

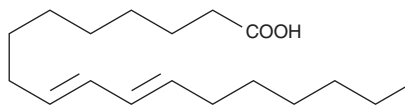
PRODUCT INFORMATION



9(E),11(E)-Conjugated Linoleic Acid

Item No. 90370

CAS Registry No.: 544-71-8
Formal Name: 9E,11E-octadecadienoic acid
Synonyms: 9E,11E-CLA, FA 18:2, Isolinoleic Acid, Mangold's acid
MF: C₁₈H₃₂O₂
FW: 280.5
Purity: ≥98%
UV/Vis.: λ_{max}: 233 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

9(E),11(E)-CLA is supplied as a crystalline solid. A stock solution may be made by dissolving the 9(E),11(E)-CLA in the solvent of choice, which should be purged with an inert gas. 9(E),11(E)-CLA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 9(E),11(E)-CLA 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. Organic solvent-free aqueous solutions of 9(E),11(E)-CLA can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 9(E),11(E)-CLA in 0.15 M Tris-HCl buffer (pH 8.5) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

9(E),11(E)-CLA is the 9,11 all-*trans* isomer of linoleic acid. It is a component of CLAs, anticarcinogenic fatty acids found predominantly in animal food sources. CLA was originally identified in ground beef, but it is also present in a variety of dairy products. CLA is effective at reducing mammary tumors in rats at levels as low as 0.1% by weight of their diet.¹

Reference

1. Ip, C., Scimeca, J.A., and Thompson, H.J. Conjugated linoleic acid. *Cancer* **74**(3), 1050-1054 (1994).

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, 02/13/2024

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