

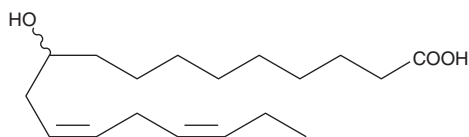
PRODUCT INFORMATION



(±)-10-hydroxy-12(Z),15(Z)-Octadecadienoic Acid

Item No. 37280

CAS Registry No.: 34932-14-4
Formal Name: 10-hydroxy-12Z,15Z-octadecadienoic acid
Synonyms: 12(Z),15(Z)-10-HODE,
(±)-10-hydroxy-12(Z),15(Z)-ODE
MF: C₁₈H₃₂O₃
FW: 296.5
Purity: ≥98%
Supplied as: A 5 mg/ml 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

(±)-10-hydroxy-12(Z),15(Z)-Octadecadienoic acid is an oxylipin. It is produced from α-linolenic acid (Item Nos. 90210 | 21910) by the gut bacteria *L. plantarum*.¹ (±)-10-hydroxy-12(Z),15(Z)-Octadecadienoic acid prevents increases in mRNA expression of the gene encoding SREBP-1c by the liver X receptor (LXR) agonist T0901317 (Item No. 71810) in HepG2 cells. It reduces LPS-induced increases in nitric oxide (NO) production in RAW 264.7 macrophages when used at a concentration of 10 μM.²

References

1. Nanthirudjanar, T., Furumoto, H., Zheng, J., *et al.* Gut microbial fatty acid metabolites reduce triacylglycerol levels in hepatocytes. *Lipids* **50(11)**, 1093-1102 (2015).
2. Yang, H.-E., Li, Y., Nishimura, A., *et al.* Synthesized enone fatty acids resembling metabolites from gut microbiota suppress macrophage-mediated inflammation in adipocytes. *Mol. Nutr. Food Res.* **61(10)**, (2017).

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, 11/06/2025

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