

Echelon Biosciences Inc.

Sheep Anti-Human DGK α Peptide Antibody

Z-DGK α

Support: echelon@echelon-inc.com

Description:

Sheep polyclonal antibody against diacylglycerol kinase α (DGK α)

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 = MAKERGLISPSDFAQL

Clonality – polyclonal

Isotype – IgG

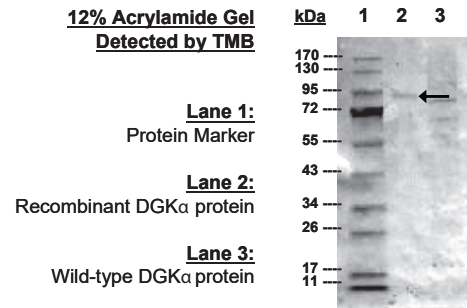
Specificity:

Shown to recognize an 80-85kDa recombinant DGK α protein with non-specific cross activity to wild-type DGK ζ 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. The alpha isoform is a type I DGKs, and has slightly different calcium-responsiveness.

Data: Western Blot



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

References:

1. Schaap D, de Widt J, van der Wal J, Vandekerckhove J, van Damme J, Gussow D, Ploegh HL, van Blitterswijk WJ, van der Bend RL. (1990) Purification, cDNA-cloning and expression of human diacylglycerol kinase. FEBS Lett. Nov 26;275(1-2):151-8.

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 ζ	Z-DGKz

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