

Anti- PLC γ 1

Rabbit Polyclonal Antibody

Catalog # P40-63R

Lot # R133-7

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the PLC γ 1 protein

Cross Reactivity

- Western blot human PLC γ 1

PLC γ 1 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 769-773 of human PLC γ 1

Formulation

PBS, pH 7.4, 0.02% NaN₃, 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

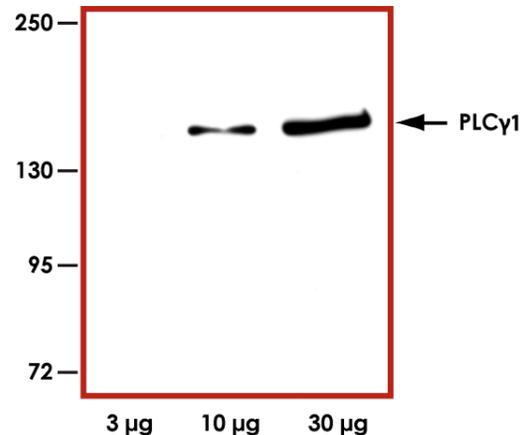
PLC γ 1 or phospholipase C gamma 1 plays an important role in the intracellular transduction of message from receptor-mediated tyrosine kinase activation. PLC γ 1 catalyzes the formation of two second-messenger molecules inositol 1, 4, 5-trisphosphate and diacylglycerol from phosphatidylinositol 4, 5-bisphosphate. PLC γ 1 is a major substrate for heparin-binding growth factor 1 activated tyrosine kinase. PLC γ 1 inhibits growth factor receptor-mediated invasion of breast and prostate tumor cells and is a physiological guanine nucleotide exchange factor for the nuclear GTPase PIKE (1).

References

1. Ye, K. et al: Phospholipase C-gamma-1 is a physiological guanine nucleotide exchange factor for the nuclear GTPase PIKE. Nature 415: 541-544, 2002.

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

Sample Data



Representative Western blot with Anti- PLC γ 1 (1:1000) using 3µg, 10µg and 30µg of Jurkat cell lysate.

Anti-PLC γ 1

Rabbit Polyclonal Antibody

Catalog Number

P40-63R

Specific Lot Number

R133-7

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.