

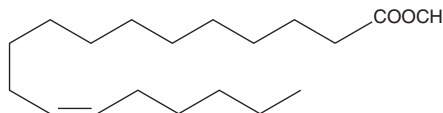
PRODUCT INFORMATION



12(Z)-Octadecenoic Acid methyl ester

Item No. 10010180

CAS Registry No.: 2733-86-0
Formal Name: 12Z-octadecenoic acid, methyl ester
Synonym: SFE 19:1
MF: C₁₉H₃₆O₂
FW: 296.5
Purity: ≥96%
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

12(Z)-Octadecenoic 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 12(Z)-octadecenoic acid methyl ester in these solvents is approximately 20 and 15 mg/ml, respectively.

12(Z)-Octadecenoic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 12(Z)-octadecenoic acid methyl ester should be diluted with the aqueous buffer of choice. The solubility of 12(Z)-octadecenoic acid methyl ester in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

cis-12-Octadecenoic acid is a monounsaturated C-18 fatty acid. It is a positional isomer of oleic acid. The methyl ester is a more lipid soluble form of the free acid that can be used as a standard for analysis.

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, 02/15/2024

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