

# Thermo Fisher S C I E N T I F I C

# Tips and Tricks for HPLC and UHPLC

Jan Pettersson Nordics Sales Support Specialist Chromatography Thermo Fisher Scientific, Hägersten/Sweden

### Overview



Thermo Scientific™ Vanquish™ UHPLC system



Thermo Scientific™ UltiMate™ 3000 UHPLC system



### Common Recommendations

## **Mobile Phase**







#### Common Problems: Mobile Phase

### Solvent compatibility

- Try to use pre-mixed solvents
  - Add 5-10% of organic eluent to the aqueous eluent
  - Add 5-10% aqueous eluent to the organic eluent
  - Avoids local crystallization in the pump (with buffers)
- Eluents with salt buffers
  - Change eluents with salt buffers regularly
  - Filtrate buffers
  - Use water with 18,2M Ohm AND <5ppb TOC</li>



- The quality of the eluent is very important to keep the noise as low as possible.
- Make sure that the eluents are good by running them without injection (Sample type "Blank").
- For MS and Thermo Scientific<sup>™</sup>
   Corona<sup>™</sup> charged aerosol
   detector only use volatile
   buffers.

UV-spectra at 200–250 nm of two methanol samples (both LC/MS grade)

Optimizing and Monitoring Solvent Quality for UV-Vis Absorption, Fluorescence and Charged Aerosol Detectors

Melanie Neubauer and Holger Franz Thermo Fisher Scientific, Germering, Germany

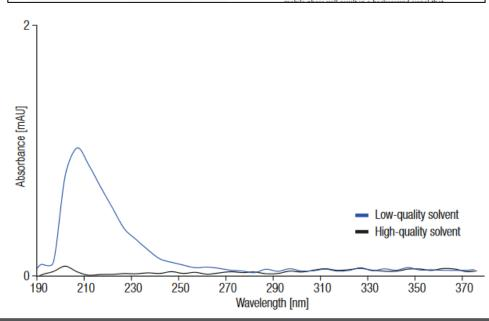
#### Key Words

Eluent Quality, Mobile Phase, UHPLC, Liquid Chromatography

#### Goal

Provide guidance on how to find out if mobile phase quality is sufficient for application specific UV-Vis, fluorescence, and charged aerosol detection requirements. Give assistance in laboratory solvent quality monitoring and solvent cost control.

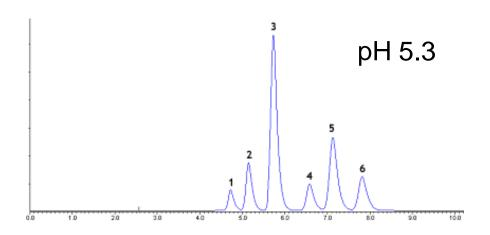
Beyond these precautions for the system mobile phase there are also detector- and application-related requirements. Optimizing the quality of mobile phase solvents can contribute to an improvement of the chromatographic or mass spectrometric properties of the analyte as well as the overall detection limits of the LC system.<sup>2</sup> To achieve lowest limits of detection (LOD) with optical detectors, the solvent should respond as little as possible to the selected wavelengths. Absorption or fluorescence of the

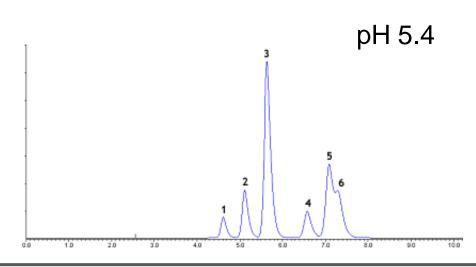




### Common Recommendations: Mobile Phase

- Symptoms
  - Peaks shift
  - Loss of resolution
- Causes
  - Mobile phase pH changes
  - Prevention
    - Use correct buffer for pH range.
    - Control pH of mobile phase.
    - Maintain buffer strength in aqueous phase.
    - Control temperature.



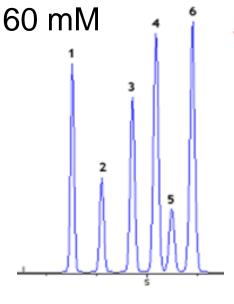




## Common Buffers and their Effective pH Range

Buffer	p <i>K</i> a	pH range		
Phosphate	2.1, 7.2, 12.3	13.1, 6.2-8.2, 11.3.13.3		
Citrate	3.1, 4.7, 6.4	2.1- 4.1, 3.7-5.7, 5.4-7.4		
Carbonate	6.1, 10.3	5.1-7.1, 9.3-11.3		
Formate	3.8	2.8-4.8		
Acetate	48	3.8-5.8		
Ammonia	9.3	8.3-10.3		
Borate	9.2	8.2-10.2		

#### Common Problems: Mobile Phase



57 mM 48.5 6

The two chromatograms illustrate the effect of a small reduction (3 mM) in buffer concentration on the separation of six food additives:

- Symptoms
  - Peaks shift
  - Loss of resolution
- Causes
  - Mobile phase buffer strength changes
  - Prevention
    - Maintain buffer strength in aqueous phase.
    - Control temperature.
    - Filter solvents rather than using vacuum degassing.

