PRODUCT DATA SHEET



Methyl 2-hydroxydodecanoate

Catalog number: 1702

Common names: 2-Hydroxy C12:0 methyl

ester

Source: synthetic

Solubility: methanol, chloroform, ethyl ether

CAS number: 51067-85-7

Molecular Formula: C₁₃H₂₆O₃

Molecular Weight: 230

Storage: -20°C

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

confirmed by MS

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

Appearance: solid

Application Notes:

alpha-Hydroxy fatty acids are abundant in nervous tissues and are components of cerebrosides and sulfatides, which are mostly found in the myelin of nervous tissues. They are common in cosmetics, skin creams, and lotions. 2-hydroxy acids display complex monolayer phase behavior due to the additional hydrogen bonding afforded by the presence of the second hydroxy group and therefore play an important role in the membrane structure.\(^1\) 2-hydroxydodecanoic acid demonstrates antifungal properties, although to a lesser extent than 3-hydroxy fatty acids.\(^2\) 2-Hydroxydodecanoic acid (along with 2-hydroxytetradecanoic acid) induced a hyphal growth response of the arbuscular mycorrhizal (AM) fungus, \(Gigaspora\) gigantean while other 2-hydroxy, 3-hydroxy, and non-hydroxy fatty acids did not. This indicates that 2-hydroxy fatty acids are a putative category of root exudate signal perceived by \(Gigaspora\) species, stimulating an increase in elongated lateral branches.\(^3\) alpha-Oxidation of 2-hydroxy fatty acids to CO2 and saturated acids occurs in the peroxisome and is unique from the \(alpha\)-Oxidation of \(beta\)-carbon branched fatty acids such as phytanic acid. Cells from Zellweger syndrome and peroxisome-deficient cells are unable to undergo \(alpha\)-oxidation although patients with other peroxisomal disorders such as X-linked adrenoleukodystrophy, Refsum disease, and rhizomelic chondrodysplasia punctata are able to.\(^4\)

Selected References:

- 1. C. Lendrum et al. "Nonequilibrium 2-Hydroxyoctadecanoic Acid Monolayers: Effect of Electrolytes" Langmuir, vol. 27 pp. 4430-4438, 2011
- 2. J. Sjogren et al. "Antifungal 3-Hydroxy Fatty Acids from Lactobacillus plantarum MiLAB 14" Applied and Environmental Microbiology, vol. 69 pp. 7554-7557, 2003
- 3. G. Nagahashi and D. Douds Jr. "The effects of hydroxy fatty acids on the hyphal branching of germinated spores of AM fungi" Fungal Biology, vol. 115 pp. 351-358, 2011
- 4. R. Sandhir, M. Khan, and I. Singh "Identification of the Pathway of α-Oxidation of Cerebronic Acid in Peroxisomes" *Lipids*, Vol. 35(10) pp. 1127-1133, 2000

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.