

Paul Altiero Applications Chemist, Agilent September 2019

Introduction: What and Why System
Configuration:
Measure
Delay Volume

Development:
Scouting
Gradient to
Analytical
Method

Transferring
Gradients to
Different
Columns

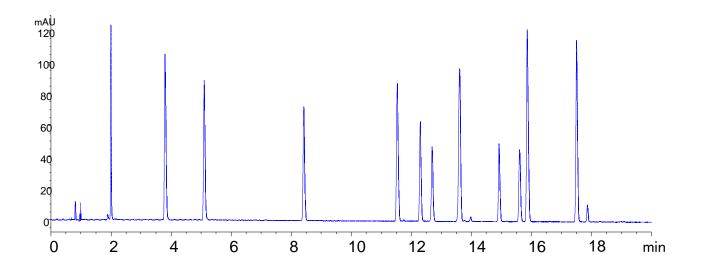
Introduction: What and Why System
Configuration:
Measure
Delay Volume

Development:
Scouting
Gradient to
Analytical
Method

Transferring
Gradients to
Different
Columns

What is Gradient Elution?

A separation that occurs by continuously increasing the solvent strength of the mobile phase



- Hydroquinone
- Resourcinol
- 3. Catechol
- 4. Phenol
- 5. 4-Nitrophenol
- 6. p-cresol
- 7. o-cresol
- 8. 2-Nitrophenol
- 9. 3,4 di methyl phenol
- 10. 2,3 di methyl phenol
- 11. 2,5 di methyl phenol
- 12. 1-napthol

Column: Poroshell 120 EC-C18, 4.6 x 150 mm, 2.7 μm

Mobile Phase: Solvent A: water with 0.1% formic acid, Solvent B: acetonitrile

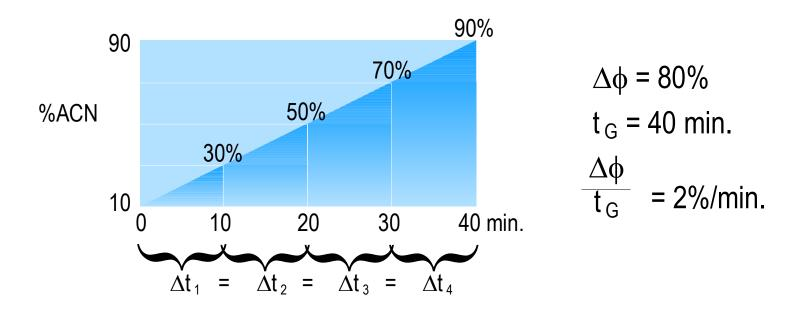
Gradient: 0–3 min 5% B, 5%-60% B over 22 minutes

Agilent 1200 SL controlled temperature at 25 °C, 2 µl flow cell



Gradient Elution for Reversed-Phase HPLC

Increasing the solvent strength = Increasing the % organic in the mobile phase Linear solvent strength gradient = % per min is a constant



For every 20% change in ACN, $\triangle t$ is 10 min.

Gradient Elution Analysis of a Complex Sample:

EU banned Azo Colorants in textiles



- 1. 1,4-phenylendiamine
- 2. 4-methoxy-m-phenylenediamine 15. o-dianisidine
- 3. 4-methyl-m-phenylenediamine
- 4. Aniline
- 5. benzidine
- 6. o-anisidine
- 7. 4,4'-oxydianiline
- 8. o-toluidine
- 9. 4-chloroaniline
- 10. 5-nitro-o-toluidine
- mAŪ 11. 4,4'-methylenedianiline
 - 12. p-cresidine
 - 13. 2,6-dimethylaniline

2

- 14. 2,4-dimethylaniline
- 16. 4,4'-bi-o-toluidine
- 17. 4,4'-thiodianiline
- 18. 2-naphthylamine
- 19. 4-chloro-o-toluidine
- 20. 2,4,5-trimethylaniline
- 21. 4,4'-methylenedi-o-toluidine
- 22. biphenyl-4-ylamine
- 23. 3.3'-dichlorobenzidine
- 24. 4-aminoazobenzene
- 25. 2,2'-dichloro-4.4'-methylene-dianiline
- 26. o-aminoazotoluene

