



Fast Analysis of Illicit Drug Residues on Currency using Agilent Poroshell 120

Application Note

Forensics and Toxicology

Authors

Anne E. Mack, James R. Evans and
William J. Long
Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808
USA

Abstract

Illicit drugs, like cocaine, are frequently found on US currency. While a more interesting perception might be that all bills were used to inhale the drug, the truth is much more mundane. Drug trafficking is thought to be the initial source of drug residues on a small percentage of bills, and because these compounds are fine powders, they are easily transferable from one surface to another. As money is processed through counting machines and automated teller machines (ATM), small amounts of drugs are readily transferred. An Agilent application note (Agilent Publication Number 5990-4254EN) details an application kit for the screening of 25 compounds considered in forensic and toxicology analyses using an Agilent 1200 Series LC system with an Agilent 6410 Triple Quadrupole LC/MS. In this work, an Agilent Poroshell 120 EC-C18 column is used to analyze 25 compounds found in the Agilent LC/MS Toxicology Test Mixture (Agilent p/n 5190-0470). This ammonium formate/acetonitrile gradient analysis is scaled using faster flow rates to shorten analysis time and exploit the low back pressure of this superficially porous column. Calibration curves for each of the 25 compounds are generated, and as a demonstration of the method a \$1 bill was extracted into methanol, analyzed and quantified.



Agilent Technologies

Introduction

The interest in superficially porous particles has led to discussions of method transfer from larger 5- μm totally porous particles, as well as from sub-2- μm totally porous particles. The high efficiency of superficially porous particles is similar to sub-2- μm totally porous particles. This is due to short mass transfer distance and substantially narrower particle size distribution.

The benefit of transferring from larger particle columns is very significant time savings, because the superficially porous particles are optimally run at faster flow rates (usually double) and are able to achieve similar resolution with a much shorter column length [1-2]. Because analysts will likely change column length and flow rate when transferring from larger totally porous particles to superficially porous columns, calculations must be performed to proportionally scale a gradient method and preserve the chromatographic selectivity (Equation 1).

Equation 1

$$t_2 = \frac{t_1 \cdot d_2^2 \cdot L_2 \cdot F_1}{d_1^2 \cdot L_1 \cdot F_2}$$

Where:

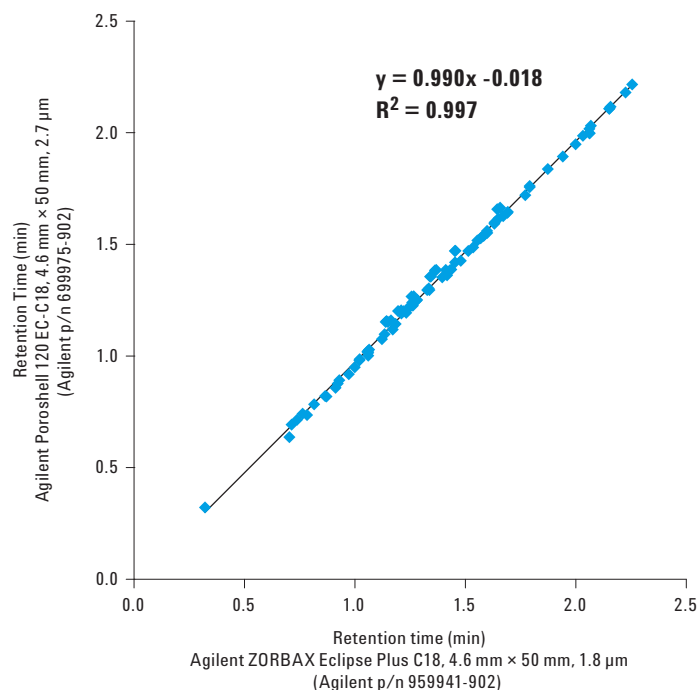
- t_1 and t_2 are the original and new gradient times (min)
- d_1 and d_2 are the original and new column internal diameters (mm)
- L_1 and L_2 are the original and new column lengths (mm)
- F_1 and F_2 are the original and new flow rates (mL/min)

In some cases, it may be useful to take advantage of the lower back pressure associated with superficially porous columns as compared to totally porous sub-2- μm columns. Depending upon operating conditions, the back pressure can be up to 50% less. This can give analysts the freedom to increase flow rates for higher throughput, or to increase column length to enhance resolution without exceeding the system pressure limits. Adjustments to flow rate and/or column length will require gradient scaling (Equation 1).

