

Anti-Phospho-APC1 (Ser377)

Rabbit Polyclonal Antibody

Catalog # A87-365R

Lot # B3216-16

Cited Applications

ELISA, IHC, WB

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

Specificity

Recognizes the APC1 protein phosphorylated at Serine 377

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antibody was produced against synthesized peptide corresponding to amino acids 373-382 of human APC1 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

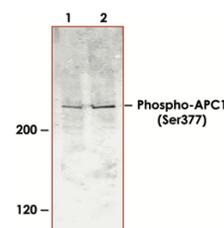
Scientific Background

APC1 (also known as Anaphase promoting complex subunit 1, Cyclosome subunit 1, Protein Tsg24, Mitotic checkpoint regulator and ANAPC1) is 1 of at least 11 subunits of the anaphase-promoting complex (APC), which functions at the metaphase-to-anaphase transition of the cell cycle and is regulated by spindle checkpoint proteins. The APC is an E3 ubiquitin ligase that targets cell cycle regulatory proteins for degradation by the proteasome, thereby allowing progression through the cell cycle.

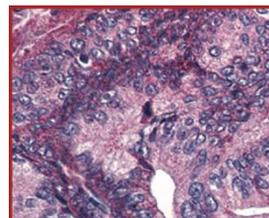
References:

1. Jorgensen, P.M., et al.: Characterization of the human APC1, the largest subunit of the anaphase-promoting complex. *Gene*. 2001;262 (1-2): 51-59.
2. Kraft C, et al.: Mitotic regulation of the human anaphase-promoting complex by phosphorylation. *Embo J*. 2003;22(24):6598-609.

Sample Data



Western blot using Anti-Phospho-APC1 (Ser377) (1:1,000) shows detection of a band ~215 kDa corresponding to phosphorylated human APC1 (arrowhead). Lane 1 shows lysate from asynchronous cells. Lane 2 shows lysate from cells treated with nocodazole. Each lane contains approximately 30 µg of HeLa whole cell lysates, separated by 4-8% SDS-PAGE followed by transfer to nitrocellulose.



Anti-Phospho-APC1 (Ser377) antibody was used at 5.0 µg/ml to detect signal in a variety of tissues including multi-human, multi-brain and multi-cancer slides at 60X. Tissue was formalin-fixed and paraffin embedded.

Anti-phospho-APC1 (Ser377)

Rabbit Polyclonal Antibody

Catalog #

A87-365R

Lot #

B3216-16

Purification

Immunoaffinity chromatography followed by solid phase adsorption.

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.