



MULTIPLE HEART-CUTTING 2D LC FOR CHALLENGING SEPARATION PROBLEMS

Grant Von Wald, Matthias Pursch¹, Binghe Gu

Analytical Sciences

The Dow Chemical Company, Midland, MI; Stade, Germany¹

gvwald@dow.com

®Trademark of The Dow Chemical Company ("Dow") or an affiliated company of Dow

- Chemical samples in certain areas are of high and increasing complexity or dimensionality
 - Polymers can have a chemical composition distribution (CCD), functional end-group distribution (FTD) and/or molecular weight distribution (MWD)
 - New chemicals are derived from natural compounds which contain a significant number of possible structures (isomers etc.)
 - Determination of additives in polymer matrices
- Very complex samples require analytical tools of more than one dimension
- 2D LC has gained increasing attention due to high separation power, but requires significant method development for each application
- Recent introduction of more dedicated 2D LC instrumentation by several vendors is facilitating a more widespread use of this technology

Agilent 1290 Infinity 2D LC Solution:

- Based on 1290 Infinity LC system (binary pump)
- 1290 Infinity quaternary pump, TCC and DAD used in 1D
- 1200 bar switching valve (8 port) for comprehensive and heart-cutting 2D analysis
- Multiple heart-cutting interface
- LCIImage Software for comprehensive 2D LC data analysis

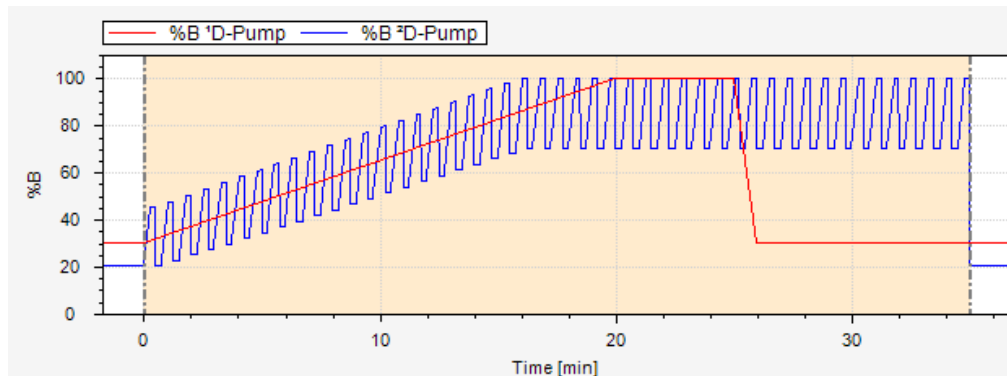


2D LC Separation Modes



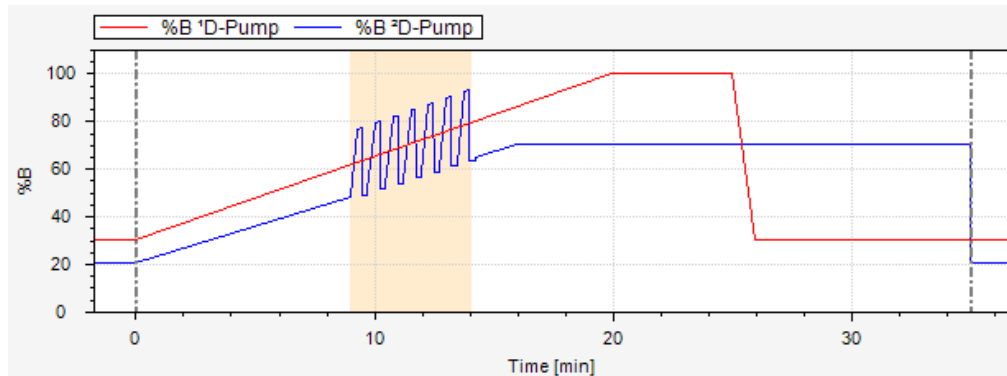
Comprehensive 2D (LCxLC):

- Complete 1st dimension effluent is sampled in discrete fractions (20-80 μL) – two loops
- Modulation time (20s – 1.5 min)



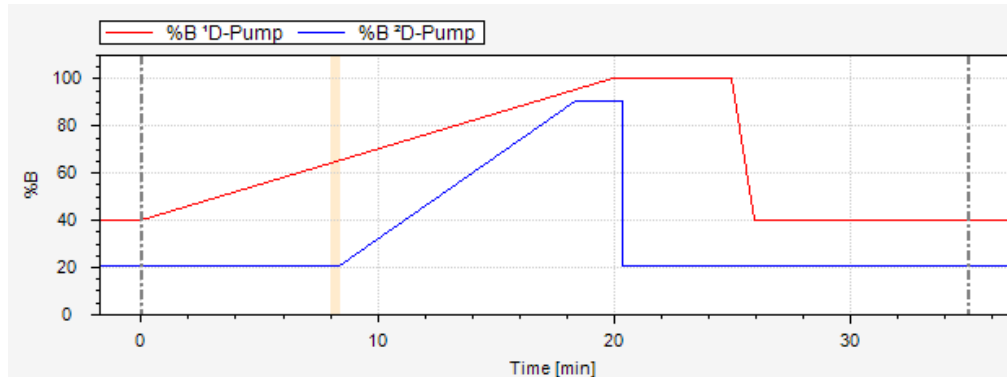
Selective Comprehensive 2D (sLCxLC):

- Defined region of interest in 1D is sampled in discrete fractions
- Transferred continuously to a second column via loops



Heart-cutting 2D LC:

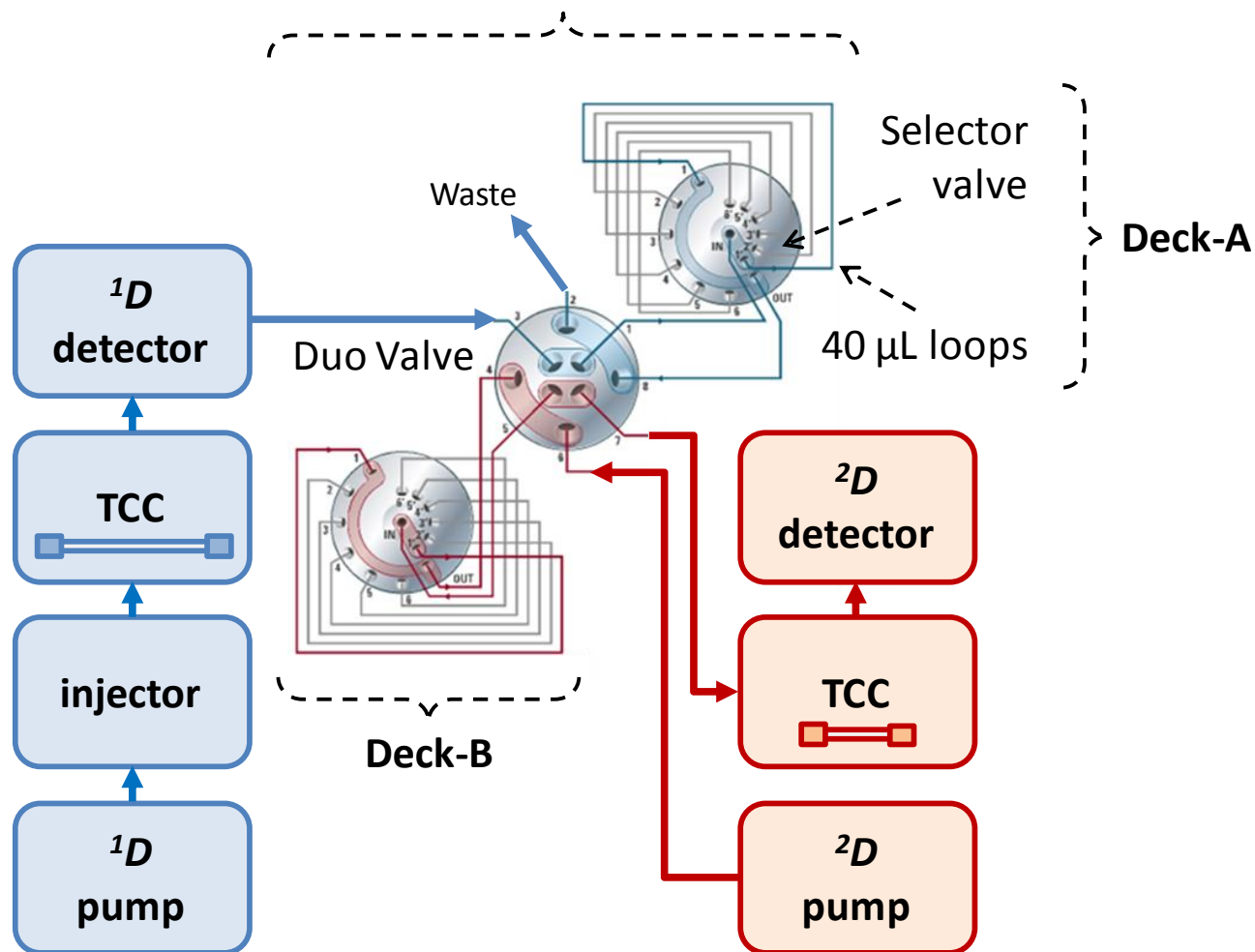
- Areas of interest are transferred from D1 to a second column via one loop (e.g., 40 μL)
- D2 analysis time is independent of D1 run time (“decoupled”), can be on order of 1.5 – 10 min



