

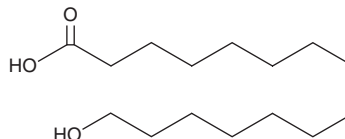
PRODUCT INFORMATION



15-hydroxy Pentadecanoic Acid

Item No. 24648

CAS Registry No.: 4617-33-8
Formal Name: 15-hydroxy-pentadecanoic acid
Synonyms: FA 15:0;O, ω-hydroxy Pentadecanoic Acid
MF: C₁₅H₃₀O₃
FW: 258.4
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

15-hydroxy Pentadecanoic acid is supplied as a solid. A stock solution may be made by dissolving the 15-hydroxy pentadecanoic acid in the solvent of choice, which should be purged with an inert gas. 15-hydroxy Pentadecanoic acid is soluble in organic solvents such as ethanol, methanol, and chloroform.

Description

15-hydroxy Pentadecanoic acid is a hydroxylated fatty acid that has been used in the synthesis of the macrocyclic lactone musk compounds exaltolide and phoracantholide I.¹

Reference

1. Cossy, J. and Pete, J.P. A three step synthesis of exaltolide and phoracantholide I. *Tetrahedron Lett.* **27(21)**, 2369-2370 (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, 02/24/2027

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