

The Agilent InfinityLab 2D-LC Solution with Active Solvent Modulation

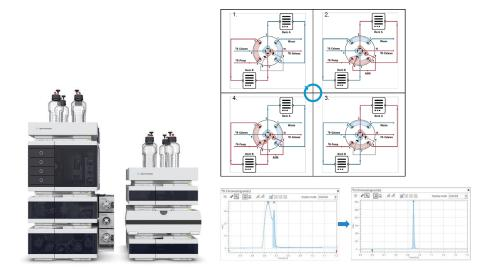
Achieving Improved Resolution and Sensitivity for Challenging Combinations of Separation Conditions

Author

Sonja Krieger Agilent Technologies, Inc.

Abstract

A major challenge during method development in two-dimensional liquid chromatography (2D-LC) can be the incompatibility between the first (¹D) and second dimension (²D) solvent systems. For challenging combinations of separation systems, volume overload effects such as breakthrough or broad and distorted peaks in the ²D can be observed. Active solvent modulation (ASM) provides a valve-based approach to address this problem, leading to improved resolution and sensitivity in the ²D. This Technical Overview explains the principle of ASM, and demonstrates its effect in the combination of two reversed-phase (RP) LC separations and a combination of hydrophilic interaction liquid chromatography (HILIC) and RPLC.



Introduction

In situations where the sample to be analyzed is complex, or contains compounds of interest that are difficult to separate, the separation that can be achieved using one-dimensional liquid chromatography (1D-LC) might not be sufficient. Two-dimensional liquid chromatography (2D-LC) is then the method of choice¹. The Agilent InfinityLab 2D-LC Solution enables comprehensive (LCxLC), multiple heart-cutting (MHC), and high-resolution sampling (HiRes) 2D-LC to be performed. Therefore, enabling either improved separation for the entire sample (LCxLC) or separation of target compounds (MHC and HiRes 2D-LC), that could not be separated using 1D-LC.

One challenge in 2D-LC method development can be the incompatibility between the first (1D) and second dimension (2D) solvent systems2. In loop-based 2D-LC, the analyte is transferred from the ¹D to the ²D in the ¹D effluent. When the volume of ¹D effluent that is transferred to the ²D becomes significant relative to the ²D column dead volume, the ¹D effluent essentially becomes the ²D mobile phase for a certain amount of time¹. For combinations of separation systems where the analytes show low retention on the ²D column in the ¹D effluent, this can lead to volume overload effects. such as breakthrough or broad and distorted peaks in the ²D^{1,2}. Active solvent modulation (ASM) addresses this problem by diluting the ¹D effluent with ²D mobile phase ahead of the ²D column. enabling focusing of the analytes at the ²D column head².

This Technical Overview explains the principle of active solvent modulation, and demonstrates its effect in the combination of two reversed-phase (RP) LC separations as well as a combination of hydrophilic interaction liquid chromatography (HILIC) and RPLC.

Experimental

Equipment

The Agilent 1290 Infinity II 2D-LC System comprised the following modules:

- Two Agilent 1290 Infinity II High Speed Pumps (G7120A)
- Agilent 1290 Infinity II Multisampler (G7167B) with cooler (option #100)
- Two Agilent 1290 Infinity II Multicolumn Thermostats (G7116B)
- Two Agilent 1290 Infinity II Diode Array Detectors (G7117B) with a 10-mm Max-Light Cartridge Cell (G4212-60008)
- Agilent 1290 Infinity Valve Drive (G1170A) with 2D-LC Valve, Active Solvent Modulation (G4243A)
- Two Agilent 1290 Infinity Valve Drives (G1170A) with Multiple heart-cutting Valves (G4242-64000) equipped with 40-µL loops

Mass selective detection (MSD) was performed using an Agilent 6150 Single Quadrupole LC/MS (G6150BA) equipped with an Agilent Jet Stream ESI source (G1958-65538).

Software

Agilent OpenLAB CDS ChemStation Edition Rev. C.01.07 SR3 [465] with 2D-LC Software version A.01.04.

