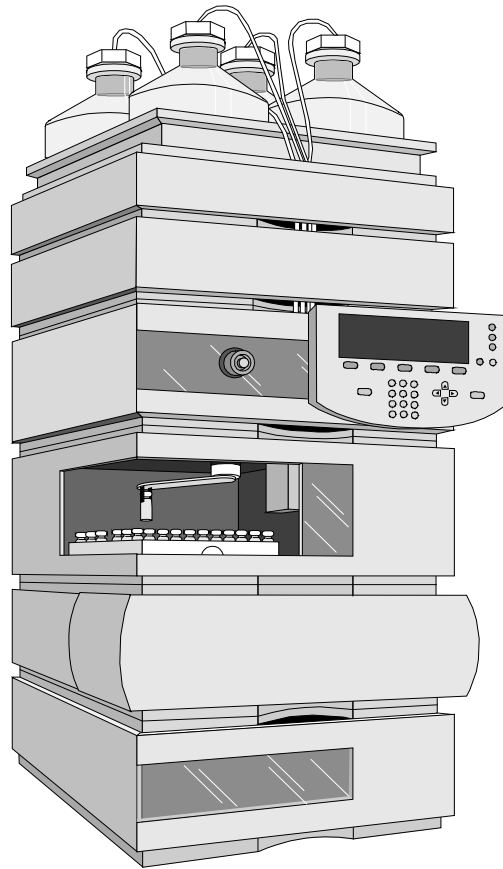


***Agilent Technologies
1200/1100 Series HPLC
Troubleshooting and Maintenance
Seminar***

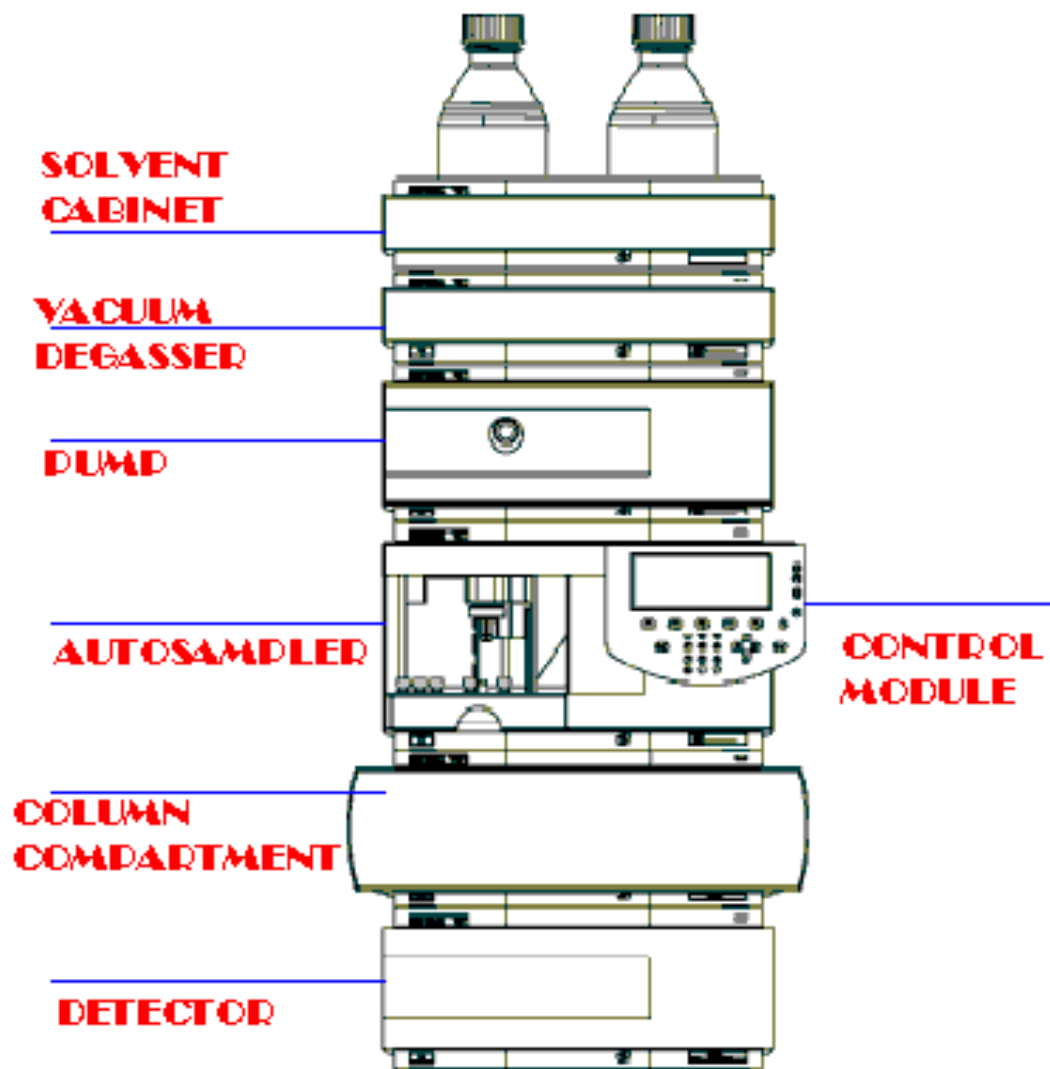


***Patrick Cronan & Sue D'Antonio
LC Applications Scientists***

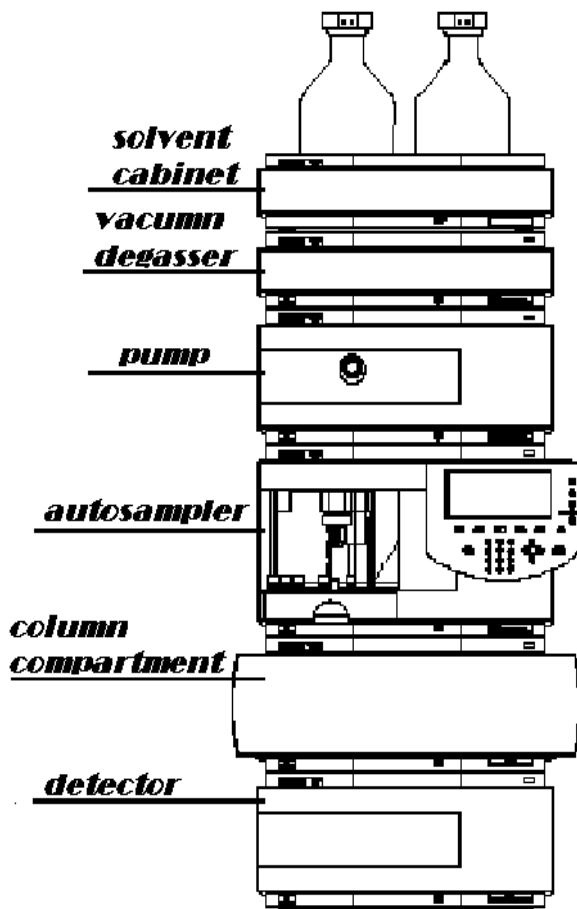
Agilent Technologies HPLC Systems



- All stackable modules for small bench space
- Single system control module
- Front access for customer maintenance
- Ergonomic tubing organization for lowest delay volume and bandspreading
- Single CAN cable connection
- All modules have RS 232, HP IB, Start/Stop, CAN



