

PRODUCT INFORMATION



Eicosapentaenoyl-L-carnitine (chloride)

Item No. 9003924

CAS Registry No.: 2133456-02-5

Formal Name: 3-carboxy-N,N,N-trimethyl-2-[[1-oxo-5Z,8Z,11Z,14Z,17Z-eicosapentaen-1-yl]oxy]-1-propanaminium, monochloride

Synonyms: CAR 20:5, C20:5 Carnitine, L-Carnitine eicosapentaenoyl ester, L-Carnitine timnodonoyl ester, L-Eicosapentaenoylcarnitine, L-Timnodonoylcarnitine

MF: $C_{27}H_{44}NO_4 \cdot Cl$

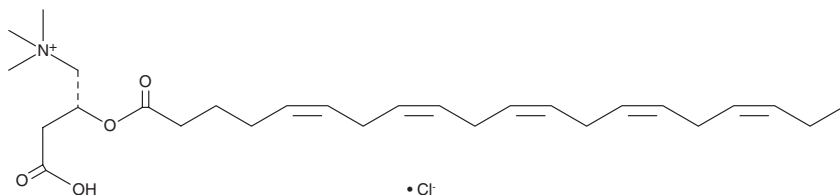
FW: 482.1

Purity: $\geq 90\%$

Supplied as: A 25 mg/ml solution in ethanol

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

Stability: ≥ 1 year



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

Description

Eicosapentaenoyl-L-carnitine is a long-chain acylcarnitine composed of eicosapentaenoic acid (EPA; Item Nos. 90110 | 90110.1 | 21908) and L-carnitine (Item No. 21489).¹ The levels of eicosapentaenoyl-L-carnitine are increased in the eyes of mice fed a diet high in *n*-3/*n*-6 polyunsaturated fatty acids (PUFAs), but not in mice fed normal diet, in a mouse model of myopia induced by out-of-focus lenses.

Reference

1. Mori, K., Kuroha, S., Hou, J., *et al.* Lipidomic analysis revealed *n*-3 polyunsaturated fatty acids suppressed choroidal thinning and myopia progression in mice. *FASEB J.* **36**(6), e22312 (2022).

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, 01/13/2026

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