Method transfer can be especially easy, when columns like the superficially porous Agilent Poroshell 120 EC-C18 and totally porous Agilent ZORBAX Eclipse Plus C18 are manufactured to have similar bonding chemistries and use similar retention mechanisms. Figure 1 shows the similar retention of 90 compounds on Poroshell 120 EC-C18 and Eclipse Plus C18 columns using a generic gradient analysis with a variety of compounds from different chemical classifications. The high correlation coefficient (R^2) indicates a high degree of similarity between the interactions involved in the separation on the two C18 columns, while the slope ≈ 1 implies similar interaction strengths [3-4]. However, while many compounds give similar selectivity, it cannot be guaranteed that every application will transfer without adjustment.

This application note shows how a Poroshell 120 column can be used in a complex analysis, previously performed on a 1.8 μm column. This separation was demonstrated on Eclipse Plus in a previous Agilent application note (Publication Number 5990-4254EN) [5]. A 25-component LC/MS Toxicology Test Mixture (Agilent p/n 5190-0470) is used to illustrate the interchangeability between the two columns. Calibration curves for each of the 25 compounds on Poroshell 120 are constructed. A \$1 bill is extracted in methanol to show significant presence of cocaine, as well as noticeable quantities of oxycodone, methamphetamine, PCP and THC. Trace amounts of several more illicit and prescription drugs can be detected also. Drug trafficking is assumed to be the cause for their initial presence on US currency, while ATM's and counting machines are likely the cause of their widespread presence [6]. Additionally, this gradient analysis is transferred to a Poroshell 120 SB-C18 column, which shows some selectivity differences; however it can be run at higher temperatures to allow for even faster flow rates and analysis times. Agilent Poroshell 120 columns are available with two different C18 phases in order to change selectivity and still have a C18 column choice. Flow rates were increased to reach 400 and 600 bar to show performance achievable on both conventional HPLC's and newer UHPLC's.

Agilent Poroshell 120 EC-C18 has Very Similar Selectivity to Agilent ZORBAX Eclipse Plus C18



Mobile phase: A: 10 mM ammonium formate, pH 3
B: Acetonitrile

Gradient: 5% B at t_0 ramp to 95% B in 2 min, hold 95% B for 1 min

Flow rate: 2 mL/min

Sample: 1 μ L of 1 mg/mL standard in H₂O

Furazolidone	Biphenyl	Acetanilide	DL phenylalanine	Oxybutynin chloride 1
Chloramphenicol	Acenaphthene	Fenoprofen	Doxepin hydrochloride	Diphenhydramine
Pyrimethamine	Methoxy naphthalene	Catechol	Ephedrine hydrochloride	Diffunisal
Sulfaquinoxaline	Anisole	Phenol	Loperamide	Nisoldipine
Sulfamonomethoxine	Dimethoxy benzene	Resorcinol	Procaine hydrochloride	Diclofenac sodium salt
Nimopidol	Corticosterone	Hydroquinone	Fenoprofen calcium salt hydrate	Hydrocortisone
Sulfadimethoxine	Alpha hydroxyprogesterone	4 nitro phenol	Erythromycin	4 hydroxybenzoic acid
Sulfamethoxazole	Porgesterone	O cresol	Econazole nitrate	Procainamide hydrochloride
Sulfachloropyridazine	Alpha hydroxyprogesterone	P cresol	Gemfibrozil	Lidocaine
Sulfamethoxy-pyridazine	Prednisolone	3,4 dimethyl phenol	Beta estradiol	Terfenadine
Sulfamethizole	Mestranol	2,3 dimethyl phenol	Metoprolol	Terfenadine
Sulfamethazine	Deoxycorticosterone	2 nitro phenol	Prednisone	Chlortetracycline hydrochloride
Sulfamerazine	Progesterone	2,4 dimethyl phenol	Protriptyline	Chlorpheniramine maleate salt
Sulfathiazole	Chlorphenamine	2,5 dimethyl phenol	2-hydroxyhippuric acid	Chloramphenicol
Sulfadiazine	Berberine	1 naphthol	Hydroxyisophthalic acid	Buspirone hydrochloride
Benzaldehyde	Impramithue	Imipramine hydrochloride	Flufenamic acid	Benzocaine
Iodobenzene	Norethindrone	D methionine	Pramoxine hydrochloride	Antipyrine
Phenanthrene	Phenacetin	3,4 dihydroxy-L-phenyl alanine	Naproxen	Acetylsalicylic acid

Figure 1. Scatter plot of retention time of 90 compounds on Agilent Poroshell 120 EC-C18 versus Agilent Eclipse Plus C18.