Multiple heart-cutting / selective comprehensive 2D



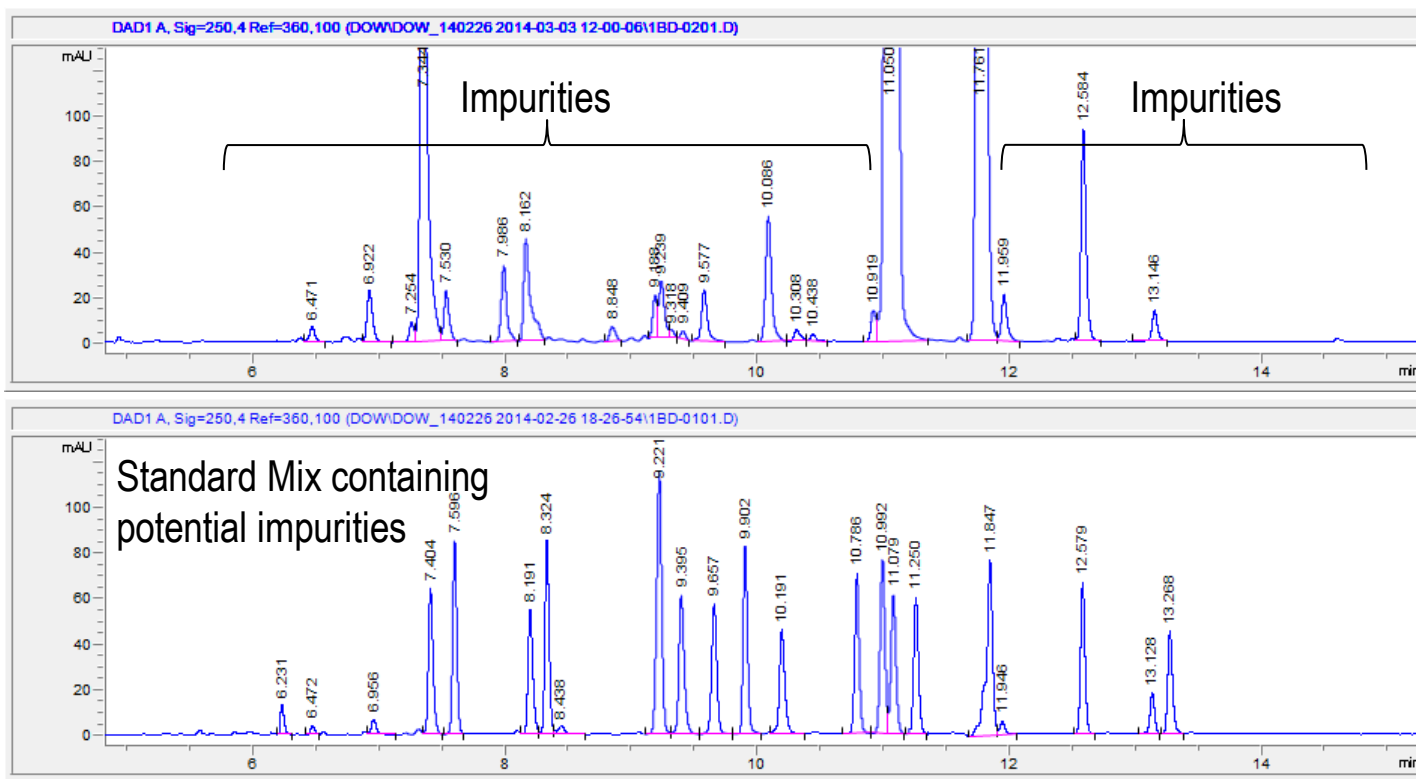
Multiple Heart-Cutting (MHC) Interface



Insecticide analysis



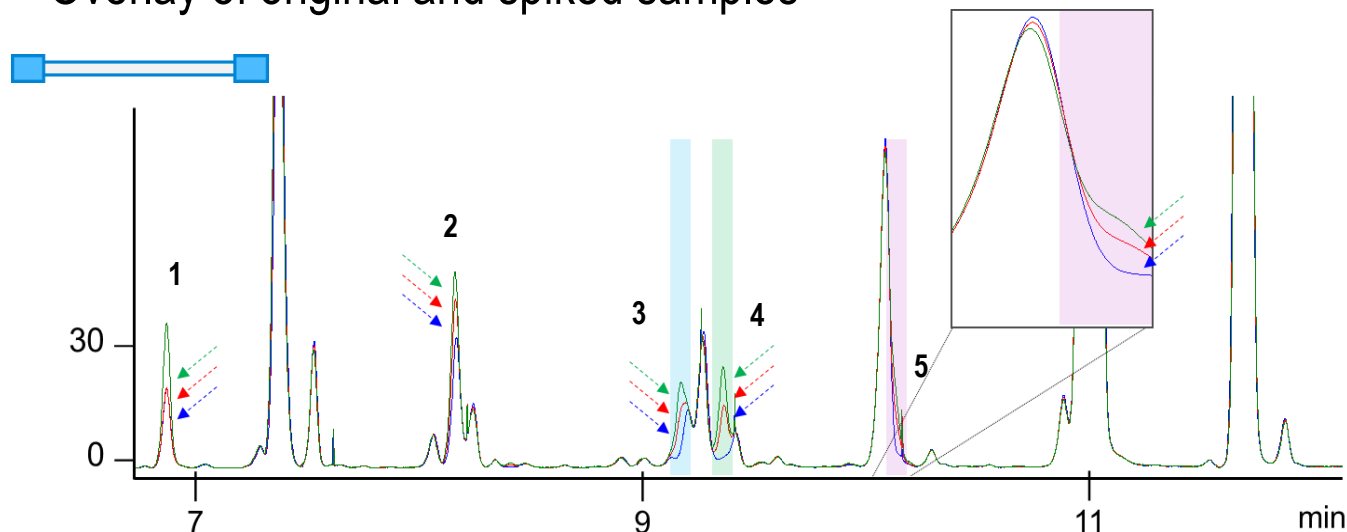
- Dow AgroSciences insecticide – contains 2 main components and 30-40 impurities
- D1: C₁₈ high resolution separation (150 mm, sub 2 µm particle column)
- Separation is good, but might be improved in certain areas



Multiple heart-cutting



Overlay of original and spiked samples



- Sample spiked w/ impurity standards.

Original sample.

Spike-1.

Spike-2.

- Integration hardly possible.