10

6

4



3.0 x 150 mm, 2.7 µm (p/n: 693975-302)

Column temperature: 40 °C

Flow rate: 0.8 mL/min

Mobile phase:

A: 0.575 g Mono ammonium phosphate, 0.7 g Disodium hydrogen phosphate in 900 mL water added with 100 mL

methanol and adjusted using phosphoric

acid to pH 6.9

B: methanol

see table

26	Gradient:
	Num-
12	min

24

22

10

Time (min)	B%	
0	14	
4.5	29	
8	29	
9	50	
10.5	65	
12	90	
14	100	
15	14	

-10

8

Three Major Reasons to Choose Gradient Elution

- 1. Faster separation of samples having components that vary in polarity.
- 2. To separate mixtures having a large number of components
- 3. To separate high molecular weight mixtures (i.e., large molecules, peptides and proteins)



Gradient Separation Reduces Analysis Time and Improves Resolution – Even with the Shortest Columns

Columns: Poroshell 120 EC-C18, 3.0 x 30 mm, 2.7 µm

Mobile Phase: A: 10 mM ammonium acetate, pH= 6.8, B: Acetonitrile. Temperature: 30 °C

Mobile Phase: A: 20 mM Phosphate buffer, B: Acetonitrile. Temperature: 30°C. Detection: UV 245 nm.

Sample: Acetaminophen impurities.

Isocratic elution:

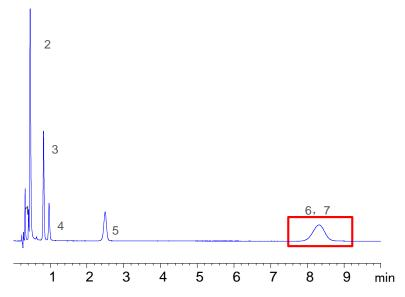
90% agueous/10% ACN

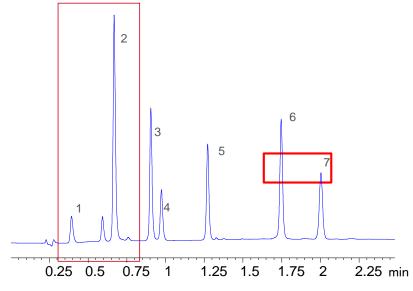
Flow Rate: 0.65 mL/min



5-50% ACN in 2 min.

Flow rate: 0.65 mL/min





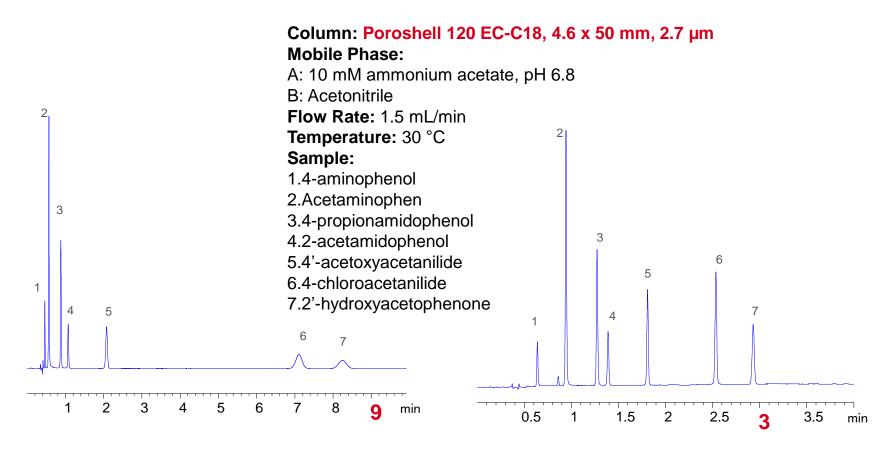
- The gradient analysis is much faster using the same 30 mm Poroshell 120 column.
- And resolution is dramatically improved for all peak pairs!

