

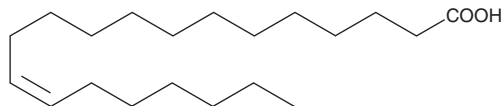
PRODUCT INFORMATION



13(Z)-Eicosenoic Acid

Item No. 26169

CAS Registry No.: 17735-94-3
Synonyms: C20:1(13Z), C20:1 *n*-7,
cis-13-Eicosenoic Acid, FA 20:1,
Paullinic Acid
MF: C₂₀H₃₈O₂
FW: 310.5
Purity: ≥95%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

13(Z)-Eicosenoic acid is supplied as a neat solid. A stock solution may be made by dissolving the 13(Z)-eicosenoic acid in the solvent of choice. 13(Z)-Eicosenoic acid is soluble in the organic solvent ethanol, which should be purged with an inert gas, at a concentration of approximately 20 mg/ml.

13(Z)-Eicosenoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 13(Z)-eicosenoic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 13(Z)-Eicosenoic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

13(Z)-Eicosenoic acid is an ω-7 fatty acid found in a variety of fish from the Indian, Atlantic, and Pacific oceans.¹ It increases triglyceride accumulation in 3T3-L1 cells when used at a concentration of 50 μM.²

References

1. Senarath, S., Yoshinaga, K., Nagai, T., *et al.* Quantitative analysis of the distribution of *cis*-eicosenoic acid positional isomers in marine fishes from the Indian ocean. *J. Oleo. Sci.* **66(2)**, 187-197 (2017).
2. Senarath, S., Yoshinaga, K., Nagai, T., *et al.* Differential effect of *cis*-eicosenoic acid positional isomers on adipogenesis and lipid accumulation in 3T3-L1 cells. *Eur. J. Lipid Sci. Technol.* **120(6)**, 1700512 (2018).

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, 03/06/2024

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