

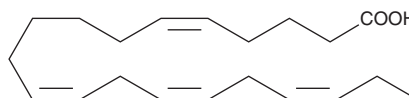
PRODUCT INFORMATION



Juniperonic Acid

Item No. 15198

CAS Registry No.: 18016-45-0
Formal Name: 5Z,11Z,14Z,17Z-eicosatetraenoic acid
Synonyms: C20:4(5Z,11Z,14Z,17Z), C20:4 n-3, FA 20:4
MF: $C_{20}H_{32}O_2$
FW: 304.5
Purity: $\geq 90\%$
Supplied as: A 1 mg/ml 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

Juniperonic acid is an ω -3 polyunsaturated fatty acid that has been found in plant oils, including those of *J. communis*.¹ Juniperonic acid can be biosynthesized from α -linolenic acid (Item Nos. 90210 | 21910) in *C. elegans* and reverses growth reductions and lifespan increases in a *C. elegans* mutant lacking Δ^6 -desaturase that cannot biosynthesize certain fatty acids, such as arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) and ω -3 arachidonic acid (Item No. 90011).² It reduces bombesin-induced proliferation of 3T3-Swiss mouse fibroblasts when used at a concentration of 50 μM .³

References

1. Güvenç, A., Küçükboyacı, N., and Gören, A.C. Fatty acid composition of *Juniperus* species (*Juniperus section*) native to Turkey. *Nat. Prod. Commun.* **7**(7), 919-922 (2012).
2. Guha, S., Calarco, S., Gachet, M.S., et al. Juniperonic acid biosynthesis is essential in *Caenorhabditis elegans* lacking Δ^6 desaturase (*fat-3*) and generates new ω -3 endocannabinoids. *Cells* **9**(9), 2127 (2020).
3. Morishige, J., Amano, N., Hirano, K., et al. Inhibitory effect of juniperonic acid (Δ -5c,11c,14c,17c-20:4, ω -3) on bombesin-induced proliferation of Swiss 3T3 cells. *Biol. Pharm. Bull.* **31**(9), 1786-1789 (2008).

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, 04/23/2026

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