

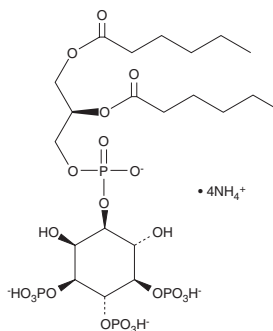
# PRODUCT INFORMATION



## PtdIns-(3,4,5)-P<sub>3</sub> (1,2-dihexanoyl) (ammonium salt)

Item No. 10008390

**CAS Registry No.:** 799268-62-5  
**Formal Name:** 1-(1,2R-dihexanoylphosphatidyl)inositol-3,4,5-triphosphate, tetraammonium salt  
**Synonyms:** DHPI-3,4,5-P<sub>3</sub>, Phosphatidylinositol-3,4,5-triphosphate C-6, PIP3C-16  
**MF:** C<sub>21</sub>H<sub>42</sub>O<sub>22</sub>P<sub>4</sub> • 4NH<sub>3</sub>  
**FW:** 838.6  
**Purity:** ≥98%  
**Supplied as:** A lyophilized powder  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

PtdIns-(3,4,5)-P<sub>3</sub> (1,2-dihexanoyl) (ammonium salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the PtdIns-(3,4,5)-P<sub>3</sub> (1,2-dihexanoyl) (ammonium salt) in water. The solubility of PtdIns-(3,4,5)-P<sub>3</sub> (1,2-dihexanoyl) (ammonium salt) in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

The phosphatidylinositol (PtdIns) phosphates represent a small percentage of total membrane phospholipids. However, they play a critical role in the generation and transmission of cellular signals.<sup>1,2</sup> PtdIns-(3,4,5)-P<sub>3</sub>, also known as PIP<sub>3</sub>, is resistant to cleavage by PI-specific phospholipase C (PLC).<sup>3</sup> Thus, it is likely to function in signal transduction as a modulator in its own right, rather than as a source of inositol tetraphosphates. PIP<sub>3</sub> can serve as an anchor for the binding of signal transduction proteins bearing pleckstrin homology (PH) domains.<sup>4,5</sup> Protein binding to PIP<sub>3</sub> is important for cytoskeletal rearrangement and membrane trafficking.<sup>6</sup> PtdIns-(3,4,5)-P<sub>3</sub> (1,2-dihexanoyl) is a synthetic analog of natural PIP<sub>3</sub> with C<sub>6</sub>:0 fatty acids at the *sn*-1 and *sn*-2 positions. The compound features the same inositol and diacylglycerol (DAG) stereochemistry as that of the natural compound. The short fatty acid chains of this analog give it different physical properties from naturally-occurring PIP<sub>3</sub>, including higher solubility in aqueous media.

### References

1. Guan, X. and Wenk, M.R. *Front. Biosci.* **13(9)**, 3239-3251 (2008).
2. Kashiwada, M., Lu, P., and Rothman, P.B. *Immunol. Res.* **39(1-3)**, 194-224 (2007).
3. Serunian, L.A., Haber, M.T., Fukui, T., et al. *J. Biol. Chem.* **264(30)**, 17809-17815 (1989).
4. Tanaka, K., Imajoh-Ohmi, S., Sawada, T., et al. *Eur. J. Biochem.* **245(2)**, 512-519 (1997).
5. Yang, X., Rudolf, M., Carew, M.A., et al. *J. Biol. Chem.* **274(27)**, 18973-18980 (1999).
6. Hamaguchi, N., Ihara, S., Ohdaira, T., et al. *Biochem. Biophys. Res. Commun.* **361(2)**, 270-275 (2007).

#### 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, 08/06/2021

#### 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