

# PRODUCT INFORMATION



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

Item No. 10007602

**CAS Registry No.:** 93772-82-8  
**Formal Name:** 2Z-eicosenoic acid  
**Synonym:** FA 20:1  
**MF:** C<sub>20</sub>H<sub>38</sub>O<sub>2</sub>  
**FW:** 310.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 209 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

$\Delta^2$ -cis Eicosenoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the  $\Delta^2$ -cis eicosenoic acid in the solvent of choice, which should be purged with an inert gas.  $\Delta^2$ -cis Eicosenoic acid is soluble in the organic solvent ethanol at a concentration of approximately 20 mg/ml

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

$\Delta^2$ -cis Eicosenoic acid is an  $\alpha,\beta$ -unsaturated fatty acid that has been extracted from fresh water clams and purified. 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>  $\Delta^2$ -cis Eicosenoic acid and its salts have potential medicinal use for treating diabetes and improving lipid metabolism.

### Reference

1. Iwamura, J., Hayashi, C., and Nozomu, T. Medicinal use of  $\alpha,\beta$ -unsaturated fatty acids. *Inabata & Co., Ltd.* 4,472,432 (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, 03/28/2024

#### 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