

PRODUCT DATA SHEET

Methyl-3-hydroxydodecanoate

Catalog number: 1732

Common Name: 3-Hydroxy C12:0 methyl ester

Source: synthetic

Solubility: chloroform, ethyl ether, ethanol

CAS number: 72864-23-4

Molecular Formula: C₁₃H₂₆O₃

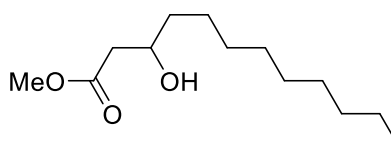
Molecular Weight: 230

Storage: -20°C

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

TLC System: hexane/ethyl ether, (85:15)

Appearance: liquid



Application Notes:

This 3-hydroxydodecanoic acid methyl ester 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-hydroxydodecanoic acid may have pharmaceutical properties. The biologically natural chiral (R)-3-hydroxydodecanoic acid is an intermediate in fatty acid biosynthesis. 3-Hydroxydodecanoic acid is the major fatty acid in some lipopolysaccharides.¹ 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.^{2,3}

Selected References:

1. K. Kunii et al. "A new extraction method for Acinetobacter species ODB-L2 rough form lipopolysaccharide from culture broth" *Microbios.*, vol. 105:412 pp. 153-61, 2001
2. 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 β -Oxidation" *Clinical Chemistry*, vol. 46, pp. 149-155, 2000
3. 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.

DS1732 Rev. #6
December 22nd, 2021