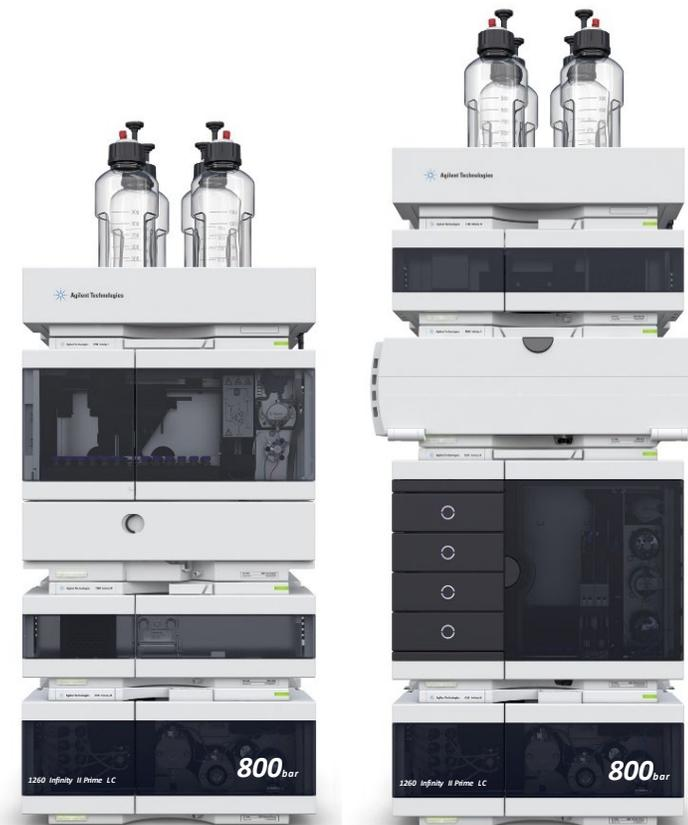


Eliminate your Application and Chromatography Challenges

LC Application Scientist Session

Information Contributed by USA HPLC Applications Scientist Team

Speaker: Lorena Lopez
LC Application Scientist
Lexington, MA



This Session will cover

Troubleshooting your Method, Application / Chromatography

- Sample prep considerations
- Troubleshooting by following the LC flow. Mobile Phase and System Hygiene
- Step through Method Setup to highlight parameters that are critical but often overlooked or misunderstood

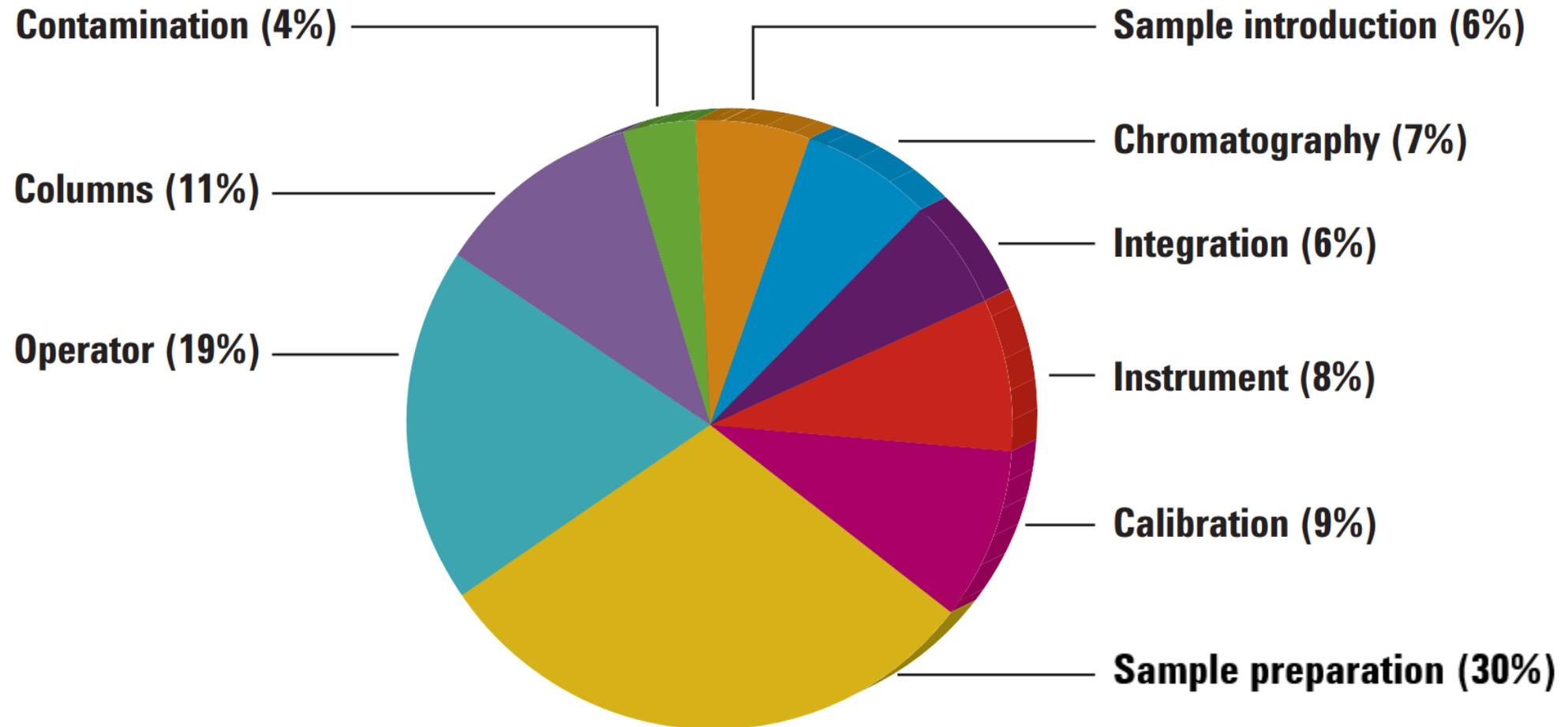
Optimizing your Application / Method Transfer considerations

- What needs to be considered when implementing App notes and transferring method between systems
- Delay volume, column void volumes
- Capillary selection and connections
- Column considerations (dimension and particle size)

Advancing your Application / Chromatography

- How to choose the appropriate LC system for the application that is going to be run.
- Where to find resources or info

Sources of Error Generated During Chromatographic Analysis



Data taken from Agilent Technologies survey

What is Method Transfer?

Different Aspects

From particle size to particle size (HPLC to UHPLC)

- ⇒ often associated with a change of instruments due to constraints in max. pressure, extra-column band broadening etc.. Some fine-tuning of method might be required due to frictional heating effects
- ⇒ available tools: Agilent Method translator, Third Party Method development SW (S-Matrix, ACD Labs, ChromSword, etc.)

From one eluent type or phase chemistry to an other

- ⇒ Method development
- ⇒ available tools: Method development SW (ACD Labs, ChromSword, S-Matrix etc.)

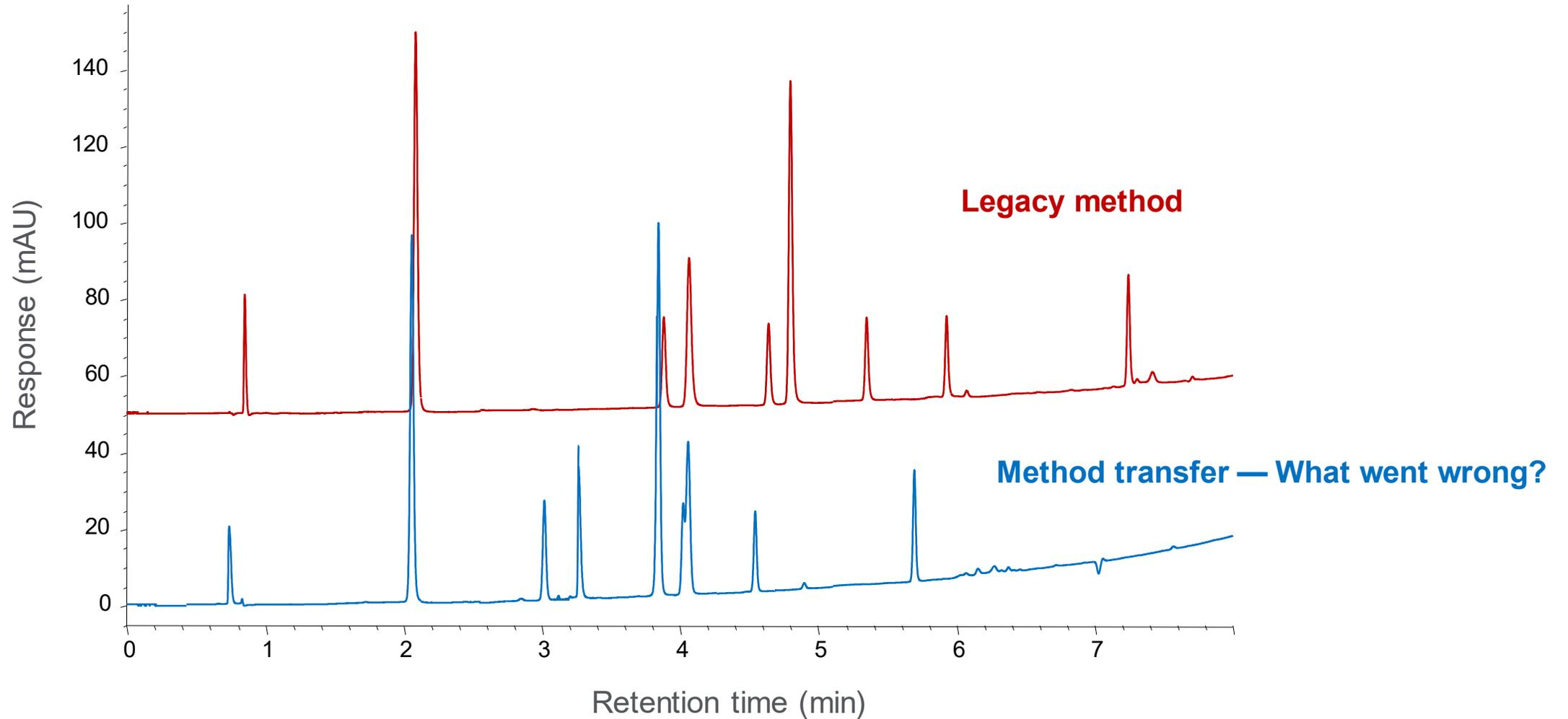
From column dimension to column dimension (e.g. 4.6 to 2.1 mm i.d., 50 mm to 100 mm)

- ⇒ recalculate flow rates, recalculate gradient times, adjust connection capillaries and flow cells, due to delay volume results may vary, method might need revalidation
- ⇒ available tools: Method Translator

From Instrument to Instrument

- ⇒ with method/instrument change: Isocratic Hold / Pre-injection, Vol. Modification
- ⇒ available tools w/o method or instrument change: none until ISET

Method Transfer Considerations & Application Optimization



Design Differences (U)HPLC Systems

- Delay volumes (gradient formation)
- Power range (flow x pressure)
- Extra column volume
- Temperature
- Data rates
- Sensitivity

Problem:

