

Anti-CDK2

Rabbit Polyclonal Antibody

Catalog # **C29-63R**

Lot # **B3216-31**

Cited Applications

ELISA, IHC, IP, WB

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

Specificity

Recognizes the 33kDa CDK2 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 CDK2 protein conjugated to KLH.

Formulation

0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

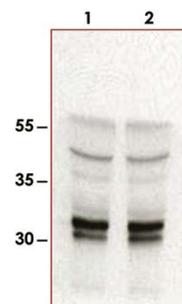
Scientific Background

Anti cdk2 antibody recognizes a protein encoded by a gene member of the Ser/Thr protein kinase family. This protein kinase is highly similar to the gene products of *S. cerevisiae* cdc28, and *S. pombe* cdc2. It is a catalytic subunit of the cyclin-dependent protein kinase complex, whose activity is restricted to the G1-S phase, and essential for cell cycle G1/S phase transition. This protein associates with and regulated by the regulatory subunits of the complex including cyclin A or E, CDK inhibitor p21Cip1 (CDKN1A) and p27Kip1 (CDKN1B). Its activity is also regulated by its protein phosphorylation. Two alternatively spliced variants and multiple transcription initiation sites of this gene have been reported.

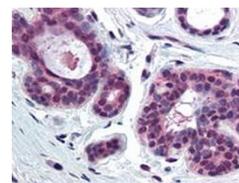
References:

1. Elledge, S.J., et al.: A new human p34 protein kinase CDK2, identified by complementation of a cdc28 mutation, is a homolog of *Xenopus* Eg1. *EMBO J.* 1991; 10; 2653-2659.
2. Giordano, A., et al.: Cell cycle regulation of histone H1 kinase activity associated with the adenoviral protein E1A. *Science.* 1991; 253; 1271-1275.
3. Tsai, L.H., et al.: Isolation of the human CDK2 gene that encodes the cyclin A and adenovirus E1A-associated p33 kinase. *Nature.* 1991; 353; 174-177.

Sample Data



Anti-CDK2 (1:200 dilution) used to detect cdk-2 in asynchronous HeLa cell lysates (run in duplicate). Cdk-2 is indicated as 32-33 kDa band. Note that multiple isoforms as well as phosphorylated forms of cdk-2 may be detected by this antibody.



Anti-CDK2 antibody was diluted 1:500 to detect CDK2 in human skin tissue. Tissue was formalin fixed and paraffin embedded. No pre-treatment of sample was required. The image shows the localization of antibody as the precipitated red signal, with a hematoxylin purple nuclear counter stain.

Anti-CDK2

Rabbit Polyclonal Antibody

Catalog #	C29-63R
Lot #	B3216-31
Purification	Delipidation and Defibrination
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°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
www.signalchem.com

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