

PRODUCT DATA SHEET

2-Fluoropalmitic acid

Catalog No: 1717

Activity: Acyl-CoA synthase inhibitor

Source: synthetic

Solubility: chloroform, methanol, ethanol

CAS No: 16518-94-8

Molecular Formula: C₁₆H₃₁FO₂

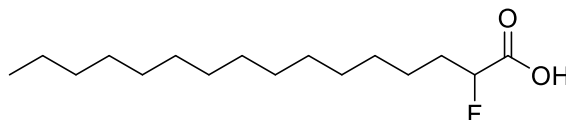
Molecular Weight: 274

Storage: -20°C

Purity: TLC > 98%, identity confirmed by MS

TLC System: hexane/ethyl ether/acetic acid
(70:30:2 by vol.)

Appearance: solid



Application Notes:

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.