

PDK2 Antibody [#C18101]

Catalog Number:	C18101
Concentration:	1mg/ml
Swiss-Prot No.:	Q15119
Other Names:	EC 2.7.11.2; kinase Pyruvate dehydrogenase kinase 2; PDK2; Pyruvate dehydrogenase [lipoamide] kinase isozyme 2, mitochondrial precursor
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 PDK2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PDK2 antibody detects endogenous levels of total PDK2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Gudi R., J. Biol. Chem. 270:28989-28994(1995). The MGC Project Team, Genome Res. 14:2121-2127(2004). Greenman C., Nature 446:153-158(2007).