1D separation:

Zorbax SB-C18 150x2.1 mm 1.8 μ m

40°C, 0.5 mL/min

A: 10 mM ammonium acetate pH=6,

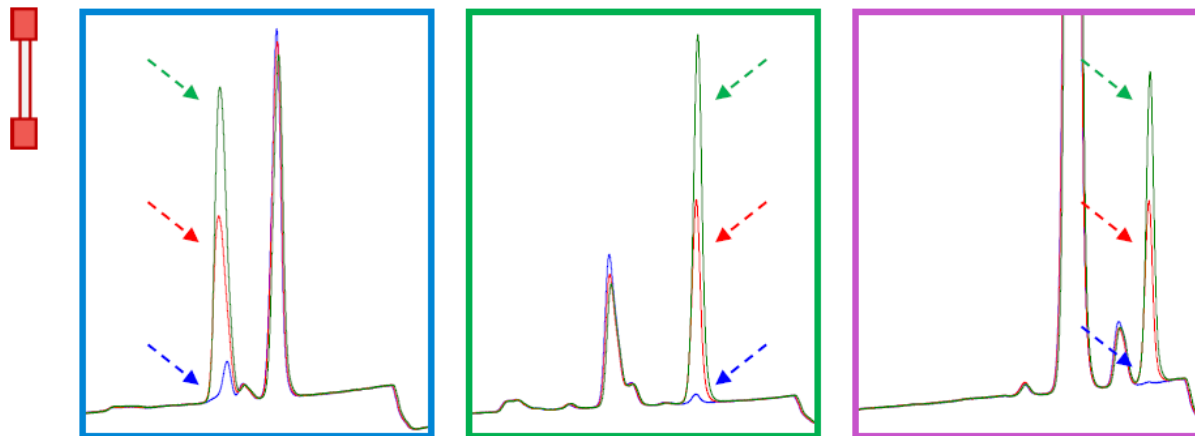
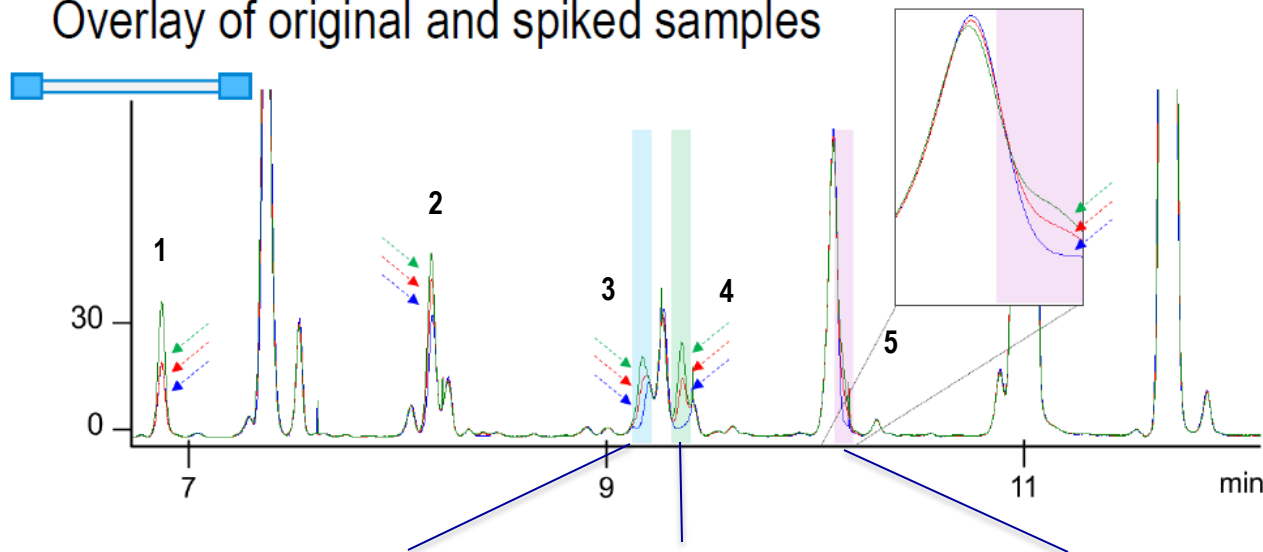
B: 20/80 methanol/acetonitrile

30-95% B/12 min

Multiple heart-cutting



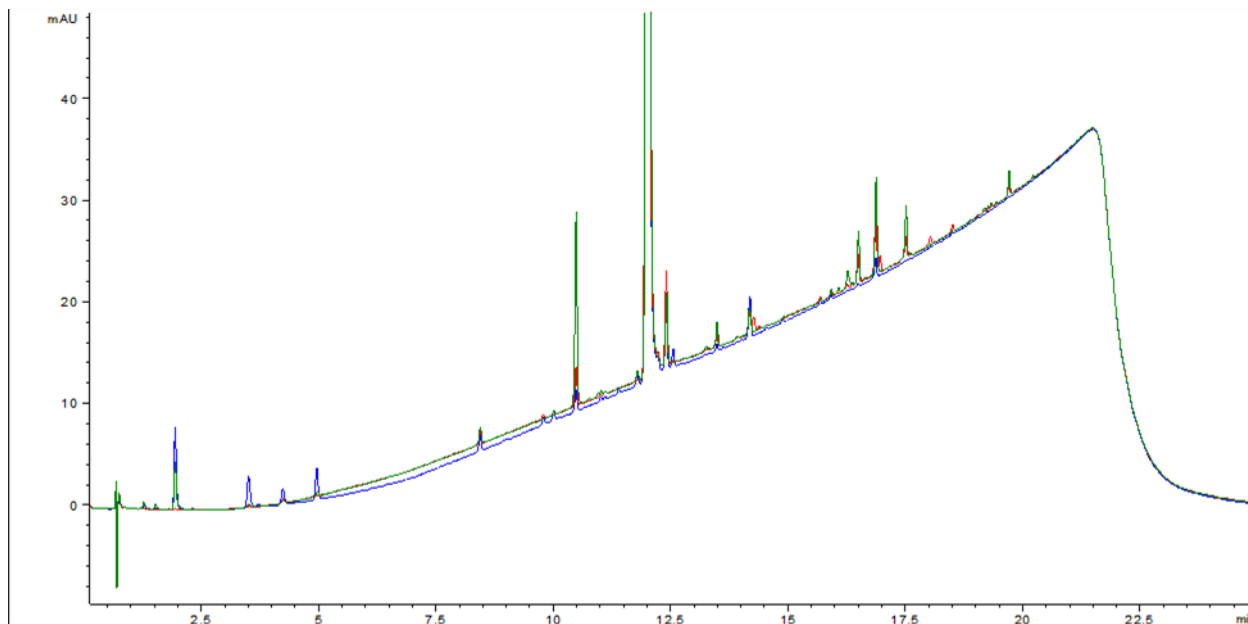
Overlay of original and spiked samples



- Multiple heart-cutting approach
- Much better separation in D2
- more accurate quantitation can be obtained with a suitable second dimension analysis

D2: Poroshell HPH-C18
50x3 mm 1.8 μ m
40°C 2 mL/min
A: 10 mM ammonium hydroxide pH=11
B: acetonitrile
55-85% B/1.3 min (1.5 min cycle time)