Experimental

An Agilent 1200 Series Rapid Resolution LC (RRLC) system with an Agilent 6410 Triple Quadrupole LC/MS system was used for this work:

- G1312B Binary Pump SL with mobile phase A: 5 mM ammonium formate with 0.01% formic acid, and B: acetonitrile with 0.01% formic acid. Gradient was 10% B at t_0 , ramp to 15% B, ramp to 50% B, then ramp to 95% B and hold 95% B. Gradient times vary depending on column dimensions and flow rate (Table 1).
- G1367C Automatic Liquid Sampler (ALS) SL. Injection volume was 1.0 μ L.
- G1316B Thermostated Column Compartment (TCC) SL with temperature set to 60 °C or 90 °C (on Poroshell 120 SB-C18 only).
- G6410A Triple Quadrupole LC/MS: electrospray AP-ESI, drying gas temperature and flow: 350 °C, 12 L/min, nebulizer gas pressure: 30 psi, capillary voltage: 2000 V, in dMRM mode, transitions found in Table 2.
- MassHunter versions B.02.01, B.02.00 and B.03.01 were used for data acquisition, qualitative and quantitative analyses respectively.

Three Agilent columns were used in this work:

- Agilent Poroshell 120 EC-C18, 2.1 mm $\times$ 100 mm, 2.7 μ m (p/n 695775-902)
- Agilent Poroshell 120 SB-C18, 2.1 mm $\times$ 100 mm, 2.7 μ m (p/n 685775-902)
- Agilent ZORBAX RRHT Eclipse Plus C18, 2.1 mm $\times$ 100 mm, 1.8 μ m (p/n 959764-902)

The compounds of interest are shown in Table 2, with their respective retention times on Poroshell 120 EC-C18 at 0.5 mL/min, and their qualitative and quantitative MRM transitions. Sample is a 1 μ g/mL standard in methanol purchased from Agilent Technologies (LC/MS Toxicology Test Mixture, Agilent p/n 5190-0470). Serial dilutions in methanol were prepared for the calibration standards. The \$1 bill sample was extracted in 7 mL of methanol and ultrasonicated for 30 min. Additionally, acetonitrile, formic acid and ammonium formate were purchased from Sigma Aldrich (Bellefont, PA). Methanol was purchased from Honeywell, Burdick and Jackson (Muskegon, MI). Water used was 18 M- Ω Milli-Q water (Bedford, MA).

Table 1. HPLC Method Parameters for Various Columns and Conditions

Gradient and method parameters	2.1 $\times$ 100 mm 1.8- μ m Agilent ZORBAX Eclipse Plus C18	2.1 $\times$ 100 mm 2.7- μ m Agilent Poroshell 120 EC-C18	2.1 $\times$ 100 mm 2.7- μ m Agilent Poroshell 120 EC-C18	2.1 $\times$ 100 mm 2.7- μ m Agilent Poroshell 120 EC-C18	2.1 $\times$ 100 mm 2.7- μ m Agilent Poroshell 120 SB-C18	2.1 $\times$ 100 mm 2.7- μ m Agilent Poroshell 120 SB-C18	2.1 $\times$ 100 mm 2.7- μ m Agilent Poroshell 120 SB-C18
Flow rate (mL/min)	0.5	0.5	0.7	1.0	0.5	0.9	1.4
10% B (min)	0.00	0.00	0.00	0.00	0.00	0.00	0.00
15% B (min)	0.50	0.50	0.36	0.25	0.50	0.28	0.18
50% B (min)	3.00	3.00	2.14	1.50	3.00	1.67	1.07
95% B (min)	4.00	4.00	2.86	2.00	4.00	2.22	1.43
95% B (min)	6.00	6.00	4.29	3.00	6.00	3.33	2.14
Stop time (min)	6.00	6.00	4.29	3.00	6.00	3.33	2.14
Post run time (min)	2.00	2.00	1.43	1.00	2.00	1.11	0.71
Overall cycle time (min)	8.00	8.00	5.71	4.00	8.00	4.44	2.86
TCC temperature (°C)	60	60	60	60	90	90	90
Injection volume (μ L)	1.0	1.0	1.0	1.0	1.0	1.0	1.0
System pressure (bar)	375	280	385	550	195	370	595

