



Solvents

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent 490 Micro GC and HayeSep A column module separates four solvents in a gas mix in 220 seconds.



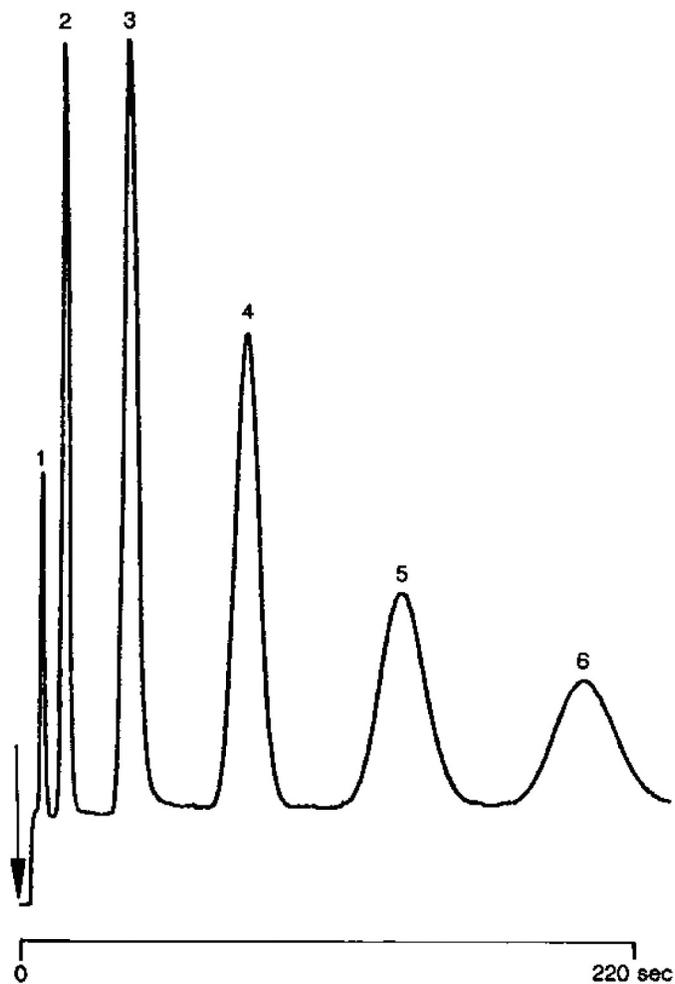
Agilent Technologies

Conditions

Technique : Micro GC
Column : Agilent CP-HayeSep A
Temperature : 120 °C
Carrier Gas : He
Detector : TCD

Peak identification

1. air
2. ethylene
3. water
4. chloroethane
5. ethanol
6. diethyl ether



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

A00888



Agilent Technologies