



Ketones, C₃ – C₈

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT U column separates seven C₃ to C₈ ketones in ten minutes.



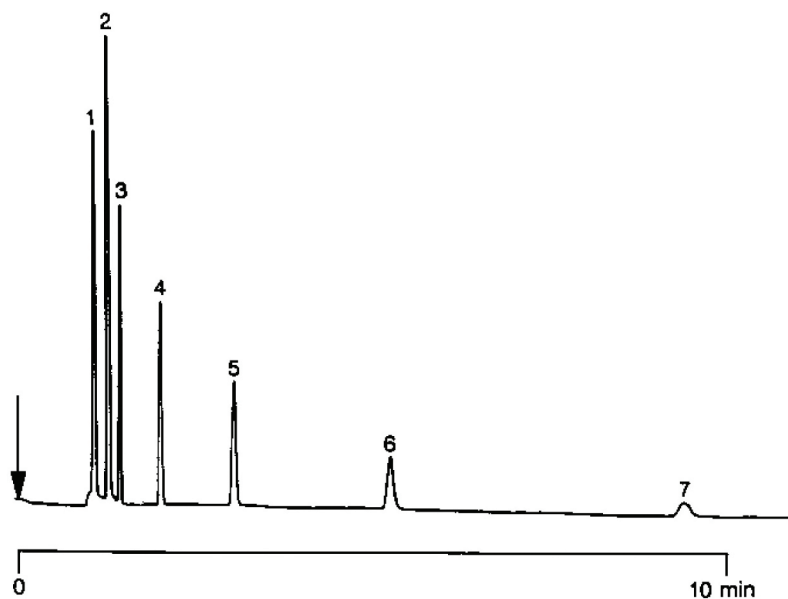
Agilent Technologies

Conditions

Technique : GC-Wide-bore
Column : Agilent PoraPLOT U, 0.53 mm x 25 m fused silica
PLOT PoraPLOT U (df = 20 μ m) (Part no. CP7584)
Temperature : 190 °C
Carrier Gas : H₂
Injector : Split, 100 mL/min
Detector : FID, 64 x 10⁻¹² Afs
Sample Size : 1% in H₂O

Peak identification

1. acetone
2. unknown
3. butanone
4. pentanone
5. hexanone
6. heptanone
7. octanone



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

A00682



Agilent Technologies