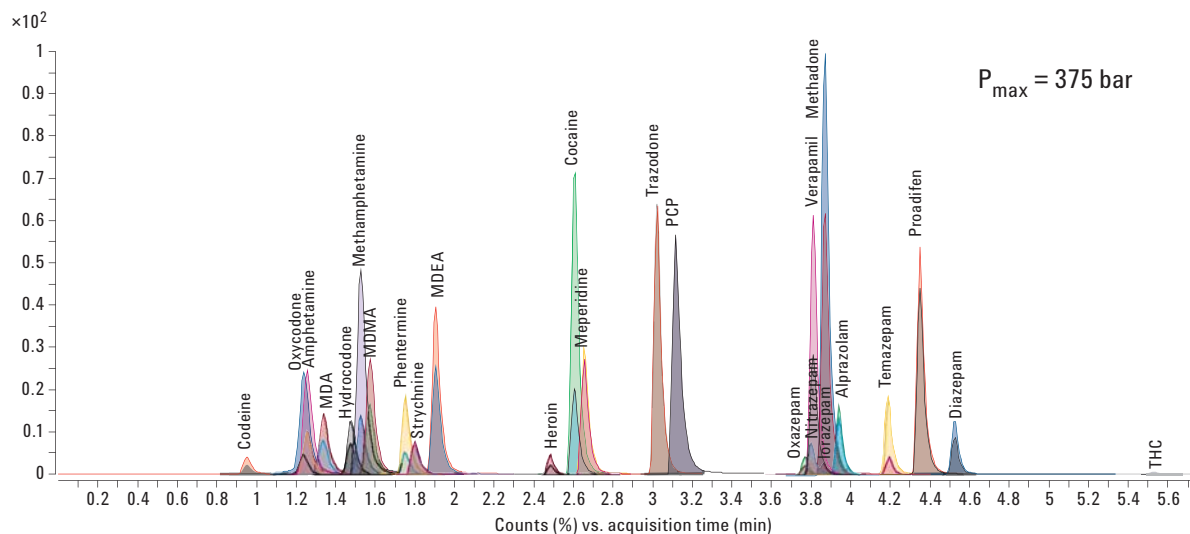
Table 2. MRM Transitions for 25 Compounds in Toxicology Test Mixture

Compound name	Precursor ion	Fragmentor voltage	Product ion 1	Collision energy 1	Product ion 2	Collision energy 2	Retention time (min)	Delta retention time
Codeine	300.2	158	165.1	45	58.1	29	0.89	0.4
Oxycodone	316.2	143	298.1	17	256.1	25	1.14	0.4
Amphetamine	136.1	66	119.1	5	91	17	1.19	0.4
MDA	180.1	61	163	5	105	21	1.25	0.4
Hydrocodone	300.2	159	199	29	128	65	1.34	0.4
Methamphetamine	150.1	92	119	5	91	17	1.43	0.4
MDMA	194.1	97	163	9	105	25	1.46	0.4
Strychnine	335.2	195	184	41	156	53	1.66	0.4
Phentermine	150	66	133	5	91	25	1.66	0.4
MDEA	208.1	107	163	9	105	25	1.8	0.4
Heroin	370.2	149	268.1	37	165	61	2.4	0.4
Cocaine	304.2	138	182.1	17	77	61	2.52	0.4
Meperidine	248.2	128	220.1	21	174.1	17	2.59	0.4
Trazodone	372.2	159	176	25	148	37	2.95	0.4
PCP	244.2	86	91	41	86.1	9	3.05	0.4
Oxazepam	287	150	269	12	241	20	3.66	0.4
Nitrazepam	282.1	148	236.1	25	180	41	3.66	0.4
Verapamil	455.3	158	165	37	150	45	3.75	0.4
Lorazepam	321	102	275	21	194	49	3.75	0.4
Methadone	310.2	112	265.1	9	105	29	3.83	0.4
Alprazolam	309.1	179	281	25	205	49	3.84	0.4
Temazepam	301.1	117	255.1	29	177	45	4.05	0.4
Proadifen	354.2	153	167	29	91.1	45	4.33	0.4
Diazepam	285.1	169	193	45	154	25	4.41	0.4
THC	315.2	150	193.2	20	123.3	30	5.4	0.4

Results and Discussion

Figure 2 shows the original method developed by P. Stone on an Agilent ZORBAX Eclipse Plus C18 2.1 mm × 100 mm, 1.8 µm column. This analysis is accomplished in 6 min with a 2-min post run time at 375 bar. Figure 3 shows the same method with an Agilent Poroshell 120 EC-C18 2.1 mm × 100 mm, 2.7 µm column. Analysis and post run time are identical to the Eclipse Plus method, while the system back pressure is reduced to 280 bar. While there are slight variations between elution patterns in Figures 2 and 3, overall selectivity is very similar, as would be predicted by Figure 1.

