

Echelon Biosciences Inc.

Rabbit Anti-Human DGK ζ Peptide Antibody

Z-DGKz

Support: echelon@echelon-inc.com

Description:

Rabbit polyclonal antibody against human diacylglycerol kinase zeta (DGK ζ)

Applications:

WB – 1:1000

Other in vitro and cellular applications are possible using this antibody, but have not been verified by Echelon Biosciences.

Properties:

Form – liquid

Storage instructions – Store at 4 °C for up to 30 days. Aliquot and store at -20 °C if longer storage is necessary. Antibody is stable for at least 2 years when stored at -20 °C. Avoid repeated freeze/thaw cycles.

Storage buffer – 0.01% thiomersal

Concentration – N/A

Purity – in serum

Immunogen – peptide = LENRQHYQMIQREDQE

Clonality – polyclonal

Isotype – IgG

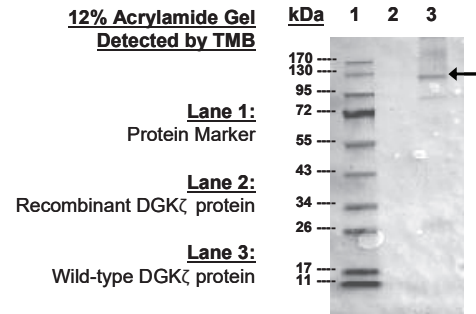
Specificity:

Shown to recognize a ~117kDa protein in wild-type DGK ζ protein with no cross activity to recombinant DGK α protein.

Background:

DGK catalyzes a reaction terminating the protein kinase C (PKC) mediated signaling by converting diacylglycerol (DAG) to phosphatidic acid (PA), a signaling lipid which is involved in DNA synthesis, phospholipid synthesis, and regulation of several enzymes. DAG is an intermediate in phospholipid synthesis which functions as an allosteric activator of PKC. Research shows the zeta isoform of DGK undergoes tissue-specific alternative splicing resulting in an elongated N terminus enzyme predominantly in muscle.

Data: Western Blot



Anti-DGK ζ peptide polyclonal antibody (1:1000)

References:

1. Luo B, Prescott SM, Topham MK. (2004) Diacylglycerol kinase zeta regulates phosphatidylinositol 4-phosphate 5-kinase I α by a novel mechanism. Cell Signal. 16(8):891-7.

Related Products:

Products	Catalog Number
Assays and Reagents	
LPA ELISA	K-2800s
Phosphatidic Acid Beads	P-B0PA
PI3K Inhibitor (LY294002)	B-0294
Lipids and Antibodies	
Biotin-phosphatidylglycerol	L-51B16
Anti-DGK α	Z-DGKa

Technical Data Sheet Rev. 5, 06-20-18 - For research use only. Not intended or approved for diagnostic or therapeutic use.



Echelon Biosciences Inc.

Ph: 866-588-0455

Fax: 801-588-0497

Echelon-inc.com