

PRODUCT INFORMATION



Linoleoyl-L-carnitine (chloride)

Item No. 26560

CAS Registry No.: 173686-75-4
Formal Name: [R-(Z,Z)]-3-carboxy-N,N,N-trimethyl-2-[(1-oxo-9,12-octadecadienyl)oxy]-1-propanaminium, monochloride

Synonyms: CAR 18:2, C18:2 Carnitine, L-Carnitine linoleoyl ester, L-Linoleoylcarnitine

MF: C₂₅H₄₆NO₄ • Cl

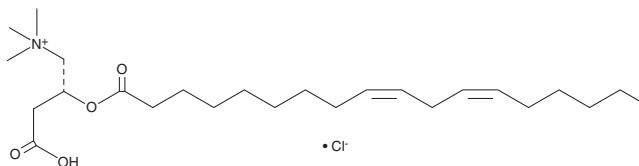
FW: 460.1

Purity: ≥98%

Supplied as: A 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

Linoleoyl-L-carnitine (chloride) is supplied as a solid. A stock solution may be made by dissolving the linoleoyl-L-carnitine (chloride) in the solvent of choice. Linoleoyl-L-carnitine (chloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of linoleoyl-L-carnitine (chloride) in ethanol and DMF is approximately 20 mg/ml and approximately 15 mg/ml in DMSO.

Description

Linoleoyl-L-carnitine is a naturally occurring long-chain acylcarnitine.¹ Hepatic levels of linoleoyl-L-carnitine are increased following high-dose (200 mg/kg) administration of isoniazid (Item No. 20378) in mice.

Reference

1. Li, F., Wang, P., Liu, K., *et al.* A high dose of isoniazid disturbs endobiotic homeostasis in mouse liver. *Drug Metab. Dispos.* **44(11)**, 1742-1751 (2016).

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