

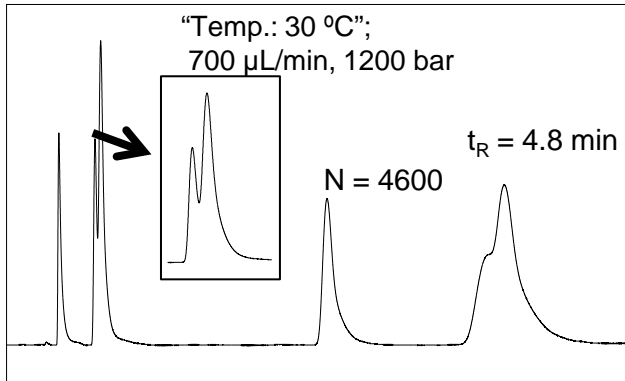


From HPLC to UHPLC: Method Transfer, What Should I Pay Attention to?

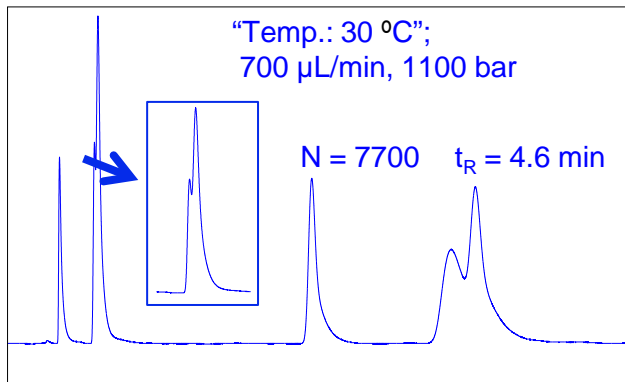
Dr. Frank Steiner
Thermo Fisher Scientific, Germering/Germany

A Well-Known Situation in the Lab

- Why does a chromatogram recorded on system X...

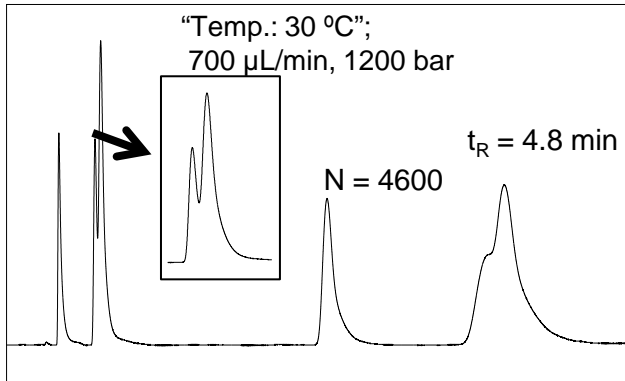


- ...look so different on system Y?

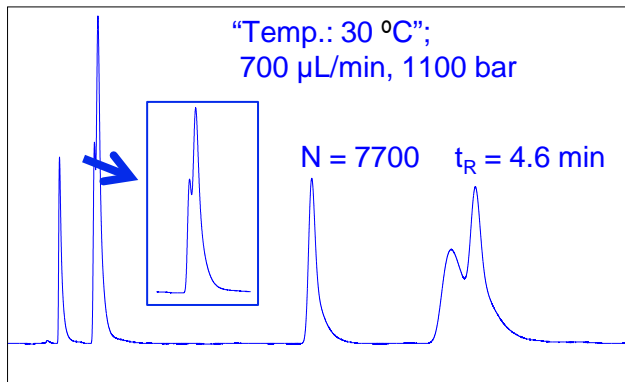


A Well-Known Situation in the Lab

- Why does a chromatogram recorded on system X...



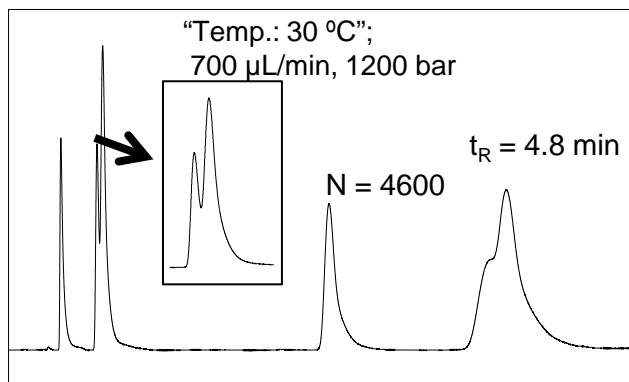
- ...look so different on system Y?



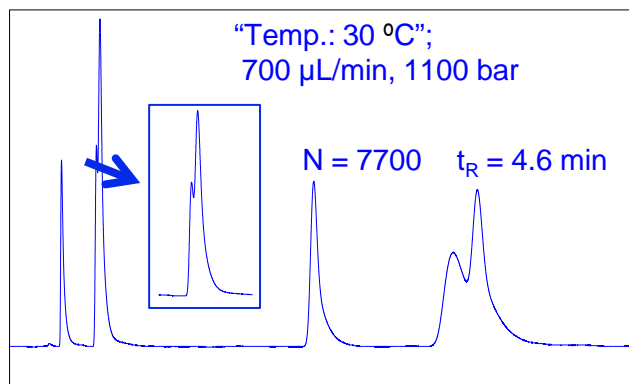
- Reasons are differences in the fluidics...
 - Gradient separations mostly affected
 - Impacts primarily retention times
 - Main reason: Gradient delay volume (GDV)

A Well-Known Situation in the Lab

- Why does a chromatogram recorded on system X...



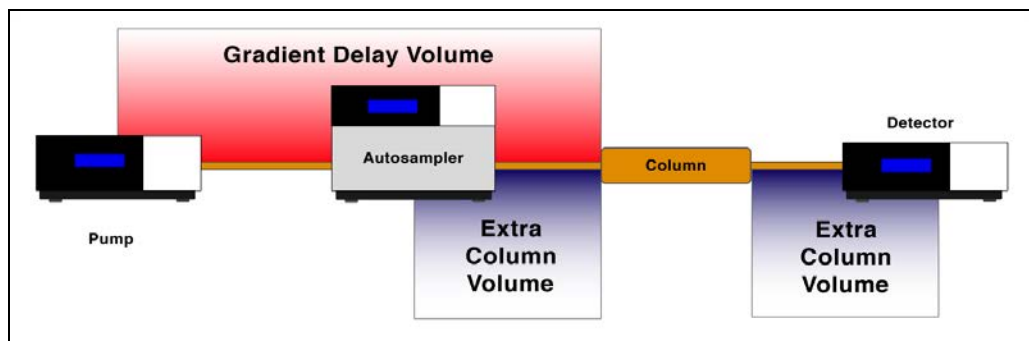
- ...look so different on system Y?



- Reasons are differences in the fluidics...
 - Gradient separations mostly affected
 - Impacts primarily retention times
 - Main reason: Gradient delay volume (GDV)
- ...and the thermostating concept
 - Affects both isocratic and gradient separations
 - Impacts retention times and peak shape (efficiency)
 - Reasons: Operational principle of the column thermostat, eluent pre-heating

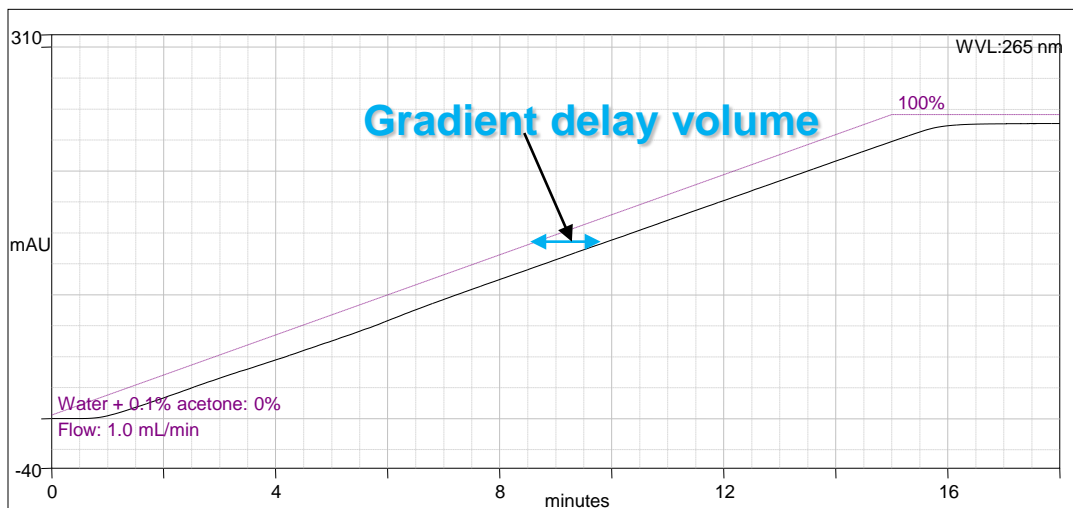
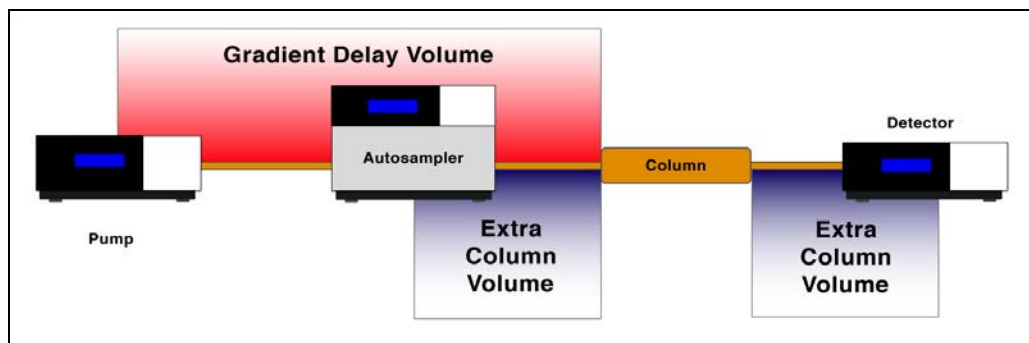
The Gradient Delay Volume or
even a Journey of a Thousand Miles
begins with a Single (Isocratic) Step

The Gradient Delay Volume (GDV)



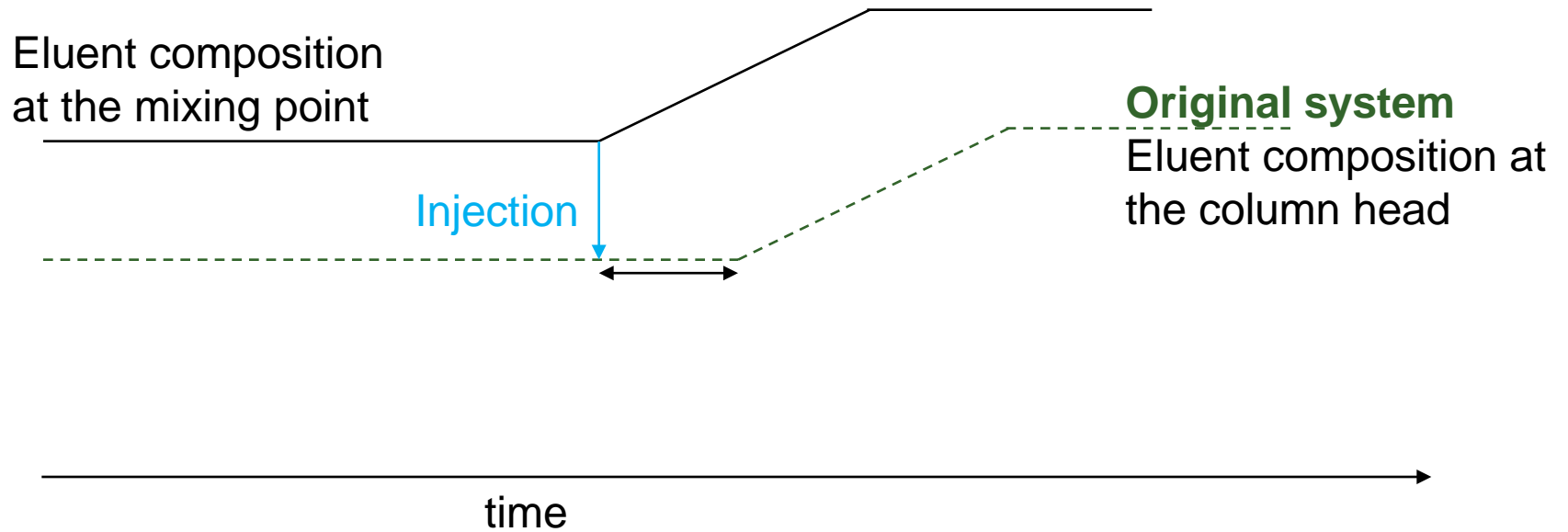
- GDV influences or generates
 - An isocratic step at the very beginning of every gradient separation
 - The gradient sharpness
 - The required equilibration time and thus the total analysis time
- Weakly retained compounds are much more affected by the GDV influence than late-eluting compounds.

