

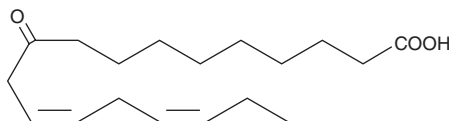
# PRODUCT INFORMATION



## 10-oxo-12(Z),15(Z)-Octadecadienoic Acid

Item No. 36976

**CAS Registry No.:** 1036406-73-1  
**Formal Name:** 10-oxo-12Z,15Z-octadecadienoic acid  
**Synonyms:** 10-oxo-12(Z),15(Z)-18:2,  
10-oxo-12(Z),15(Z)-ODE  
**MF:** C<sub>18</sub>H<sub>30</sub>O<sub>3</sub>  
**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

10-oxo-12(Z),15(Z)-Octadecadienoic acid is an oxylipin. It is produced from  $\alpha$ -linolenic acid (Item Nos. 90210 | 21910) by the gut bacteria *L. plantarum*.<sup>1</sup> Octadecadienoic acid reduces the increase in mRNA expression levels of the gene encoding SREBP-1c induced by the liver X receptor agonist T0901317 (Item No. 71810) in HepG2 cells.

### Reference

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).

#### 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, 01/16/2024

#### 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](http://WWW.CAYMANCHEM.COM)