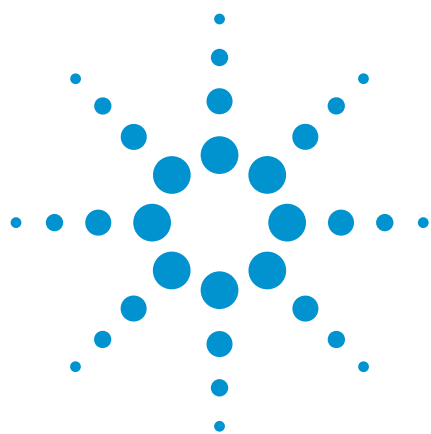




# Ketones, C<sub>3</sub> – C<sub>10</sub>

## Application Note

Energy & Fuels

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates 15 C<sub>3</sub> to C<sub>10</sub> ketones in ten minutes.



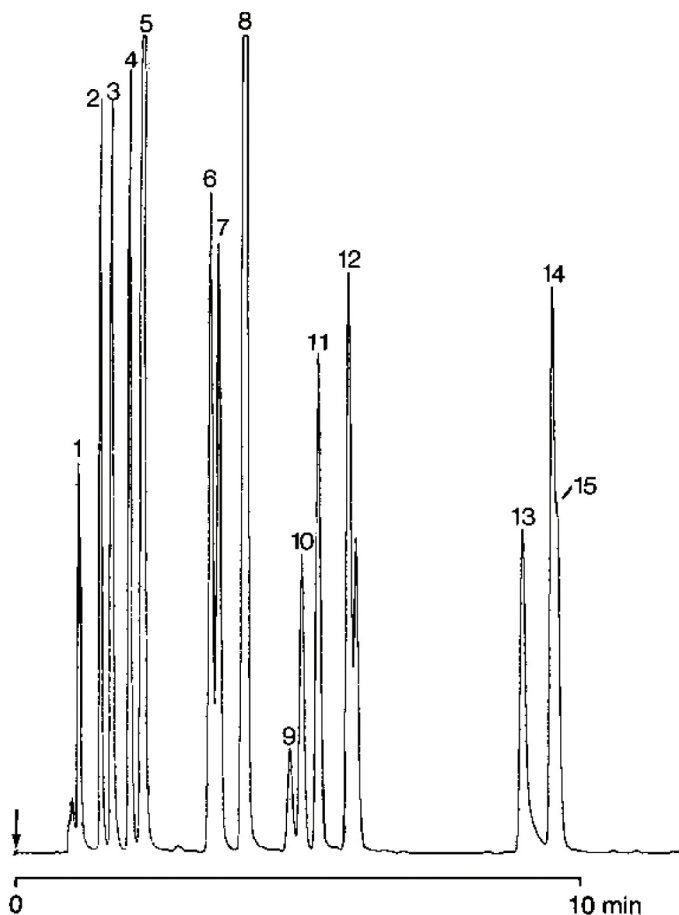
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Wax 52 CB, 0.53 mm x 25 m fused silica  
WCOT CP-Wax 52 CB (2.0  $\mu$ m) (Part no. CP7658)  
Temperature : 50 °C  $\rightarrow$  200 °C, 10 °C/min  
Carrier Gas : N<sub>2</sub>, 47 cm/s (10 mL/min)  
Injector : Direct  
T = 250 °C  
Detector : FID, 100 x 10<sup>-12</sup> Afs  
T = 275 °C  
Sample Size : 0.2  $\mu$ L  
Concentration range : 1 %

## Peak identification

1. acetone
2. butanone
3. methylisopropylketone
4. diethylketone
5. diisopropylketone
6. dipropylketone
7. methylisoamylketone
- B. diisobutylketone + 2-heptanone + cyclopentanone
9. 3-octanone (ethylamylketone)
10. pentoxone (4-methoxy-4-methylpentanone-2)
11. cyclohexanone
12. dibutylketone
13. isophorone (3,5,5-trimethyl-2-cyclohexen-1-one)
14. acetophenone
15. methylpyrrolidone



[www.agilent.com/chem](http://www.agilent.com/chem)

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

Published in UK, August 03, 2010

A00178



**Agilent Technologies**