

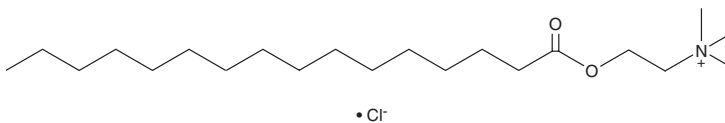
PRODUCT INFORMATION



Palmitoylcholine (chloride)

Item No. 9003456

CAS Registry No.: 2932-74-3
Formal Name: N,N,N-trimethyl-2-[(1-oxohexadecyl)oxy]-ethanaminium, monochloride
MF: C₂₁H₄₄NO₂ • Cl
FW: 378.0
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Palmitoylcholine (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the palmitoylcholine (chloride) in the solvent of choice, which should be purged with an inert gas. Palmitoylcholine (chloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of palmitoylcholine (chloride) in ethanol and DMF is approximately 20 mg/ml and approximately 14 mg/ml in DMSO.

Description

Palmitoylcholine is an acyl choline.¹ It inhibits protein kinase C activity when used at a concentration of 100 μM.² Palmitoylcholine induces hemolysis in rat erythrocytes.³ Plasma levels of palmitoylcholine are decreased in female patients with myalgic encephalomyelitis/chronic fatigue syndrome (ME/CFS).²

References

1. Germain, A., Barupal, D.K., Levine, S.M., *et al.* Comprehensive circulatory metabolomics in ME/CFS reveals disrupted metabolism of acyl lipids and steroids. *Metabolites* **10(1)**, 34 (2020).
2. Nakadate, T. and Blumberg, P.M. Modulation by palmitoylcarnitine of protein kinase C activation. *Cancer Res.* **47(24 Pt 1)**, 6537-6542 (1987).
3. Cho, K.S. and Proulx, P. Interactions of acyl carnitines and other lysins with erythrocytes and reconstituted erythrocyte lipoproteins. *Biochim. Biophys. Acta* **318(1)**, 50-60 (1973).

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, 07/07/2020

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