Crop Protection Chemical Application



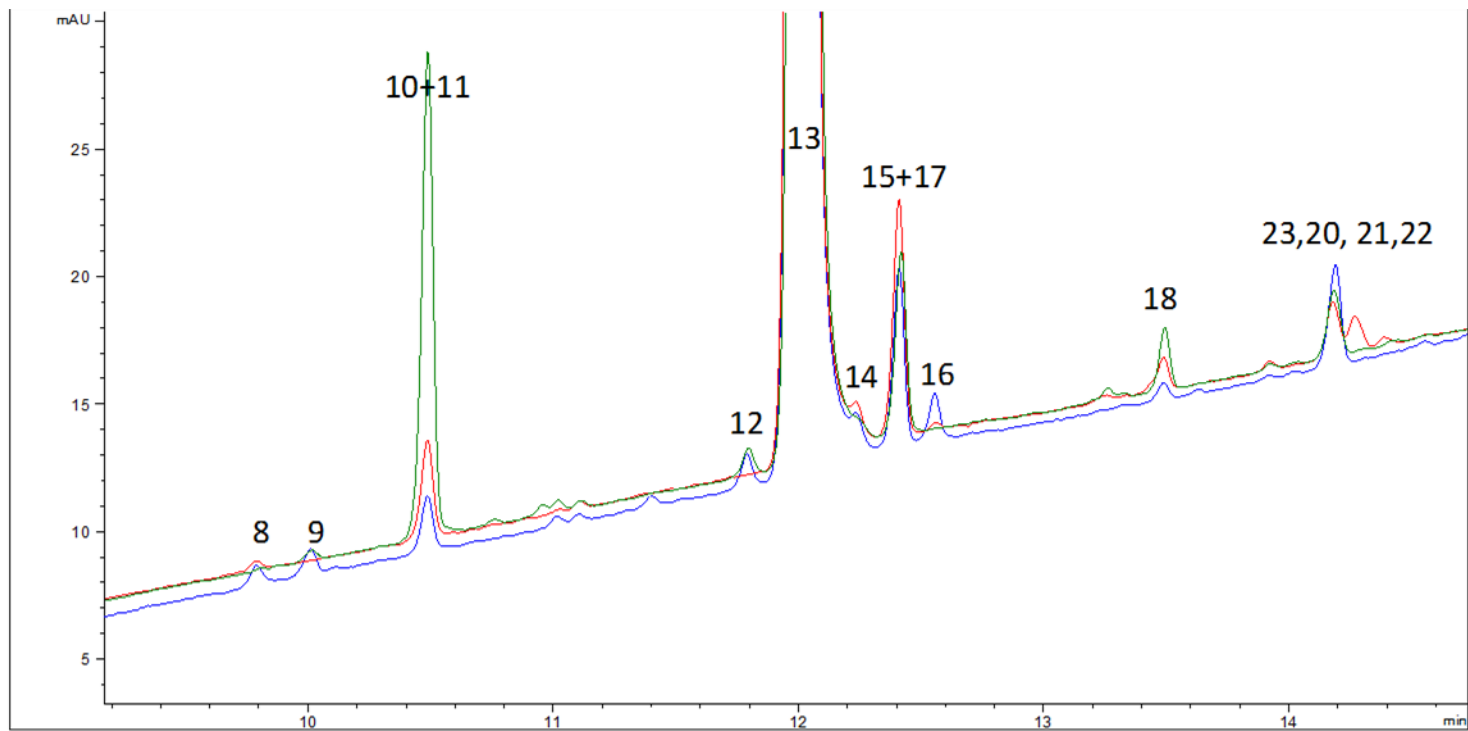
Poroshell 120 PhenylHexyl, 2.7 μm , 2.1x150mm

0.5 mL/min; 1 μL injection ; 35 $^{\circ}\text{C}$ column temp.; 245 nm;

Gradient: A: 0.1% formic acid/water B: 0.1% formic acid/acetonitrile

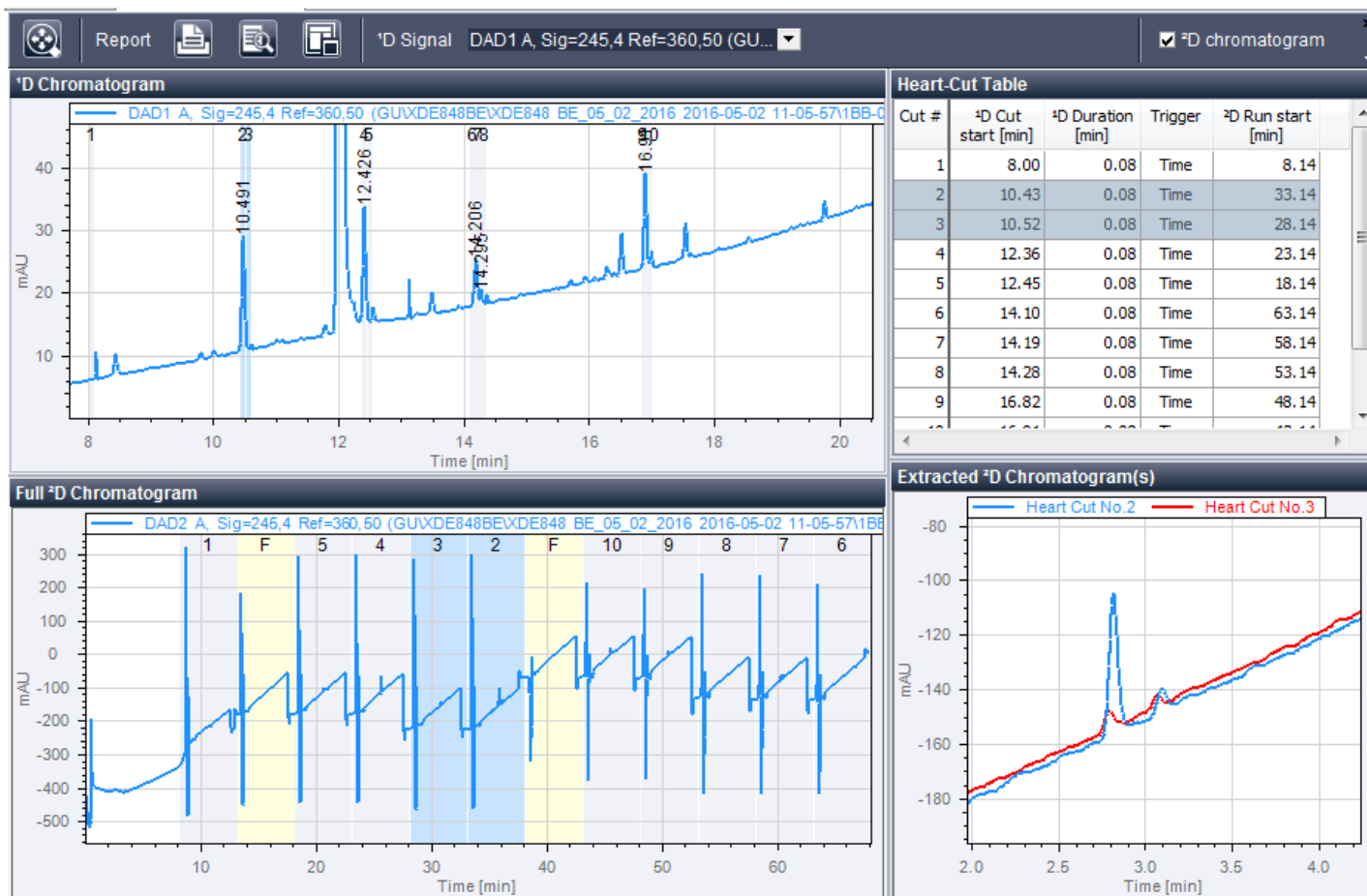
0-1 min 30% B, 1-20 min 30-100% B, 20.1-25 min 30% B

Expanded View



Co-elutions: 10/11; 15/17; 23/20;

