

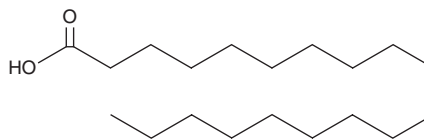
PRODUCT INFORMATION



Nonadecanoic Acid

Item No. 19723

CAS Registry No.: 646-30-0
Formal Name: nonadecanoic acid
Synonyms: FA 19:0, Nonadecanoate, n-Nonadecanoic Acid, Nonadecylic Acid, NSC 11914
MF: C₁₉H₃₈O₂
FW: 298.5
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 is supplied as a crystalline solid. A stock solution may be made by dissolving the nonadecanoic acid in the solvent of choice, which should be purged with an inert gas. Nonadecanoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of nonadecanoic acid in ethanol and DMF is approximately 25 mg/ml and approximately 10 mg/ml in DMSO.

Nonadecanoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nonadecanoic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Nonadecanoic acid 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 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, 02/08/2024

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