

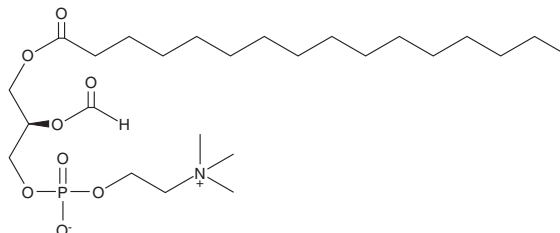
PRODUCT INFORMATION



1-Palmitoyl-2-formyl PC

Item No. 9002943

CAS Registry No.: 2194571-25-8
Formal Name: (R)-2-(formyloxy)-3-(palmitoyloxy)propyl
(2-(trimethylammonio)ethyl) phosphate
Synonyms: 1-Palmitoyl-2-formyl
Phosphatidylcholine, PFPC
MF: C₂₅H₅₀NO₈P
FW: 523.6
Purity: ≥95%
UV/Vis.: λ_{max}: 278 nm
Supplied as: A crystalline 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

1-Palmitoyl-2-formyl PC is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-palmitoyl-2-formyl PC in the solvent of choice. 1-Palmitoyl-2-formyl PC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 1-palmitoyl-2-formyl PC in ethanol is approximately 30 mg/ml and approximately 0.25 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 1-palmitoyl-2-formyl PC can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 1-palmitoyl-2-formyl PC in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

1-Palmitoyl-2-formyl PC is a formate-containing lysophospholipid.¹ Its biological significance is unknown.

Reference

1. Owen, J.S., Wykle, R.L., Samuel, M.P., *et al.* An improved assay for platelet-activating factor using HPLC-tandem mass spectrometry. *J. Lipid. Res.* **46**(2), 373-382 (2005).

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, 05/01/2024

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