

## Anti-CDK1

Rabbit Polyclonal Antibody

Catalog # C22-63R

Lot # B3216-27

### Cited Applications

ELISA, IHC, IP, WB

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

### Specificity

Recognizes the 34kDa cdc2 protein

### Cross Reactivity

Human, Rat and Mouse

### Host/Isotype/Clone#

Rabbit, Antiserum

### Immunogen

The antibody was produced against synthesized peptide corresponding to the C-terminus of the human protein conjugated to KLH.

### Formulation

0.01% (w/v) Sodium Azide

### Stability

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

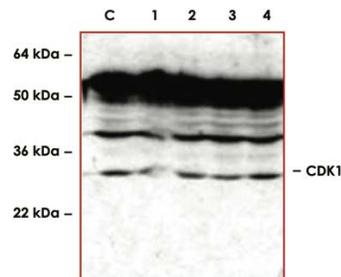
### Scientific Background

CDK1 or Cell Division Control protein 1 is essential for the completion of START, the controlling event in the cell cycle that is required to initiate mitosis. CDK1 is a catalytic subunit of a protein kinase complex, called the M-Phase Promoting Factor that induces entry into mitosis and is universal among eukaryotes (1). Phosphorylation of Bcl-2 in G2/M phase-arrested cells following photodynamic therapy with hypericin involves a CDK1-mediated signal and delays the onset of apoptosis. Therapeutic potential of CDK inhibitor NU2058 in androgen-independent prostate cancer has also been demonstrated (2)

### References:

1. Vantieghem, A. et al: Phosphorylation of Bcl-2 in G2/M phase-arrested cells following photodynamic therapy with hypericin involves a CDK1-mediated signal and delays the onset of apoptosis. J. Biol. Chem. 2002; 277(40):37718-31.
2. Rigas, A.C. et al: Therapeutic potential of CDK inhibitor NU2058 in androgen-independent prostate cancer. Oncogene. 2007; 18.

### Sample Data



Western blot using Anti-CDK1 antibody to detect CDK1 in MCF-7 cell lysates.

## Anti-CDK1

Rabbit Polyclonal Antibody

Catalog #	C22-63R
Lot #	B3216-27
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.**