

Cyclin B1 (Ab-147) Antibody [#B0635]

Catalog Number:	B0635
Concentration:	1mg/ml
Swiss-Prot No.:	P14635
Other Names:	CCN-2, CCNB, CCNB1, CCNB1-RS1, CYCB, CYCB1, G2/mitotic-specific cyclin B1
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 B1 around the phosphorylation site of serine 147 (A-F-S ^P -D-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Cyclin B1 (Ab-147) antibody detects endogenous levels of total Cyclin B1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Catherine A. Rugg, AACR Meeting Abstracts, Apr 2005; 2005: 1047. Jianping Gong, Cancer Res., Aug 1994; 54: 4285 - 4288. Lourdes A. Gomez, Mol. Cancer Ther., May 2007; 6: 1534 - 1543.