

DGKI Antibody [#C10197]

Catalog Number:	C10197
Concentration:	1mg/ml
Swiss-Prot No.:	O75912
Other Names:	DAG kinase iota, DGK-IOTA, DGK-iota, DGKI, EC 2.7.1.107, diacylglycerol kinase, iota, diglyceride kinase
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 C-terminal of human DGKI.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	DGKI antibody detects endogenous levels of total DGKI protein.
Reactivity:	Human, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Ding L., J. Biol. Chem. 273:32746-32752(1998). Bowne S.J., Mol. Vis. 6:6-9(2000).