

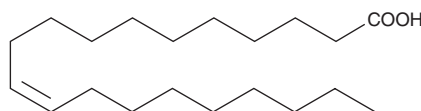
PRODUCT INFORMATION



11(Z)-Eicosenoic Acid

Item No. 20606

CAS Registry No.: 5561-99-9
Formal Name: 11Z-eicosenoic acid
Synonyms: cis-11-Eicosenoic Acid, FA 20:1
MF: C₂₀H₃₈O₂
FW: 310.5
Purity: ≥95%
Supplied as: A 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.

Laboratory Procedures

11(Z)-Eicosenoic 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 11(Z)-eicosenoic 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 11(Z)-eicosenoic acid is needed, it can be prepared by evaporating the 11(Z)-eicosenoic acid and directly dissolving the neat oil in aqueous buffers. The solubility of 11(Z)-eicosenoic acid in PBS (pH 7.2) and 0.15 M tris-HCl (pH 8.5) is approximately 0.15 and 1 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

11(Z)-Eicosenoic acid is one of several monounsaturated 20-carbon fatty acids. The combined C20:1 isomers constitute 70% of the total fatty acid pool in jojoba seed oil isolated from plants in the Arizona desert.¹

Reference

1. Miwa, T.K. Jojoba oil wax esters and derived fatty acids and alcohols: Gas chromatographic analyses. *J. Am. Chem. Soc.* **48(6)**, 259-264 (1971).

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, 01/24/2024

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