

CKS2 Antibody [#C10543]

Catalog Number:	C10543
Concentration:	1mg/ml
Swiss-Prot No.:	P33552
Other Names:	Cyclin-dependent kinases regulatory subunit 2, CKS-2, CKS2
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 CKS2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CKS2 antibody detects endogenous levels of total CKS2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Richardson H.E., Genes Dev. 4:1332-1344(1990). The MGC Project Team, Genome Res. 14:2121-2127(2004). Parge H.E., Science 262:387-395(1993).