

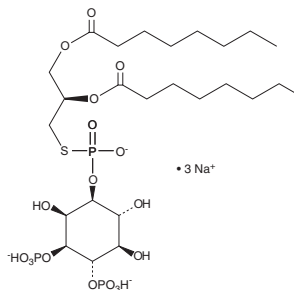
PRODUCT INFORMATION



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

Item No. 10010181

Formal Name: 1-(1-2R-phosphothiotidyl)inositol-3,4-bisphosphate, trisodium salt
Synonym: DOsPI-3,4-P₂
MF: C₂₅H₄₆O₁₈P₃S • 3Na
FW: 828.6
Purity: ≥95%
Supplied as: A lyophilized powder
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Ptd(S)Ins-(3,4)-P₂ (1,2-dioctanoyl) (sodium salt) is supplied as a lyophilized powder. Ptd(S)Ins-(3,4)-P₂ (1,2-dioctanoyl) (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-dioctanoyl) (sodium salt) be prepared by directly dissolving the lyophilized powder in aqueous buffers. The solubility of Ptd(S)Ins-(3,4)-P₂ (1,2-dioctanoyl) (sodium salt) in water is approximately 10 mg/ml. We do not recommend storing the aqueous solutions for more than one day.

Description

Ptd(S)Ins-(3,4)-P₂ (1,2-dioctanoyl) 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-P₂ (1,2-dioctanoyl) or its use as a chromogenic substrate for PLC.

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, 11/14/2017

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