

## PRODUCT DATA SHEET

## 14-Methylhexadecanoic acid

**CAS number:** 5918-29-6

Catalog number: 1616 Molecular Formula: C<sub>17</sub>H<sub>34</sub>O<sub>2</sub>

Common Name: anteiso-Heptadecanoic acid; Molecular Weight: 270

anteiso C17 Fatty acid **Storage:** -20°C

**Source:** synthetic **Purity:** TLC >98%, GC >98%

Solubility: chloroform, ethyl ether, ethanol TLC System: hexane/ethyl ether/acetic acid,

(80:20:1 by vol.)

Appearance: solid

## **Application Notes:**

This high purity anteiso-heptadecanoic acid is ideal for use as a standard <sup>1</sup> and in biological studies. Many types of bacteria and marine organisms produce anteiso-fatty acids which can be used for their characterization. <sup>2</sup> 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. <sup>3</sup> 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. This anteiso-14-methylhexadecanoic acid has been found in pine seed oil, a major exception to branched chain fatty acids which are found almost exclusively in animals and microorganisms. <sup>4</sup>

## **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. 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
- 3. 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
- 4. R. Wolff, W. Christie and D. Coakley "The unusual occurrence of 14-methylhexadecanoic acid in pinaceae seed oils among plants" *Lipids*, vol. 32 pp, 971-973, 1997

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.