

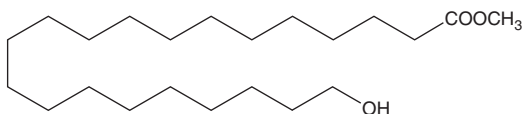
PRODUCT INFORMATION



21-hydroxy Heneicosanoic Acid methyl ester

Item No. 24654

CAS Registry No.: 94035-98-0
Synonyms: Methyl
21-hydroxyheneicosanoate,
SFE 22:0;O
MF: C₂₂H₄₄O₃
FW: 356.6
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

21-hydroxy Heneicosanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 21-hydroxy heneicosanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 21-hydroxy Heneicosanoic acid methyl ester is soluble in chloroform, warm ethanol, and ethyl ether.

Description

21-hydroxy Heneicosanoic acid methyl ester is a hydroxylated fatty acid methyl ester and substrate for lactonizing lipase from *Pseudomonas* nov. sp. 109, a lipase that catalyzes the synthesis of macrocyclic lactones.¹

Reference

1. Ihara, F., Kageyama, Y., Hirata, M., *et al.* Purification, characterization, and molecular cloning of lactonizing lipase from *Pseudomonas* species. *J. Biol. Chem.* **266(27)**, 18135-18140 (1991).

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