

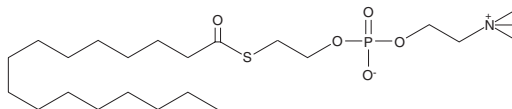
Product Information



HEPC

Catalog No. 10006695

CAS Registry No.: 60793-01-3
Formal Name: 4-oxide-4-hydroxy-N,N,N-trimethyl-9-oxo-3,5-dioxo-8-thia-4-phosphatetracosan-1-aminium, inner salt
Synonym: 2-Hexadecanoylthio-1-Ethylphosphorylcholine
MF: C₂₃H₄₈NO₅PS
FW: 481.7
Purity: ≥95%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that HEPC be stored as supplied at -20°C. It should be stable for at least one year.

HEPC is sparingly soluble in organic solvents such as DMSO and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of HEPC be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of HEPC in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Thioester analogs of glycerophospholipids, in combination with Ellman's reagent, are convenient colorimetric substrates for the measurement of phospholipase activity. HEPC is a truncated analog of phosphatidylcholine with the entire *sn*-1 substituent deleted, and with a hexadecylthioester in the *sn*-2 position.¹ HEPC is a substrate for Type II phospholipase A₂ (PLA₂) enzymes, such as porcine pancreatic, bee venom, and snake venom PLA₂. The intrinsic biological activity of HEPC has not been actively investigated.

Reference

1. Bhat, M.K., Mueller-Harvey, I., Sumner, I.G., *et al.* Simplified methods for the synthesis of 2-hexadecanoylthio-1-ethylphosphorylcholine and for the determination of phospholipase A₂ activity. *Biochem. Biophys. Acta* **1166**, 244-250 (1993).

Related Products

1,2-*bis*(heptanoylthio) Glycerophosphocholine - Cat. No. 62235 • Arachidonoyl Thio-PC - Cat. No. 62240 • 10-Pyrene-PC - Cat. No. 62245

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 02/12/2010