

Anti-PKCδ

Rabbit Polyclonal Antibody

Catalog # P64-63R

Lot # H270-2

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 PKCδ

PKCδ from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 643-647 of human PKCδ

Formulation

PBS pH 7.4, 0.02% sodium azide, 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.

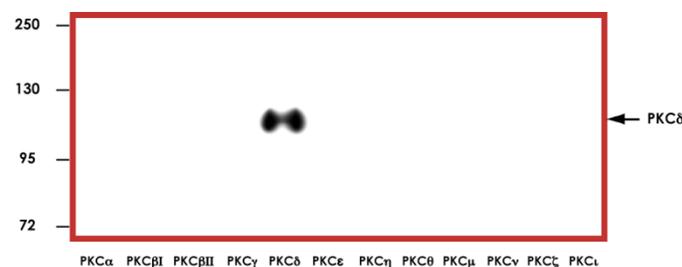
Scientific Background

Protein kinase C delta (PKCδ) is a member of the protein kinase C (PKC) family of serine-threonine kinases. It is a 79 kd 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 PKCdelta is widely distributed in almost all the tissues and is a major isoform of PKC expressed in hemopoietic cells (2). PKCdelta is involved in fundamental cellular functions regulated by diacylglycerols and mimicked by phorbol esters.

References

1. 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.

Sample Data



Representative western blot with Anti- PKCδ (1:1000) using 20ng of GST-tagged human recombinant PKCα, PKCβI, PKCβII, PKCγ, PKCδ, PKCε, PKCη, PKCθ, PKCμ, PKCν, PKCζ, and PKCι.

2. 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.

Anti-PKCδ

Rabbit Polyclonal Antibody

Catalog Number	P64-63R
Specific Lot Number	H270-2
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.