

PLCD4 Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog # P442-30DG

Lot # U1812-3

Product Description

Recombinant full-length human PLCD4 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_032726](#).

Gene Aliases

EIEE12; PI-PLC; PLC-154; PLC-I; PLC154; PLCD4A; PLCD4B

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

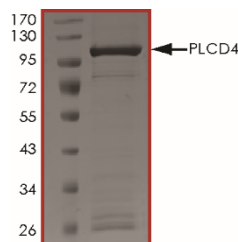
Scientific Background

1-phosphatidylinositol 4, 5-bisphosphate phosphodiesterase delta-4 (PLCD4) is one of the subtypes of phosphoinositide-specific phospholipase C (PI-PLC). Hydrolyzes the phosphatidylinositol 4, 5-bisphosphate (PIP2) to generate 2 second messenger molecules diacylglycerol (DAG) and inositol 1, 4, 5-trisphosphate (IP3). PLCD4 has been implicated to have a key role in cell proliferation. It was found that aberrant expression of PLCD4 may cause oncogenesis in certain tissues through upregulation of ErbB and activation of ERK pathway.

References

1. Leung DW, et al: Phospholipase C delta-4 overexpression upregulates ErbB1/2 expression, Erk signaling pathway, and proliferation in MCF-7 cells. Mol Cancer. 3:15, 2004.
2. Liu N, et al: A new phospholipase C delta 4 is induced at S-phase of the cell cycle and appears in the nucleus. J Biol Chem. 271(1): 355-60, 1996.

Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **110 kDa**.

PLCD4 Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog #	P442-30DG
Lot #	U1812-3
Purity	>90%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.