

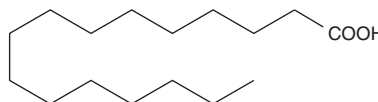
PRODUCT INFORMATION



Palmitic Acid

Item No. 10006627

CAS Registry No.: 57-10-3
Formal Name: hexadecanoic acid
Synonyms: C16:0, FA 16:0, Cetylic Acid, Hexadecanoic Acid, NSC 5030
MF: C₁₆H₃₂O₂
FW: 256.4
Purity: ≥98%
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

Palmitic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the palmitic acid in the solvent of choice, which should be purged with an inert gas. Palmitic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of palmitic acid in ethanol is approximately 30 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Palmitic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, palmitic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Palmitic acid has a solubility of approximately 0.25 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Palmitic acid is a common 16-carbon saturated fat that represents 10-20% of human dietary fat intake and comprises approximately 25 and 65% of human total plasma lipids and saturated fatty acids, respectively.^{1,2} Acylation of palmitic acid to proteins facilitates anchoring of membrane-bound proteins to the lipid bilayer and trafficking of intracellular proteins, promotes protein-vesicle interactions, and regulates various G protein-coupled receptor functions.¹ Palmitic acid (200 μM) increases NF-κB p65 levels, matrix metalloproteinase-9 (MMP-9) activity, and production of reactive oxygen species (ROS) in AsPC-1 pancreatic cancer cells, as well as increases migration of AsPC-1 cells.³ It increases COX-2 levels in RAW 264.7 cells and increases LPS-induced IL-1β levels and caspase-1 activity in isolated mouse peritoneal macrophages.^{4,5} Dietary administration of palmitic acid (2.2% w/w for 12 weeks) increases mouse hippocampal β-secretase 1 (BACE1) activity and amyloid β (1-42) (Aβ42; Item No. 20574) levels.⁶ It also induces systolic contractile dysfunction in isolated mouse hearts.⁷ Red blood cell palmitic acid levels are increased in patients with metabolic syndrome compared to patients without metabolic syndrome and are also increased in the plasma of patients with type 2 diabetes compared to individuals without diabetes.^{8,9}

References

1. Fatima, S., Hu, X., Gong, R.-H., et al. *Cell Mol. Life Sci.* **76(13)**, 2547-2557 (2019).
2. Santos, M.J., López-Jurado, M., Llopis, J., et al. *Ann. Nutr. Metab.* **39(1)**, 52-62 (1995).
3. Binker-Cosen, M.J., Richards, D., Oliver, B., et al. *Biochem. Biophys. Res. Commun.* **484(1)**, 152-158 (2017).
4. Lee, J.Y., Sohn, K.H., Rhee, S.H., et al. *J. Biol. Chem.* **276(20)**, 16683-16689 (2001).
5. Karasawa, T., Kawashima, A., Usui-Kawanishi, F., et al. *Arterioscler. Thromb. Vasc. Biol.* **38(4)**, 744-756 (2018).
6. Marwarha, G., Rostad, S., Lilek, J., et al. *J. Alzheimers Dis.* **57(3)**, 907-925 (2017).
7. Knowles, C.J., Cebova, M., and Pinz, I.M. *J. PLoS One* **8(4)**, e61369 (2013).
8. Yi, L.-Z., He, J., Liang, Y.-Z., et al. *FEBES J.* **580(3)**, 6837-6845 (2006).
9. Kabagambe, E.K., Tsai, M.Y., Hopkins, P.N., et al. *Clin. Chem.* **54(1)**, 154-162 (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, 03/11/2024

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