

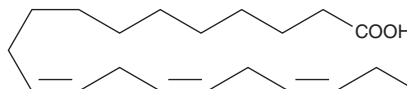
PRODUCT INFORMATION



11(Z),14(Z),17(Z)-Eicosatrienoic Acid

Item No. 90192

CAS Registry No.: 17046-59-2
Formal Name: 11Z,14Z,17Z-eicosatrienoic acid
Synonyms: Dihomo- α -Linolenic Acid, FA 20:3
MF: C₂₀H₃₄O₂
FW: 306.5
Purity: \geq 98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

11(Z),14(Z),17(Z)-Eicosatrienoic 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 11(Z),14(Z),17(Z)-eicosatrienoic acid in these solvents is approximately 100 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 11(Z),14(Z),17(Z)-eicosatrienoic acid is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. For maximum aqueous solubility, 11(Z),14(Z),17(Z)-eicosatrienoic acid can be directly dissolved in 0.1 M Na₂CO₃ (1.7 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. We do not recommend storing the aqueous solution for more than one day.

Description

Eicosatrienoic Acid (20:3 ω -3) is a rare polyunsaturated fatty acid of the ω -3 series. In normal humans, it represents less than 0.25% of serum phospholipid fatty acids. However, it is one of the most active essential fatty acids when assayed for the inhibition of fatty acid elongation/desaturation reactions which convert dietary C-18 fatty acids to C-20 eicosanoid precursors.¹

References

1. Holman, R.T. Control of polyunsaturated acids in tissue lipids. *J. Am. Coll. Nutr.* **5**, 183-211 (1986).

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, 02/12/2024

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