

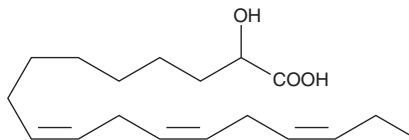
PRODUCT INFORMATION



(±)2-HOTrE

Item No. 39100

CAS Registry No.: 22324-71-6
Formal Name: 2-hydroxy-9Z,12Z,15Z-octadecatrienoic acid
Synonyms: FA 18:3;O, (±)2-9(Z),12(Z),15(Z)-HOTrE, α-Hydroxylinolenic Acid, 2-Hydroxylinolenic Acid, (±)-2-hydroxy-9(Z),12(Z),15(Z)-OTrE
MF: C₁₈H₃₀O₃
FW: 294.4
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

(±)2-HOTrE is an oxylipin.¹ It is active against the plant pathogenic fungi *A. niger*, *S. cerevisiae*, and *P. membranae-faciens* (MICs = 1.5, 3, and 1 g/L, respectively), as well as *P. roqueforti*, *C. albicans*, and *C. valida* (MICs = 2.33, 3, and 2 g/L, respectively). Hepatic levels of 2-HOTrE are increased in a mouse model of non-alcoholic fatty liver disease (NAFLD) induced by a high-fat diet.²

References

- Liang, N., Dacko, A., Tan, A.K., *et al.* Structure-function relationships of antifungal monohydroxy unsaturated fatty acids (HUFA) of plant and bacterial origin. *Food Res. Int.* **134**, 109237 (2020).
- Wu, Z., Xiao, H., Rao, D., *et al.* Analytical strategy for oxylipin annotation by combining chemical derivatization-based retention index algorithm and feature tandem mass spectrometric fragmentation as a biomarker discovery tool. *Anal. Chem.* **95(43)**, 15933-15942 (2023).

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