Instrument to Instrument Method Transferability

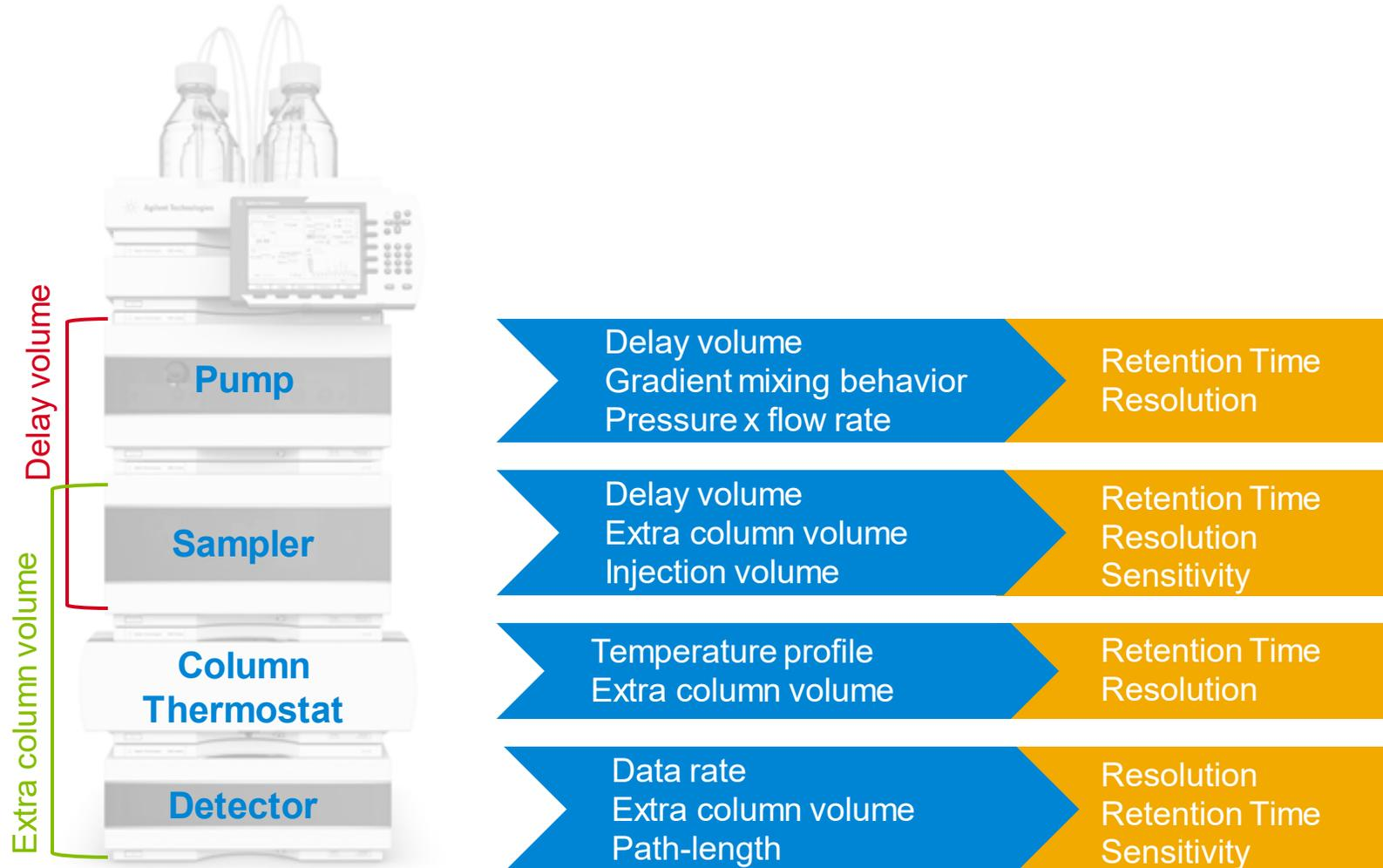
- Not possible
- Requires Re-Development
- Requires Re-Validation

Resulting in huge additional cost factor



Instrument to Instrument Method Transferability

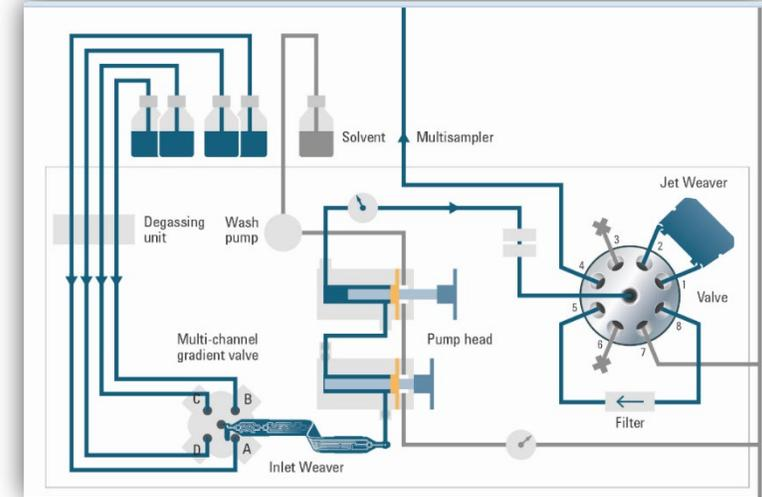
- Important Parameters



Comparison of Gradient Delay Volume (Dwell Volume)

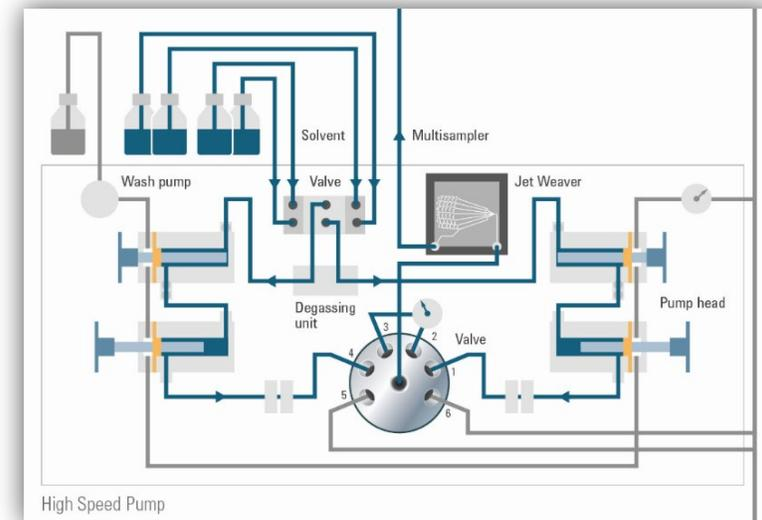
1290 Infinity II Flexible Pump (Quaternary)

- Integrated Degasser
- 4 solvent channels with concurrent mixing of all 4 channels
- Lower in price, typically, than binary pump



1290 Infinity II High Speed Pump (Binary)

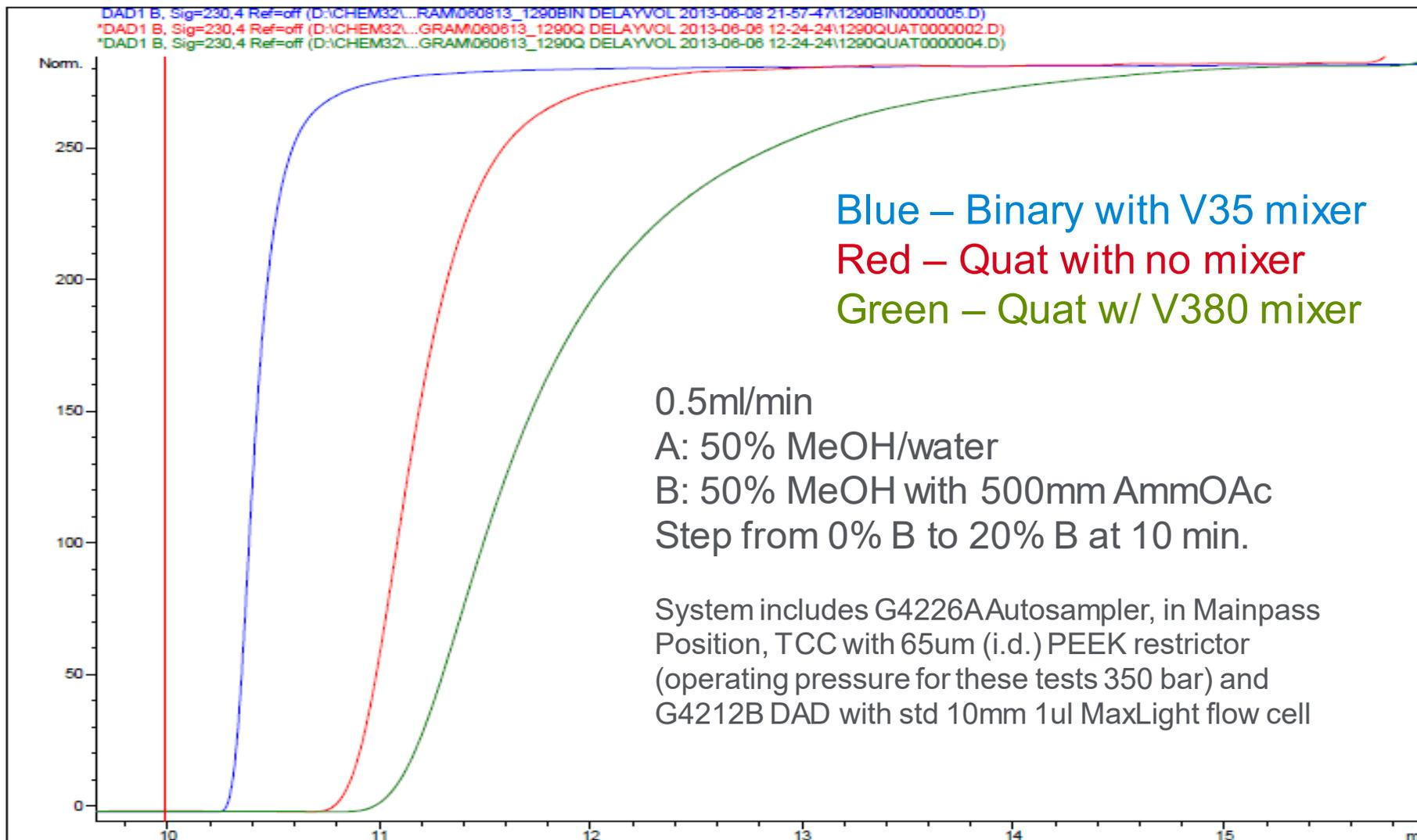
- Integrated Degasser
- 4 solvent channels available, mixing of 2 channels possible
- Better performance concept is widely accepted
- Greater control over dwell volume vs. Quaternary pump



