

# Phosphoinositide Kinases

## Key References

- Alessi, D.R., Downes, C.P. "The role of PI 3-kinase in insulin action." *Biochim. Biophys. Acta* **1436**, 151-164 (1998).
- Anderson, R.A. et al. "Phosphatidylinositol phosphate kinases, a multifaceted family of signaling enzymes." *J. Biol. Chem.* **274**, 9907-9910 (1999).
- Ciruela, A. et al. "Nuclear targeting of the beta isoform of type II phosphatidylinositol phosphate kinase (phosphatidylinositol 5-phosphate 4-kinase) by its alpha-helix 7." *Biochem. J.* **346**, 587-591 (2000).
- De Camilli, P. et al. "Phosphoinositides as regulators in membrane traffic." *Science* **271**, 1533-1539 (1996).
- Gehrmann, T., Heilmeyer, L.M. "Phosphatidylinositol 4-kinases." *Eur. J. Biochem.* **253**, 357-370 (1998).
- Hinchliffe, K.A. et al. "PIPKins1, their substrates and their products: new functions for old enzymes." *Biochim. Biophys. Acta* **1436**, 87-104 (1998).
- Martin, T.F.J. "Phosphoinositide lipids as signalling molecules: Common themes for signal transduction, cytoskeletal regulation and membrane trafficking." *Annu. Rev. Cell Dev. Biol.* **14**, 231-264 (1998).
- McEwen, R.K. et al. "Complementation analysis in PtdInsP kinase-deficient yeast mutants demonstrates that *S. pombe* and murine Fab1p homologues are phosphatidylinositol 3-phosphate 5-kinases." *J. Biol. Chem.* **274**, 33905-33912 (1999).
- Pacold, M.E. et al. "Crystal structure and functional analysis of ras binding to its effector phosphoinositide 3-kinase gamma." *Cell* **103**, 931-943 (2000).
- Rao, V.D. et al. "Structure of type II beta phosphatidylinositol phosphate kinase: A protein kinase fold flattened for interfacial phosphorylation." *Cell* **94**, 829-839 (1998).
- Tolias, K.F., Cantley, L.C. "Pathways for phosphoinositide synthesis." *Chem. Phys. Lipids* **98**, 69-77 (1999).
- Vanhaesebroeck, B., Waterfield, M.D. "Signaling by distinct classes of phosphoinositide 3-kinases." *Exptl. Cell Res.* **253**, 239-54 (1999).

## Overview

The membranes of mammalian cells contain substantial quantities of phosphatidylinositol (PtdIns) and small amounts of at least seven phosphorylated PtdIns derivatives (PPI). Phosphoinositide kinases synthesize PPI by adding phosphate groups to pre-existing inositol glycerophospholipids (see scheme below).

The kinase domains of the two subfamilies of PtdIns 4-kinases and all of the phosphoinositide 3-kinase (PI3K) subfamilies are closely related to each other and also to protein kinases. However, the three types of phosphatidylinositol phosphate (PIP) kinases (PIPKin), kinases which donate a phosphate to one of the three isomeric PtdInsPs, constitute a kinase family that shares some kinase domain motifs with other kinases, but are otherwise unique.

Four of the kinases are responsible for biosynthesis of PtdIns(4,5) $P_2$ , the key phosphoinositide that is the shared substrate of two widespread receptor-activated signaling pathways: the phospholipase C-catalyzed formation of Ins(1,4,5) $P_3$  and 1,2-diacylglycerol (DAG); and synthesis by Type I PI3Ks of the membrane-associated messenger molecule PtdIns(3,4,5) $P_3$ . PtdIns(4,5) $P_2$  is also essential for a substantial number of other cell functions, including actin-based motility, secretory vesicle trafficking, exocytosis and phospholipase D activation.

PtdIns(4,5) $P_2$  is synthesized, via PtdIns4P, in the plasma membrane, secretory vesicles, Golgi and nuclei, and probably elsewhere. Why cells use an uncharacterized PtdIns 5-kinase and a PIPkin-II to make a small proportion of cellular PtdIns(4,5) $P_2$  by an alternative pathway, via PtdIns5P, is not

known. Some PtdIns(4,5) $P_2$  is made and undoubtedly functions in the nucleus, but in what manner is not known. There is evidence for partial nuclear localization of PIPkin-I and of PIPkin-II.

