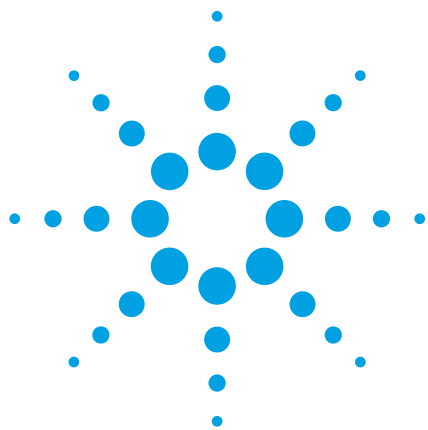




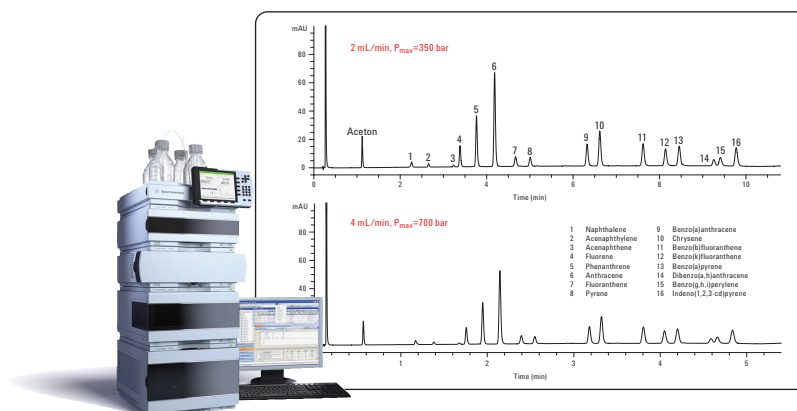
Fast analysis of polyaromatic hydrocarbons using the Agilent 1290 Infinity LC and Eclipse PAH columns

Application Note

Environmental

Author

Gerd Vanhoenacker
Research Institute for Chromatography
Kennedypark 26, B-8500 Kortrijk,
Belgium



The Agilent 1290 Infinity LC has a broader power range (the combination of pressure and flow capabilities) than any other commercially available system. This is extremely useful for method transfer from one (U)HPLC to the Agilent 1290 Infinity LC system and allows the analyst to develop methods that are impossible to run on these other systems.

The flow and pressure capabilities are illustrated by a separation of 16 polyaromatic hydrocarbons (PAHs) at high pressure and flow rate. At 2 mL/min, the analysis time is approximately 11 min. Doubling the flow rate and gradient speed allows the sample to be analyzed in 5.5 min with a maximum pressure of 700 bar. The combination of high flow (4 mL/min) and pressure is useful in this case to increase the sample throughput. The separation of the PAHs is shown in Figure 1.



Agilent Technologies

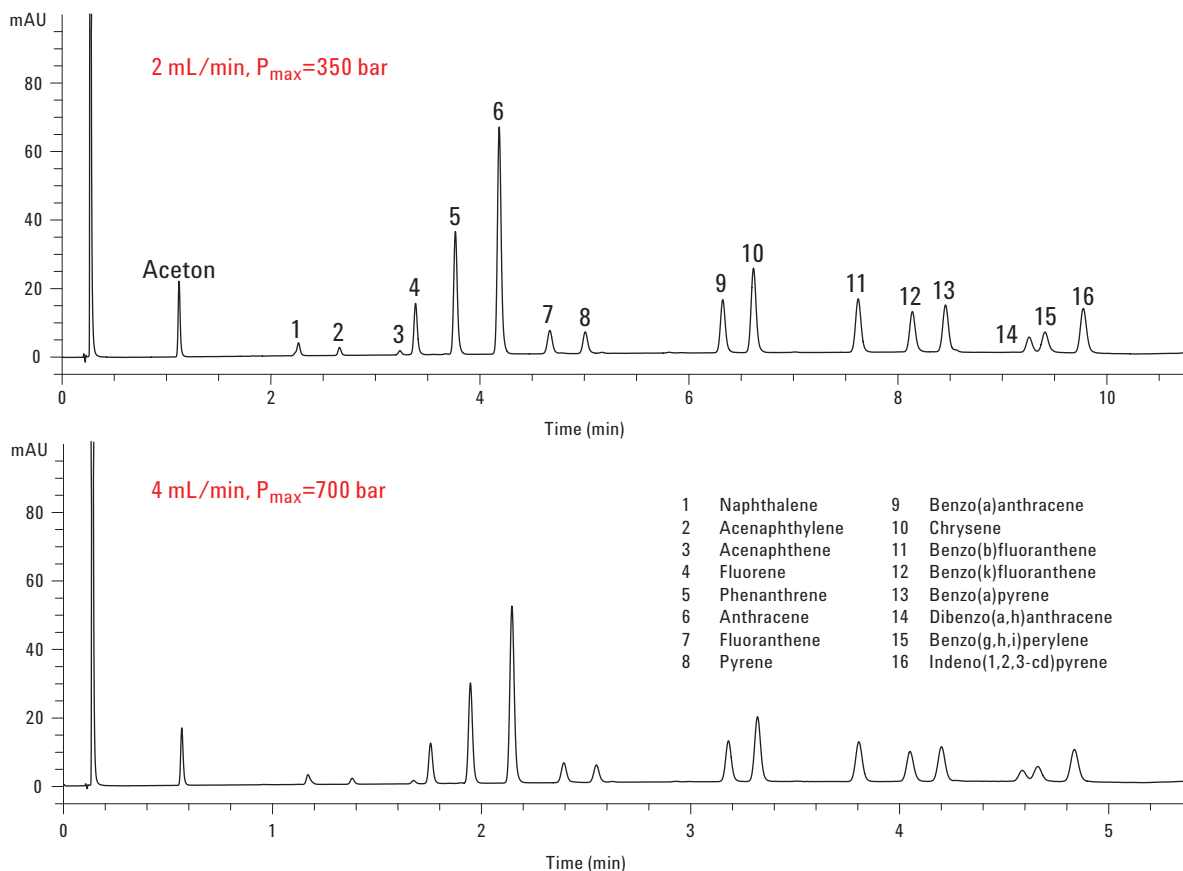


Figure 1

Analysis of 16 PAHs on the 1290 Infinity LC. Sample: standard solution of 16 PAHs, 50 µg/mL each.

Configuration:

- G4220A 1290 Infinity Binary Pump with Integrated Vacuum Degasser
- G4226A 1290 Infinity Autosampler
- G1316C 1290 Infinity Thermostatted Column Compartment
- G4212A 1290 Infinity Diode Array Detector

Method:

Column: ZORBAX Eclipse PAH 4.6 mm × 50 mm, 1.8 µm
 Mobile phase: A = water, B = acetonitrile
 Flow rate and gradient: 2 mL/min 0–0.33 min 40% B
 0.33–10 min 40–100% B
 4 mL/min 0–0.17 min 40% B
 0.17–5 min 40–100% B
 Injection volume: 0.2 µL
 Detector: Sig = 254/10 nm, Ref = off, 40 Hz
 Temperature: 25 °C

www.agilent.com/chem

© Agilent Technologies, Inc., 2009
 December 1, 2009
 Publication Number 5990-4934EN



Agilent Technologies