

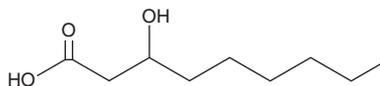
# PRODUCT INFORMATION



## 3-hydroxy Nonanoic Acid

Item No. 24611

**CAS Registry No.:** 40165-87-5  
**Formal Name:** 3-hydroxy-nonanoic acid  
**Synonyms:** FA 9:0;O,  
 $\beta$ -hydroxy Nonanoic acid  
**MF:** C<sub>9</sub>H<sub>18</sub>O<sub>3</sub>  
**FW:** 174.2  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

3-hydroxy Nonanoic acid is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxy nonanoic acid in the solvent of choice, which should be purged with an inert gas. 3-hydroxy Nonanoic acid is soluble in organic solvents such as chloroform, ethanol, and methanol.

### Description

3-hydroxy Nonanoic acid is a hydroxylated fatty acid that has been found in young peridioles of *Pisolithus* fungi, LPS from *P. aeruginosa*, bovine milk, and the methyl-branched poly(3-hydroxyalkanoate) (PHA) polymers produced by *P. oleovorans*.<sup>1-4</sup>

### References

1. Campos, A.N.R., Costa, M.D., Tótoia, M.R., *et al.* Total lipid and fatty acid accumulation during basidiospore formation in the ectomycorrhizal fungus *Pisolithus sp. R. Bras. Ci. Solo* **32(4)**, 1531-1540 (2008).
2. Uhlig, S., Negård, M., Heldal, K.K., *et al.* Profiling of 3-hydroxy fatty acids as environmental markers of endotoxin using liquid chromatography coupled to tandem mass spectrometry. *J. Chromatogr. A* **1434**, 119-236 (2016).
3. Parks, O.W. Isolation and characterization of nonesterified 3-hydroxy acids in milk. *J. Dairy Sci.* **60(5)**, 718-720 (1977).
4. Hazer, B., Lenz, R.W., and Fuller, R.C. Biosynthesis of methyl-branched poly( $\beta$ -hydroxyalkanoate)s by *Pseudomonas oleovorans*. *Macromol.* **27(1)**, 45-49 (1994).

#### 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, 02/27/2024

#### 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