



Flavors and aromas

Fingerprint analysis of deodorant

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column produces a deodorant fingerprint in 40 minutes.



Agilent Technologies

Conditions

Technique : GC-PTI

Column : Agilent CP-Wax 52 CB, 0.25 mm x 50 m fused silica
WCOT CP-Wax 52 CB (df = 0.2 µm)
(Part no. CP7723)

Temperature : 40 °C (5 min) → 210 °C, 5 °C/min

Injector : PTI

Cold trap : CP-Sil 8 CB
0.53 mm; df = 5 µm

Precool temp. : -130 °C

Precool time : 3 min

Purge time : 5 min

Purge flow : 12 mL/min

Injection temp. : 200 °C

Injection time : 1 min

Backflush flow : 50 mL/min

Detector : FID
T = 250 °C

Sample Size : 2 gr

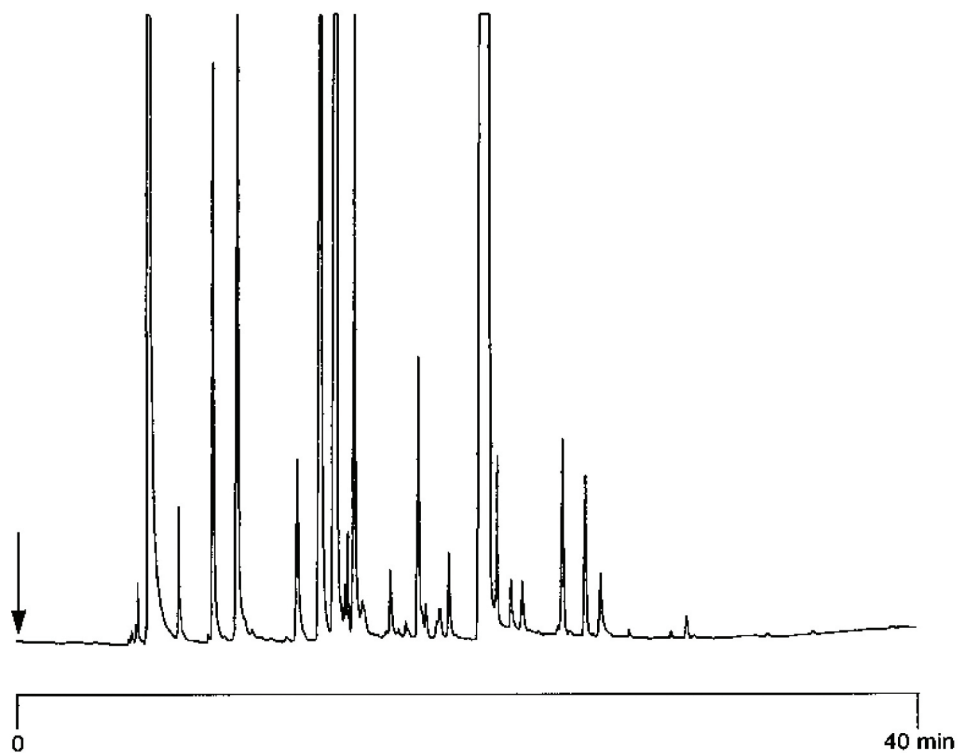
Sample Temperature : ambient

Concentration Range : ppb/ppm

Courtesy : GL Sciences, Japan

Peak identification

Peaks not identified



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

A01036



Agilent Technologies