

Anti-AKT1

Mouse Monoclonal Antibody

Catalog # **A16-60M**

Lot # V058-2

Cited Applications

- Western blot (1:2500)

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

0.1M TRIS pH 8.0, 0.1M glycine

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

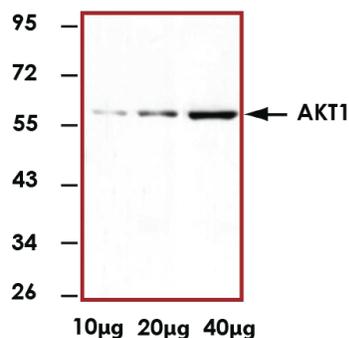
Scientific Background

AKT1/PKB α 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



Representative western blot with Anti-AKT1 (1:2500) using 10 µg, 20 µg, and 40 µg of HEK-293 cell lysate.

Anti-AKT1

Mouse Monoclonal Antibody

Catalog Number

A16-60M

Specific Lot Number

V058-2

Purification

Affinity purified using Protein G

Concentration

1.0 µg/µL

Stability

1 yr 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.