Gradient Separation is Faster than Isocratic

Separation of acetaminophen impurities on Poroshell 120 EC-C18

Isocratic elution: 85% aqueous/15% ACN

Gradient elution: 5–50% ACN in 5 min



Introduction: What and Why

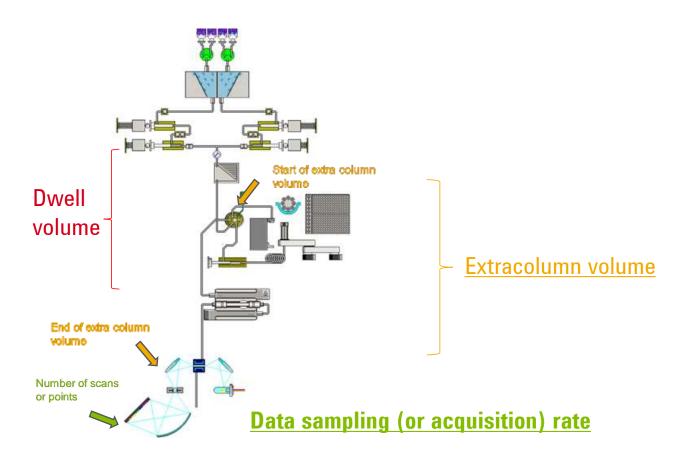
System
Configuration:
Measure
Delay Volume

Development:
Scouting
Gradient to
Analytical
Method

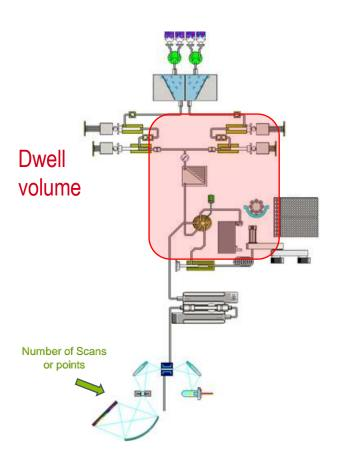
Transferring
Gradients to
Different
Columns

Gradient Separations

Instrument Impact on Column Performance



Instrument Impact on Column Performance Dwell volume



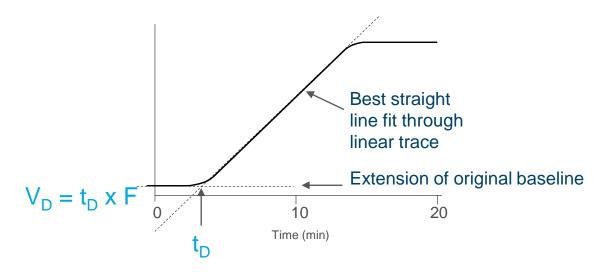
 Dwell volume = volume from formation of gradient to the column

Determining the Dwell Volume of Your System

- Look it up in the LC manual or follow the procedure below
- Replace column with short piece of HPLC stainless steel tubing
- Prepare mobile phase components
 A. Water UV-transparent
 B. Water with 0.2% acetone UV-absorbing
- Monitor at 265 nm
- Run gradient profile 0–100% B/10 min at 1.0 mL/min
- Record
- Expected dwell volume in UHPLCs μL range!



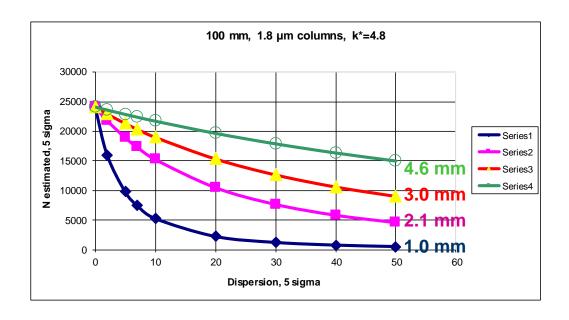
Measuring Dwell Volume (V_D)



- Intersection of the two lines identifies dwell time (t_D)
- Dwell volume is equal to product of the flow rate and the dwell time.



Efficiency Yield vs. Dispersion in LC Systems



Use the largest columns suitable for the application requirements -- they are less affected by extracolumn contributions. At the same time, consider solvent consumption and detector (ELSD, MS, etc.) compatibility.

