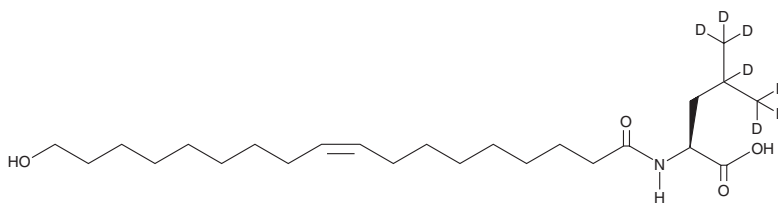


PRODUCT INFORMATION



18-hydroxy Oleoyl Leucine-d₇ Item No. 9004331

MF: C₂₄H₃₈D₇NO₄
FW: 418.7
Chemical Purity: ≥95% (18-hydroxy Oleoyl Leucine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 years



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

Laboratory Procedures

18-hydroxy Oleoyl leucine-d₇ is intended for use as an internal standard for the quantification of 18-hydroxy oleoyl leucine by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Description

18-hydroxy oleoyl leucine is a ω-hydroxy metabolite of the N-acyl amide mitochondrial uncoupling agent N-oleoyl leucine (Item No. 20064).¹ It is formed from N-oleoyl leucine by the cytochrome P450 (CYP) isoform CYP4F2.

Reference

1. Tanzo, J.T., Li, V.L., Wiggenhorn, A.L., *et al.* CYP4F2 is a human-specific determinant of circulating N-acyl amino acid levels. *J. Biol. Chem.* **299**(6), 104764 (2023).

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, 01/24/2024

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