

CDC6 Antibody [#C12172]

Catalog Number:	C12172
Concentration:	1mg/ml
Swiss-Prot No.:	Q99741
Other Names:	CDC18, CDC6 cell division cycle 6 homolog, CDC6-related protein
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 CDC6.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CDC6 antibody detects endogenous levels of total CDC6 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Williams R.S., Proc. Natl. Acad. Sci. U.S.A. 94:142-147(1997). Saha P., Mol. Cell. Biol. 18:2758-2767(1998). Yu L.-R., J. Proteome Res. 6:4150-4162(2007).