

## Anti-AKT1

Mouse Monoclonal Antibody

Catalog # **A16-60M**

Lot # K3337-1

### Cited Applications

- Western blot (1:1,000)

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

### Specificity

Recognizes the AKT1 protein

### Cross Reactivity

- Western blot AKT1 from human and mouse cells

AKT1 from other species may also be detectable

### Host/Isotype/Clone#

Mouse, IgG1, 9A4

### Immunogen

Recombinant full length human AKT1 protein

### Formulation

PBS

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. 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.

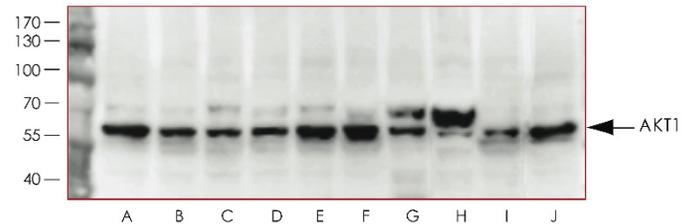
### Scientific Background

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. 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 (1).

### References

1. Anderson, KE. 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 with Anti-AKT1 (1:1,000) in (A) HEK293, (B) MDA-MB-231 kidney, (C) A549, (D) A2058, (E) K562, (F) Jurkat, (G) MCF7, (H) OVCAR-3, (I) PC3 and (J) THP1 cell lysates at 1 µg/mL.

## Anti-AKT1

Mouse Monoclonal Antibody

Catalog #	A16-60M
Lot #	K3337-1
Purification	Affinity purified using Protein G
Concentration	1.0 µg/µL
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.**