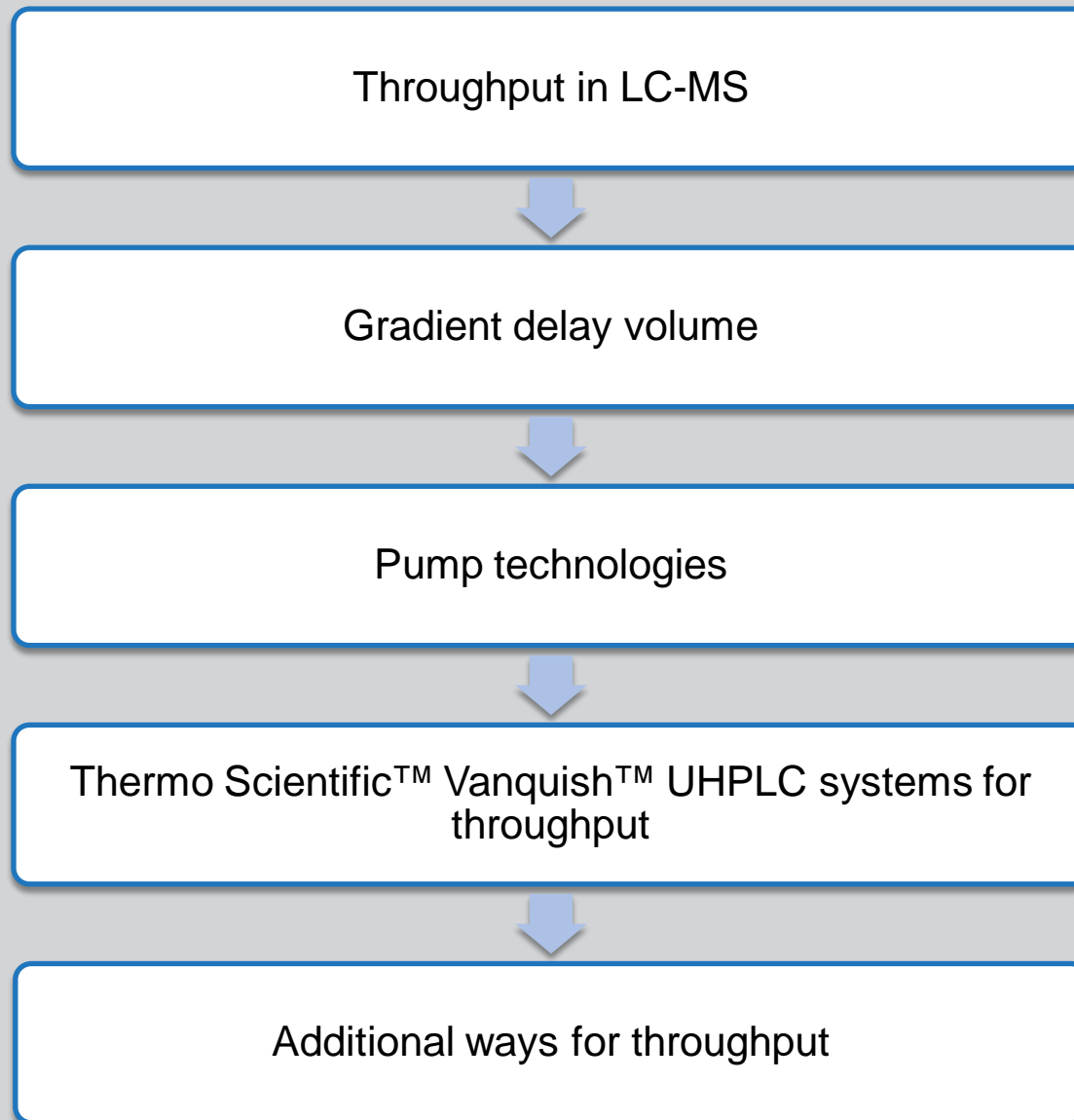




Which UHPLC for LC-MS?

Dr. Giorgia Greco
Product Manager HPLC for MS Systems
Thermo Fisher Scientific, Germering, Germany

Content of the Webinar



What is UHPLC?

- **System back pressure above 600 bar**
- **Analytical column with sub-2 micron particles**



To achieve:

- **Faster separation**
- **Better resolution**
- **Less waste**



Vanquish Flex systems

Quaternary system

- Quaternary solvent blending
- Pressures up to 1000 bar
- Flow rates of up to 8 mL/min
- Biocompatible

Binary system

- Binary high pressure solvent mixing
- Pressures up to 1000 bar
- Flow rates of up to 8 mL/min
- Biocompatible



Up to 1000 bar

Vanquish Horizon systems

- Highest pressure capability up to 1500 bar
- Flow rates up to 5 mL/min
- Lowest system dispersion and GDV
- Unmatched detection sensitivity and linearity
- Biocompatible



1500 bar

Why is Throughput in LC-MS Important?

We require sophisticated LC-MS technology.



We must manage regulatory demands.



We need systems with high-resolution, selectivity, sensitivity, robustness, large number of molecules to be discovered, identified and quantified in one single run.



We require high throughput capabilities at the same time.



Which UHPLC system is suitable for our high throughput LC-MS analysis?

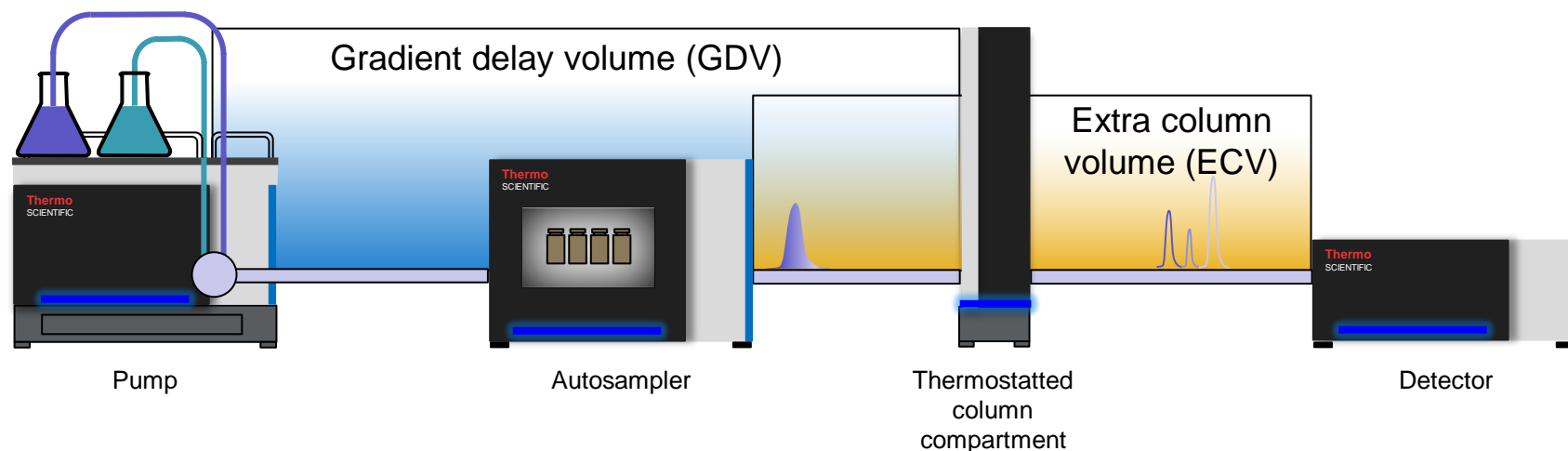


Vanquish UHPLC system
Thermo Scientific™ Q Exactive™ hybrid quadrupole Orbitrap™ MS

