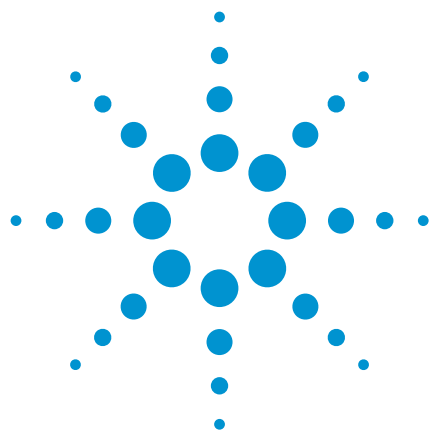




Hydrocarbons, $C_2 - C_4$

Determination of impurities in vinyl chloride

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Al₂O₃/KCl column identifies six C_2 to C_4 hydrocarbon impurities in vinyl chloride in 15 minutes.



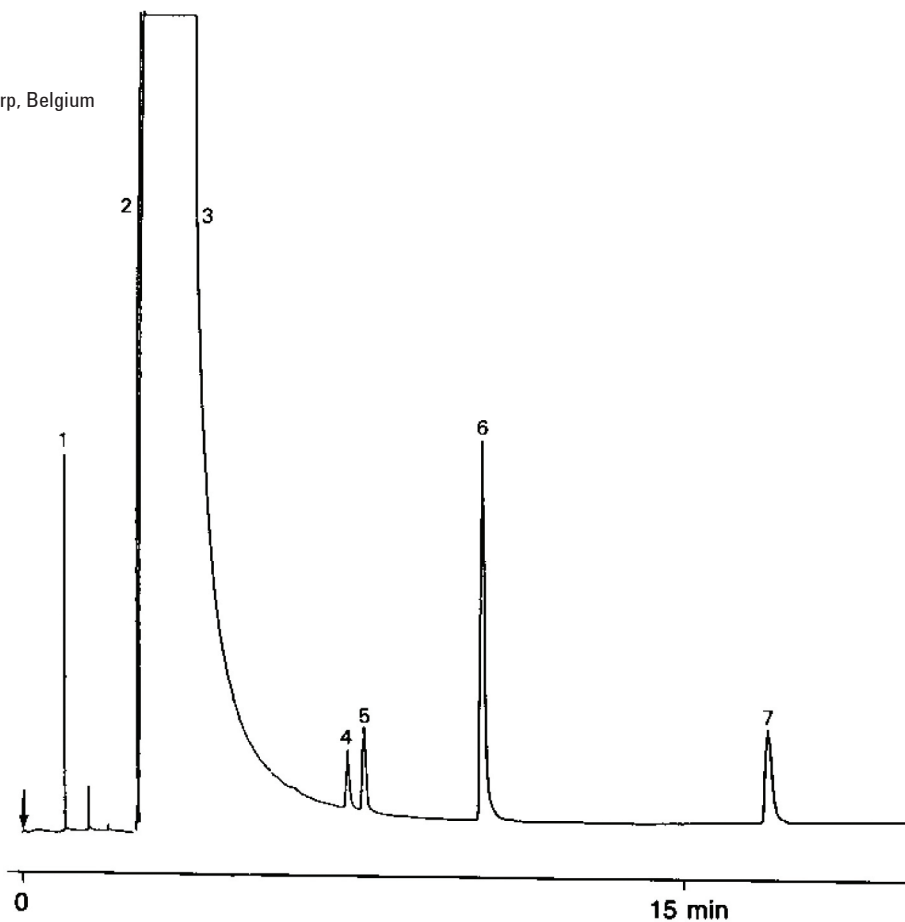
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Al₂O₃/KCl, 0.32 mm x 50 m fused silica
PLOT Al₂O₃/KCl (Part no. CP7515)
Temperature : 75 °C
Carrier Gas : H₂, 3 mL/min
Injector : 100 µL gas injection valve, Splitter
Detector : FID
Detection limit : ca. 0.1 ppm
Courtesy : Dr. R. Smits, BASF, Antwerp, Belgium

Peak identification

1. ethene
2. chloromethane
3. vinylchloride
4. vinylbromide
5. 1,3-butadiene
6. chloroethane
7. mono-vinylacetylene



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

A00270



Agilent Technologies