Introduction: What and Why

System
Configuration:
Measure
Delay Volume

Development:
Scouting
Gradient to
Analytical
Method

Transferring
Gradients to
Different
Columns

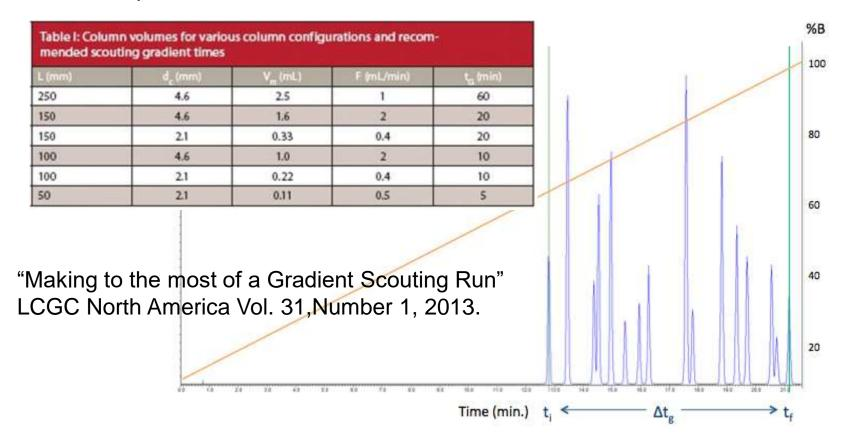
Starting Point Scouting Gradient

A good starting point for work is a scouting gradient.

The conditions recommended by John Dolan are 5–95% acetonitrile, low pH, and are dependent on the column length.

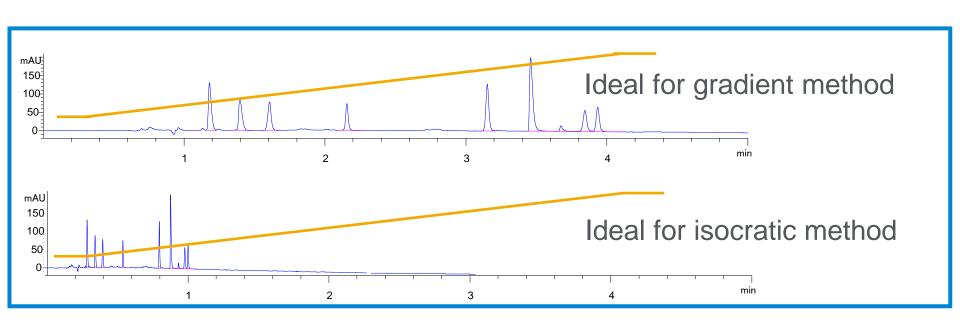
Where 10 cm columns are chosen, use a 10 minute gradient.

This example shows a 150 mm column.



Gradients are Critical Tools for Faster Methods

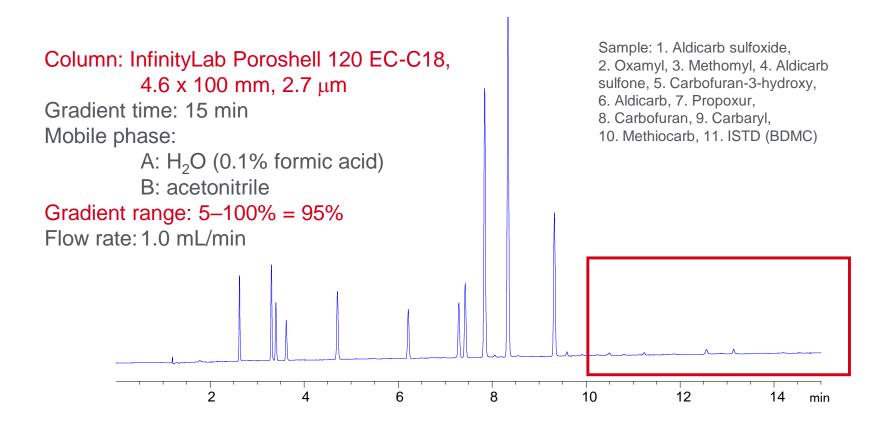
- Run a scouting method 5% to 95% organic (reversed phase)
- · Quick evaluation: how much of the gradient is occupied
 - $-\frac{\Delta t_G}{t_G} \le 25\%$ isocratic is recommended
 - $-\frac{\Delta t_G}{t_G} \ge 40\%$ gradient is recommended





