

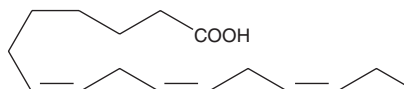
PRODUCT INFORMATION



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

Item No. 10005156

CAS Registry No.: 7561-64-0
Formal Name: (7Z,10Z,13Z)-hexadecatrienoic acid
Synonyms: C16:3 *n*-3, C16:3(7Z,10Z,13Z), FA 16:0, all-*cis*-7,10,13-Hexadecatrienoic Acid, Roughanic Acid
MF: C₁₆H₂₆O₂
FW: 250.4
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 is an ω-3 polyunsaturated fatty acid (PUFA) that has been found in *S. olusatrum* leaves.¹ It is a precursor in the biosynthesis of jasmonic acid (Item No. 88300) in *A. thaliana* leaves.²

References

1. Caprioli, G., Fiorini, D., Maggi, F., *et al.* Ascorbic acid content, fatty acid composition and nutritional value of the neglected vegetable Alexanders (*Smyrniolus olusatrum* L., Apiaceae). *J. Food Compos. Anal.* **35**(1), 30-36 (2014).
2. Snoeren, T.A.L., Van Poecke, R.M.P., and Dicke, M. Multidisciplinary approach to unravelling the relative contribution of different oxylipins in indirect defense of *Arabidopsis thaliana*. *J. Chem. Ecol.* **35**(9), 1021-1031 (2009).

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, 03/08/2024

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