

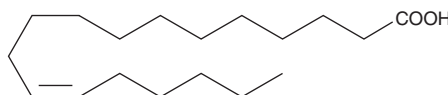
# PRODUCT INFORMATION



## cis-12-Octadecenoic Acid

Item No. 43580

**CAS Registry No.:** 13126-37-9  
**Formal Name:** 12Z-octadecenoic acid  
**Synonyms:** 18:1(12Z), FA 18:1  
**MF:** C<sub>18</sub>H<sub>34</sub>O<sub>2</sub>  
**FW:** 282.5  
**Purity:** ≥98%  
**Supplied as:** A 50 mg/ml solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Description

cis-12-Octadecenoic acid is acid is a monounsaturated fatty acid that has been found in bovine milk.<sup>1</sup>

### Reference

1. Antunes, I., Bexiga, R., Pinto, C., *et al.* Lipid profile of plant-based milk alternatives (PBMA) and cow's milk: A comparison. *J. Agric. Food Chem.* **72(32)**, 18110-18120 (2024).

#### 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, 11/21/2025

#### 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