PRODUCT DATA SHEET



Methyl 3-hydroxyheptadecanoate

Catalog number: 1742

Common names: 3-Hydroxy C17:0 methyl ester

Source: synthetic

Solubility: ethanol, methanol **CAS number:** 112538-92-8 **Molecular Formula:** C₁₈H₃₆O₃

Molecular Weight: 300

Storage: -20°C

Purity: TLC: >98%, GC: >98%; identity

confirmed by MS

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

Appearance: solid

Application Notes:

This product is a high purity 3-hydroxy fatty acid methyl ester that is ideal as a standard and for biological systems. This odd numbered fatty acid is unusual in many biological systems and can therefore often be used as a biomarker. 3-Hydroxyheptadecanoic acid has been used to characterize certain bacteria such as some gliding bacteria. 3-Hydroxy fatty acids have been found to be converted to the *omega*-fatty acid by the enzyme CYP4F11 and then into dicarboxylic acids *in vivo*. 3-hydroxy fatty acids are used as biomarkers for fatty acid oxidative disorders of both the long- and short-chain 3-hydroxy-acyl-CoA dehydrogenases. 3.4

Selected References:

- 1. P. Srisukchayakul et al. "Rapidithrix thailandica gen. nov., sp. nov., a marine gliding bacterium isolated from samples collected from the Andaman sea, along the southern coastline of Thailand" IJSEM, vol. 57 pp. 2275-2279, 2007
- 2. M. Dhar et al. "Omega oxidation of 3-hydroxy fatty acids by the human CYP4F gene subfamily enzyme CYP4F11" *Journal of Lipid Research*, vol. 49, pp. 612-624, 2008
- 3. P. Jones et al. "Improved Stable Isotope Dilution-Gas Chromatography-Mass Spectrometry Method for Serum or Plasma Free 3-Hydroxy-Fatty Acids and Its Utility for the Study of Disorders of Mitochondrial Fatty Acid beta-Oxidation" Clinical Chemistry, vol. 46, pp. 149-155, 2000
- 4. P. Jones et al. "Accumulation of free 3-hydroxy fatty acids in the culture media of fibroblasts from patients deficient in long-chain 1-3-hydroxyacyl-CoA dehydrogenase: a useful diagnostic aid" Clinical Chemistry, vol. 47(7) pp. 1190-1194, 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.