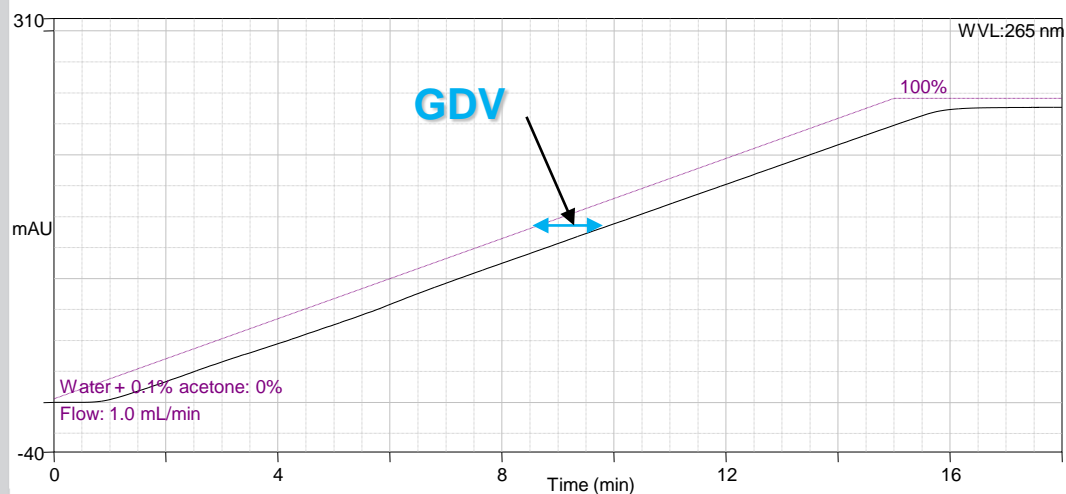


Vanquish UHPLC system
Thermo Scientific™ TSQ Quantiva™ triple quadrupole MS

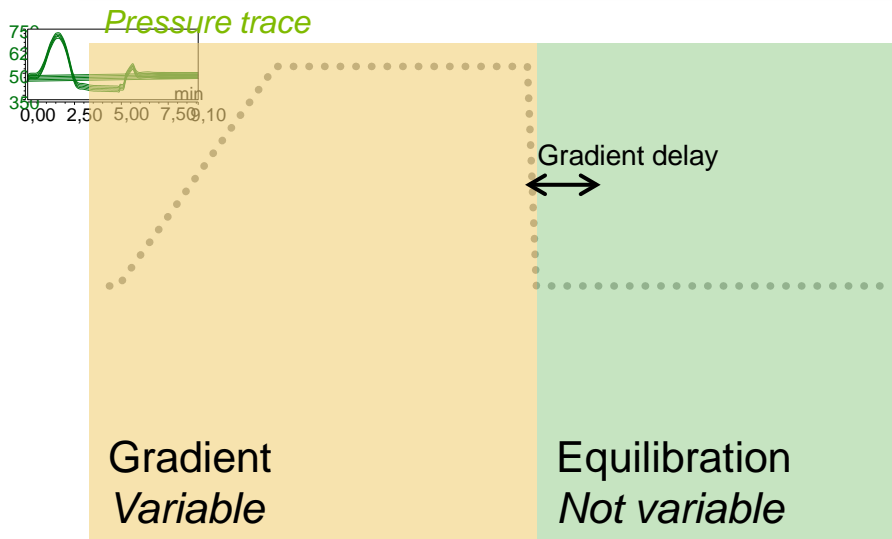
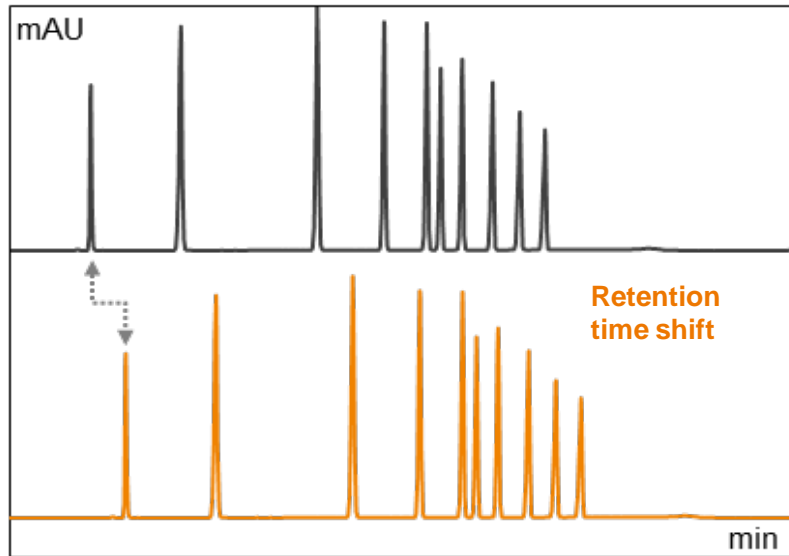
Why is the GDV of your “Front-End” Important for Throughput?



- **Gradient delay volume (GDV):**
Volume of fluid between mixing point of the gradient and column head
- **Extra column volume (ECV):**
Volume of fluid between sample injection point and midpoint of the detector's flow cell.



Why is the GDV of your “Front-End” Important for Throughput?

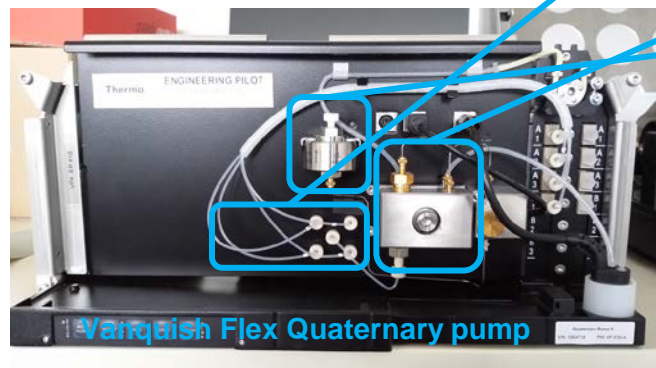
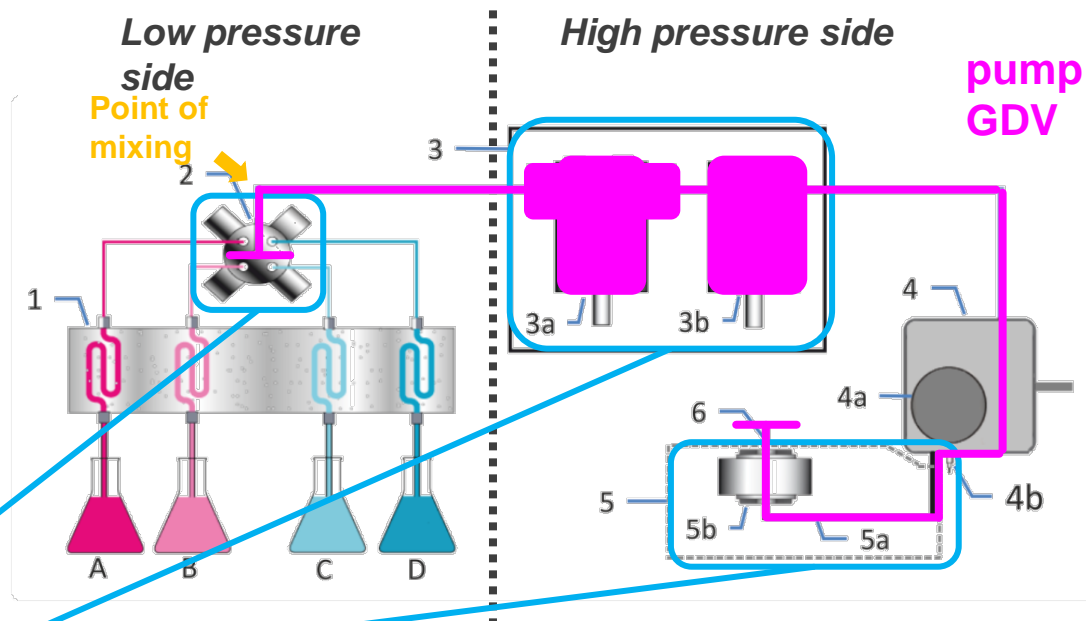


Short runtimes and higher throughput

- Column equilibration time is related to the GDV.
- Small GDV - shorter equilibration times
- Contribution of equilibration time is relevant for short runs.

