

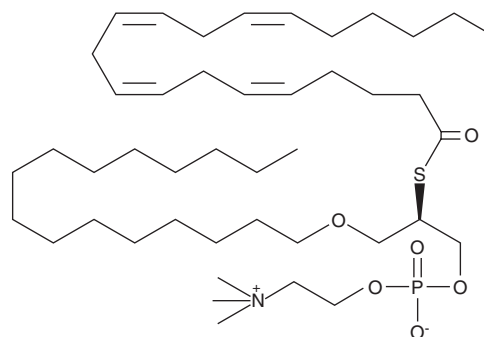
PRODUCT INFORMATION



Arachidonoyl thio-PC

Item No. 62240

CAS Registry No.: 146797-82-2
Formal Name: 1-O-hexadecyl-2-deoxy-2-thio-R-(5Z,8Z,11Z,14Z-eicosatetraenoyl)-sn-glycerol-3-phosphorylcholine
Synonym: 2-deoxy-2-thio Arachidonoyl PC
MF: C₄₄H₈₂NO₆PS
FW: 784.2
Purity: ≥98%
Supplied as: A solution in ethanol containing 0.1% BHT
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Arachidonoyl thio-PC is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of arachidonoyl thio-PC in these solvents is approximately 25 mg/ml. Arachidonoyl thio-PC is soluble in water as a concentration of approximately 0.5 mg/ml.

Description

Arachidonoyl thio-PC is a substrate for many phospholipase A₂s (PLA₂s) including sPLA₂, cPLA₂, and iPLA₂. Cleavage of the sn-2 fatty acid by PLA₂ results in generation of a free thiol which reacts with chromogenic reagents such as DTNB (Ellman's reagent) and DTP to allow quantitation of PLA₂ activity.¹ Isozyme-specific cPLA₂ activity can be measured by excluding or inhibiting sPLA₂ and iPLA₂ activities in the assay.

Reference

1. Yu, L. and Dennis, E.A. Thio-based phospholipase assay. *Methods Enzymol.* **197**, 65-75 (1991).

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, 02/15/2024

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