

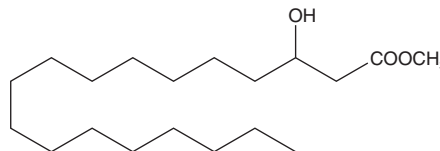
PRODUCT INFORMATION



3-hydroxy Stearic Acid methyl ester

Item No. 24647

CAS Registry No.: 2420-36-2
Formal Name: 3-hydroxy-octadecanoic acid, methyl ester
Synonyms: 3-Hydroxy C18:0 methyl ester,
Methyl 3-hydroxyoctadecanoate, SFE 19:0;O
MF: C₁₉H₃₈O₃
FW: 314.5
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 Stearic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxy stearic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 3-hydroxy Stearic acid methyl ester is soluble in ethanol and methanol.

Description

3-hydroxy Stearic acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in methyl-branched poly(3-hydroxyalkanoate) (PHA) polymers produced by *Streptomyces* sp. JM3 and in trace amounts in *M. bijuga* fruit.^{1,2} It is a chemical scent constituent in feces of adult wild Iberian wolves (*C. lupus signatus*).³

References

1. Allen, A.D., Daley, P., Ayorinde, F.O., et al. Characterization of medium chain length (R)-3-hydroxycarboxylic acids produced by *Streptomyces* sp. JM3 and the evaluation of their antimicrobial properties. *World J. Microbiol. Biotechnol.* **28(9)**, 2791-2800 (2012).
2. Puino, J.A., Marbot, R., and Agüero, J. Volatile components of geneps (*Melicocca bijuga* L.) fruit. *J. Essent. Oil Res.* **14(4)**, 247-248 (2002).
3. Martín, J., Barja, J., and López, P. Chemical scent constituents in feces of wild Iberian wolves (*Canis lupus signatus*). *Biochem. System. Ecol.* **38(6)**, 1096-1102 (2010).

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, 03/04/2024

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