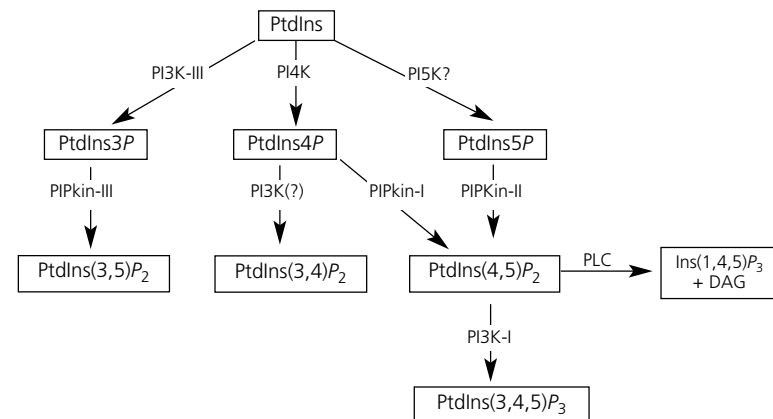
The individual PI3K subfamilies phosphorylate different phosphoinositides. Type I enzymes phosphorylate PtdIns(4,5) $P_2$  to PtdIns(3,4,5) $P_3$  and thereby transmit signals from numerous receptor tyrosine kinases (PI3K-IA: activated by the insulin receptor and many growth factor receptors) and from some G protein-coupled receptors (PI3K-IB: particularly in hemopoietic cells, by receptors that influence motility and the bactericidal oxidative burst of neutrophils).

PtdIns(3,4,5) $P_3$  is a remarkable membrane-associated second messenger molecule. It appears to have many direct target proteins, each of which it interacts with through highly PtdIns(3,4,5) $P_3$ -selective pleckstrin homology (PH) domains. Some of these are protein kinases (e.g. PDK1, PKB, Btk, Itk), while others include regula-

tors of small GTPases (GTP/GDP exchange factors and GTPase-activating factors). The PtdIns(3,4,5) $P_3$ -activated protein kinases have numerous substrates, and sometimes initiate downstream protein kinase cascades.

Type III PI3Ks convert PtdIns to PtdIns3P. This lipid is essential for vesicle trafficking from the trans-Golgi to vacuole/lysosome compartments. It appears to exert its effects through interactions with target proteins that include PtdIns3P-specific FYVE domains.

The functions of Type II PI3Ks remain uncertain. In addition, there is a diverse group of 'Type IV PI3K-related kinases' – comprising Tor (target of rapamycin) proteins, DNA-dependent protein kinase (DNA-PK) and ATM (Ataxia Telangiectasia Mutated). These have kinase domains that are related to those of *bona fide* PI3Ks, but are probably protein kinases rather than phosphoinositide kinases.



