

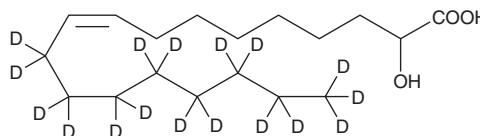
PRODUCT INFORMATION



2-hydroxy Oleic Acid-d₁₇

Item No. 9004307

CAS Registry No.: 1346606-40-3
Formal Name: (Z)-2-hydroxyoctadec-9-enoic-11,11,12,12,13,13,14,14,15,15,16,16,17,17,18,18,18-d₁₇ acid
Synonyms: FA 18:1;O-d₁₇, 2-HOME-d₁₇, 2-hydroxy-*cis*-9-Octadecenoic Acid-d₁₇, α-hydroxy-*cis*-9-Octadecenoic Acid-d₁₇, 2-hydroxy-9(Z)-OME-d₁₇, 2-Hydroxyoctadecenoic Acid-d₁₇, 2-OHOA-d₁₇
MF: C₁₈H₁₇D₁₇O₃
FW: 315.6
Chemical Purity: ≥95% (2-hydroxy Oleic Acid)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₁₇); ≤1% d₀
Supplied as: A 1 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

2-hydroxy Oleic acid-d₁₇ is intended for use as an internal standard for the quantification of 2-hydroxy oleic acid (Item No. 36044) 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).

2-hydroxy Oleic acid-d₁₇ is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy Oleic acid-d₁₇ in the solvent of choice, which should be purged with an inert gas. 2-hydroxy Oleic acid-d₁₇ is soluble in chloroform.

Description

2-hydroxy Oleic acid is a synthetic monounsaturated hydroxylated fatty acid.¹ It decreases the lamellar-to-hexagonal transition temperature of bovine liver phosphatidylethanolamine (PE) membranes by approximately 12.5°C when included at a 20:1 molar ratio of PE:2-hydroxy oleic acid. 2-hydroxy Oleic acid increases levels of protein kinase C α (PKCα) in and decreases proliferation of A549 non-small cell lung cancer (NSCLC) cells in a concentration-dependent manner. It reduces tumor growth in an A549 mouse xenograft model when administered at doses of 75 and 200 mg/kg. 2-hydroxy Oleic acid has been used in the generation of liposomes.²

References

1. Martínez, J., Vögler, O., Casas, J., *et al.* Membrane structure modulation, protein kinase Cα activation, and anticancer activity of minerval. *Mol. Pharmacol.* **67**(2), 531-540 (2004).
2. Olechowska, K., Mach, M., Hąc-Wydro, K., *et al.* Studies on the interactions of 2-hydroxyoleic acid with monolayers and bilayers containing cationic lipid: Searching for the formulations for more efficient drug delivery to cancer cells. *Langmuir* **35**(27), 9084-9092 (2019).

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, 07/14/2026

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