



Free fatty acids

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

The trifluoropropyl phase of Agilent VF-200ms is very stable, which allows the analysis of aggressive compounds. Loadability of the thin film columns, as shown here, is quite limited especially for C_3 acids and a thicker film would be preferred. For acetic acid solubility is relatively high which results in a narrow peak.



Agilent Technologies

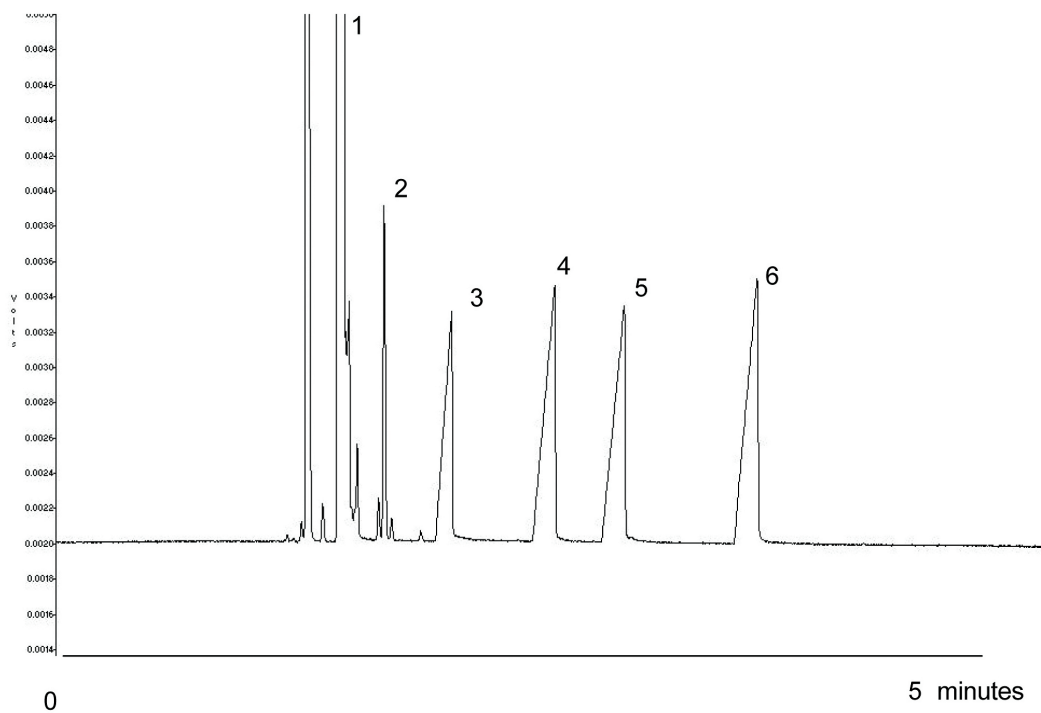
Conditions

Technique : GC
Column : Agilent FactorFour VF-200ms, 0.25 mm x 30 m
(df = 0.25 μ m) (Part no. CP8858)
Temperature : 45 °C, 10 °C/min \rightarrow 325 °C
Carrier Gas : Helium, ca. 1.0 mL/min
Pressure program : 60 kPa
Injector : Split/Splitless, in split mode, 1:100
Detector : FID
Sample Size : 1 μ L
Solvent : cyclohexane, 0.1% w/v

Courtesy : Jane Peene, Agilent Application Laboratory,
Middelburg, The Netherlands

Peak identification

1. cyclohexane
2. acetic acid
3. propionic acid
4. iso-butyric acid
5. butyric acid
6. iso-valeric 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

A02203



Agilent Technologies