

# PRODUCT INFORMATION



## 13(Z)-Docosenoic Acid (sodium salt)

Item No. 38975

CAS Registry No.: 14488-85-8

Synonyms: C22:1(13Z), C22:1 n-9,  
cis-13-Docosenoic Acid,  
Erucic Acid, FA 22:1,  
Sodium Erucate

MF:  $C_{22}H_{41}O_2 \cdot Na$

FW: 360.6

Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

13(Z)-Docosenoic acid (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 13(Z)-docosenoic acid (sodium salt) in the solvent of choice, which should be purged with an inert gas. 13(Z)-Docosenoic acid (sodium salt) is slightly soluble in ethanol.

### Description

13(Z)-Docosenoic acid is a 22-carbon monounsaturated fatty acid. Dietary administration of 13(Z)-docosenoic acid (6.6% w/w) increases weight gain and heart triglyceride levels in calcium-deprived weanling rats.<sup>1</sup>

### Reference

1. Hulan, H.W., Kramer, J.K.G., Mahadevan, S., *et al.* Relationship between erucic acid and myocardial changes in male rats. *Lipids* **11**(1), 9-15 (1975).

#### 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, 02/13/2025

#### 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