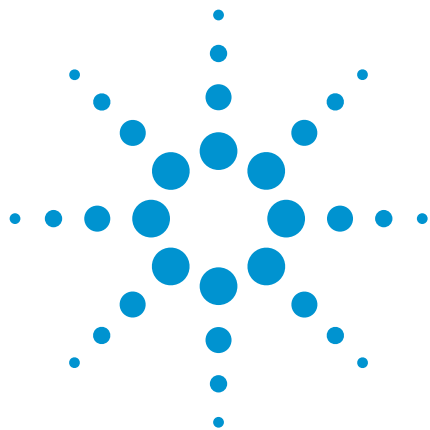




Waxes

Analysis of paraffin wax

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-SimDist Ultimetall column separates five components in a paraffin wax in 50 minutes.

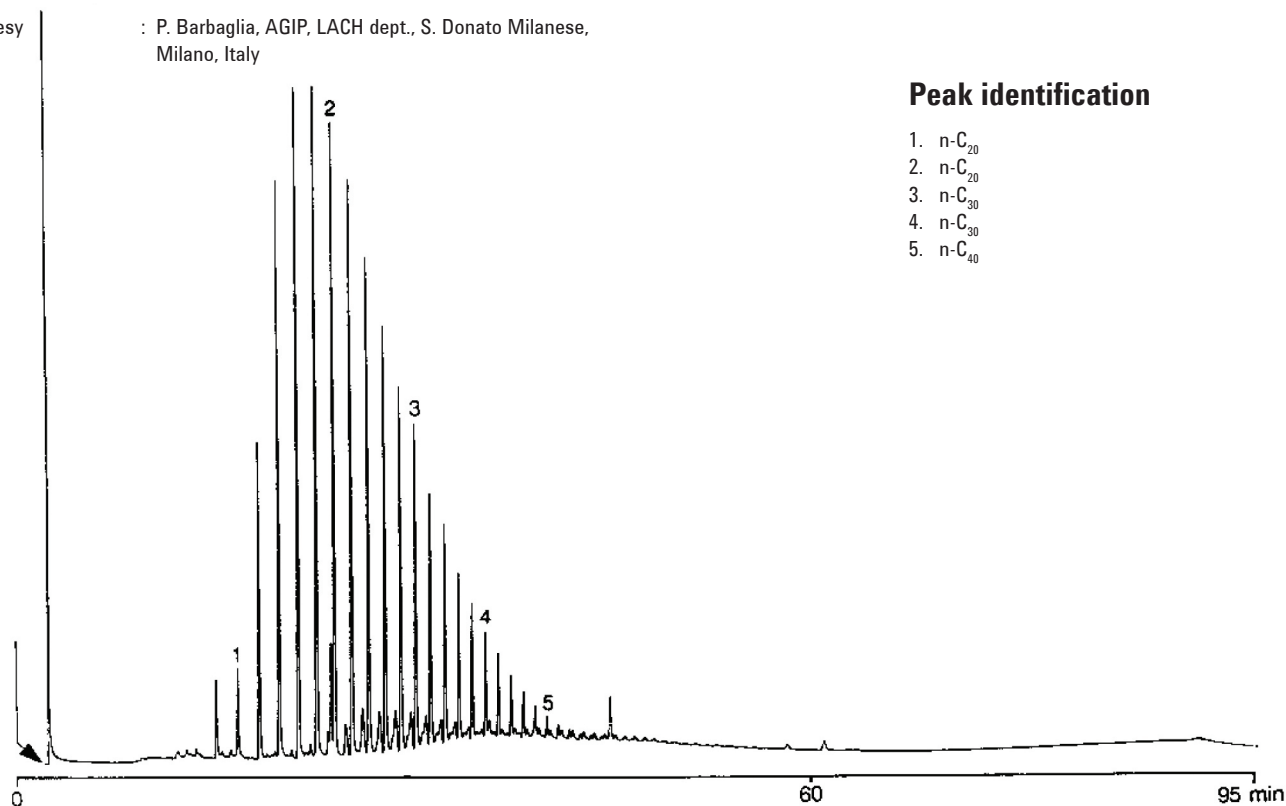


Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-SimDist CB UltiMetal, 0.5 mm x 10 m
WCOT CP-SimDist CB UltiMetal (df = 0.15 µm)
(Part no. CP7542)
with WCOT metal retention gap
0.55 mm x 75 cm methyl deactivated
(Part no. CP8067)
Temperature : 50 °C (2 min) → 100 °C, 10 °C/min;
100 °C → 300 °C, 5 °C/min;
300 °C → 430 °C, 3 °C/min; 430 °C (5 min)
Carrier Gas : He, 21 kPa (0.21 bar, 3 psi), 9 mL/min
Injector : cool on-column
Detector : FID; T = 430 °C
Sample Size : 0.5 µL
Solvent Sample : carbon disulfide; 10 mg/mL

Courtesy : P. Barbaglia, AGIP, LACH dept., S. Donato Milanese,
Milano, Italy



Peak identification

1. n-C₂₀
2. n-C₂₀
3. n-C₃₀
4. n-C₃₀
5. n-C₄₀

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

A00773



Agilent Technologies