

# Solvents, C<sub>3</sub>

## Determination of water in acetone

### Application Note

Materials Testing & Research

#### Authors

Agilent Technologies, Inc.

#### Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates water from acetone in six minutes.



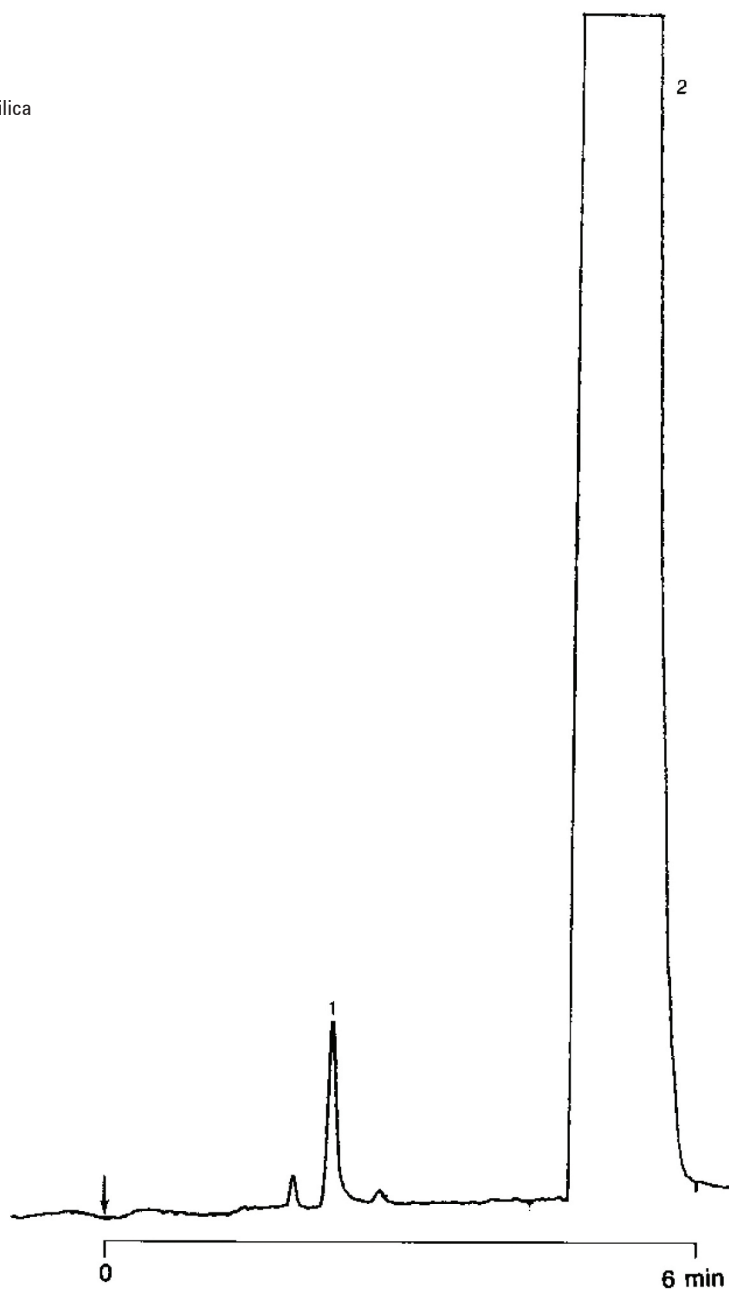
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent Pora-PLOT Q, 0.32 mm x 10 m fused silica  
PoraPLOT Q (10  $\mu$ m) (Part no. CP7550)  
Temperature : 175  $^{\circ}$ C  
Carrier Gas : H<sub>2</sub>, 50 kPa (0.5 bar, 7 psi)  
Injector : Splitter  
Detector :  $\mu$  TDC

## Peak identification

1. water (1000 ppm)
2. acetone



[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

A00286



**Agilent Technologies**