

PRODUCT DATA SHEET

<u>KEL-FIM-FAME-5 Mixture</u> (quantitative)		Carbon	Component	
		Number	Name n	ng/ml
		C8:0	Methyl octanoate (caprylate)	0.3
Catalog No:	4210	C10:0	Methyl decanoate (caprate)	0.5
_	-20°C	C12:0	Methyl dodecanoate (laurate)	1.0
Storage:		C13:0	Methyl tridecanoate	0.5
Concentration:	Various	C14:0	Methyl tetradecanoate (myristate)	0.5
Quantity:	1ml	C14:1	Methyl tetradecenoate (cis-9), (myristoleate)	0.3
-		C15:0	Methyl pentadecanoate	0.3
		C16:0	Methyl hexadecanoate (palmitate)	2.0
		C16:1	Methyl hexadecenoate (cis-9), (palmitoleate)	1.0
		C17:0	Methyl heptadecanoate (margarate)	0.5
		C18:0	Methyl octadecanoate (stearate)	1.0
		C18:1	Methyl octadecenoate (trans-9), (elaidate)	0.4
		C18:1	Methyl octadecenoate (cis-9), (oleate)	3.0
		C18:2	Methyl octadecadienoate (all cis-9,12), (linoleate)	2.0
		C20:0	Methyl eicosanoate (arachidate)	0.3
		C18:3	Methyl octadecatrienoate (all cis-9,12,15), (linolenate)	
		C20:1	Methyl eicosenoate (cis-11), (gondoate)	0.3
		C22:0	Methyl docosanoate (behenate)	0.3
		C22:1	Methyl docosenoate (cis-13), (erucate)	0.3
			Column: Rtx-2330 30m x 0.25mm x 0.2u Oven: 170°C (3 min.), 4°C/min. to 230°C (Inj.: 250°C Det.: 250°C Carrier gas: helium, 20cm/sec. Injection: 1.0uL, 100:1 split	(4min.)

Application Notes:

This fatty acid methyl ester mixture contains 19 fatty acids for the identification and quantification of unknowns. ¹ It is prepared from high purity stock material and contains saturated and unsaturated fatty acids. This mixture is very useful for bacterial identification, ² triglyceride determination and the analysis of various plant and animal lipids. ³ Understanding the role of fatty acids and fatty acid metabolism in plants and animals is important in drug development.

Selected References:

- 1. T. Murata "Analysis of fatty acid methyl esters by a gas-liquid chromatography-chemical ionization mass spectrometry computer system" *Journal of Lipid Research*, Vol. 19:166, 1978
- 2. N. Rozès et al. "A rapid method for the determination of bacterial fatty acid composition" Applied Microbiology, Vol. 3:17 pp.126, 1993
- 3. D. Welch "Applications of cellular fatty acid analysis" Clinical Microbiology Reviews, Vol. 4:4 pp. 422, 1991

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.