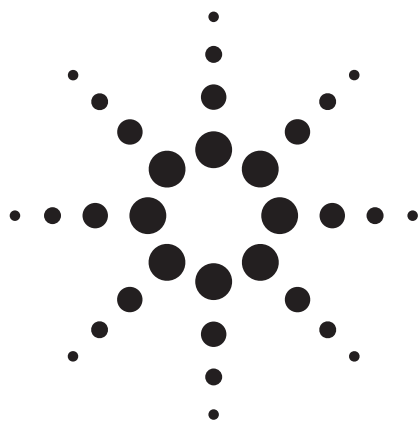




使用 QuEChERS/dSPE 样品制备方法和高效 DB-5ms 超高惰性 GC 色谱柱通过 GC/MS 分析鱼体内的多环芳烃 (PAH)

Application Note

Food Testing and Hydrocarbon Processing

Authors

Doris Smith and Ken Lynam
Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19809-1610
USA

Abstract

本应用摘要详细说明了一种快捷有效的分析方法，可用于检测鱼样品中少量及痕量的多环芳烃 (PAH)，该方法可替代之前耗时更长、技术更复杂的程序。简化的 QuEChERS（快速、简便、经济、高效、耐用和安全）方法与分散固相萃取 (dSPE) 相结合，可以在保持低浓度分析物检测能力的同时充分净化样品基质。柱后安装有压力控制三通以实现自动反吹，从而减少样品残留并且降低仪器循环次数。

通过安捷伦 J&W DB-5ms 超高惰性 (UI) 色谱柱 (20 m×0.18 mm, 0.18 μm) 成功分离出 16 种目标 PAH。实验在 10、25、50、100、250、500 和 1000 ng/mL 的 PAH 水平上对使用选择性离子监测 (SIM) 的 GC/MS 方法进行了校准，所得线性和重现性均非常出色。回收率测定的加标水平为 25、250 和 500 ng/mL。回收率范围为 80% 至 139%，RSD 低于 6%。



Agilent Technologies

Introduction

Oil spills can introduce massive amounts of crude oil and the myriad of components found in the oil, including Polycyclic Aromatic Hydrocarbons (PAHs). PAHs are a broad class of molecules that are well known, well characterized, and a worldwide focal point for regulations as both food contaminants and environmental pollutants [1-2]. PAHs have known toxicity and in some cases have been implicated as carcinogens. These molecules are persistent in the environment and bio-accumulate in fatty tissue of fish species throughout the food chain [3]. Fortunately, from an analyst's perspective, there are historical accounts of large oil spills and their impact on aquatic species and the environment that serve as case studies for investigation of future spills [4].

When large scale oil spills occur, regulatory agencies spring to action to minimize the risk of human exposure, contain the leak at its source when possible, and protect the environment as much as feasible. In the US, the National Oceanic and Atmospheric Administration (NOAA) in coordination with the US Food and Drug Administration (FDA) and other federal and state agencies set policies for both fishery closure and the reopening of affected fisheries. As part of the reopening process for a closed fishery NOAA and the FDA have established testing procedures for PAH analyses as a requirement to reopen a fishery. Levels of concern for eight targeted PAHs and their alkylated homologues have been defined for finfish, oysters, shrimp and crab species. The established levels of concern range from a high of 233 ppm for anthracene/phenanthrene in shrimp and crabs to a low of 0.03 ppm in finfish for benzo(a)pyrene [5].

Analysis at these levels can be reliably achieved using a capillary GC/MS approach. Resolution of anthracene and phenanthrene can also be routinely achieved using an Agilent J&W DB-5ms column and appropriate instrument conditions. Use of the Ultra Inert column in this case assures less interaction between the column and the sample matrix.

Sample preparation to execute the NOAA method has historically been labor intensive involving two successive steps of preparative chromatography. In this note the use of a QuEChERS (Quick Easy Cheap Effective and Rugged) approach to sample preparation is presented for use as a screening method for PAHs. Typically sample preparation using the NOAA method takes 12-14 hours per sample. Using a QuEChERS sample preparation approach, up to 60 samples can be processed in one eight-hour shift [6]. Sample preparation time savings are substantial for the QuEChERS approach enabling higher laboratory throughput for the screening of PAHs in seafood.

