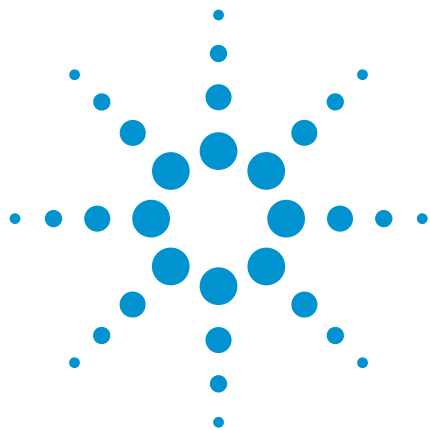




Herbicides

Separation of triazine herbicides

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 13 CB column separates 11 triazine herbicides in 15 minutes.



Agilent Technologies

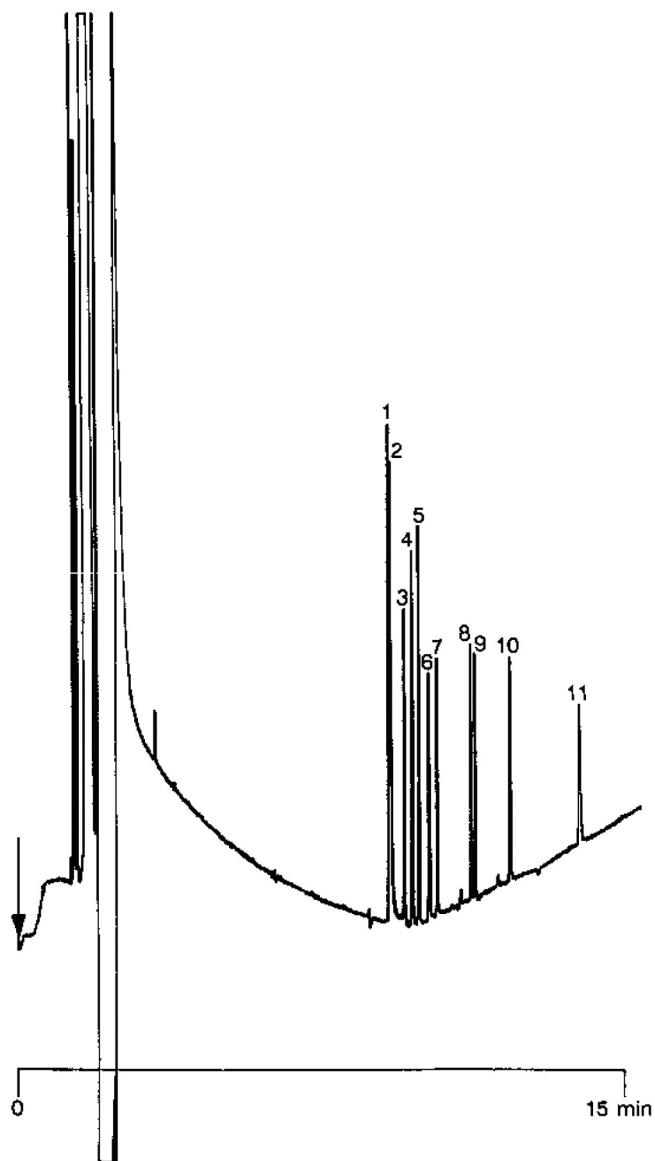
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 13 CB, 0.32 mm x 50 m fused silica
WCOT CP-Sil 13 CB (df = 0.4 µm) (Part no. CP7937)
Temperature : 80 °C → 270 °C, 20 °C/min
Carrier Gas : H₂, 100 kPa (1.0 bar, 14 psi)
Injector : On-column
Detector : NPD
T = 300 °C
Sample Size : 1 µL
Concentration Range : 100 pg/µL (100 ppb) per compound

Courtesy : Mr Lembacher,
Hipp K. G., Pfaffenhofen,
Germany

Peak identification

1. desethylatrazine
2. desisopropylatrazine
3. atraton
4. atrazine
5. trietazine
6. secbumeton
7. sebuthylazine
8. simetryn
9. dipropetryn
10. dimethametryn
11. methoprotryn



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

A00613



Agilent Technologies