



Solvents

Trace analysis of water in acetone

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT U column identifies trace water in acetone in five minutes.



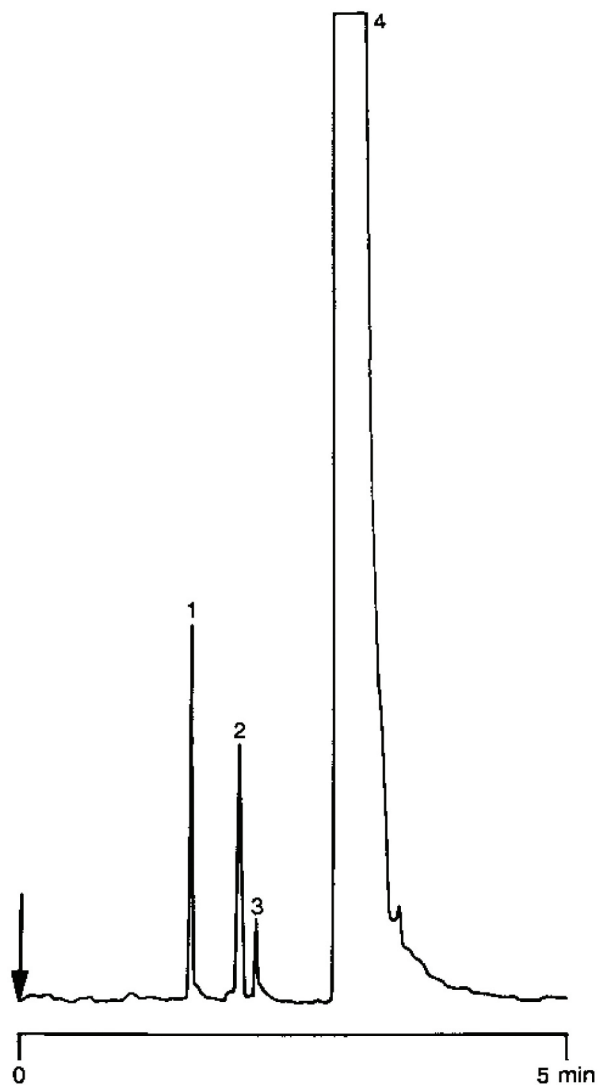
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT U, 0.32 mm x 25 m fused silica
PLOT PoraPLOT U (df = 10 μ m) (Part no. CP7581)
Temperature : 175 °C
Carrier Gas : H₂, 40 kPa (0.4 bar, 5.72 psi)
Injector : Splitter, 100 mL/min;
T = 200 °C
Detector : μ -TCD, 4 x 10
T = 100/180 °C
Sample Size : 0.5 μ L
Concentration Range : 500 ppm

Peak identification

| | |
|------------|---------|
| 1. air | |
| 2. water | 500 ppm |
| 3. unknown | |
| 4. acetone | (99.9%) |



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

A00669



Agilent Technologies