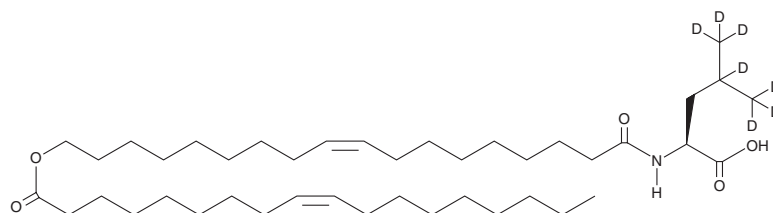


PRODUCT INFORMATION



Oleic Acid-hydroxy Oleoyl Leucine-d₇ Item No. 9004332

Synonyms:	C18:1-hydroxy Oleoyl Leucine-d ₇ , OAHOL-d ₇
MF:	C ₄₂ H ₇₀ D ₇ NO ₅
FW:	683.1
Chemical Purity:	≥95% (Oleic acid-hydroxy oleoyl leucine)
Deuterium	
Incorporation:	≥99% deuterated forms (d ₁ -d ₇); ≤1% d ₀
Supplied as:	A solution in ethanol
Storage:	-20°C
Stability:	≥1 year



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

Laboratory Procedures

Oleic acid-hydroxy oleoyl leucine-d₇ is intended for use as an internal standard for the quantification of oleic acid-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

Oleic acid-hydroxy oleoyl leucine is a very long-chain fatty acid composed of oleic acid (Item Nos. 90260 | 24659) connected to hydroxy oleoyl leucine.¹

Reference

1. Tanzo, J.T., Li, V.L., Wiggernhorn, 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, 01/26/2024

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