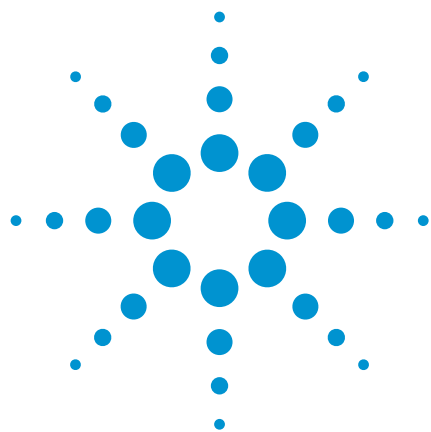




Chlorine/gases

Analysis of chlorine using its reaction with acetone

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB column is used to analyze chlorine by employing its reaction with acetone, giving results in ten minutes.



Agilent Technologies

Conditio

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 8 CB (df = 0.4 μ m) (Part no. CP7789)
Temperature : 60 °C $\rightarrow$ 150 °C, 10 °C/min
Carrier Gas : H₂
Injector : Splitter, 120 mL/min
Detector : FID
Sample Size : 0.1 μ L

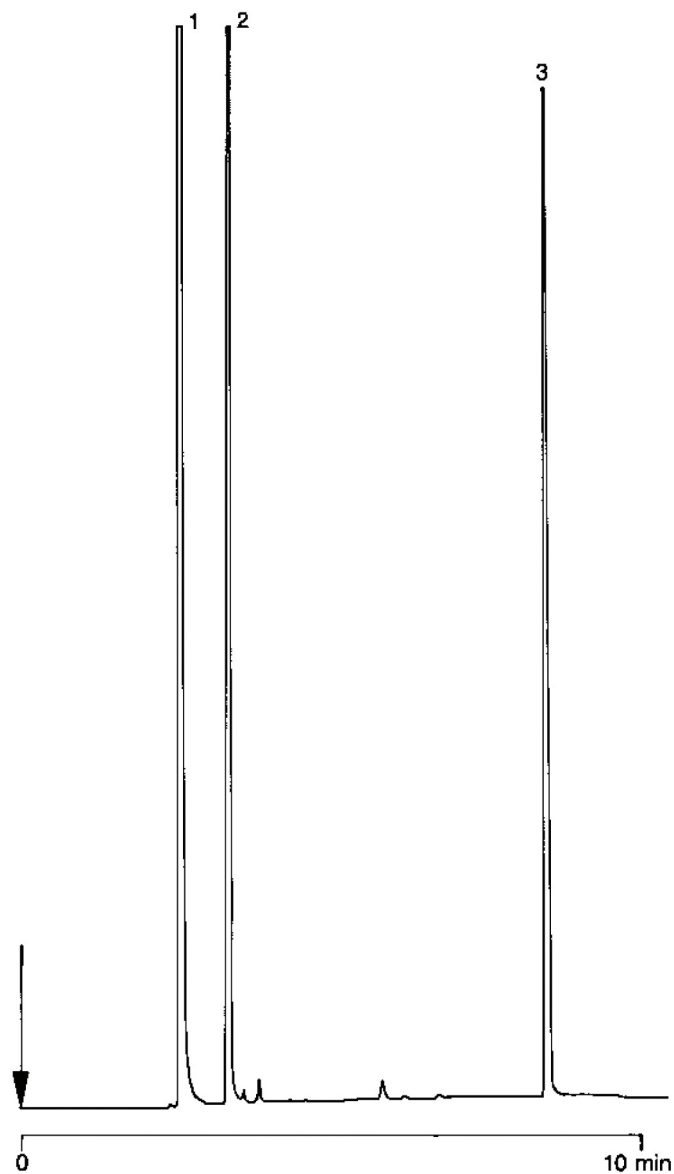
Sample preparation

Injection of 1 μ L of a sample of around
10 mL of acetone after reaction of 0.78 g
of chlorine and addition of 1 mL of n-decane
added as internal standard

Note: The reaction of chlorine with acetone
can be explosive so take care that you
don't mix too much chlorine with
acetone at the same time.

Peak identification

1. acetone
2. mono-chloro-acetone
3. decane



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

A00897



Agilent Technologies