

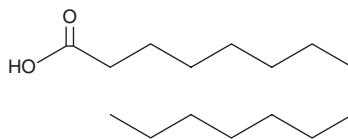
PRODUCT INFORMATION



Pentadecanoic Acid

Item No. 17399

CAS Registry No.: 1002-84-2
Formal Name: Pentadecanoic acid
Synonyms: C15:0, FA 15:0, NSC 28486, Pentadecylic Acid
MF: $C_{15}H_{30}O_2$
FW: 242.4
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

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

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

Pentadecanoic acid is a 15:0 saturated fatty acid found in esterified form in the lipids of many bacterial species and in the milk fat of ruminants. It has been used as a biological marker for the intake of dairy fat in the assessment of metabolic risk factors.^{1,2}

References

1. Smedman, A.E.M., Gustafsson, I.-B., Berglund, L.G.T., *et al.* Pentadecanoic acid in serum as a marker for intake of milk fat: Relations between intake of milk fat and metabolic risk factors. *Am. J. Clin. Nutr.* **69**(1), 22-29 (1999).
2. Wolk, A., Vessby, B., Ljung, H., *et al.* Evaluation of a biological marker of dairy fat intake. *Am. J. Clin. Nutr.* **68**(2), 291-295 (1998).

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