

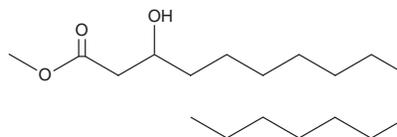
PRODUCT INFORMATION



3-hydroxy Heptadecanoic Acid methyl ester

Item No. 24646

CAS Registry No.: 112538-92-8
Formal Name: 3-hydroxy-heptadecanoic acid, methyl ester
Synonyms: Methyl 3-hydroxyheptadecanoate, SFE 18:0;O
MF: C₁₈H₃₆O₃
FW: 300.5
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-hydroxy Heptadecanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxy heptadecanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 3-hydroxy Heptadecanoic acid methyl ester is soluble in organic solvents such as ethanol and methanol.

Description

3-hydroxy Heptadecanoic acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in *B. thetaiotaomicron*.^{1,2}

References

1. Vannieuwenhuyze, F. and Sandra, P. Selectivity optimization for the capillary gas chromatographic analysis of bacterial cellular fatty acids. *Chromatographia* **23(11)**, 850-855 (1987).
2. Husmann, H., Schomburg, G., Müller, K.-D., *et al.* Gas chromatographic FAME analysis for characterization of bacteria: Formation of aldehydes from 3-hydroxy fatty acid methyl esters during sample introduction. *J. Sep. Sci.* **13(11)**, 780-782 (1990).

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, 04/02/2024

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