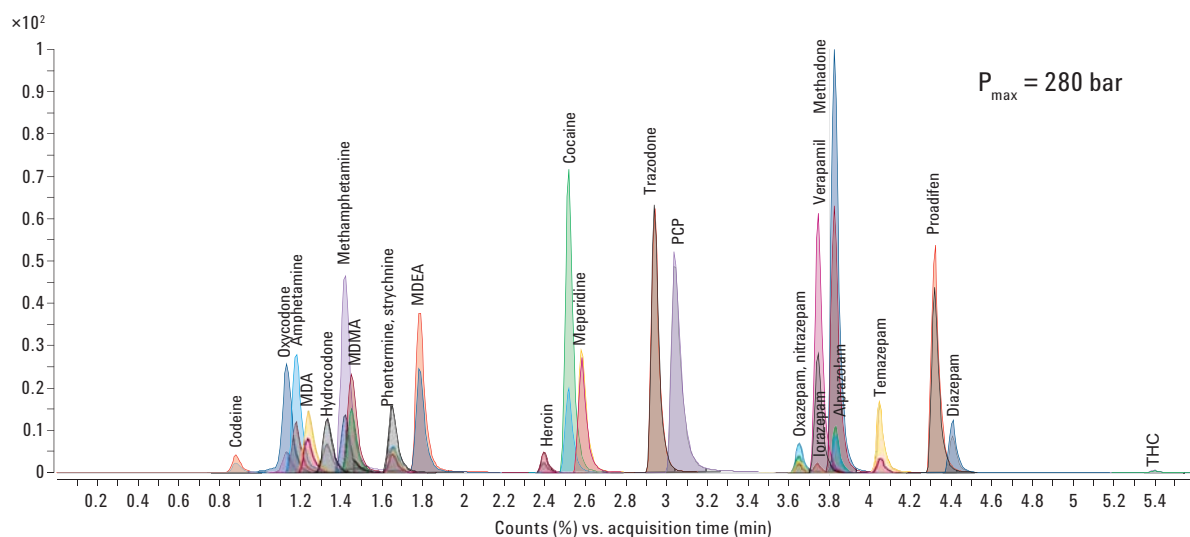
Original Toxicology Method on Agilent ZORBAX Eclipse Plus C18 2.1 mm × 100 mm, 1.8 µm (Agilent p/n 959764-902)



A: 5 mM ammonium formate w/ 0.01% formic acid (1 L water + 0.3153 g ammonium formate + 0.1 mL formic acid),
 B: acetonitrile w/ 0.01% formic acid (1 L acetonitrile + 0.1 mL formic acid); 0.5 mL/min; 10% B at t_0 , ramp to 15% B in 0.5 min,
 ramp to 50% B in 2.5 min, ramp to 95% B in 1 min, hold 95% B for 2 min; stop time 6 min, post run 2 min;
 Sample: injector program: draw 5 µL water, draw 1 µL LC/MS Toxicology Test Mixture (p/n 5190-0470), inject; TCC = 60 °C
 MS Source: electrospray AP-ESI, drying gas temperature and flow: 350 °C, 12 L/min, nebulizer gas pressure: 30 psi, capillary voltage: 2000V;
 MS Acquisition: dynamic MRM (see Table 2 for MRM transitions), positive ionization polarity

Figure 2. Agilent LC/MS Toxicology Test Mixture (Agilent p/n 5190-0470) analyzed on Agilent ZORBAX Eclipse Plus C18 via an Agilent 1200 Series LC system with detection by an Agilent 6410 Triple Quadrupole LC/MS.

Original Toxicology Method on Agilent Poroshell 120 EC-C18 2.1 mm × 100 mm, 2.7 µm (Agilent p/n 695775-902)



A: 5 mM ammonium formate w/ 0.01% formic acid (1 L water + 0.3153 g ammonium formate + 0.1 mL formic acid),
 B: acetonitrile w/ 0.01% formic acid (1 L acetonitrile + 0.1 mL formic acid); 0.5 mL/min; 10% B at t_0 , ramp to 15% B in 0.5 min,
 ramp to 50% B in 2.5 min, ramp to 95% B in 1 min, hold 95% B for 2 min; stop time 6 min, post run 2 min;
 Sample: injector program: draw 5 µL water, draw 1 µL LC/MS Toxicology Test Mixture (p/n 5190-0470), inject; TCC = 60 °C
 MS Source: electrospray AP-ESI, drying gas temperature and flow: 350 °C, 12 L/min, nebulizer gas pressure: 30 psi, capillary voltage: 2000V;
 MS Acquisition: dynamic MRM (see Table 2 for MRM transitions), positive ionization polarity

Figure 3. Agilent LC/MS Toxicology Test Mixture (Agilent p/n 5190-0470) analyzed on Agilent Poroshell 120 EC-C18 via an Agilent 1200 Series LC system with detection by an Agilent 6410 Triple Quadrupole LC/MS.

Table 3 shows calibration data for all 25 compounds found in the Agilent LC/MS Toxicology Test Mixture on Poroshell 120. All compounds exhibit strong linear correlations, with $R^2 > 0.9979$. Calibration data was used to quantify a methanol-extracted US \$1 bill sample; chromatographic and quantitative results are shown in Figure 4. A significant amount of cocaine

was found on the dollar bill. Oxycodone, methamphetamine, PCP and THC were also detected. Smaller quantities of amphetamine, hydrocodone, MDMA, heroin, methadone and diazepam were also found. Quantities of these substances on US currency are consistent with previous findings [6-8].

