

## Anti-Cyclin B1

Rabbit Polyclonal Antibody

Catalog # C79-362R

Lot # B3216-38

### Cited Applications

ELISA, IP, WB

Ideal working dilutions for each application should be empirically determined by the investigator.

### Specificity

Recognizes the 55-66kDa Cyclin B1 protein

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, Antiserum

### Immunogen

The antibody was produced against full length fusion protein corresponding to the human Cyclin B1 protein.

### Formulation

0.01% (w/v) Sodium Azide

### Stability

1yr at  $-20^{\circ}$ C from date of shipment

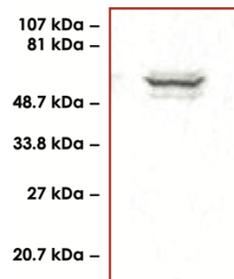
### Scientific Background

The protein encoded by this gene is a regulatory protein involved in mitosis. The gene product forms a complex with p34 (cdc2) to form the maturation-promoting factor (MPF). Two alternative transcripts have been found, a constitutively expressed transcript and a cell cycle-regulated transcript that is expressed predominantly during G2/M phase. The different transcripts result from the use of alternate transcription initiation sites.

### References:

1. Giordano, A., et al.: A 60 kd cdc-2 associated polypeptide complexes with the E1A proteins of adenovirus-infected cells. *Cell*. 1989; 58: 981-990.
2. Pines, J., et al.: Human cyclin A is adenovirus E1A-associated protein p60 and behaves differently from cyclin B. *Nature*. 1990; 346: 760-763.

### Sample Data



Western blot analysis using Anti-Cyclin B1 antibody (1:500) shows detection of  $\sim$ 62kDa human Cyclin B<sub>1</sub> present in asynchronous HN30 cell lysates. HN30 cells are from head and neck cancer tumors that over express cyclin B<sub>1</sub> and D<sub>1</sub>.

## Anti-Cyclin B1

Rabbit Polyclonal Antibody

Catalog #	C79-362R
Lot #	B3216-38
Purification	Delipidation and Defibrination
Stability	1yr at $-20^{\circ}$ C from date of shipment
Storage & Shipping	Store product at $-20^{\circ}$ C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**