

Anti-APC3 (CDC27)

Rabbit Polyclonal Antibody

Catalog # **A89-363BR**

Lot # B3216-28

Cited Applications

ELISA, IF, WB

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

Specificity

Recognizes the human Cdc27 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

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

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

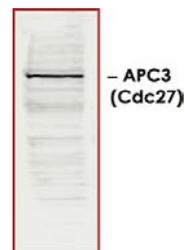
Scientific Background

Cdc27 (a.k.a. cell division cycle protein 27, CDC27Hs and H-NUC) is a component of anaphase-promoting complex (APC), which is composed of eight protein subunits. APC catalyzes the formation of cyclin B-ubiquitin conjugate that controls the ubiquitin-mediated proteolysis of B-type cyclins. Cdc27 contains a TPR (tetra-ricopeptide repeat) domain, which facilitates protein-protein interactions. This protein interacts with mitotic checkpoint proteins including Mad2, p55CDC and BUBR1, pointing to a role in mitotic timing.

References:

1. 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.
2. 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.

Sample Data



Western blot using Anti-APC3 (CDC27) antibody (1.0 µg/ml) shows detection of a band ~90 kDa corresponding to human cdc27 (arrowhead) in ~35 µg of HeLa whole cell lysate.

Anti-APC3 (CDC27)

Rabbit Polyclonal Antibody

Catalog #	A89-363BR
Lot #	B3216-28
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
www.signalchem.com

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