

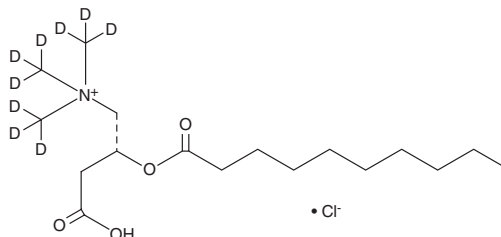
PRODUCT INFORMATION



Decanoyl-L-carnitine-d₉ (chloride)

Item No. 40352

CAS Registry No.: 2847775-92-0
Formal Name: 3-carboxy-N,N,N-tri(methyl-d₃)-2-[(1-oxodecyl)oxy]-1-propanaminium, monochloride
Synonyms: CAR 10:0-d₉, C10:0 Carnitine-d₉, L-Carnitine decanoyl ester-d₉, L-Decanoylcarnitine-d₉, (-)-Decanoylcarnitine-d₉
MF: C₁₇H₂₅D₉NO₄ • Cl
FW: 361.0
Chemical Purity: ≥98% (Decanoyl-L-carnitine (chloride))
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Decanoyl-L-carnitine-d₉ (chloride) is intended for use as an internal standard for the quantification of decanoyl-L-carnitine (Item Nos. 26549 | 26909) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Decanoyl-L-carnitine-d₉ (chloride) is supplied as a solid. A stock solution may be made by dissolving the decanoyl-L-carnitine-d₉ (chloride) in the solvent of choice, which should be purged with an inert gas. Decanoyl-L-carnitine-d₉ (chloride) is sparingly soluble (1-10 mg/ml) in ethanol and 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.¹

Reference

1. Tran, T.N., Retterstøl, K. and Christophersen, B.O. Differences in the conversion of the polyunsaturated fatty acids [1-¹⁴C]22:4(n-6) and [1-¹⁴C]22:5(n-3) to [¹⁴C]22:5(n-6) and [¹⁴C]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, 11/12/2025

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