

KAD5 Antibody [#C11988]

Catalog Number:	C11988
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y6K8
Other Names:	Adenylate kinase isoenzyme 5, EC 2.7.4.3, ATP-AMP transphosphorylase, AK5
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 KAD5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KAD5 antibody detects endogenous levels of total KAD5 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Van Rompay A.R., Eur. J. Biochem. 261:509-517(1999). The MGC Project Team; Genome Res. 14:2121-2127(2004).