Table 3. Calibration Data for 25 Toxicology Compounds on Poroshell 120

Compound name	Linear calibration curve	Correlation coefficient, R^2
Codeine	$y = 25.4023x + 3.1628$	0.99990276
Oxycodone	$y = 138.9535x - 0.6269$	0.99938632
Amphetamine	$y = 196.3425x + 50.1606$	0.99987385
MDA	$y = 121.2945x + 180.2165$	0.99945701
Hydrocodone	$y = 72.1351x - 8.1010$	0.99964622
Methamphetamine	$y = 286.7936x + 429.4970$	0.99789141
MDMA	$y = 121.4217x - 55.0435$	0.99874569
Phentermine	$y = 110.8083x - 65.1028$	0.99914972
Strychnine	$y = 39.3465x - 9.5339$	0.99964358
MDEA	$y = 200.4804x - 14.2886$	0.99980092
Heroin	$y = 18.2969x + 0.4442$	0.99987634
Cocaine	$y = 295.8654x - 5.6261$	0.99963342
Meperidine	$y = 145.0367x + 17.2273$	0.99986118
Trazodone	$y = 286.1986x - 12.4408$	0.99969366
PCP	$y = 287.4395x - 24.8090$	0.99989199
Oxazepam	$y = 14.7883x - 0.4919$	0.99900677
Nitrazepam	$y = 49.1750x + 69.2747$	0.99876656
Verapamil	$y = 273.3001x + 17.3890$	0.99986678
Lorazepam	$y = 11.2911x + 6.0687$	0.99896851
Methadone	$y = 439.7238x - 6.7890$	0.9997511
Alprazolam	$y = 80.2721x + 18.5435$	0.99969734
Temazepam	$y = 70.9899x + 15.5246$	0.99976598
Proadifen	$y = 243.9474x - 13.0696$	0.99990655
Diazepam	$y = 68.9622x + 26.0608$	0.99948978
THC	$y = 3.1838x - 2.7072$	0.99801611

Oxycodone, Amphetamine, Hydrocodone, Methamphetamine, MDMA, Heroin, Cocaine, PCP, Methadone, Diazepam and THC are Extracted from a US \$1 Bill and Quantified

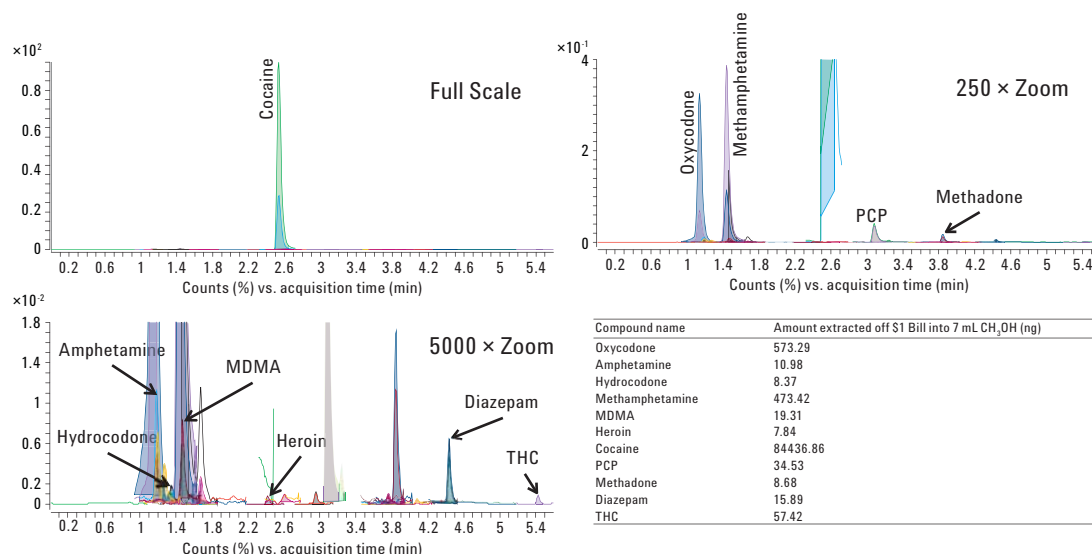


Figure 4. Chromatographic and quantitative results from a random US \$1 bill sample extracted with 7 mL of methanol and ultrasonicated for 30 minutes.