Columns

- Agilent ZORBAX Eclipse Plus PAH, 2.1 × 100 mm, 1.8 μm (p/n 959764-918)
- Agilent ZORBAX Eclipse Plus C8 RRHD, 2.1 × 50 mm, 1.8 μm (p/n 959757-906)
- Agilent ZORBAX HILIC Plus, 2.1 × 100 mm, 3.5 μm (p/n 959793-901)
- Agilent AdvanceBio Peptide Mapping, 2.1 × 100 mm, 2.7 μm (p/n 655750-902)

Chemicals

All solvents were LC grade. Acetonitrile and methanol were purchased from Merck (Darmstadt, Germany). Fresh ultrapure water was obtained from a Milli-Q Integral system equipped with a 0.22-µm membrane point-of-use cartridge (Millipak, EMD Millipore, Billerica, MA, USA). Ammonium formate and formic acid were purchased from Sigma-Aldrich (Steinheim, Germany) and VWR (Darmstadt, Germany), respectively.

Samples and Methods RPLC-RPLC Analysis of Polycyclic Aromatic Hydrocarbons (PAHs)

A mixture of 16 polycyclic aromatic hydrocarbons (PAHs) (PAH mix 25 from Dr. Ehrenstorfer, Augsburg, Germany (YA20950025AB)) was diluted with acetonitrile to obtain a concentration of 10 μ g/mL of each compound. Table 1 shows the instrument parameters.

HILIC-RPLC Analysis of Peptides

A peptide standard (Agilent 10 Peptide Standard, lyophilized 71 μg , p/n 5190-0583) was dissolved in 100 μL water/acetonitrile (50/50, v/v) containing 0.1 % formic acid. Table 2 shows the instrument parameters.

The following peptides are contained in the peptide standard:

- Bradykinin frag 1-7 (Peptide 1)
- Bradykinin (Peptide 2)
- Angiotensin II (Peptide 3)
- Neurotensin (Peptide 4)
- Angiotensin I (Peptide 5)
- Renin substrate porcine (Peptide 6)
- [Ace-F-3,-2 H-1] Angiotensinogen (1-14) (Peptide 7)
- Ser/Thr Protein Phosphatase (15-31) (Peptide 8)
- [F14] Ser/Thr Protein Phosphatase (15-31) (Peptide 9)
- Melittin (Peptide 10)

 $\textbf{Table 1.} \ \mathsf{RPLC}\text{-}\mathsf{RPLC} \ \mathsf{parameters}.$

First dimension						
Column	Agilent ZORBAX Eclipse Plus PAH, 2.1 × 100 mm, 1.8 μm					
Solvent	A) Water B) Methanol					
Gradient	0 minutes – 50 %B 15 minutes – 100 %B Stop time: 22 minutes Post time: 5 minutes Stop time and post time were set in the ² D pump to allow extension of the run time until completion of ² D analysis of all cuts.					
Flow rate	0.200 mL/min					
Temperature	40 °C					
	DAD; 230 nm/4 nm, reference 390 nm/20 nm; 20 Hz					
Detection	To protect the flow cell from pressure pulses originating from valve switches, a pressure release kit (G4236-60010) was installed between the ¹D DAD and the 2D-LC valve.					
Injection	Injection volume: 2 µL Sample temperature: 8 °C Needle wash: 3 seconds in water/acetonitrile (70/30)					
Second dimension						
Column	Agilent ZORBAX Eclipse Plus C8, 2.1 × 50 mm, 1.8 μm					
Solvent	A) Water B) Acetonitrile					
Temperature	40 °C					
Detection	DAD; 230 nm/4 nm, reference 390 nm/20 nm; 80 Hz					
	2D-LC					
2D-LC Mode	Heart-cutting					
Flow rate	1.000 mL/min					
Sampling table	Sampling table Time based heart-cutting with a sampling time of 0.20 minutes was set up using the ¹D chromatogram of the PAH mixture as reference chromatogram.					
	ASM and ² D gradient – Method without the use of ASM					
ASM	ASM disabled					
² D Gradient	0.00 minutes - 50 %B 1.00 minutes - 100 %B ² D gradient stop time: 1.20 minutes ² D cycle time: 1.70 minutes					
Α	SM and ² D gradient – Method with the use of ASM using an ASM factor of 1.5					
ASM	ASM capillary: 5500-1303 (0.12 × 680 mm) ASM enabled (ASM factor: 1.5) Flush sample loop 3.0 times (0.18 minutes)					
² D Gradient	0.00 minutes – 5 %B 0.18 minutes – 5 %B 0.19 minutes – 50 %B 1.19 minutes – 100 %B 2D gradient stop time: 1.39 minutes 2D cycle time: 1.89 minutes					
A	SM and ² D gradient – Method with the use of ASM using an ASM factor of 2.0					
ASM	ASM capillary: 5500-1302 (0.12 × 340 mm) ASM enabled (ASM factor: 2.0) Flush sample loop 3.0 times (0.24 minutes)					
² D Gradient	0.00 minutes - 5 %B 0.24 minutes - 5 %B 0.25 minutes - 50 %B 1.25 minutes - 100 %B ² D gradient stop time: 1.45 minutes ² D cycle time: 1.95 minutes					