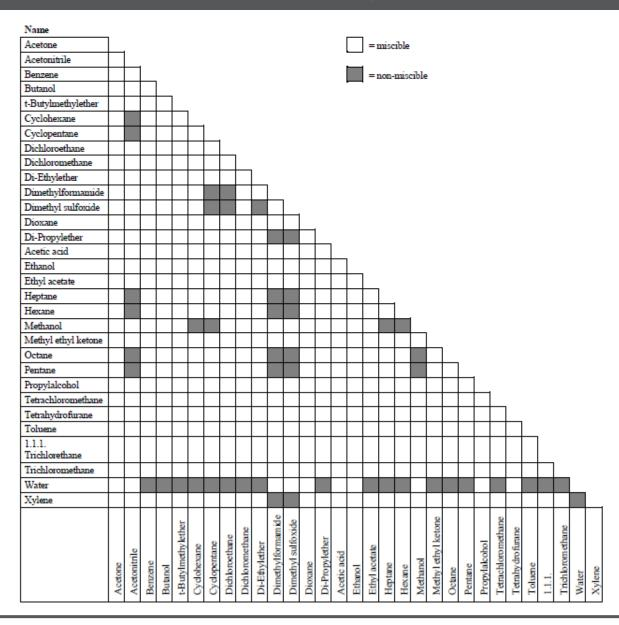
### Solvent Effects – Why Water and AcN are Popular

Solvent (nm)	Minimum wavelength
Acetonitrile	190
Water	191
Cyclohexane	195
Hexane	201
Methanol	203
Ethanol	204
Ethoxyethane	215
Dichloromethane	220
Trichloromethane	237
Tetrachloromethane	257

 Inappropriate solvent choice can cause issues with reduced sensitivity, noise and rising baselines in gradient analysis.



### Common Problems: Solvent Mixing





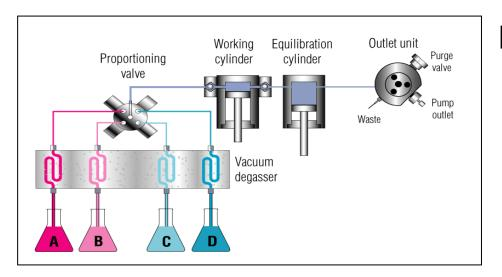
### Common Recommendations

## **The Pump**



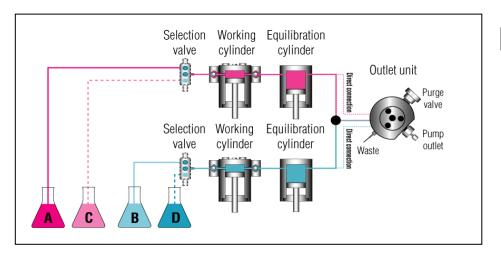


### Pump Types



### LPG pump type

- Eluent composition at the low pressure side
  Before the pump head
- Eluent is composed through proportioning valves
- Eluent segments pass working and equilibration cylinder
- Up to four solvents can be mixed



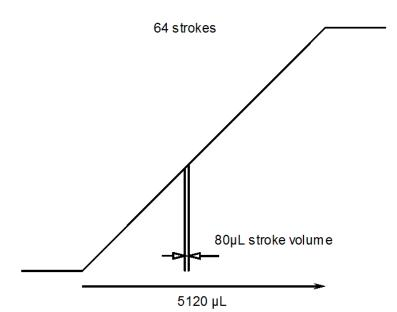
### HPG pump type

- Eluent composition at the high pressure side
  => Behind the pump head
- Pure solvents pass working and equilibration cylinder
- Eluent mixture is prepared with two pump blocks
  - => Binary eluent mixture only

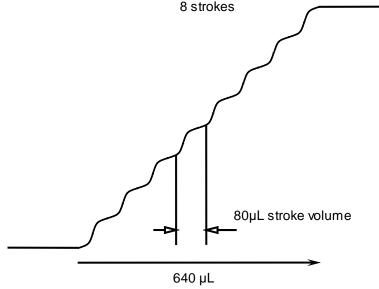


#### LPG and Ballistic Gradients

 Speed of ballistic gradients is limited by the composition change per stroke. This should not higher be than 2.0% per stroke.