The DB-5ms column has been widely used for US-EPA PAHs and offers suitable resolution of the sixteen targeted PAHs investigated by this method. The high efficiency DB-5ms Ultra Inert column was chosen for this application because of its added benefits for trace level analysis in more complex samples matrixes while retaining the same selectivity [7]. A 0.18 mm id or high efficiency GC column can also help by providing faster sample analysis than a 0.25 mm id column typically used for GC/MS analysis [8-9].

The GC/MS system used was also equipped with backflush capability. This capability enables faster instrument cycle time by backflushing late eluting matrix components back through the inlet purge valve. Long bake out times between injections are avoided by using this technique. Backflushing has the additional benefit of increasing the time intervals for source cleaning effectively clearing deleterious matrix components from the system [10-11].

Experimental

An Agilent 7890 GC and Agilent 5975B GC/MS System equipped with a multimode inlet (MMI) and Agilent 7693 automatic liquid sampler were used for this series of experiments. The GC was also fitted with a pressure controlled tee (PCT) post-column for automated backflush. Table 1 lists the chromatographic conditions used for these analyses. Table 2 lists flow path consumable supplies used in these experiments.

Table 1. Chromatographic Conditions

GC/MSD:	Agilent 7890 GC/Agilent 5975B GC/MS System
Sampler:	Agilent 7693 automatic liquid sampler, 5.0 µL syringe (Agilent p/n 5181-1273)
PCT Device:	Purged Ultimate Union (Agilent p/n G3186-60580)
Carrier:	Helium, constant flow 1.7 mL/min
Restrictor:	0.7 m × 0.15 mm id deactivated silica tubing
PCM 1:	3.8 psi constant pressure
MMI:	0.5 µL splitless; 320 °C, Purge flow 50 mL/min at 0.8 min Gas saver 30 mL/min at 2 min
Column:	Agilent J&W DB-5msUI 20 m × 0.18 mm, 0.18 µm (Agilent p/n 122-5522UI)
Oven:	50 °C (0.4 min), 25 °C/min to 195 °C (1.5 min), 8 °C/min to 265 °C, 20 °C/min to 315 °C (1.25 min)
Postrun Backflush:	7 min at 315 °C, backflush pressure 70 psi, 2 psi inlet pressure during backflush
MSD:	340 °C transfer line, 340 °C source, 150 °C quad

Table 2 Flow Path Supplies

Vials:	Amber screw top glass vials (Agilent p/n 5183-2072)
Vial Caps:	Blue screw caps (Agilent p/n 5182-0717)
Vial inserts:	100 μ L glass/polymer feet (Agilent p/n 5181-8872)
Syringe:	5 μ L (Agilent p/n 5181-1273)
Septum:	Advanced green (Agilent p/n 5183-4759)
Inlet liners:	Deactivated dual taper Helix liner (Agilent p/n G5188-5398)
Ferrules:	0.4 mm id short; 85/15 vespel/graphite (Agilent p/n 5181-3323)
PCT fittings:	Internal nut (Agilent p/n G2855-20530)
PCT ferrules:	SilTite ferrules, 0.25 mm id (Agilent p/n 5188-5361)
20x magnifier :	20x Magnifier loop (Agilent p/n 430-1020)

Reagents and Chemicals

All reagents and solvents were HPLC or Ultra Resi grade. Acetonitrile (ACN) was from Honeywell (Muskegon, MI, USA), and acetone was from VWR International (West Chester, PA, USA). The 16-component PAH standard used was obtained from Agilent (p/n 8500-6035).

Solutions and Standards

The PAH stock standard solution (502 μ g/mL of 16 polycyclic aromatic hydrocarbons) was diluted in acetone to yield spiking solutions of 0.5, 1, 5, 10 and 50 μ g/mL. The spiking solutions were used to prepare the calibration curves in the matrix blank extract by appropriate dilution.

Sample Preparation

A red snapper fish sample was purchased from a local grocery store. The fish was chopped into small cubes and frozen at -80°C overnight. The samples were then comminuted thoroughly to achieve sample homogeneity. The sample extraction method used the QuEChERS method followed by dSPE [5]. Figure 1 illustrates the sample preparation procedure graphically in a flow chart.

Samples containing 3.0 grams of fish were weighed into centrifuge tubes. QC samples were spiked with an appropriate amount of PAH spiking solution to yield QC samples with concentrations of 25, 250, and 500 ng/mL. Each sample received 12.0 mL aliquot of deionized water and 15 mL aliquot of ACN. Two ceramic bars (Agilent p/n 5982-9313) were added to each sample to aid in sample extraction. The samples were vortexed for 1 minute. An Agilent original QuEChERS extraction salt packet (Agilent p/n 5982-5555) containing 6 grams of MgSO_4 and 1.5 grams sodium chloride was added to each centrifuge tube. The capped tubes were shaken on a Geno

Grinder at 1500 rpm for 1 minute. The samples were centrifuged at 4000 rpm for 5 min.

