

Cyclin B1 (Ab-126) Antibody [#B0068]

- Catalog Number:** BOO68-1, BOO68-2, BOO68
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P14635
- All Names:** CCN-2, CCNB, CCNB1, CCNB1-RS1, CYCB, CYCB1, G2/mitotic-specific cyclin B1
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human Cyclin B1 around the phosphorylation site of serine 126 (T-A-S-P-S).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Cyclin B1 (Ab-126) antibody detects endogenous levels of total Cyclin B1 protein.
- Reactivity:** Human, Mouse
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
IF: 1:500~1:1000 ELISA: 1:20000
- References:** Jing Yang, J. Biol. Chem., Jan 2001; 276: 3604 - 3609.
Susan Walsh, Mol. Cancer Res., Feb 2003; 1: 280.
Florian Bassermann, J. Biol. Chem., Mar 2007; 10.1074.
Jianke Li, PNAS, Jan 1997; 94: 502.