

PRODUCT INFORMATION



Nonadecanoyl-L-carnitine (chloride)

Item No. 9003265

Formal Name: (R)-3-carboxy-N,N,N-trimethyl-2-(nonadecanoyloxy)propan-1-aminium, monochloride

Synonyms: CAR 19:0, C19:0 Carnitine, L-Carnitine nonadecanoyl ester, L-Carnitine pristanoyl ester, L-Nonadecanoylcarnitine, Pristanoylcarnitine

MF: $C_{26}H_{52}NO_4 \cdot Cl$

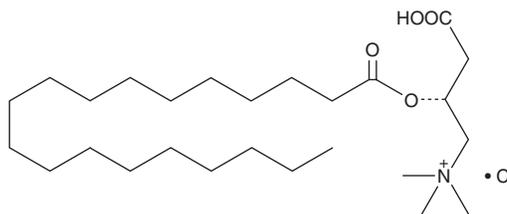
FW: 478.2

Purity: $\geq 95\%$

Supplied as: A crystalline 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

Nonadecanoyl-L-carnitine (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the nonadecanoyl-L-carnitine (chloride) in the solvent of choice, which should be purged with an inert gas. Nonadecanoyl-L-carnitine (chloride) is soluble in methanol.

Description

Nonadecanoyl-L-carnitine is a long-chain acylcarnitine composed of nonadecanoic acid (Item No. 19723) and L-carnitine (Item No. 21489). Serum levels of nonadecanoyl-L-carnitine are decreased in children and adolescents chronically exposed to industrial carcinogenic pollutants.¹

Reference

1. Chen, C.-H.S., Kuo, T.-C., Kuo, H.-C., *et al.* Metabolomics of children and adolescents exposed to industrial carcinogenic pollutants. *Environ. Sci. Technol.* **53(9)**, 5454-5465 (2019).

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, 12/09/2022

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