Step 1: Choose, Shorter Efficient Column

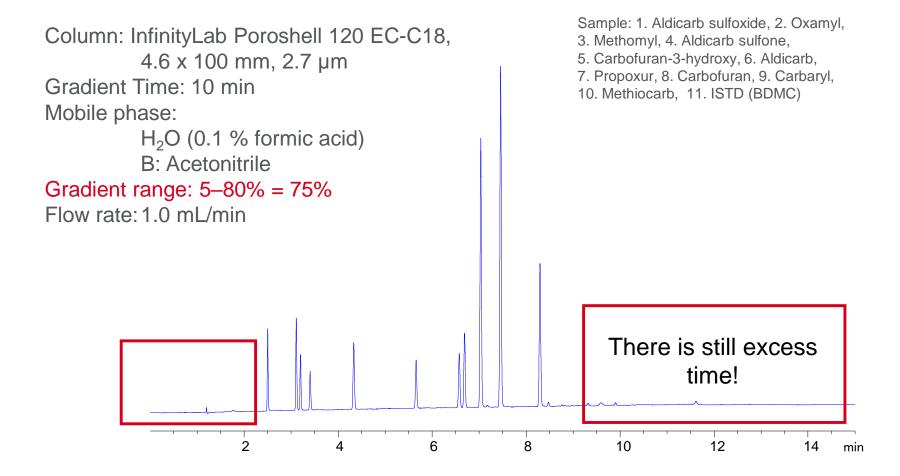
Perform Gradient Scouting from 5%B-100% in 15 min



The scouting shows that there is wasted time in this chromatogram and resolution of all components can be achieved. Optimization possible!

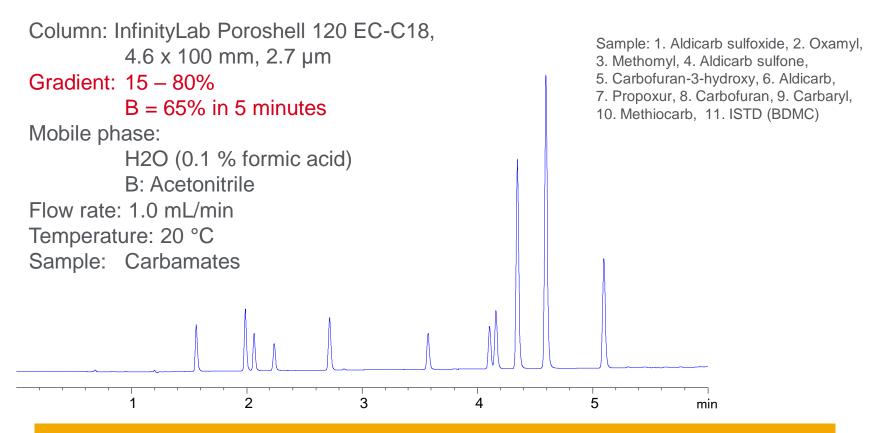


Step 2: Reduce Gradient Range to Minimize Time Adjust gradient from 5%B-80% in 10 min



