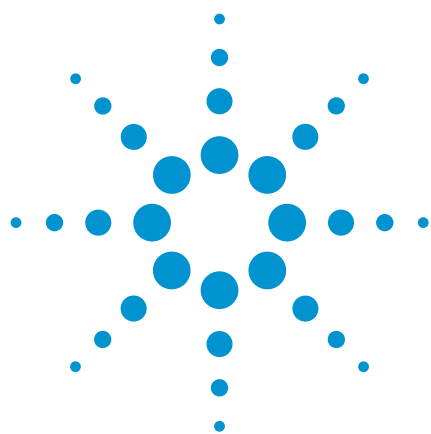




Hydrocarbons, $C_6 - C_7$

Separation of impurities in cyclohexane

Application Note

Materials Testing & Research

Authors

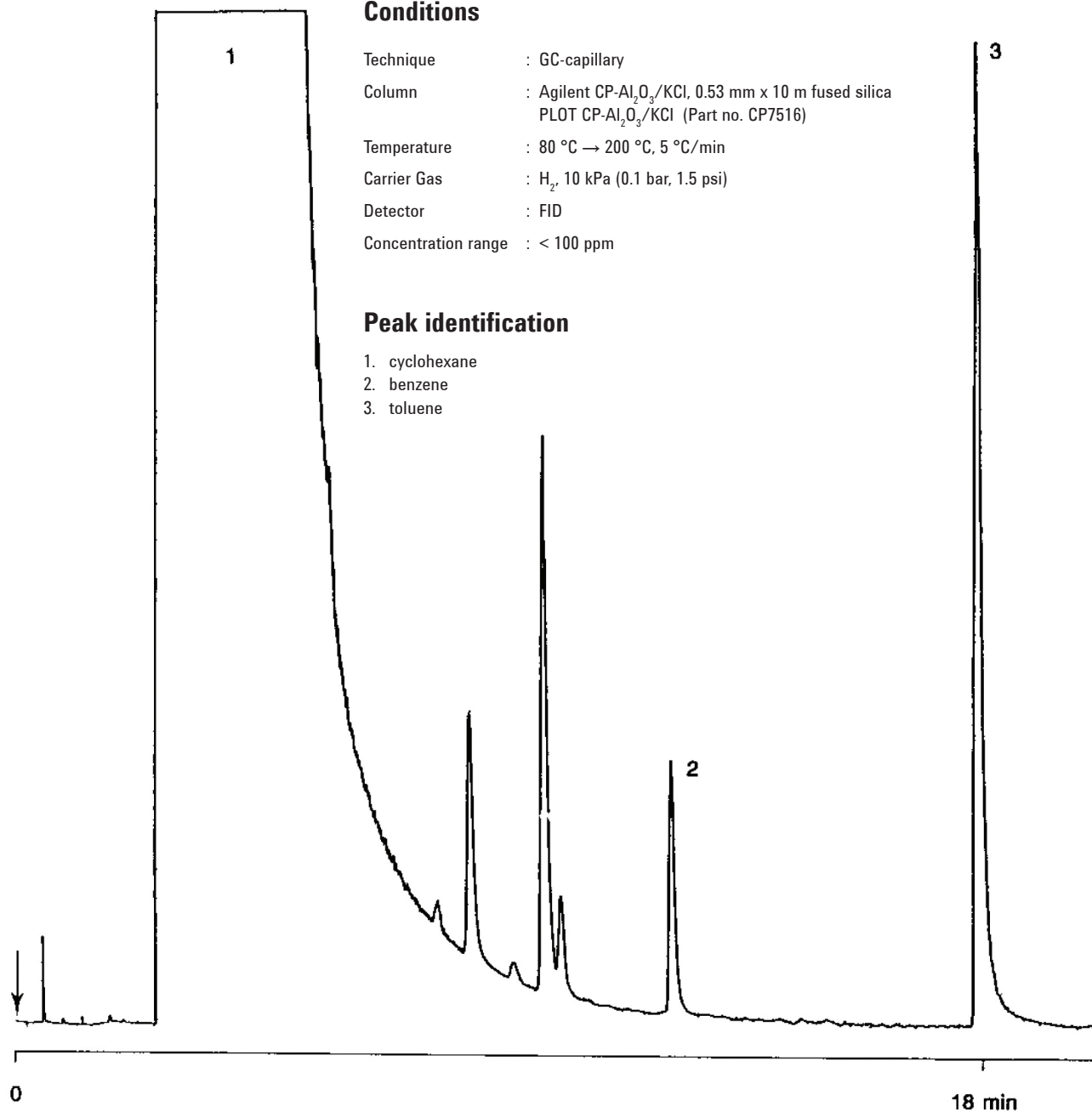
Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Al₂O₃/KCl column separates three C_6 to C_7 aromatic impurities in cyclohexane in 18 minutes.



Agilent Technologies



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

A00260



Agilent Technologies