

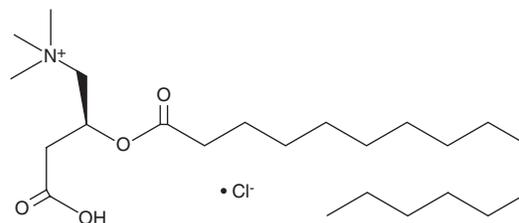
PRODUCT INFORMATION



Palmitoyl-D-carnitine (chloride)

Item No. 26552

CAS Registry No.: 28330-02-1
Formal Name: (S)-3-carboxy-N,N,N-trimethyl-2-(palmitoyloxy)propan-1-aminium, monochloride
Synonyms: CAR 16:0, C16:0 Carnitine, D-Carnitine hexadecanoyl ester, D-Carnitine palmitoyl ester, D-Hexadecanoylcarnitine, Hexadecanoyl-D-carnitine, D-Palmitoylcarnitine
MF: C₂₃H₄₆NO₄ • Cl
FW: 436.1
Purity: ≥98%
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-D-carnitine (chloride) is supplied as a solid. A stock solution may be made by dissolving the palmitoyl-D-carnitine (chloride) in the solvent of choice. Palmitoyl-D-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 palmitoyl-D-carnitine (chloride) in ethanol and DMF is approximately 20 mg/ml and approximately 14 mg/ml in DMSO.

Description

Palmitoyl-D-carnitine is a long-chain acylcarnitine, an isomer of palmitoyl-L-carnitine (Item No. 26553), and the D enantiomer of palmitoyl-DL-carnitine (Item No. 11095).¹ It inhibits carnitine palmitoyltransferase with a K_i value of 2.1 mM for ¹⁴C-palmitoylcarnitine synthesis by erythrocyte membranes.

Reference

1. Wittels, B. and Hochstein, P. The identification of carnitine palmitoyltransferase in erythrocyte membranes. *J. Biol. Chem.* **242**(1), 126-130 (1967).

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