 Table 2. HILIC-RPLC parameters.

First dimension						
Column	Agilent ZORBAX HILIC Plus, 2.1 × 100 mm, 3.5 μm					
Solvent	A) 50 mM ammonium formate in water adjusted to pH 4.0 B) Acetonitrile					
O minutes – 85 %B 10 minutes – 60 %B 11 minutes – 60 %B 12 minutes – 85 %B Stop time: 12 minutes Post time: 15 minutes Stop time and post time were set in the ² D pump to allo extension of the run time until completion of ² D analysicuts.						
Flow rate	0.500 mL/min					
Temperature	40 °C					
	DAD; 220 nm/4 nm, reference 360 nm/100 nm; 20 Hz					
Detection	To protect the flow cell from pressure pulses originating from valve switches, a pressure release kit (G4236-60010) was installed between the ¹ D DAD and the 2D-LC valve.					
Injection	Injection volume: 5 µL Sample temperature: 8 °C Needle wash: 3 seconds in water/acetonitrile (10/90)					
	Second dimension					
Column	AdvanceBio Peptide Mapping, 2.1 × 100 mm, 2.7 μm					
Solvent	A) Water + 0.1 % formic acid B) Acetonitrile + 0.1 % formic acid					
Temperature 40 °C						
Detection	DAD and mass selective detection (MSD)					
DAD	220 nm/4 nm, reference 360 nm/100 nm; 40 Hz					
	Mass selective detection (MSD)					
MSD Signals	Polarity: Positive Mode: Scan Mass range: 100–1,350 m/z Fragmentor: 100 V Gain: 1.00 Threshold: 150 Step size: 0.10 Peak width: 0.10 minutes Data storage: Full Stop time: No Limit The stop time was set to No Limit to allow extension of the run time until completion of ² D analysis of all cuts.					
Drying gas flow: 10.0 L/min Nebulizer: 35 psig Drying gas temperature: 200 °C Sheath gas temperature: 300 °C Sheath gas flow: 10.0 L/min Capillary voltage: 2,500 V Nozzle voltage: 300 V						

	2D-LC				
2D-LC Mode	Heart-cutting				
Flow rate	0.500 mL/min				
Sampling table	1.00 minute time based with 0.08 minutes sampling time 4.00 minutes peak based with 0.15 minutes sampling time 6.50 minutes peak based with 0.25 minutes sampling time 8.00 minutes peak based with 0.15 minutes sampling time 8.50 minutes peak based with 0.45 minutes sampling time 10.00 minutes off				
Peak detection	Mode: slope Upslope: 3 mAU/s Downslope: 1 mAU/s				
ASM and ² D gradient – Method without the use of ASM					
ASM	ASM disabled				
² D Gradient	0.00 minutes – 15 %B 3.00 minutes – 60 %B ² D gradient stop time: 3.00 minutes ² D cycle time: 5 minutes				
ASM and ² D gradi	ent - Method with the use of ASM using an ASM factor of 3.0				
ASM	ASM capillary: 5500-1301 (0.12 × 170 mm) ASM enabled (ASM factor: 3.0) Flush sample loop 3.0 times (0.74 minutes)				
² D Gradient	0.00 minutes – 3 %B 0.74 minutes – 3 %B 0.75 minutes – 15 %B 3.75 minutes – 60 %B ² D gradient stop time: 3.75 minutes ² D cycle time: 5.75 minutes				
ASM and ² D gradi	ent - Method with the use of ASM using an ASM factor of 5.1				
ASM capillary: 5500-1300 (0.12 × 85 mm) ASM enabled (ASM factor: 5.1) Flush sample loop 3.0 times (1.23 min)					
² D Gradient	0.00 minutes - 3 %B 1.23 minutes - 3 %B 1.24 minutes - 15 %B 4.24 minutes - 60 %B ² D gradient stop time: 4.24 minutes ² D cycle time: 6.24 minutes				