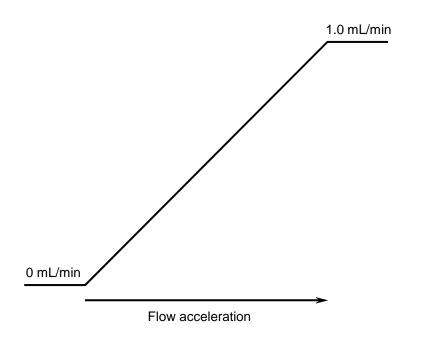
5.0 min gradient time @ 1 mL/min Gradient volume =  $5120 \,\mu$ L Piston strokes for gradient = 64 For 0-50% gradient the composition changer per stroke is 0.8%



38 s gradient time @ 1 mL/min Gradient volume = 640  $\mu$ L Piston strokes for gradient = 8 For 0-50% gradient the composition changer per stroke is 6.3%

#### **HPG** and Ballistic Gradients

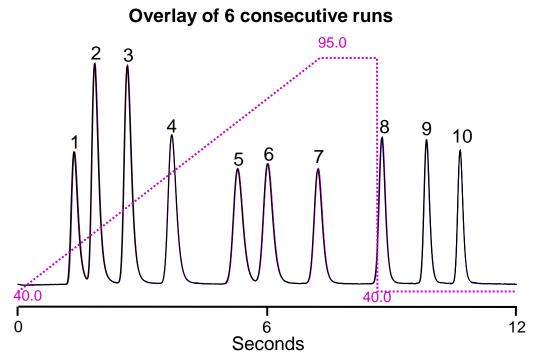
 Speed of ballistic gradients limited by the flow acceleration and deceleration of the pump



Ballistic gradient formation via ultra-precise acceleration and deceleration with an internal resolution of 125 Hz

→ Ballistic gradients faster than a second are theoretically possible

#### 10 Peaks in 10 Seconds with a Ballistic Gradient



#### **Test conditions**

Column: C18, 30 x 2.1 mm,

 $1.8 \mu m$ 

Eluents: A: Water

B: Acetonitrile

Flow: 3.70 mL/min @ 725 bar

(10,500 psi)

Temperature: 100 ° C

Inj. volume: 1 μL

Test mixture: Uracil and

9 alkylphenones

Resolution (Critical peak pair): 1.7

	Peak number									
	1	2	3	4	5	6	7	8	9	10
Retention time RSD [%]	0.76	0.41	0.29	0.14	0.15	0.14	0.11	0.09	0.08	0.05
Retention time SD [ms]	8.54	8.81	8.61	9.14	17.00	18.34	16.26	15.25	11.09	10.97





### The Comprehensive SpinFlow Mixer Portfolio

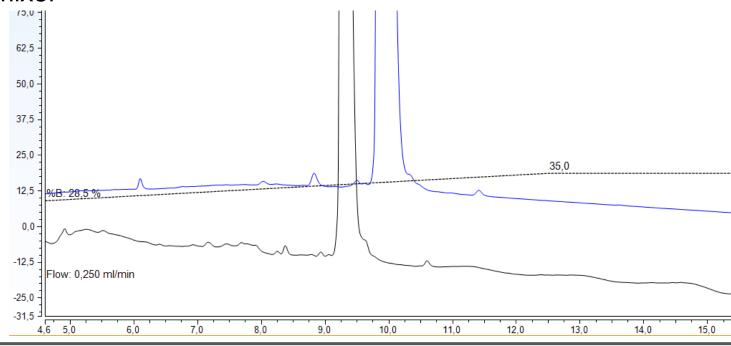
 Range of static mixers suitable from low-GDV LC/MS application to high-sensitive TFA application



Thermo Scientific™ SpinFlow™ technology

#### **Effects of Different Mixers**

- These samples were run with the same Vanquish Horizon system, the only change was the size of the mixer
- Flow 0,250µl/min
- A: (0.1% TFA, 50 mM NaCl):MeCN 95:5
- B: (0.1% TFA, 50 mM NaCl):MeCN 30:70
  - Black 10µl mixer
  - Blue 350µl mixer



### Easy Tunable for Optimum Application Performance

- For fast separations where the mixing ripple does not interfere with the detection (e.g., Corona charged aerosol detector or MS detectors), use the low mixer volumes (35 μL, 100 μL).
- Use the medium sized mixers (200 μL, 400 μL) as the best balance between fast separation and low mixing ripple in UV detection.
- For highest sensitivity and when mixing ripples interfere with the detection (e.g., due to use of UV-absorbing solvents), use a larger mixer volume (400 μL, 800 μL).
- For UV-absorbing solvent additives that amplify the mixing ripples by interaction with the stationary phase (e.g., TFA application), use for highest sensitivity the largest mixer volumes (800 μL, 1550 μL).



