

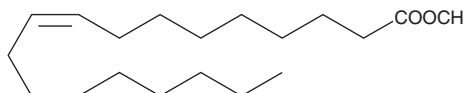
PRODUCT INFORMATION



Oleic Acid methyl ester

Item No. 20604

CAS Registry No.: 112-62-9
Formal Name: 9Z-octadecenoic acid, methyl ester
Synonyms: Methyl *cis*-9-Octadecenoate,
Methyl Oleate,
cis-9-Octadecenoic Acid methyl ester,



MF: C₁₉H₃₆O₂

FW: 296.5

Purity: ≥98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: As supplied, 1 year from the QC date provided on the Certificate of Analysis, when stored properly

Laboratory Procedures

Oleic acid methyl ester is supplied as a solution in ethanol. To change the solvent, evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO or dimethyl formamide purged with an inert gas can be used. The solubility of oleic acid methyl ester in these solvents is at least 100 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of oleic acid methyl ester is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of oleic acid methyl ester in PBS, pH 7.2 and 0.15 M Tris-HCl, pH 8.5, is less than 0.1 and at least 1 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Oleic acid methyl ester is an esterified version of the free acid which is less water soluble but more amenable for the formulation of oleate-containing diets and dietary supplements. Oleic acid (Item No. 90260) is a monounsaturated fatty acid and is one of the major components of membrane phospholipids. Oleic acid contributes about 17% of the total fatty acids esterified to phosphatidylcholine, the major phospholipid class in porcine platelets.¹

Reference

1. Wahle, K.W.J. and Peacock, L.I.L. Effects of isomeric *cis* and *trans* eighteen carbon monounsaturated fatty acids on porcine platelet function. *Biochim. Biophys. Acta* **1301**, 141-149 (1996).

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, 04/12/2017

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