

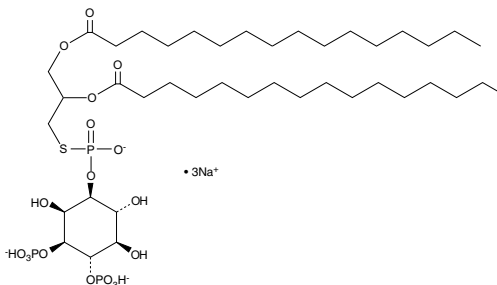
Product Information



Ptd(S)Ins-(3,4)-P₂ (1,2-dipalmitoyl) (sodium salt)

Item No. 10010112

Formal Name: 1-(1,2R-dihexadecanoylphosphothiotidyl) inositol-3,4-bisphosphate, trisodium salt
Synonym: DOsPI-3,4-P₂
MF: C₄₁H₇₈O₁₈P₃S • 3Na
FW: 1,053.0
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A lyophilized powder



Laboratory Procedures

For long term storage, we suggest that Ptd(S)Ins-(3,4)-P₂ (1,2-dipalmitoyl) (sodium salt) be stored as supplied at -20°C. It should be stable for at least one year.

Ptd(S)Ins-(3,4)-P₂ (1,2-dipalmitoyl) (sodium salt) is supplied as a lyophilized powder. Ptd(S)Ins-(3,4)-P₂ (1,2-dipalmitoyl) (sodium salt) is sparingly soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of Ptd(S)Ins-(3,4)-P₂ (1,2-dipalmitoyl) (sodium salt) be prepared by directly dissolving the lyophilized powder in aqueous buffers. The solubility of Ptd(S)Ins-(3,4)-P₂ (1,2-dipalmitoyl) (sodium salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solutions for more than one day.

Ptd(S)Ins-(3,4)-P₂ (1,2-dipalmitoyl) is an analog of naturally occurring PtdIns-(3,4)-P₂ that contains sulfur rather than oxygen at the *sn*-3 position of the glycerol backbone. Potential hydrolysis by phospholipase C (PLC) would yield a free thiol which could then react with chromogenic reagents such as DTNB (Ellman's reagent). There are no published reports on the biological activity of Ptd(S)Ins-(3,4)-P₂ (1,2-dipalmitoyl) or its use as a chromogenic substrate for PLC.

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10010112

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

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