The Gradient Delay Volume (GDV)

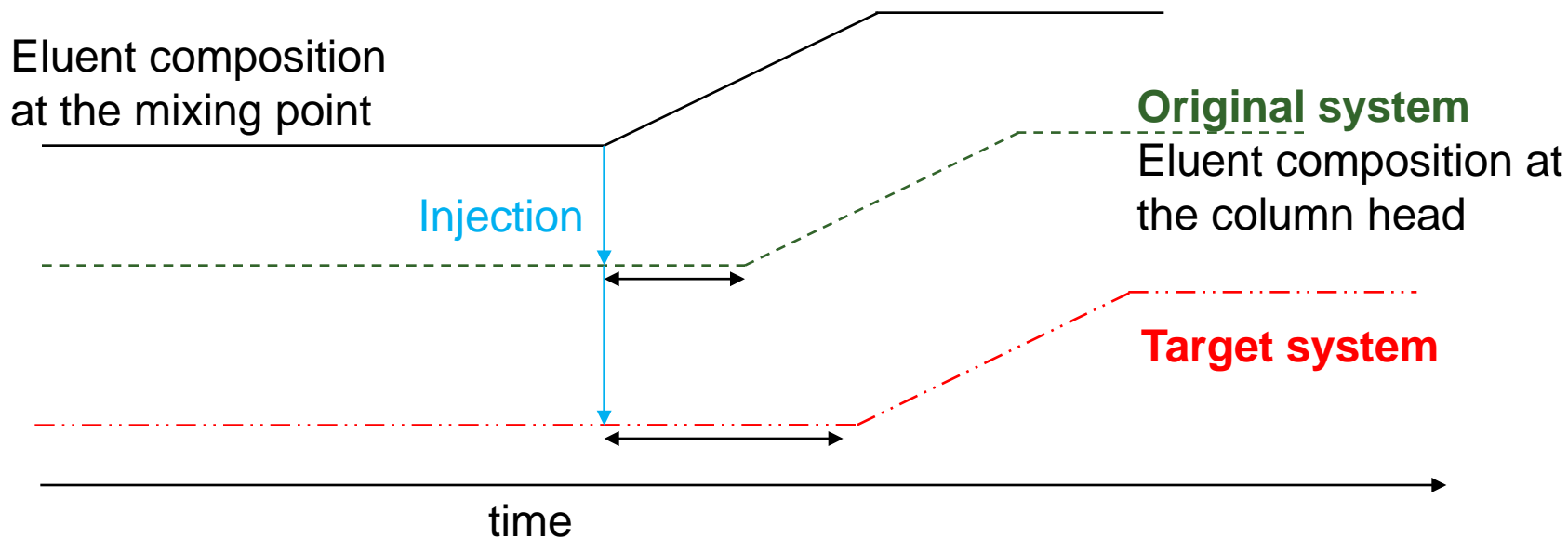


- GDV influences or generates
 - An isocratic step at the very beginning of every gradient separation
 - The gradient sharpness
 - The required equilibration time and thus the total analysis time
 - Weakly retained compounds are much more affected by the GDV influence than late-eluting compounds.
- ⇒ Needs special attention with steep/ballistic gradients and/or low flow rates

What Does the GDV Do with My Method Transfer?

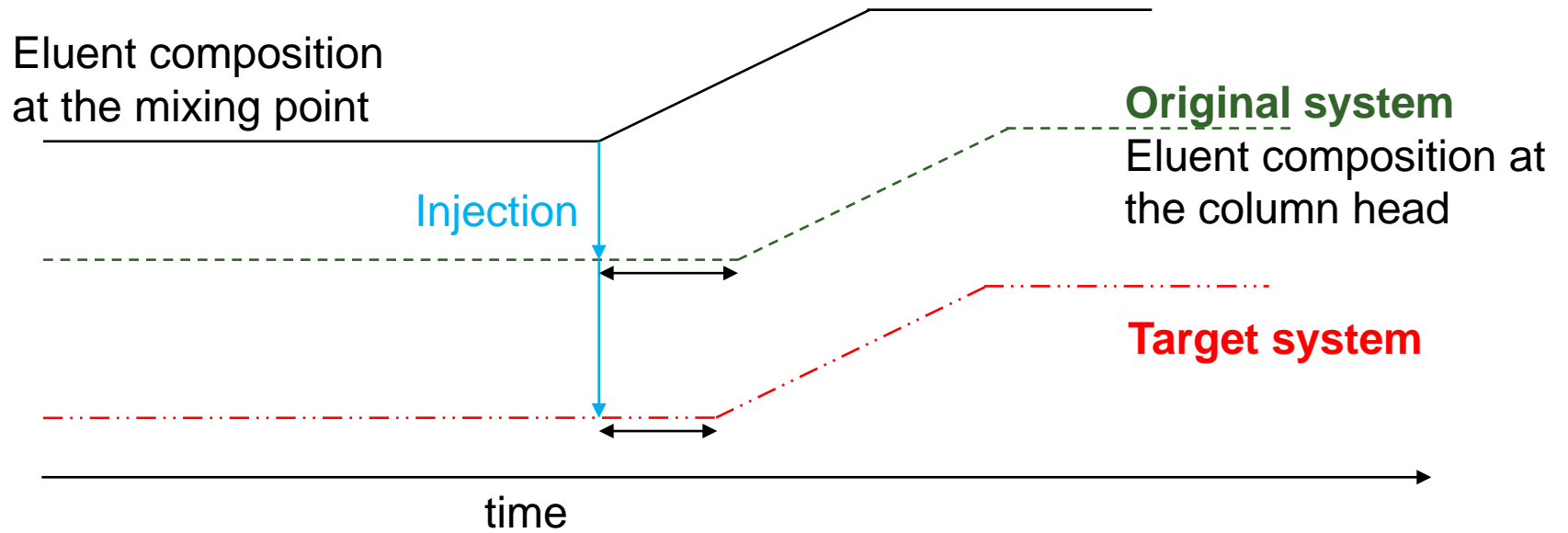


What Does the GDV Do with My Method Transfer?



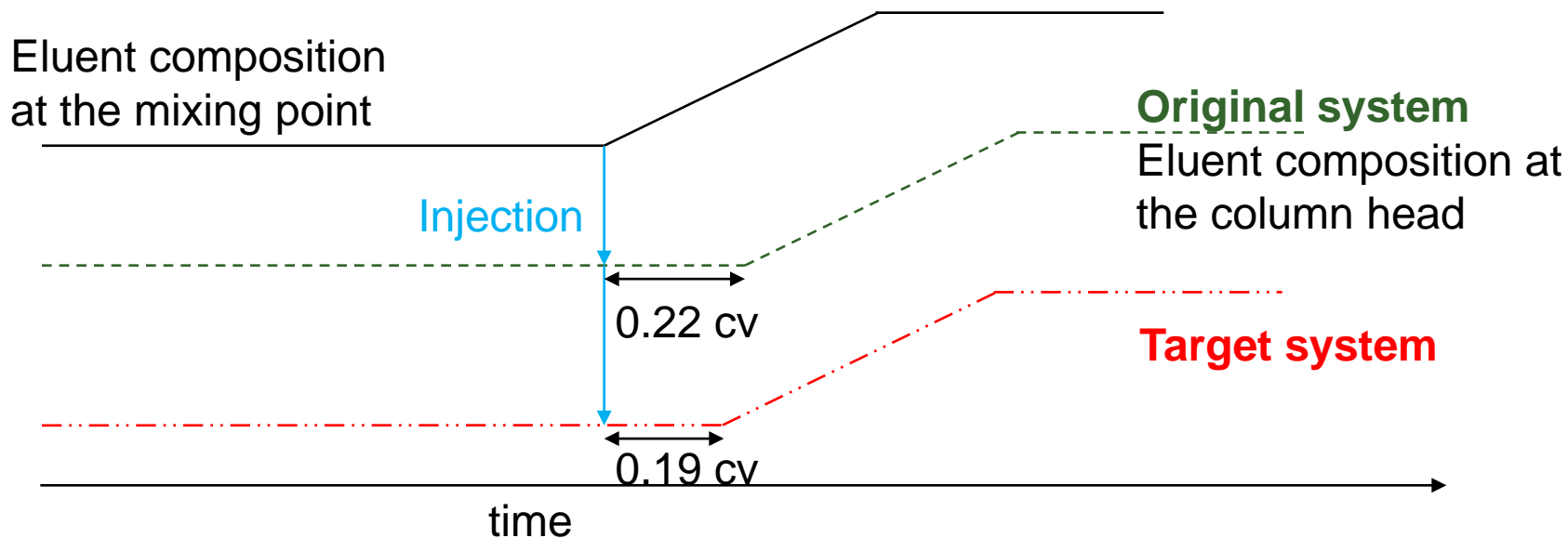
- Compare the GDVs (Represented as GDV per column volume)
- If the target system gives you a smaller isocratic segment:
 - Add an additional hold-up step before the gradient start to ensure identical isocratic start conditions
- If the target system gives you a larger isocratic segment:
 - Minimize the isocratic segment (E.g. with microflow kits, smaller tubing I.D., autosampler bypass)

What Does the GDV Do with My Method Transfer?



$$GDV_{per}(CV) = \frac{GDV}{CV}$$

What Does the GDV Do with My Method Transfer?