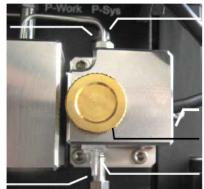
### Practical Use of a HPLC Pump

Prime the pump – but be sure that there is eluent in the pump

 Manually or with the autosampler

Pump Type	Purge Flow	Purge Time	
Analytical pump	3 mL/min	5 min	
Micro pump	2 mL/min	5 min	
Semipreparative pump	30 mL/min	5 min	

Connection port for the right pump head (if available)



Connection port for the left pump head

Pressure transducer for the system pressure

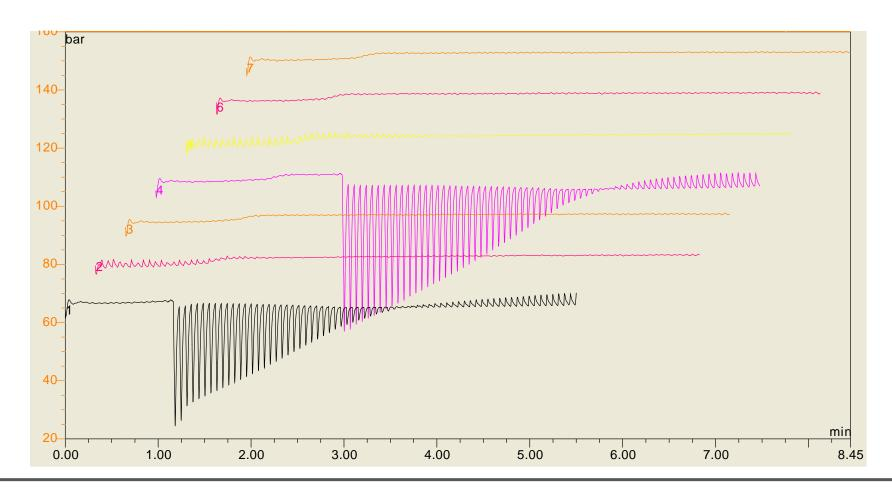
Purge valve knob

Connection port (capillary mixer/inline filter/pulse damper)

Purge outlet nozzle

### Prime and Degas

- Air bubble stuck in the pump
- Prime the pump
- Degas the eluents 5 min in an ultrasonic water bath

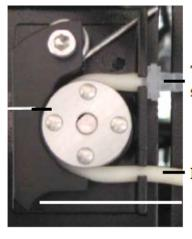


### Practical Use of a HPLC Pump

- Use eluent filters if necessary
- Use a degasser it might be a good idea to degas in an ultrasonic water bath for 5 min

Detector of the seal wash system

- Seal wash helps the pump seal to survive
- If the pump is out of wash solution it should not start
- If the pump starts to leak the Thermo Scientific<sup>™</sup> Chromeleon<sup>™</sup> chromatography data software will give a warning



Tubing from seal wash reservoir

PharMed tubing Lever





Securing clip

Rotor

### Common Recommendations

# Autosampler







### Autosamplers

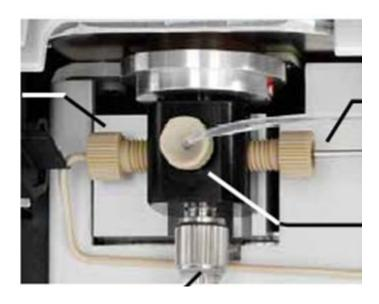
- Modern autosamplers should be considered as pipetting robots
- Other than injecting the sample they can:
  - Dilute samples
  - Do pre-column derivatization
  - Spike the samples

#### Autosamplers

- Before you run samples flush the syringe
  - Even a tiny air bubble ruins the performance
  - Autosamplers need a transport liquid
  - Usually eluent A is connected to the sampler and the pump
  - It is possible to use a separate bottle of transport liquid
  - The transport liquid is used to wash the needle







#### Important Parameters

#### DrawSpeed

- Defines the speed at which the sample is drawn by the syringe
- In analytical range (5 100 μL) a draw cycle should normally take
  3 4 seconds.

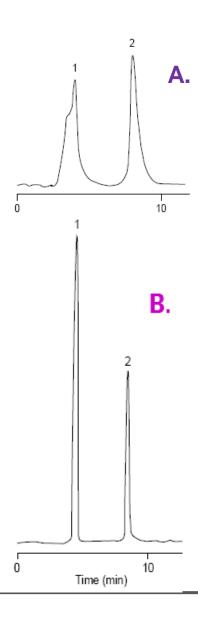
At lower volumes, a draw cycle should take approximately 10 times longer

**Examples** (normal HPLC eluents and samples dissolved in starting eluent)

- => 10 μL injection volume -> Recommended draw speed: 2 3 μL/sec
- => 2 μL injection volume -> Recommended draw speed: 0.2 0.3 μL/sec
- Draw speed has to be adapted in case of samples and eluents with higher viscosity.
- Incorrect setting is frequently responsible for area precision problems.



