

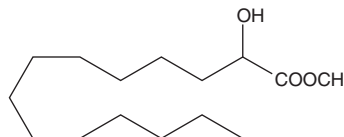
PRODUCT INFORMATION



2-hydroxy Myristic Acid methyl ester

Item No. 24588

CAS Registry No.: 56009-40-6
Formal Name: 2-hydroxy-tetradecanoic acid, methyl ester
Synonym: SFE 15:0;O
MF: C₁₅H₃₀O₃
FW: 258.4
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

2-hydroxy Myristic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy myristic acid methyl ester in the solvent of choice. 2-hydroxy Myristic acid methyl ester is soluble in organic solvents such as methanol, chloroform, and ethyl ether, which should be purged with an inert gas.

Description

2-hydroxy Myristic acid methyl ester is a rare hydroxylated fatty acid methyl ester that has been isolated from aquatic microbial sources.¹ Levels of 2-hydroxy myristic acid methyl ester decrease to <1% in aquatic microbial communities after administration of the pesticide Atradox at a concentration of 245 mg/L.

Reference

1. Littlefield-Wyer, J.G., Brooks, P., and Katouli, M. Application of biochemical fingerprinting and fatty acid methyl ester profiling to assess the effect of the pesticide Atradox on aquatic microbial communities. *Environ. Pollut.* **153**(2), 393-400 (2008).

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, 11/08/2022

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