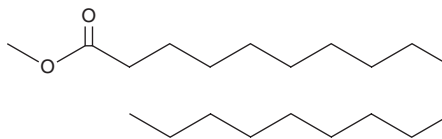


PRODUCT INFORMATION

Nonadecanoic Acid methyl ester

Item No. 20607

CAS Registry No.: 1731-94-8
Formal Name: nonadecanoic acid, methyl ester
Synonyms: Methyl Nonadecanoate, n-Nonadecanoic Acid methyl ester, SFE 20:0
MF: C₂₀H₄₀O₂
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

Nonadecanoic acid methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the nonadecanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Nonadecanoic acid methyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of nonadecanoic acid methyl ester in ethanol and DMF is approximately 25 mg/ml and approximately 10 mg/ml in DMSO.

Nonadecanoic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nonadecanoic acid methyl ester should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Nonadecanoic acid methyl ester has a solubility of approximately 0.25 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Nonadecanoic acid methyl ester is an esterified version of the free acid which is less water soluble but more amenable for the formulation of nonadecanoate-containing diets and dietary supplements. Nonadecanoic acid (Item No. 19723) is a 19-carbon saturated fatty acid that is found in fats and vegetable oils. It has been shown to inhibit HL-60 cancer cell proliferation with an IC₅₀ value of 68 μM.¹

Reference

1. Fukuzawa, M., Yamaguchi, R., Hide, I., et al. Possible involvement of long chain fatty acids in the spores of *Ganoderma lucidum* (Reishi Houshi) to its anti-tumor activity. *Biol. Pharm. Bull.* **31(10)**, 1933-1937 (2008).

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, 03/24/2024

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