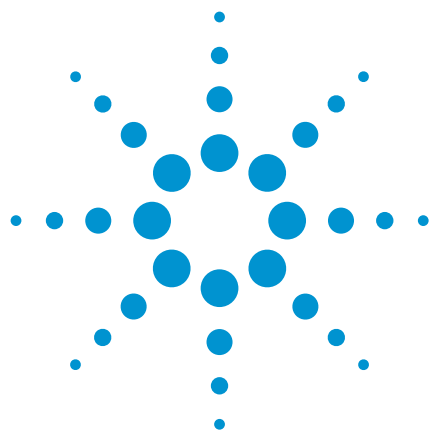




FAMES and FABEs

Comparison of fatty acid methyl esters and fatty acid iso-butyl esters

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 88 for FAME column separates four components in a comparison of fatty acid methyl esters and fatty acid –isobutyl esters in 27 minutes.



Agilent Technologies

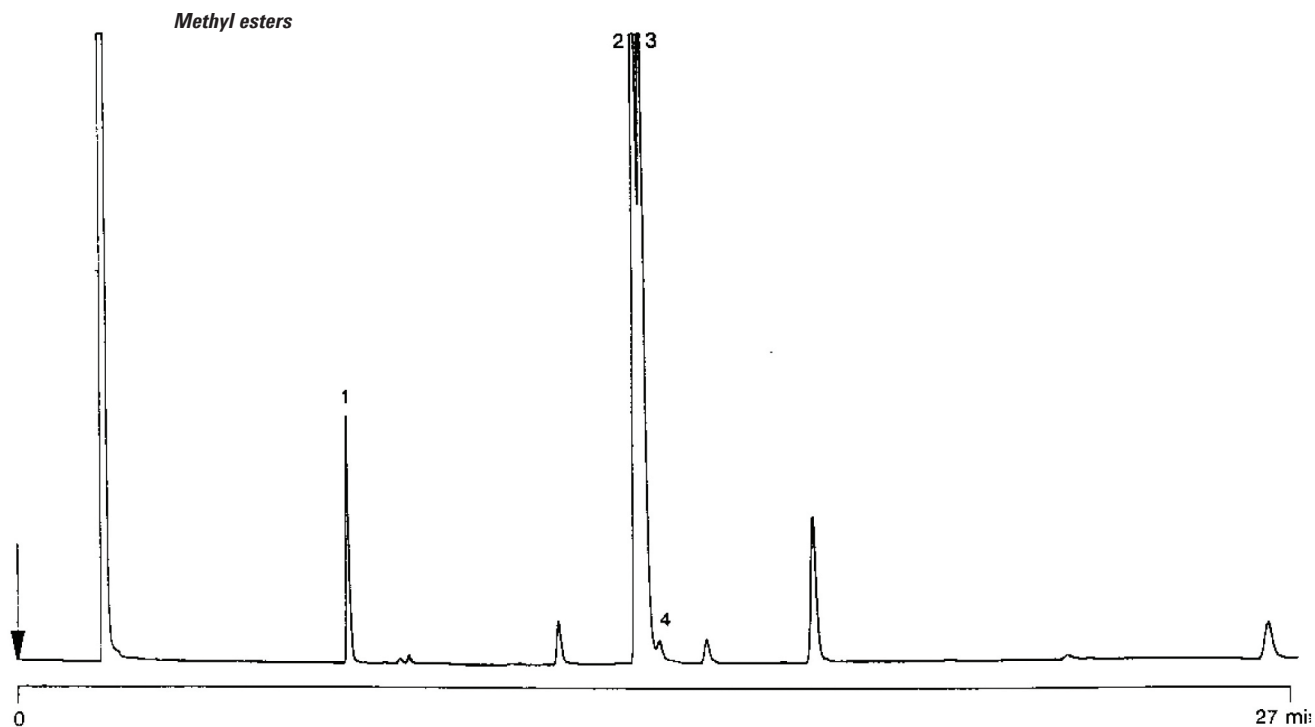
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 88 for FAME, 0.25 mm x 50 m fused silica WCOT CP-Sil 88 (df = 0.2 µm)
(Part no. CP7488)
Temperature : 170 °C
Carrier Gas : H₂
Injector : Splitter, 2 mL/min
T = 250 °C
Detector : FID
T = 250 °C
Sample Size : 1 µL
Solvent Sample : heptane

