



Halogenated hydrocarbons

Anesthetic gases by headspace analysis

Application Note

BioPharma

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent PoraPLOT Q column separates four anesthetic gases in 18 minutes.



Agilent Technologies

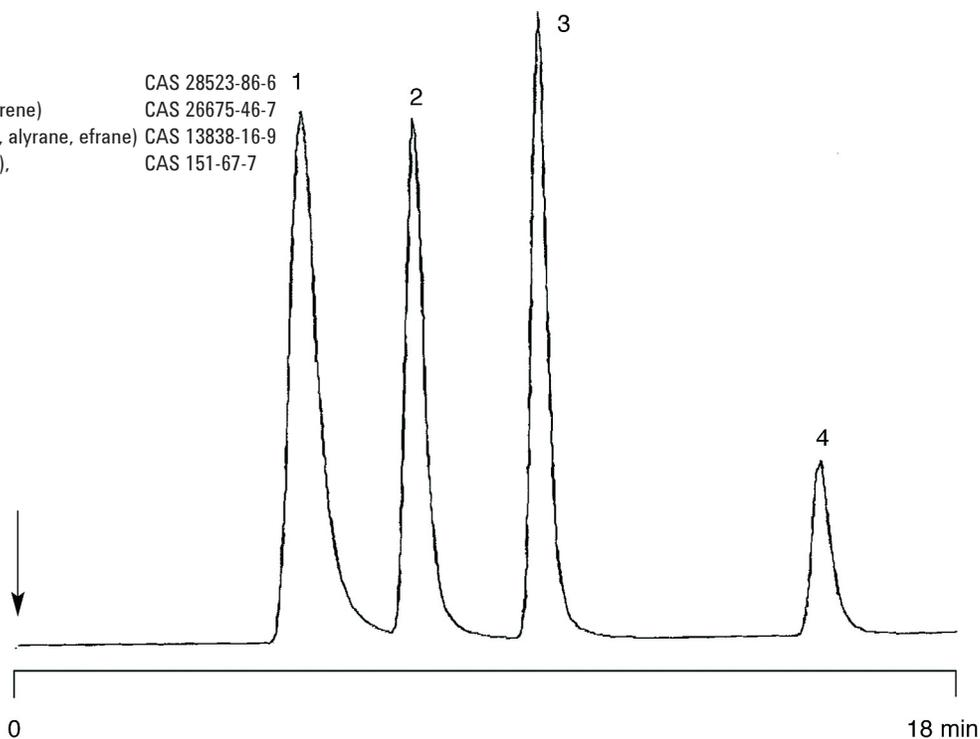
Conditions

Technique : GC-capillary
Column : Agilent PoraPLOT Q, 032 mm x 25 m fused silica
PLOT (df = 10 μ m) (Part no. CP7551)
Temperature : 40 °C (2 min) \rightarrow 150 °C, 10 °C/min :
150 °C (0 min) \rightarrow 200 °C, 6 °C/min
Carrier Gas : He, 10 mL/min, 180 kPa (1.8 bar, 24 psi)
Injector : Headspace
T = 150 °C
Detector : MS, TIC,
T = 250 °C
Sample Size : 1 mL

Courtesy : Dr. Laura Zaratina, Fondazione S. Maugeri, Padova,
Italy

Peak identification

- | | | | | |
|---------------|---|--|----------------|---|
| 1. sevofrane | (C ₄ H ₃ F ₇ O) | | CAS 28523-86-6 | 1 |
| 2. isoflurane | (C ₃ H ₂ ClF ₅ O) (forane, forene) | | CAS 26675-46-7 | 2 |
| 3. ethrane | (C ₃ H ₂ ClF ₅ O) (enflurane, alyrane, efrane) | | CAS 13838-16-9 | 3 |
| 4. halothane | (C ₂ HBrClF ₃) (fluothane), | | CAS 151-67-7 | 4 |



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

A01627



Agilent Technologies