

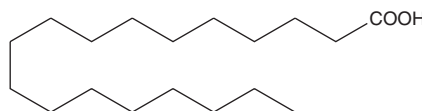
PRODUCT INFORMATION



Stearic Acid

Item No. 10011298

CAS Registry No.: 57-11-4
Formal Name: octadecanoic acid
Synonyms: C18:0, Octadecanoic Acid
MF: C₁₈H₃₆O₂
FW: 284.5
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Stearic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the stearic acid in an organic solvent purged with an inert gas. Stearic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of stearic acid in these solvents is approximately 20, 10, and 30 mg/ml, respectively.

Description

Stearic acid is a long-chain saturated fatty acid. It is a major component of cocoa butter and has also been found in beef fat and vegetable oils.¹⁻³ Unlike many long-chain saturated fatty acids, dietary stearic acid does not induce hypercholesterolemia or raise LDL-cholesterol.⁴

References

1. Chuparova, E., Chobanov, D., and Popov, A. Quantitative analysis of fatty acids by liquid-partition chromatography. *Izv. Inst. Org. Khim. Bulgar. Akad. Nauk* **2**, 31-35 (1965).
2. Westerling, D.B. and Hedrick, H.B. Fatty acid composition of bovine lipids as influenced by diet, sex and anatomical location and relationship to sensory characteristics. *J. Anim. Sci.* **48(6)**, 1343-1348 (1979).
3. Demirbaş, A. Chemical and fuel properties of seventeen vegetable oils. *Energy Sources* **25(7)**, 721-728 (2003).
4. Grundy, S.M. Influence of stearic acid on cholesterol metabolism relative to other long-chain fatty acids. *Am. J. Clin. Nutr.* **60(6 Suppl.)**, 986S-990S (1994).

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, 07/12/2019

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