Courtesy : C. Mariani, Stazione sperimentale per le industrie degli oli e dei grassi, Milano, Italy

Peak identification

1. C16:0 (palmitic acid)
2. C16:1ω6, cis (petroselinic acid)
3. C16:1ω9, cis (oleic acid)
4. C18:1ω11, trans (vaccenic acid)

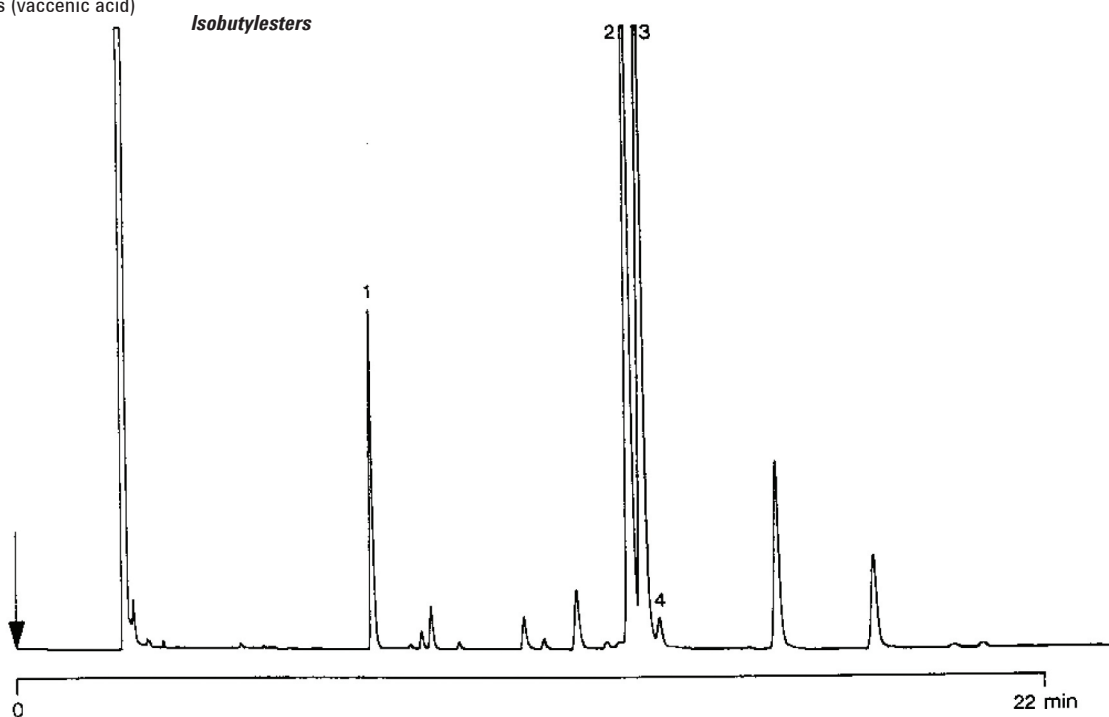


Conditions

Technique	: GC-capillary
Column	: Agilent CP-Sil 88 for FAME, 0.25 mm x 50 m fused silica WCOT CP-Sil 88 (df = 0.2 µm) (Part no. CP7488)
Temperature	: 170 °C
Carrier Gas	: H ₂
Injector	: Splitter, 2 mL/min T = 250 °C
Detector	: FID T = 250 °C
Sample Size	: 1 µL
Solvent Sample	: heptane
Courtesy	: C. Mariani, Stazione sperimentale per le industrie degli oli e dei grassi, Milano, Italy

Peak identification

1. C16:0 (palmitic acid)
2. C16:1 ω 6, cis (petroselinic acid)
3. C16:1 ω 9, cis (oleic acid)
4. C18:1 ω 11, trans (vaccenic acid)



www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A00769



Agilent Technologies