

PRODUCT INFORMATION



Docosahexaenoyl-L-carnitine (chloride)

Item No. 38350

Formal Name: 3-carboxy-N,N,N-trimethyl-2R-[[1-oxo-4Z,7Z,10Z,13Z,16Z,19Z-docosahexaen-1-yl]oxy]-1-propanaminium, inner salt, monochloride

Synonyms: CAR 22:6, C22:6 Carnitine, L-Cervonoylcarnitine, L-Carnitine cervonoyl ester

MF: $C_{29}H_{46}NO_4 \cdot Cl$

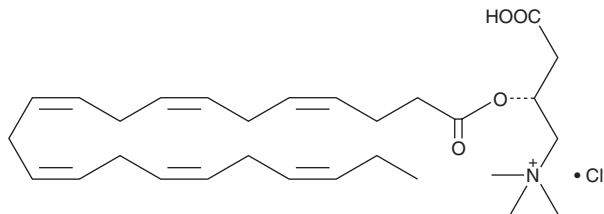
FW: 508.1

Purity: $\geq 95\%$

Supplied as: A 10 mg/ml solution in ethanol

Storage: $-80^{\circ}C$

Stability: ≥ 2 years



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

Description

Docosahexaenoyl-L-carnitine is a long-chain acylcarnitine composed of docosahexaenoic acid (DHA; Item No. 90310) and L-carnitine (Item No. 21489). Decreased venous cord plasma levels of docosahexaenoyl-L-carnitine are associated with babies born small for gestational age.¹

Reference

1. Horgan, R.P., Broadhurst, D.I., Walsh, S.K., *et al.* Metabolic profiling uncovers a phenotypic signature of small for gestational age in early pregnancy. *J. Proteome Res.* **10(8)**, 3660-3673 (2011).

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