

Anti-PKCδ

Rabbit Polyclonal Antibody

Catalog # P64-63CR

Lot # M191-2

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the PKCδ protein

Cross Reactivity

- Western blot human and mouse PKCδ

PKCδ from other species may also be detectable

Host

Rabbit

Immunogen

Synthesized peptide derived from C-terminus of human PKCδ protein

Formulation

TBS, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) 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.

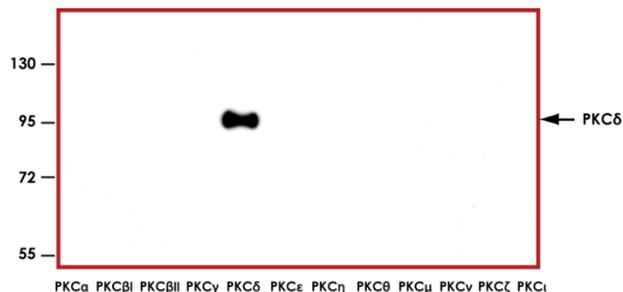
Scientific Background

Protein kinase C delta (PKCδ) is a member of the protein kinase C (PKC) family of serine-threonine kinases. It is a 79 kDa protein kinase that shows strict dependence on the presence of phospholipids, but shows no activation by Ca²⁺ (1). Phosphatidylinositol is the most potent activator of PKC delta. Northern blot analysis indicated that PKC delta is widely distributed in almost all the tissues and is a major isoform of PKC expressed in hemopoietic cells (2). PKC delta is involved in fundamental cellular functions regulated by diacylglycerols and mimicked by phorbol esters.

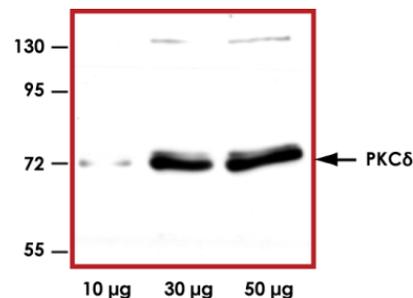
References

- Leibersperger, H. et al: Immunological demonstration of a calcium-unresponsive protein kinase C of the delta-type in different species and murine tissues. Predominance in epidermis. *J Biol Chem.* 1991 Aug 5; 266(22):14778-84.
- Mischak, H. et al: Mouse protein kinase C-delta, the major isoform expressed in mouse hemopoietic cells: sequence of the cDNA, expression patterns, and characterization of the protein. *Biochemistry.* 1991 Aug 13; 30(32):7925-31.

Sample Data



Representative Western blot with Anti- PKCδ (1:1000) using 15ng of GST-tagged human recombinant PKCα, PKCβI, PKCβII, PKCγ, PKCδ, PKCε, PKCη, PKCθ, PKCμ, PKCν, PKCξ, and PKCι.



Representative Western blot (1:1000) with PKCδ using 10 µg, 30µg, 50µg of 293 cell lysates.

Anti-PKCδ

Rabbit Polyclonal Antibody

Catalog Number

P64-63CR

Specific Lot Number

M191-2

Purification

Affinity chromatography

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
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.