

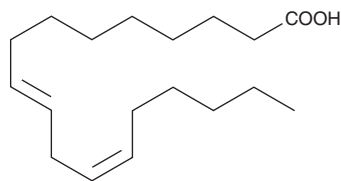
# PRODUCT INFORMATION



## 9(E),12(Z)-Octadecadienoic Acid

Item No. 10005146

CAS Registry No.: 2420-42-0  
Synonyms: C18:2(9E,12Z), C18:2 n-6, FA 18:2  
MF:  $C_{18}H_{32}O_2$   
FW: 280.5  
Purity:  $\geq 95\%$   
Supplied as: A solution in ethanol  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 2$  years



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

### Description

9(E),12(Z)-Octadecadienoic acid is an  $\omega$ -6 polyunsaturated fatty acid and an isomer of linoleic acid (Item Nos. 90150 | 90150.1 | 21909) that contains a *trans* double bond at the C9 position. It has been found as a minor component of bovine milk fat and in partially hydrogenated vegetable oils.<sup>1</sup> 9(E),12(Z)-Octadecadienoic acid levels increase in rabbit meat following supplementation with heated sunflower oil,  $\alpha$ -tocopheryl acetate, and zinc.<sup>2</sup>

### References

1. Precht, D. and Molckentin, J. Trans-geometrical and positional isomers of linoleic acid including conjugated linoleic acid (CLA) in German milk and vegetable fats. *Eur. J. Lipid Sci. Technol.* **99(9)**, 319-326 (1997).
2. Tres, A., Bou, R., Codony, R., *et al.* Oxidized oils and dietary zinc and  $\alpha$ -tocopheryl acetate supplementation: effects on rabbit plasma, liver and meat fatty acid composition and meat Zn, Cu, Fe and Se content. *Animal* **4(11)**, 1929-1939 (2010).

#### 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