Original system with 0.9 mL GDV and 4.15 mL CV:

0.22 GDV per column volume

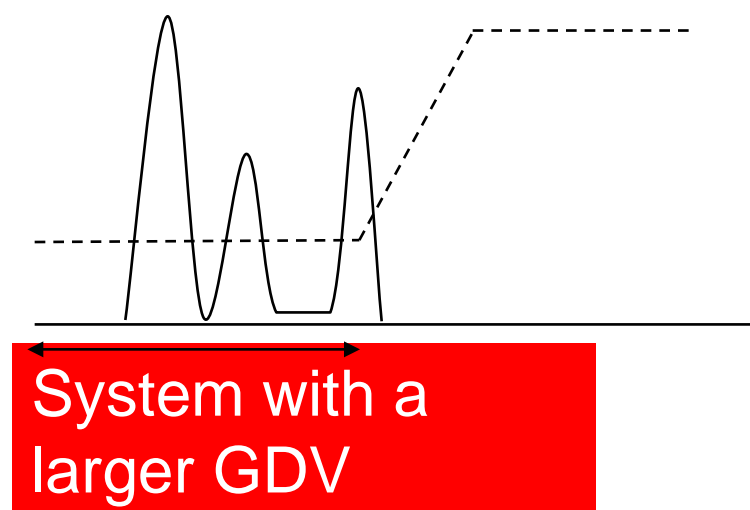
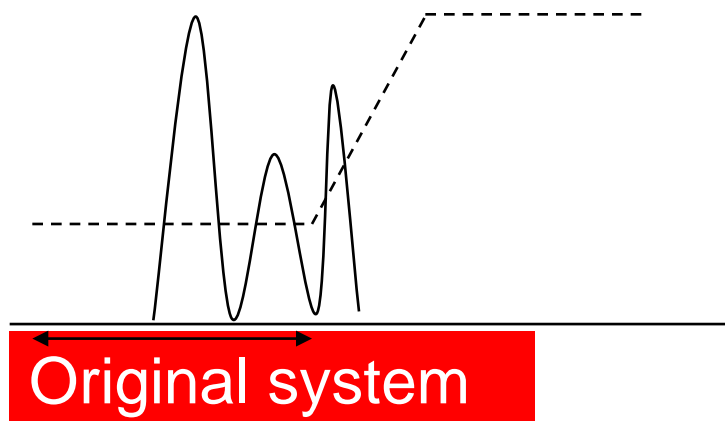
$$GDV_{per}(CV) = \frac{GDV}{CV}$$

Target system with 0.1 mL GDV and 0.52 mL CV:

0.19 GDV per column volume

- Change the GDV to column volume ratio as little as possible!

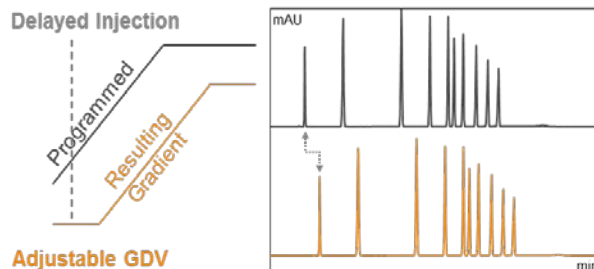
Different GDVs and their Influence on Early-Eluting Compounds



How a UHPLC System can Help to Handle GDV Challenges

- **Larger GDV required**

- Add additional volumes to the flow path:
 - Individually settable metering device in the autosampler to fine-tune GDV
 - Adjust mixer and (Thermo Scientific™ Viper™) capillaries
- Add isocratic hold-up step at the beginning of the separation



Chromleon



Pump H and F



Sampler HT and FT

Thermo Scientific™
Chromleon™ software



Thermo Scientific™
Vanquish™
Horizon
system



Thermo Scientific™
Vanquish™
Flex
system

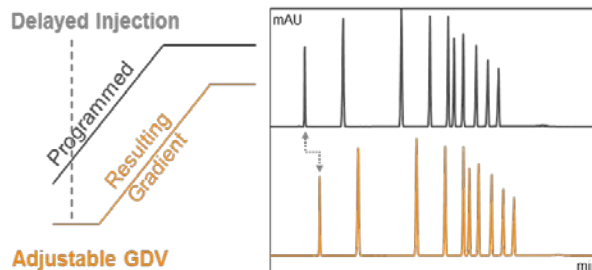


Thermo Scientific™
UltiMate™ 3000
(Bio)RS system

How a UHPLC system can help to handle GDV challenges

- **Larger GDV required**

- Add additional volumes to the flow path:
 - Individually settable metering device in the autosampler to fine-tune GDV
 - Adjust mixer and (Viper) capillaries
- Add isocratic hold-up step at the beginning of the separation



- **Smaller GDV required**

- Reduce volumes in the flow path:
 - Individually settable Metering Device in the autosampler to fine-tune GDV
 - Adjust mixer and (Viper) capillaries
- Delayed Injection at the program start



Chromleon



Pump H and F



Sampler HT and FT



Thermo Scientific™
Vanquish™
Flex
system



Thermo Scientific™
Vanquish™
Horizon
system



Thermo Scientific™
UltiMate™ 3000
(Bio)RS system

Thermostating with UHPLC or “Some like it Hot”

The Role of Temperature in UHPLC Separations

- Temperature is a key factor for selectivity optimization.
- Retention depends inversely on temperature:

$$\ln k = -\frac{\Delta H^0}{R} \cdot \frac{1}{T} + \frac{\Delta S^0}{R} + \ln \beta \quad (\text{Van't Hoff equation})$$

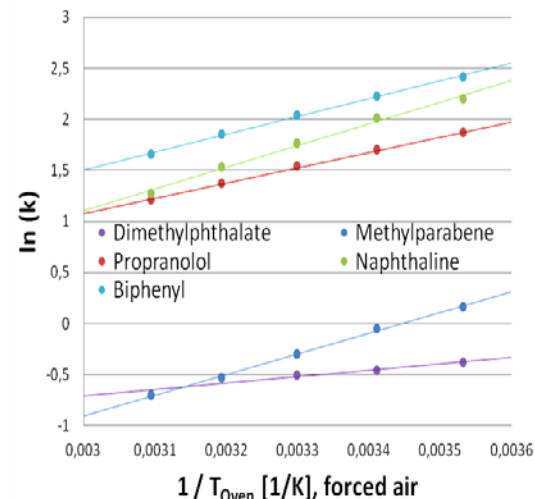
The Role of Temperature in UHPLC Separations

- Temperature is a key factor for selectivity optimization.
- Retention depends inversely on temperature:

$$\ln k = -\frac{\Delta H^0}{R} \cdot \frac{1}{T} + \frac{\Delta S^0}{R} + \ln \beta \quad (\text{Van't Hoff equation})$$

$$\ln k \sim \frac{1}{T}$$

- Even minor temperature variations can lead to changes in the elution order.



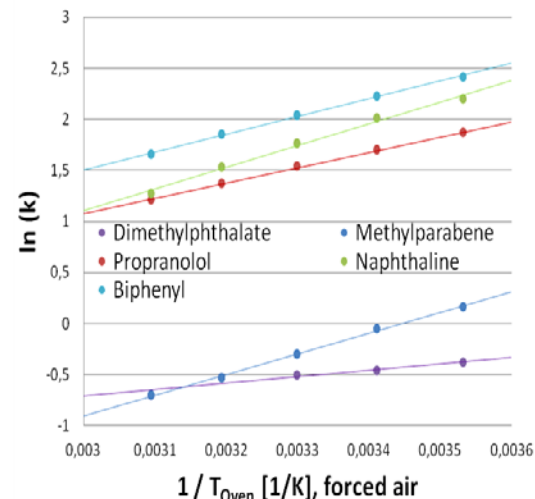
The Role of Temperature in UHPLC Separations

- Temperature is a key factor for selectivity optimization.
- Retention depends inversely on temperature:

$$\ln k = -\frac{\Delta H^0}{R} \cdot \frac{1}{T} + \frac{\Delta S^0}{R} + \ln \beta \quad (\text{Van't Hoff equation})$$

$$\ln k \sim \frac{1}{T}$$

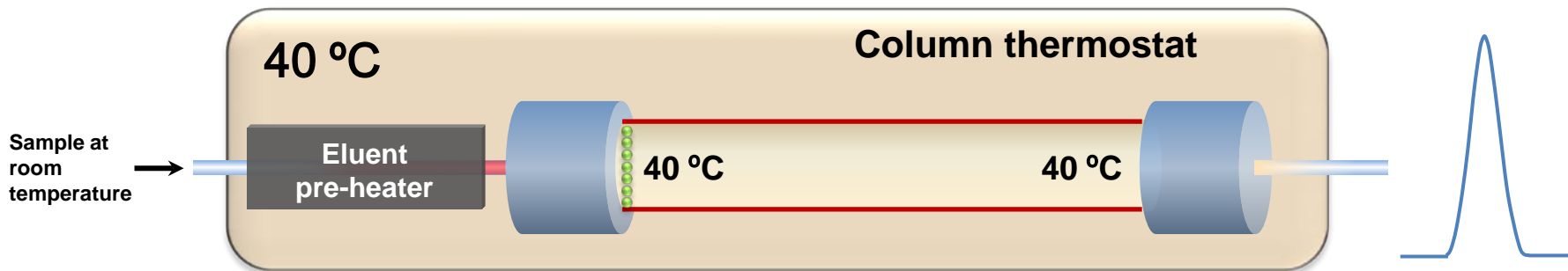
- Even minor temperature variations can lead to changes in the elution order.
- High pressure in UHPLC creates frictional heat inside the column.
- Two design concepts for column compartments:
 - Forced air thermostats – Continuous air circulation inside the column chamber
 - Still air thermostats – No air circulation with the goal to thermally insulate the separation column



- Still air thermostating always leads to better chromatographic resolution due to higher efficiencies / plate counts.
- Therefore, UHPLC methods which measurably increase the temperature inside the column by viscous friction always benefit from still air thermostating.
- Yes, but wait...:
 - This *can* be true, but it doesn't have to, if the rising internal column temperature results in a lower selectivity of critical peak pairs.

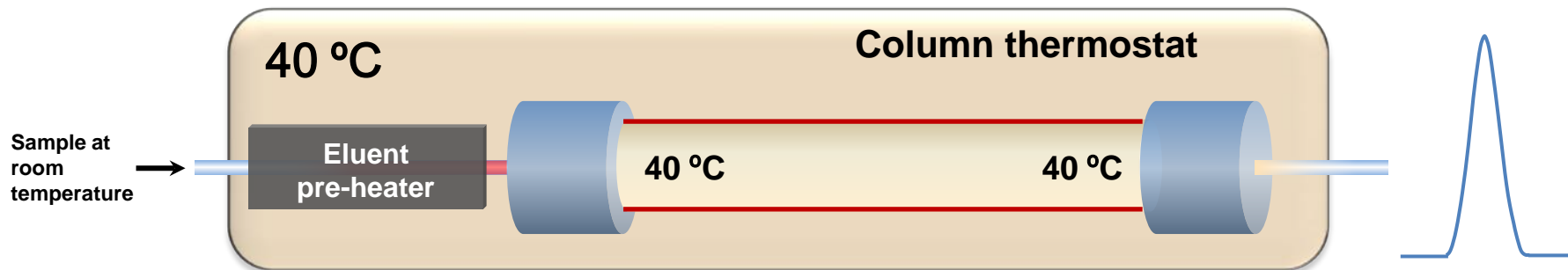
Temperature Effect with Viscous Friction Inside the Column

HPLC column, run without notable frictional heating:

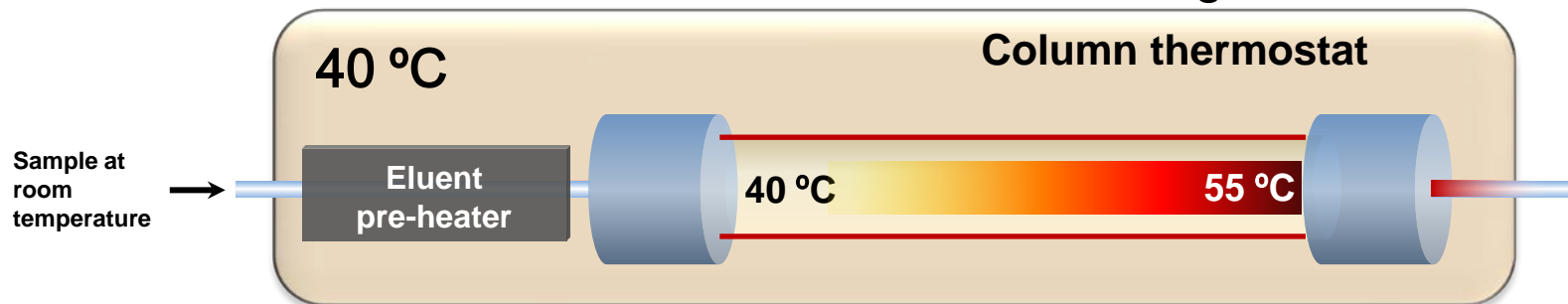


Temperature Effect with Viscous Friction Inside the Column

HPLC column, run without notable frictional heating:

