

Anti-APC1 (Ser355)

Rabbit Polyclonal Antibody

Catalog # A87-363BR

Lot # B3216-17

Cited Applications

ELISA, IP, WB

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

Specificity

Recognizes the human AC1 protein phosphorylated at Serine 355

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antibody was produced against synthesized peptide corresponding to amino acids 351-359 of Human APC1.

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

Anti-APC1 (Ser355)

Rabbit Polyclonal Antibody

| | |
|--------------------|--|
| Catalog # | A87-363BR |
| Lot # | B3216-17 |
| 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.