Results and Discussion

The incompatibility between the ¹D and ²D solvent systems is often referred to as one of the major challenges in 2D-LC method development². ASM provides a valve-based approach to address this problem by diluting the ¹D effluent with ²D mobile phase ahead of the ²D column².

To enable ASM, a valve topology consisting of the 2D-LC Valve Active Solvent Modulation (G4243A) and two MHC valves (G4242-64000) is used. Figure 1 shows the valve configuration.

Figure 2 explains the principle of active solvent modulation by showing the valve switching cycle.

In position 1, a fraction from the ¹D is sampled in Deck A (MHC valve A). Meanwhile, the flow from the ²D pump flows through Deck B (MHC valve B). Once the analysis of the fraction sampled in Deck A starts, the ASM valve switches to position 2. The flow coming from the ²D pump enters the ASM valve at port 5, and is split inside the valve between Deck A (port 8) and the ASM capillary (port 6). At port 10, the 2D flows through Deck A and the ASM capillary are reunited and led towards the ²D column. In this valve position, the fraction of ¹D effluent that was sampled in Deck A is diluted with ²D solvent before it reaches the ²D column. Using weak ²D solvent, focusing of the analytes at the head of the ²D column is achieved. The ASM valve stays in position 2 until the fraction of ¹D effluent is entirely transferred to the ²D (ASM phase). Once the transfer of the fraction is finished, the ASM valve switches to position 3. Now the entire flow coming from the ²D pump flows through Deck A, and the ASM capillary is disconnected from the flow. The ASM valve stays in position 3 until the end of the ²D analysis of the transferred fraction. Meanwhile, fractions from the ¹D effluent can be sampled on Deck B.

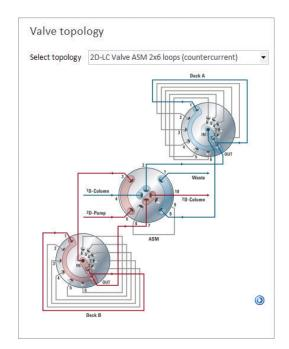


Figure 1. Valve topology for ASM, countercurrent configuration.

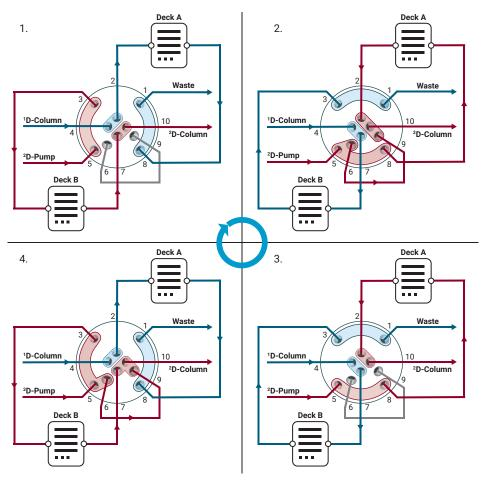


Figure 2. Valve switching cycle for ASM, countercurrent configuration.

Once the analysis of a fraction of ¹D effluent sampled in Deck B starts, the ASM valve switches to position 4 during the ASM phase to enable dilution of the ¹D effluent with ²D solvent. After the ASM phase, the ASM valve switches back to position 1, and the ASM capillary is again disconnected from flow.

