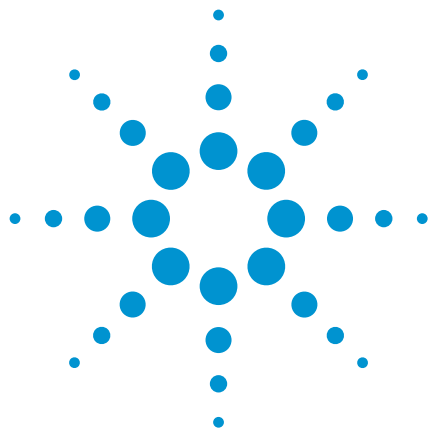




Residual solvents

Analysis of styrene in yoghurt

Application Note

Food Testing & Agriculture

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Wax 52 CB column separates residual solvents in yoghurt in 16 minutes.



Agilent Technologies

Conditions

Technique : GC-PTI

Column : Agilent CP-Wax 52 CB, 0.32 mm x 25 m fused silica
WCOT CP-Wax 52 CB (df = 1.2 µm)
(Part no. CP7763)

Temperature : 80 °C (5 min) → 210 °C, 15 °C/min; 210 °C (7 min)

Carrier gas : He, 50 kPa (0.5 bar, 7.2 psi)

Injector : TCT

Cold trap : CP-Sil 8 CB
0.53 mm; df = 5 µm

Precool temp. : -100 °C

Precool time : 3 min

Purge time : 10 min

Purge flow : 10 mL/min

Injection temp. : 200 °C

Injection time : 1 min

Backflush flow : 50 mL/min

Detector : FID
T = 250 °C

Sample Size : 25 mL, 1:1 dilution with water and
addition of anti-foaming agent

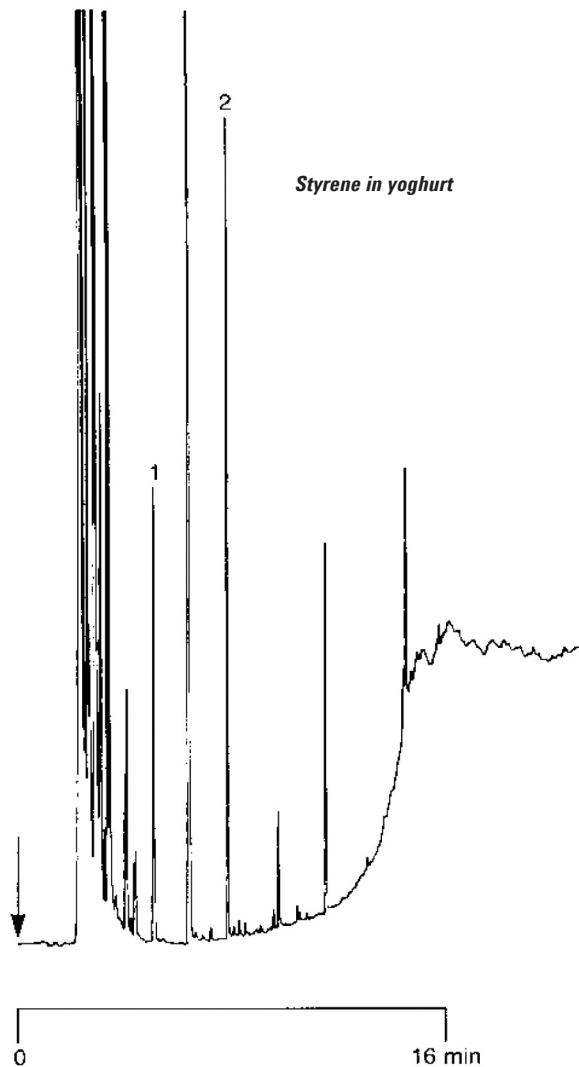
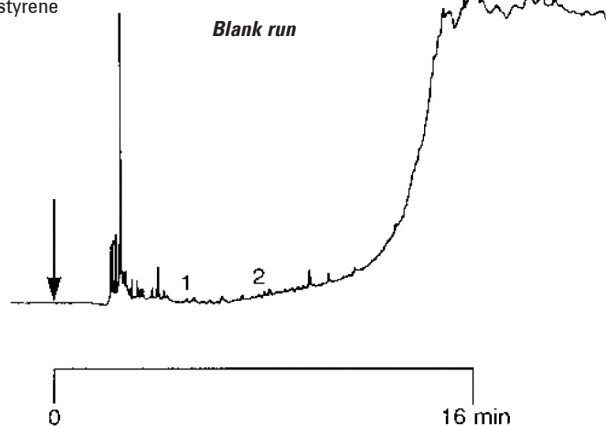
Sample Temperature : ambient

Concentration Range : ppb/ppm

Courtesy : DOW Benelux, Temeuzen, The Netherlands

Peak identification

1. ethylbenzene
2. styrene



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

A01046



Agilent Technologies