THE STACK



Flow connections in the stack:
Example setup with **0.17mm ID green capillaries**

Solvent bottles - degasser:

G1311-60003 (bottle-head assembly, PTFE-tubings)

Degasser - pump:

G1322-67300 (PTFE-tubings)

Pump - autosampler:

G1312-67305 (SST, green)

**control
module**

Autosampler - column compartment:

G1313-87305 (SST, green)

Column compartment - column:

G1316-87300 (SST, green)

Column - detector:

DAD G1315-87311 (SST, coated)

VWD 5062-8522 (PEEK)

Detector - waste:

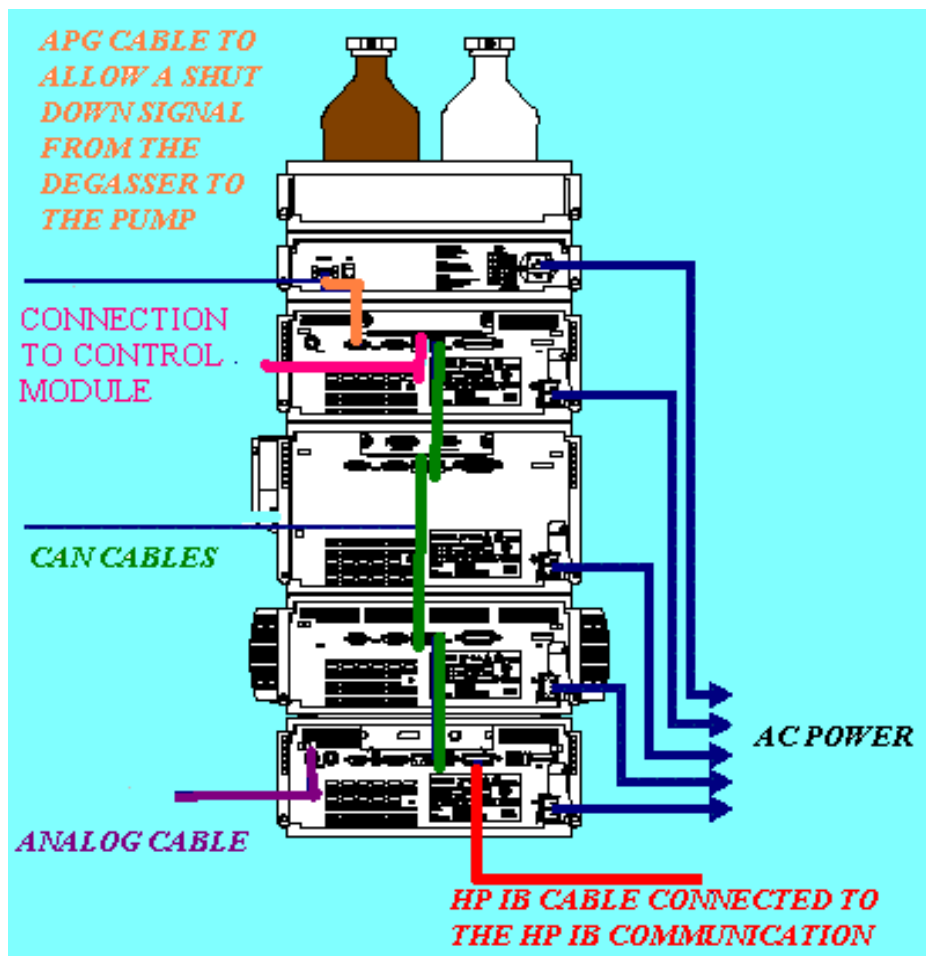
DAD 0890-1713 (PTFE, wide bore)

