

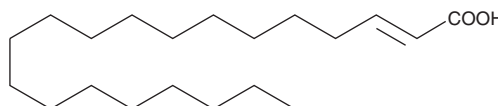
# PRODUCT INFORMATION



## $\Delta^2$ -trans Eicosenoic Acid

Item No. 10007622

CAS Registry No.: 2127095-72-9  
Formal Name: 2E-eicosenoic acid  
Synonym: FA 20:1  
MF:  $C_{20}H_{38}O_2$   
FW: 310.5  
Purity:  $\geq 98\%$   
Supplied as: A crystalline solid  
Storage:  $-20^\circ\text{C}$   
Stability:  $\geq 2$  years



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

### Laboratory Procedures

$\Delta^2$ -trans Eicosenoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the  $\Delta^2$ -trans eicosenoic acid in an organic solvent purged with an inert gas.  $\Delta^2$ -trans Eicosenoic acid is soluble in organic solvents such as ethanol. The solubility of  $\Delta^2$ -trans eicosenoic acid in this solvent is approximately 20 mg/ml.

$\Delta^2$ -trans Eicosenoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers,  $\Delta^2$ -trans eicosenoic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice.  $\Delta^2$ -trans 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

$\Delta^2$ -trans Eicosenoic acid is an  $\alpha,\beta$ -unsaturated fatty acid that is a bi-product of the synthesis of  $\Delta^2$ -cis eicosenoic acid.  $\Delta^2$ -cis Eicosenoic acid and its salts have potential medical use for treating diabetes and improving lipid metabolism. A related compound, 2-octadecenoic acid, has been shown to improve liver function and decrease blood sugar in streptozocin-induced diabetic rats.<sup>1</sup>

### Reference

1. Iwamura, J., Hayashi, C., and Nozomu, T. 1984. Medicinal use of  $\alpha,\beta$ -unsaturated fatty acids. U.S. Patent 4,472,432, filed May 12, 1982, and issued Sept. 18, 1984.

#### 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, 03/13/2024

#### 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