

## Anti-Phospho-APC3 (Ser427)

Rabbit Polyclonal Antibody

**Catalog # A89-365BR**

Lot # B3216-29

### Cited Applications

ELISA, WB

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

### Specificity

Recognizes the human Cdc27 protein phosphorylated at Serine 427.

### Cross Reactivity

Human, Rat, Dog, Chicken and Chimpanzee

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

The antibody was produced against synthesized peptide corresponding to amino acids 422-430 of human Cdc27 protein.

### Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide

### Stability

1yr at -20°C from date of shipment

### Scientific Background

Cdc27 (also known as cell division cycle protein 27 homologue, CDC27Hs and H-NUC) shares strong similarity with *Saccharomyces cerevisiae* protein Cdc27, and the gene product of *Schizosaccharomyces pombe* nuc 2. Cdc27 is a component of anaphase-promoting complex (APC), which is composed of eight protein subunits and highly conserved in eukaryotic cells. APC catalyzes the formation of cyclin B-ubiquitin conjugate that is responsible for the ubiquitin-mediated proteolysis of B-type cyclins. This protein and 3 other members of the APC complex contain the TPR (tetratricopeptide repeat), a protein domain important for protein-protein interaction. This protein was shown to interact with mitotic checkpoint proteins including Mad2, p55CDC and BUBR1, and thus may be involved in controlling the timing of mitosis. Cdc27 has a nuclear localization.

### References:

1. Gabellini, D., et al.: Early mitotic degradation of the homeoprotein HOXC10 is potentially linked to cell cycle progression. *EMBO J.* 2003; 22 (14); 3715-3724.
2. Topper, L.M., et al. The dephosphorylated form of the anaphase-promoting complex protein Cdc27/Apc3 concentrates on kinetochores and chromosome arms in mitosis. *Cell Cycle.* 2002; 1 (4); 282-292.
3. Wassmann, K., et al.: Mad2 transiently associates with an APC/p55Cdc complex during mitosis. *Proc. Natl. Acad. Sci. U.S.A.* 1988; 95 (19); 11193-11198.

## Anti-Phospho-APC3 (Ser427)

Rabbit Polyclonal Antibody

|                    |  |
|--------------------|--|
| Catalog #          | A89-365BR  |
| Lot #              | B3216-29   |
| Purification       | Immunoaffinity chromatography  |
| 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](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.**