Why do UHPLC Systems have Different System GDVs?

UHPLC systems can use different pump technologies: *Low pressure gradient (LPG)*



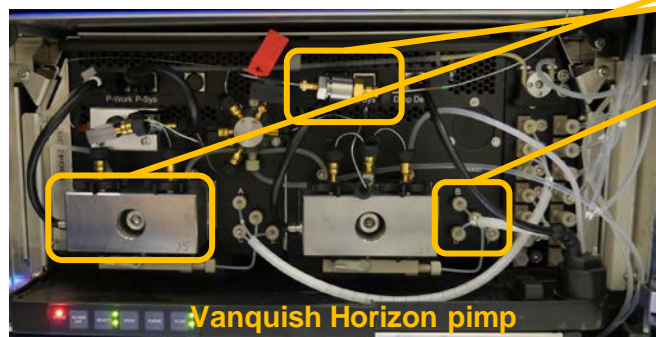
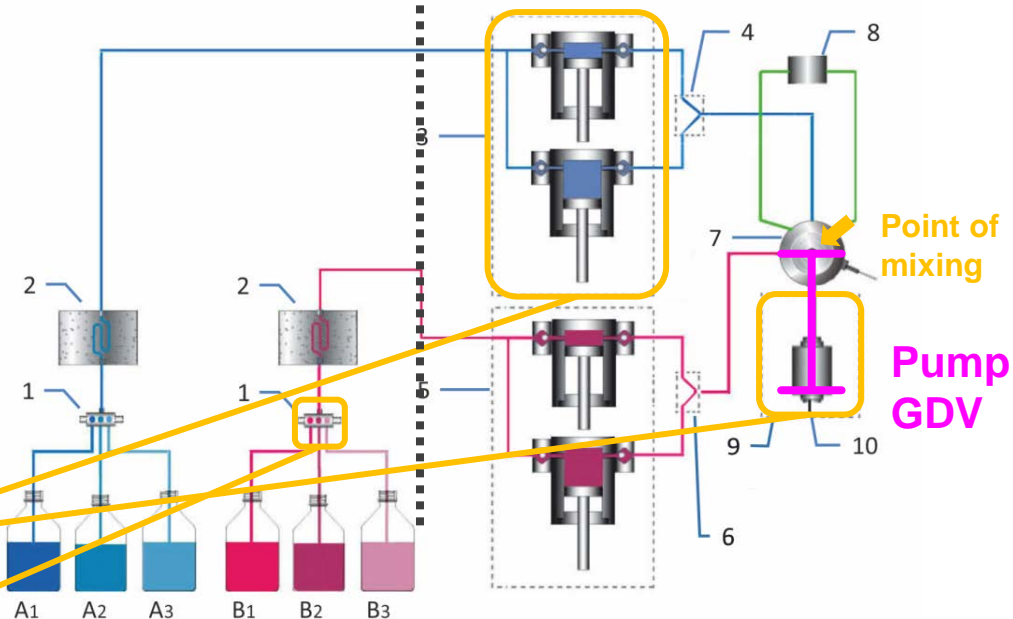
By default, LPG pump systems have large GDVs.

Why do UHPLC Systems have Different System GDVs?

UHPLC systems can use different pump technologies: *High pressure gradient (HPG)*

Low pressure side

High pressure side



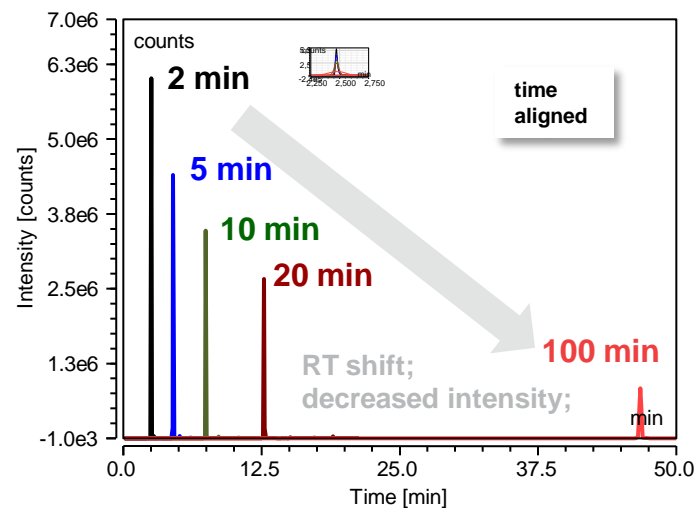
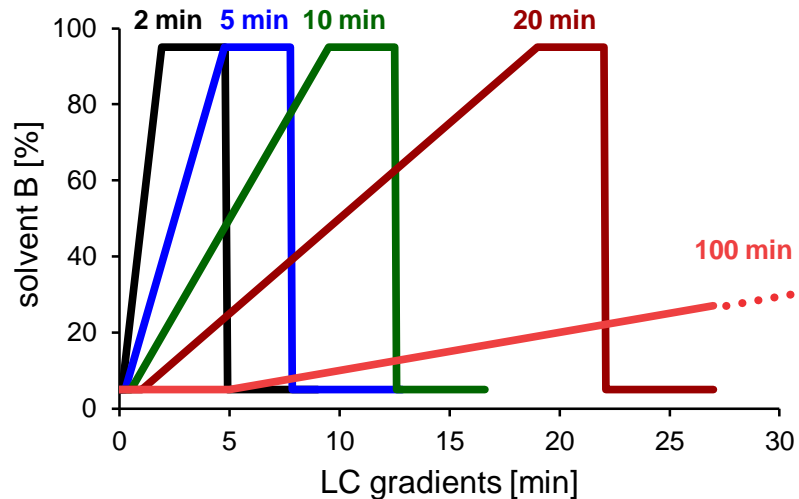
By default, HPG pump systems have small GDVs.

What UHPLC for LC-MS?

