

PRODUCT INFORMATION



Palmitoyl-L-carnitine-d₉

Item No. 40353

Formal Name: 2R-3-carboxy-N,N,N-tri(methyl-d₃)-2-[(1-oxohexadecyl)oxy]-1-propanaminium, inner salt, monochloride

Synonyms: CAR 16:0-d₉, C16:0 Carnitine-d₉, L-Carnitine hexadecanoyl ester-d₉, L-Carnitine palmitoyl ester-d₉, Hexadecanoyl-L-carnitine-d₉, L-Hexadecanoylcarnitine-d₉, L-Palmitoylcarnitine-d₉

MF: C₂₃H₃₇D₉NO₄ • Cl
FW: 445.1

Chemical Purity: ≥98% (Palmitoyl-L-carnitine)

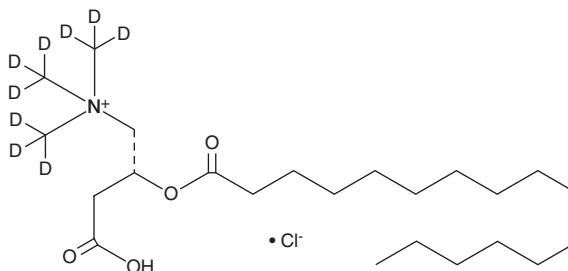
Deuterium

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

Palmitoyl-L-carnitine-d₉ (chloride) is intended for use as an internal standard for the quantification of palmitoyl-L-carnitine 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).

Palmitoyl-L-carnitine-d₉ (chloride) is supplied as a solid. A stock solution may be made by dissolving the palmitoyl-L-carnitine-d₉ (chloride) in the solvent of choice, which should be purged with an inert gas. Palmitoyl-L-carnitine-d₉ (chloride) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Palmitoyl-L-carnitine is a long-chain acylcarnitine and an isomer of palmitoyl-DL-carnitine (Item No. 11095) and palmitoyl-D-carnitine (Item No. 26552).¹ Palmitoyl-L-carnitine is transported into mitochondria via carnitine palmitoyl transferase II to deliver palmitate for fatty acid oxidation and energy production.² It inhibits lecithin:cholesterol acyltransferase activity in isolated rat, but not human, plasma when used at a concentration of 500 μM.³ Serum and hepatic levels of palmitoyl-L-carnitine are increased in mice during cold exposure, and it is taken up by brown adipose tissue.⁴ Palmitoyl-L-carnitine also protects against age-induced cold sensitivity in mice.

References

1. Bezaire, V., Bruce, C.R., Heigenhauser, G.J.F., *et al.* *Am. J. Physiol. Endocrinol. Metab.* **290**(3), E509-E515 (2006).
2. El-Hayek, R., Valdivia, C., Valdivia, H.H., *et al.* *Biophys. J.* **65**(2), 779-789 (1993).
3. Bell, F.P. *Int. J. Biochem.* **15**(2), 133-136 (1983).
4. Simcox, J., Geoghegan, G., Maschek, J.A., *et al.* *Cell Metab.* **26**(3), 509-522 (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, 11/01/2024

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