

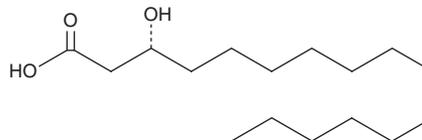
PRODUCT INFORMATION



3(R)-hydroxy Palmitic Acid

Item No. 45136

CAS Registry No.: 20595-04-4
Formal Name: (3R)-3-hydroxy-hexadecanoic acid
Synonyms: FA 16:0;O, D-3-hydroxy Hexadecanoic Acid, NSC 179484, D-3-hydroxy Palmitic Acid
MF: C₁₆H₃₂O₃
FW: 272.4
Purity: ≥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

3(R)-hydroxy Palmitic acid is supplied as a solid. A stock solution may be made by dissolving the 3(R)-hydroxy palmitic acid in the solvent of choice, which should be purged with an inert gas. 3(R)-hydroxy Palmitic acid is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

3(R)-hydroxy Palmitic acid is a hydroxy fatty acid and a component of lipid A and plipastatin A1 (Item No. 38916).^{1,2}

References

- Holst, O., Borowiak, D., Weckesser, J., et al. Structural studies on the phosphate-free lipid A of *Rhodocrobium vannielii* ATCC 17100. *Eur. J. Biochem.* **137(1-2)**, 325-333 (1983).
- Umezawa, H., Ayoagi, T., Nishikiori, T., et al. Plipastatins: New inhibitors of phospholipase A₂, produced by *Bacillus cereus* BMG302-ff67. I. Taxonomy, production, isolation and preliminary characterization. *J. Antibiot. (Tokyo)* **39(6)**, 737-744 (1986).

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, 02/17/2026

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