

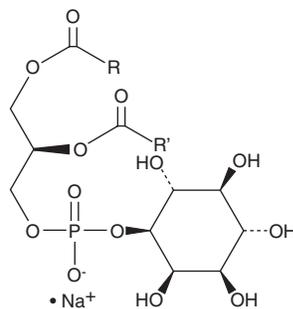
# PRODUCT INFORMATION



## Phosphatidylinositol (plant, wheat germ) (sodium salt)

Item No. 33671

**CAS Registry No.:** 84776-78-3  
**Formal Name:** (R)-2,3-bis(((9Z,12Z)-octadeca-9,12-dienoyl)oxy)propyl ((1S,2R,3R,4S,5S,6R)-2,3,4,5,6-pentahydroxycyclohexyl) phosphate  
**MF:** C<sub>45</sub>H<sub>78</sub>O<sub>13</sub>P • Na (for linoleoyl)  
**FW:** 881.1  
**Purity:** ≥98%  
**Supplied as:** A solution in chloroform  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Phosphatidylinositol (plant, wheat germ) (sodium salt) is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice. Phosphatidylinositol (plant, wheat germ) (sodium salt) is soluble in ethyl ether.

### Description

Phosphatidylinositols are glycerophospholipids that contain a glycerol backbone, two non-polar fatty acid tails, and a polar inositol head group.<sup>1,2</sup> They are synthesized from cytidine diphosphate diacylglycerol (CPD-DAG) and myo-inositol by phosphoinositide synthase and represent approximately 10% of total cellular phospholipids. Phosphatidylinositols can be phosphorylated on their inositol rings to produce phosphoinositides, which have been implicated in calcium regulation, vesicle trafficking, mitogenesis, cell survival, and rearrangement of actin. This product contains phosphatidylinositol molecular species with primarily C16:0 and C18:0 fatty acyl chain lengths.

### References

1. De Craene, J.-O., Bertazzi, D.L., Bär, S., *et al.* Phosphoinositides, major actors in membrane trafficking and lipid signaling pathways. *Int. J. Mol. Sci.* **18(3)**, E634 (2017).
2. Toliás, K.F. and Cantley, L.C. Pathways for phosphoinositide synthesis. *Chem. Phys. Lipids* **98(1-2)**, 69-77 (1999).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/16/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM