

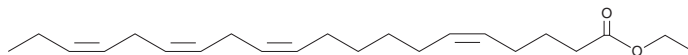
PRODUCT INFORMATION



Juniperonic Acid ethyl ester

Item No. 45623

CAS Registry No.: 143202-57-7
Formal Name: (all-Z)-5,11,14,17-eicosatetraenoic acid, ethyl ester
Synonyms: C20:4 (cis-5,cis-11,cis-14,cis-17) ethyl ester, Ethyl eicosa-5Z,11Z,14Z,17Z-tetraenoate, SFE 20:4(5Z,11Z,14Z,17Z)
MF: C₂₂H₃₆O₂
FW: 332.5
Purity: ≥95%
Supplied as: A 5 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Juniperus communis*



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

Description

Juniperonic acid ethyl ester is a polyunsaturated fatty acid ethyl ester and the esterified form of juniperonic acid. It is formed during the purification process of juniperonic acid from *B. orientalis* seed oil.¹ Juniperonic acid ethyl ester is a neutral, more lipid soluble form of the free acid.

Reference

1. Montañés, F., Tallon, S., and Catchole, O. Isolation of non-methylene interrupted or acetylenic fatty acids from seed oils using semi-preparative supercritical chromatography. *J. Am. Oil. Chem. Soc.* **94**, 981-991 (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, 04/03/2026

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