VWD 5062-8535 (PEEK)

5062-2463 (corrugated waste tubing, reorder pack)



REAR CONNECTIONS



CAN CABLES BETWEEN
MODULES

*CAN CONNECTION TO
THE CONTROLLER*

ANALOG OUT PUT FOR
DATA
COMMUNICATION TO
A NON HP DATA
SYSTEM

*HP IB CABLE TO
COMPUTER*

*LAN Connections on
Newer Systems*

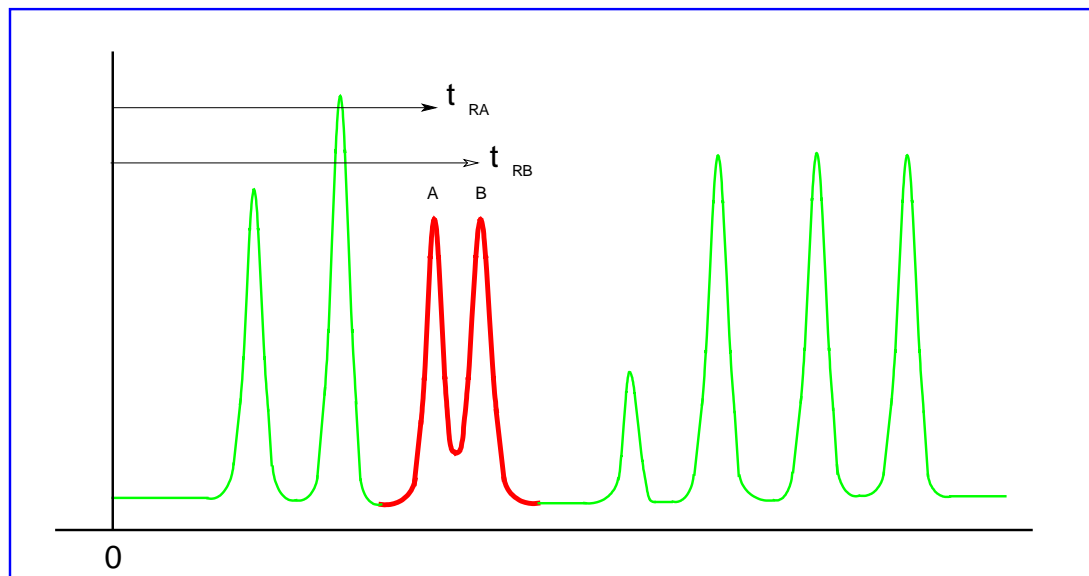
What are Chromatographers Looking for?

Better performance

- ▶ Baseline separation in shortest time
- ▶ Repeatability of results
- ▶ Accuracy of results
- ▶ Sensitive detection
- ▶ Standard, narrow bore and capillary column capability

