

PRODUCT DATA SHEET

Methyl 2-fluoropalmitate

Catalog number: 1718

Activity: Inactive ester of 2-fluoropalmitic acid

Source: synthetic

Solubility: chloroform, ethanol, methanol

CAS number: 137676-82-5

Molecular Formula: C₁₇H₃₃FO₂

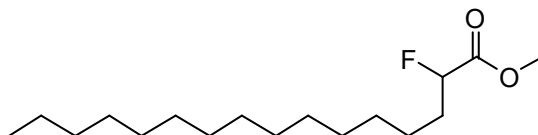
Molecular Weight: 288

Storage: -20C

Purity: TLC: >98%; identity confirmed by MS

TLC System: hexane/ethyl ether (70:30)

Appearance: solid



Application Notes:

This product is the inactive methyl ester of the acyl-CoA synthase inhibitor 2-fluoropalmitic acid. 2-Fluoropalmitic acid is a synthetic inhibitor of palmitoyl-CoA formation by long chain acyl-CoA synthetase thereby inhibiting sphingosine biosynthesis and protein palmitoylation. This product is very useful in the investigation of sphingosine synthesis¹, protein acylation², and other biological mechanisms³. Other *alpha*-fluoro-fatty acids also inhibit their respective acyl-CoAs.⁴

Selected References:

1. R. Soltysiak et al. "D,L-*alpha*-Fluoropalmitic acid inhibits sphingosine base formation and accumulates in membrane lipids of cultured mammalian cells" *Biochim Biophys Acta*, Vol. 792(2) pp. 214-226, 1984
2. J. Zhang et al. "Novel bimodal effects of the G-protein tissue transglutaminase on adrenoreceptor signaling" *Journal of Biochemistry*, Vol. 343, 541-549, 1999
3. G. DeJesus and O. Bizzozero "Effect of 2-Fluoropalmitate, Cerulenin and Tunicamycin on the Palmitoylation and Intracellular Translocation of Myelin Proteolipid Protein" *Neurochemical Research*, Vol. 27(12) pp. 1669-1675, 2002
4. M. Grillo et al. "Effect of *alpha*-Fluorination of Valproic Acid on Valproyl-S-Acyl-CoA Formation in Vivo in Rats" *Drug Metabolism and Disposition*, Vol. 29(9) pp. 1210-1215, 2001

This product is to be used for research only. It is not intended for drug or diagnostic use, human consumption or to be used in food or food additives. Matreya assumes no liability for any use of this product by the end user. We believe the information, offered in good faith, is accurate.