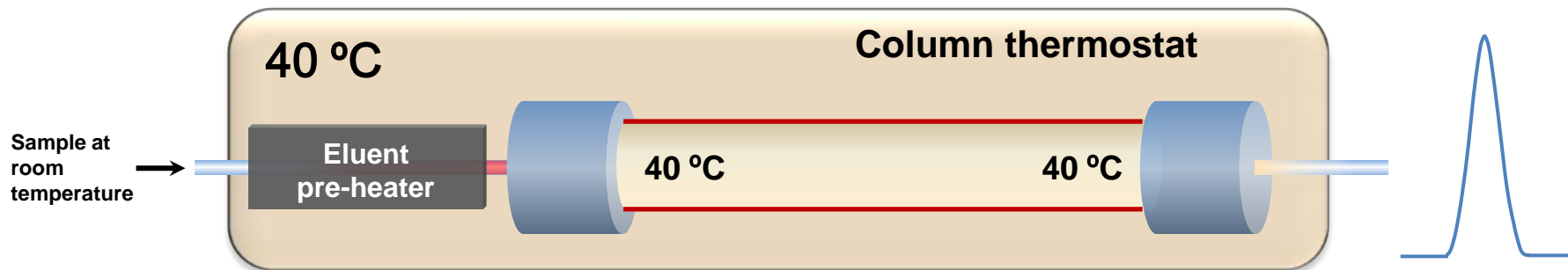


UHPLC column, with substantial frictional heating:

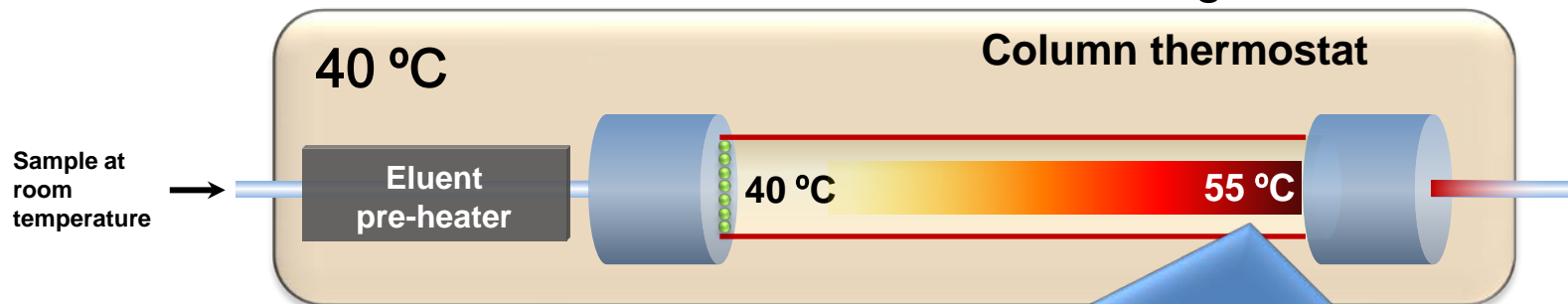


Temperature Effect with Viscous Friction Inside the Column

HPLC column, run without notable frictional heating:



UHPLC column, with substantial frictional heating:

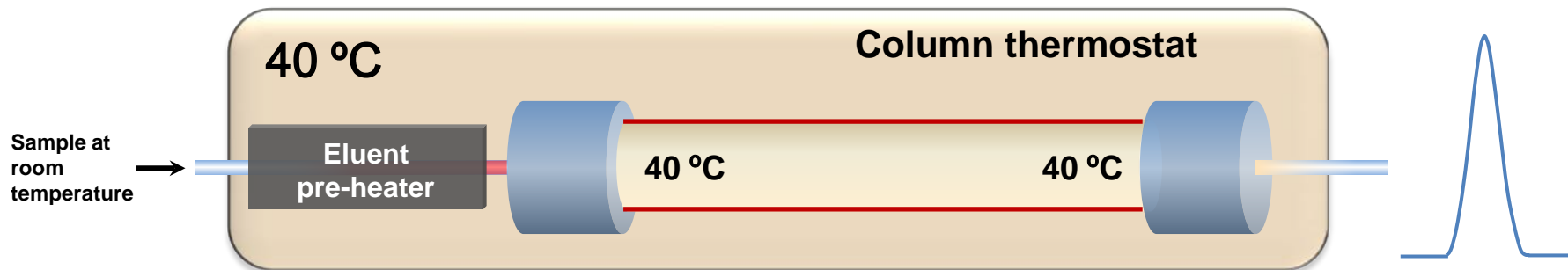


Mismatch with radial temperature gradients:

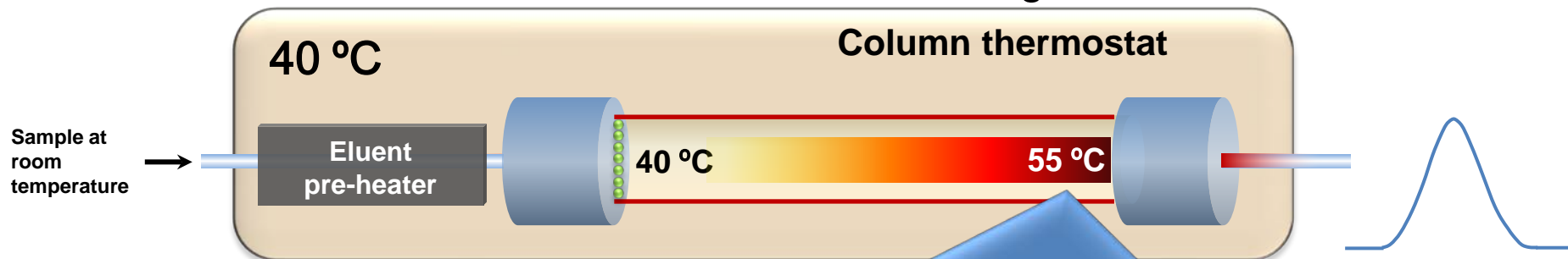
- Column heats up substantially beyond the compartment temperature
 - Lower viscosity leads to higher linear velocity
 - Lower retention

Temperature Effect with Viscous Friction Inside the Column

HPLC column, run without notable frictional heating:



UHPLC column, with substantial frictional heating:

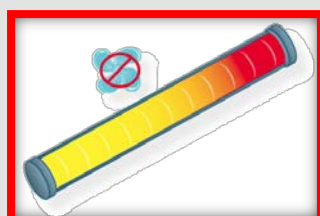
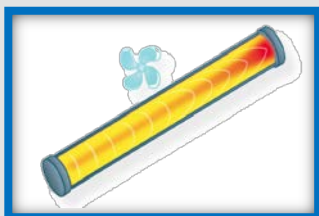
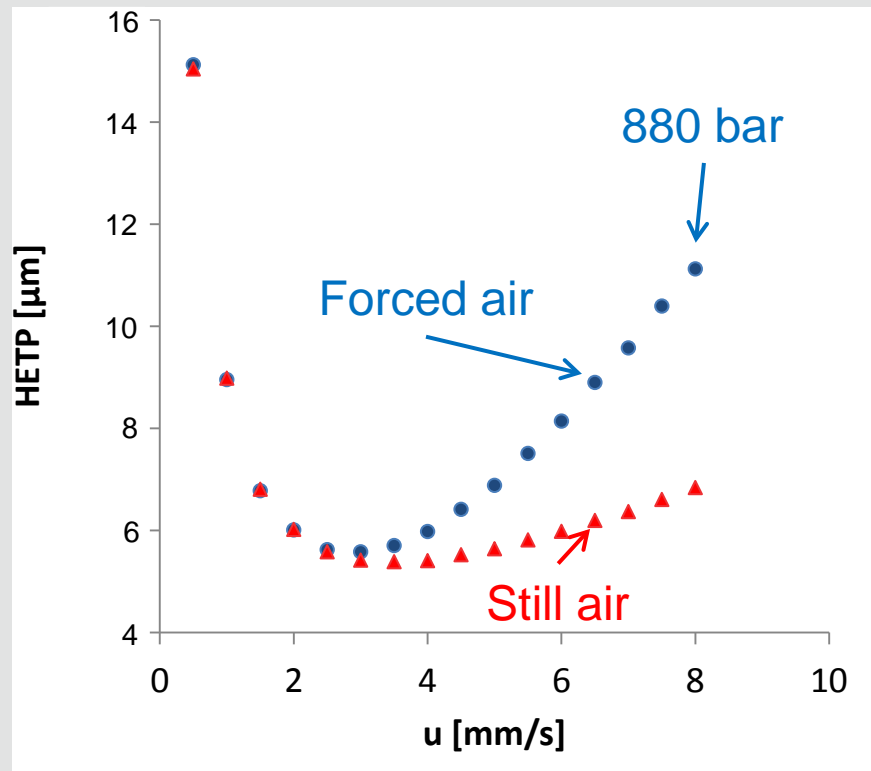


Mismatch with radial temperature gradients:

- Column heats up substantially beyond the compartment temperature
 - Lower viscosity leads to higher linear velocity
 - Lower retention

Frictional Heat Impacting Efficiency and Retention

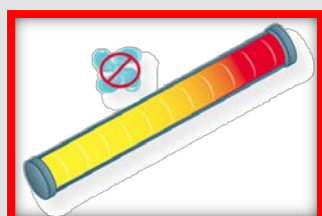
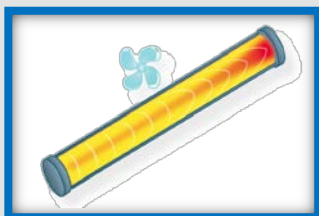
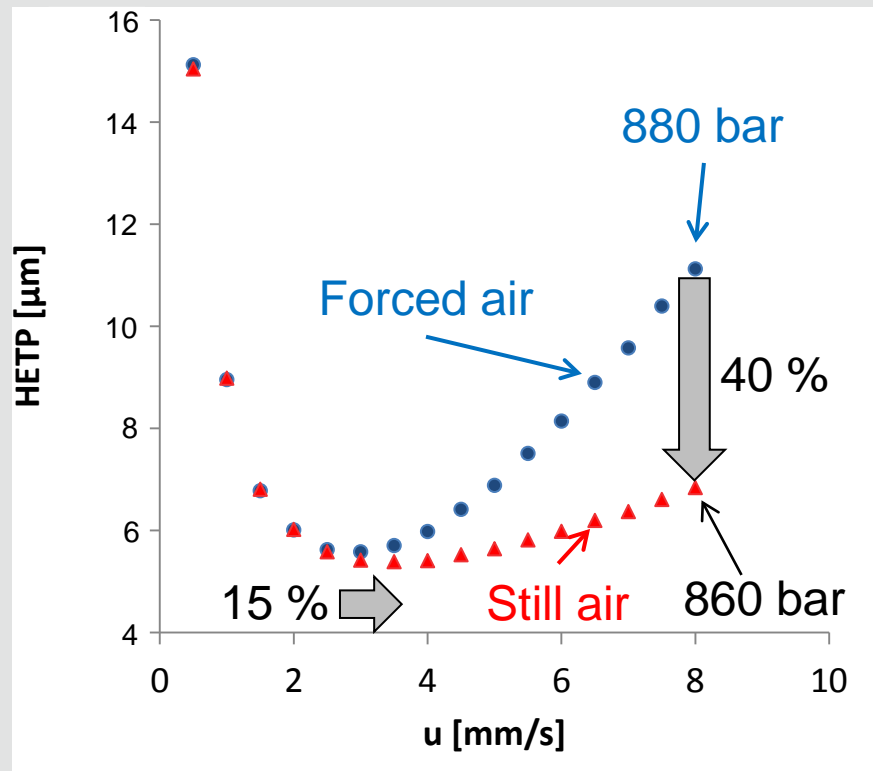
Efficiency:



Column: Thermo Scientific™ Hypersil GOLD™, 1.9 μm, 2.1 x 100 mm
Mobile phase: 50% water/ 50% acetonitrile
Sample: Hexanophenone
Temperature: CC and pre-heater 30 °C, forced air and still air, resp.

