

PRODUCT INFORMATION



Decanoyl-L-carnitine (hydrate)

Item No. 26909

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

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

MF: $C_{17}H_{33}NO_4 \cdot XH_2O$

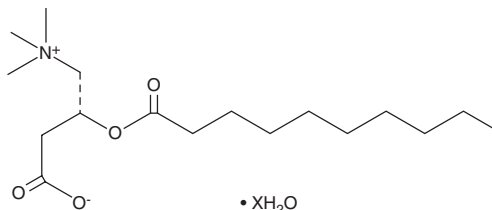
FW: 315.5

Purity: $\geq 95\%$

Supplied as: A solid

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

Stability: ≥ 2 years



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

Laboratory Procedures

Decanoyl-L-carnitine (hydrate) is supplied as a solid. A stock solution may be made by dissolving the decanoyl-L-carnitine (hydrate) in water. We do not recommend storing the aqueous solution for more than one day.

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.¹

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, 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