The Goal of Separation

-Resolution Between Sample Components



$$R=2\left(\frac{t_{RB}-t_{RA}}{W_A+W_B}\right)$$

$$R=1.176\left(\frac{t_{RB}-t_{RA}}{W_{1/2A}+W_{1/2B}}\right)$$

R - resolution

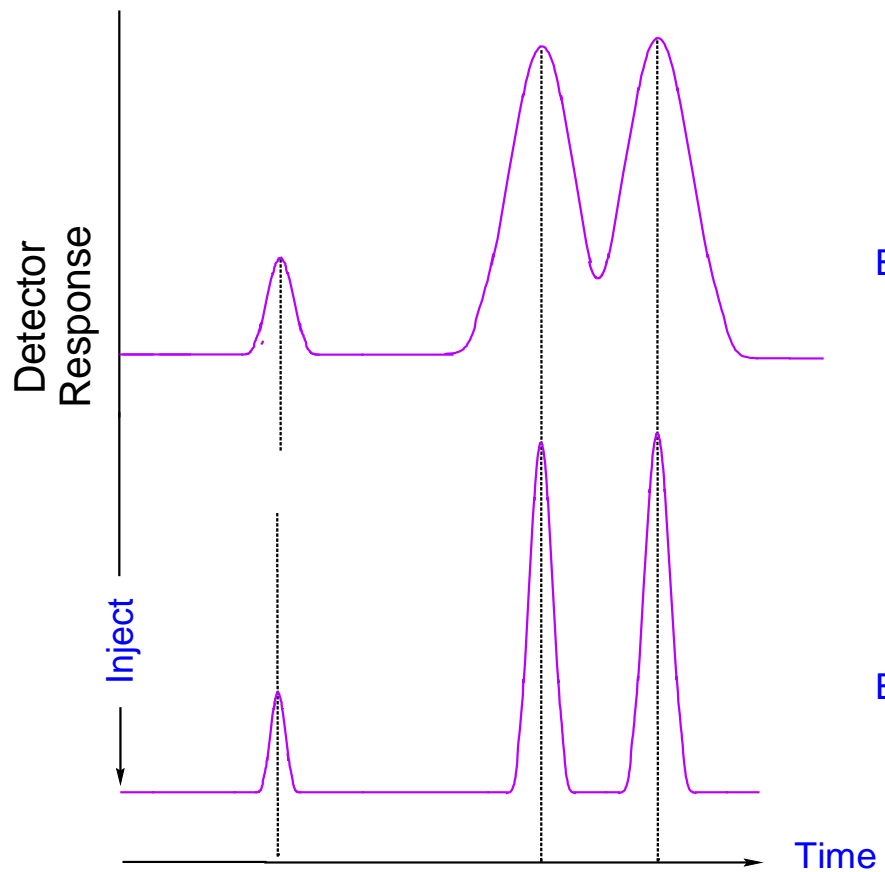
t - retention time of component B

t - retention time of component A

w - width at base of peak

w - width at half-height

The Goal of Separation -Resolution Between Sample Components



$$R = \underbrace{1/4 \sqrt{N}}_{\text{Efficiency}} \times \underbrace{\frac{\alpha - 1}{\alpha}}_{\text{Selectivity}} \times \underbrace{\frac{k'}{1 + k'}}_{\text{Capacity}}$$

Resolution

R_s , defined as the amount of separation between two adjacent peaks, is given by:

$$R_s = (1/4) (\alpha - 1) (N)^{1/2} [k/(1 + k)]$$

where **k** is the average value for the two peaks.

Performance Characteristics of an HPLC System

Influenced by one module...

Flow: accuracy, precision
Composition: accuracy, precision

Injection volume precision
Linearity, dynamic range
Carry over

Column temperature accuracy
Column temperature precision

Wavelength: accuracy, precision
Signal linearity

Spectral resolution (DAD only)

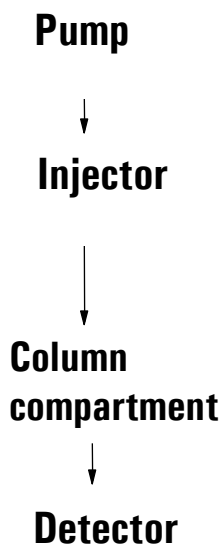
Influenced by several modules...

Repeatability of retention times
Delay volume

Repeatability of peak areas
Dead volume

Peak elution order

Baseline: noise, drift and wander



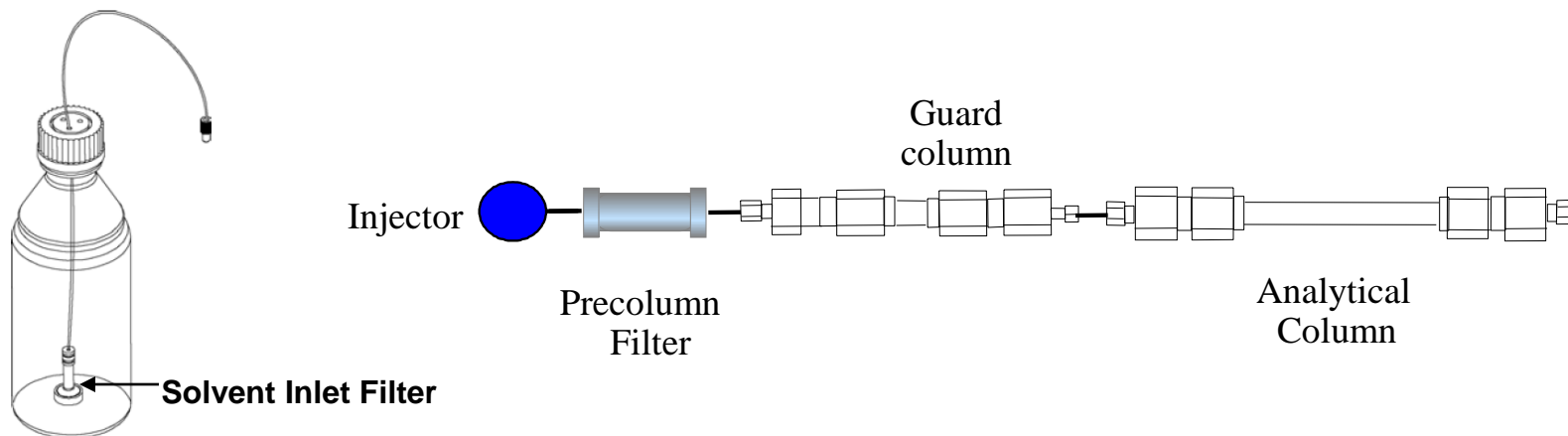
Instrument Status Indicator

The instrument status indicator indicates one of four possible instrument conditions:

- **When the status indicator is OFF** (and power switch light is on), the quaternary pump is in a prerun condition, and is ready to begin an analysis.
- **A green status indicator**, indicates the quaternary pump is performing an analysis (run mode).
- **A yellow indicator** indicates a not-ready condition. The quaternary pump is in a not-ready state when it is waiting for a specific condition to be reached or completed (for example, immediately after changing a setpoint), or while a self-test procedure is running.
- **An error condition is indicated when the status indicator is red.** An error condition indicates the quaternary pump has detected an internal problem which affects correct operation of the quaternary pump. Usually, an error condition requires attention (for example, leak, defective internal components). An error condition always interrupts the analysis.



