

DAK Antibody [#C11155]

Catalog Number:	C11155
Concentration:	1mg/ml
Swiss-Prot No.:	Q3LXA3
Other Names:	Dihydroxyacetone kinase, EC 2.7.1.29, Glycerone kinase, DHA kinase, DAK
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 DAK.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	DAK antibody detects endogenous levels of total DAK protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Ota T., Nat. Genet. 36:40-45(2004). Totoki Y., Submitted (APR-2005) to the EMBL/GenBank/DDBJ databases. The MGC Project Team, Genome Res. 14:2121-2127(2004).