The ASM factor describes the dilution of the sampled fraction of ¹D effluent, which results from the split ratio between the deck and the ASM capillary. This split ratio depends on the backpressures generated at the ASM capillary and the flowpath through the deck. In the flowpath through the deck, the backpressure is mainly generated by the transfer capillaries between the ASM valve and the MHC valve. Different ASM capillaries are available to adjust the split ratio and the ASM factor (Table 3). If, for example, ASM capillary 5500-1302 (0.12 × 340 mm) is installed, it will generate the same backpressure as the transfer capillaries between the ASM valve and the MHC valve (two capillaries of 0.12 × 170 mm). Therefore, the flow from the ²D pump is split with a ratio of 1:1 between the deck and the ASM capillary. This dilutes the sampled fraction of ¹D effluent by a factor of approximately 2 (ASM factor 2).

To use ASM, the ASM valve needs to be installed and configured in the 2D-LC configuration, as shown in Figure 1. The installed ASM capillary as well as transfer capillaries need to be defined in the 2D-LC configuration, as shown in Figure 3.

In the 2D-LC method, the use of ASM is activated and the length of the ASM phase is defined by specifying how many sample loop volumes are flushed through the loop to the ²D in the valve position where dilution takes place (Figure 4). The resulting duration of the ASM phase is displayed. Flushing the sample loop three times is typically enough to achieve quantitative transfer and the recommended default.

Table 3. Available ASM capillaries.

ASM Capillary	Length (mm)	ID (mm)	Volume (µL)	ASM factor*	Split ratio* (loop: ASM)
5500-1300	85	0.12	0.96	5	1:4
5500-1301	170	0.12	1.9	3	1:2
5500-1302	340	0.12	3.8	2	1:1
5500-1303	680	0.12	7.7	1.5	1:0.5

^{*} ASM factor and split ratio are based on transfer capillaries 5500-1270 (0.12 × 170 mm) and depend on solvent viscosity in the flowpaths through the ASM and transfer capillaries.

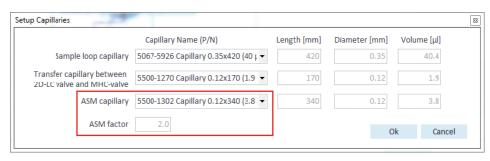


Figure 3. Configuration of the ASM capillary in the 2D-LC configuration.

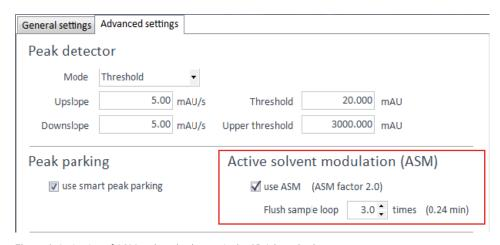


Figure 4. Activation of ASM and method setup in the 2D-LC method.

In the 2D-LC method, ²D gradients that were initially programmed without the use of ASM need to be delayed by the duration of the ASM phase. The ²D gradient stop time and ²D cycle time need to be increased accordingly.

For effective use of ASM, a solvent composition with low strength is chosen during the ASM phase. Figure 5 demonstrates the change from a ²D gradient initially programmed without the use of ASM (Figure 5A) to a

²D gradient with the use of ASM using an ASM phase of 0.24 minutes duration (Figure 5B).

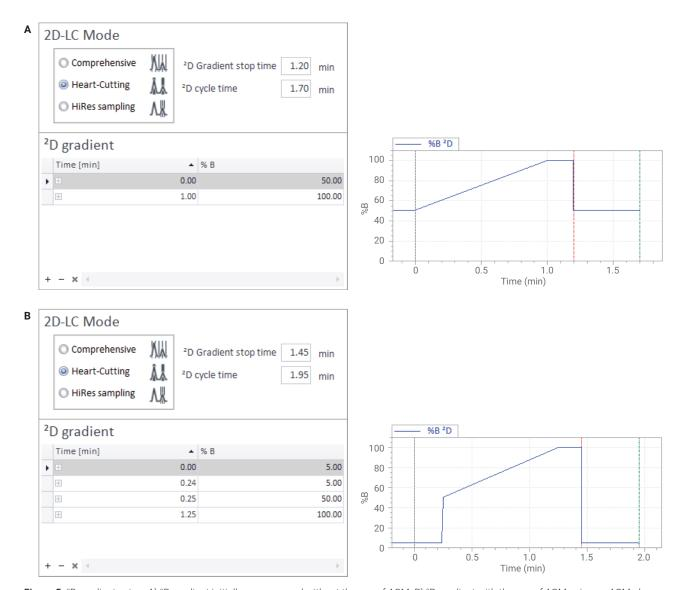


Figure 5. ²D gradient setup. A) ²D gradient initially programmed without the use of ASM. B) ²D gradient with the use of ASM using an ASM phase of 0.24 minutes duration.

