

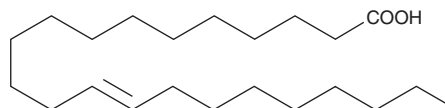
PRODUCT INFORMATION



13(E)-Docosenoic Acid

Item No. 22580

CAS Registry No.: 506-33-2
Formal Name: 13E-docosenoic acid
Synonyms: Brassidic Acid, *trans*-13-Docosenoic Acid, FA 22:1, NSC 59686
MF: C₂₂H₄₂O₂
FW: 338.6
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

13(E)-Docosenoic acid is supplied as a neat oil. A stock solution may be made by dissolving the 13(E)-docosenoic acid in the solvent of choice, which should be purged with an inert gas. 13(E)-Docosenoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 13(E)-docosenoic acid in ethanol is approximately 10 mg/ml and approximately 100 mg/ml in DMSO and DMF.

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 13(E)-docosenoic acid can be prepared by directly dissolving the neat oil in aqueous buffers. The solubility of 13(E)-docosenoic acid in PBS (pH 7.2) is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

13(E)-Docosenoic acid is a 22-carbon monounsaturated fatty acid. It is the *trans* isomer of 13(Z)-docosenoic acid (Item No. 90175).

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, 01/25/2024

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