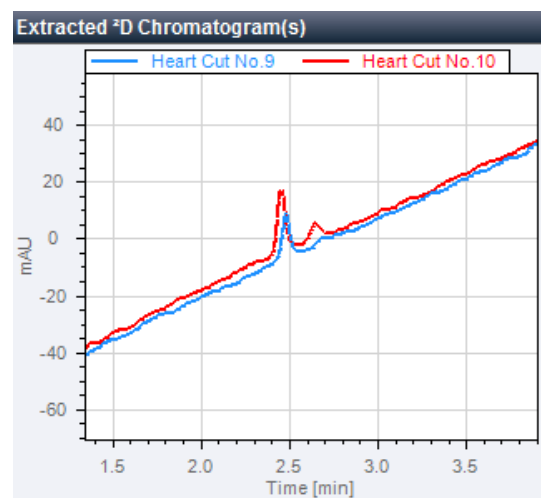
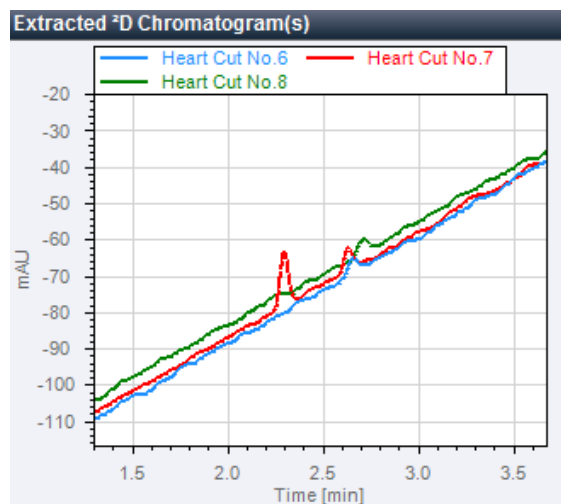
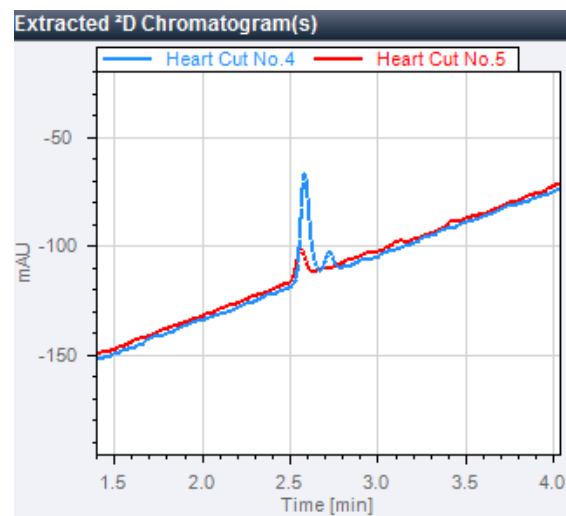
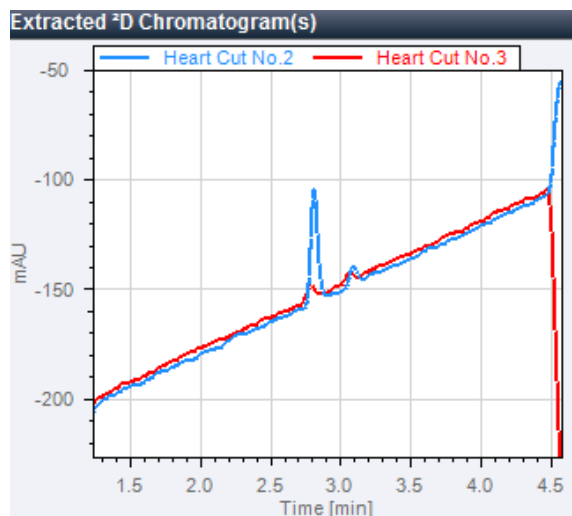
2-D MHC with Shifted Gradient



2D gradient:
 0 min 20% B
 (start % to 60%
 over 20 min)
 4 min 30% B
 (final % to 70%
 over 20 min)

1st dimension: Agilent Poroshell 120 Phenylhexyl, 2.7 μ m, 2.1x150 mm; 0.5 mL/min; 20 min from 30-100% ACN containing 0.1% FA;
 2nd dimension: Agilent Poroshell 120 EC-C18, 2.7 μ m, 4.6x50 mm; 0.5 mL/min; 4 min 20-30% THF containing 0.1% FA; cycle time, 5min.

Views of Each Cut





Reproducibility of Peak Areas

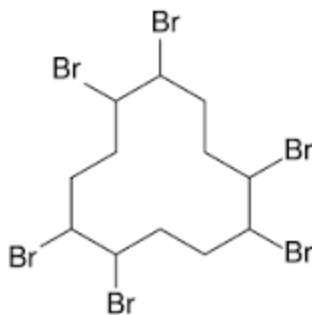
| Component | Average Area | R.S.D. (%) |
|----------------|--------------|------------|
| A (cut 5 only) | 66.3 | 24 |
| A (cut 4 only) | 184 | 13 |
| A (sum) | 250 | 4.5 |
| B* | 39.0 | 11 |
| C | 216 | 2.1 |
| D | 56.9 | 7.4 |
| E | 134 | 2.8 |
| F* | 26.6 | 19 |

*S/N for peaks B and F are <10. S/N of >10 for the remainder of the peaks.

HBCD analysis in Polystyrene

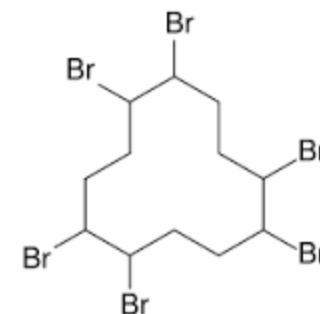
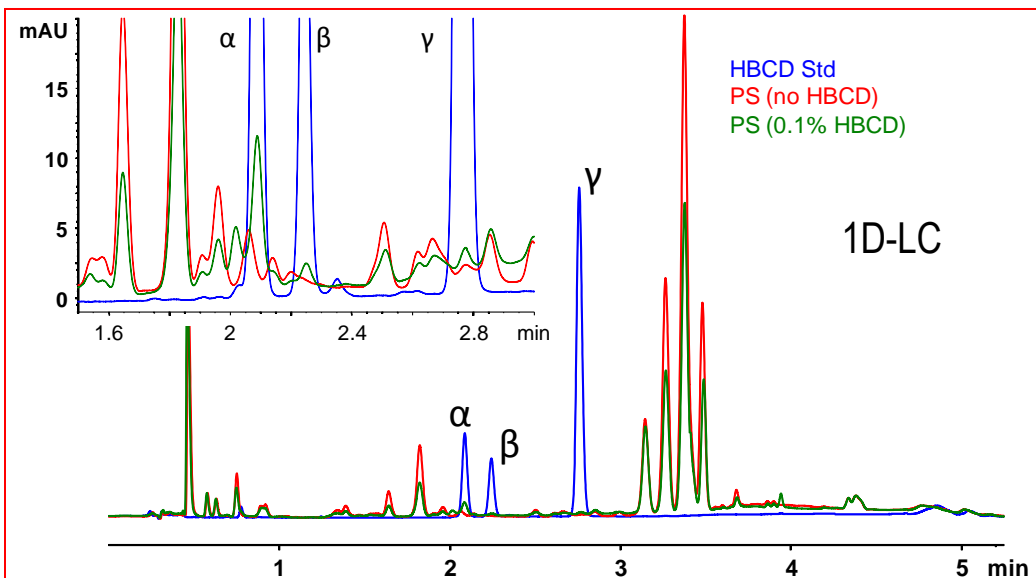


- Many polymers contain additives for improved UV-stability, flexibility and/or flame retardant properties
- In Polystyrene (PS) foam brominated compounds such as hexabromocyclododecane (HBCD) are used as flame retardant
- Need to measure HBCD flame retardant levels in PS foam at low level (0.1% and lower)
- Sensitive and selective method is needed

