



Free fatty acids

Analysis of free fatty acids, C₂-C₄

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-FFAP CB column separates free fatty acids in five minutes.



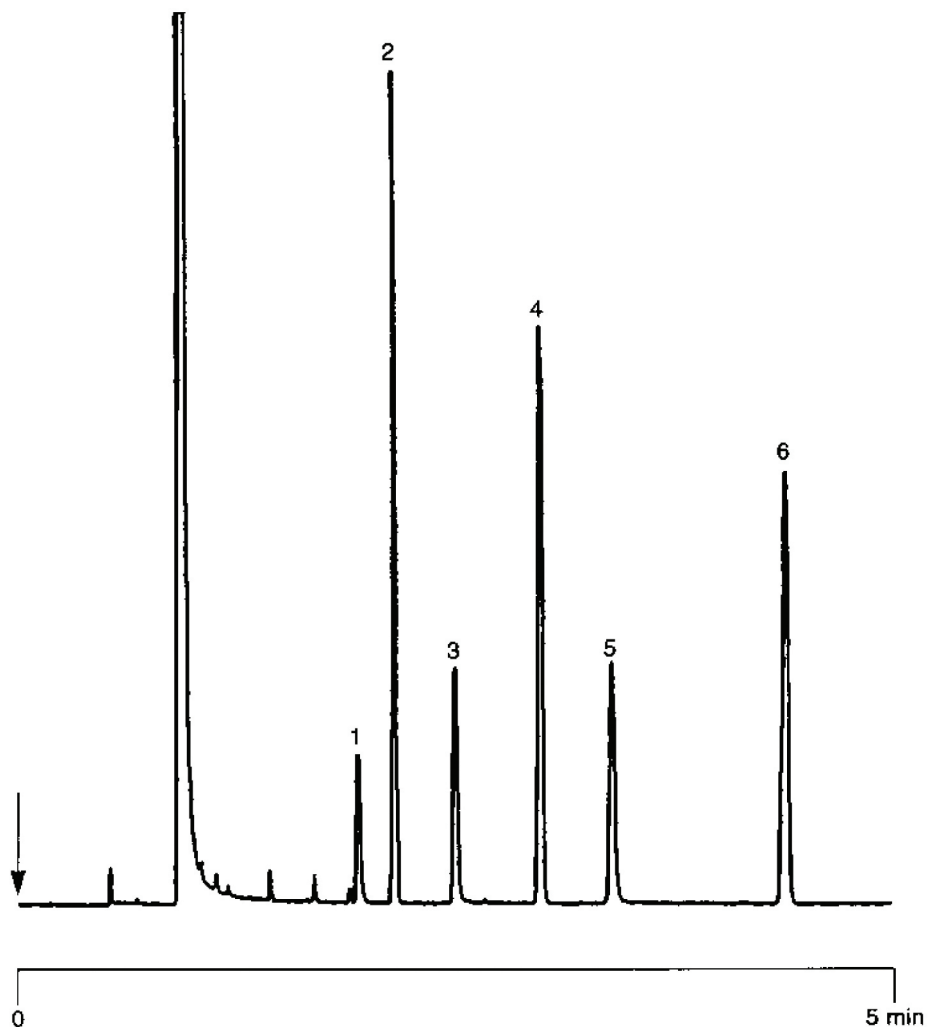
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-FFAP CB, 0.15 mm x 25 m fused silica
WCOT CP-FFAP CB (df = 0.25 µm)
(Part no. CP7686)
Temperature : 130 °C
Carrier Gas : H₂, 200 kPa (2.0 bar, 24 psi)
Injector : Splitter
Detector : FID
Sample Size : 0.3 µL
Sample : 0.1% in cyclohexane

Peak identification

1. acetic acid
2. n-pentadecane
3. propionic acid
4. n-hexadecane
5. butyric acid
6. n-heptadecane



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

A01093



Agilent Technologies