

PRODUCT DATA SHEET

12-Methyltetradecanoic acid

Catalog number: 1615

Common Name: anteiso-Pentadecanoic acid;
anteiso C15 Fatty acid

Source: synthetic

Solubility: chloroform, ethyl ether, ethanol

CAS number: 5502-94-3

Molecular Formula: C₁₅H₃₀O₂

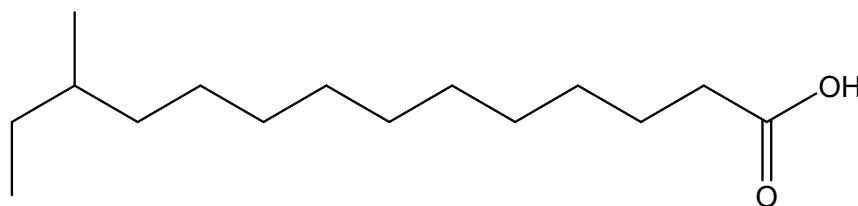
Molecular Weight: 242

Storage: -20°C

Purity: TLC > 98%, GC > 98%

TLC System: hexane/ethyl ether/acetic acid,
(80:20:1 by vol.)

Appearance: solid



Application Notes:

This high purity anteiso-pentadecanoic acid is ideal for use as a standard in the verification of samples¹ and for use in biological studies. anteiso-Pentadecanoic acid is one of the most abundant fatty acids in members of the genus *Bacillus*.² Many types of bacteria and marine organisms produce anteiso-fatty acids which can be used for their characterization.³ Some bacteria have iso- but not anteiso-fatty acids while others have anteiso- but not iso-fatty acids. anteiso-Fatty acids are also found in ruminant animals (from the rumen microorganisms) and in ruminant milk mainly from the food chain but also due to some *de novo* synthesis. Branched chain fatty acids have been found in the gut of newborn animals where they have a role in the microorganism gut colonization.⁴ Some bacteria handle stress (such as heat and toxicity) by changing the ratio of anteiso/iso-fatty acids in the cell membrane. Although anteiso-fatty acids are not usually found in plant oils the waxy surface of leaves can contain significant amounts of these fatty acids. Branched chain fatty acids are critical for the regulation of fluidity in membranes and in membrane transport for many types of bacteria due to their having a significantly lower transition temperature than straight chain fatty acids.

Selected References:

1. N. Jensen and M. Gross "Fast Atom Bombardment and Tandem Mass Spectrometry for Determining Iso- and Anteiso- Fatty Acids" *Lipids*, Vol. 21(5) pp. 362-365, 1986
2. H. Daron "Nutritional Alteration of the Fatty Acid Composition of a Thermophilic Bacillus Species" *Journal of Bacteriology*, vol. 116 pp. 1096-1099, 1973
3. E. Kim et al. "Fatty Acid Profiles Associated with Microbial Colonization of Freshly Ingested Grass and Rumen Biohydrogenation" *Journal of Dairy Science*, Vol. 88 pp. 3220-3230, 2005
4. R. Ran-Ressler et al. "Branched Chain Fatty Acids Are Constituents of the Normal Healthy Newborn Gastrointestinal Tract" *Pediatric Research*, Vol. 64(6) pp. 605-609, 2008

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.