



Alcohols, C₂ – C₃

Application Note

Materials Testing & Research

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column determines 0.1% iso-propanol in ethanol in three minutes.



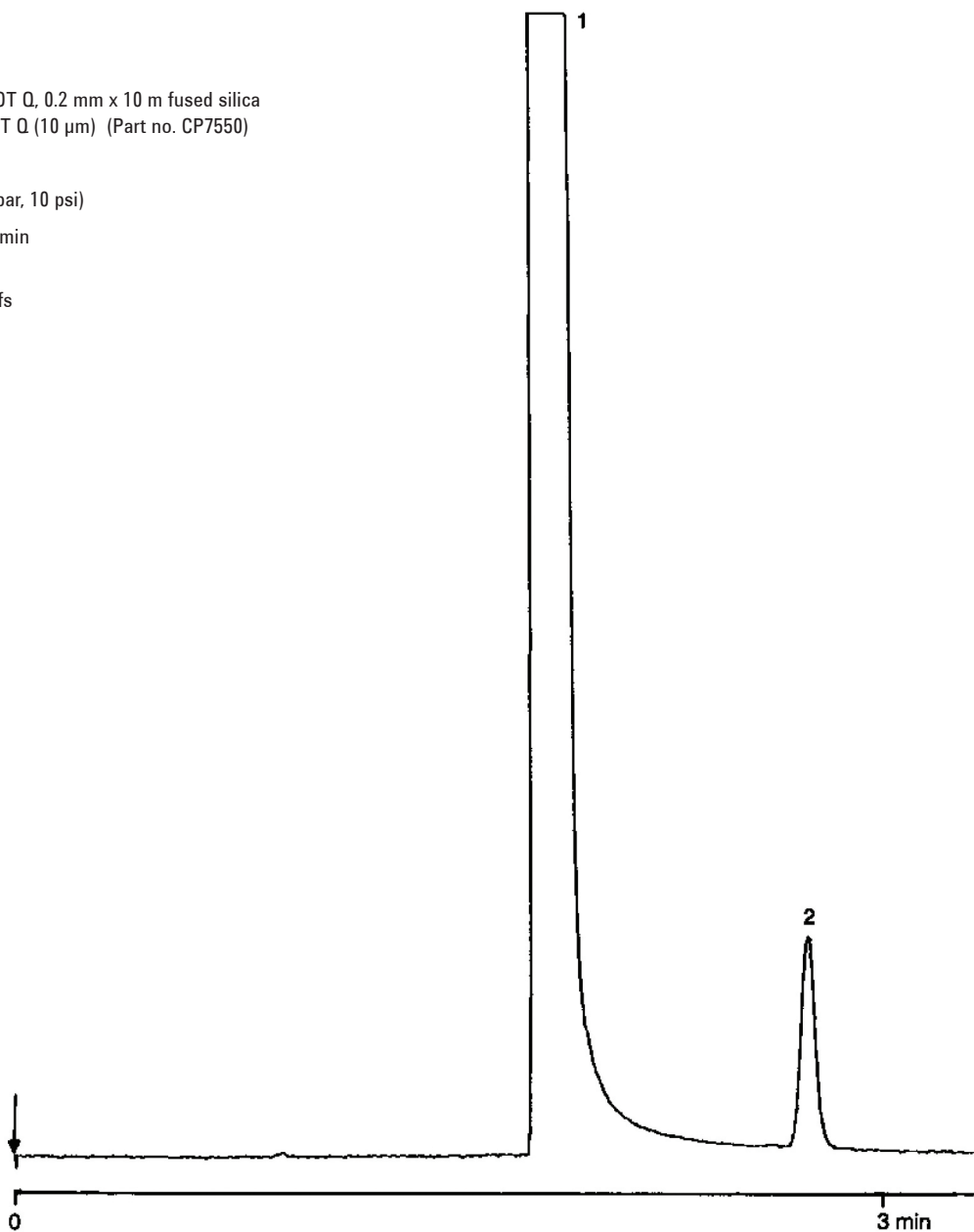
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 0.2 mm x 10 m fused silica
WCOT PoraPLOT Q (10 μ m) (Part no. CP7550)
Temperature : 150 °C
Carrier Gas : H₂, 70 kPa (0.7 bar, 10 psi)
Injector : Splitter, 50 mL/min
T = 250 °C
Detector : FID, 16 x 10⁻¹² Afs
T = 290 °C
Sample Size : 0.1 μ L
Concentration range : 0.1%

Peak identification

1. ethanol
2. iso-propanol



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

A00350



Agilent Technologies