

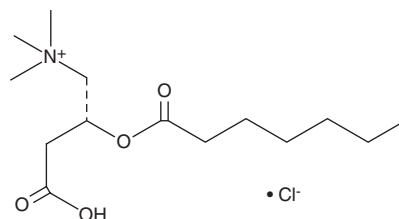
# PRODUCT INFORMATION



## Heptanoyl-L-carnitine (chloride)

Item No. 26551

**CAS Registry No.:** 162041-08-9  
**Formal Name:** (R)-3-carboxy-N,N,N-trimethyl-2-[(1-oxoheptyl)oxy]-1-propanaminium, monochloride  
**Synonyms:** C7:0 Carnitine, CAR 7:0, L-Carnitine heptanoyl ester, L-Heptanoylcarnitine  
**MF:**  $C_{14}H_{26}NO_4 \cdot Cl$   
**FW:** 309.8  
**Purity:**  $\geq 95\%$   
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Heptanoyl-L-carnitine (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the heptanoyl-L-carnitine (chloride) in the solvent of choice. Heptanoyl-L-carnitine (chloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of heptanoyl-L-carnitine (chloride) in ethanol and DMF is approximately 20 mg/ml and approximately 10 mg/ml in DMSO.

### Description

Heptanoyl-L-carnitine is a medium-chain acylcarnitine and derivative of L-carnitine (Item No. 21489).

**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, 02/16/2024

### 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