### Sample Solvent



- Matching injection volume with solvent strength
  - Ideally the sample solvent will not affect chromatographic separation
  - If a stronger sample solvent is required, injection volumes should be kept to a minimum
  - In a strong solvent, the sample moves more quickly through the mobile phase and often split or distorted peak shapes are observed as in chromatogram A
  - The syringe is usually connected to eluent A

### Vials and Septa

- Do not overfill the vial the septum is used to clean the needle. Fill 2/3 of the volume
- If you use inserts make sure there are no airbubbles in them
- Autosamplers with PEEK needle must have caps with slit septa
- Never shake the vial. If you do there is going to be sample on the underside of the septum which gives carry over
- · Adjust the needle height according to the vial dimensions
- Rubber septas get dry and can block the injector

... And use the vial and septum only once!







### **Common Recommendations**

# **Column Compartment**

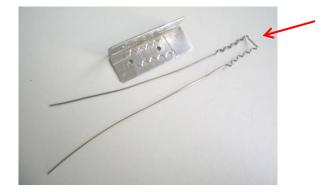


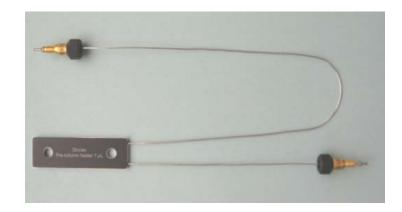


#### Pre-Column Heater

- Serpentineshaped capillary embedded in aluminum block
- Different types available

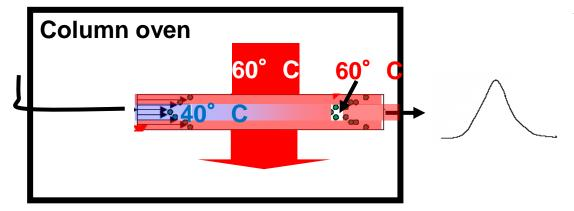
2μL, 0.12mm ID 7μL, 0.18mm ID 11μL, 0.25mm ID





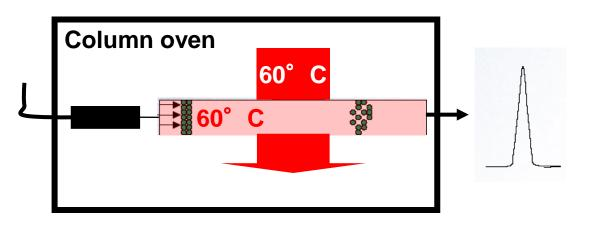


#### Pre-Column Heater

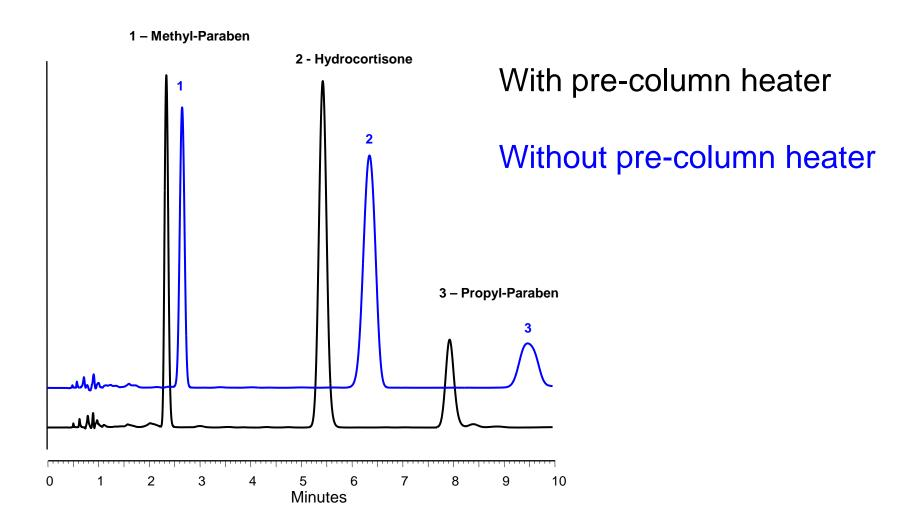


#### Without pre-heating:

- Poor peak resolution
- Peak broadening
- Peak splitting
- Extended analysis time



#### Pre-Column Heater





### Common Recommendations

### **UV-Detectors**



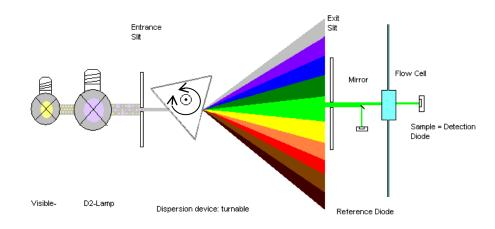




### Operating Principle Variable Wavelength Detector (VWD)

#### Forward optics design

- Only the selected wavelength passes the flow cell and is detected by the sample diode
- A part of the light beam is redirected to the reference diode
- Reference signal is not influenced by the content of the flow cell
  - 'True' reference