Due to the low system back pressure generated with the Poroshell 120 column, the flow rate can be increased from 0.5 mL/min to 0.7 mL/min without exceeding 400 bar for use on a standard HPLC, or it can be increased to 1 mL/min without exceeding 600 bar for use on a UHPLC, as shown in Figure 5. The increased flow rate may be desirable when high throughput is important and when a UHPLC is available for use. Overall cycle time can be decreased by 2.3 minutes while keeping pressure below 400 bar, or by 4 minutes while keeping pressure below 600 bar (a 50% reduction in cycle time). Increasing the flow rate to this degree does cause some loss in resolution, but with MS detection this is not critical.

Significant Time Savings are Possible by Increasing Flow Rate with Agilent Poroshell 120 EC-C18 to LC System Pressure Limits, whether 400 or 600 bar

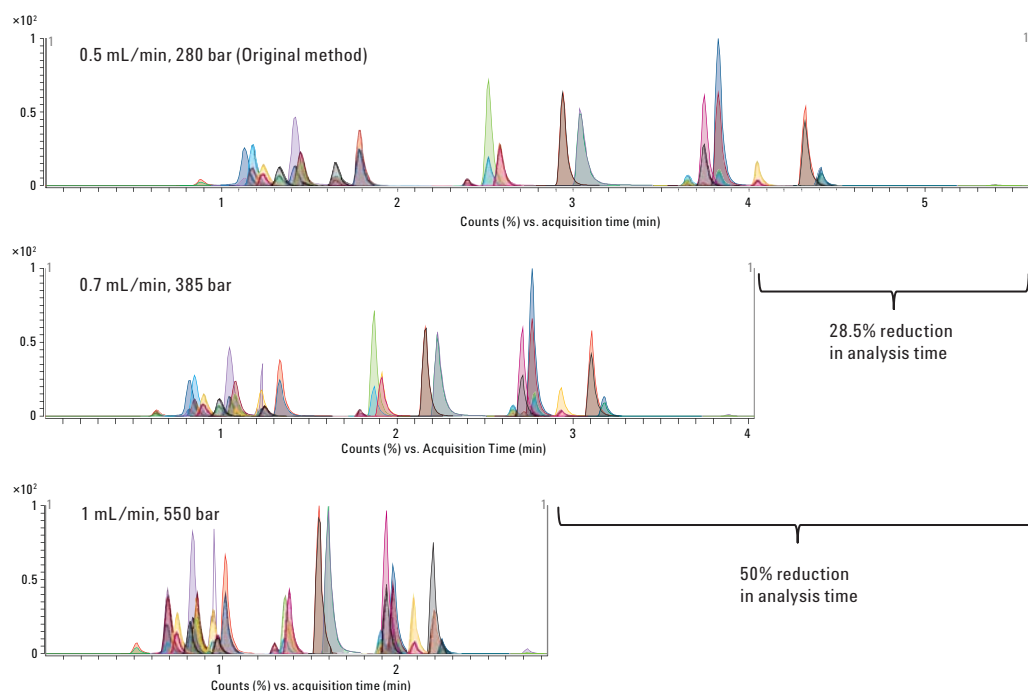


Figure 5. Overlay of Agilent Poroshell 120 EC-C18 toxicology analysis showing time savings by increasing flow rate to reach a 400 or 600 bar system limit.

Flow rate can be further increased by elevating temperature, thereby reducing mobile phase viscosity. The original method however was run at 60 °C, which is the maximum operating temperature for both Eclipse Plus C18 and Poroshell 120 EC-C18. In order to perform this analysis at a higher temperature, the column must be replaced with a Poroshell 120 SB-C18, which has a maximum operating temperature of 90 °C. Figure 6 shows the fast chromatography possible with Poroshell 120 SB-C18. With a 600 bar system pressure limit, it is possible to reduce run time by 64.3%, however this comes

at the cost of reduced resolution. For an analysis as complex as this toxicology method, this loss of resolution and significant coelution will cost the analysts a reduction in data points across all peaks, therefore reducing the quality of the results. A simple solution may be to increase column length. A slight increase in column length from 100 mm to 150 mm will increase the resolution of all compounds. While the longer column cannot be run at quite as fast flow rates the analyst can still glean significant time savings by running it at its respective highest flow rate without exceeding system limitations.

