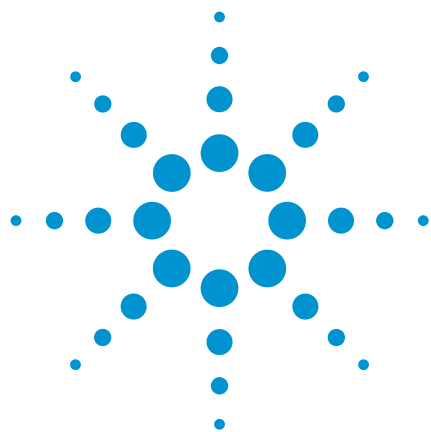




Chlorinated herbicides

Analysis of chlorinated herbicides to EPA 515 and EPA 8150

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates ten chlorinated herbicides according to EPA 515 and EPA 8150 in 25 minutes.



Agilent Technologies

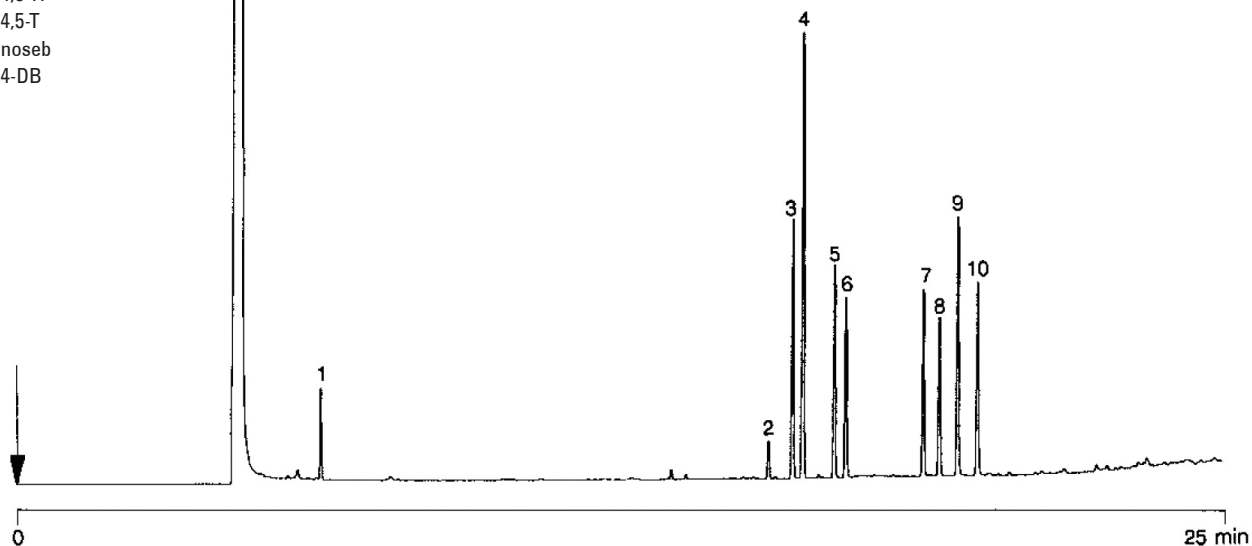
Conditions

Technique	: GC-capillary
Column	: Agilent CP-Sil 8 CB for pesticides, 0.25 mm x 50 m fused silica WCOT CP-Sil 8 CB (df = 0.12 µm) with retention gap (Part no. CP7481)
Temperature	: 60 °C → 300 °C, 10 °C/min
Carrier Gas	: He, 150 kPa (1.5 bar, 21 psi)
Injector	: Split, 100 mL/min T = 275 °C
Detector	: FID T = 300 °C

Peak identification

Methyl esters of

1. dalapon
2. dicamba
3. MCPP
4. MCPA
5. dichlorprop
6. 2,4-D
7. 2,4,5-TP
8. 2,4,5-T
9. dinoseb
10. 2,4-DB



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

A00803



Agilent Technologies