RPLC-RPLC Analysis of Polycyclic Aromatic Hydrocarbons (PAHs)

Figure 6 shows the MHC 2D-LC analysis of PAHs using RPLC in both dimensions. For this analysis, an Agilent ZORBAX Eclipse Plus PAH column was chosen as the ¹D column, and a ZORBAX Eclipse Plus C8 column was used in the ²D. The retention of the PAHs is higher on the ZORBAX Eclipse Plus PAH column compared to the ZORBAX Eclipse Plus

C8 column. Generally, it is beneficial for an RPLC-RPLC analysis to place the more retentive column in the ²D to achieve focusing at the head of the ²D column^{2,3}. Here, the more retentive column was deliberately placed in the ¹D to show the challenge of this column order, and demonstrate that ASM enables focusing at the head of the ²D column. ASM therefore improves ²D resolution and sensitivity for this challenging combination.

As shown in Figure 6B, without the use of ASM, the ²D PAH peaks are distorted due to the lack of focusing at the ²D column head. This lack of focusing is because of the high percentage of organic solvent contained in the fractions of ¹D effluent transferred to the ²D. Figures 6C and 6D demonstrate that ASM enables focusing at the ²D column head, resulting in symmetrical and narrow ²D peaks.

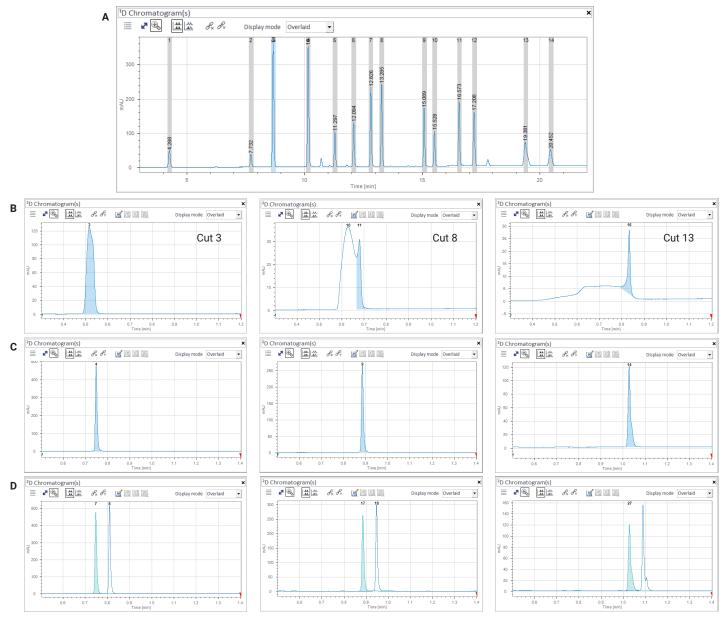


Figure 6. Multiple heart-cutting 2D-LC analyses of PAHs. A) ¹D chromatogram. B) ²D Chromatograms from the analysis without the use of ASM. C) ²D Chromatograms from the analysis with the use of ASM using an ASM factor of 1.5. D) ²D Chromatograms from the analysis with the use of ASM using an ASM factor of 1.5 (green trace) and using an ASM factor of 2 (blue trace).

In Figure 6C, ASM was used with an ASM factor of 1.5, which leads to symmetrical and narrow ²D peaks. Figure 6D compares the ²D separations obtained using ASM with an ASM factor of 1.5 (green trace) and an ASM factor of 2 (blue trace). Due to the higher dilution of ¹D effluent enabled using an ASM factor of 2, a further reduction in ²D peak width and increased peak height and sensitivity was obtained. In cut 13, a coelution that was already visible as a shoulder using an ASM factor of 1.5 was partially resolved using an ASM factor of 2.

