



Solvents

Analysis of USP 467 solvents

Application Note

BioPharma

Authors

Agilent Technologies, Inc.

Introduction

GC analysis of 14 USP 467 solvents on an Agilent FactorFour VF-200ms column takes 15 minutes.



Agilent Technologies

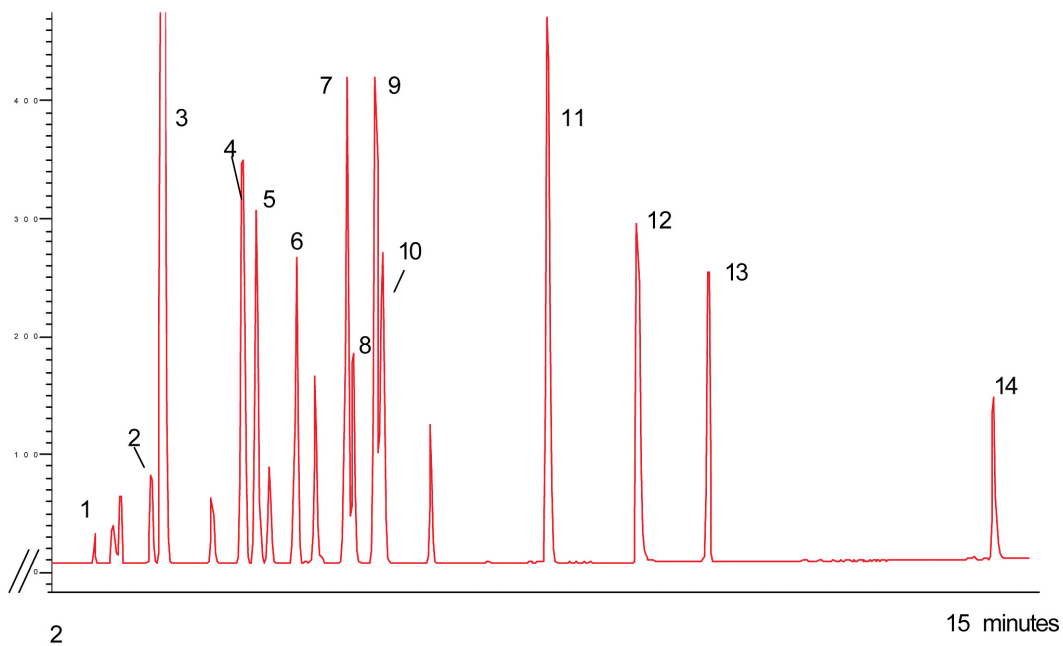
Conditions

Technique : GC
Column : Agilent FactorFour VF-200ms, 0.25 mm x 30 m
(df = 0.25 µm) (Part no. CP8858)
Temperature : 45 °C, 10 °C/min → 325 °C
Carrier Gas : Helium, ca. 1.0 mL/min
Pressure program : 60 kPa
Injector : Split/Splitless, in split mode, 1:100
Detector : FID
Sample Size : 1 µL
Solvent : DMSO, 0.1% w/v

Courtesy : Jane Peene, Agilent Application laboratory,
Middelburg, The Netherlands

Peak identification

1. pentane
2. 2-propanol
3. hexane
4. chloroform
5. cyclohexane
6. heptane
7. benzene
8. tetrahydrofuran
9. trichloroethylene
10. ethylacetate
11. toluene
12. pyridine
13. MEK methyl isobutylketone
14. n,n-dimethylformamide



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

A02204



Agilent Technologies