

PIP5K3 Protein

Recombinant protein expressed in Sf9 cells

Catalog # P17-31G

Lot # X429-2

Product Description

Recombinant human PIP5K3 (1493-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_015040](#).

Gene Aliases

PIKFYVE; CFD; FAB1; PIP5K; FLJ37746; KIAA0981; MGC40423

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

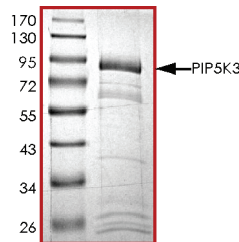
Scientific Background

PIP5K3 (also known as phosphatidylinositol-3-phosphate 5-kinase type III or PIPKIII) encodes an enzyme that phosphorylates the D-5 position in PtdIns and phosphatidylinositol-3-phosphate (PtdIns3P) to make PtdIns5P and PtdIns(3,5)biphosphate. PIP5K3 belongs to a large family of lipid kinases that alter the phosphorylation status of intracellular phosphatidylinositol but also have protein kinase activity (1). PIP5K3 regulate cytoskeletal functions, membrane trafficking, and receptor signaling by recruiting protein complexes to cell- and endosomal-membranes. The inhibition of PIP5K3 renders the late endosome/lysosome compartment refractory to fusion with autophagosomes and multivesicular bodies (2).

References

1. Shisheva, A..et.al: Cloning, characterization, and expression of a novel Zn(2+)-binding FYVE finger-containing phosphoinositide kinase in insulin-sensitive cells. *Molec. Cell Biol.* 19: 623-634, 1999.
2. de Lartigue, J.et.al: PIKfyve regulation of endosome-linked pathways. *Traffic* 10: 883-893, 2009.

Purity



The purity of PIP5K3 was determined to be **>75%** by densitometry.
Approx. MW **93kDa**.

PIP5K3 Protein

Recombinant protein expressed in Sf9 cells

Catalog #	P17-31G
Lot #	X429-2
Purity	>75%
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.