Tested Vanquish systems

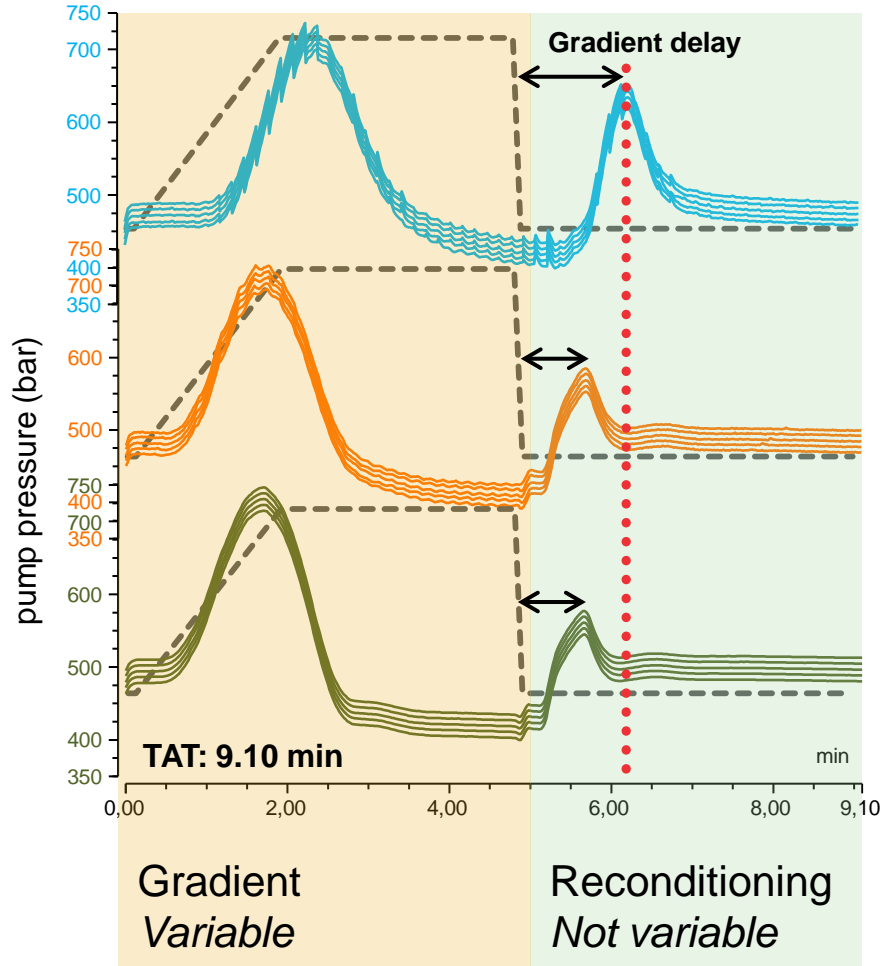
Vanquish Flex Quaternary	Vanquish Flex Binary	Vanquish Horizon
Low pressure gradient (LPG)	High pressure gradient (HPG)	High pressure gradient (HPG)
used mixers: 25 + 10 = 35 μ L	used mixer: 25 + 10 = 35 μ L	used mixer: 25 + 10 = 35 μ L
1034 bar	1034 bar	1500 bar

LC-MS analysis of pesticides



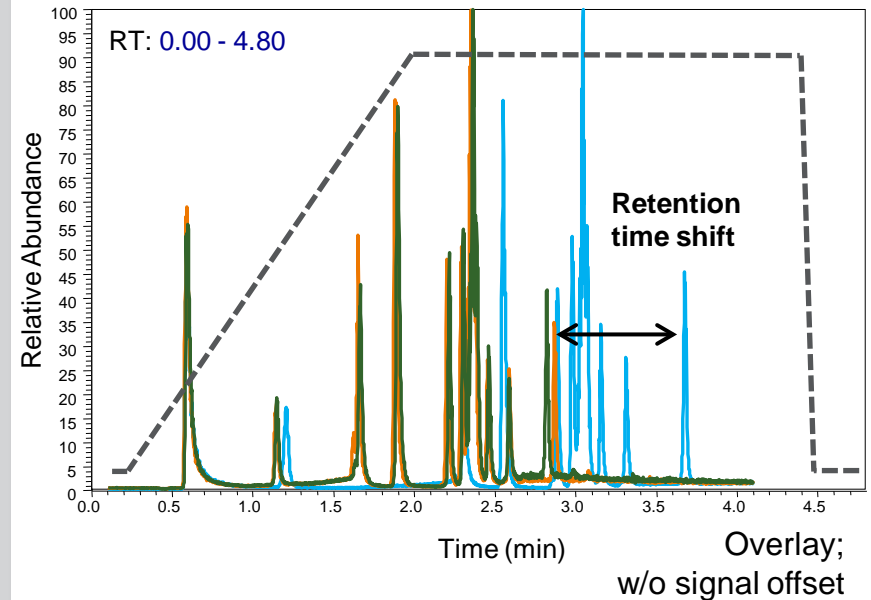
Total Analysis Time (TAT) of Vanquish UHPLC Systems

Pressure traces (2 min gradient)