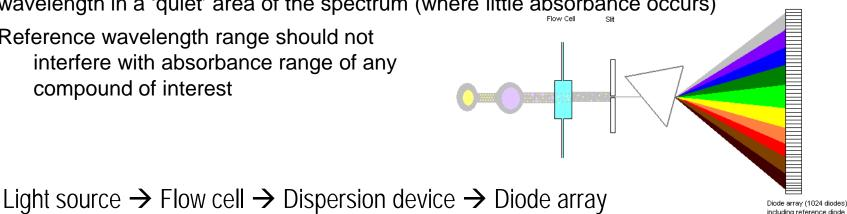


Light source → Dispersion device → Flow cell → Sample diode

#### Recommended Parameters: Reference Wavelength DAD

#### Reversed optics design

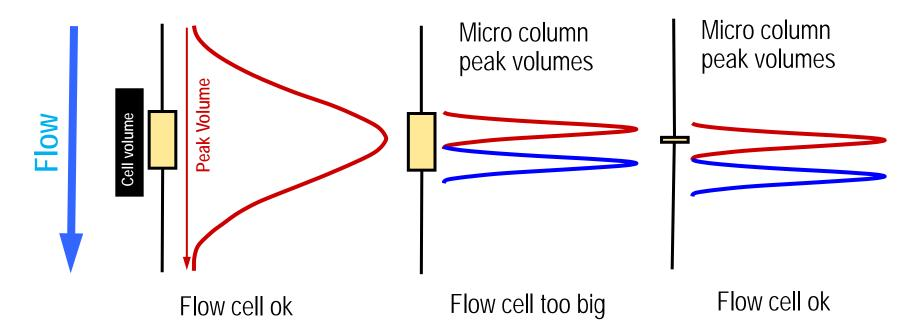
- Light beam passes the flow cell before being diffracted
  - No true reference signal can be obtained
- Instead -and with limitations-, any diode or bunch of diodes can be selected as a reference
  - If selected reference and acquisition wavelength are the same, the resulting signal would be zero (0)
  - As a consequence either don't use a reference (preferred) or select a reference wavelength in a 'quiet' area of the spectrum (where little absorbance occurs)
  - Reference wavelength range should not interfere with absorbance range of any compound of interest





#### Flow Cell Volume

- Flow cell volume depends on peak volume
  - Rule: Flow cell volume should not exceed 10% of the peak volume

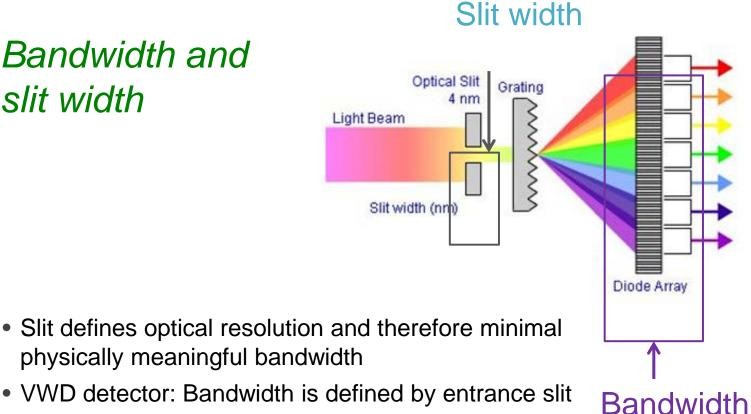


Note: Besides lamp age the light intensity is highly dependent on the installed flow cell

Smaller cell volume → Less light is passing through the flow cell

### **Diode Array Detector**

## Bandwidth and slit width



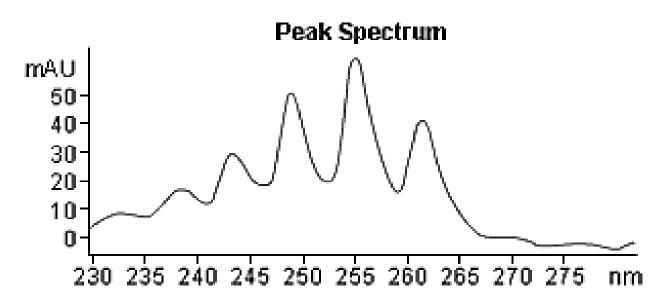
physically meaningful bandwidth

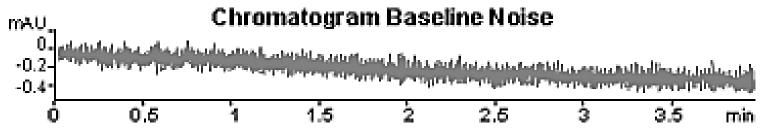
VWD detector: Bandwidth is defined by entrance slit