Solvent Filters



Solvent Inlet Filter

Stainless Steel or glass with 10 micron porosity.

Removes particulates from solvent.

Precolumn Filter

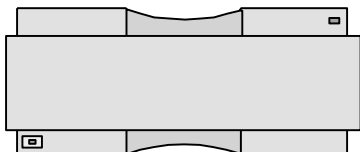
Used between the injector and guard column.

2 to 0.5 micron

Removes particulates from sample and autosampler wear debris.

Must be well designed to prevent dispersion.

Agilent 1200 Series Vacuum Degasser



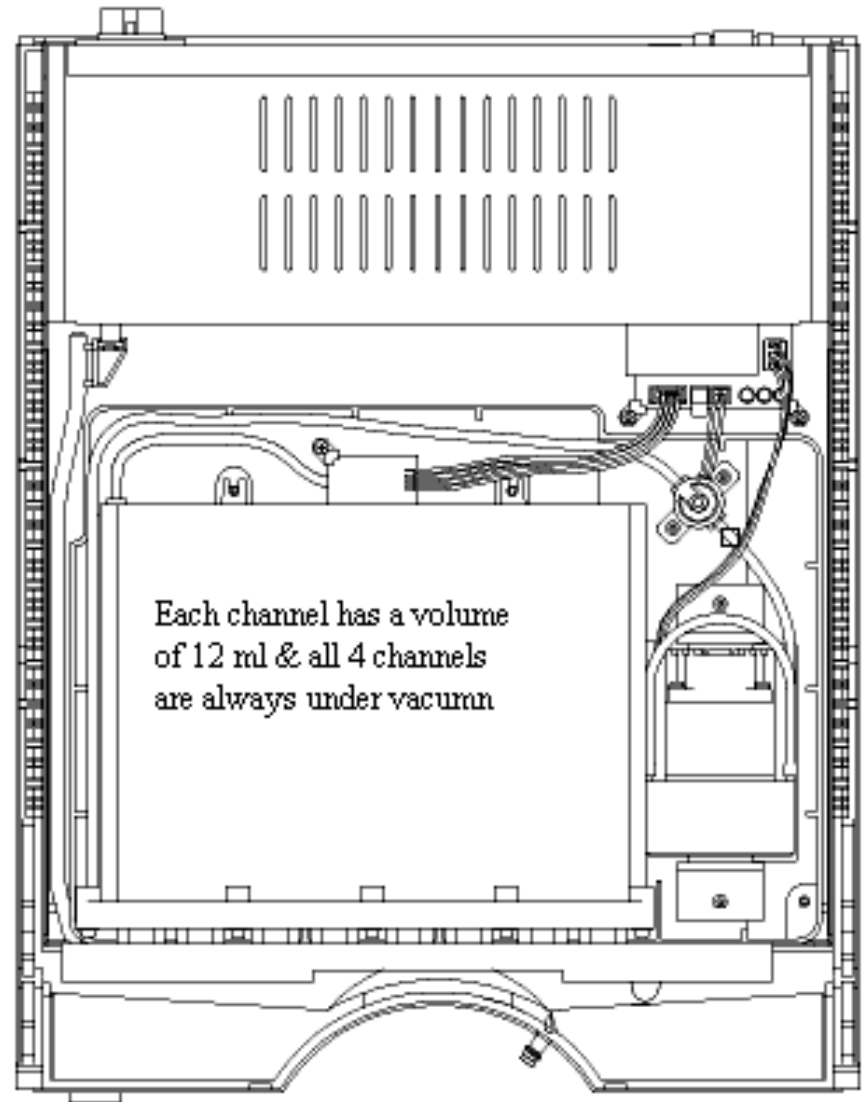
Recommended for ...

- reliable performance with gradient pumps using low pressure mixing
- improved performance with all pump designs at low flowrates
- use with detectors requiring oxygen-free mobile phase

Features/Benefits

- Convenient and cost effective alternative to helium sparging
- High degassing efficiency for trouble-free system operation (<1.5 ppm oxygen at 10 ml/min)
- Low internal volume (12 ml) for fast solvent changeover
- Up to 4 channels for highest versatility
- Stackable with other Agilent 1200 Series modules

