

Clock Antibody [#C10086]

Catalog Number:	C10086
Concentration:	1mg/ml
Swiss-Prot No.:	O15516
Other Names:	circadian locomotor output cycles kaput, hCLOCK, CLOCK, KIAA0334
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 Clock.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Clock antibody detects endogenous levels of total Clock protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Steeves T.D.L., Genomics 57:189-200(1999). Nagase T., DNA Res. 4:141-150(1997). Sato T.K., Nat. Genet. 38:312-319(2006).