Step 3: Finalize Your Results

Increase starting % organic and reduce time

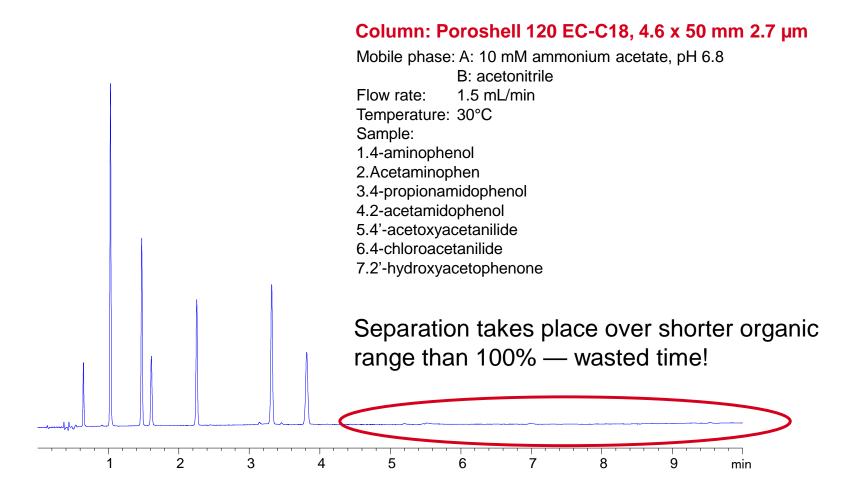


Saved 50% of the time with method optimization. Used Poroshell 120 for high efficiency and resolution.



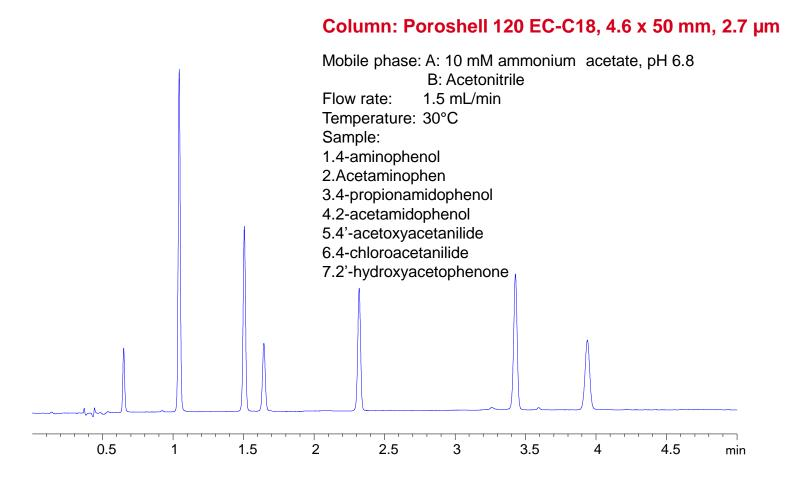
Gradient Scouting Works for Any Sample

Gradient from 5%-100% in 10 min for acetaminophen





Optimizing Gradient from 5%–50% in 5 min to Reduce Wasted Time

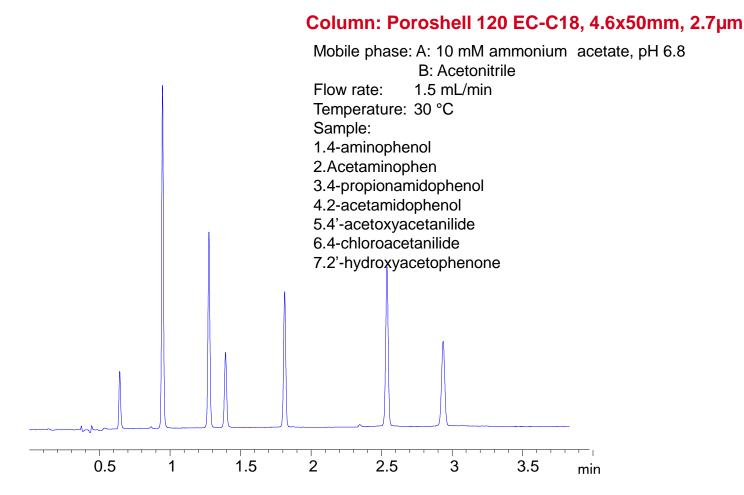


Excellent resolution and distribution of peaks in the gradient – within 5 minutes.



