C14383
Concentration:	1mg/ml
Swiss-Prot No.:	P40394
Other Names:	ADH-4; alcohol dehydrogenase class 4 mu/sigma; alcohol dehydrogenase class IV mu/sigma; EC 1.1.1.1; gastric alcohol dehydrogenase; retinol dehydrogenase
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 ADH7.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ADH7 antibody detects endogenous levels of total ADH7 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Satre M.A., J. Biol. Chem. 269:15606-15612(1994). Zgombic-Knight M., J. Biol. Chem. 270:4305-4311(1995). Farres J., Eur. J. Biochem. 224:549-557(1994).