To demonstrate that focusing at the ²D column head depends on the dilution of ¹D effluent before the ²D column and not solely on the ²D gradient used, the analysis was performed using the same ²D gradient with an initial hold at 5 %B (²D gradient used during the analysis with an ASM factor of 2) with the use of ASM (Figure 7, green trace) and without the use of ASM (Figure 7, blue trace). The fact that symmetrical and narrow ²D peaks are only obtained with the use of ASM clearly demonstrates the necessity of ASM to achieve focusing at the ²D column head.

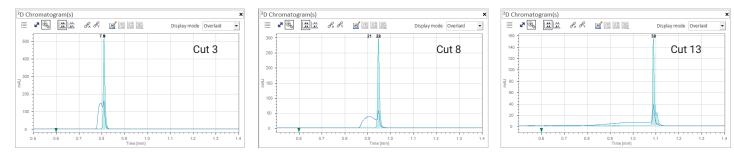


Figure 7. ²D Chromatograms from the multiple heart-cutting 2D-LC analysis of PAHs using the same ²D gradient with the use of ASM (green trace) and without the use of ASM (blue trace).

HILIC-RPLC Analysis of Peptides

Figure 8 shows the MHC 2D-LC analysis of peptides using HILIC in the ¹D and RPLC in the ²D. In the ¹D chromatogram (Figure 8A), coelution of peptides is already visible for the later-eluting peaks. In Figures 8B–D, the ²D extracted ion chromatograms (EICs) from the detected peptides are shown. Here, the most abundant ion detected for each

peptide was extracted (for example, for peptide 9, EIC 1,050 was extracted, which corresponds to the [M+2H]²⁺ ion of mol. wt. 2,099).

Figure 8B shows the analysis without the use of ASM. The peptides react differently to the high amount of organic solvent contained in the fractions of ¹D effluent transferred to the ²D compared to the PAHs. The peptides

show a partial breakthrough with the plug of ¹D effluent that was transferred to the ²D column. The percentage of peptide retained on the ²D column compared to the percentage that shows breakthrough is different for the different peptides analyzed. To confirm that the unretained peaks observed in the EICs originate from partial breakthrough of the peptides, the mass spectra of the

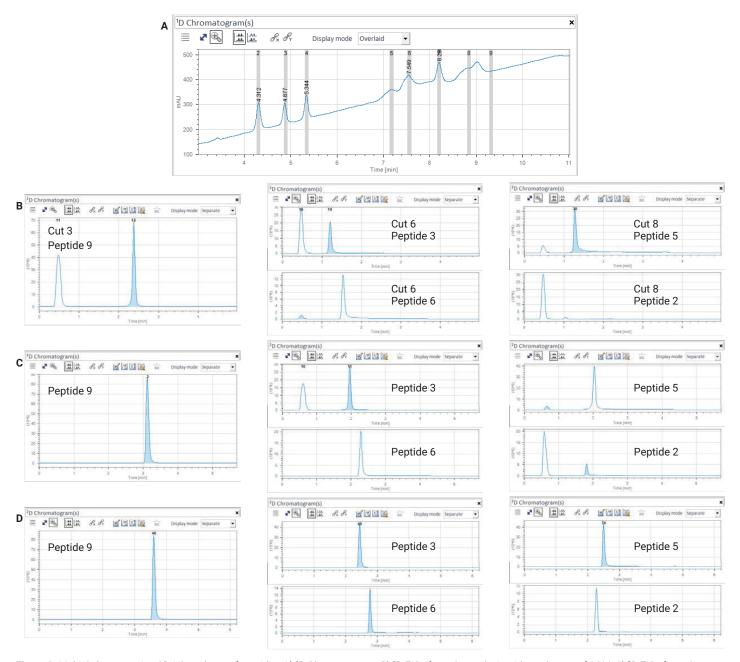


Figure 8. Multiple heart-cutting 2D-LC analyses of peptides: A) 1 D Chromatogram. B) 2 D EICs from the analysis without the use of ASM. C) 2 D EICs from the analysis with the use of ASM using an ASM factor of 3. D) 2 D EICs from the analysis with the use of ASM using an ASM factor of 5.

retained and unretained peaks were compared. The similarity of these mass spectra confirms the partial breakthrough. As an example, this is shown for cut 3 in Figure 9.

