

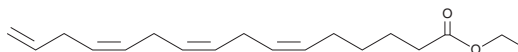
# PRODUCT INFORMATION



## 6(Z),9(Z),12(Z),15(Z)-Hexadecatetraenoic Acid ethyl ester

Item No. 36842

**CAS Registry No.:** 845779-89-7  
**Formal Name:** 6Z,9Z,12Z,15-hexadecatetraenoic acid, ethyl ester  
**Synonyms:** Ethyl 6(Z),9(Z),12(Z)15(Z)-Hexadecatetraenoate, SFE 18:4  
**MF:** C<sub>18</sub>H<sub>28</sub>O<sub>2</sub>  
**FW:** 276.4  
**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

6(Z),9(Z),12(Z)15(Z)-Hexadecatetraenoic Acid ethyl ester is a polyunsaturated fatty ester.<sup>1</sup> Dietary administration of 6(Z),9(Z),12(Z)15(Z)-hexadecatetraenoic acid ethyl ester reduces plasma triglyceride content in mice.

### Reference

1. Hosomi, R., Tanizaki, T., Ikawa, S., *et al.* Effect of 6,9,12,15-hexadecatetraenoic acid (C16:4n-1)-ethyl ester on lipid content and fatty acid composition in the blood and organs of mice. *J. Oleo. Sci.* **70(5)**, 703-712 (2021).

#### 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, 02/13/2024

#### 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