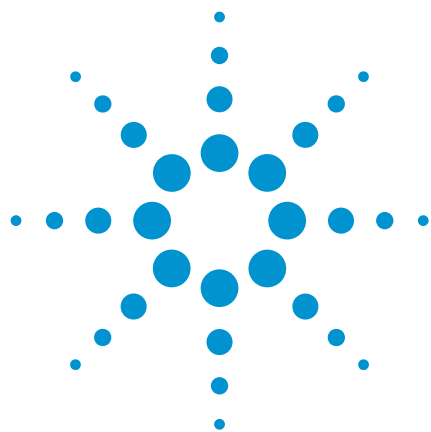




# Chlorinated hydrocarbons, C<sub>2</sub>

## Determination of impurities in vinyl chloride

### Application Note

Materials Testing & Research

#### Authors

Agilent Technologies, Inc.

#### Introduction

Gas chromatography with an Agilent CP-PoraPLOT Q column separates six impurities in vinyl chloride in two minutes.



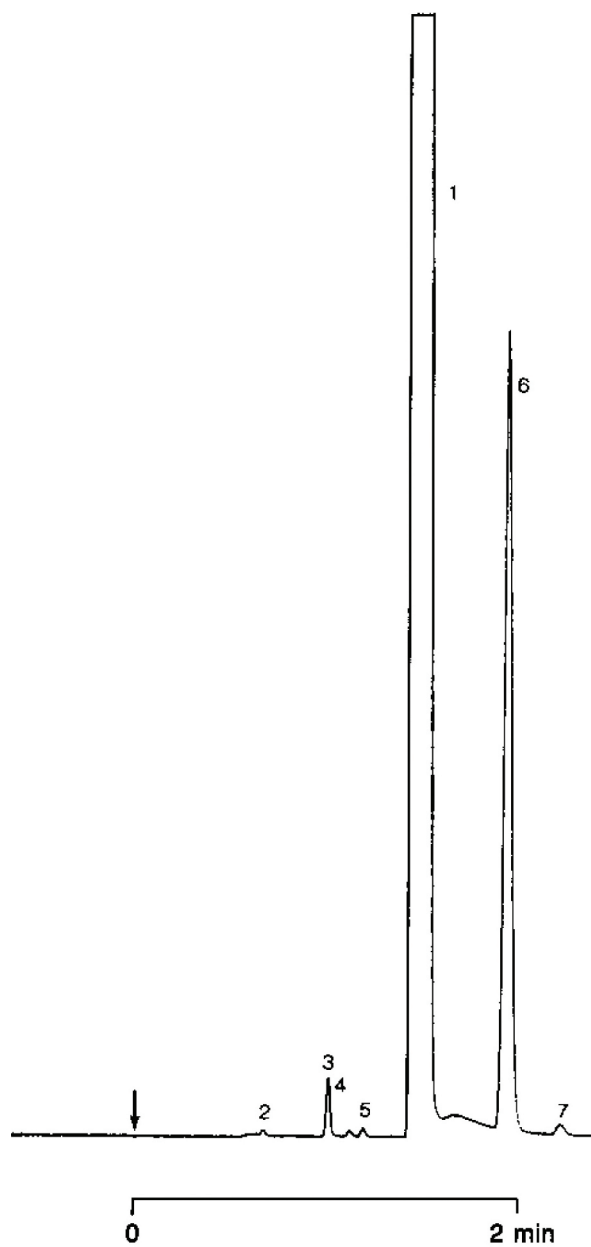
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-PoraPLOT Q, 0.32 mm x 10 m fused silica  
PoraPLOT Q (10 µm) (Part no. CP7550)  
Temperature : 120 °C  
Carrier Gas : H<sub>2</sub>, 40 kPa (0.4 bar, 5.8 psi)  
Injector : Splitter  
Detector : FID

## Peak identification

1. vinylchloride  
2-7. impurities



[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

First published prior to 11 May, 2010

A00275



**Agilent Technologies**