The Vacuum Degasser

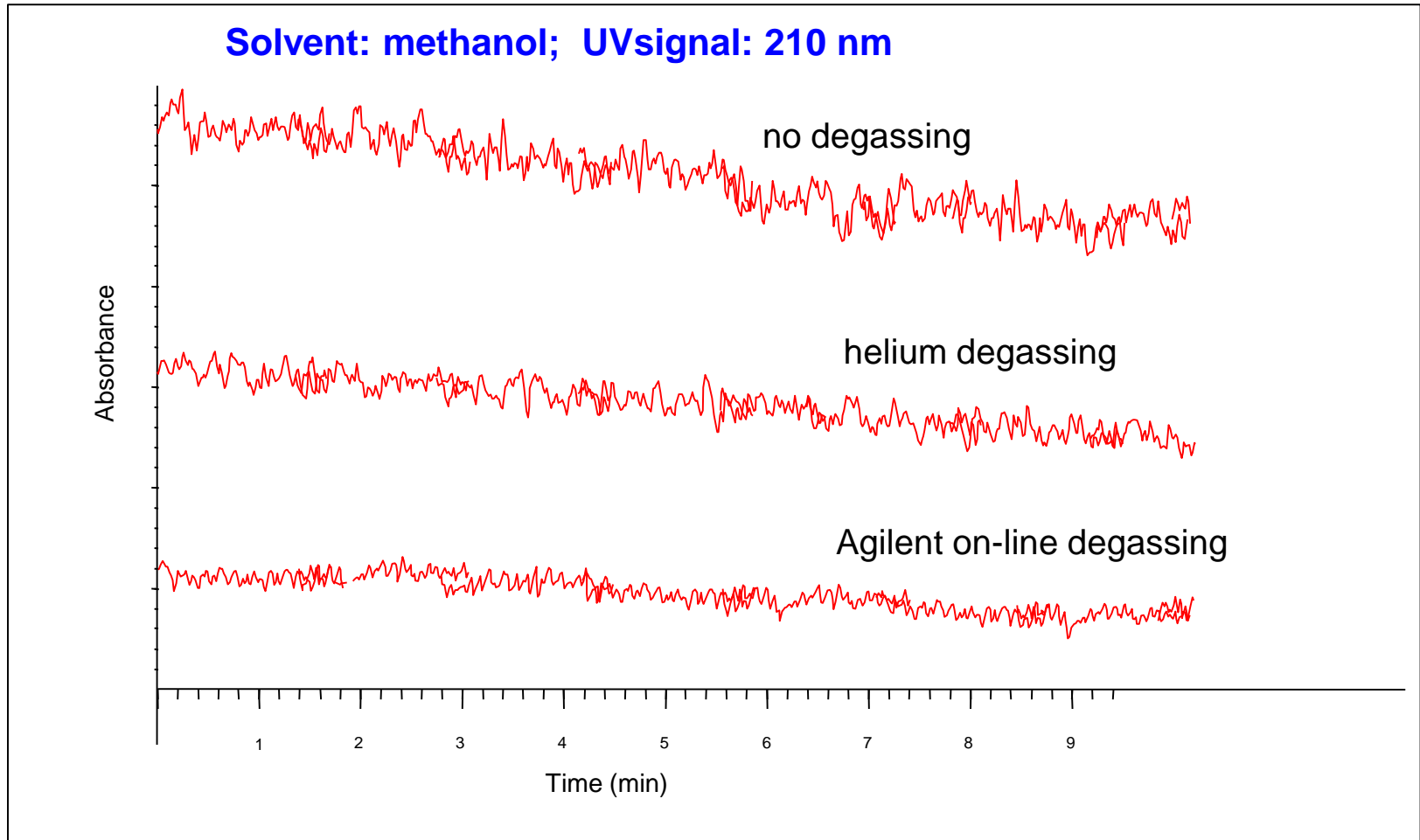


Priming and Purging the System

- The system can be primed either by drawing solvent through the degasser with a syringe or by pumping with the pump.
- Priming the system with a syringe is recommended, when:
 - vacuum degasser or connected tubings are used for the first time or vacuum tubes are empty or
 - changing to solvents that are immiscible with the solvent currently in the vacuum tubes.
- Priming the system by using the pump at high flow rate (3–5 ml/min) is recommended, when:
 - pumping system was turned off for a length of time (for example, overnight) and if volatile solvent mixtures are used, or
 - solvents have been changed.

