

## PLCB1 Protein

Full length human recombinant protein expressed in Sf9 cells

**Catalog # P441-30G**

Lot # U1800-4

### Product Description

Recombinant full-length human PLCB1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_015192](#).

### Gene Aliases

EIEE12; PI-PLC; PLC-154; PLC-I; PLC154; PLCB1A; PLCB1B

### 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.

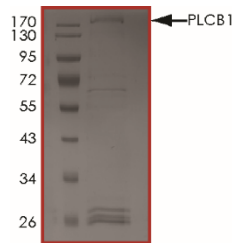
### Scientific Background

1-phosphatidylinositol 4, 5-bisphosphate phosphodiesterase beta-1 (PLCB1) is one of the best characterized isoenzymes in the inositide specific phospholipase C (PI-PLC) family. Responsible for the hydrolysis of the phosphatidylino-4, 5-bisphosphate (PIP<sub>2</sub>), generating the second messengers inositol-1, 4, 5-triphosphate and diacylglycerol. It is a key molecule for nuclear inositide signaling, where it plays a role in cell cycle progression, proliferation and differentiation.

### References

1. Faenza I, et al: A role for PLCbeta1 in myotonic dystrophies type 1 and 2. *Faseb J.* 26:3042-3048, 2012.
2. Faenza I, et al: Nuclear inositide specific phospholipase C signalling - interactions and activity. *FEBS J.* 280(24):6311-21, 2013.

### Purity



The purity was determined to be **>70%** by densitometry. Approx. MW **180 kDa**.

## PLCB1 Protein

Full length human recombinant protein expressed in Sf9 cells

|                    |   |
|--------------------|---|
| Catalog #          | P441-30G  |
| Lot #              | U1800-4   |
| Purity             | >70%  |
| Concentration      | 0.05 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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