An 8 mL aliquot of the upper layer was transferred to an Agilent QuEChERS fatty sample dispersive SPE 15 mL tube (Agilent p/n 5982-5158). The dSPE tube was vortexed for 1 minute and then centrifuged at 4000 rpm for 5 minutes to complete the sample extraction. The liquid from the dSPE tube was transferred to a GC vial and analyzed by SIM GC/MS using the chromatographic conditions listed in Table 1.

Extractions of water and acetonitrile aliquots were prepared in the same manner as the samples and served as a reagent blanks.

QuEChERS/dSPE Sample Preparation Workflow

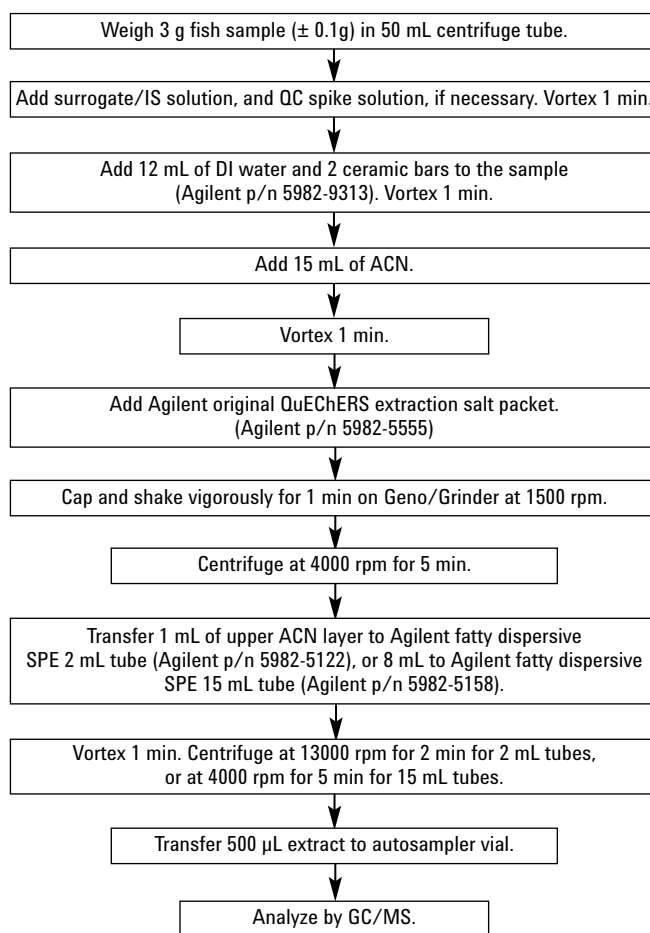


Figure 1. Flow chart of the Agilent SampliQ QuEChERS modified extraction procedure for fish sample.

Discussion of Results

The sixteen PAH analytes were resolved on the Agilent J&W DB-5msUI 20 m × 0.18 mm, 0.18 µm analysis column in less than 20 minutes. Figure 2 shows the separation of a 500 ng/mL PAH standard solution. Quantitative GC/MS analysis of the target PAHs was performed using selective ion monitoring (SIM), which gave excellent sensitivity down to 10 ppb for all the analytes.

The performance of the DB-5msUI high efficiency column yielded excellent linearity and recovery over the calibration range of this study. The linearity of the column as defined by the r^2 values of the PAH standard curve ranged from 0.9988-0.9999. The individual PAH analyte values are shown in Table 3. Excellent signal-to-noise ratios were also seen for the lowest calibration standard on the column as shown in Figure 3. The method limit of quantitation (MLQ) of 10 ppb for benzo[a]pyrene is well below the level of concern of 30 ppb set by the NOAA and FDA.