Overlay (n=5); w/ signal offset

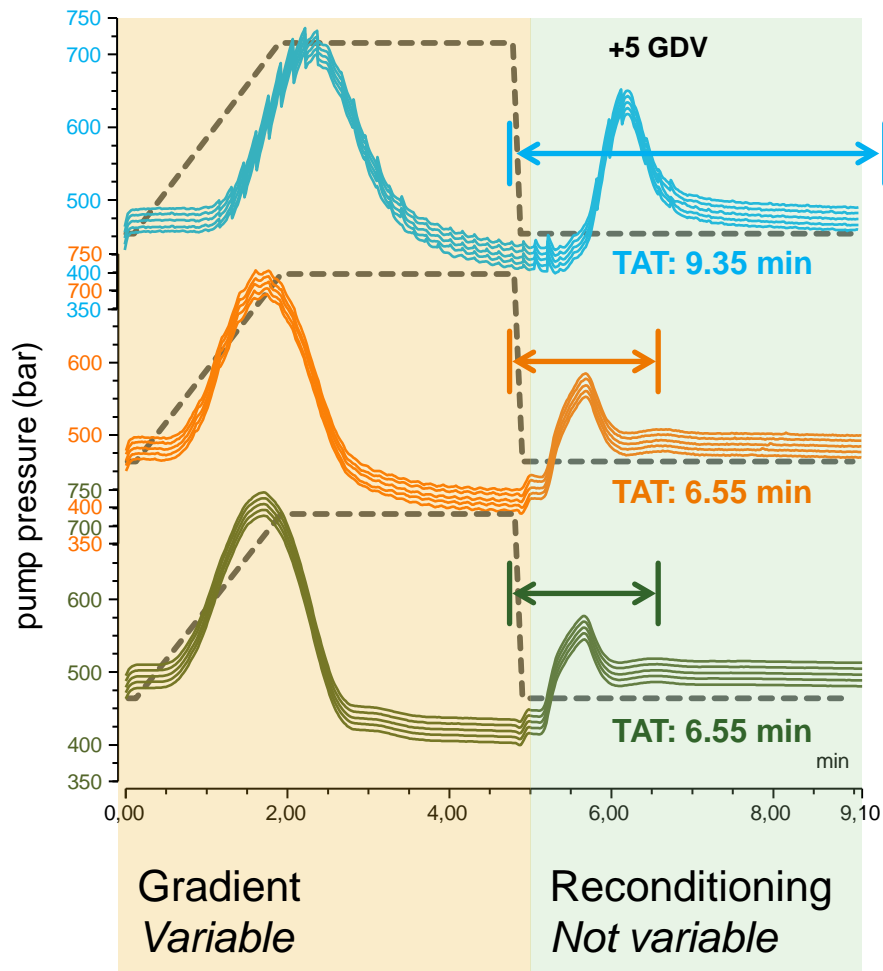
Extracted ion chromatograms



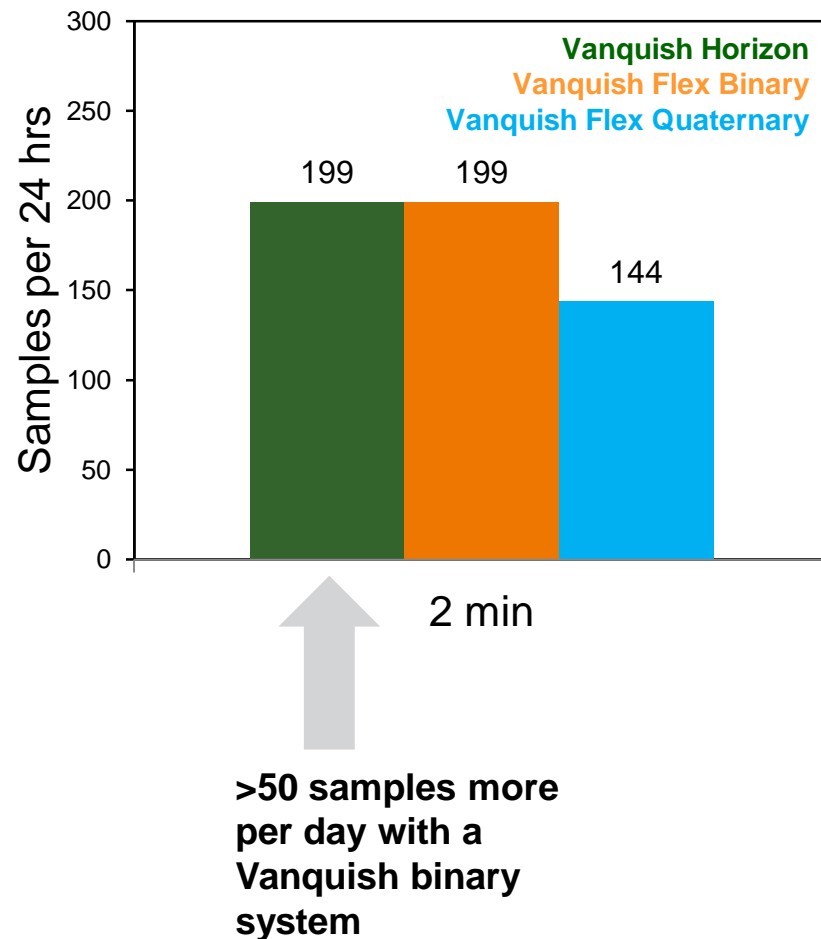
Vanquish Flex Quaternary
Vanquish Flex Binary
Vanquish Horizon

TAT Optimization of Vanquish UHPLC Systems

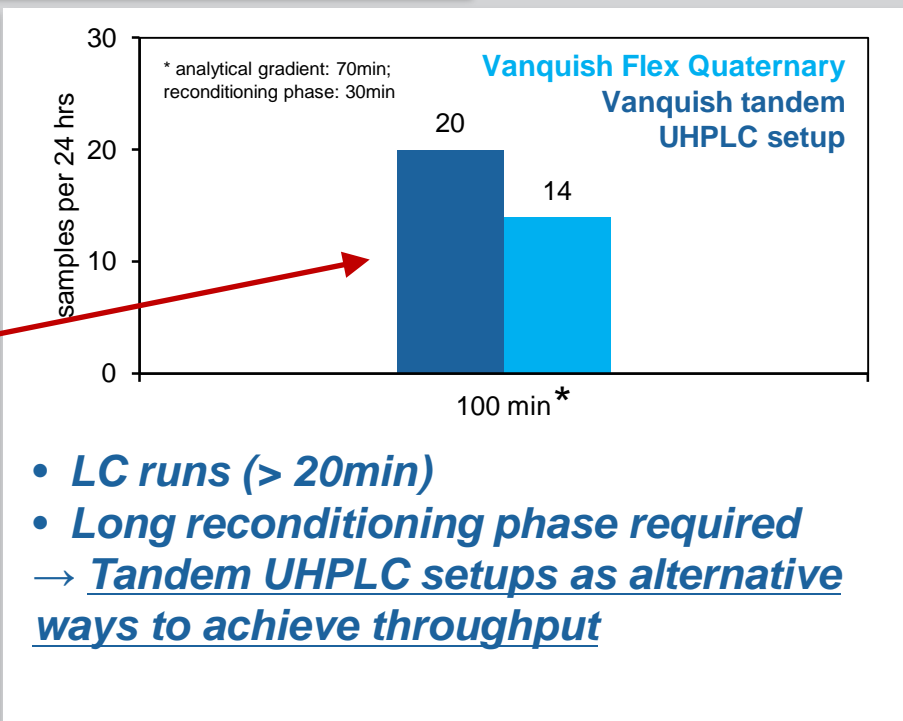
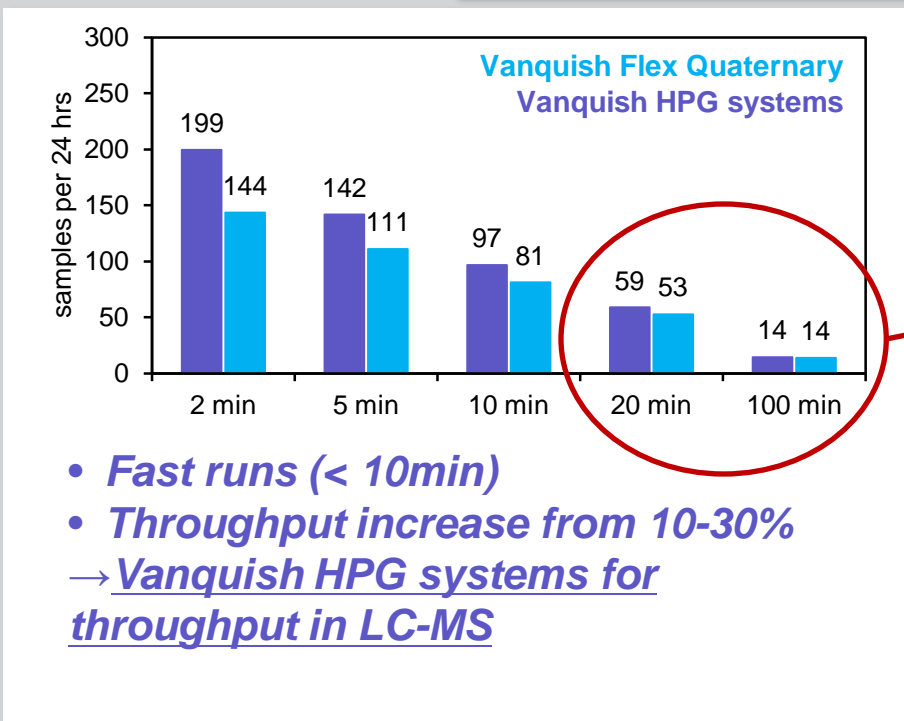
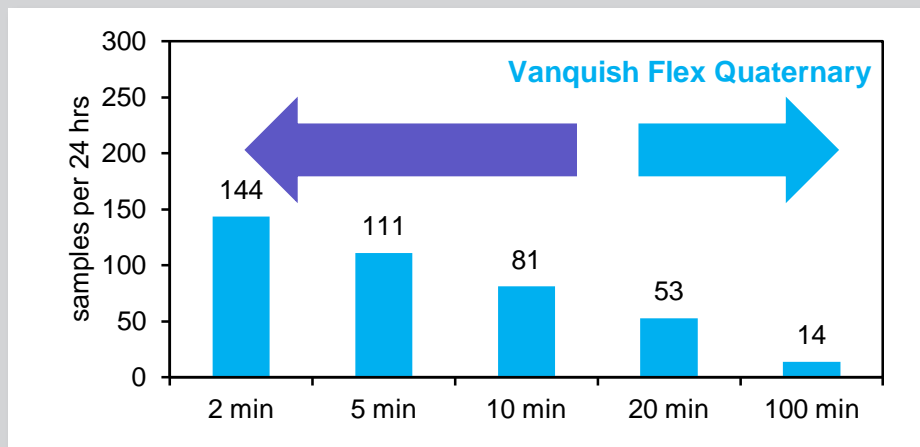
Pressure traces (2 min gradient)



