

# PRODUCT DATA SHEET

## Methyl 3-hydroxynonanoate

**Catalog number:** 1726

**Common names:** 3-Hydroxy C9:0 methyl ester

**Source:** synthetic

**Solubility:** chloroform, ethanol, ethyl ether

**CAS number:** 83968-06-3

**Molecular Formula:** C<sub>10</sub>H<sub>20</sub>O<sub>3</sub>

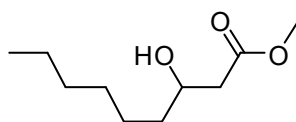
**Molecular Weight:** 188

**Storage:** -20°C

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

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

**Appearance:** liquid



### Application Notes:

This 3-hydroxynonanoic acid methyl ester is a high purity standard that is ideal for analysis and biological systems. 3-Hydroxy fatty acids are intermediates in fatty acid biosynthesis and have been found to be converted to the *omega*-fatty acid by the enzyme CYP4F11 and then into dicarboxylic acids *in vivo*.<sup>1</sup> 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.<sup>2</sup> 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 and are used as biodegradable plastics.<sup>3</sup> Medium chain-length polyhydroxyalkenoate monomers may have pharmaceutical properties.

### Selected References:

1. 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
2. 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
3. J. Gangotri et al. "Production of Chiral (*R*)-3-Hydroxyoctanoic Acid Monomers, Catalyzed by *Pseudomonas fluorescens* GK13 Poly(3-Hydroxyoctanoic Acid) Depolymerase" *Applied and Environmental Microbiology*, vol. 76 pp. 3554-3560, 2010

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.