

# PRODUCT INFORMATION



## Decanoyl-L-carnitine (chloride)

Item No. 26549

CAS Registry No.: 369651-88-7

Formal Name: (2R)-3-carboxy-N,N,N-trimethyl-2-[(1-oxodecyl)oxy]-1-propanaminium, monochloride

Synonyms: CAR 10:0, C10:0 Carnitine, L-Carnitine decanoyl ester, L-Decanoylcarnitine, (-)-Decanoylcarnitine

MF:  $C_{17}H_{34}NO_4 \cdot Cl$

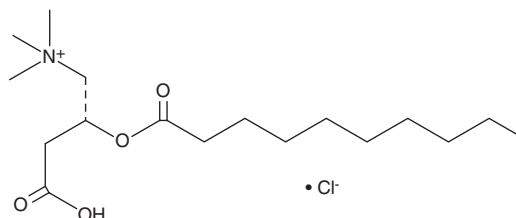
FW: 351.9

Purity:  $\geq 98\%$

Supplied as: A crystalline solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

Decanoyl-L-carnitine (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the decanoyl-L-carnitine (chloride) in the solvent of choice. Decanoyl-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 decanoyl-L-carnitine (chloride) in ethanol and DMF is approximately 20 mg/ml and approximately 14 mg/ml in DMSO.

### Description

Decanoyl-L-carnitine is an ester derivative of L-carnitine (Item No. 21489). It increases the formation of C24 fatty acid intermediates, as well as docosapentaenoic and docosahexaenoic acid (Item No. 90310) in rat hepatocytes.<sup>1</sup>

### Reference

1. Tran, T.N., Retterstøl, K., and Christophersen, B.O. Differences in the conversion of the polyunsaturated fatty acids  $[1-^{14}C]22:4(n-6)$  and  $[1-^{14}C]22:5(n-3)$  to  $[^{14}C]22:5(n-6)$  and  $[^{14}C]22:6(n-3)$  in isolated rat hepatocytes. *Biochim Biophys. Acta.* **1532**(1-2), 137-147 (2001).

#### 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, 02/04/2025

#### 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