

PRODUCT INFORMATION



POV-PC

Item No. 10031

CAS Registry No.: 121324-31-0

Formal Name: 1-palmitoyl-2-(5-oxovaleroyl)-sn-glycero-3-phosphatidylcholine

Synonym: 2-(5-oxovaleryl) Phosphatidylcholine

MF: C₂₉H₅₆NO₉P

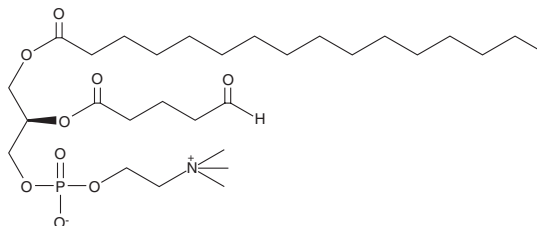
FW: 593.7

Purity: ≥98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: As supplied, 1 year from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

POV-PC is supplied as a solution in ethanol. To change the solvent, simply evaporate the POV-PC under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of POV-PC in ethanol is approximately 30 mg/ml and approximately 10 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. If an organic solvent-free solution of POV-PC is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of POV-PC in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Oxidized low-density lipoprotein (oxLDL) particles contain low molecular weight species which are cytotoxic and pro-atherogenic.¹ Many of these substances were recently isolated and purified from oxLDL, and identified as phosphatidylcholine species containing a fragmented, oxidized short-chain fatty acid remnant at the sn-2 position.² 1-(Palmitoyl)-2-(5-oxovaleroyl)-phosphatidylcholine, or POV-PC, is one of the oxLDL species derived from 2-arachidonoyl or eicosapentanoyl phospholipids.³ POV-PC confers CD36 scavenger receptor binding affinity more potently than any hydroperoxy PC species, and may be one of the more important structural determinants of oxLDL. Treatment of cultured endothelial cells with POV-PC stimulates monocyte binding, stimulates intracellular cAMP production, and strongly inhibits the LPS-induced binding of neutrophils.⁴

References

1. Podrez, E.A., Febbraio, M., Sheibani, N., *et al.* *J. Clin. Invest.* **105(8)**, 1095-1108 (2000).
2. Podrez, E.A., Batyreva, E., Shen, Z., *et al.* *J. Biol. Chem.* **277(41)**, 38517-38523 (2002).
3. Podrez, E.A., Batyreva, B., Shen, Z., *et al.* *J. Biol. Chem.* **277(41)**, 38503-38516 (2002).
4. Leitinger, N., Tyner, T.R., Oslund, L., *et al.* *Proc. Natl. Acad. Sci. USA* **96(21)**, 12010-12015 (1999).

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, 03/06/2017

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