Table 3. The r^2 Values for the PAH Calibration Standards Over the 10 ng/mL to 1000 ng/mL Range of this Study

Analyte	R^2
Naphthalene	0.9997
Acenaphthylene	0.9999
Acenaphthalene	0.9999
Fluorene	0.9999
Phenanthrene	0.9999
Anthracene	0.9999
Fluoranthene	0.9999
Pyrene	0.9998
Benzo[a]anthracene	0.9993
Chrysene	0.9994
Benzo[b]fluoranthene	0.9990
Benzo[k]fluoranthene	0.9988
Benzo[a]pyrene	0.9992
Indeno[1,2,3-c,d]pyrene	0.9992
Dibenz[a,h]anthracene	0.9991
Benzo[g,h,i]perylene	0.9995

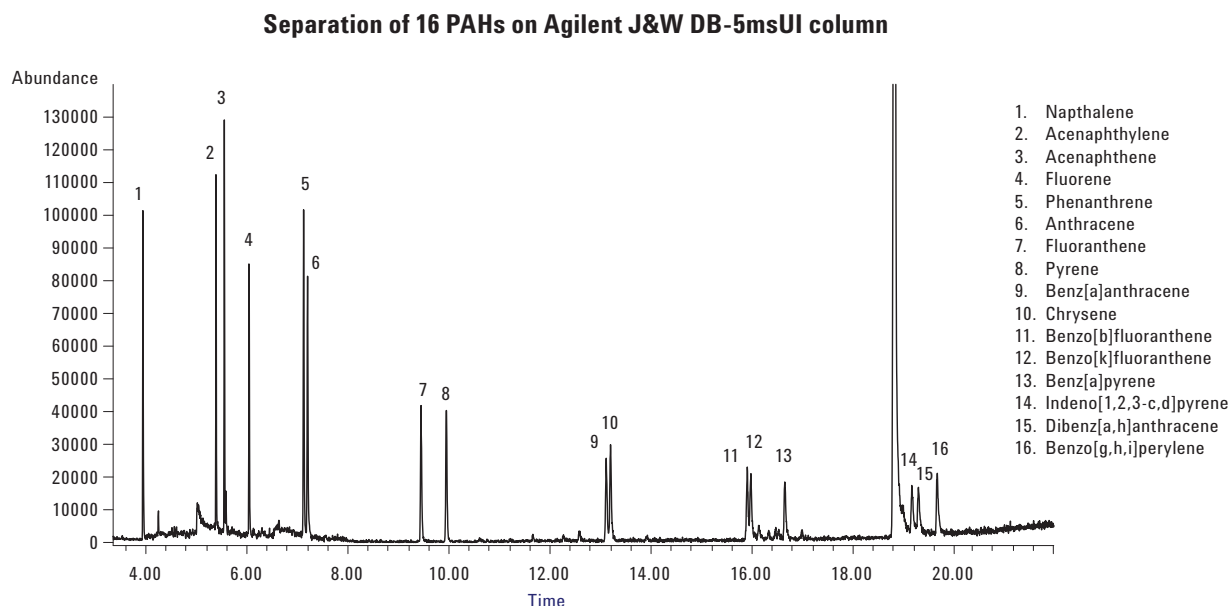


Figure 2. GC/MS chromatogram of the 500 ng/mL PAH standard prepared in sample matrix analyzed on an Agilent J&W DB-5msUI 20 m × 0.18 mm, 0.18 µm capillary GC column (Agilent p/n 122-5522UI). Chromatographic conditions are listed in Table 1.

Excellent Signal-to-Noise Ratios at Trace Levels

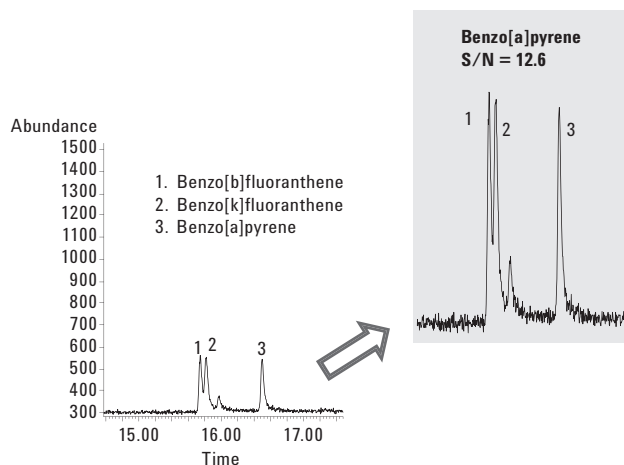


Figure 3. Enlarged view chromatogram of the benzo[a]pyrene peak in the 10 ng/mL PAH calibration standard prepared in sample matrix analyzed on the Agilent J&W DB-5msUI capillary column (Agilent p/n 122-5522UI). Chromatographic conditions are listed in Table 1.

