

Product Information



PtdIns-(4,5)-P₂ (1,2-dipalmitoyl)-d₆₂ (sodium salt)

Item No. 10005615

Formal Name: 1-(1,2-dihexadecanoyl(2,2',3,3',4,4',5,5',6,6',7,7',8,8',9,9',10,10',11,11',12,12',13,13',14,14',15,16,16-d₃₁)phosphatidyl)inositol-4,5-bisphosphate, trisodium salt

Synonyms: DPPI-4,5-P₂-d₆₂, Phosphatidylinositol-4,5-diphosphate C16-d₆₂

MF: C₄₁H₁₆D₆₂O₁₉P₃ • 3Na

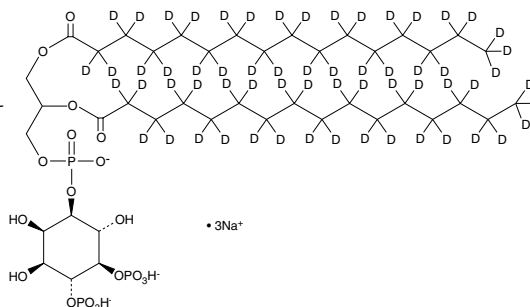
FW: 1,099.3

Chemical Purity: ≥98% PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt)

Deuterium Incorporation: ≥99% deuterated forms (d₁-d₆₂); ≤1% d₀

Stability: ≥1 year at -20°C

Supplied as: A lyophilized powder



Laboratory Procedures

PtdIns-(4,5)-P₂ (1,2-dipalmitoyl)-d₆₂ (sodium salt) contains 62 deuterium atoms at the 2, 2', 3, 3', 4, 4', 5, 5', 6, 6', 7, 7', 8, 8', 9, 9', 10, 10', 11, 11', 12, 12', 13, 13', 14, 14', 15, 15', 16, 16, and 16 positions of each fatty acyl chain. It is intended for use as an internal standard for the quantification of PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt) by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt) be stored as supplied at -20°C. It should be stable for at least one year.

PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt) in an organic solvent purged with an inert gas. PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt) is soluble in organic solvent such as chloroform:methanol:water (4:3:1). The solubility of PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt) in this solvent is at least 1 mg/ml. PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt) will not be stable in aqueous solutions for more than 24 hours.

PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt) is used as an internal standard for the quantification of PtdIns-(4,5)-P₂ (1,2-dipalmitoyl) (sodium salt) by stable isotope dilution MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

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.^{1,2} PtdIns-4,5-P₂-(1,2-dipalmitoyl)-d₆₂ is a synthetic analog of natural PtdIns containing deuterated C16:0 fatty acids at the *sn*-1 and *sn*-2 positions. This synthetic standard features the same inositol and diacyl glycerol (DAG) stereochemistry as that of the natural compound. The natural compound is the product of phosphatidylinositol 4-phosphate 5-kinase acting on PtdIns-(4)-P₁. Hydrolysis of PtdIns-(4,5)-P₂ by phosphoinositide (PI)-specific phospholipase C generates inositol triphosphate (IP₃) and DAG which are key second messengers in an intricate biochemical signal transduction cascade.

References

1. Exton, J.H. *Annu. Rev. Pharmacol. Toxicol.* **36**, 481-509 (1996).
2. Majerus, P.W. *Annu. Rev. Biochem.* **61**, 225-250 (1992).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10005615

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 02/14/2012

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com