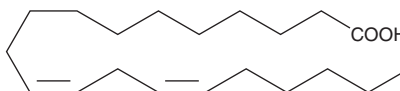


PRODUCT INFORMATION

11(Z),14(Z)-Eicosadienoic Acid

Item No. 90330

CAS Registry No.: 5598-38-9
 Formal Name: 11Z,14Z-eicosadienoic acid
 MF: $C_{20}H_{36}O_2$
 FW: 308.5
 Purity: $\geq 98\%$
 Supplied as: A solution in ethanol
 Storage: -20°C
 Stability: ≥ 2 years
 Special Conditions: Oxygen and light sensitive



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

Laboratory Procedures

11(Z),14(Z)-Eicosadienoic 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 11(Z),14(Z)-Eicosadienoic acid in these solvents is approximately 100 mg/ml.

11(Z),14(Z)-Eicosadienoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 11(Z),14(Z)-eicosadienoic acid should be diluted with the aqueous buffer of choice. The solubility of 11(Z),14(Z)-eicosadienoic acid in PBS (pH 7.2) is approximately 100 $\mu\text{g/ml}$. We do not recommend storing the aqueous solution for more than one day.

Description

11(Z),14(Z)-Eicosadienoic acid is an uncommon, naturally occurring PUFA. It inhibits the binding of $[^3\text{H}]\text{-LTB}_4$ to pig neutrophil membranes with a K_i of 3 μM .¹

Reference

1. Yagaloff, K.A., Franco, L., Simko, B., *et al.* Essential fatty acids are antagonists of the leukotriene B_4 receptor. *Prostaglandins, Leukot. Essent. Fatty Acids* **52(5)**, 293-297 (1995).

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, 06/13/2023

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