Very Significant Time Savings are Possible by Increasing Temperature and Flow Rate with Agilent Poroshell 120 SB-C18 to LC System Pressure Limits, whether 400 or 600 bar

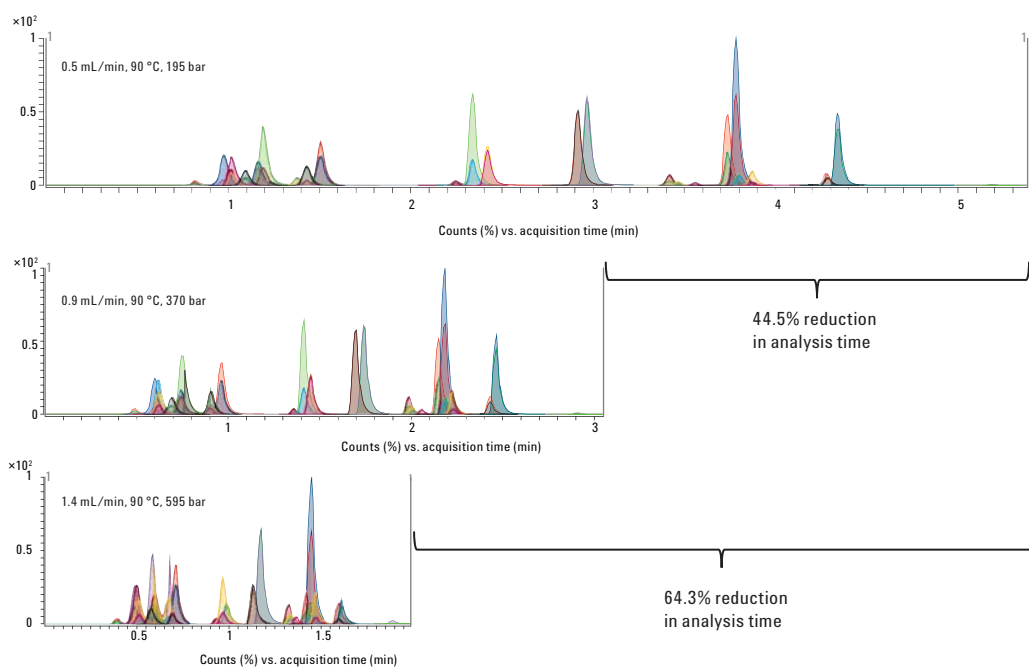


Figure 6. Overlay of Agilent Poroshell 120 SB-C18 toxicology analysis showing time savings by increasing temperature and flow rate to reach a 400 or 600 bar system limit.

Conclusion

A complex analysis of 25 toxicology compounds, that was originally performed on an Agilent ZORBAX Eclipse Plus C18 column, was easily carried out on a superficially porous Agilent Poroshell 120 EC-C18 column with high quality results and substantial time savings. Other complex analyses can likely be transferred from 1.8- μ m Eclipse Plus C18 to Poroshell 120 EC-C18 of the same dimensions without method modification, due to very similar selectivity and efficiency. The lower back pressure of Poroshell 120's 2.7- μ m particles can be exploited for productivity gains; faster flow rates may be used to shorten analysis time without exceeding system pressure limits for 400 bar HPLC's or higher pressure UHPLC's. This method was used to detect and quantify several drugs of abuse found on a \$1 bill, including: cocaine, oxycodone, methamphetamine, PCP and THC.

References

1. A. Gratzfeld-Hüsgen, E. Naegele, "Maximizing efficiency using Agilent Poroshell 120 columns," Agilent Technologies publication 5990-5602EN, 2010.
2. V. Meyer, "Practical High Performance Liquid Chromatography," Fourth Ed., p 34, Wiley, 2004.
3. K. Croes, A. Steffens, D. Marchand, L. Snyder, "Relevance of π - π and dipole-dipole interactions for retention on cyano and phenyl columns in reversed-phase liquid chromatography", *Journal of Chromatography A*, Volume 1098, Issues 1-2, 9 December 2005, Pages 123-130.
4. W. Long, A. Mack, "Comparison of Selectivity Differences Among Different Agilent ZORBAX Phenyl Columns using Acetonitrile or Methanol," Agilent Technologies publication 5990-4711EN, 2009.
5. P. Stone, "An Application Kit for the Screening of Samples for Analytes of Forensic and Toxicological Interest using LC/QQQ MS/MS with a Dynamic MRM Transition Database," Agilent Technologies publication 5990-4254EN, 2009.
6. J. Oyler, W. Darwin, E. Cone, "Cocaine Contamination of United States Paper Currency," *Journal of Analytical Toxicology*, Volume 20, Number 4, July 1996, Pages 213-216.
7. A. Jenkins, "Drug contamination of US paper currency, Forensic Science International," Volume 121, Issue 3, 1 October 2001, Pages 189-193.
8. Y. Zuo, K. Zhang, J. Wu, C. Rego, J. Fritz, "An accurate and nondestructive GC method for determination of cocaine on US paper currency," *Journal of Separation Science*, Volume 31, Issue 13, 21 July 2008, Pages 2444-2450.

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
September 2, 2010
5990-6345EN



Agilent Technologies