Frictional Heat Impacting Efficiency and Retention

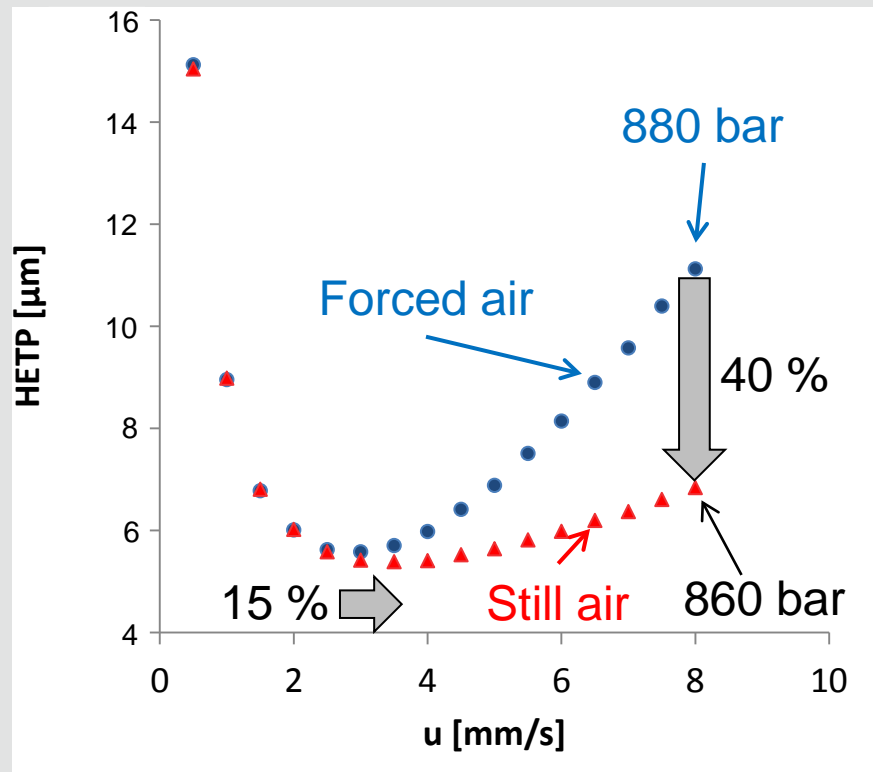
Efficiency:



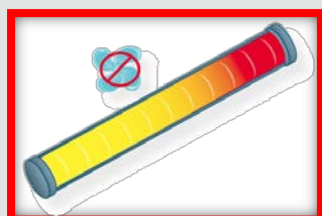
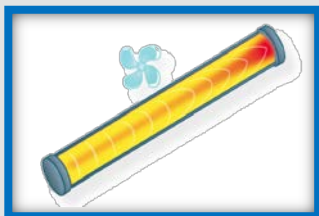
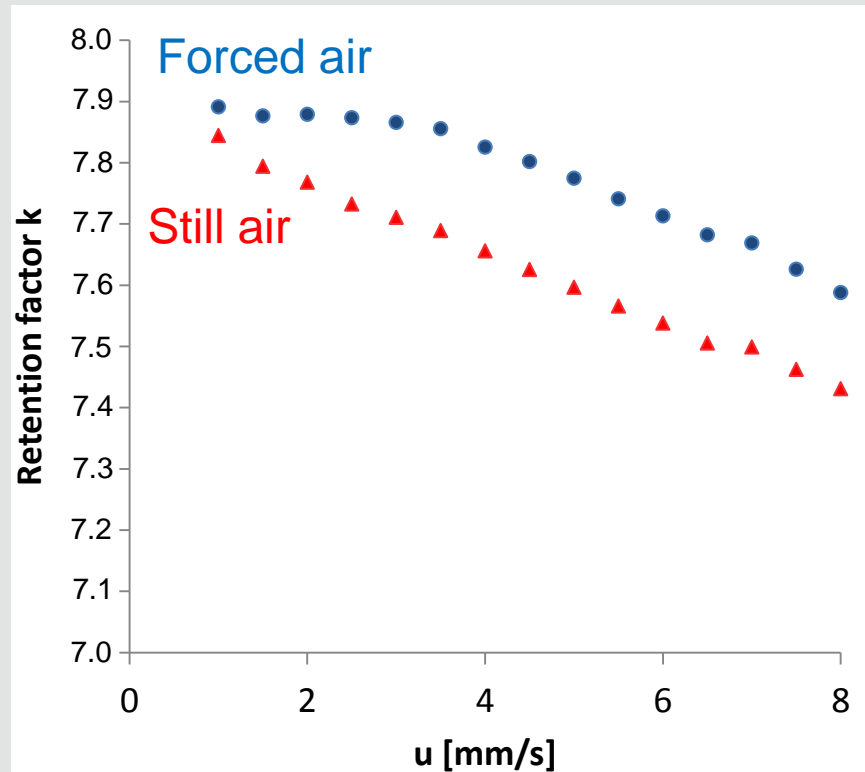
Column: Thermo Scientific™ Hypersil GOLD™, 1.9 μm, 2.1 x 100 mm
Mobile phase: 50% water/ 50% acetonitrile
Sample: Hexanophenone
Temperature: CC and pre-heater 30 °C, forced air and still air, resp.

Frictional Heat Impacting Efficiency and Retention

Efficiency:



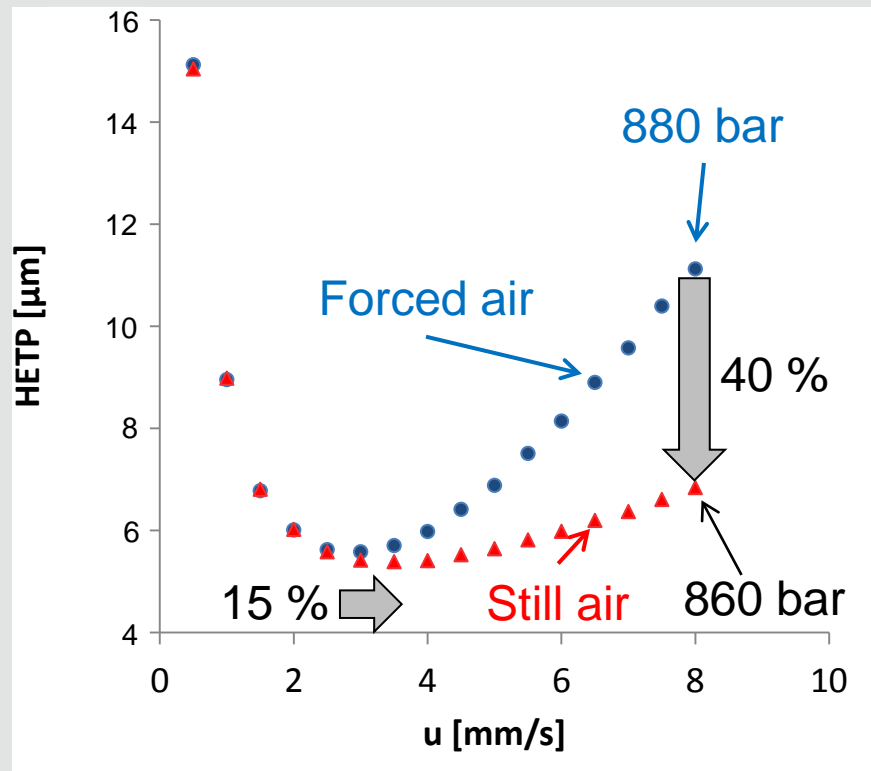
Retention:



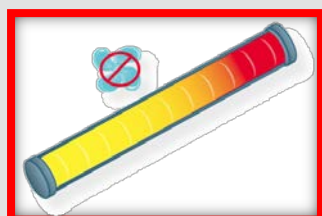
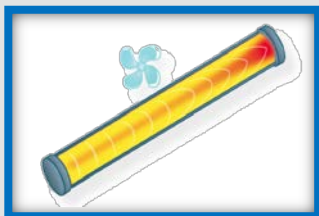
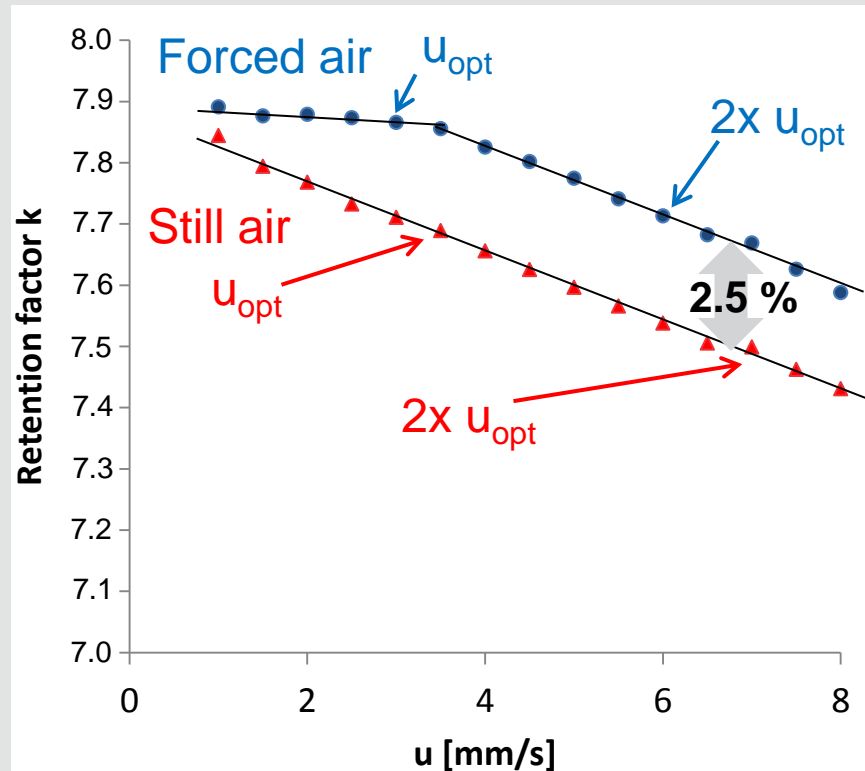
Column: Thermo Scientific™ Hypersil GOLD™, 1.9 μm , 2.1 x 100 mm
Mobile phase: 50% water/ 50% acetonitrile
Sample: Hexanophenone
Temperature: CC and pre-heater 30 $^{\circ}\text{C}$, forced air and still air, resp.

