



15-Methylhexadecanoic acid

Catalog number: 1606 Synonyms: iso-Heptadecanoic acid; *iso* C17 Fatty acid; 15-Methylpalmitic acid Source: synthetic Solubility: chloroform, ethyl ether, ethanol CAS number: 1603-03-8 Molecular Formula: C₁₇H₃₄O₂ Molecular Weight: 270 Storage: -20°C Purity: TLC > 98%; GC > 98% TLC System: hexane/ethyl ether (80:20) Appearance: solid



Application Notes:

15-Methylhexadecanoic acid is present in many plants, animals, and other organisms and has a central role in the biosynthesis of other *iso*-fatty acids in *Stigmatella aurantiaca*.¹ In bacteria, the *iso*-fatty acid² content and composition can often be used as a taxonomic marker³ because *iso*-fatty acids are often found in bacteria but not commonly in other microorganisms. Some bacteria have *iso*- but not *anteiso*-fatty acids while others have *anteiso*- but not *iso*-fatty acids. *Iso*-fatty acids are found in small amounts in marine organisms and ruminants, mainly due to the food chain but also due to some *de novo* synthesis. *Iso*-fatty acids with a total even number of carbons are more common than a total odd number. Some waxy materials such as lanolin, as well as secretions near animal eyes, contain an unusually high amount of *iso*-fatty acids being employed for their lubricating effect. Branched chain fatty acids are critical for the regulation of fluidity in membranes and in membrane transport for many types of bacteria due to their having a significantly lower transition temperature than straight chain fatty acids. Some bacteria have been found to be activators for various enzymes and systems and are used as protein modifiers. Although *iso*-even chain fatty acids via α-oxidation of *iso*-fatty acids.⁴ While *iso*-fatty acids are not usually found in plant oils the waxy surface of leaves can contain significant amounts of these fatty acids.

Selected References:

- 1.J. Dickschat et al. "Biosynthesis of iso-fatty acids in myxobacteria" Organic and Biomolecular Chemistry, Vol. 3 pp. 2824-2831, 2005
- T. Kaneda "Iso- and Anteiso-Fatty Acids in Bacteria: Biosynthesis, Function, and Taxonomic Significancet" American Society for Microbiology, Vol. 55(2) pp. 288-302, 1991
- 3. N. Jensen and M. Gross "Fast Atom Bombardment and Tandem Mass Spectrometry for Determining Iso- and Anteiso-Fatty Acids" Lipids, Vol. 21(5) pp. 362-365, 1986
- 4. H. Bode et al. "Biosynthesis of *Iso*-Fatty Acids in Myxobacteria: *Iso*-Even Fatty Acids Are Derived by α-Oxidation from *Iso*-Odd Fatty Acids" *J. Am. Chem. Soc.*, Vol. 127(2) pp 532–533, 2005

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.