

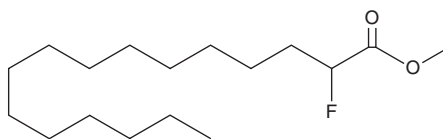
PRODUCT INFORMATION



2-fluoro Palmitic Acid methyl ester

Item No. 24526

CAS Registry No.: 137676-82-5
Formal Name: 2-fluoro-hexadecanoic acid, methyl ester
Synonyms: Methyl 2-fluoro Palmitate,
Methyl 2-fluoro Hexadecanoate
MF: C₁₇H₃₃FO₂
FW: 288.4
Purity: ≥98%
Supplied as: A 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

2-fluoro Palmitic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 2-fluoro palmitic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 2-fluoro Palmitic acid methyl ester is soluble in organic solvents such as chloroform, methanol, and ethanol.

Description

2-fluoro Palmitic acid methyl ester is a monofluorinated form of palmitic acid methyl ester (Item No. 10007358). It is an intermediate in the synthesis of 2-iodoaldehydes, monofluorinated fatty acids, and polyfluoro ketones.¹⁻³

References

1. Panneels, V., Van den Bergen, H., Jacoby, C., *et al.* Inhibition of H₂O₂ production by iodoaldehydes in cultured dog thyroid cells. *Mol. Cell Endocrinol.* **102(1-2)**, 167-176 (1994).
2. Delgado, A., Ruiz, M.A., Camps, F., *et al.* Synthesis of potential inhibitors of the biosynthesis of the sex pheromone of *Spodoptera littoralis*. Part I: Monofluorinated fatty acids. *Chem. Phys. Lipids* **59(2)**, 127-135 (1991).
3. Baskakis, C., Magrioti, V., Cotton, N., *et al.* Synthesis of polyfluoro ketones for selective inhibition of human phospholipase A₂ enzymes. *J. Med. Chem.* **51(24)**, 8027-8037 (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, 11/14/2022

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