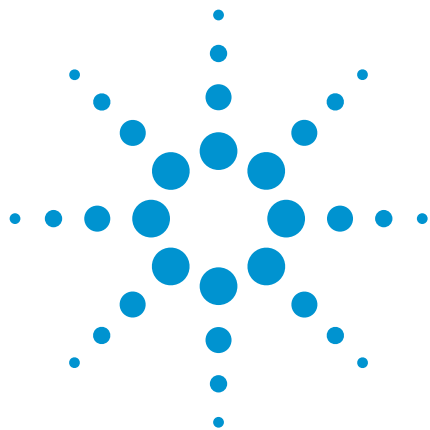




# Herbicides

## Application Note

Environmental

### Authors

Agilent Technologies, Inc.

### Introduction

Gas chromatography using an Agilent CP-Sil 24 CB column separates ten phenoxy herbicides in 34 minutes.



**Agilent Technologies**

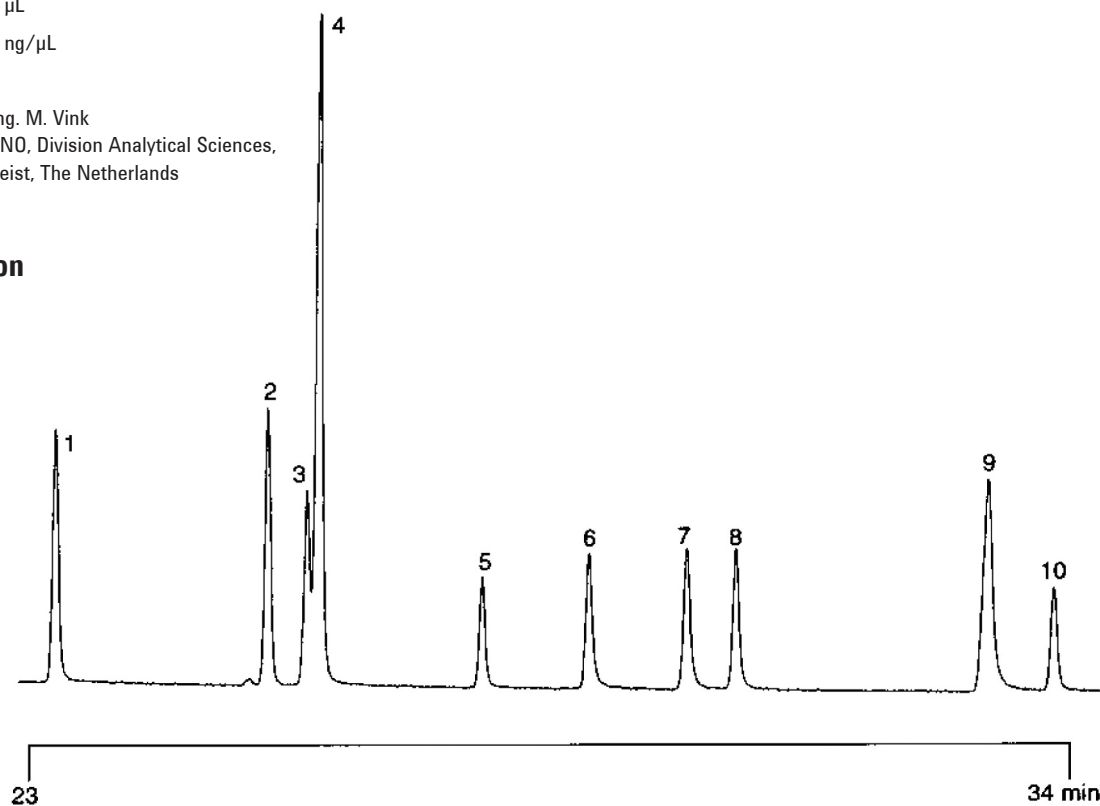
## Conditions

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) → 150 °C, 15 °C/min; 150 °C (0 min)  
→ 270 °C, (30 min)  
Carrier Gas : He, 60 kPa (0.6 bar, 8.6 psi)  
Injector : Splitless  
T = 220 °C  
Detector : NPD  
T = 300 °C  
Sample Size : 1 µL  
Concentration Range : 2 ng/µL

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

## Peak identification

1. MCPP
2. dicamba
3. MCPA
4. dichlorprop (2,4-DP)
5. dikegulac
6. 2,4-D
7. 2,4,5-TP
8. 2,4,5-T
9. MCPB
10. bentazon



[www.agilent.com/chem](http://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

A01230



**Agilent Technologies**