
PLCZ1 Polyclonal Antibody

(Catalog #A65778)

Background

The production of the second messenger molecules diacylglycerol (DAG) and inositol 1,4,5-trisphosphate (IP3) is mediated by activated phosphatidylinositol-specific phospholipase C enzymes. In vitro, hydrolyzes PtdIns(4,5)P2 in a Ca(2+)-dependent manner. Triggers intracellular Ca(2+) oscillations in oocytes solely during M phase and is involved in inducing oocyte activation and initiating embryonic development up to the blastocyst stage. Is therefore a strong candidate for the egg-activating soluble sperm factor that is transferred from the sperm into the egg cytoplasm following gamete membrane fusion. May exert an inhibitory effect on phospholipase-C-coupled processes that depend on calcium ions and protein kinase C, including CFTR trafficking and function.

Description

PLCZ1 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

Q86YW0

Purification

>95%, Protein G purified

Immunogen

Recombinant Human 1-phosphatidylinositol 4,5-bisphosphate phosphodiesterase zeta-1 protein (1-415AA)

Storage

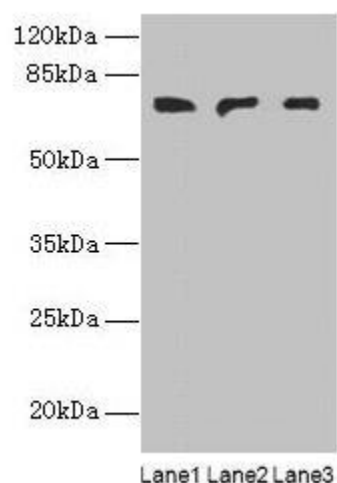
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

1-phosphatidylinositol 4, 5-bisphosphate phosphodiesterase zeta-1, Phosphoinositide phospholipase C-zeta-1, Phospholipase C-zeta-11, PLCZ1

Application

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



Western blot

All lanes: PLCZ1 antibody at 3ug/ml

Lane 1: Mouse liver tissue

Lane 2: Mouse kidney tissue

Lane 3: Human placenta tissue

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 71, 48, 58 kDa

Observed band size: 71 kDa