

# PRODUCT INFORMATION



## Heptadecanoic Acid

Item No. 19722

CAS Registry No.: 506-12-7

Synonyms: C17:0, FA 17:0, n-Heptadecoic Acid, Heptadecylic Acid, Margaric Acid, NSC 3743

MF:  $C_{17}H_{34}O_2$

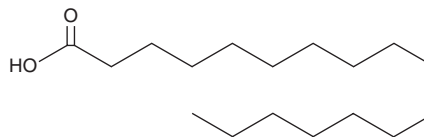
FW: 270.5

Purity:  $\geq 95\%$

Supplied as: A crystalline solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

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

Heptadecanoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, heptadecanoic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Heptadecanoic 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

Heptadecanoic acid is an odd-chain saturated fatty acid that contains seventeen carbons and has been found in milk fat.<sup>1</sup> Heptadecanoic acid has been used as an internal standard for the quantification of fatty acids in human plasma by LC- and GC-MS and as a biomarker for dairy fat intake.<sup>1,2</sup>

### References

1. Jenkins, B., West, J.A., and Koulman, A. A review of odd-chain fatty acid metabolism and the role of pentadecanoic acid (c15:0) and heptadecanoic acid (c17:0) in health and disease. *Molecules* **20**(2), 2425-2444 (2015).
2. Yakoob, M.Y., Shi, P., Hu, F.B., et al. Circulating biomarkers of dairy fat and risk of incident stroke in U.S. men and women in 2 large prospective cohorts. *Am. J. Clin. Nutr.* **100**(6), 1437-1447 (2014).

#### 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