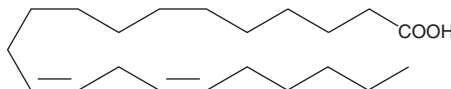


PRODUCT INFORMATION

12(Z),15(Z)-Heneicosadienoic Acid

Item No. 22594

CAS Registry No.: 191545-08-1
 Formal Name: 12Z,15Z-heneicosadienoic acid
 Synonyms: C21:2, C21:2n-6,9, FA 21:2
 MF: $C_{21}H_{38}O_2$
 FW: 322.5
 Purity: $\geq 98\%$
 Supplied as: A 10 mg/ml solution in ethanol
 Storage: -20°C
 Stability: ≥ 4 years



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

Laboratory Procedures

12(Z),15(Z)-Heneicosadienoic acid is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 12(Z),15(Z)-heneicosadienoic acid in these solvents is approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of 12(Z),15(Z)-heneicosadienoic acid is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 12(Z),15(Z)-heneicosadienoic acid in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

12(Z),15(Z)-Heneicosadienoic acid is an ω -6 very long-chain polyunsaturated fatty acid. It is a positional isomer of heneicosadienoic acid (Item No. 22593) that is rare in living organisms.¹

Reference

1. Kawashima, H., Ohnishi, M., Ogawa, S., *et al.* Unusual fatty acid isomers of triacylglycerols and polar lipids in female limpet gonads of *Cellana grata*. *Lipids* **43**(6), 559-567 (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, 03/25/2026

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