

PRODUCT INFORMATION



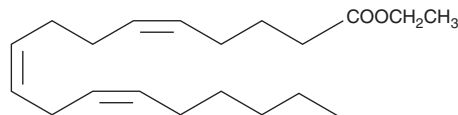
Pinolenic Acid ethyl ester (technical grade)

Item No. 9004488

CAS Registry No.: 493015-74-0

Formal Name: 5Z,9Z,12Z-octadecatrienoic acid, ethyl ester

Synonyms: C18:3 (cis-5,9,12) ethyl ester,
Ethyl all-cis-5,9,12-Octadecatrienoate,
Ethyl Pinolenate,
all-cis-5,9,12-Octadecatrienoic Acid ethyl ester,
all-Z-5,9,12-Octadecatrienoic Acid ethyl ester,
PLA ethyl ester, SFE 20:3



MF: $C_{20}H_{34}O_2$

FW: 306.5

Purity: $\geq 80\%$

Supplied as: A 500 mg/ml solution in ethanol with 0.2% vitamin E stabilizer

Storage: -20°C

Stability: ≥ 1 year

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

Description

Pinolenic acid ethyl ester is a polyunsaturated fatty acid ethyl ester.¹ Oral administration of pinolenic acid ethyl ester (100 mg/kg) decreases blood glucose levels in an oral glucose tolerance test in mice.

Reference

1. Christiansen, E., Watterson, K.R., Stocker, C.J., *et al.* Activity of dietary fatty acids on FFA1 and FFA4 and characterisation of pinolenic acid as a dual FFA1/FFA4 agonist with potential effect against metabolic diseases. *Br. J. Nutr.* **113**(11), 1677-1688 (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, 11/24/2025

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