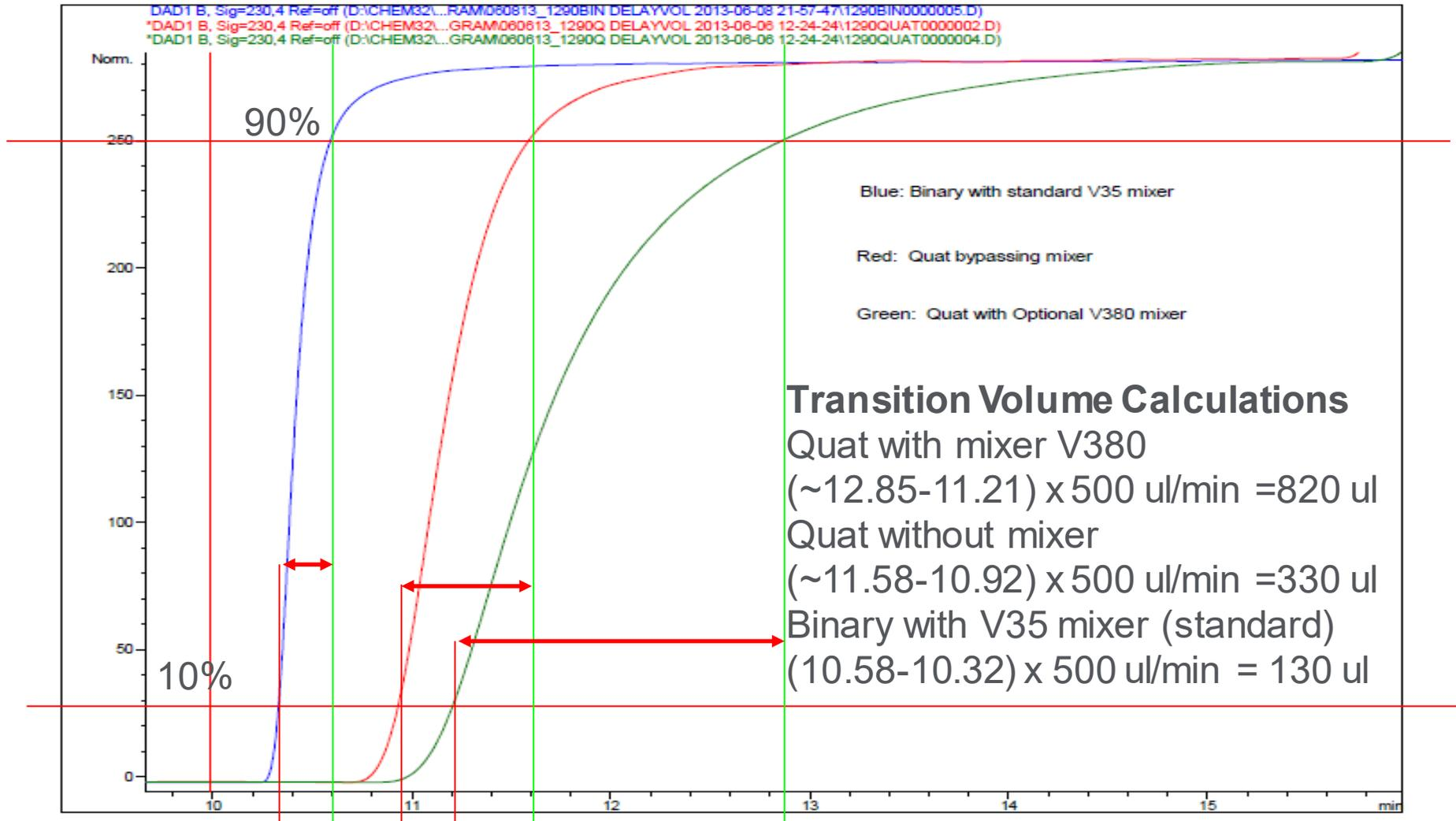
*Incl. Degasser for both pumps

**with solvent selection valve

1290 Family Delay Volume Profiles

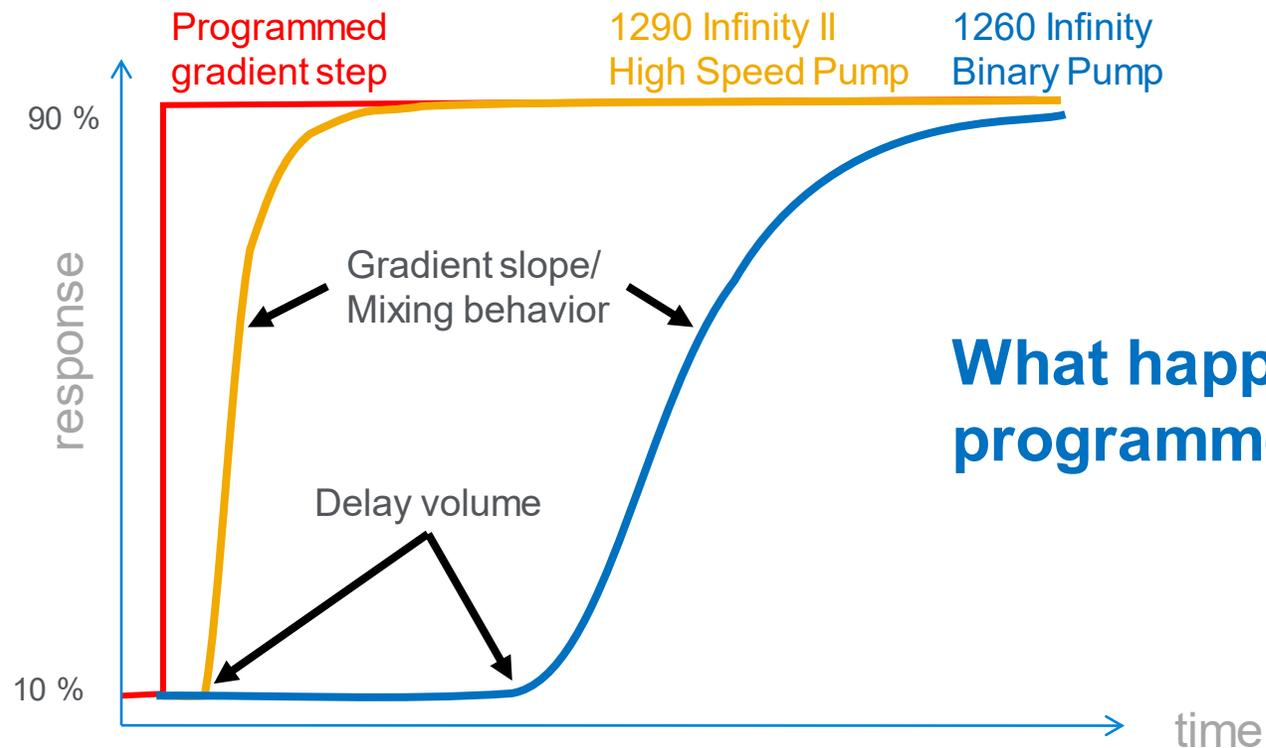


1290 Delay Volume – 10-90% Transition



Method Transfer to UHPLC instruments

Impact of delay volume and mixing behavior...



What happens with a programmed gradient?

