

PRODUCT INFORMATION



Heptadecanoyl-L-carnitine (chloride)

Item No. 37934

CAS Registry No.: 173686-69-6

Formal Name: 3-carboxy-N,N,N-trimethyl-

Synonyms: 2-[(1-oxoheptadecyl)oxy]-1-propanaminium, monochloride
CAR 17:0, C17:0 Carnitine,
L-Carnitine heptadecanoyl ester,
L-Heptadecanoylcarnitine

MF: $C_{24}H_{48}NO_4 \cdot Cl$

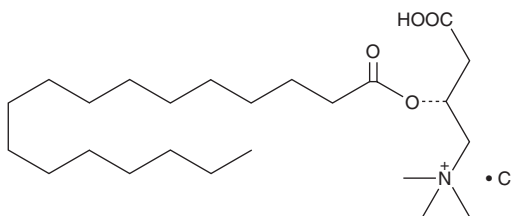
FW: 450.1

Purity: $\geq 95\%$

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Heptadecanoyl-L-carnitine (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the heptadecanoyl-L-carnitine (chloride) in the solvent of choice, which should be purged with an inert gas. Heptadecanoyl-L-carnitine (chloride) is soluble in chloroform and methanol. The solubility of heptadecanoyl-L-carnitine (chloride) in these solvents is approximately 30 mg/ml.

Description

Heptadecanoyl-L-carnitine is a long-chain acylcarnitine composed of heptadecanoic acid (Item No. 19722) and L-carnitine (Item No. 21489). Levels of heptadecanoyl-L-carnitine are increased in dried blood spots from neonates with propionic or methylmalonic acidemias, inborn errors of metabolism characterized by a deficiency in the activity of propionyl-CoA carboxylase or L-methylmalonyl-CoA mutase or its cofactor 5-deoxyadenosylcobalamin, respectively.¹

Reference

1. Malvagia, S., Haynes, C.A., Grisotto, L., *et al.* Heptadecanoylcarnitine (C17) a novel candidate biomarker for propionic and methylmalonic acidemias during expanded newborn screening. *Clin. Chim. Acta* 450, 342-348 (2015).

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