Frictional Heat Impacting Efficiency and Retention

Efficiency:



Retention:

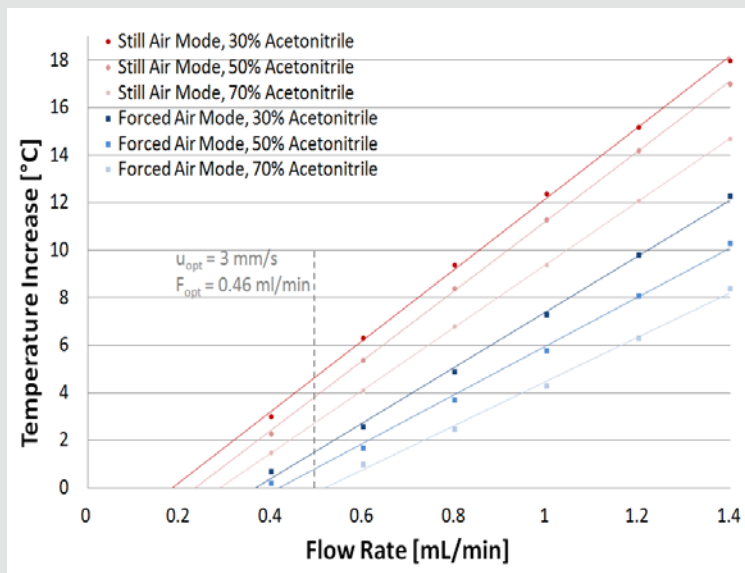


Column: Thermo Scientific™ Hypersil GOLD™, 1.9 μm, 2.1 x 100 mm
Mobile phase: 50% water/ 50% acetonitrile
Sample: Hexanophenone
Temperature: CC and pre-heater 30 °C, forced air and still air, resp.

Measured Temperature Increase in Both Modes

Experimental conditions:

Instrument: Vanquish Horizon UHPLC System
 Column: Hypersil GOLD, 1.9 μm , 2.1 x 100 mm
 Eluent: H₂O/ACN
 Flow rate: up to 1.40 mL/min
 Temperature: 30 °C (Column compartment and eluent pre-heater)



| 50% acetonitrile | Still air mode | | Forced air mode | | |
|------------------|--------------------|----------------|---------------------------|----------------|---------------------------|
| | Flow rate [mL/min] | Pressure [bar] | Temperature increase [°C] | Pressure [bar] | Temperature increase [°C] |
| | 0.2 | 173 | 0.0 | 177 | 0.0 |
| | 0.4 | 343 | 2.3 | 356 | 0.2 |
| | 0.6 | 506 | 5.4 | 527 | 1.7 |
| | 0.8 | 659 | 8.4 | 685 | 3.7 |
| | 1.0 | 809 | 11.3 | 830 | 5.8 |
| | 1.2 | 958 | 14.2 | 975 | 8.1 |
| | 1.4 | 1079 | 17.0 | 1099 | 10.3 |

Effects Resulting from Viscous Friction at 1400 bar

Experimental conditions:

Column: Thermo Scientific™ Accucore™ Vanquish™ C18, 1.5 μm, 2.1 x 100 mm

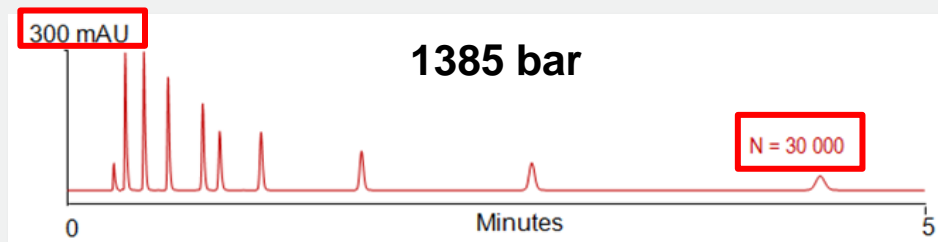
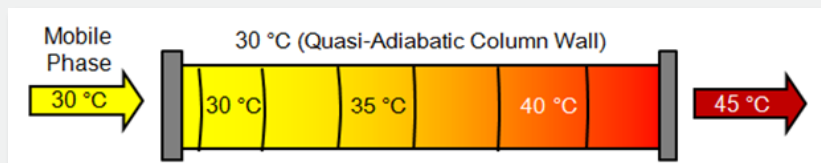
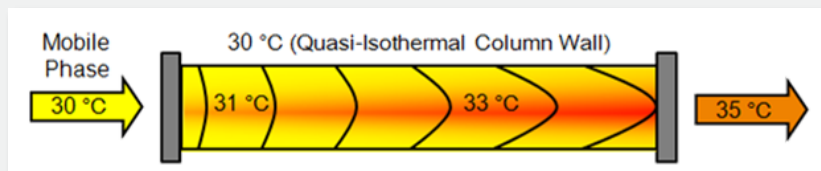
Sample: Uracil, acetanilide and 8 alkylphenones

Inj. volume: 1 μL

Eluent: 45:55 H₂O:ACN (v/v)

Flow rate: 0.65 mL/min

Thermostatting: Compartment and active pre-heater at 30 °C, fan speed 5 (forced air) or fan speed 0 (still air)



Effects Resulting from Viscous Friction at 1400 bar

Experimental conditions:

Column: Thermo Scientific™ Accucore™ Vanquish™ C18, 1.5 μm, 2.1 x 100 mm

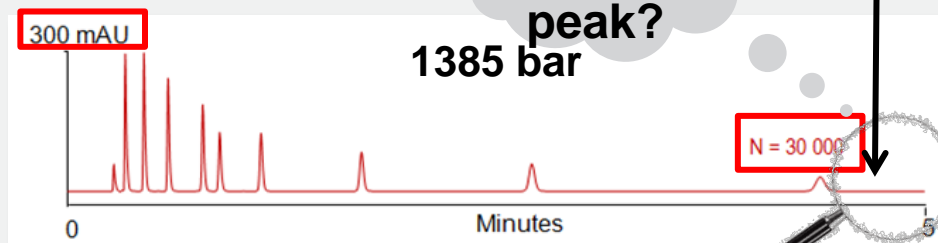
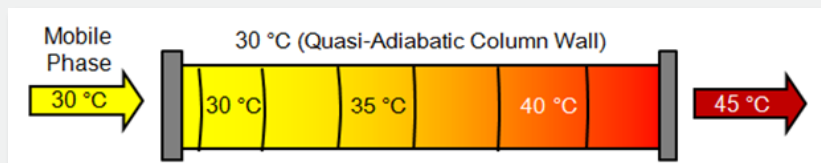
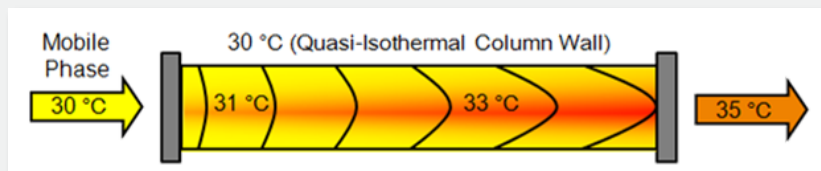
Sample: Uracil, acetanilide and 8 alkylphenones

Inj. volume: 1 μL

Eluent: 45:55 H₂O:ACN (v/v)

Flow rate: 0.65 mL/min

Thermostatting: Compartment and active pre-heater at 30 °C, fan speed 5 (forced air) or fan speed 0 (still air)



Where is my peak?



Effects Resulting from Viscous Friction at 1400 bar

Experimental conditions:

Column: Thermo Scientific™ Accucore™ Vanquish™ C18, 1.5 μm, 2.1 x 100 mm

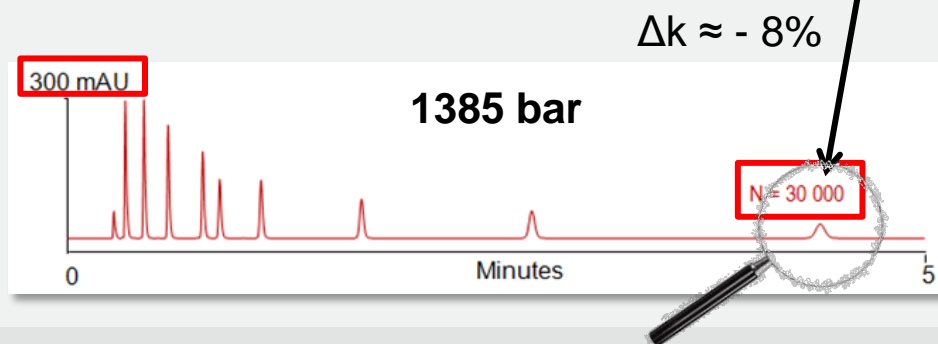
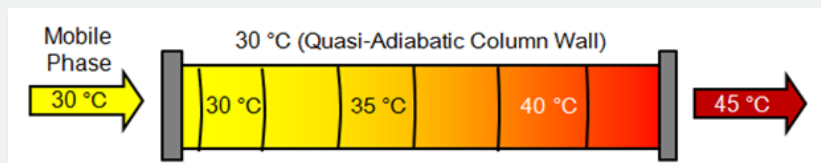
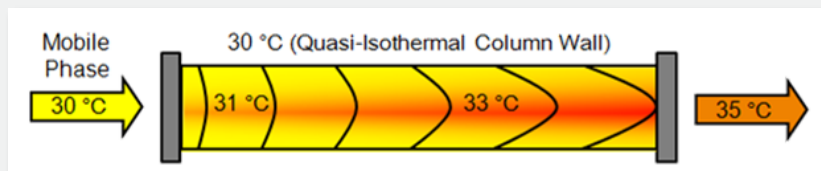
Sample: Uracil, acetanilide and 8 alkylphenones

Inj. volume: 1 μL

Eluent: 45:55 H₂O:ACN (v/v)

Flow rate: 0.65 mL/min

Thermostatting: Compartment and active pre-heater at 30 °C, fan speed 5 (forced air) or fan speed 0 (still air)



Effects Resulting from Viscous Friction at 1400 bar

Experimental conditions:

Column: Thermo Scientific™ Accucore™ Vanquish™ C18, 1.5 μm, 2.1 x 100 mm

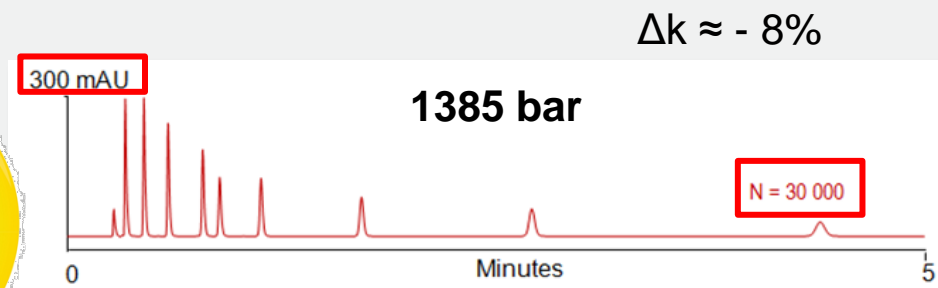
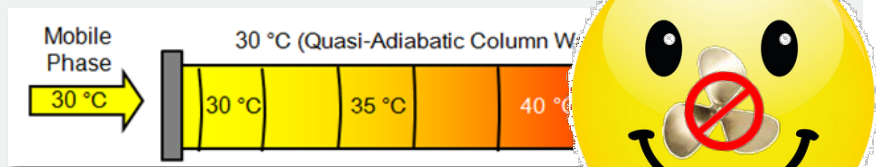
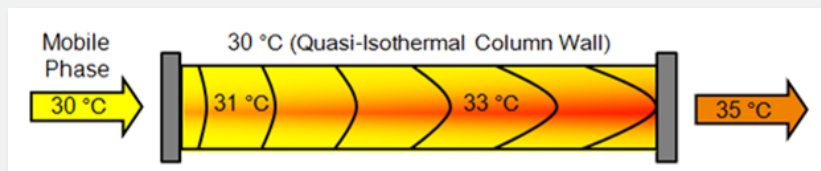
Sample: Uracil, acetanilide and 8 alkylphenones