1200 Series

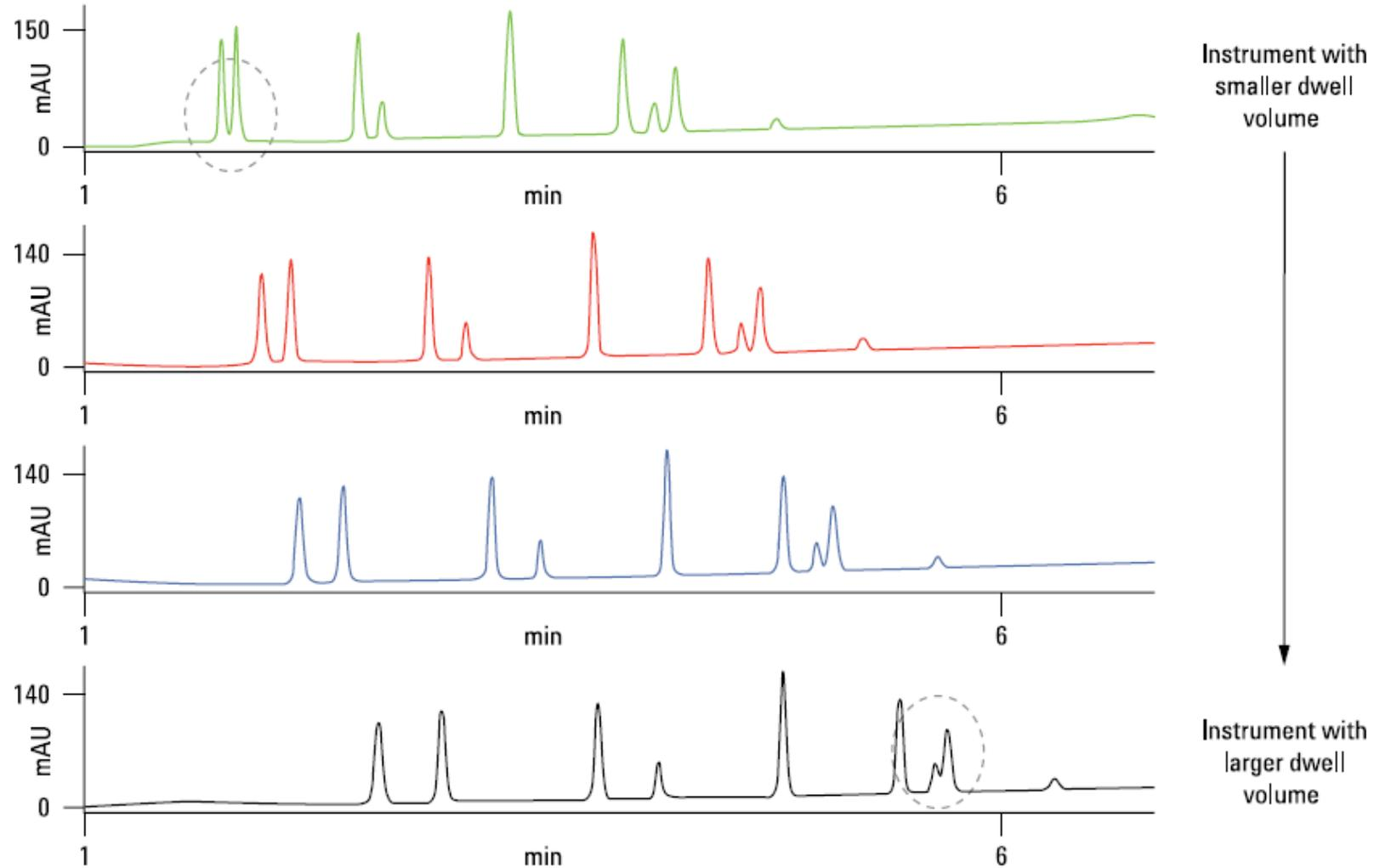


1290 Infinity II

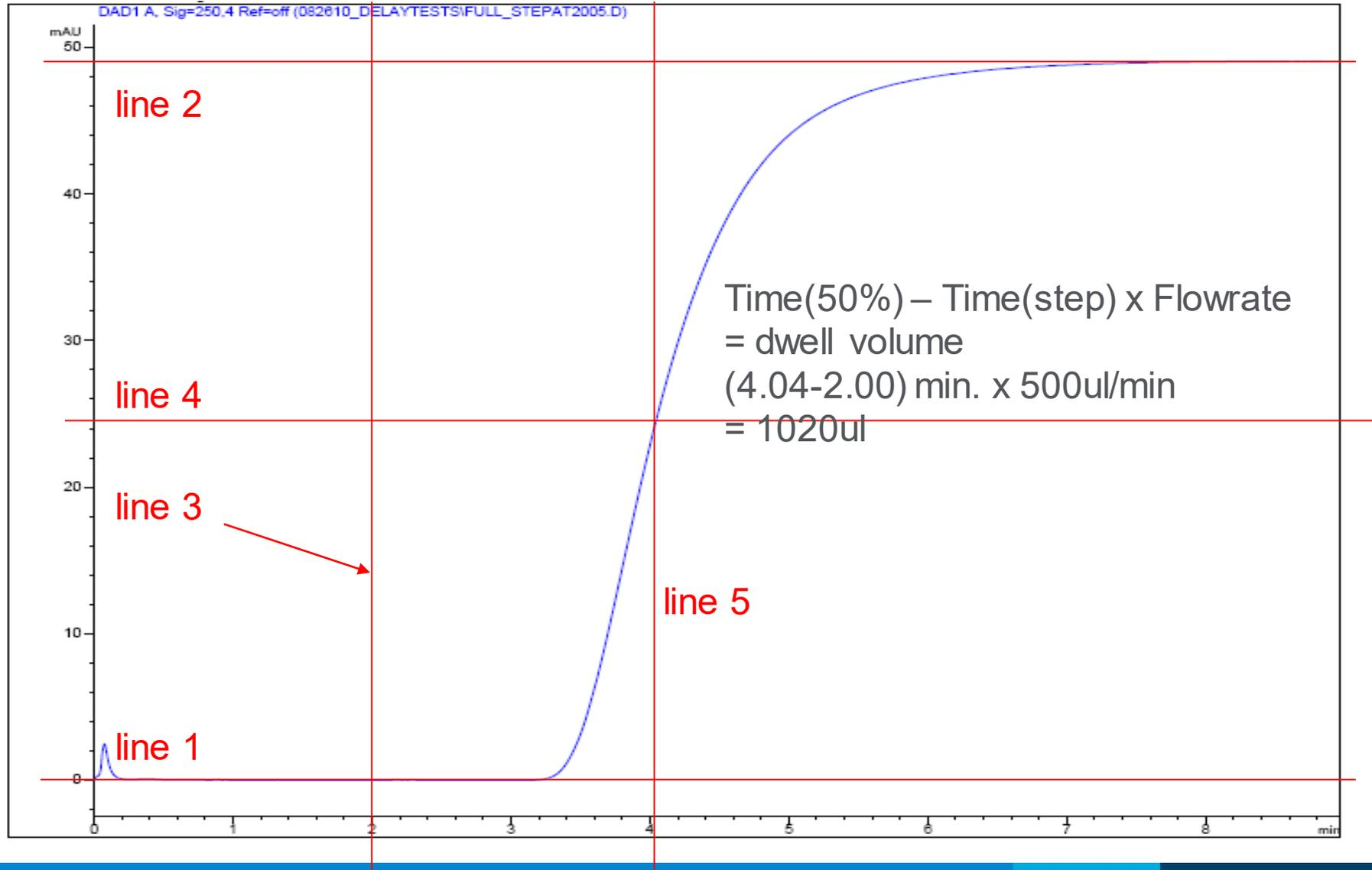


Total delay volume of the system (sum of capillaries, mixer, cells, valves..)

Chromatographic Test Result; Different Delay Volumes



Example -- How to Calculate Delay Volume



Summary of Available Data

Delay Volume Summary for Agilent Infinity 1290 Systems

Pump	Configuration	Sampler	Liftoff	Midpoint	Transition
1290 bin @ 350 Bar	35ul JetWeaver	G4226A	125	225	130
1290 quat @ 350 Bar	no mixer	G4226A	375	570	330
1290 quat @ 350 Bar	380ul JetWeaver	G4226A	450	820	820
For Reference...					
1260 bin @ 350 Bar	as shipped	G1367E	720	1105	545

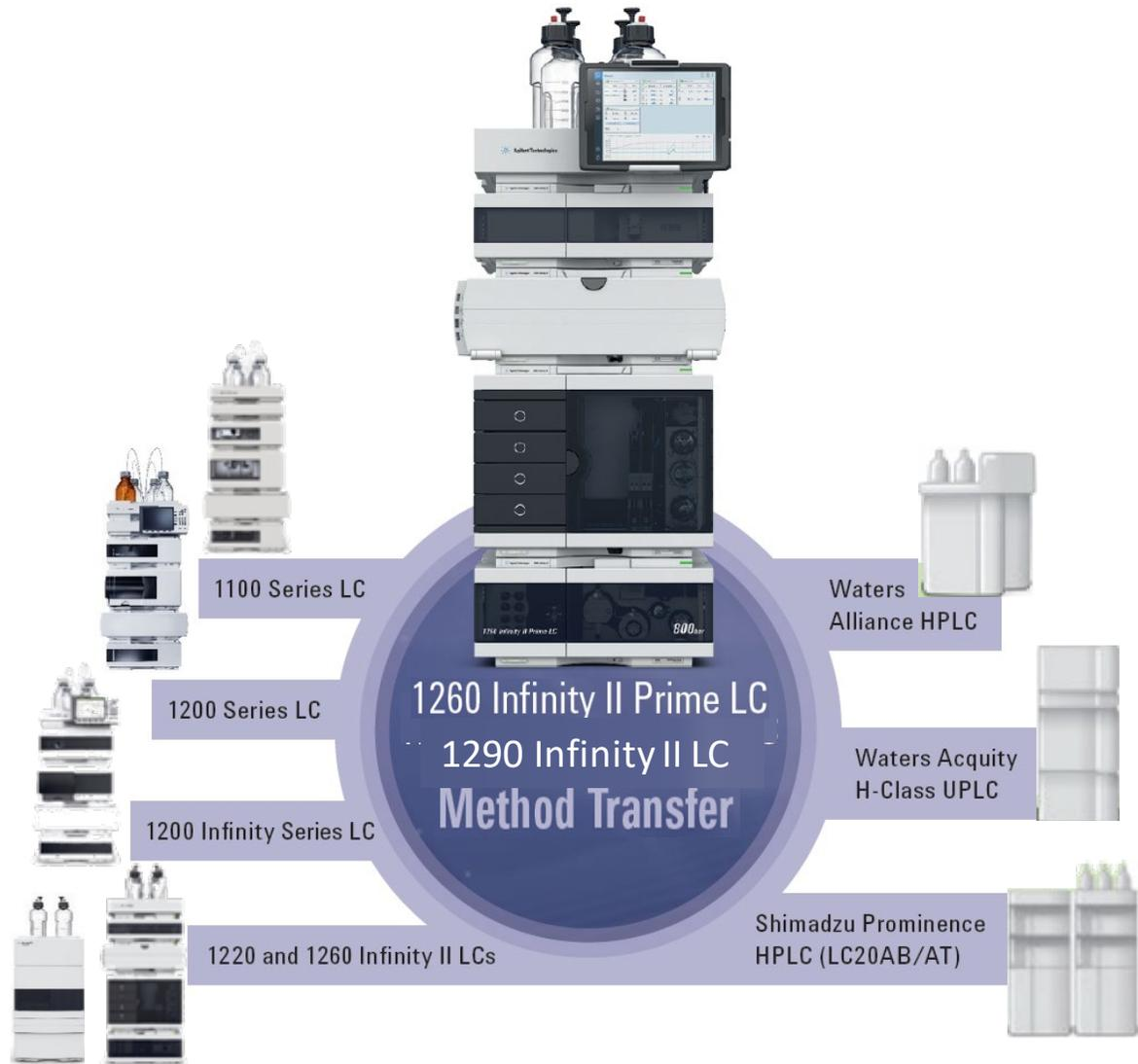
Summary of Available Data

Delay Volume Summary for Binary 1260 Systems

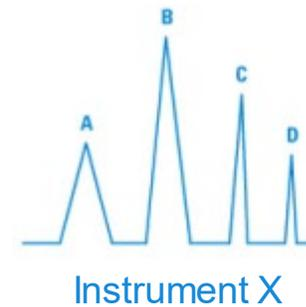
Pump	Configuration	Sampler	Liftoff	Midpoint	Transition
1260 bin @ 350 Bar	as shipped	G1367E	720	1105	545
1260 bin @ 350 Bar	minus damper	G1367E	635	955	460
1260 bin @ 350 Bar	minus damper with purge valve in channel A	G1367E	550	835	425
1260 bin @ 350 Bar	minus mixer/damper with purge valve in channel A	G1367E	270	375	175
1260 bin @ 350 Bar	minus mixer/damper with purge valve in channel A	G1367E with bypass (ADVR)		110	
For Reference					
1290 bin @ 350 Bar	35ul JetWeaver	G4226A	125	225	130

Intelligent System Emulation Technology (ISET)

One-Click LC System Emulation

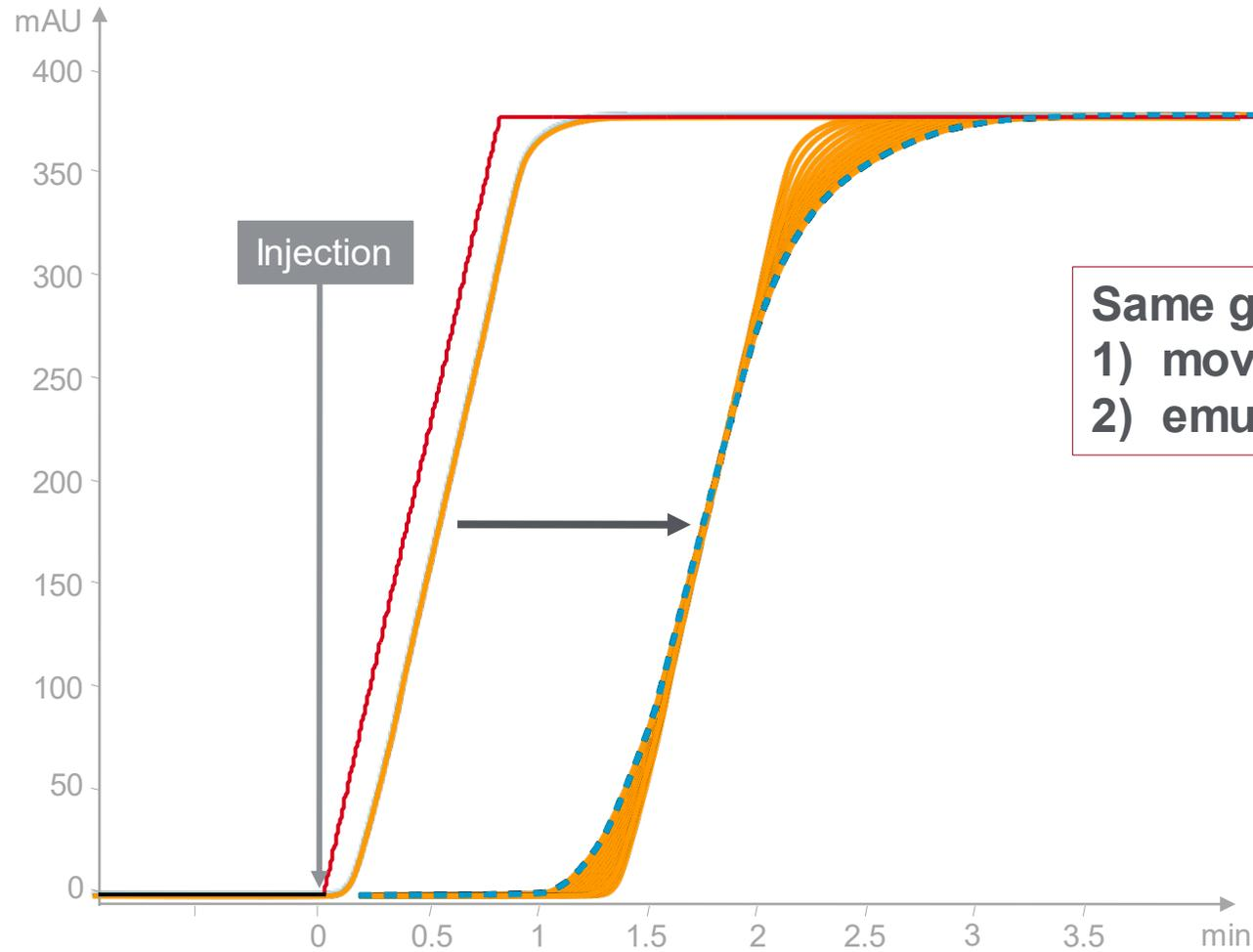


- Emulates other (U)HPLC instruments – by a simple mouse click
- Runs existing (U)HPLC methods – without modifying method or system
- Delivers same retention times and peak resolution – for infinitely better method transfer



