



Herbicides

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 24 CB column separates 11 urea herbicides, derivatized with methyl iodide, in 50 minutes.



Agilent Technologies

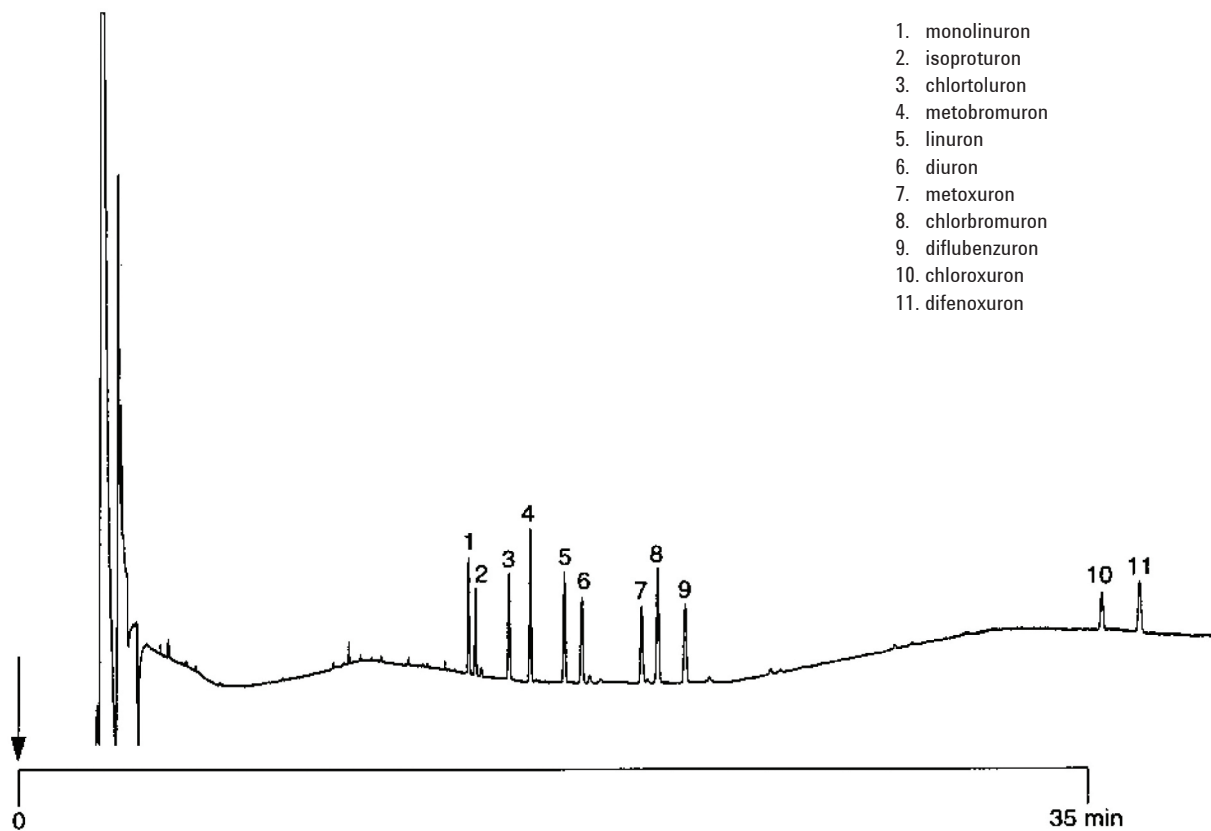
Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 24 CB, 0.32 mm x 30 m fused silica
WCOT CP-Sil 24 CB (df = 0.25 µm) (Part no. CP7831)
Temperature : 60 °C (1 min) → 205 °C, 15 °C/min
Carrier Gas : He, 50 kPa (0.5 bar, 7.2 psi)
Injector : Splitless
T = 220 °C
Detector : NPD
T = 300 °C
Sample Size : 3 µL

Courtesy : Ing. M. Vink
TNO, Division Analytical Sciences,
Zeist, The Netherlands

Peak identification

1. monolinuron
2. isoproturon
3. chlortoluron
4. metobromuron
5. linuron
6. diuron
7. metoxuron
8. chlorbromuron
9. diflubenzuron
10. chloroxuron
11. difenoxuron



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

A01228



Agilent Technologies