Agilent 1200 Series On-line Degasser

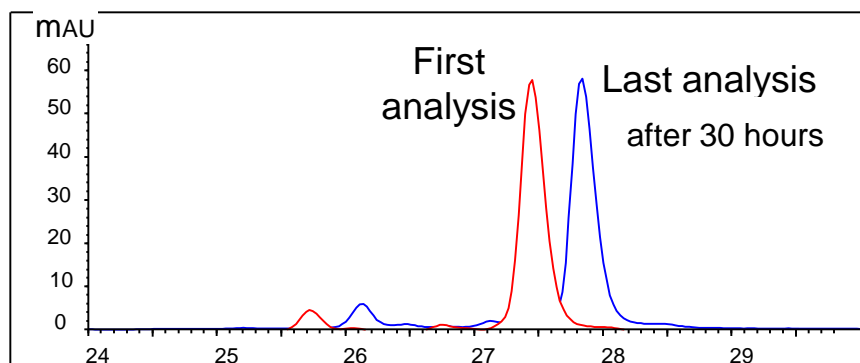
-Influence on Detector Baseline



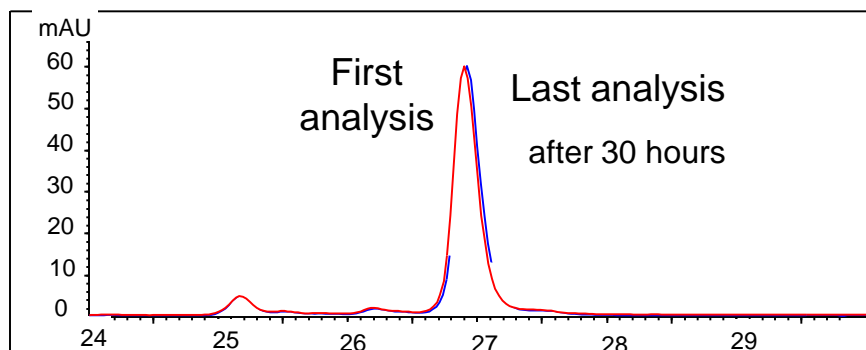
Agilent 1200 Series On-line Degasser

-Influence on Reproducibility

Lysozyme Analysis



**Helium
Degassing**



**On-line Vacuum
Degassing**

Pump

- Important characteristics

- ▶ Common to isocratic and gradient pumps

- Flow accuracy
- Flow precision
- Pressure pulsation

- ▶ Gradient pumps only

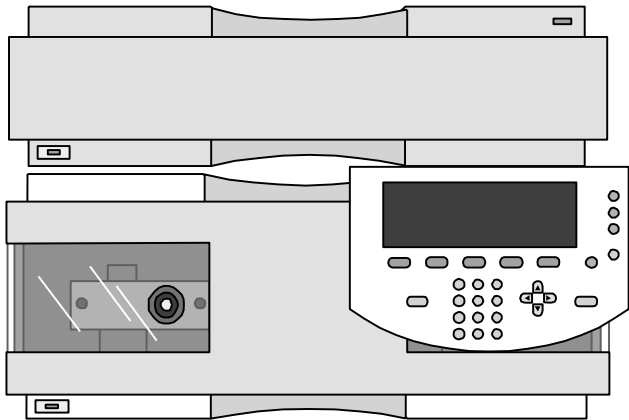
- Delay volume in low and high pressure mixing
- Composition accuracy
- Composition precision

- Influence on...

- Retention time and peak area precision (system to system)
- Retention time and peak area precision (within one system)
- Baseline noise

- Gradient shape and precision
- Retention time and peak area precision (system to system)
- Retention time and peak area precision (within one system)

