

CDK8 Antibody [#C10643]

Catalog Number:	C10643
Concentration:	1mg/ml
Swiss-Prot No.:	P49336
Other Names:	Cell division protein kinase 8, EC 2.7.11.22, EC 2.7.11.23, Protein kinase K35, CDK8
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 N-terminal of human CDK8.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CDK8 antibody detects endogenous levels of total CDK8 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Tassan J.-P., Proc. Natl. Acad. Sci. U.S.A. 92:8871-8875(1995). Fryer C.J., Mol. Cell 16:509-520(2004).