

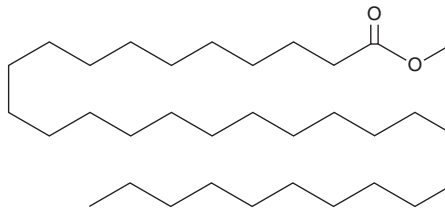
PRODUCT INFORMATION



Dotriacontanoic Acid methyl ester

Item No. 26728

CAS Registry No.: 41755-79-7
Formal Name: dotriacontanoic acid, methyl ester
Synonyms: C32:0 FAME, Lacceric Acid methyl ester, Methyl Dotriacontanoate, SFE 33:0
MF: C₃₃H₆₆O₂
FW: 494.9
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

Dotriacontanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the dotriacontanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Dotriacontanoic acid methyl ester is soluble in chloroform and methylene chloride.

Description

Dotriacontanoic acid methyl ester is a naturally occurring fatty acid methyl ester that has been found in the cuticular wax of *P. abies* needles.¹ It has also been found in sediment samples from the Harney River in Florida and Lake Kivu in the East African rift valley.^{2,3}

References

1. Sümmechen, P., Markstädter, C., and Wienhaus, O. Composition of the epicuticular wax esters of *Picea abies* (L.) Karst. *Zeitschrift für Naturforschung* **50(1-2)**, 11-14 (2014).
2. 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).
3. Al-Mutlaq, K., Standley, L.J., and Simoneit, B.R. Composition and sources of extractable organic matter from a sediment core in Lake Kivu, East African rift valley. *Appl. Geochem.* **23(5)**, 1023-1040 (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, 02/13/2024

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