

DGKG Polyclonal Antibody

(Catalog # A62406)

Background

Reverses the normal flow of glycerolipid biosynthesis by phosphorylating diacylglycerol back to phosphatidic acid.

Description

DGKG Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P49619

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Diacylglycerol kinase gamma protein (1-255AA)

Storage

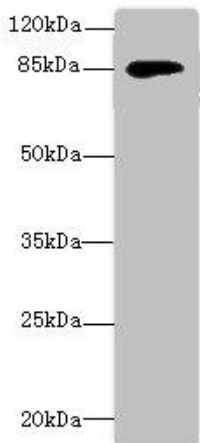
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Diacylglycerol kinase gamma, DAG kinase gamma, Diglyceride kinase gamma, DGK-gamma, DGKG, DAGK3

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: DGKG Polyclonal Antibody at
4ug/ml+Mouse brain tissue
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 90, 87, 85 kDa
Observed band size: 89 kDa