

Period Circadian Protein 2 Antibody [#B1223]

Catalog Number:	B1223
Concentration:	1mg/ml
Swiss-Prot No.:	O15055
Other Names:	KIAA0347, Period circadian protein 2, Mper2, Per2
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 Period Circadian Protein 2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Period Circadian Protein 2 antibody detects endogenous levels of total Period Circadian Protein 2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 90%, Positives = 90%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Ikeda M., Submitted (MAR-1998) to the EMBL/GenBank/DDBJ databases. Camacho F., FEBS Lett. 489:159-165(2001). Nagase T., DNA Res. 4:141-150(1997).