

ADCK2 Antibody [#C11266]

Catalog Number:	C11266
Concentration:	1mg/ml
Swiss-Prot No.:	Q7Z695
Other Names:	Uncharacterized aarF domain-containing protein kinase 2, EC 2.7.11.-, ADCK2, AARF
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 ADCK2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ADCK2 antibody detects endogenous levels of total ADCK2 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Huang C.Q., Submitted (MAY-2003) to the EMBL/GenBank/DDBJ databases. Hillier L.W., Nature 424:157-164(2003). The MGC Project Team, Genome Res. 14:2121-2127(2004).