

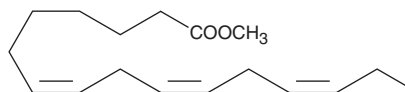
PRODUCT INFORMATION



7(Z),10(Z),13(Z)-Hexadecatrienoic Acid methyl ester

Item No. 10005161

CAS Registry No.: 77388-16-0
Synonyms: C16:3 (*cis*-7,10,13) methyl ester,
Methyl *cis*-7,10,13-Hexadecatrienoate,
Roughanic Acid methyl ester
MF: C₁₇H₂₈O₂
FW: 264.4
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Description

7(Z),10(Z),13(Z)-Hexadecatrienoic acid methyl ester is a fatty ester that has been found in *A. fumigatus* and a methyl ester form of 7(Z),10(Z),13(Z)-hexadecatrienoic acid (Item No. 10005156).¹

Reference

1. Mohamed, A.F., Abuamara, T.M.M., Amer, M.E., *et al.* Genetic and histopathological alterations in Caco-2 and HuH-7 cells treated with secondary metabolites of marine fungi. *J. Gastrointest. Cancer.* **53(2)**, 480-495 (2022).

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, 02/12/2024

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