

# 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<sub>29</sub>H<sub>46</sub>NO<sub>4</sub> • Cl

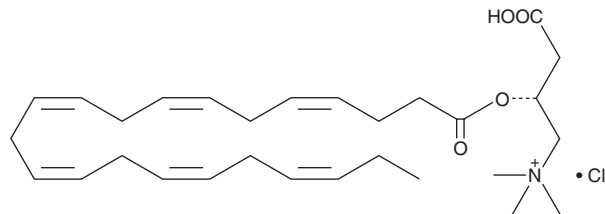
**FW:** 508.1

**Purity:** ≥95%

**Supplied as:** A solution in ethanol

**Storage:** -80°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.<sup>1</sup>

### 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, 08/14/2023

#### 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