

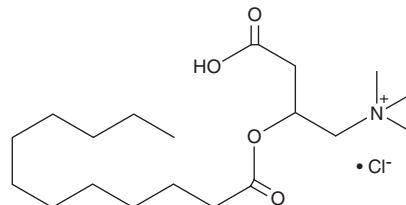
PRODUCT INFORMATION



Lauroyl-DL-Carnitine (chloride)

Item No. 16006

CAS Registry No.: 14919-37-0
Formal Name: 3-carboxy-N,N,N-trimethyl-2-[(1-oxododecyl)oxy]-1-propanaminium, monochloride
Synonyms: CAR 12:0, C12:0 Carnitine, DL-Carnitine lauroyl ester, DL-Lauroylcarnitine
MF: C₁₉H₃₈NO₄ • Cl
FW: 380.0
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

Lauroyl-DL-carnitine (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the lauroyl-DL-carnitine (chloride) in the solvent of choice, which should be purged with an inert gas. Lauroyl-DL-carnitine (chloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of lauroyl-DL-carnitine (chloride) in ethanol and DMF is approximately 20 mg/ml and approximately 10 mg/ml in DMSO.

Description

Lauroyl-DL-carnitine is a zwitterionic, long-chain acylcarnitine used to improve *in vivo* absorption of certain hydrophilic compounds, especially through mucosal membranes.¹⁻³

References

1. Lecluyse, E.L., Sutton, S.C., and Fix, J.A. *In vitro* effects of long-chain acylcarnitines on the permeability, transepithelial electrical resistance and morphology of rat colonic mucosa. *J. Pharmacol. Exp. Ther.* **265**(2), 955-962 (1993).
2. Hayashi, M., Sakai, T., Hasegawa, Y., et al. Physiological mechanism for enhancement of paracellular drug transport. *J. Control. Release* **62**(1-2), 141-148 (1999).
3. Kagatani, S., Inaba, N., Fukui, M., et al. Nasal absorption kinetics of human growth hormone enhanced by acylcarnitines in rats. *Int. J. Pharm.* **169**, 245-253 (1997).

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, 12/12/2022

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