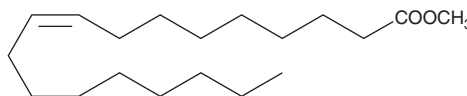


PRODUCT INFORMATION

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,
 SFE 19:1
MF: C₁₉H₃₆O₂
FW: 296.5
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

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.

Oleic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of oleic acid methyl ester should be diluted with the aqueous buffer of choice. The solubility of oleic acid methyl ester in PBS, pH 7.2 and 0.15 M Tris-HCl, pH 8.5, is approximately 0.1 and 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, 04/02/2024

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