Sample throughput



Throughput Considerations

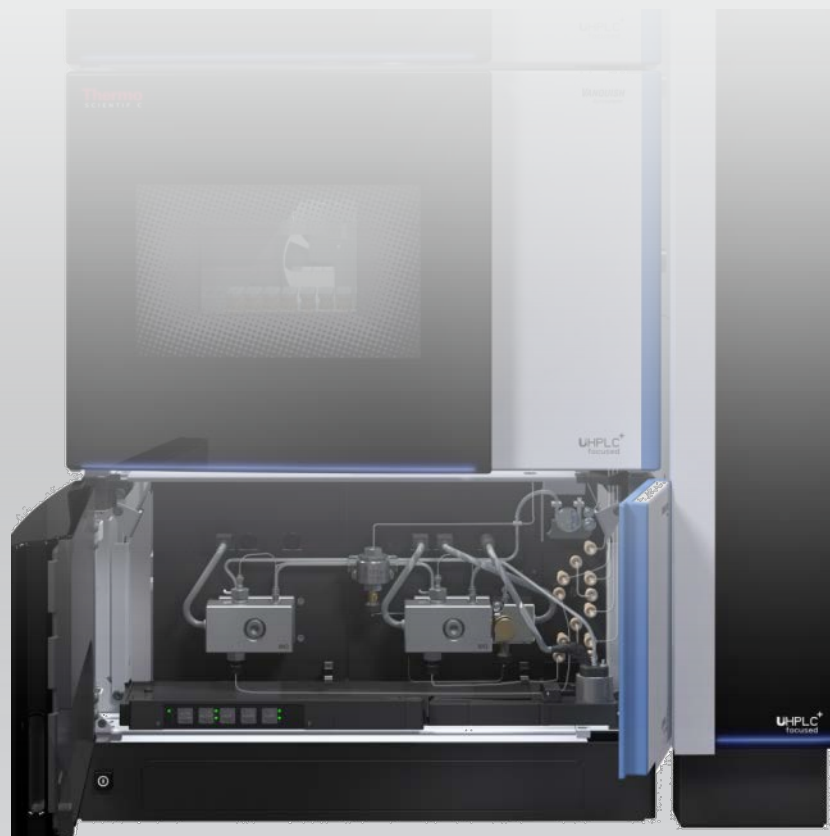


- **Fast runs (< 10min)**
- **Throughput increase from 10-30%**
→ Vanquish HPG systems for throughput in LC-MS

- **LC runs (> 20min)**
- **Long reconditioning phase required**
→ Tandem UHPLC setups as alternative ways to achieve throughput

Vanquish Flex Binary and Horizon Systems – What's the difference?

Vanquish Flex Binary



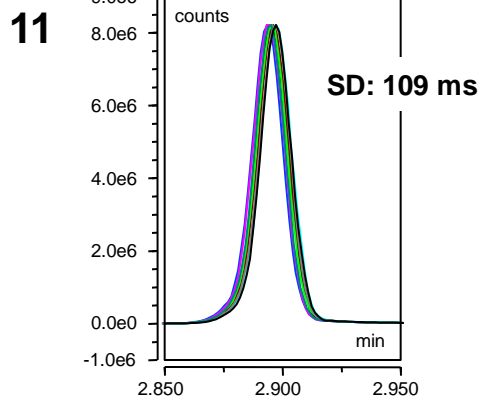
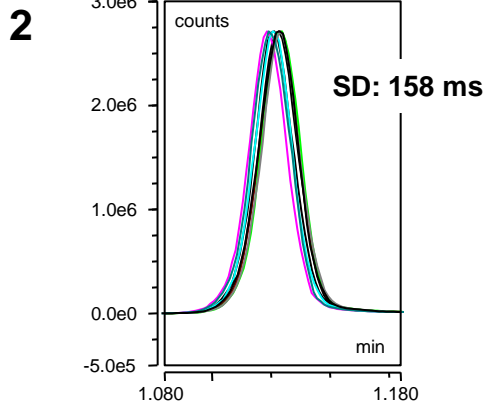
Vanquish Horizon



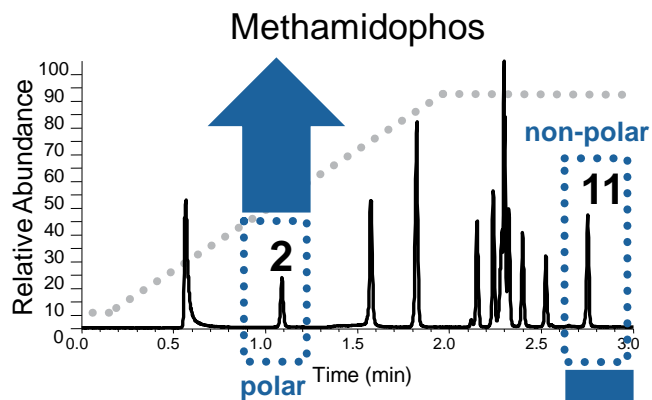
Vanquish Horizon system operates with the superior pump technology for critical and challenging applications.

Vanquish Flex Binary and Horizon Systems – What's the difference?

Vanquish Flex Binary



60 ms \equiv 0.001 min

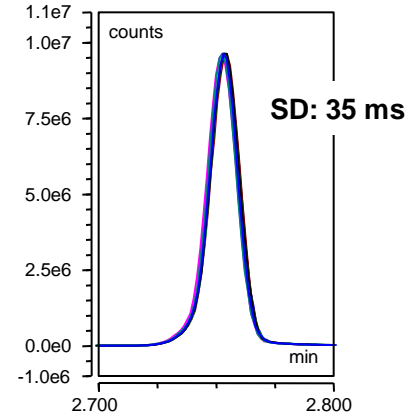
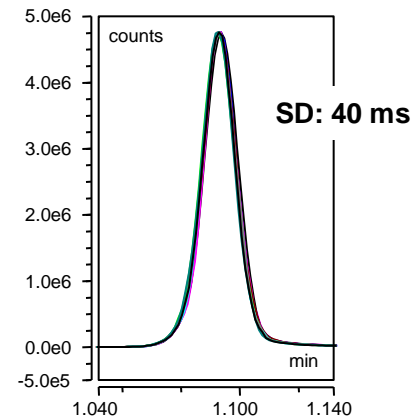


2 min

Quinoxifen

RT precision for 199 consecutive injections

Vanquish Horizon



Higher RT precision of the Vanquish Horizon system can reduce your number of technical replicates, one important aspect to be considered for high-throughput.