The extraction process using the QuEChERS method followed by dispersive SPE was effective in retaining the PAHs in the spiked fish sample and providing sufficient cleanup of the sample matrix for GC/MS analysis. Figure 4 shows the separation of the extracted PAHs in a spiked fish sample on the DB-5ms UI column.

The recoveries were determined at the 25, 250, and 500 ng/mL PAH levels. Recoveries for the individual PAHs are shown in Table 4. The recoveries ranges (80% to 139%) and RSDs were excellent with the DB-5ms UI column for all PAHs investigated.

GC/MS SIM Chromatogram of Red Snapper Fish Extracts Blank Relative to Spiked Sample after Agilent's QuEChERS Extraction and Dispersive SPE

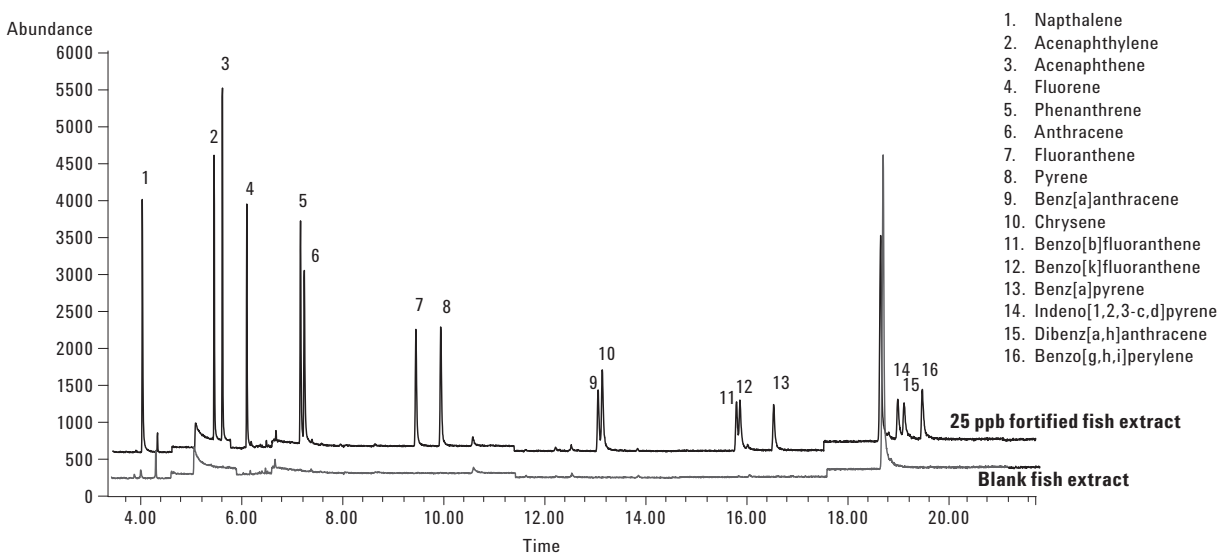


Figure 4. GC/MS SIM chromatogram of the fish extract blank and the 25 ng/mL spiked fish extract analyzed on Agilent J&W DB-5msUI capillary column (Agilent p/n 122-5522UI). Chromatographic conditions are listed in Table 1.

Table 4. Recovery and Repeatability of PAHs in Spiked Red Snapper Fish with Agilent J&W DB-5msUI Column (Agilent p/n 122-5522UI)

Analyte	25 ng/mL fortifield QC		250 ng/mL fortifield QC		500 ng/mL fortifield QC	
	% Recovery	RSD (n=6)	% Recovery	RSD (n=6)	% Recovery	RSD (n=6)
Naphthalene	80.35	3.29	96.77	4.23	98.64	1.88
Acenaphthylene	95.28	2.30	103.36	2.80	101.02	2.27
Acenaphthalene	92.28	2.51	101.18	2.87	100.69	2.34
Fluorene	95.98	2.99	105.94	2.82	105.00	1.28
Phenanthrene	100.51	3.46	104.93	2.71	103.25	1.70
Anthracene	107.38	3.51	105.95	3.45	105.38	1.74
Fluoranthene	113.27	3.87	105.76	3.33	103.64	1.81
Pyrene	113.55	3.51	103.99	3.24	102.29	1.94
Benz[a]anthracene	129.79	3.41	101.45	3.91	100.61	3.24
Chrysene	116.75	4.01	98.55	4.17	95.95	5.61
Benzo[b]fluoranthene	131.20	3.70	98.77	4.08	98.08	3.24
Benzo[k]fluoranthene	139.45	2.52	99.13	3.98	95.31	4.54
Benzo[a]pyrene	125.30	3.68	95.33	3.89	96.82	1.80
Indeno[1,2,3-c,d]pyrene	119.51	3.47	94.57	3.23	93.71	2.55
Dibenz[a,h]anthracene	126.35	3.54	98.55	3.50	98.85	2.24
Benzo[g,h,i]perylene	114.91	4.93	97.30	3.37	95.63	1.83

