

PRODUCT INFORMATION



Palmitoleic Acid methyl ester

Item No. 20605

CAS Registry No.: 1120-25-8

Formal Name: (9Z)-hexadecenoic acid, methyl ester

Synonyms: Methyl *cis*-9-Hexadecenoate,
Methyl Palmitoleate

MF: C₁₇H₃₂O₂

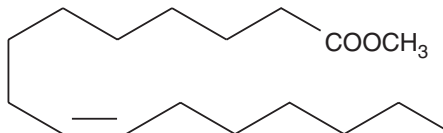
FW: 268.4

Purity: ≥98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

Palmitoleic acid methyl ester 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 ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of palmitoleic acid methyl ester in ethanol is approximately 50 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

Palmitoleic acid methyl ester is an ester version of palmitoleic acid (Item No. 10009871), an ω-7 monounsaturated fatty acid that is a common constituent of the triglycerides of human adipose tissue. Palmitoleic acid-based diets raise low-density lipoprotein (LDL) cholesterol and lower high-density lipoprotein (HDL) cholesterol, even when dietary intake of cholesterol is maintained at a low level.¹ Methyl esters of fatty acids are commonly used for the formulation of fatty acid-containing diets and dietary supplements.

Reference

1. Nestel, P., Clifton, P., and Noakes, M. Effects of increasing dietary palmitoleic acid compared with palmitic and oleic acids on plasma lipids of hypercholesterolemic men. *J. Lipid Res.* **35**, 656-662 (1994).

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, 03/07/2017

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