Slit width	Baseline noise	Spectral resolution	Bandwidth	S/N ratio	Spectral resolution	
$\downarrow$	1	<b>↑</b>	$\uparrow$	<b>↑</b>	$\downarrow$	
<b>↑</b>	$\downarrow$	$\downarrow$	$\downarrow$	$\downarrow$	<b>↑</b>	

#### Effects of Slit Width

#### Slit width - 1nm



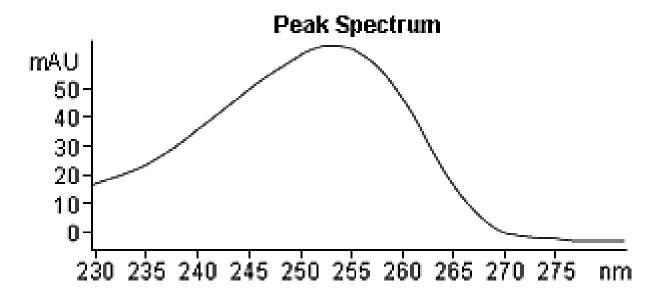


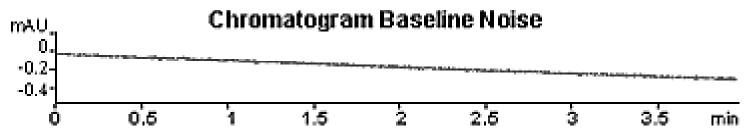
 $4 \times light > 0.5 \times noise$ 



#### Effects of Slit Width

#### Slit width - 16nm

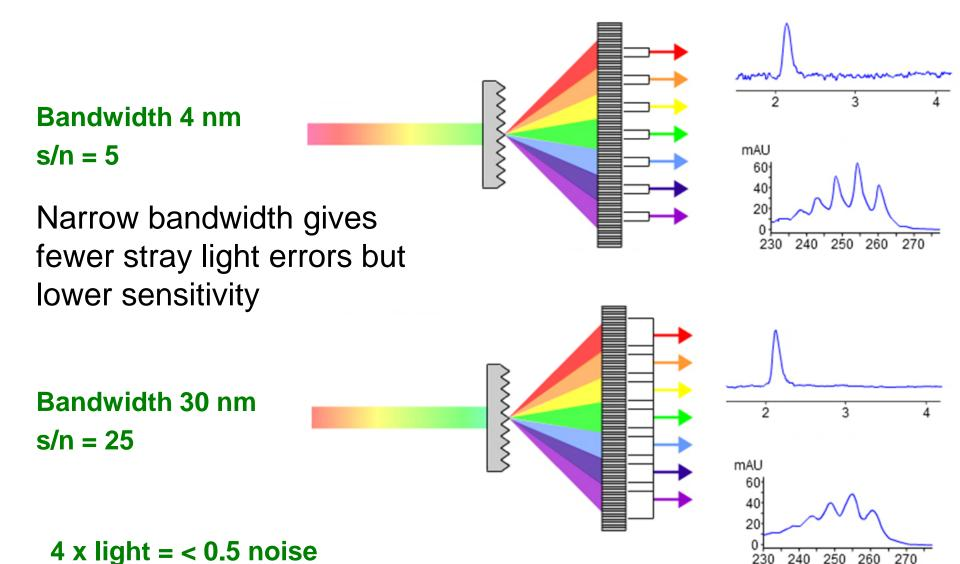




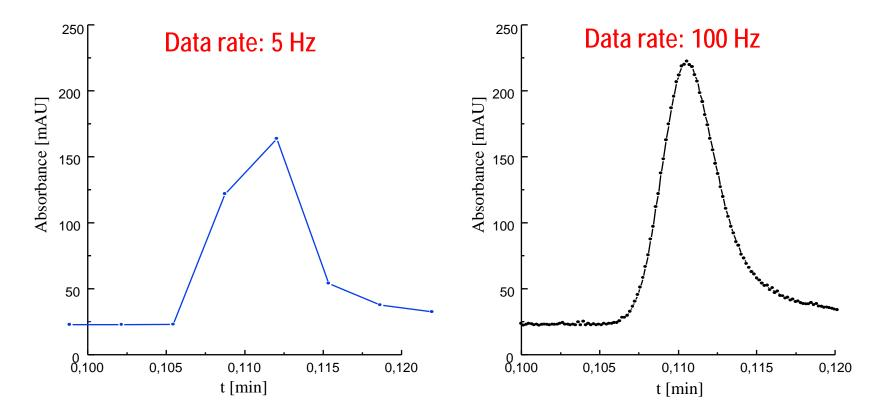
 $4 \times light > 0.5 \times noise$ 



### **Setting Bandwidth**



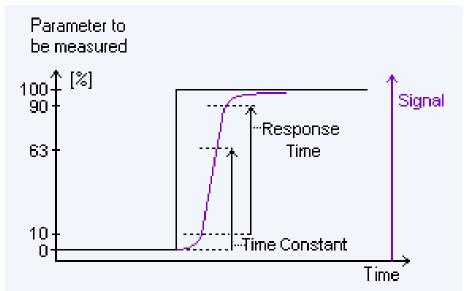
#### Recommended Parameters: Data Acquisition



