



Ketones

Analysis of impurities in methyl ethyl ketone

Application Note

Energy & Fuels

Authors

Agilent Technologies, Inc.

Introduction

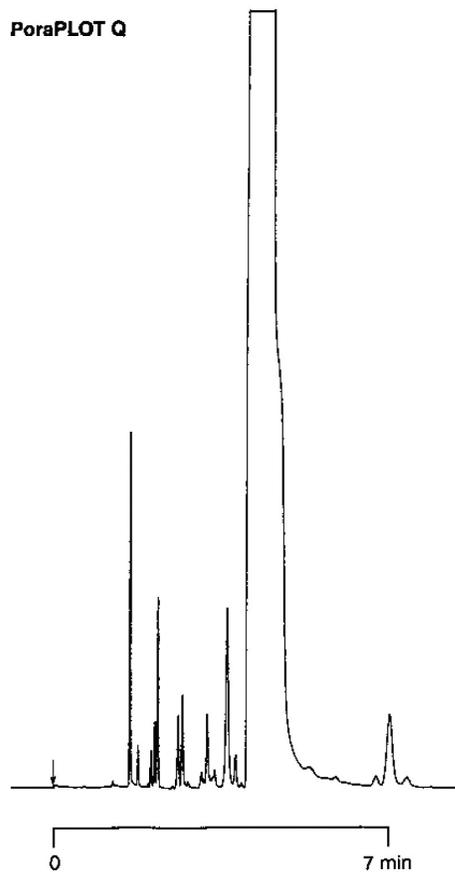
Gas chromatography with Agilent PoraPLOT Q and PoraPLOT S columns separate impurities in methyl ethyl ketone in seven minutes.



Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q and S, 0.53 mm x 25 m (ID x L)
fused silica PoraPLOT Q and S (df = 20 μ m)
(Part no. CP7554 and CP7574)
Temperature : 175 $^{\circ}$ C
Carrier Gas : H₂, 30 kPa (0.3 bar, 4.2 psi)
Injector : Split, 100 mL/min
Detector : FIO, 64 x 10⁻¹² Afs
Sample : pure MEK



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

A00454



Agilent Technologies