

PRODUCT DATA SHEET

Methyl 14-methylhexadecanoate

Catalog number: 1614 Synonyms: anteiso-Heptadecanoic methyl ester; anteiso C17:0 methyl ester Source: synthetic Solubility: chloroform, ethyl ether, ethanol CAS number: 2490-49-5 Molecular Formula: C₁₈H₃₆O₂ Molecular Weight: 284 Storage: -20°C Purity: TLC > 98%, GC > 98% TLC System: hexane/ethyl ether (85:15) Appearance: liquid



Application Notes:

This anteiso-heptadecanoic methyl ester is ideal for use as a standard in the verification of samples.¹ Many types of bacteria and marine organisms produce anteiso fatty acids which can be used to characterize the various organisms.² anteiso-Fatty acids, especially anteiso-heptadecanoic acid, are also found in ruminant animals (from the rumen microorganisms) and in ruminant milk. Branched chain fatty acids have been found in the gut of newborn animals where they have a role in the microorganism gut colonization.³ anteiso-Heptadecanoic acid is found in a very few plant species such as in the pine seed oil of *Pinaceae*.

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*

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

2. 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.