



Herbicides

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

GC/MS with an Agilent CP-Sil 8 CB Low Bleed/MS column separates seven polar N-herbicides in 25 minutes.



Agilent Technologies

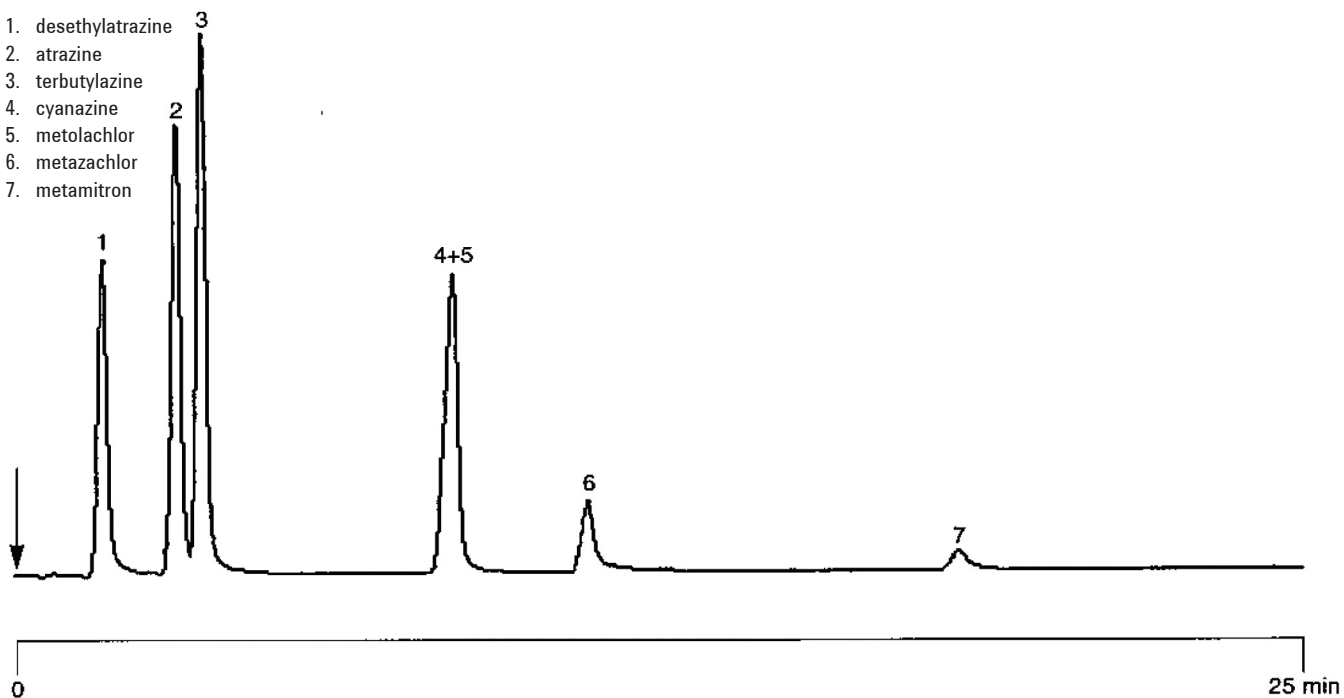
Conditions

Technique : GC-capillary
Column : CP-Sil 8 CB low bleed/MS, 0.25 mm x 60 m fused silica WCOT CP-Sil 8 CB (df = 0.251 μm) (Part no. CP5861)
Temperature : 30 °C \rightarrow 270 °C, 25 °C/min; 270 °C \rightarrow 300 °C, 25 °C/min; 300 °C (13 min)
Carrier Gas : He, 100 kPa (1.0 bar, 14 psi)
Injector : On-column
Detector : MS
Concentration Range : 5 ng
Solvent Sample : diethyl ether

Courtesy : H. v. Cauwenbergh,
Vlaamse Maatschappij voor Watervoorziening,
Haasrode., Belgium

Peak identification

1. desethylatrazine
2. atrazine
3. terbutylazine
4. cyanazine
5. metolachlor
6. metazachlor
7. metamitron



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

A01111



Agilent Technologies