

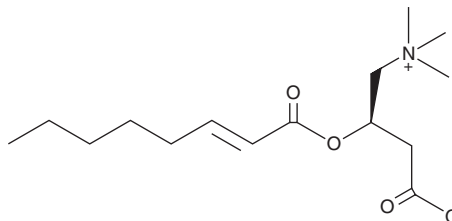
PRODUCT INFORMATION



trans-2-Octenoyl-L-carnitine

Item No. 30995

CAS Registry No.: 152064-94-3
Formal Name: (2R)-3-carboxy-N,N,N-trimethyl-2-[[[(2E)-1-oxo-2-octen-1-yl]oxy]-1-propanaminium, inner salt
Synonyms: CAR 8:1, C8:1 Carnitine, L-Carnitine *trans*-2-octenoyl ester, L-*trans*-2-Octenoylcarnitine
MF: C₁₅H₂₇NO₄
FW: 285.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

trans-2-Octenoyl-L-carnitine is supplied as a solid. A stock solution may be made by dissolving the *trans*-2-octenoyl-L-carnitine in the solvent of choice, which should be purged with an inert gas. *trans*-2-Octenoyl-L-carnitine is slightly soluble in DMSO (warmed) and sparingly soluble in methanol.

Description

trans-2-Octenoyl-L-carnitine is a synthetic carnitine.¹ It has been used as a standard for the detection of carnitines in human urine.

Reference

1. Yang, S., Minkler, P., and Hoppel, C. *cis*-3,4-methylene-heptanoylcarnitine: Characterization and verification of the C8:1 acylcarnitine in human urine. *J. Chromatogr. B. Analyt. Technol. Biomed. Life Sci.* **857**(2), 251-258 (2007).

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