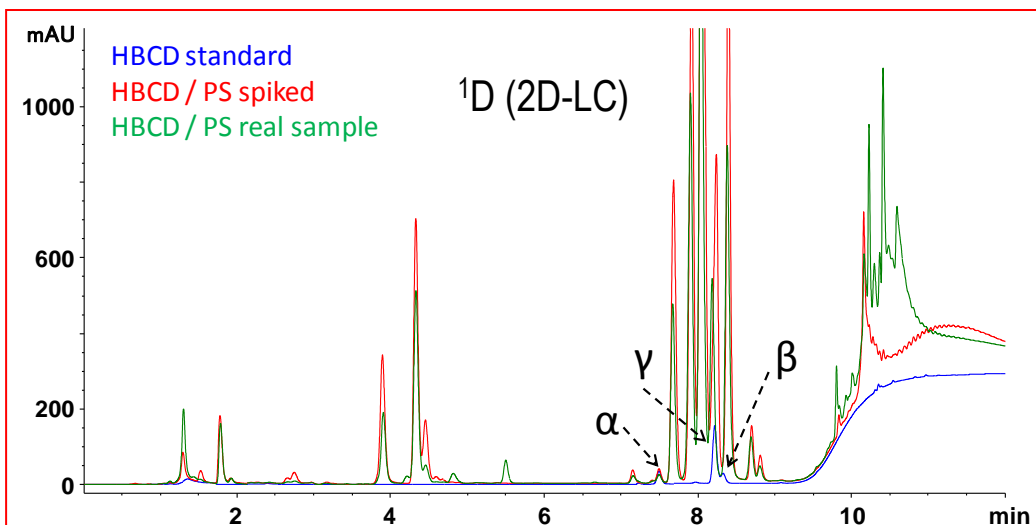


Hexabromocyclododecane (HBCD) –
exists in several isomeric forms

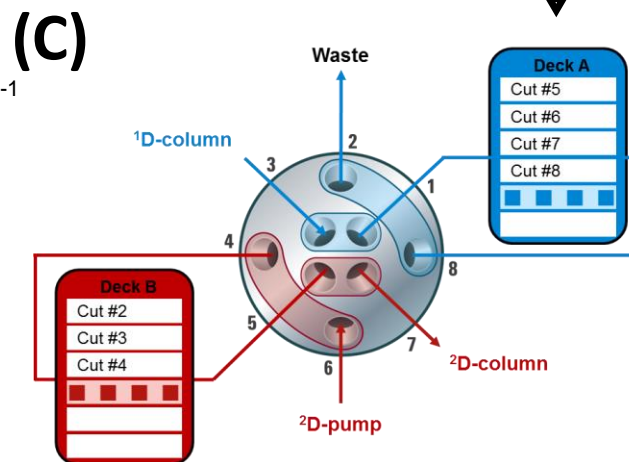
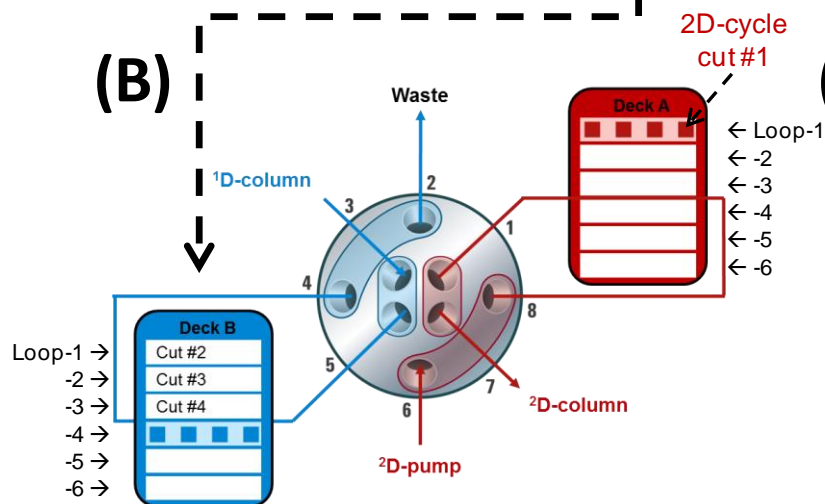
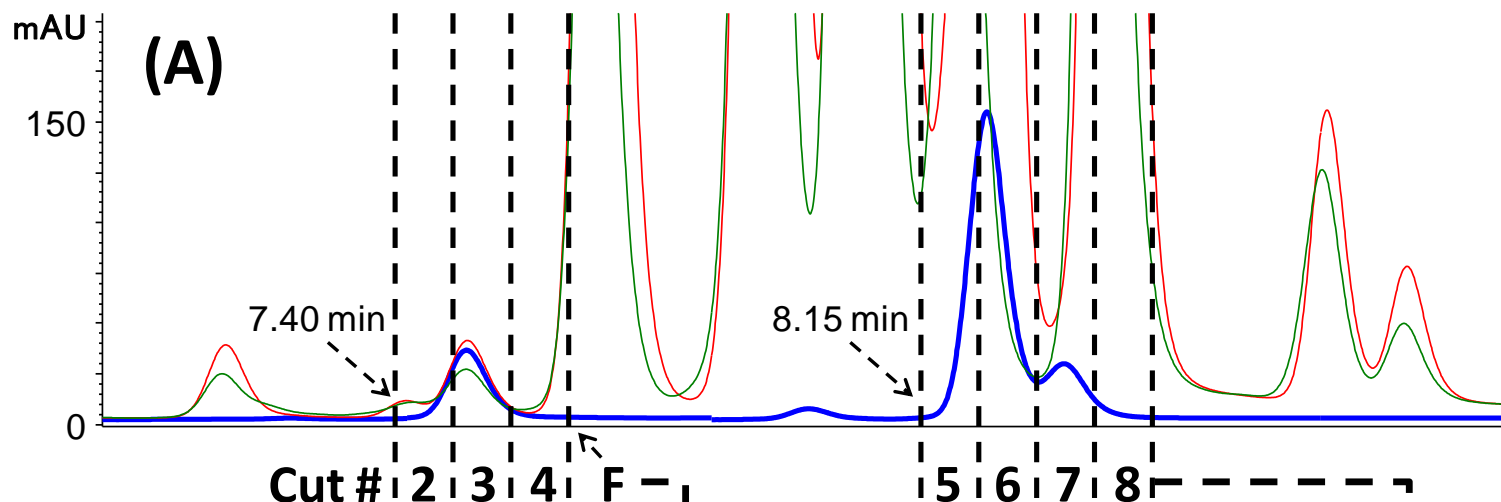
Multiple heart-cutting 2D LC



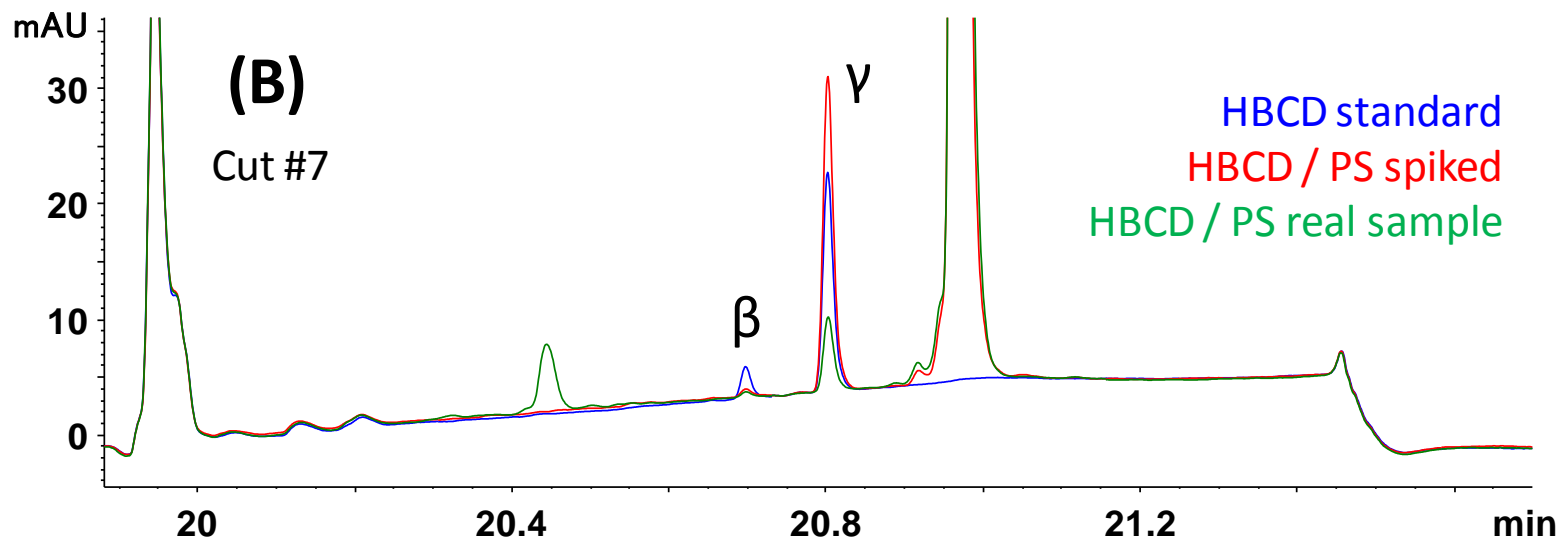
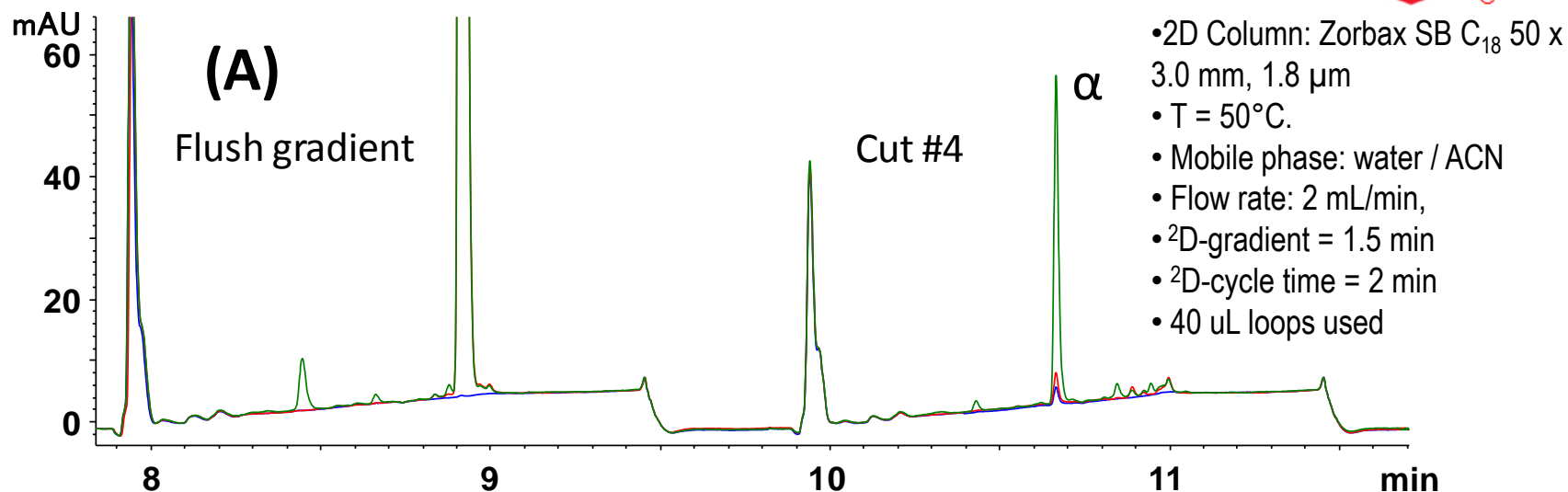
Hexabromocyclododecane (HBCD)



