

PRODUCT INFORMATION



3-hydroxy Heptadecanoic Acid

Item No. 28735

CAS Registry No.: 40165-89-7

Synonyms: β -hydroxy Heptadecanoic Acid,
3-hydroxy Heptadecylic Acid,
3-hydroxy Margaric Acid, FA 17:0;O

MF: $C_{17}H_{34}O_3$

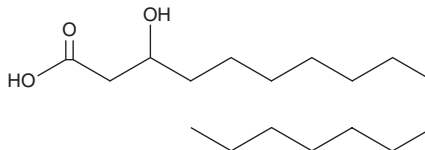
FW: 286.5

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

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

Description

3-hydroxy Heptadecanoic acid is a hydroxylated fatty acid. It is a component of LPS in several species of Gram-negative bacteria.^{1,2} Levels of 3-hydroxy heptadecanoic acid are decreased in the feces of patients with Crohn's disease.³

References

1. Weintraub, A., Larsson, B.E., and Lindberg, A.A. Chemical and immunochemical analyses of *Bacteroides fragilis* lipopolysaccharides. *Infect. Immun.* **49**(1), 197-201 (1985).
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. De La Cochetiere, M.F., Rouge, C., Szponar, B., et al. 3-Hydroxy fatty acids in faeces of patients with Crohn's disease or ulcerative colitis as a non-invasive diagnostic approach. *Microb. Ecol. Health Dis.* **19**(1), 1-6 (2007).

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, 03/26/2024

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