Intelligent System Emulation Technology (ISET)

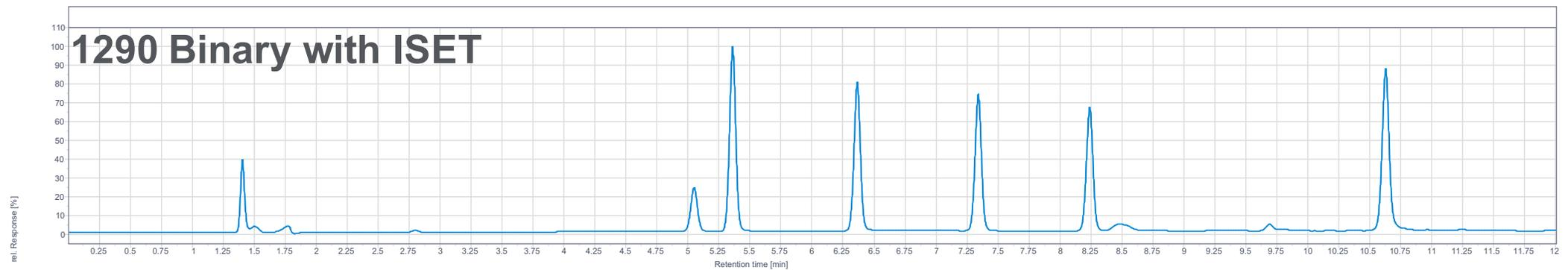
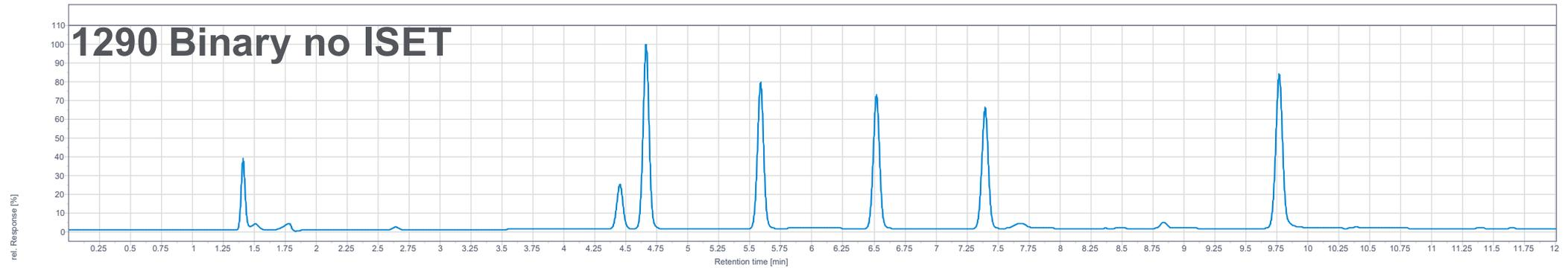
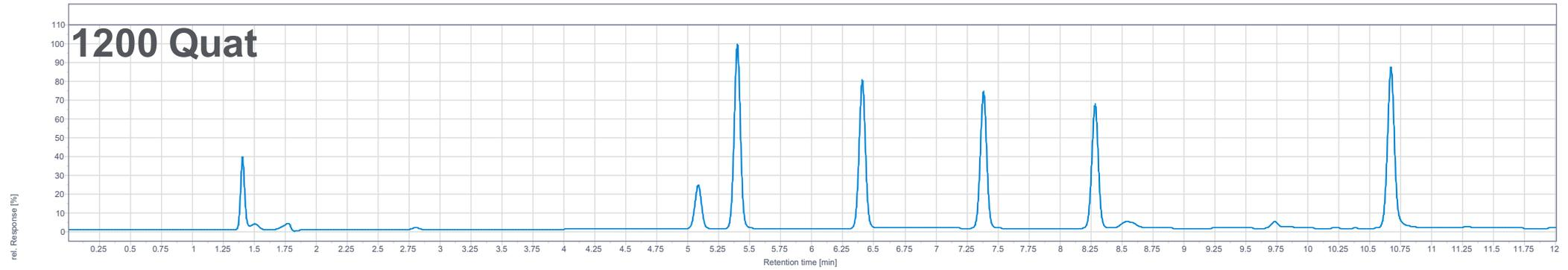
How does it work?

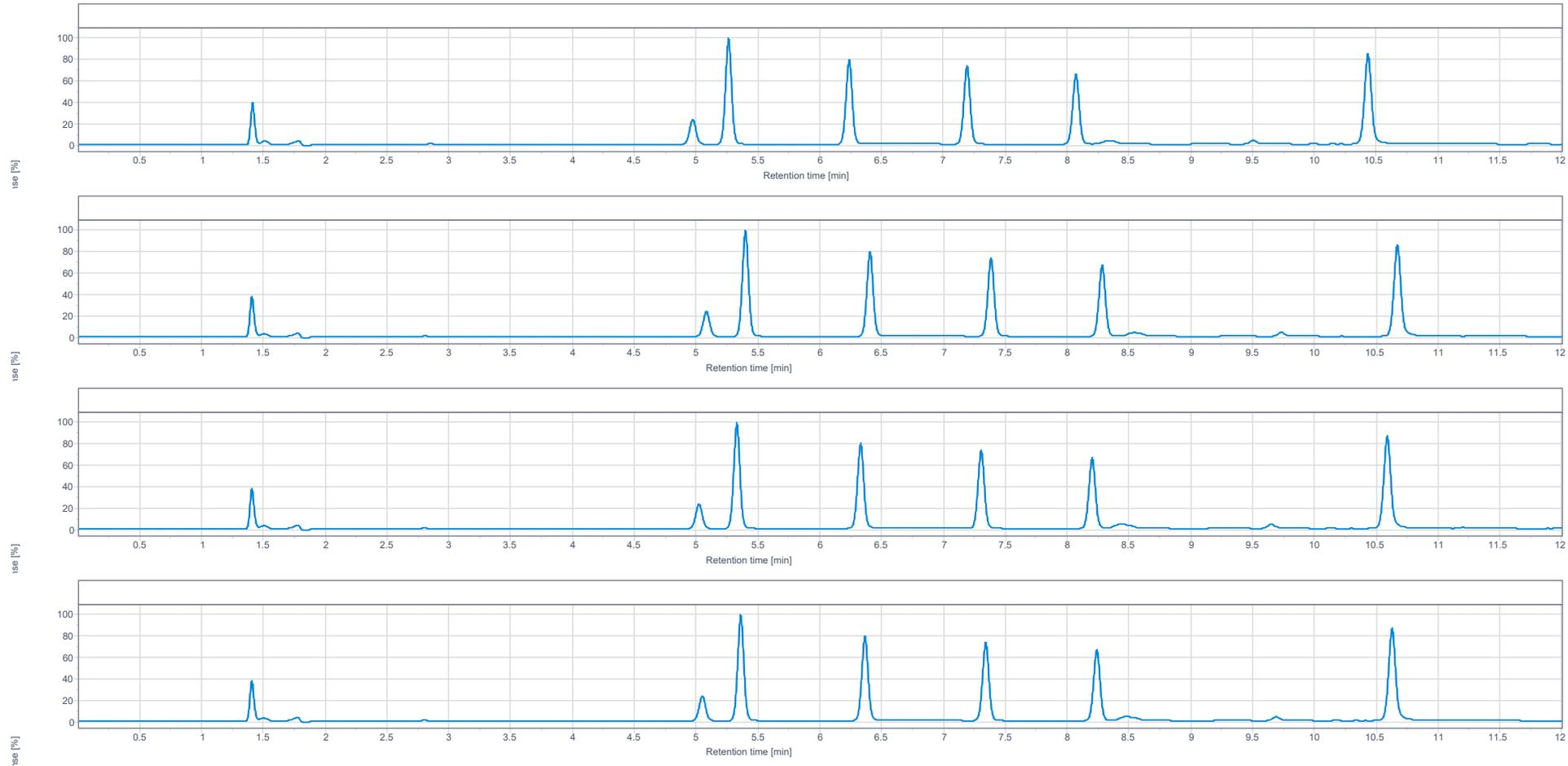


Same gradient conditions by
1) moving the gradient
2) emulating the gradient mixing behavior



Intelligent System Emulation Technology (ISET)





Can you tell which is which?
 You can run the fastest UHPLC methods, and still run your legacy methods

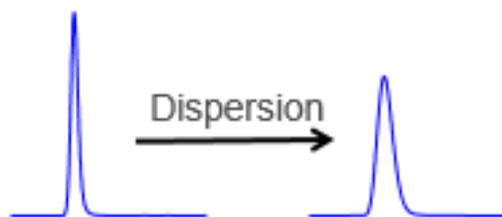
Lowest dispersion for highest resolution

Agilent **A-Line Quick Connect UHPLC column fittings** for truly dead-volume-free fluidic connections

- Tool-free connection up to 1300 bar
- Spring loaded design for easy and **truly zero-dead volume** connection
- Removable and reusable for all column types

System dispersion

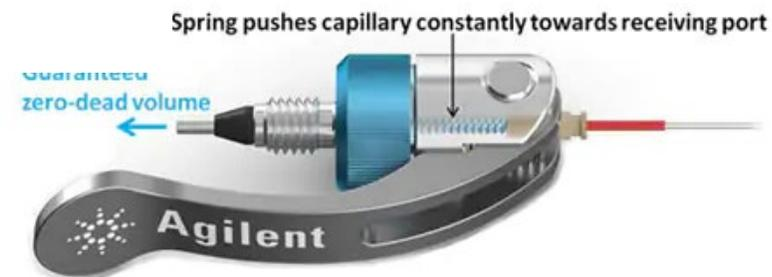
- *“Dispersion is the sample bandspreading or dilution which occurs in connecting tubing, sample valves, flow cells and in column end-fittings.”*



Peak height: Loss of sensitivity

Peak width: Loss of resolution

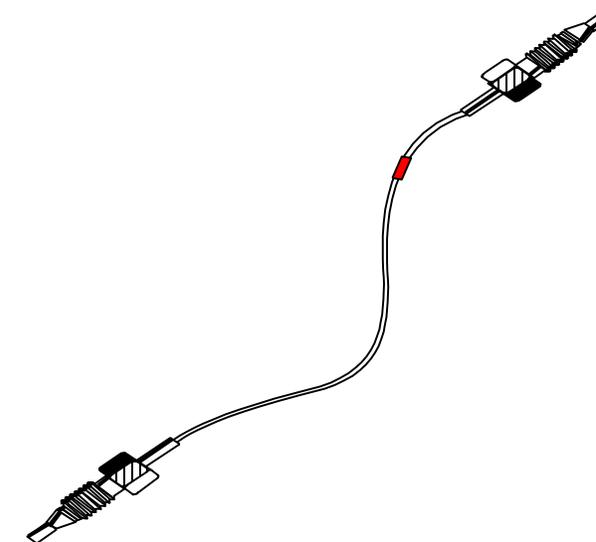
- Capillaries (Inner diameter, length)



Tubing Pressure and Internal Volume

Parameter	Black	Red	Green	Blue
Tubing Description	1	2	3	4
i.d. (mm)	0.075	0.127	0.178	0.254
Total Length (mm)	500	500	500	500
Flow rate (ml/min)	1.000	1.00	1.00	1.00
Solvent Name	ACN/wa	ACN/wa	ACN/wa	ACN/wa
Viscosity (cP)	1.2	1.2	1.2	1.2
Internal Volume (ul)	2.209	6.334	12.4	25.3
Expected pressure (bar)	129.1	15.7	4.1	1.0