Changing from C18 (top) to phenyl (bottom) column in 1D provides a significant change in separation selectivity



2D MHC chromatograms

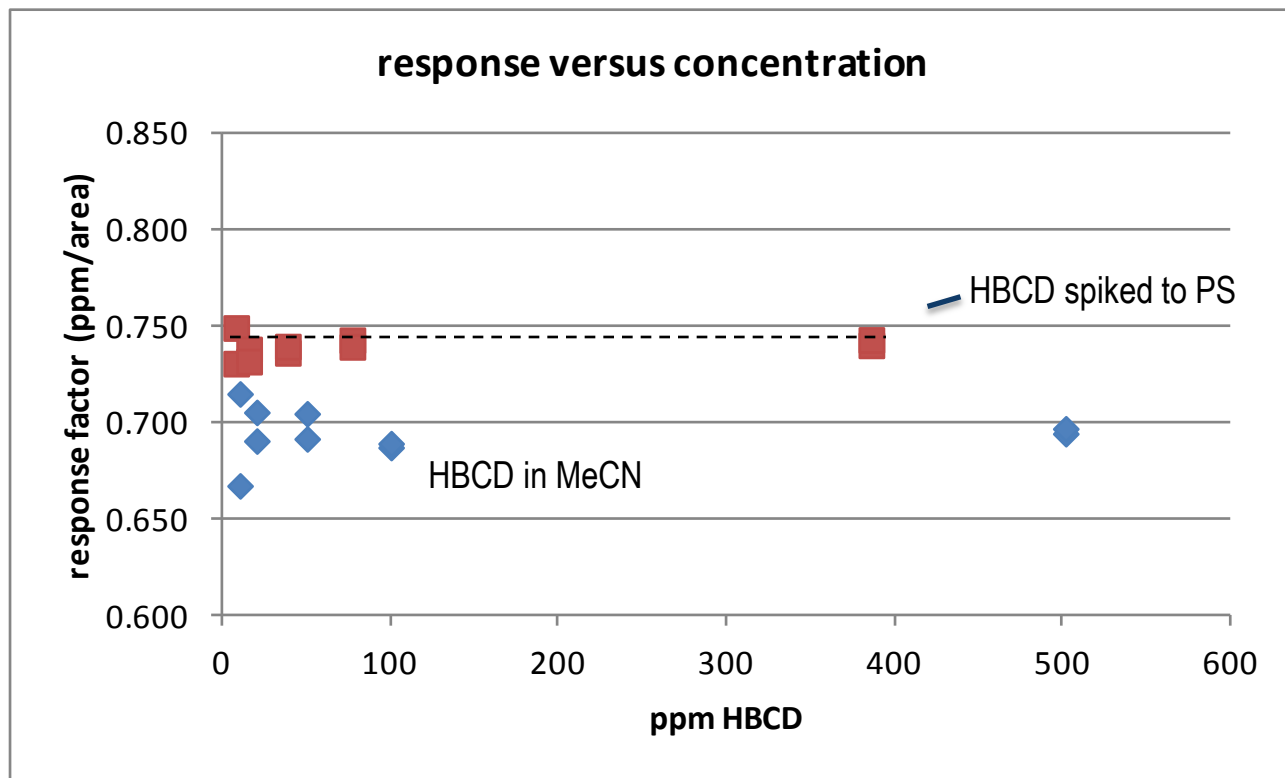


- Target analytes are fully separated from polymer components

Quantitation via 2D LC



- very good linearity for additive in MeCN and additive spiked to PS (response factors plotted against concentration)
- difference between standard and spiked polymer due to loss of additive in polymer upon precipitation



