

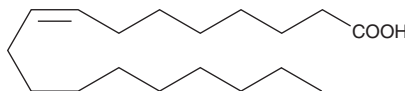
PRODUCT INFORMATION



cis-8-Octadecenoic Acid

Item No. 27447

CAS Registry No.: 5684-71-9
Formal Name: 8Z-octadecenoic acid
Synonyms: C18:1(8Z), *cis*-8-Octadecenoate, Δ^8 -*cis*-Octadecenoic Acid, FA 18:1
MF: C₁₈H₃₄O₂
FW: 282.5
Purity: ≥98%
Supplied as: A 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.

Laboratory Procedures

cis-8-Octadecenoic 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 *cis*-8-octadecenoic acid in these solvents is approximately 30 mg/ml.

cis-8-Octadecenoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of *cis*-8-octadecenoic acid should be diluted with the aqueous buffer of choice. *cis*-8-Octadecenoic acid has a solubility of approximately 0.25 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method.

Description

cis-8-Octadecenoic acid is a monounsaturated fatty acid and an isomer of oleic acid (Item Nos. 90260 | 24659), *trans*-vaccenic acid (Item No. 15301), *trans*-petroselinic acid (Item No. 20026), and *cis*-petroselinic acid (Item No. 20024). It has been found in partially hydrogenated vegetable oil and milk fat.¹

Reference

1. Yoshinaga, K., Asanuma, M., Mizobe, H., *et al.* Characterization of *cis*- and *trans*-octadecenoic acid positional isomers in edible fat and oil using gas chromatography-flame ionisation detector equipped with highly polar ionic liquid capillary column. *Food Chem.* **160**, 39-45 (2014).

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, 03/12/2024

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