<u>Color</u>	<u>i.d.</u>
Black	0.075 mm (0.003 inches)
Red	0.12 mm (0.005 inches)
Green	0.17 mm (0.007 inches)
Blue	0.25 mm (0.01 inches)
Clear	0.50 mm (0.02 inches)



500mm – typical sample flow path with UV detn. LC/MS -- near 1000mm

Tubing Dimensions and System Dispersion

	HPLC 1260		UHPLC 1290	
Column to Detector	18-30cm 0.17 18-30cm 0.12		18-30cm 0.12 25cm 0.075	
TCC outlet to column	10cm 0.17 10cm 0.12		10cm 0.12 10cm 0.075	
ALS to TCC inlet	18-30cm 0.17 18-30cm 0.12		18-30cm 0.12 45cm 0.075	
Needle Seat	Green 0.17 - 2.3uL Red 0.12 - 1.7uL		Red 0.12 - 1.7uL Black 0.075 - 0.9uL	
Pump to ALS	20-40cm 0.17 20-40cm 0.12		20-40cm 0.12 20cm 0.075	

https://www.agilent.com/cs/library/usermanuals/public/Copy%20of%20UltraLowDispersionKit_TN_EN.pdf

Optimize Dispersion for Small Columns

2.1 x 50 mm Poroshell 120 1.9 μ m

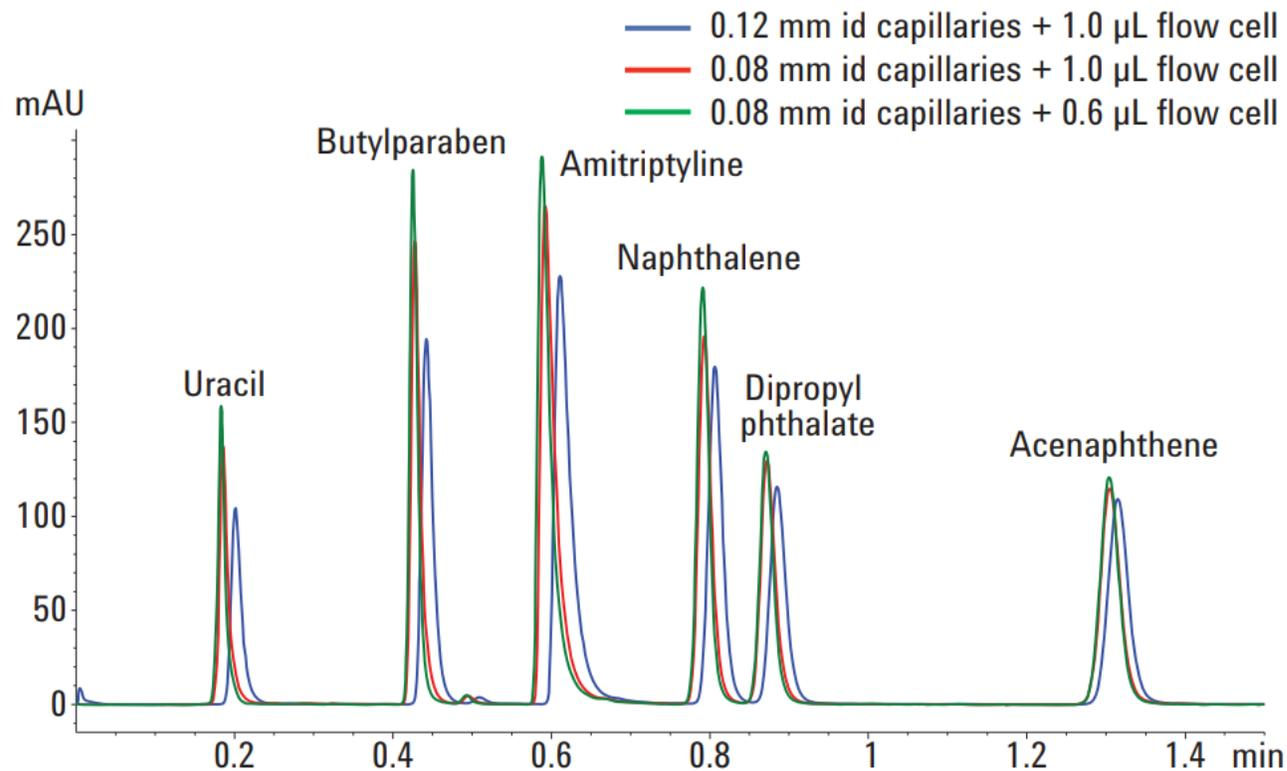


Figure 1A. The performance of an Agilent InfinityLab Poroshell 1.9 μ m column is improved when LC system volume is reduced by using smaller internal diameter capillaries and a smaller volume detector flow cell.

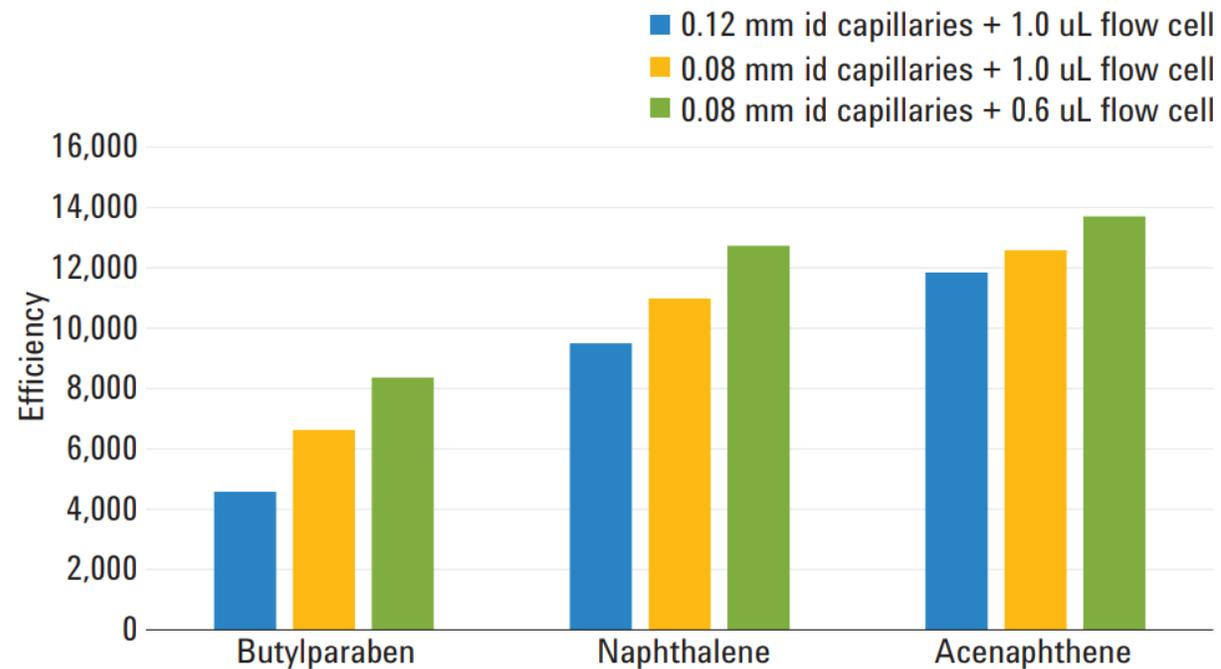
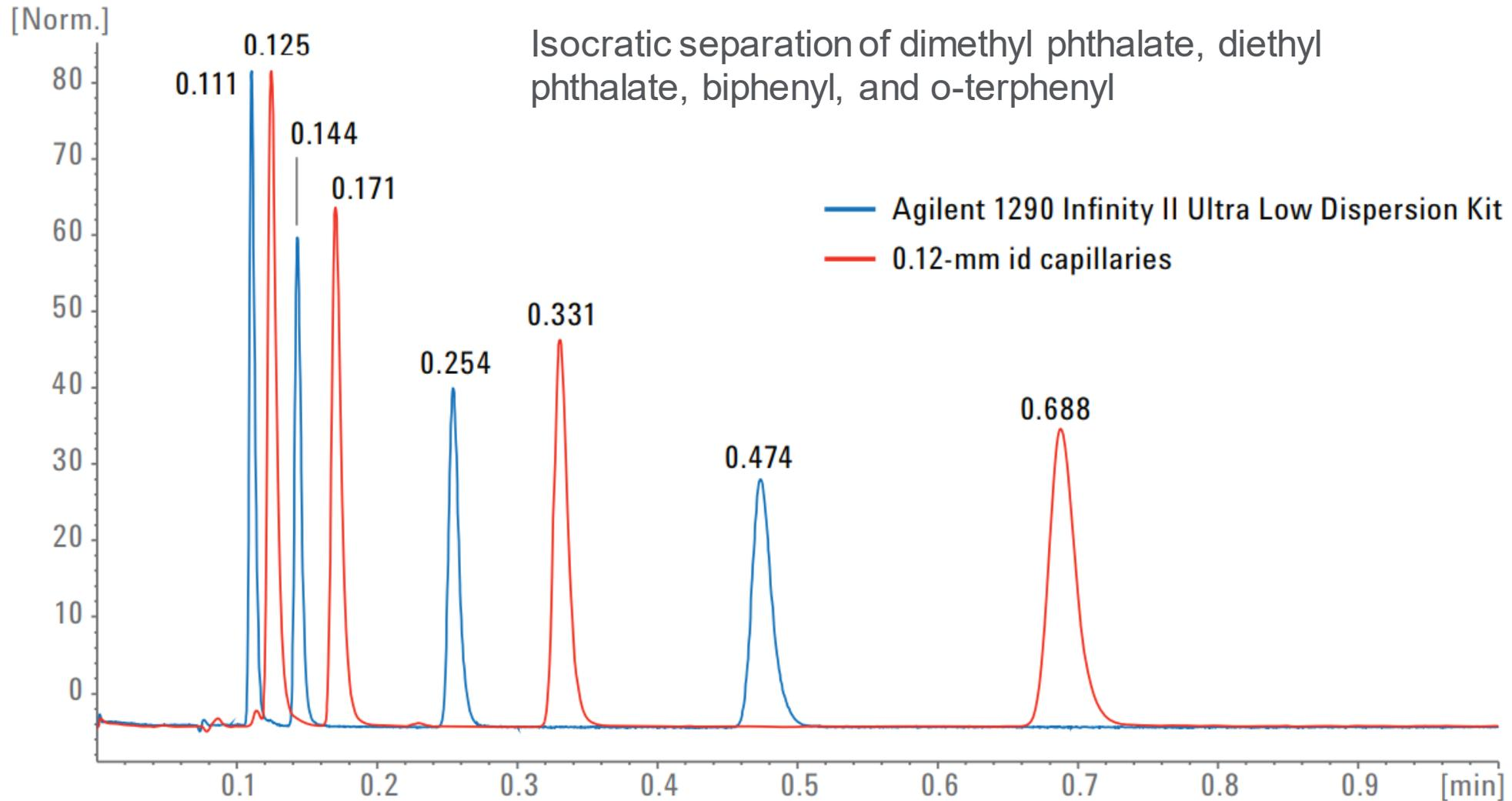


Figure 1B. The performance of an Agilent InfinityLab Poroshell 1.9 μ m column is improved when LC system volume is reduced by using smaller internal diameter capillaries and a smaller volume detector flow cell.

