



Chlorophenoxycarboxylic acids

Application Note

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 24 CB column resolves 22 methylated chlorophenoxycarboxylic acids in 35 minutes.



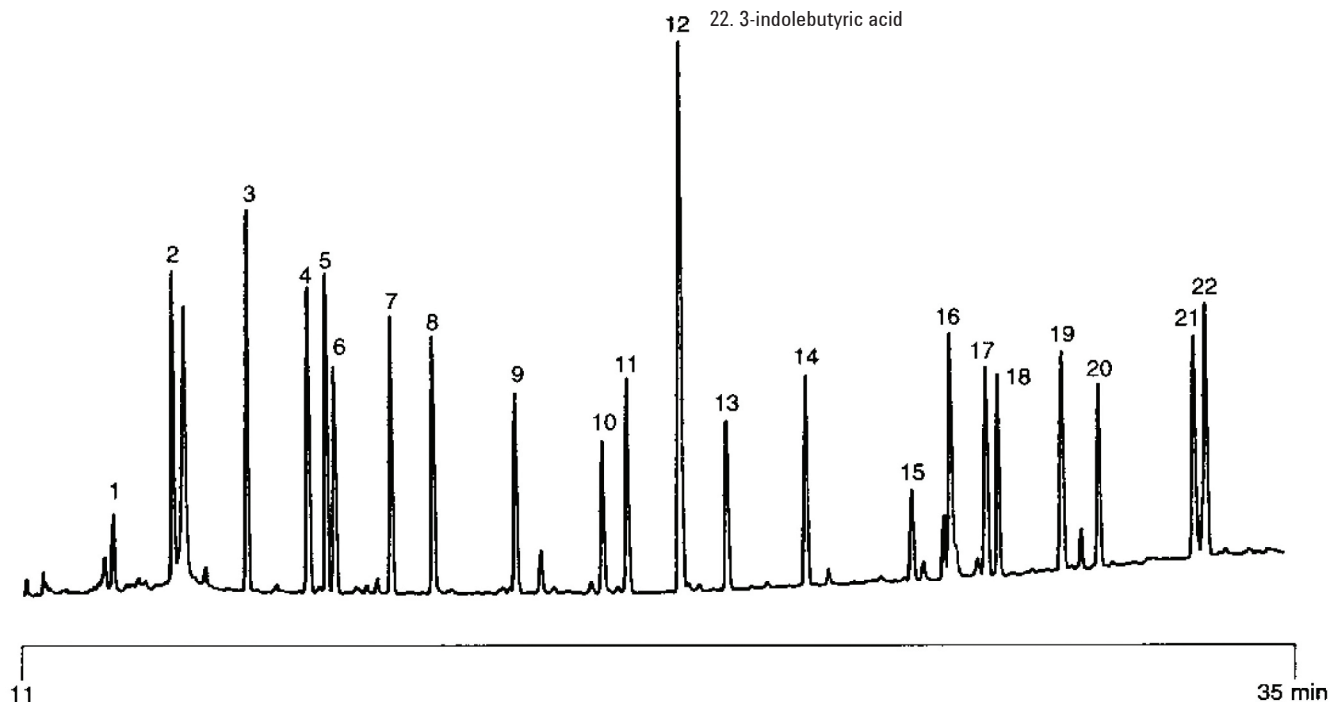
Agilent Technologies

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 (2 min) → 150 °C, 15 °C/min;
150 °C → 260 °C, 3.5 °C/min
Carrier Gas : H₂
Injector : 60 s Splitless, then with a Split of 50 mL/min
T = 275 °C
Detector : FID
T = 325 °C
Injection volume : < 1 ng per compound

Peak identification

1. dibromobenzoic acid
2. 2,4-dichlorobenzoic acid
3. clofibric acid
4. 4-chlorophenoxyacetic acid
5. MCPP
6. dicamba
7. MCPA
8. 2,4-DP (dichlorprop)
9. 2,4-D
10. triclopyr
11. 2,4,5-TP
12. MCPB + naphthylacetic acid
13. 2,4,5-T
14. 2,4-DB
15. fluroxypur
16. indoleacetic acid
17. bentazon
18. fluazitop
19. flurenol
20. haloxyfop
21. benazolin
22. 3-indolebutyric acid



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

A01090



Agilent Technologies