Figures 8C and 8D demonstrate that ASM enables complete retention of the peptides transferred to the ²D, and avoids the partial breakthrough observed without the use of ASM. Because of the focusing at the ²D column head, symmetrical ²D peaks were obtained. In Figure 8C, ASM was used with an ASM factor of 3, which avoids breakthrough for certain peptides (for example, peptide 9 in cut 3). For other peptides (for example, peptide 3 in cut 6), an ASM factor of 3 did not provide sufficient dilution of the ¹D effluent with weak ²D solvent to allow complete retention of the peptide. For these peptides, an ASM factor of 5 was necessary to completely avoid breakthrough, as shown in Figure 8D.

To demonstrate that avoiding breakthrough of the peptides depends on the dilution of ¹D effluent before the ²D column and not solely on the ²D gradient used, the analysis was performed using the same ²D gradient with an initial hold at 3 %B (²D gradient used during the analysis with an

ASM factor of 5) with the use of ASM (Figure 10, green trace) and without the use of ASM (Figure 10, blue trace). Without the use of ASM, breakthrough of the peptides was observed, confirming the necessity of ASM to achieve retention of the peptides.

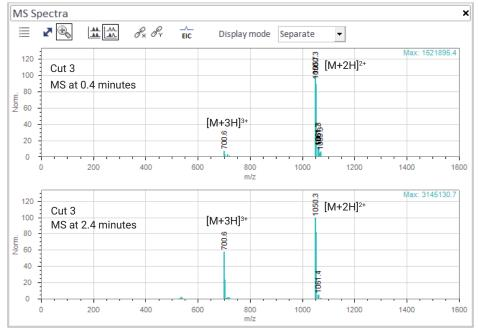
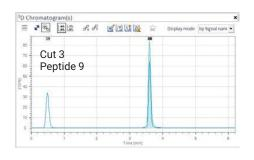
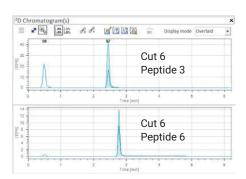


Figure 9. Mass spectra of the unretained (top) and retained (bottom) peak observed in cut 3 of the multiple heart-cutting 2D-LC analysis of peptides without the use of ASM.





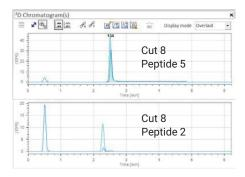


Figure 10. ²D EICs from the multiple heart-cutting 2D-LC analysis of peptides using the same ²D gradient with the use of ASM (green trace) and without the use of ASM (blue trace).

Conclusion

One major challenge in 2D-LC method development can be the incompatibility between the ¹D and ²D solvent systems. For combinations of separation systems where the analytes show low retention on the ²D column in the ¹D effluent, volume overload effects such as breakthrough or broad and distorted peaks in the ²D can be observed. Active solvent modulation provides a valve-based approach to address this problem by diluting the ¹D effluent with ²D mobile phase ahead of the ²D column, thereby enabling focusing at the ²D column head.

For an RPLC-RPLC analysis of PAHs, broad and distorted ²D peaks were observed without the use of ASM. Using ASM with an ASM factor of 2 enabled focusing at the ²D column head, leading to symmetrical and narrow ²D peaks. During an HILIC-RPLC analysis of peptides, partial breakthrough of the peptides was observed in the ²D without the use of ASM. Complete retention of the peptides was achieved using ASM with an ASM factor of 5.

References

- Stoll, D. R.; Carr, P. W.
 Two-Dimensional Liquid
 Chromatography: A State of the Art
 Tutorial. Analytical Chemistry 2017,
 89, 519-531.
- Stoll, D. R.; et al. Active Solvent Modulation: A Valve-Based Approach To Improve Separation Compatibility in Two-Dimensional Liquid Chromatography. Analytical Chemistry 2017, 89, 9260-9267.
- 3. Li, D.; Jakob, C.; Schmitz, O. Practical considerations in comprehensive two-dimensional liquid chromatography systems (LCxLC) with reversed-phases in both dimensions. Analytical and *Bioanalytical Chemistry* **2015**, *407*, 153-167.

www.agilent.com/chem