Optimizing System Dispersion on the Agilent 1290 Infinity II LC



Increased Pressure Requirements (column particle size)

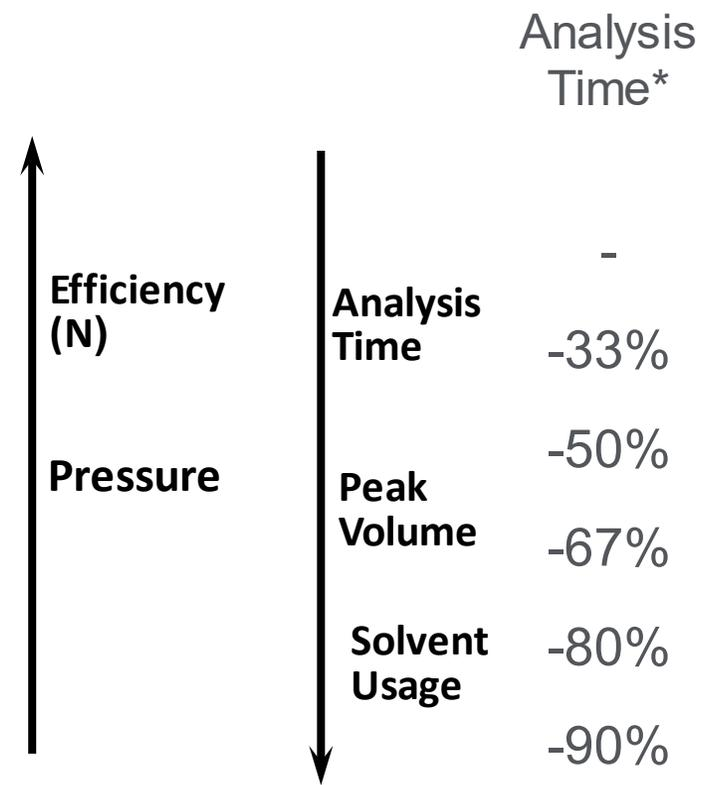
d_p (μm)	ΔP (bar)	N
5.0	14.5	25,000
3.0	66.9	41,000
1.5	531	83,000
1.0	1800	125,000
0.75	4270	166,000
0.50	14400	250,000

For a 30 cm column, 5.0 to 0.50 μm particle size reduction:

- N (theoretical plates) increases 10-fold
- ΔP (change in pressure) increases 1000-fold

Column Length and Particle Size

Column Length (mm)	Column Efficiency N(5 µm)	Column Efficiency N(3.5 µm)	Column Efficiency N(1.8 µm)
150	12,500	21,000	35,000
100	8,500	14,000	23,250
75	6000	10,500	17,500
50	4,200	7,000	12,000
30	N.A.	4,200	6,500
15	N.A.	2,100	2,500



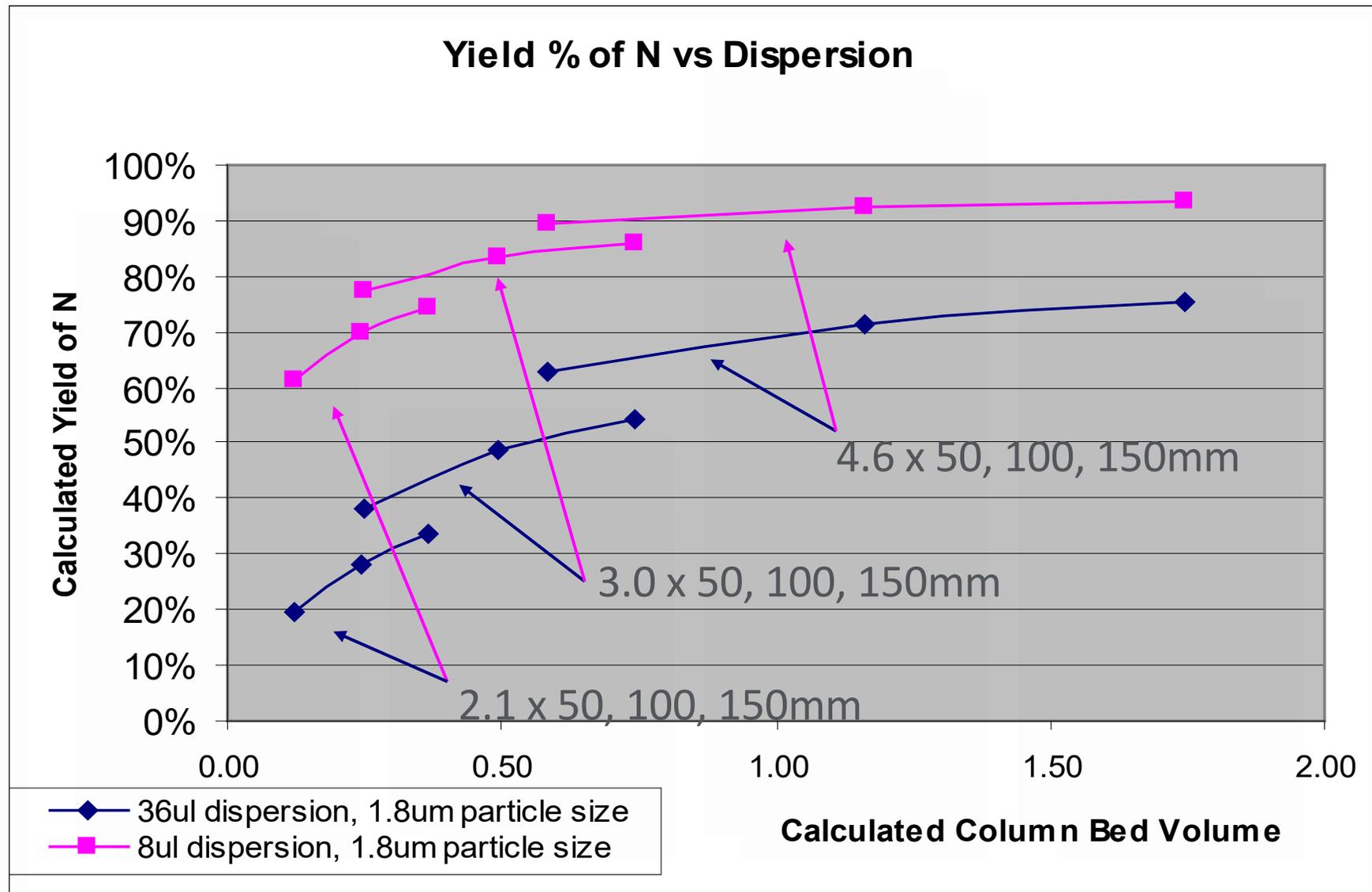
* Reduction in analysis time compared to 150 mm column; all columns 4.6-mm i.d.

Predicting Column Pressure and Performance

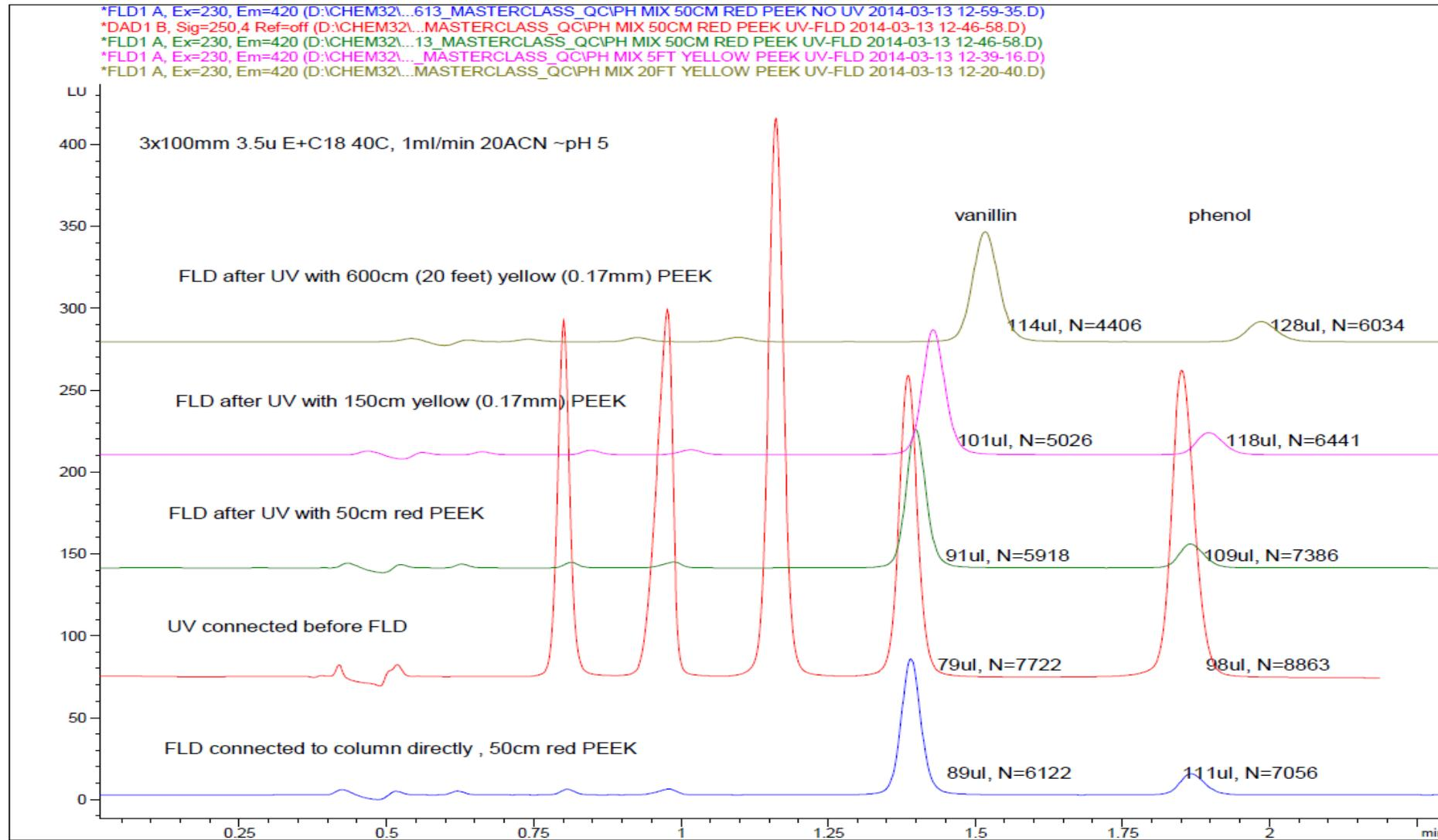
Parameter		Media	ZORBAX			Poroshell 120	
i.d. (mm)			2.100	3.000	3.000	3.000	3.000
Length (mm)			50.0	50.0	100.0	100.0	100.0
Particle Size (µM)			1.8	1.8	1.8	1.9	2.7
Column void fraction			0.60	0.60	0.60	0.60	0.60
Permeability (PF)			900	900	900	900	900
Reduced plate height value			2.30	2.30	2.30	1.70	1.70
Flow rate (ml/min)			0.5000	1.0000	1.000	1.000	1.000
Linear Velocity (mm/sec)			4.01	3.93	3.93	3.93	3.93
Viscosity (cP) 25C ACN/Water			1.2	1.2	1.2	1.2	1.2
Column Volume (ml)			0.104	0.212	0.424	0.424	0.424
Expected pressure (bar)			401.0	393.0	785.9	705.4	349.3
Estimated Efficiency** (Neue)			12077	12077	24155	30960	21786

**Estimates do not consider system dispersion effects. Actual results will vary.

Yield estimates – 1.8um particles 2.1-4.6mm i.d.



