

PRODUCT INFORMATION



3-hydroxy Stearic Acid

Item No. 22692

CAS Registry No.: 17773-30-7
Formal Name: 3-hydroxy-octadecanoic Acid
Synonyms: 3-hydroxy Octadecanoic Acid,
 β -hydroxy Octadecanoic Acid,
 β -hydroxy Stearic acid

MF: $C_{18}H_{36}O_3$

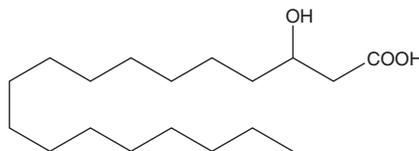
FW: 300.5

Purity: $\geq 95\%$

Supplied as: A crystalline 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-hydroxy Stearic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-hydroxy stearic acid in the solvent of choice, which should be purged with an inert gas. 3-hydroxy Stearic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 3-hydroxy stearic acid in these solvents is approximately 20, 10, and 30 mg/ml, respectively.

Description

3-hydroxy Stearic acid is a long-chain saturated fatty acid and intermediate in the production of 3-hydroxy octadecanedioic acid from stearic acid (Item No. 10011298).¹ 3-hydroxy Stearic acid is found in *C. pyloridis*, a bacteria implicated in the etiology of peptic ulcers.²

References

1. Dhar, M., Sepkovic, D.W., Hirani, V., *et al.* Omega oxidation of 3-hydroxy fatty acids by the human CYP4F gene subfamily enzyme CYP4F11. *J. Lipid Res.* **49(3)**, 612-624 (2008).
2. Lambert, M.A., Patton, C.M., Barrett, T.J., *et al.* Differentiation of *Campylobacter* and *Campylobacter*-like organisms by cellular fatty acid composition. *J. Clin. Microbiol.* **25(4)**, 706-713 (1987).

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, 01/04/2023

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