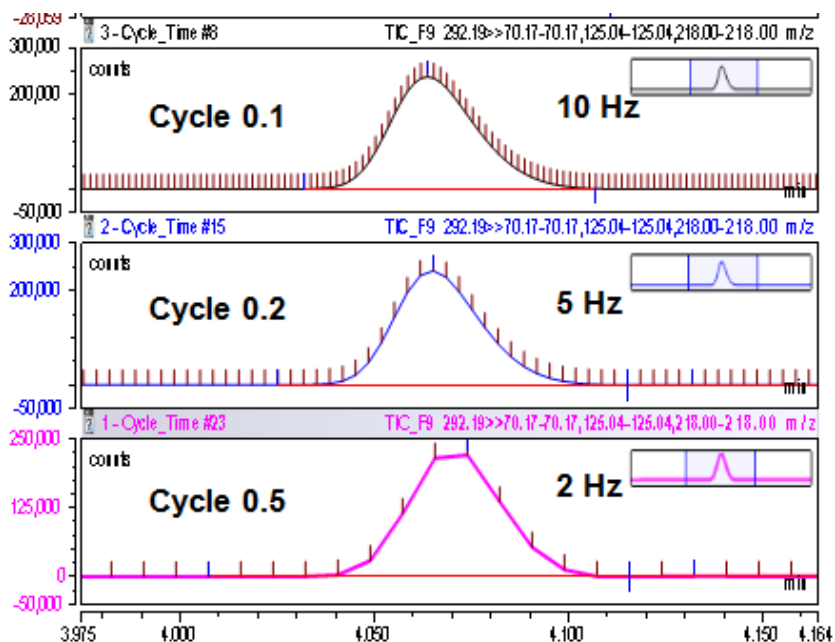
What Level of Throughput is Possible with the Vanquish Horizon?

LC-MS analysis of pesticides

LC conditions

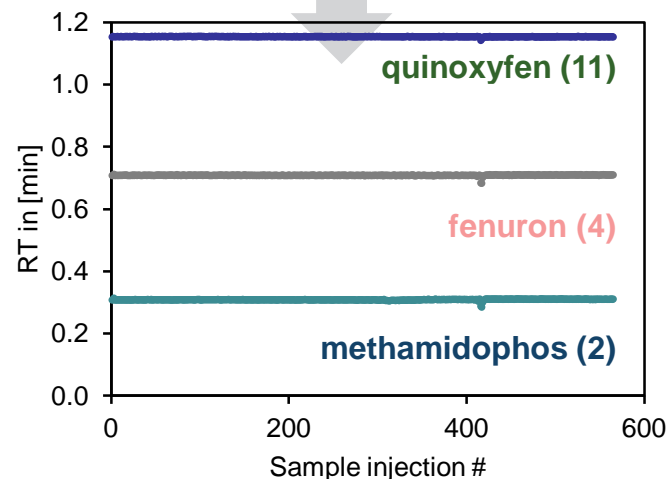
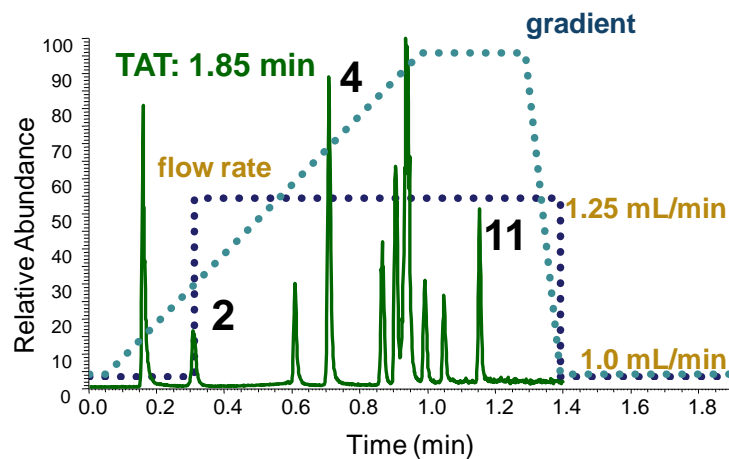
Column Thermo Scientific™ Hypersil GOLD™
50 x 2.1 mm, 1.9 μm

567 samples per day



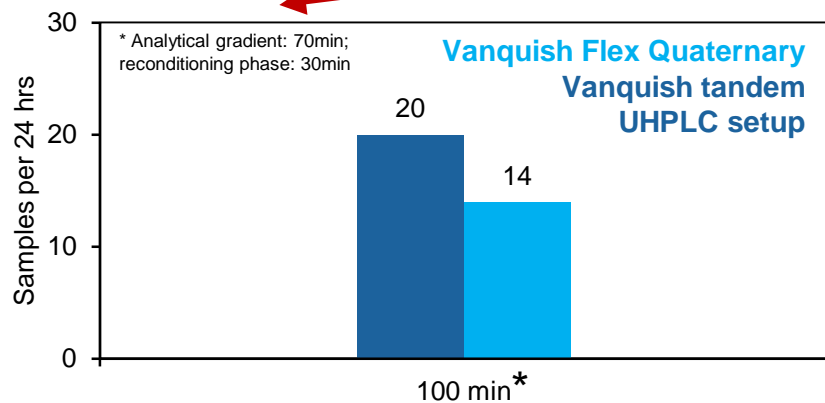
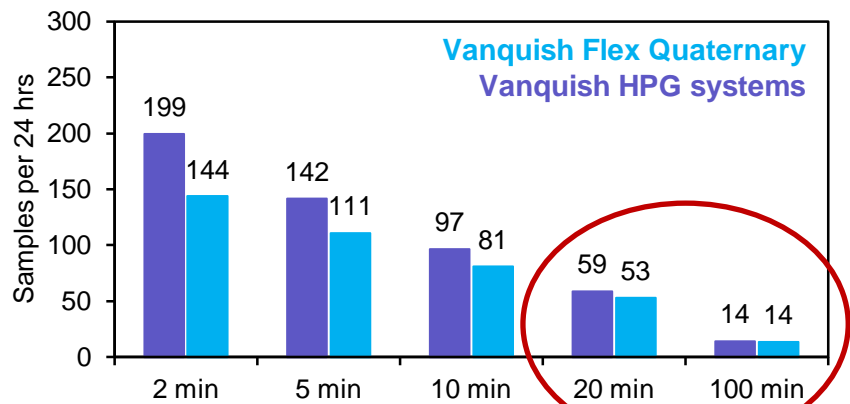
At least 10 Hz are required to obtain RT SD = 0.001 min for peak width at 50% of 0.025 min (1.5 s).

