

## Anti-AKT1

Rabbit Polyclonal Antibody

**Catalog # A16-63CR**

Lot # O2109-41

### Cited Applications

WB, ELISA, IF, IHC

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

### Specificity

Recognizes the AKT1 protein

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

AKT1 antibody was raised against a 16 amino acid synthetic peptide from near the amino-terminus of human AKT1

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

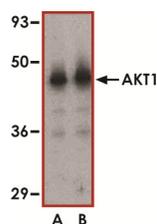
### Scientific Background

AKT1/PKBα is a serine/threonine kinase that belongs to the AKT family. AKT1 is activated in cells in response to diverse stimuli such as hormones, growth factors and extracellular matrix components and is involved in glucose metabolism, transcription, survival, cell proliferation, angiogenesis, and cell motility (1). AKT1 is frequently overexpressed and active in many types of human cancers including cancers of colon, breast, brain, pancreas and prostate as well as lymphomas and leukemias (2).

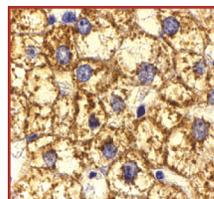
### References

1. Coffey, P.J. et al: Protein kinase B (c-Akt): a multifunctional mediator of phosphatidylinositol 3-kinase activation. *Biochem J.* 1998 Oct 1; 335 (Pt 1): 1-13.
2. Anderson, K.E. et al: Translocation of PDK-1 to the plasma membrane is important in allowing PDK-1 to activate protein kinase B. *Curr Biol.* 1998 Jun 4;8(12): 684-91.

### Sample Data



Western blot analysis of AKT1 in human liver cell lysate with AKT1 antibody at (A) 1 and (B) 2 µg/ml.



Immunohistochemistry of AKT1 in human liver cells with AKT1 antibody at 2 µg/ml.

## Anti-AKT1

Rabbit Polyclonal Antibody

Catalog Number

A16-63CR

Specific Lot Number

O2109-41

Purification

Affinity 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.**