

PRODUCT DATA SHEET

17-Hydroxyheptadecanoic acid

Catalog number: 1760

Common names: *omega*-Hydroxy C17:0 fatty acid

Source: synthetic

Solubility: chloroform, warm ethanol, ethyl ether

CAS number: 13099-34-8

Molecular Formula: C₁₇H₃₄O₃

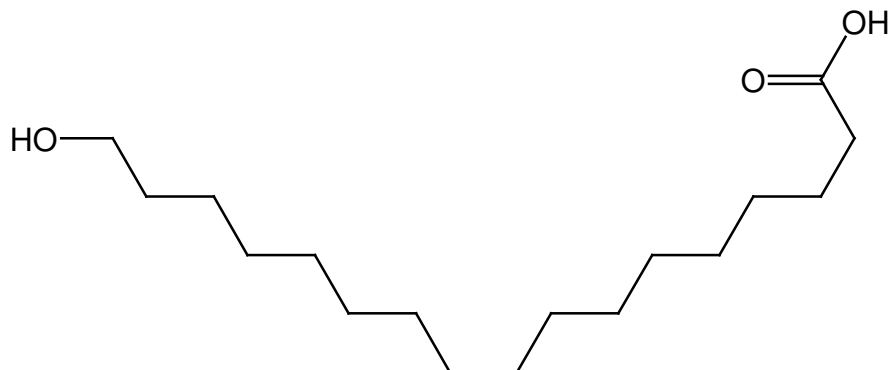
Molecular Weight: 286

Storage: room temperature

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

TLC System: chloroform/methanol/acetic acid (90:10:2)

Appearance: solid



Application Notes:

This odd numbered *omega*-hydroxy heptadecanoic acid is ideal as an internal standard for studies involving medium to long-chain *omega*-hydroxy fatty acids.¹ *omega*-Hydroxy C17:0 fatty acid and other *omega*-hydroxy fatty acids can be lactonized by certain enzymes into mono- and oligolactones.² *omega*-Oxidation is a minor fatty acid pathway used for fatty acid metabolism and usually occurs in the smooth endoplasmic reticulum. Stimulation of *omega*-hydroxylation has been proposed as a method for treating X-linked adrenoleukodystrophy, a disease that is characterized by elevated levels of very long chain fatty acids.³ *omega*-Hydroxy fatty acids have an important role while acylated to various lipids. *omega*-Hydroxylated very long chain fatty acid (VLCFA) ceramides are vital to skin barrier functions and a deficiency of these lipids can cause death from water loss through the skin. In atopic dermatitis, a common skin disease, there are considerable deficiencies of *omega*-hydroxy long chain fatty acids acylated to ceramides and this may contribute to the severely damaged permeability barrier found in this disease.⁴

Selected References:

1. M. Nakano, E. Kelly, and A. Rettie "Expression and Characterization of CYP4V2 as a Fatty Acid ω -Hydroxylase" *Drug Metabolism and Disposition*, vol. 37 pp. 2119-2122, 2009
2. U. Antczak et al. "Enzymatic lactonization of 15-hydroxypentadecanoic and 16-hydroxyhexadecanoic acids to macrocyclic lactones" *Enzyme and Microbial Technology*, vol. 13 pp. 589-593, 1991
3. R. Sanders et al. "Omega-Oxidation of Very Long-Chain Fatty Acids in Human Liver Microsomes: Implications for X-Linked Adrenoleukodystrophy" *Journal of Biological Chemistry*, Vol. 281 pp. 13180-13187, 2006
4. O. Macheleidt, H. Kaiser, K. Sandhoff "Deficiency of epidermal protein-bound omega-hydroxyceramides in atopic dermatitis" *J Invest Dermatol.*, vol. 119 pp. 166-173, 2002

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.