

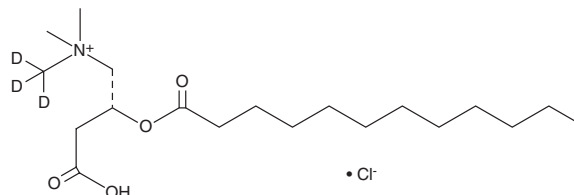
PRODUCT INFORMATION



Lauroyl-L-carnitine-d₃ (chloride)

Item No. 26571

CAS Registry No.: 2687960-76-3
Formal Name: (2R)-3-carboxy-N,N-dimethyl-N-(methyl-d₃)-2-[(1-oxododecyl)oxy]-1-propanaminium, monochloride
Synonyms: C12:0 Carnitine-d₃, CAR 12:0-d₃, L-Carnitine lauroyl ester-d₃, L-Lauroylcarnitine-d₃
MF: C₁₉H₃₅D₃NO₄ • Cl
FW: 383.0
Chemical Purity: ≥98% (Lauroyl-L-carnitine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
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

Lauroyl-L-carnitine-d₃ (chloride) is intended for use as an internal standard for the quantification of lauroyl-L-carnitine (Item No. 26550) 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).

Lauroyl-L-carnitine-d₃ (chloride) is supplied as a solid. A stock solution may be made by dissolving the lauroyl-L-carnitine-d₃ (chloride) in the solvent of choice. Lauroyl-L-carnitine-d₃ (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 lauroyl-L-carnitine-d₃ (chloride) in DMSO is approximately 10 mg/ml and approximately 20 mg/ml in ethanol and DMF.

Description

Lauroyl-L-carnitine is an acylcarnitine and a quaternary ammonium-containing cationic surfactant.¹ It has been used to permeabilize porcine enterocytes for delivery of the polar fluorescent probe lucifer yellow (Item No. 25573).

Reference

1. Danielsen, E.M. and Hansen, G.H. Intestinal surfactant permeation enhancers and their interaction with enterocyte cell membranes in a mucosal explant system. *Tissue Barriers* **5**(3), e1361900 (2017).

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, 03/27/2023

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