

PRODUCT DATA SHEET

2-Hydroxy Methyl Ester Mixture (quantitative)

Catalog No: 1722 Solvent: chloroform Storage: -20°C Concentration: 10mg/ml Volume: 1ml

GC Conditions:

T1 4.

Column: SPD-1 or RTX-1 Carrier Gas: helium Make-up Gas: nitrogen Split Ratio: 100:1 Oven Initial: 280°C Oven Final: 280°C Detector: FID. 320 °C

Linear Velocity: 20cm/sec Flow Rate: 40ml/min Vent Flow: 70ml/min Program Rate: isothermal Hold Time: Injector: 300°C

Elution	Carbon		
Order	Number	Component Name	% Conc. by weight
1	C14:0	Methyl 2-hydroxytetracosanoate	20.0
2	C16:0	Methyl 2-hydroxyhexadecanoate	20.0
3	C18:0	Methyl 2-hydroxyoctadecanoate	15.0
4	C20:0	Methyl 2-hydroxyeicosanoate	15.0
5	C22:0	Methyl 2-hydroxydocosanoate	10.0
6	C23:0	Methyl 2-hydroxytricosanoate	10.0
7	C24:0	Methyl 2-hydroxytetracosanoate	10.0

Composition in weight percent determined by synthesis, not by analysis.

Application Notes:

This fatty acid mixture contains common *alpha*-hydroxy fatty acids in methylene chloride for the quantitative identification and quantitation of unknowns. Microbial fatty acid profiles are unique from one species to another and can therefore be used in the determination of bacterial identity. All materials are analyzed to verify their identity and to determine their purity. All analytes are $98^+\%$ pure. This standard is accurately prepared by gravimetric technique (+/- 0.5%) and all glassware is class A rated. Ampules are purged with nitrogen/argon before and after filling and chilled before being flame sealed. Store ampules sealed as received and process without delay immediately after opening the ampule.

Selected References:

1. M. Or-Rashid, N. Odongo and B. McBride, "Fatty acid composition of ruminal bacteria and protozoa, with emphasis on conjugated linoleic acid, vaccenic acid, and odd-chain and branched-chain fatty acids" *Journal of Animal Science*, Vol. 85 pp. 1228, 2007

2. Y.Zhang, S. White, and C. Rock "Inhibiting Bacterial Fatty Acid Synthesis" *The Journal of Biological Chemistry*, Vol. 281(26) pp. 17541, 2006 3. N. Rozès, S. Garbay, M. Denayrolles, A. Lonvaud-Funel "A rapid method for the determination of bacterial fatty acid composition" *Applied*

Microbiology, Vol. 3(17) pp. 126 1993

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.