

ME2 Antibody [#C16813]

Catalog Number:	C16813
Concentration:	1mg/ml
Swiss-Prot No.:	P23368
Other Names:	EC 1.1.1.38; malate dehydrogenase; malic enzyme 2; Malic enzyme, mitochondrial; MAOM; NAD-dependent malic enzyme, mitochondrial; NAD-ME; pyruvic-malic carboxylase
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 ME2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ME2 antibody detects endogenous levels of total ME2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Loeber G., J. Biol. Chem. 266:3016-3021(1991). Yanaihara N., Genomics 72:169-179(2001). Abts H.; Thesis (1995), University of Cologne, Germany.