

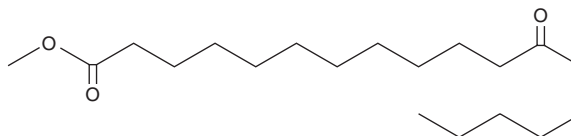
PRODUCT INFORMATION



12-oxo Stearic Acid methyl ester

Item No. 10007012

CAS Registry No.: 2380-27-0
Formal Name: 12-oxo-octadecanoic acid, methyl ester
Synonyms: Methyl 12-oxo Octadecanoate,
Methyl 12-keto Stearate,
Methyl 12-oxo Stearate
MF: C₁₉H₃₆O₃
FW: 312.5
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

12-oxo Stearic acid methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the 12-oxo stearic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 12-oxo Stearic acid methyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 12-oxo stearic acid methyl ester in these solvents is approximately 50 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. Organic solvent-free aqueous solutions of 12-oxo stearic acid methyl ester can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 12-oxo stearic acid methyl ester in PBS, pH 7.2, is approximately 0.15 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

12-oxo Stearic acid methyl ester is an esterified form of 12-oxo stearic acid. It has been found in esterified *L. chinensis* and *L. decipiens* pulp oils and serves as a measure of 12-oxo stearic acid content.¹

Reference

1. Kadry, H., Shoala, S., El Gindi, O., *et al.* Chemical characterization of the lipophilic fraction of *Livistona decipiens* and *Livistona Chinensis* fruit Pulps (palmae) and assessment of their anti-hyperlipidemic and anti-ulcer activities. *Nat. Prod. Commun.* **4(2)**, 265-270 (2009).

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, 10/24/2022

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