

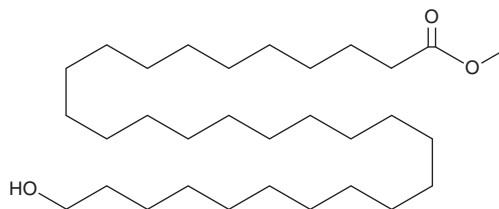
PRODUCT INFORMATION



30-hydroxy Triacontanoic Acid methyl ester

Item No. 24808

CAS Registry No.: 79162-70-2
Formal Name: 30-hydroxy-triacontanoic acid, methyl ester
MF: $C_{31}H_{62}O_3$
FW: 482.8
Purity: $\geq 97\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

30-hydroxy Triacontanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 30-hydroxy triacontanoic acid methyl ester in the solvent of choice. 30-hydroxy Triacontanoic acid methyl ester is soluble in the organic solvent chloroform, which should be purged with an inert gas.

Description

30-hydroxy Triacontanoic acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in sediment samples from the Harney River.¹

Reference

1. Jaffé, R., Rushdi, A.I., Medeiros, P.M., *et al.* Natural product biomarkers as indicators of sources and transport of sedimentary organic matter in a subtropical river. *Chemosphere* **64(11)**, 1870-1884 (2006).

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, 06/04/2018

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