

CSK Antibody [#C10598]

Catalog Number:	C10598
Concentration:	1mg/ml
Swiss-Prot No.:	P41240
Other Names:	Tyrosine-protein kinase CSK, EC 2.7.10.2, C-SRC kinase, Protein-tyrosine kinase CYL, CSK
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 CSK.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CSK antibody detects endogenous levels of total CSK protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Partanen J., Oncogene 6:2013-2018(1991). Braeuninger A., Proc. Natl. Acad. Sci. U.S.A. 88:10411-10415(1991). Brauninger A., Gene 110:205-211(1992).