Inj. volume: 1 μL

Eluent: 45:55 H₂O:ACN (v/v)

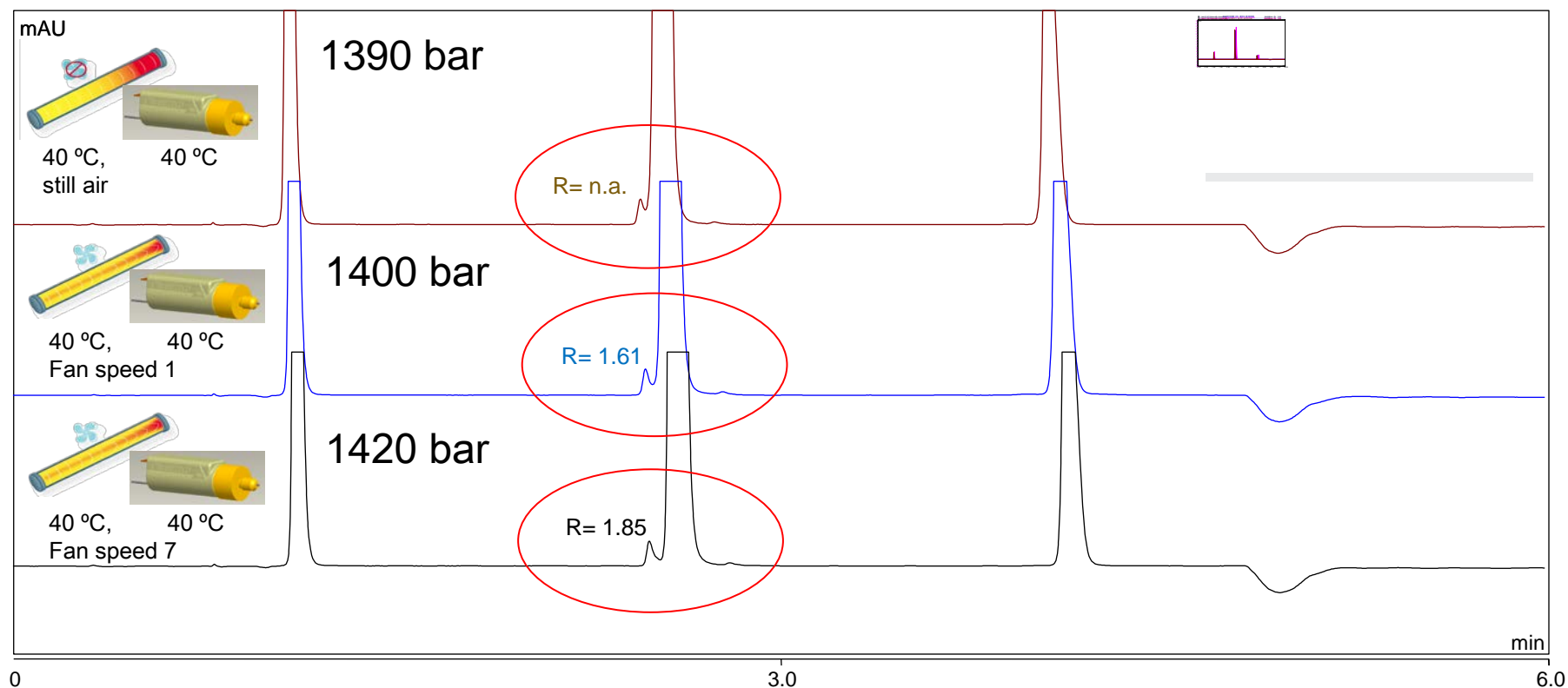
Flow rate: 0.65 mL/min

Thermostatting: Compartment and active pre-heater at 30 °C, fan speed 5 (forced air) or fan speed 0 (still air)



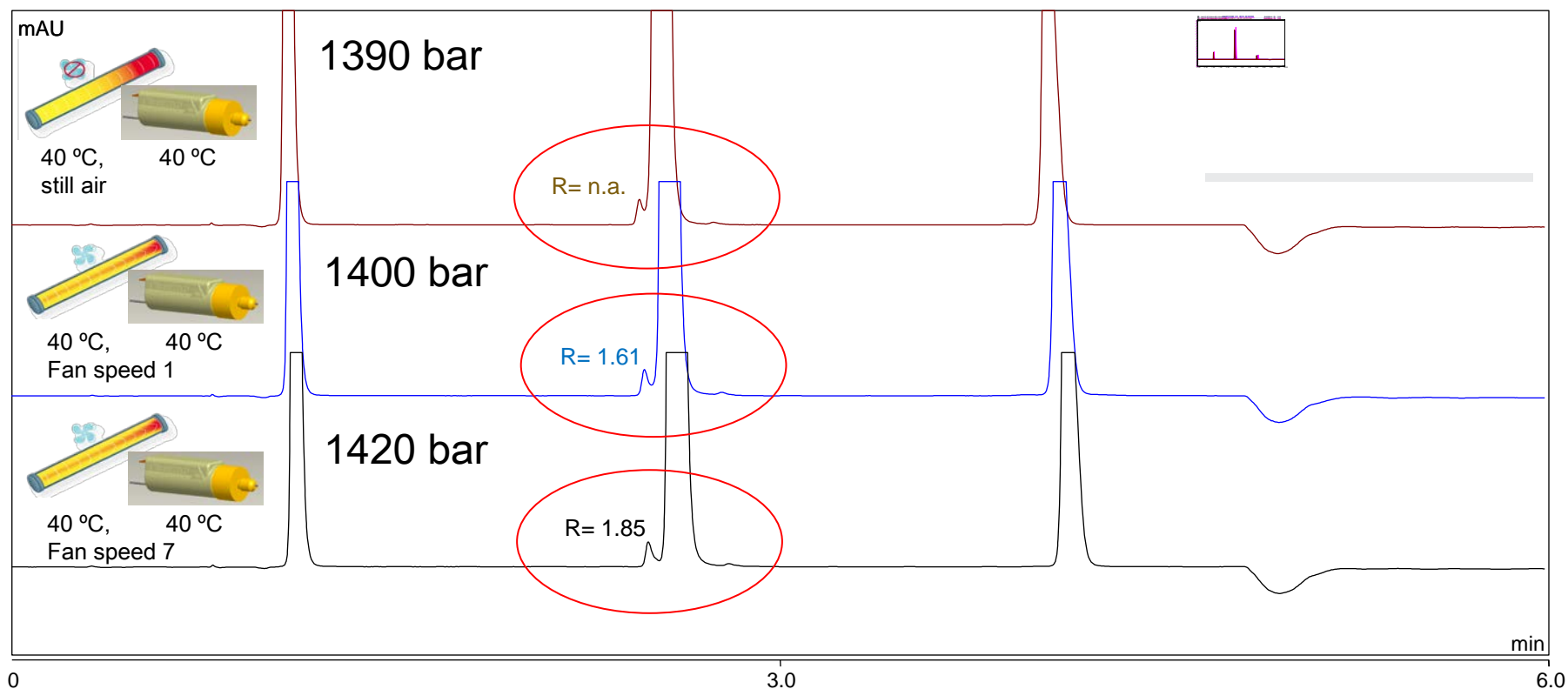
Still air is the better mode

Fan Speed Variation with Another Separation Example



Instrument: Vanquish Horizon UHPLC system
Gradient: From 5% to 30% B in 4 min,
A: 0.1% formic acid in H₂O, B: ACN
Column: Hypersil GOLD Vanquish, 1.9 μm, 2.1 x 200 mm
Sample: Mix of gallic acid, caffeic acid, and salicylic acid
Detection: UV 300 nm, 20 Hz, 0.2 s response time
Flow rate: 0.8 mL/min

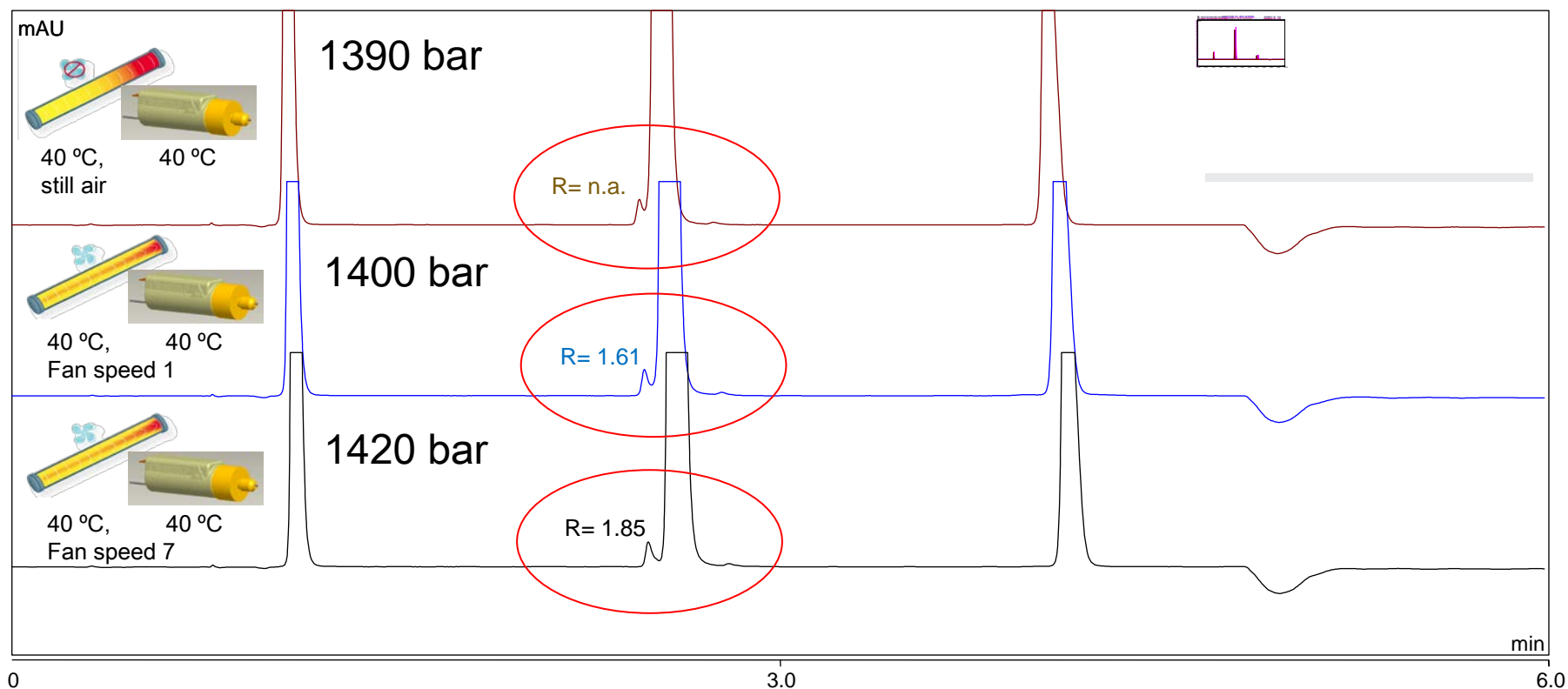
Fan Speed Variation with Another Separation Example



- Enhanced air circulation lowers the column temperature.
- The selectivity for the impurity peak improves.

Instrument: Vanquish Horizon UHPLC system
Gradient: From 5% to 30% B in 4 min,
A: 0.1% formic acid in H₂O, B: ACN
Column: Hypersil GOLD Vanquish, 1.9 μm, 2.1 x 200 mm
Sample: Mix of gallic acid, caffeic acid, and salicylic acid
Detection: UV 300 nm, 20 Hz, 0.2 s response time
Flow rate: 0.8 mL/min

Fan Speed Variation with Another Separation Example



- Enhanced air circulation lowers the column temperature.
- The selectivity for the impurity peak improves.