Throughput maximization

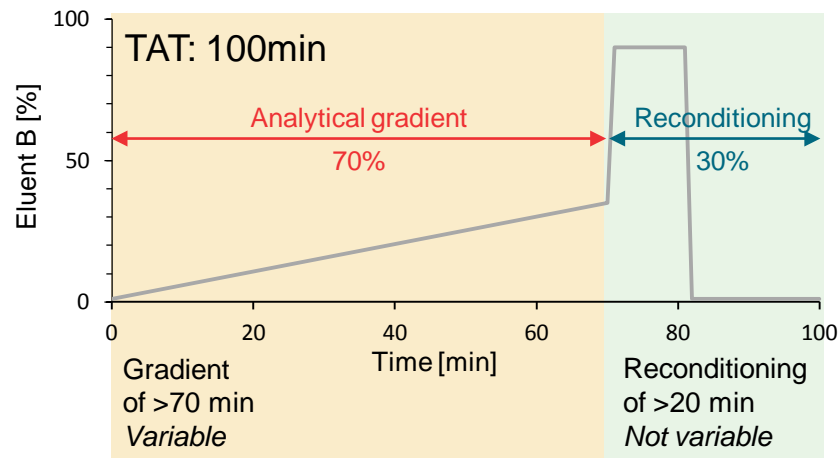


Additional Ways For Throughput

Discussion starter



Shallow LC gradient



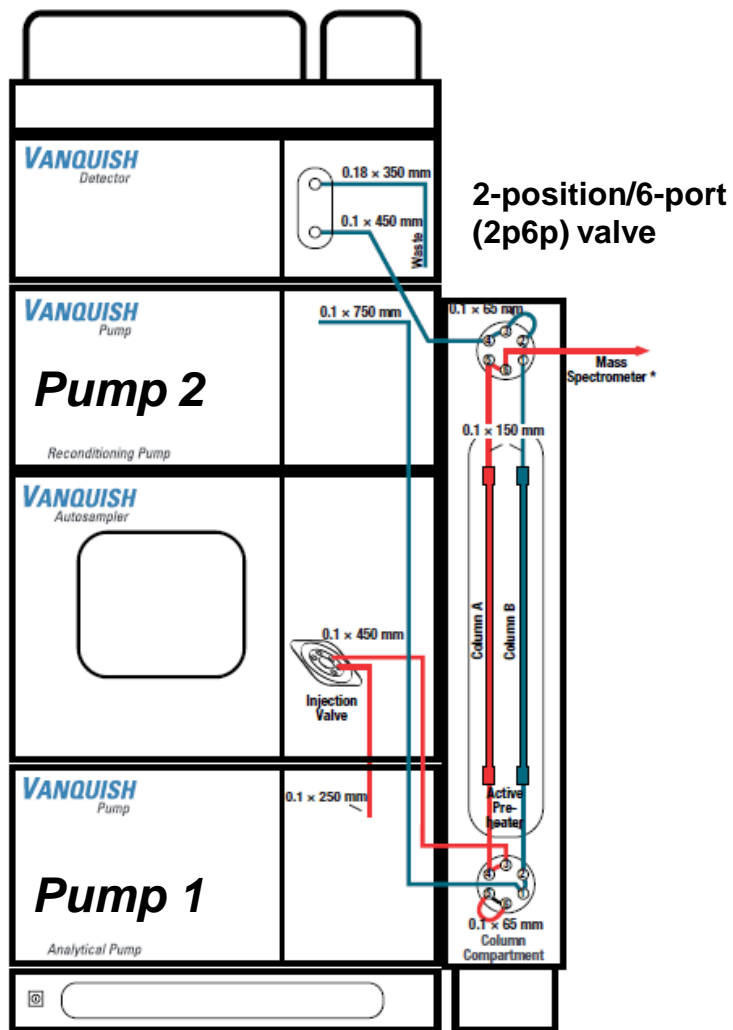
Reconditioning time may take more than 20 min until next sample injection.



Target of tandem UHPLC setups!!!

Vanquish UHPLC Tandem LC Setup to Increase Productivity

Vanquish tandem LC and LC-MS workflow



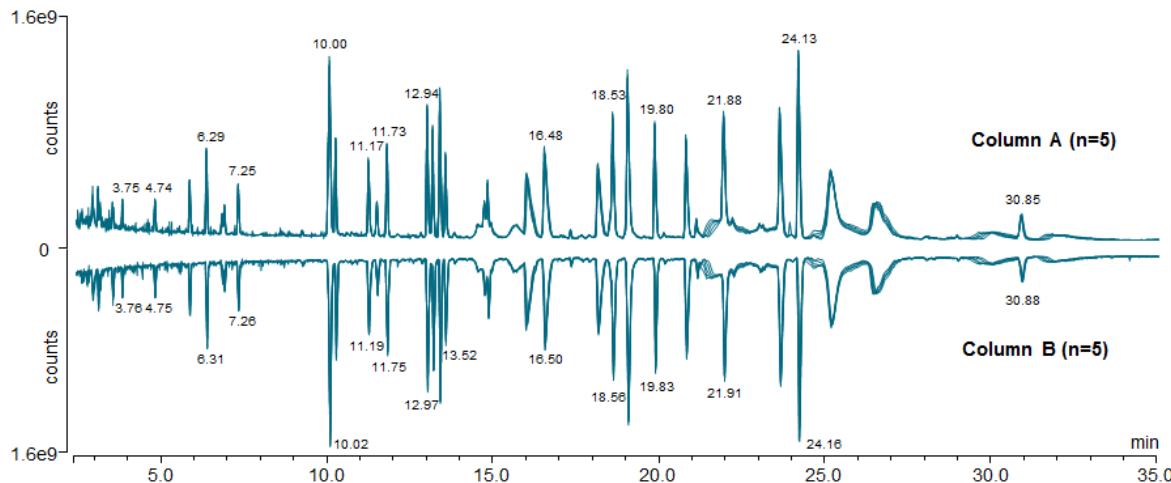
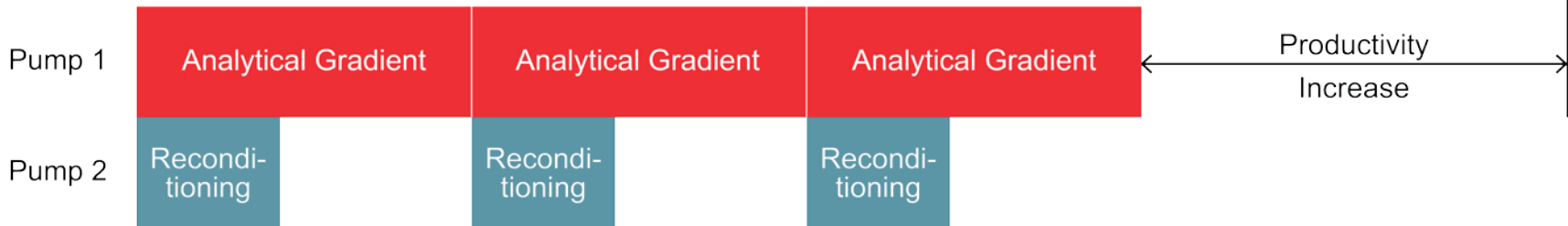
Vanquish UHPLC Tandem LC Setup to Increase Productivity

Vanquish tandem LC-MS workflow

Standard LC-MS Method



Tandem LC-MS Method



Overlay of 5 total ion current chromatograms Vanquish tandem setup:
 2x Thermo Scientific™ Acclaim™ VANQUISH™ column C18 2.1x250 mm, 2.2 μm, Lot No. 1425071; **A:** 0.1/100 FA/Water (v/v); **B:** 0.1/100 FA/Acetonitrile (v/v)
Gradient: 1-45% B in 40 min, 60 °C; 400 μL/min; Injection volume 5 μL, 1 μg Infliximab SMART digest; **Reconditioning:** Two times 90% B for 5 min, 24 min equilibration at 1% B;
Detection: Q Exactive HF MS, R=15k, mass range 140-2000; UV detection@214 nm

Summary of the Webinar

1. With an UHPLC system you can focus on peak capacity and sample throughput.
2. The system gradient delay volume (GDV) is the crucial parameter for your sample throughput.
3. The operation principle of the pump and the pump mixer configurations contribute significantly to the GDV of your UHPLC system.
4. Low pressure gradient (LPG) or high pressure gradient (HPG) pump technology is chosen in alignment with your application requirements.
5. With HPG pump technology you can achieve maximum sample throughput.
6. With LPG pump technology you can use ternary or quaternary gradients and advanced column chemistries for e.g. method development.
7. For LC-MS applications with very long run times, you can increase your productivity with tandem UHPLC system setups.

Thank you very much for your attention!



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.