

PRODUCT DATA SHEET

FIM-FAME-8 Mixture (quantitative)

Catalog No: 2012

Solvent: Methylene chloride

Storage: -20°C

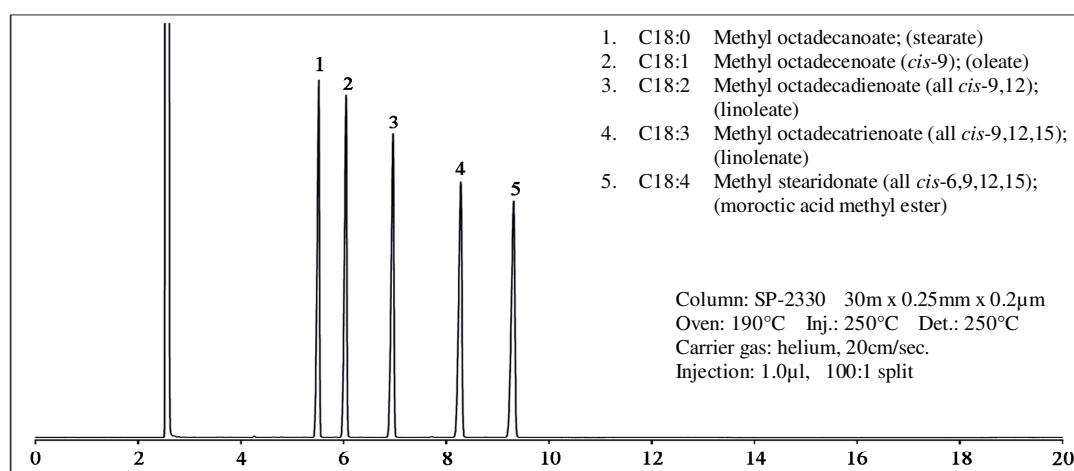
Column: SP-2330, 30m x 0.25mm x 0.2µm

Oven: 190°C

Carrier: Helium @ 20cm/sec.

Detector: FID, 250 °C

Injector: 250°C



Application notes:

This FAME mixture contains 5 fatty acid methyl esters for the identification and quantification of unknowns.¹ It is prepared from high purity stock material and contains saturated and unsaturated fatty acids in the 18 carbon series. 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.