Conclusions

This application note successfully shows a quick and efficient analytical method for monitoring low and trace level PAHs in fish samples suitable to address current food safety concerns. This method demonstrates the feasibility of a simplified approach for routine fish screening as an alternative to more time consuming and complex techniques.

The Agilent QuEChERS method followed by dSPE for fatty samples was effective at providing enough sample cleanup to avoid matrix interferences while still maintaining low level analyte detection. The simple QuEChERS extraction method allows faster sample prep, facilitating higher sample throughput. Any residual sample matrix carryover is removed through use of backflush, which eliminates the need for a bakeout cycle, and significantly reduces analytical run times.

The Agilent J&W DB-5ms UI High Efficiency column was effective at analyzing 16 PAHs in a fish matrix following sample matrix cleanup by QuEChERS and dSPE. The DB-5ms UI provided satisfactory resolution of the four known critical pairs; phenanthrene/anthracene, benz[a]anthracene/chrysene, benzo[b]/[k]fluoranthene, and indeno[1,2,3-c,d]pyrene/dibenz[a,h]anthracene. The performance of the DB-5ms UI yielded excellent linearity over the range of concentrations studied with r^2 values between 0.9988-0.9999 for the PAH compounds. Recovery and reproducibility was shown to be greater than 80% with an RSD below 6.0. The method limit of quantitation of 10 ppb for the benz[a]pyrene using this approach was significantly lower than current levels of concern.

Acknowledgements

The authors wish to thank Joan Stevens for her help and suggestions with the Agilent QuEChERS extraction procedure.

References

1. Environmental Protection Agency (EPA) of the United States of America, Compendium Method TO-13A, EPA, Cincinnati, OH, USA, 1999.
2. European Union, Commission Recommendation 2005/108/EC, Off. J. Eur. Comm. L34 (2005) 433.
3. Haruhiko Nakata, Yasufumi Sakai, Takashi Miyawaki, and Akira Takemura "Bioaccumulation and Toxic Potencies of Polychlorinated Biphenyls and Polycyclic Aromatic Hydrocarbons in Tidal Flat and Coastal Ecosystems of the Ariake Sea, Japan," Environmental Science and Technology, 2003,37, 3513-3521.
4. Zhendi Wand and Scott A. Stout, "Oil Spill Environmental Forensics: Finger Printing and Source Identification," Academic Press, Burlington, MA 01803, USA, 2007.
5. NOAA Cleared Document Dated 5/29/2010 "Protocol for Interpretation and Use of Sensory Testing and Analytical Chemistry Results from Re-Opening Oil-Impacted Areas Closed to Seafood Harvesting."
6. AOAC Method 2007.01
7. Kenneth Lynam and Doris Smith, "Polycyclic Aromatic Hydrocarbon (PAH) Analysis Using an Agilent J&W DB-5ms Ultra Inert Capillary GC Column," Agilent Technologies publication 5989-9181EN
8. Kenneth Lynam and Mike Szelewski, "Analysis of Semivolatiles using High Efficiency Capillary GC Columns," Agilent Technologies publication 5989-7500EN
9. To download Agilent Method Translation software please visit the link below:
<http://www.chem.agilent.com/cag/servsup/usersoft/files/GCTS.htm>
10. Chin Kai Meng, "Improving Productivity and Extending Column Life with Backflush," Agilent Technologies publication 5989-6018EN
11. Harry Prest, "User Quick Guide to Pressure Controlled Tee (PCT) Operation – Post Run Backflushing," Agilent Technologies publication 5990-5484EN

For More Information

For more information on our products and services, visit our Web site at www.agilent.com/chem.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2010
Printed in the USA
November 2, 2010
Publication Number 5990-6668EN



Agilent Technologies