

PRODUCT DATA SHEET

Methyl 3-hydroxyhexadecanoate

Catalog number: 1740

Common names: 3-Hydroxy C16:0 methyl ester; Methyl 3-hydroxypalmitate

Source: synthetic

Solubility: ethanol, methanol

CAS number: 51883-36-4

Molecular Formula: C₁₇H₃₄O₃

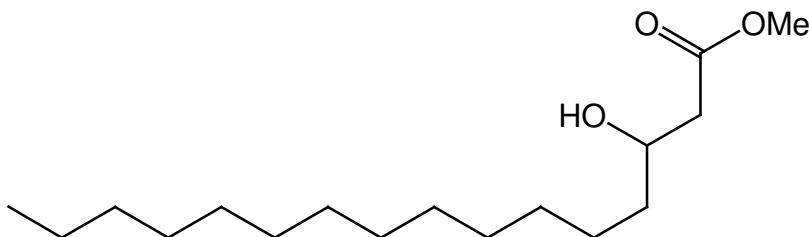
Molecular Weight: 286

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 standard that is useful for the investigation of disorders and diseases. Polyhydroxyalkenoates, polyesters produced by bacteria fermentation, are used for carbon and energy storage and are of interest in studies regarding their synthesis, properties and mechanisms. Medium chain-length polyhydroxyalkenoate monomers such as 3-hydroxyhexadecanoic acid may have pharmaceutical properties. The biologically natural chiral (R)-3-hydroxyhexadecanoic acid is an intermediate in fatty acid biosynthesis. It has been suggested that the methyl ester of 3-hydroxyhexadecanoic acid is an intercellular signal that autoregulates virulence gene expression¹ and that it modulates production of PhcA-regulated virulence factors in *Ralstonia (Pseudomonas) solanacearum*. 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*.² 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. A. Flavier et al. "Identification of 3-hydroxypalmitic acid methyl ester as a novel autoregulator controlling virulence in *Ralstonia solanacearum*" *Mol Microbiol*, vol. 26 pp. 251-259, 1997
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.