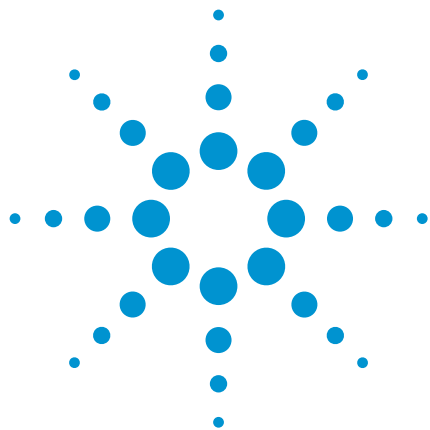




Hydrocarbons, C₂₄

Separation of squalane and squalene

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 5 CB column separates squalane and squalene in ten minutes.



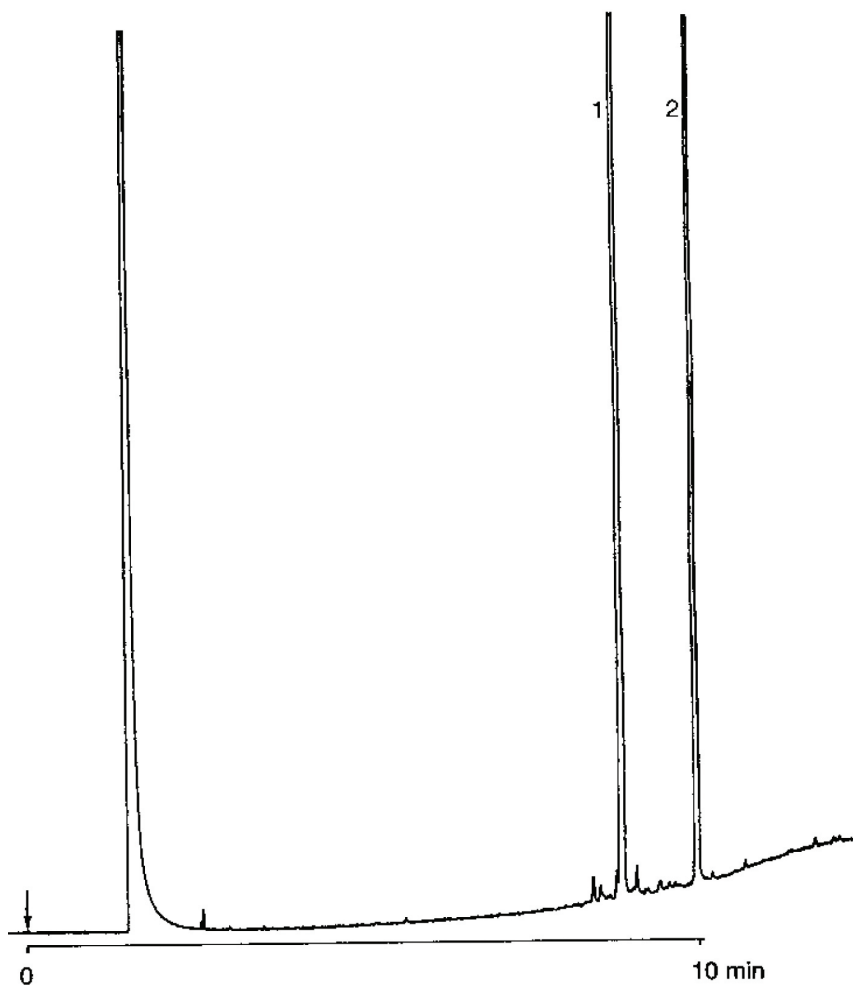
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 5 CB, 0.22 mm x 25 m fused silica
WCOT CP-Sil 5 CB (0.12 μ m) (Part no. CP7710)
Temperature : 200 °C (1 min) $\rightarrow$ 300 °C, 10 °C/min
Carrier Gas : He, 110 kPa (1.1 bar, 16 psi), 28 cm/s
Injector : Splitter, 82 mL/min
T = 300 °C
Detector : FID, 8×10^{-12} Afs
T = 300 °C
Sample Size : 0.3 μ L
Concentration Range : 0.5 % w/v/component
Solvent Sample : dichloromethane

Peak identification

1. squalane (0.5%)
2. squalene (0.5%)



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

A00156



Agilent Technologies