Multi-Detector Connections Bandspreading



Use The Right Tool for the Job



We might get this to work, but the results may not be what we needed

(U)HPLC Systems

- Each component of the system has settable values, however not all settable values can be achieved
- Before using or purchasing a system, the user must be certain that the desired parameters can be run accurately and reproducibly
- Instrument specifications should be obtained from the appropriate manuals
- From these specifications, it can be determined if that component may be used

Pump Parameters

The screenshot displays the 'Flo' (Flow) parameters section, including solvent composition (A: 95.0% Water, B: 5.0% Acetonitrile), pressure limits (Min: 0.00 ba, Max: 1300.00 ba), and injection options. The 'Advanced' section shows a 'Timetable (2/100 events)' table with the following data:

Time [min]	A [%]	B [%]	Flow [mL/min]	Max. Pressur Limit [bar]
0.00	95.00	5.00	0.200	1300.00
3.00	95.00	5.00	---	---
100.00	40.00	60.00	---	---

Annotations in the image include:

- A red box around the flow rate '0.200 mL/min' with an arrow pointing to the text 'Note flow rate 0.2 ml/min'.
- A red box around the '40.00' value in the B [%] column of the timetable, with an arrow pointing to the text 'Note gradient change (%B/min)'.

- Low flow rate (<0.5 ml/min) start to move to binary pump due lower delay volumes
- Shallow gradient (<1%B/min) requires binary pump due to lower delay volume and more precise changes

AutoSampler Parameters

Injecti

Injection volu 0.50 µ

Needle W

Multi-wash

Stopti Postti

As Pump/No Li Of

1.00 mi 1.00 mi

Advanced

Injection Path Cleaning

Standard W

Mode: Wash Vial

Time: 3 s

Locati

Repea 3

Multi-wa

Step	Solvent	Time [s]	Seat Back FI	Needle W	Comment
1	S1	0	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	
2	S2	0	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	
3	S3	0	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	
Start C...	S1		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	

Newer, UHPLC methods may have lower injection volumes

'Sticky samples or more sensitive detectors may require aggressive cleaning

- Shorter (<150mm) and narrower columns (<3mm) may have lower injection volumes. Make sure the sampler precision and accuracy are correct for low volumes
- More sensitive detectors (or higher salt concentrations in the buffers) may require additional wash steps for both the needle and needle seat

Detector Parameters

Signal

Acquire	Wavelength	Bandwidth	Reference Wavelength	Reference Bandwidth
Signal A	254.0	4.0	360.0	100.0
Signal B	210.0	4.0	360.0	100.0
Signal C	214.0	4.0	360.0	100.0
Signal D	230.0	4.0	360.0	100.0
Signal E	260.0	4.0	360.0	100.0
Signal F	273.0	4.0	360.0	100.0
Signal G	280.0	4.0	360.0	100.0
Signal H	250.0	4.0	360.0	100.0

Advanced

Spectru

Store: All

Range from: 190.0 t 600.0 n

Step: 20 n

Analog Out

Zero Offset: 5 %

Attenuation: 1000 mA

Margin for negative Abso

100 mA

Autobala

Lamps on required for acq

Prerun UV Lamp

Postrun

Stopti

- > 0.013 min (0.25 s response time) (20)
- < 0.0016 min (0.016 s response time) (12...
- > 0.0016 min (0.031 s response time) (12...
- > 0.0031 min (0.063 s response time) (80...
- > 0.0063 min (0.13 s response time) (40...
- > 0.013 min (0.25 s response time) (20 H...
- > 0.025 min (0.5 s response time) (10 Hz)
- > 0.05 min (1 s response time) (5 Hz)
- > 0.1 min (2 s response time) (2.5 Hz)
- > 0.2 min (4 s response time) (1.25 Hz)
- > 0.4 min (8 s response time) (0.62 Hz)
- > 0.85 min (16 s response time) (0.31 Hz)

As columns become shorter peaks elute more quickly requiring detector data rates to increase.

Resources — Primers

[5990-7595EN](#)

The LC Handbook

Guide to LC Columns and Method Development

[5991-2359EN](#)

Two Dimensional Liquid Chromatography

[5990-3777EN](#)

High Performance Capillary Electrophoresis

[5991-5509EN](#)

Supercritical Fluid Chromatography

[5989-6639EN](#)

Principles in Preparative HPLC

[5991-3326EN](#)

Sample Preparation Fundamentals for Chromatography

[5980-1397EN](#)

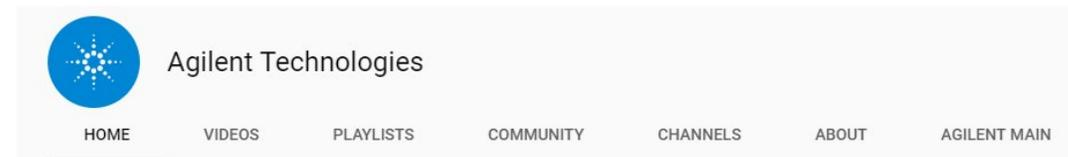
Fundamentals of UV-visible Spectroscopy



Resources for Support

- Collection of LC resources:
https://community.agilent.com/docs/DOC-1852-lc-insights-to-go#jive_content_id_LC_Troubleshooting
- Agilent support resources:
<https://community.agilent.com/community/resources>
- Agilent University: <http://www.agilent.com/crosslab/university>
- Agilent resource center:
<http://www.agilent.com/chem/agilentresources>
- InfinityLab Supplies Catalog ([5991-8031EN](#))
- Your local FSE and Specialists
- Youtube – [Agilent Channel](#)

- Sales and support phone assistance (US and Canada):
1-800-227-9770 [Phone Tree Navigation Assistance](#)



The Chemical Standards Song

6,089 views • 3 v

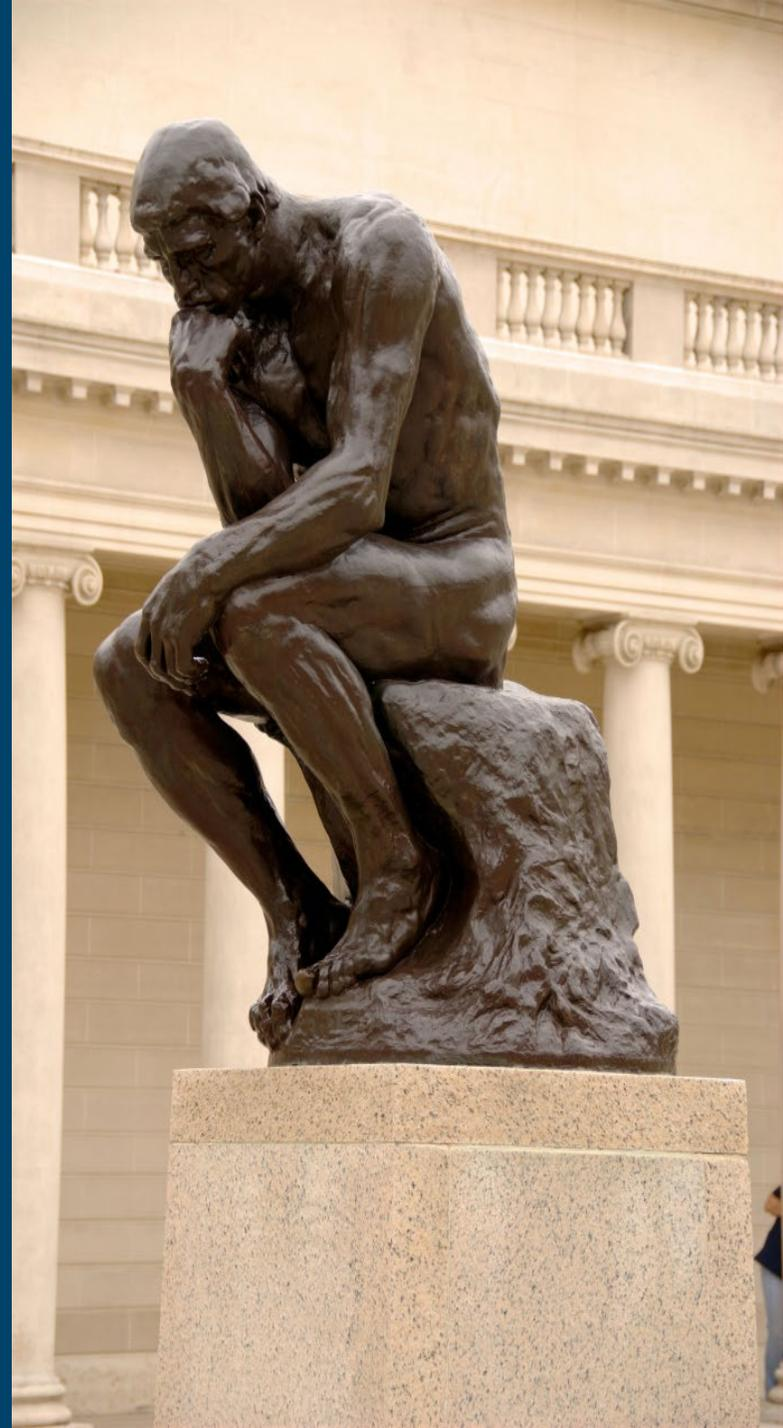
<http://www.agilent.com>
When you calibrate
great.
#standardsrule



gc-column-support@agilent.com
lc-column-support@agilent.com
spp-support@agilent.com
spectro-supplies-support@agilent.com

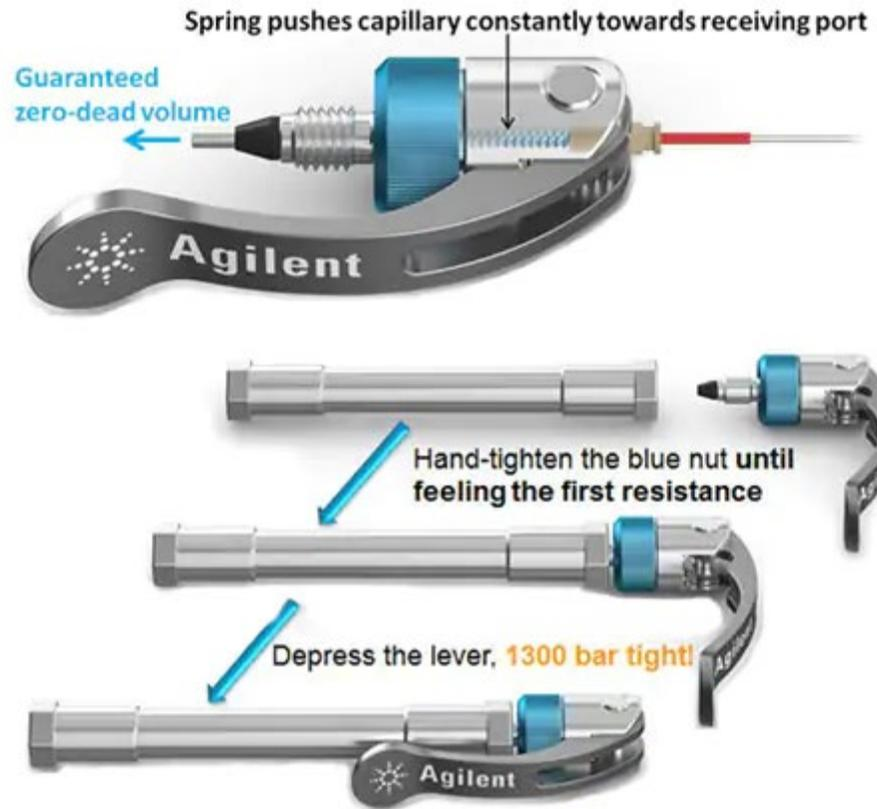


Thanks for your
attention!



Universal Low Dispersion Fittings

Quick Connect



Quick Turn



Agilent application note 5991-5525EN