## Phosphoinositide Kinases

| KINASE                               | PtdIns 4-kinase (PI4K)   | PtdIns4P 5-kinase (PIPkin-I)   | PtdIns5P 4-kinase (PIPkin-II)  | PtdIns3P 5-kinase (PIPkin-III)  |
|--------------------------------------|--|--|--|---|
| <b>REACTION <i>In Vivo</i></b>       | PtdIns ( <a href="#">P 8443</a> , <a href="#">P 2517</a> ) > PtdIns4P ( <a href="#">P 9638</a> )                                     | PtdIns4P ( <a href="#">P 9638</a> ) > PtdIns(4,5)P <sub>2</sub> ( <a href="#">P 9763</a> ) | PtdIns5P > PtdIns(4,5)P <sub>2</sub> ( <a href="#">P 9763</a> )                          | PtdIns3P > PtdIns(4,5)P <sub>2</sub> ( <a href="#">P 9763</a> )                       |
| <b>STRUCTURE</b>                     | Two subfamilies:<br>I: ~100 kDa – prototype is yeast Pik1p<br>II: ~200 kDa (some have SH3 and PH domains) – prototype is yeast Stt2p | ~60-90 kDa: multiple forms<br>- prototype is yeast Mss4p                                   | ~45 kDa: multiple forms  | ~200-250 kDa: includes FYVE, CCP1-chaperone-like domains<br>-prototype is yeast Fab1p |
| <b>LOCALIZATION/CONTROL/COMMENTS</b> | Some types may be regulated by phosphorylation   | Major route for PtdIns(4,5)P <sub>2</sub> synthesis  | Minor route for PtdIns(4,5)P <sub>2</sub> synthesis. May be regulated by phosphorylation | Hyperosmotic stress in yeast, unknown in animal cells                                 |
| <b>INHIBITORS</b>                    | Adenosine ( <a href="#">A 9251</a> )<br>Wortmannin for some forms ( <a href="#">W 1628</a> )   | Stimulated by phosphatidate ( <a href="#">P 9511</a> )                                     | Heparin ( <a href="#">H 3393 (p)</a> , <a href="#">H 0777 (b)</a> )                      |   |

## Phosphoinositide Kinases (continued)

| KINASE                               | Type IA PI3K  | Type IB PI3K   | Type II PI3K  | Type III PI3K   |
|--------------------------------------|---|--|---|---|
| <b>REACTION <i>In Vivo</i></b>       | PtdIns(4,5)P <sub>2</sub> ( <a href="#">P 9763</a> ) > PtdIns(3,4,5)P <sub>3</sub> (also has protein kinase activity) | PtdIns(4,5)P <sub>2</sub> ( <a href="#">P 9763</a> ) > PtdIns(3,4,5)P <sub>3</sub>           | PtdIns4P ( <a href="#">P 9638</a> ) > PtdIns(3,4)P <sub>2</sub> | PtdIns ( <a href="#">P 8443</a> , <a href="#">P 2517</a> ) > PtdIns3P |
| <b>STRUCTURE</b>                     | p85(α,β) or p55γ regulatory subunit:<br>p110 (α, β or δ) catalytic subunit  | p101 regulatory subunit:<br>p110γ catalytic subunit  | >170 kDa, C-terminal C2 domain                                  | ~80-110 kDa<br>- prototype is yeast Vps34p                            |
| <b>LOCALIZATION/CONTROL/COMMENTS</b> | p85 SH2 domain interacts with P-Tyr residues. p110 activated by p85* and by activated Ras                             | p101 activated by βγ-complexes liberated from activated G <sub>i</sub> and/or G <sub>o</sub> | May be activated by some tyrosine kinases                       | Putative 150 kDa regulatory partner in mammalian cells                |
| <b>INHIBITORS</b>                    | Wortmannin ( <a href="#">W 1628</a> )<br>LY-294002 ( <a href="#">L 9908</a> )<br>Dominant-negative p85 constructs     | Wortmannin ( <a href="#">W 1628</a> )<br>LY-294002 ( <a href="#">L 9908</a> )                |   |   |

### ABBREVIATIONS

**PI3K:** phosphoinositide 3-kinase

**PIK:** PtdIns kinase

**PIP:** Phosphatidylinositol phosphate

**PIPkin:** PtdInsP kinase

**PPI:** Polyphosphoinositides

**PtdIns:** Phosphatidylinositol

**PtdIns3P:** PtdIns 3-phosphate

**PtdIns4P:** PtdIns 4-phosphate

**PtdIns5P:** PtdIns 5-phosphate

**PtdIns(3,4)P<sub>2</sub>:** PtdIns 3,4-bisphosphate

**PtdIns(3,5)P<sub>2</sub>:** PtdIns 3,5-bisphosphate

**PtdIns(4,5)P<sub>2</sub>:** PtdIns 4,5-bisphosphate

**PtdIns(3,4,5)P<sub>3</sub>:** PtdIns 3,4,5-trisphosphate

**Ins(1,4,5)P<sub>3</sub>:** Inositol 1,4,5-trisphosphate