

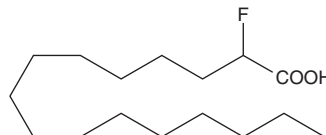
PRODUCT INFORMATION



2-fluoro Palmitic Acid

Item No. 90380

CAS Registry No.: 16518-94-8
Formal Name: 2-fluorohexadecanoic acid
Synonyms: 2-fluoro Hexadecanoate,
2-fluoro Palmitate
MF: C₁₆H₃₁FO₂
FW: 274.4
Purity: ≥95%
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

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

Description

2-fluoro Palmitic acid inhibits sphingosine biosynthesis and long-chain acyl-CoA synthetase with an IC₅₀ value of 0.2 mM.¹

Reference

1. Soltysiak, R.M., Matsuura, F., Bloomer, D., *et al.* D,L- α -Fluoropalmitic acid inhibits sphingosine base formation and accumulates in membrane lipids of cultured mammalian cells. *Biochimica et Biophysica Acta* **792**, 214-226 (1984).

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, 05/02/2024

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