- Too few data points effect peak form, reproducibility and area precision
  - A minimum of 20, ideally 30-40 data points/peak is required

#### **Time Constant**

- The parameter is a measure of how quickly the detector responds to a change in signal
- Defined as the time it takes the detectors output signal to rise from 10% of its final value to 90%

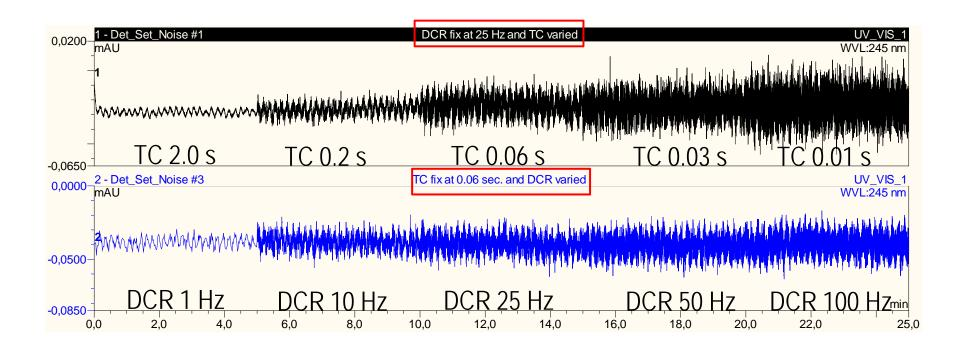


The Rise Time (Response Time) is closely releated to the time constant:

Rise time =  $2.2 \times \text{Time constant}$ 



### Recommended Parameters: Data Acquisition

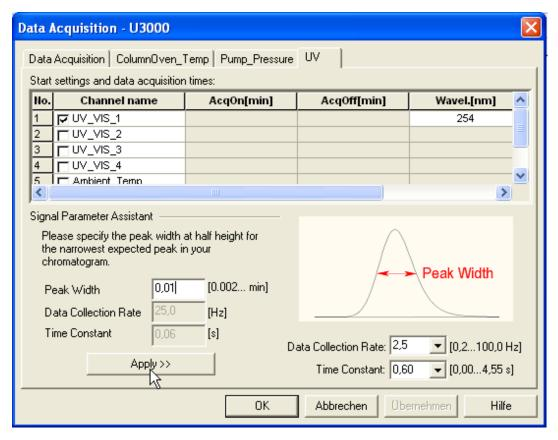


 Noise is much more influenced by time constant (TC) than by data collection rate (DCR)



#### Recommended Parameters: Data Acquisition

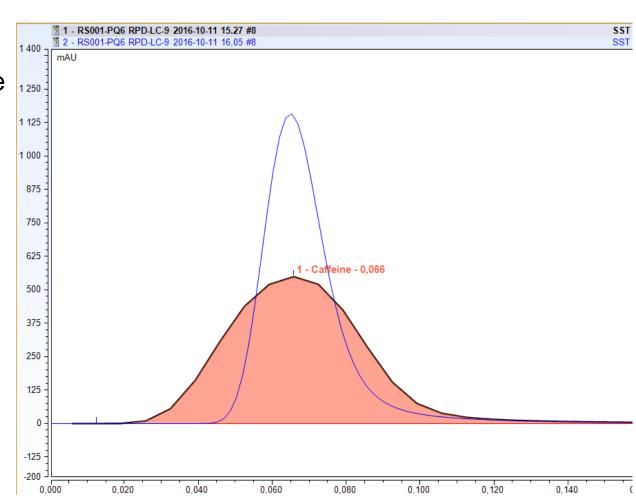
- The Program Wizard of Chromeleon has a dedicated step for setting the correct 'Data Collection Rate' and 'Time Constant'
- The internal calculation is based on the peak width at half peak height of the slimmest peak in the chromatogram





### Sampling and Rise Time

- The same instrument, back pressure loop, eluent and sample. The area is the same the peakshape is very different.
- 2,5 Hz 2s response time
- 10 Hz 0,5s response time





#### THANK YOU!

 Technical Support for Chromatography Columns and Consumables <u>www.thermoscientific.com/chromexpert</u>

Applications Library Resource <u>www.thermoscientific.com/AppsLab</u>



# 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.