

ADCK1 Antibody [#C11274]

Catalog Number:	C11274
Concentration:	1mg/ml
Swiss-Prot No.:	Q86TW2
Other Names:	Uncharacterized aarF domain-containing protein kinase 1 [Precursor], EC 2.7.11.-, ADCK1
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 ADCK1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ADCK1 antibody detects endogenous levels of total ADCK1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	The MGC Project Team, Genome Res. 14:2121-2127(2004). Li W.B., Submitted (FEB-2003) to the EMBL/GenBank/DDBJ databases.