

ALDOA Antibody [#C10268]

Catalog Number:	C10268
Concentration:	1mg/ml
Swiss-Prot No.:	P04075
Other Names:	ALDA, ALDOA, ALFA, EC 4.1.2.13, Fructose-bisphosphate aldolase A, Lung cancer antigen NY-LU-1, Muscle-type aldolase
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 ALDOA.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ALDOA antibody detects endogenous levels of total ALDOA protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Sakakibara M., Biochem. Biophys. Res. Commun. 131:413-420(1985). Izzo P., Eur. J. Biochem. 164:9-13(1987). Mukai T., Eur. J. Biochem. 195:781-787(1991)