Final Optimization to Reduce Time

Gradient from 5%-50% in 3 min



Introduction: What and Why

System
Configuration:
Measure
Delay Volume

Development:
Scouting
Gradient to
Analytical
Method

Transferring
Gradients to
Different
Columns

Resolution Relationship for Gradient Elution

$$R \approx \frac{\sqrt{N}}{4} \alpha^* k^*$$

k* - represents the fact that k changes constantly during a gradient

$$\textbf{K*} = \begin{array}{c} 87 \ t_g \, \textbf{F} \\ \hline S \ (\Delta \% B) \ V_m \end{array} \begin{array}{c} \Delta \% B = \\ S = \\ F = \\ V_m = \end{array} \begin{array}{c} \text{difference between initial and final \% B values} \\ \text{constant (\approx 4 for 100 - 500 Da)} \\ \text{flow rate (mL/min.)} \\ \text{gradient time (min.)} \\ \text{column void volume (mL)} \\ \end{array}$$



Maintaining k*

To keep relative peak position unchanged while changing analysis parameters

Any decrease in

Column length

Can be offset by a proportional

Decrease in t_G or F

Increase in ∆%B

Column volume (i.d.)

Decrease in t_G or F

Increase in ∆%B

∆%B (same column)

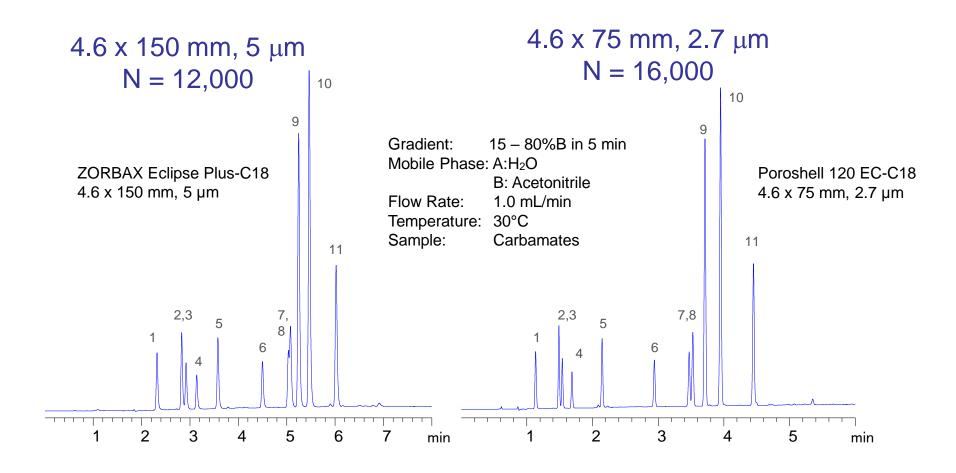


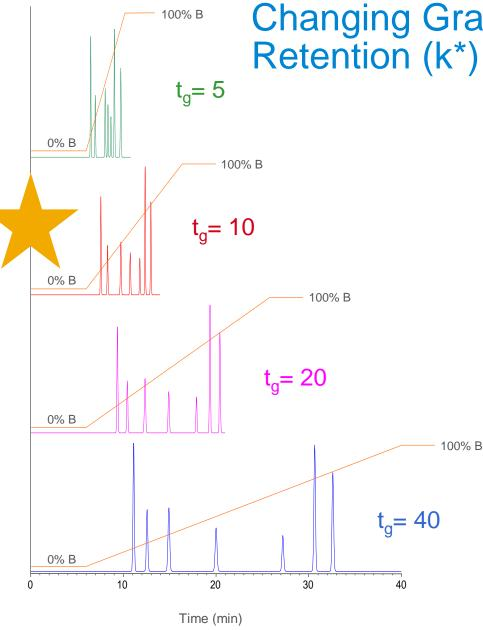
Decrease in t_G or F

$$k^* \propto \frac{t_G \bullet F}{S \bullet \Delta \Phi \bullet Vm}$$

A Shorter Column (smaller Vm)

Reduce run-time and improve resolution while maintaining constant N





Changing Gradient Time to Affect Retention (k*) and Resolution

$$k^* = \frac{t_g F}{S \Delta \% B V_m}$$

1/k* = gradient steepness = b

 $\Delta\Phi$ = change in volume fraction of B solvent

S = constant

F = flow rate (mL/min.)

 t_{o} = gradient time (min.)

