

ALDH1B1 Antibody [#C14388]

Catalog Number:	C14388
Concentration:	1mg/ml
Swiss-Prot No.:	P30837
Other Names:	acetaldehyde dehydrogenase 5; AL1B1; aldehyde dehydrogenase X, mitochondrial; ALDH class 2; ALDH5; ALDHX; EC 1.2.1.3
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 ALDH1B1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ALDH1B1 antibody detects endogenous levels of total ALDH1B1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Hsu L.C., J. Biol. Chem. 266:12257-12265(1991). Kalnine N., Submitted (MAY-2003) to the EMBL/GenBank/DDBJ databases. Humphray S.J., Nature 429:369-374(2004).