



Herbicides

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 24 CB column separates 11 triazine herbicides in 30 minutes.



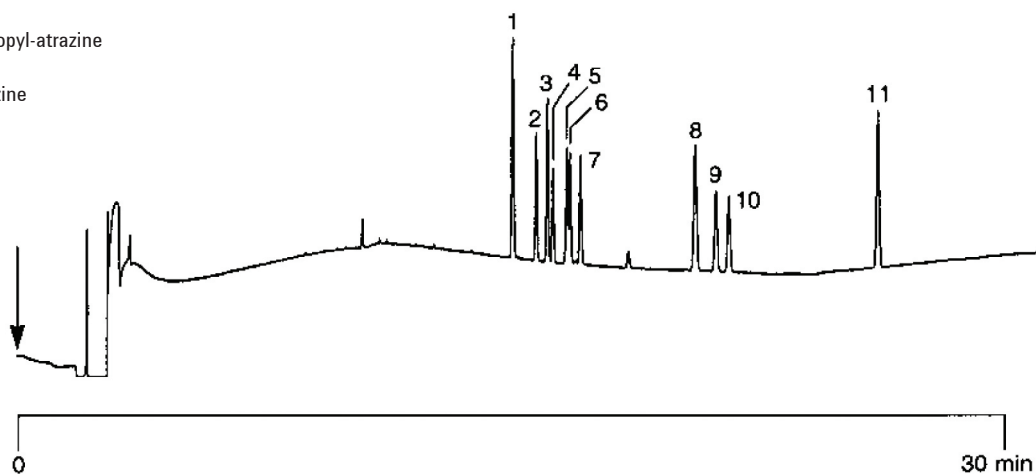
Agilent Technologies

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. DEDIPA | desethyl-desisopropyl-atrazine |
| 2. DEA | desethyl-atrazine |
| 3. DIPA | desisopropyl-atrazine |
| 4. propazine | |
| 5. atrazine | |
| 6. simazine | |
| 7. terbutylazine | |
| 8. prometryn | |
| 9. ametryn | |
| 10. terbutryn | |
| 11. cyanazine | |



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

A01227



Agilent Technologies