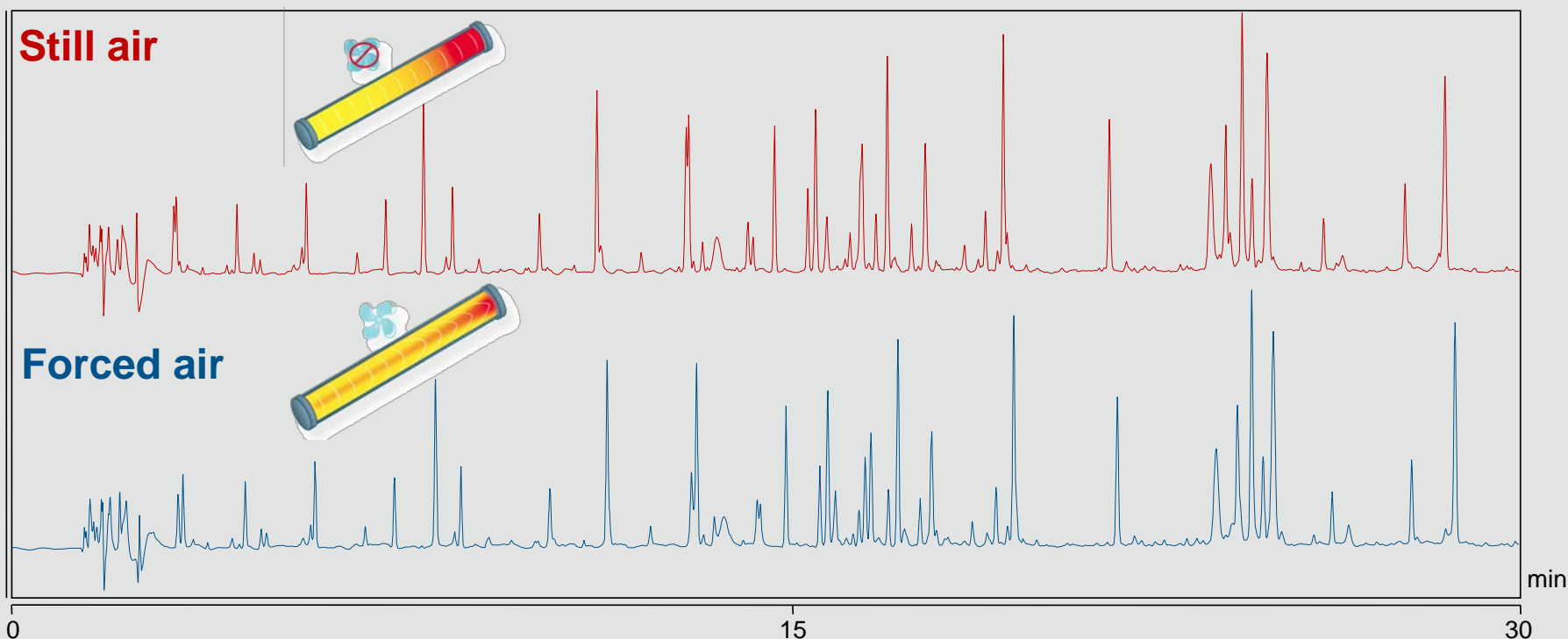


Instrument: Vanquish Horizon UHPLC system
Gradient: From 5% to 30% B in 4 min,
A: 0.1% formic acid in H₂O, B: ACN
Column: Hypersil GOLD Vanquish, 1.9 μm, 2.1 x 200 mm
Sample: Mix of gallic acid, caffeic acid, and salicylic acid
Detection: UV 300 nm, 20 Hz, 0.2 s response time
Flow rate: 0.8 mL/min



Forced air is the better mode

Column Thermostating for Peptide Mapping



Vanquish Horizon UHPLC system with 10 mm Thermo Scientific™ Lightpipe™ flow cell

Column: Thermo Scientific™ Acclaim™ RSLC 120 C18, 2.2 μm, 2.1 x 250 mm

Sample: 1 μL Monoclonal IgG Tryptic Digest (2 mg/mL)

Detection: 214 nm, 100 Hz

Temperature: CC and pre-heater 80 °C, **still air** and **forced air**, respectively

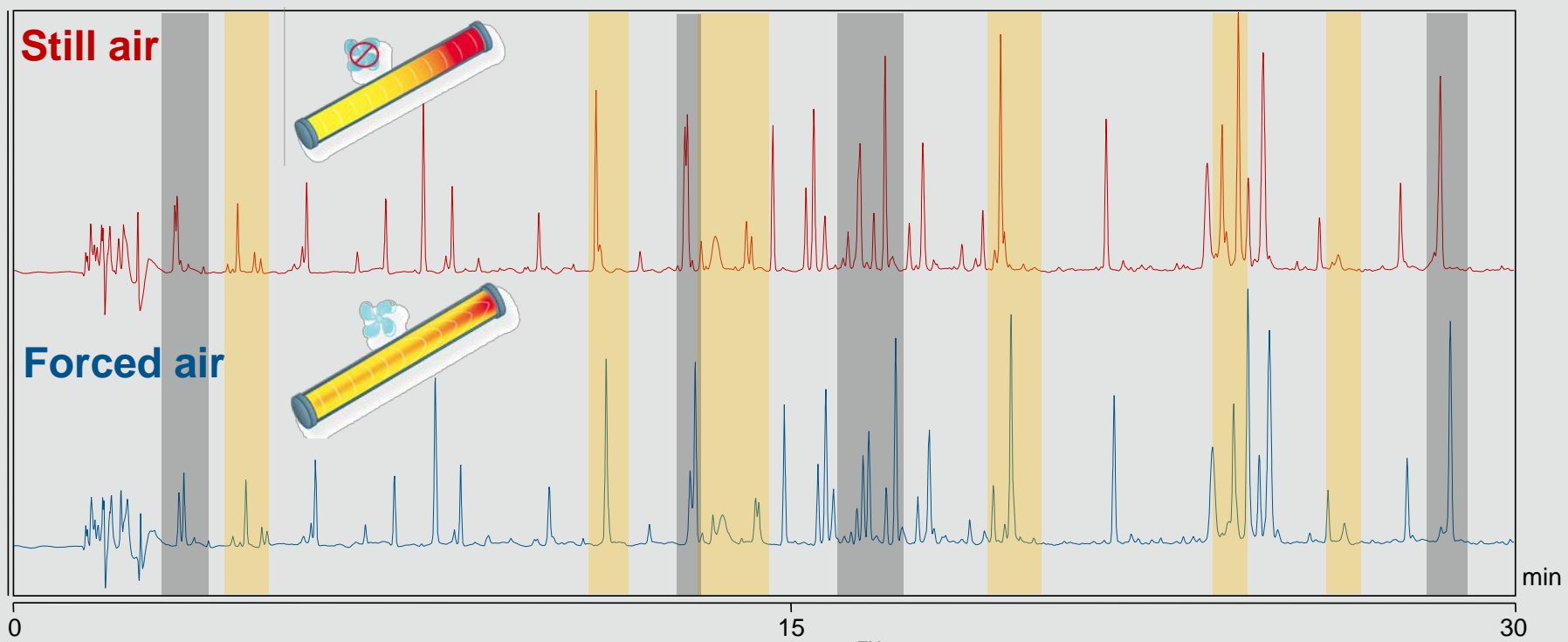
Mobile phase A: TFA 0.05% in water

Mobile phase B: TFA 0.05% in 2/8 water/ACN

Flow rate: 400 μL/min


Gradient: From 4% to 55% B in 30 min


Column Thermostatting for Peptide Mapping



Vanquish Horizon UHPLC system with 10 mm Thermo Scientific™ Lightpipe™ flow cell
Column: Thermo Scientific™ Acclaim™ RSLC 120 C18, 2.2 μm, 2.1 x 250 mm
Sample: 1 μL Monoclonal IgG Tryptic Digest (2 mg/mL)
Detection: 214 nm, 100 Hz
Temperature: CC and pre-heater 80 °C, **still air** and **forced air**, respectively

Mobile phase A: TFA 0.05% in water
Mobile phase B: TFA 0.05% in 2/8 water/ACN
Flow rate: 400 μL/min
Gradient: From 4% to 55% B in 30 min

 Still air mode beneficial → Partially better resolution due to enhanced efficiency

 Forced air mode beneficial → Partially better resolution due to enhanced selectivity

 **Both modes are complementary**

Summary: Variation of Frictional Heat Dissipation

These effects are relevant with methods beyond 500 bar:

- Air circulation in the column thermostat is hardware-depending.
- Still air thermostats typically provide better resolution with frictional heat.
- Different thermostats can lead to different chromatograms even with the same effective temperature.

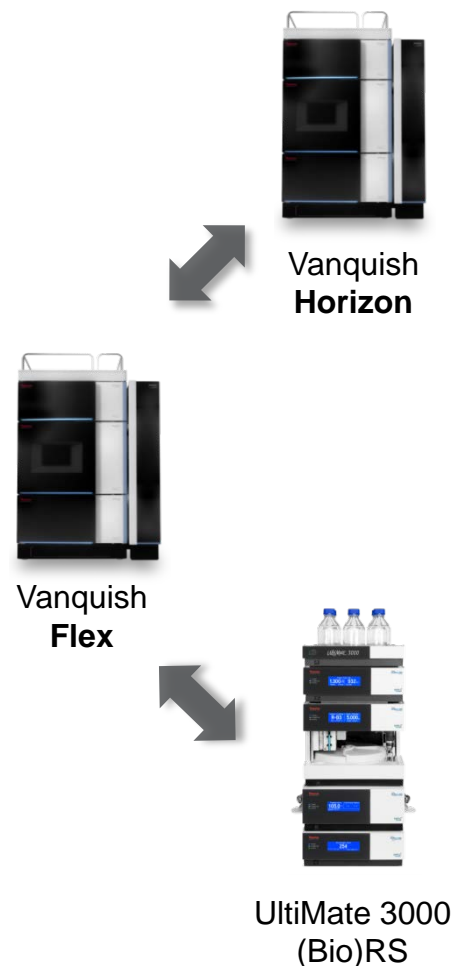
Hints:

- Vary the forced air ventilation if feasible (Vanquish UHPLC systems).
- Try to insulate the column inside a forced-air thermostat.
- Try to vary the position of the column inside the thermostat.

Conclusions

Gradient delay volume (GDV)

- The GDV decides on how long the sample is separated isocratically by the hold-up at the beginning of the gradient elution.
- Knowing and adjusting the GDV are decisive for the successful method transfer.
- Different GDVs between two systems can be adjusted
 - Either by the fluidics (e.g. by changing fluidic parts).
 - Or by changing the gradient program.
- An individually adjustable metering device in the autosampler of the Vanquish UHPLC systems allows an easy and automated GDV adjustment.



Conclusions

Thermostating

- Column thermostats are frequently underrated – “no oven is like the other”
- The operation principle of the thermostat has a substantial impact on the success of a method transfer
- **Still air thermostats**
 - Lock the frictional heat in the column
 - Thus lead to better efficiencies and sharper peaks
 - But internal frictional heating has the potential to change selectivities
- **Forced air thermostats**
 - Dissipate frictional heat from the column surface
 - Which may result in peak distortion due to a radial temperature gradient
- Both modes are complementary and can be used for fine-tuning of a separation



Vanquish
Horizon



Vanquish
Flex



UltiMate 3000
(Bio)RS

Any questions?



**Do you have additional questions
or do you want to talk to an expert from
Thermo Fisher Scientific?**

**Please send an E-Mail to
analyze.eu@thermofisher.com
and we will get back to you.**