

Analysis of Ylang Oil using Agilent J&W FactorFour VF-WAXms GC Columns

Application Note

Author

Helena Jacobse
Agilent Technologies, Inc.

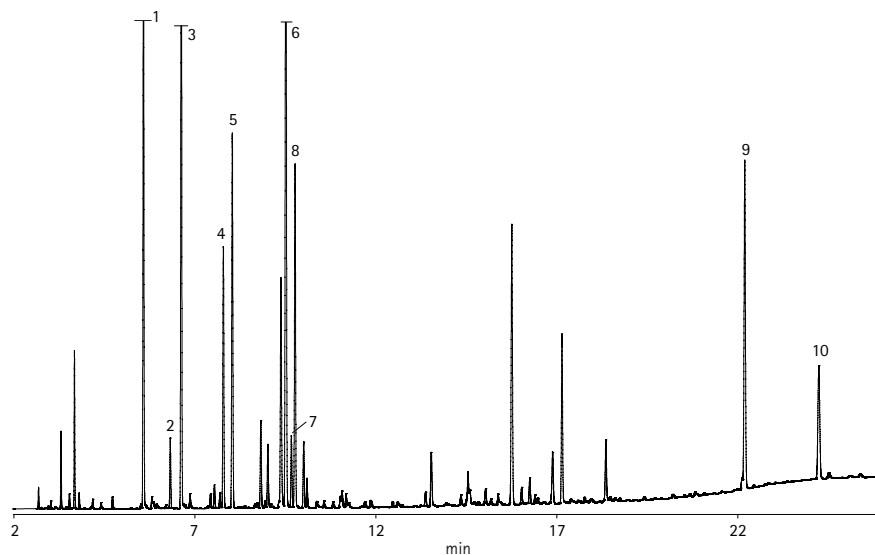
Conditions

Technique: GC-FID
Sample: Ylang oil, 0.1 % (acetone)
Sample Size: 1.0 µL
Column: VF-WAXms 30 m x 0.25 mm df =
0.25 µm (part number CP9205)

Carrier Gas: Helium, 1.0 bar
Injector: 250 °C, split 75 mL/min
Temperature: 100 °C to 260 °C with 7 °C/min
Detection: 275 °C, FID

Peak Identification:

1. Benzene, 1-methoxy-4-methyl-
2. *a*-Copaene
3. Linalool
4. *b*-Caryophyllene
5. Methyl benzoate
6. Benzyl acetate
7. *a*-Farnesene
8. Geranyl acetate
9. Benzyl benzoate
10. Benzyl salicylate



Analysis of ylang oil using VF-WAXms GC columns

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2010

Published in UK, October 08, 2010

SI-00945



Agilent Technologies