 $V_m = \text{column void volume (mL)}$

- S ≈ 4–5 for small molecules
- 10 < S < 1000 for peptides and proteins

Adapting Gradient Methods to Different Column Dimensions

To adjust gradient methods to different column dimensions keep gradient steepness (b) the same.

gradient
$$S \bullet \Delta \Phi \bullet Vm$$
1/k* ∞ Steepness = $b = \frac{t_G \bullet F}{}$

S = constant

 $\Delta\Phi$ = change in % organic during the gradient run

Vm = void volume of column

F = flow rate

 t_G = gradient time

k* = k of solute at mid point of column

If "b" is kept constant from run-to-run peaks will elute in the same relative pattern.



Adjusting a Gradient from a 4.6 x 150 mm Column to a 2.1 x 100 mm Column

4.6 x 150 mm

$$\Delta \Phi = 40 (20\% - 60\%)$$

Vm = 1.5 mL

$$F = 1.0 \text{ mL/min}$$

$$t_G = 15 \text{ min}$$

$$b = \frac{\Delta \Phi_1 \cdot V_{m_1}}{F_1 \cdot t_{G_1}} = \frac{\Delta \Phi_2 \cdot V_{m_2}}{F_2 \cdot t_{G_2}}$$

2.1 x 100 mm

$$\Delta \Phi = 40 (20\% - 60\%)$$

$$Vm = 0.2 mL$$

$$F = 0.2 \text{ mL/min}$$

$$t_G = ? (10 min)$$

$$t_{G_2} = t_{G_1} \cdot \frac{\Delta \Phi_2}{\Delta \Phi_1} \cdot \frac{V_{m_2}}{V_{m_1}} \cdot \frac{F_1}{F_2}$$

$$t_{G_2} = 15 \cdot \frac{40}{40} \cdot \frac{0.2}{1.5} \cdot \frac{1}{1.5} = 10$$



Adjusting a Gradient from a 4.6 x 150 mm Column to a 2.1 x 100 mm Column for constant %B

$$t_{G_2} = t_{G_1} \cdot \frac{F_1}{F_2} \cdot \left(\frac{d_2}{d_1}\right)^2 \frac{L_2}{L_1} \cdot \frac{\Delta \Phi_2}{\Delta \Phi_1}$$

for constant %B

$$t_{G_2} = t_{G_1} \cdot \frac{F_1}{F_2} \cdot \left(\frac{d_2}{d_1}\right)^2 \frac{L_2}{L_1}$$

<u>4.6 x 150 m</u>

F = 1.0 mL

2.1 x 100 mm

 $F = 0.2 \, mL$

%A	%B	t_{old}	t_{new}
95	5	0	0
95	5	1	0.7
35	65	13	9.0
35	65	14	9.7
0	100	14.5	10.0
0	100	16.5	11.5
95	5	17	11.8
95	5	20	13.9

Adjusting a Gradient from a 4.6 x 150 mm Column to a 2.1 x 100 mm Column for constant %B

$$t_{G_2} = t_{G_1} \cdot \frac{F_1}{F_2} \cdot \left(\frac{d_2}{d_1}\right)^2 \frac{L_2}{L_1} \cdot \frac{\Delta \Phi_2}{\Delta \Phi_1}$$

for constant %B

$$t_{G_2} = t_{G_1} \cdot \frac{F_1}{F_2} \cdot \left(\frac{d_2}{d_1}\right)^2 \frac{L_2}{L_1}$$

<u>4.6 x 150 m</u>

F = 1.0 mL

2.1 x 100 mm

 $F = 0.2 \, mL$

%A	%B	t_{old}	t_{new}
95	5	0	0
95	5	1	0.7
35	65	13	9.0
35	65	14	9.7
0	100	14.5	10.0
0	100	16.5	11.5
95	5	17	11.8
95	5	20	13.9

Contact Agilent Chemistries and Supplies Technical Support



1-800-227-9770 Option 3, Option 3:

Option 1 for GC/GCMS Columns and Supplies

Option 2 for LC/LCMS Columns and Supplies

Option 3 for Sample Preparation, Filtration and QuEChERS

Option 4 for Spectroscopy Supplies

Available in the USA & Canada 8-5 all time zones



gc-column-support@Agilent.com

<u>lc-column-support@agilent.com</u>

spp-support@agilent.com

spectro-supplies-support@agilent.com