

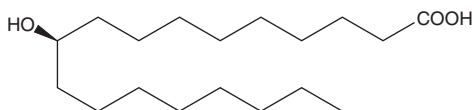
PRODUCT INFORMATION



10(R)-hydroxy Stearic Acid

Item No. 42929

CAS Registry No.: 5856-32-6
Formal Name: 10R-hydroxy-octadecanoic acid
Synonyms: FA 18:0;O, 10(R)-HSA, 10(R)-hydroxy Octadecanoic Acid, 10(R)-hydroxy ODA
MF: $C_{18}H_{36}O_3$
FW: 300.5
Purity: $\geq 98\%$
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

10(R)-hydroxy Stearic acid is supplied as a solid. A stock solution may be made by dissolving the 10(R)-hydroxy stearic acid in the solvent of choice, which should be purged with an inert gas. 10(R)-hydroxy Stearic acid is soluble (≥ 10 mg/ml) in DMSO and ethanol.

Description

10(R)-hydroxy Stearic acid is a hydroxy fatty acid and hydroxylation product of oleic acid (Item Nos. 90260 | 24659) formed by gut microbiota.^{1,2}

References

1. Prem, S., Helmer, C.P.O., Dimos, N., *et al.* Towards an understanding of oleate hydratases and their application in industrial processes. *Microb. Cell Fact.* **21(1)**, 58 (2022).
2. Serra, S. and De Simeis, D. Use of *Lactobacillus rhamnosus* (ATCC 53103) as whole-cell biocatalyst for the regio- and stereoselective hydration of oleic, linoleic, and linolenic acid. *Catalysts* **8(3)**, 109 (2018).

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, 02/06/2025

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