

PRODUCT INFORMATION



PtdIns-(4,5)-P₂ (1,2-dioctanoyl)

Item No. 39270

CAS Registry No.: 204858-53-7

Formal Name: 1-[(2R)-2,3-bis[(1-oxooctyl)oxy]propyl hydrogen phosphate] 4,5-bis(dihydrogen phosphate) D-myo-inositol

Synonyms: DOPI-4,5-P₂, L-α-Phosphatidyl-D-myo-inositol 4,5-diphosphate, dioctanoyl, PI(4,5)P₂ (8:0/8:0), PIP2[4',5'] (8:0/8:0)

MF: C₂₅H₄₉O₁₉P₃

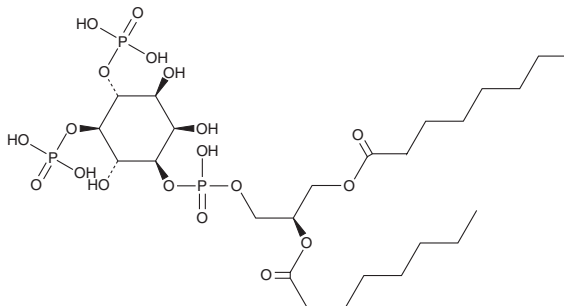
FW: 746.6

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

PtdIns-(4,5)-P₂ (1,2-dioctanoyl) is supplied as a solid. A stock solution may be made by dissolving the PtdIns-(4,5)-P₂ (1,2-dioctanoyl) in water. We do not recommend storing the aqueous solution for more than one day.

Description

PtdIns-(4,5)-P₂ (1,2-dioctanoyl) is a synthetic derivative of phosphatidylinositol.¹ It has been used to study the substrate specificity profile of inositol polyphosphate-5-phosphatase A (5ptase).

Reference

1. Bru, M., Kotkar, S.P., Kar, N., *et al.* Development of a solid phase synthesis strategy for soluble phosphoinositide analogues. *Chem. Sci.* **3**(6), 1893-1902 (2012).

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, 09/21/2023

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