Repeatability



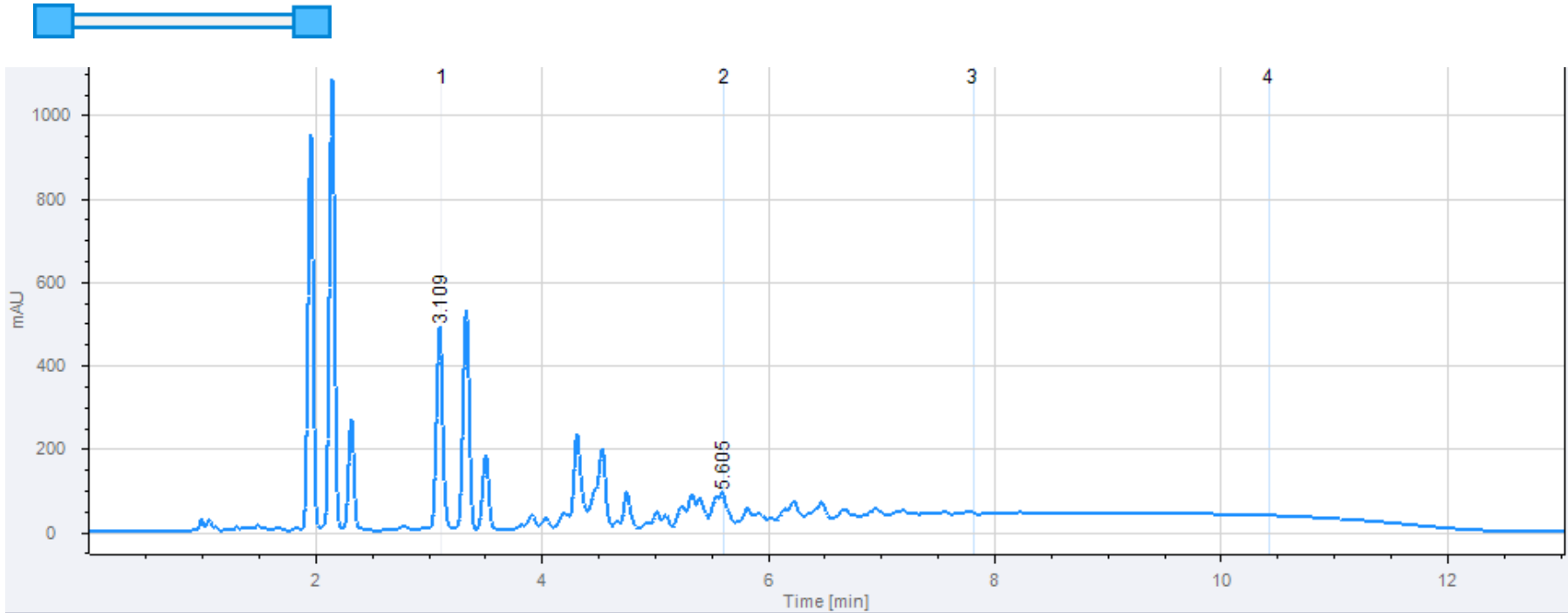
| | Peak Area | | | | | | | | | | | |
|---------|----------------------|--------|--------|-------|---------------------|--------|-------|----------------------|--------|--------|-------|------------|
| | HBCD isomer α | | | | HBCD isomer β | | | HBCD isomer γ | | | | Total HBCD |
| Run # | Cut #2 | Cut #3 | Cut #4 | Total | Cut #7 | Cut #8 | Total | Cut #5 | Cut #6 | Cut #7 | Total | |
| 1 | 21.90 | 46.30 | 1.20 | 69.40 | 0.75 | 10.00 | 10.75 | 1.20 | 1.60 | 5.90 | 8.70 | 88.85 |
| 2 | 20.00 | 48.10 | 1.40 | 69.50 | 0.78 | 10.20 | 10.98 | 1.20 | 1.70 | 6.10 | 9.00 | 89.48 |
| 3 | 16.90 | 51.00 | 1.60 | 69.50 | 0.53 | 10.10 | 10.63 | 1.10 | 1.30 | 6.40 | 8.80 | 88.93 |
| 4 | 16.60 | 51.30 | 1.60 | 69.50 | 0.46 | 10.00 | 10.46 | 1.20 | 1.00 | 6.40 | 8.60 | 88.56 |
| 5 | 16.30 | 51.60 | 1.60 | 69.50 | 0.34 | 10.10 | 10.44 | 1.40 | 1.20 | 6.20 | 8.80 | 88.74 |
| 6 | 12.60 | 54.90 | 1.80 | 69.30 | 0.37 | 10.00 | 10.37 | 1.40 | 1.10 | 6.50 | 9.00 | 88.3 |
| 7 | 18.10 | 49.90 | 1.50 | 69.50 | 0.48 | 10.10 | 10.58 | 1.20 | 1.40 | 6.10 | 8.70 | 88.78 |
| 8 | 11.20 | 56.10 | 2.10 | 69.40 | 0.33 | 9.80 | 10.13 | 1.50 | 0.74 | 6.60 | 8.84 | 88.04 |
| 9 | 21.50 | 46.90 | 1.40 | 69.80 | 0.70 | 10.20 | 10.90 | 1.20 | 1.60 | 6.10 | 8.90 | 89.6 |
| 10 | 28.20 | 40.70 | 1.10 | 70.00 | 1.30 | 9.90 | 11.20 | 1.10 | 2.30 | 5.50 | 8.90 | 90.1 |
| Average | 18.33 | 49.68 | 1.53 | 69.54 | 0.60 | 10.04 | 10.64 | 1.25 | 1.39 | 6.18 | 8.82 | 88.94 |
| StdDev | 4.90 | 4.46 | 0.29 | 0.21 | 0.30 | 0.13 | 0.32 | 0.14 | 0.44 | 0.32 | 0.13 | 0.62 |
| RSD | 26.72 | 8.97 | 18.75 | 0.30 | 48.96 | 1.26 | 3.00 | 10.83 | 31.32 | 5.22 | 1.49 | 0.70 |

MHC 2D LC of oligomers/polymers

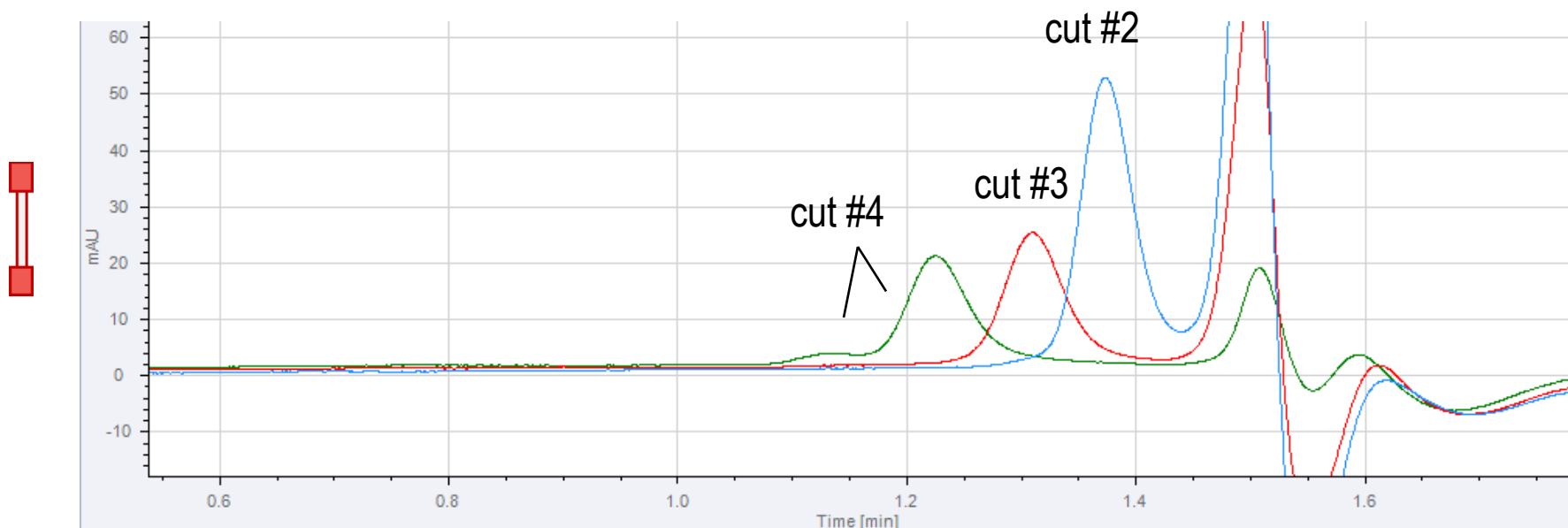


- Correlation of polarity with molecular weight (hydrodynamic volume)
- Epoxy resin used as example
- D1: Reversed phase separation – Core-shell C18 (MeCN/water) gradient
- D2: Size Exclusion Chromatography – THF

MHC 2D LC of oligomers/polymers



- D1: Oligomer and isomer separation according to differences in polarity & molecular weight
- Initial peaks are well resolved
- Resolution deteriorates toward the end of the separation
- Cuts are made at several areas of the 1D separation



- D2: SEC analysis
- Information on molecular weight can be obtained (calibration with standards or sample of known composition)
- Two species of different molecular weight are seen in cut #4 (two components of similar polarity but different chemical structure)
- MHC 2D LC provides additional structure information

- 2D LC is a valuable technique for challenging separation problems, including qualitative and quantitative target analysis
- Allows to differentiate complex samples in a way which would not be possible with 1D methods
- Depending on the problem either full comprehensive mode or heart-cutting/selective comprehensive modes can be applied
- Heart-cutting methods reduce dependence of 2D on 1D and simplify method development. Crucial parameters include:
 - To reduce peak broadening in the 2D:
 - » 1D column should be smaller I.D., less retentive
 - » 2D column should be larger I.D., more retentive
 - Transfer volume, measured to be 19 μL
 - Cuts from the same peak need to be sent to the same deck
 - Shifted gradient is an asset, but the appearance of the shift is misleading in the method editor
- Using accurate mass MS detection, thorough structural analysis can be performed

Acknowledgment



Marian Rinken, Paul Lewer, Andreas Schweizer-Theobaldt, Antje Wegener

Agilent Technologies (Stephan Buckenmaier, Edgar Naegele, Jens Trafkowski)

HBCD results were previously reported in *Anal. Chem.* **2015**, 87, 5310-5317

and C&EN Webinar September 30, 2015 “Applications of 2D-LC and 2D-LC/MS for Polymer and Crop Protection Chemical Analysis”