

Cyclin H (Ab-315) Antibody [#B0881]

Catalog Number:	B0881
Concentration:	1mg/ml
Swiss-Prot No.:	P51946
Other Names:	Cyclin-H, MO15-associated protein, p37, p34, CCNH
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 non-phosphopeptide derived from human Cyclin H around the phosphorylation site of threonine 315 (E-W-T ^P -D-D).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Cyclin H (Ab-315) antibody detects endogenous levels of total Cyclin H protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Maekelae T.P., Nature 371:254-257(1994). Fisher R.P., Cell 78:713-724(1994). Shiekhattar R., Nature 374:283-287(1995).