

GAK Antibody [#C10064]

Catalog Number:	C10064
Concentration:	1mg/ml
Swiss-Prot No.:	O14976
Other Names:	Cyclin G-associated kinase, EC 2.7.11.1, kinase GAK
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 GAK.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GAK antibody detects endogenous levels of total GAK protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Kimura S.H., Genomics 44:179-187(1997). The MGC Project Team; Genome Res. 14:2121-2127(2004). Greener T., J. Biol. Chem. 275:1365-1370(2000).