The Agilent 1260/1200/1100 Series Pumps

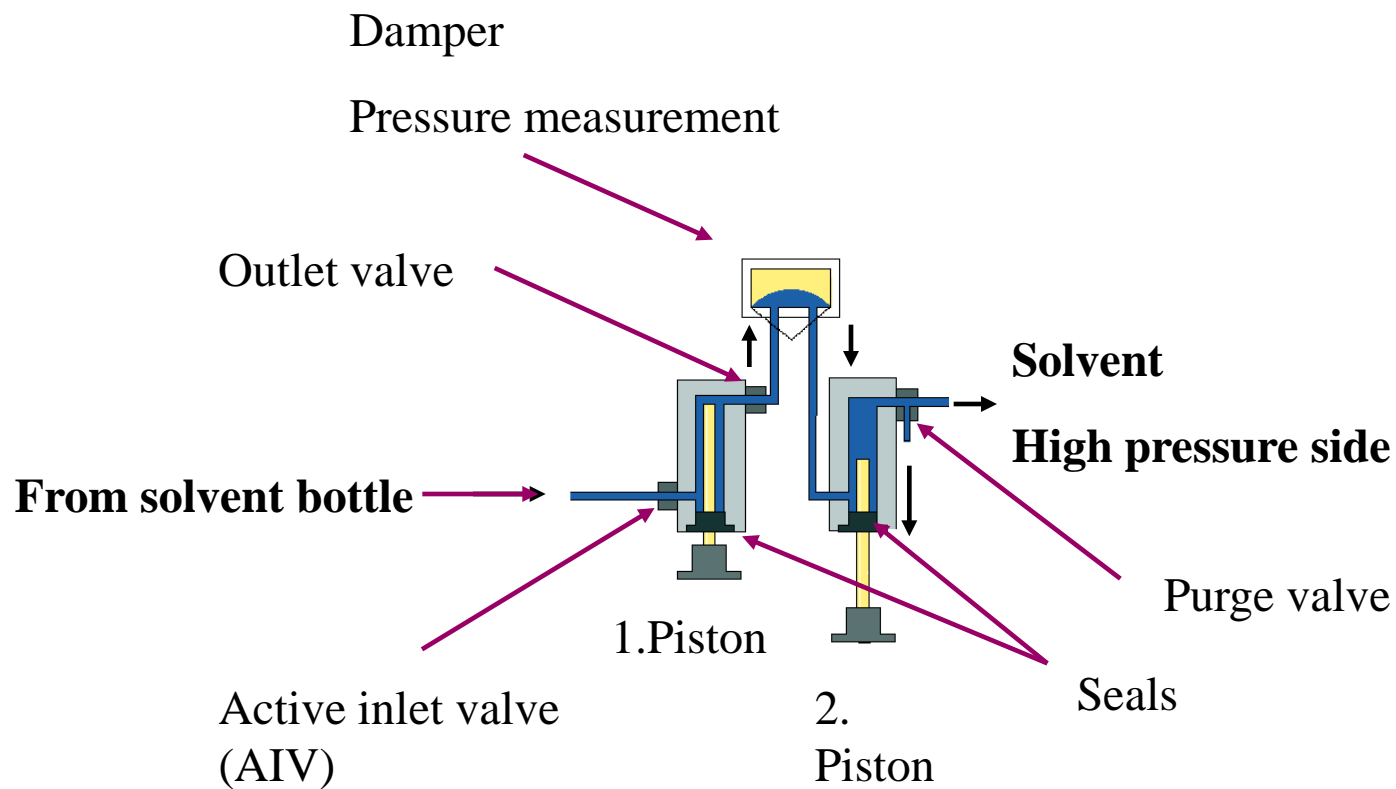


Common Features

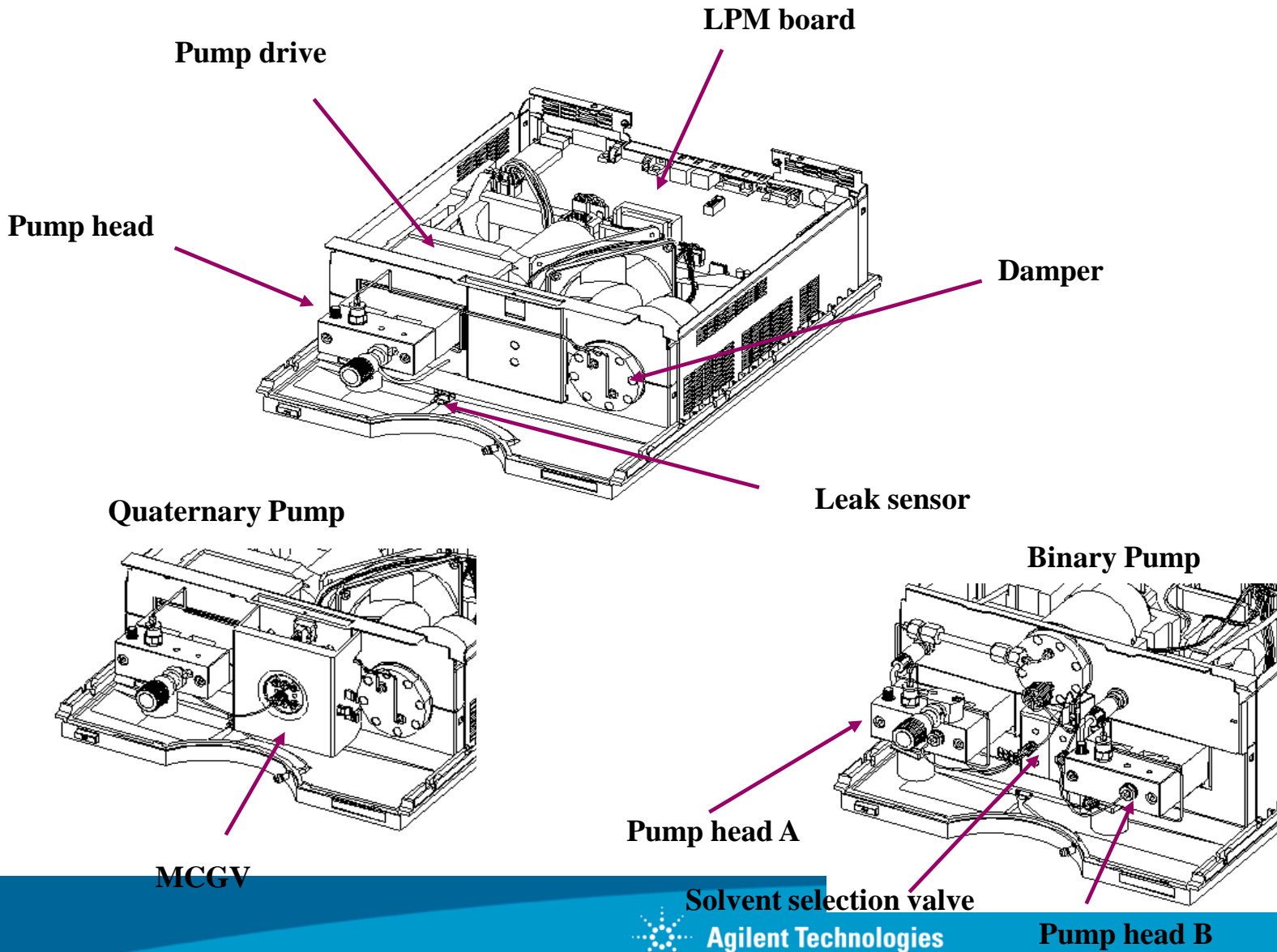
- Complete with solvent bottle, filter, cabinet, purge valve and tubing's for fast start up
- Dual-piston, variable stroke volume, and pulse dampener for pulseless flow
- Automatic stroke volume adjustment for excellent mixing and noise-free baselines
- Improved valves for longer life and lower replacement cost
- Optional seal wash for trouble-free operation with high salt mobile phases
- Easy to maintain and repair



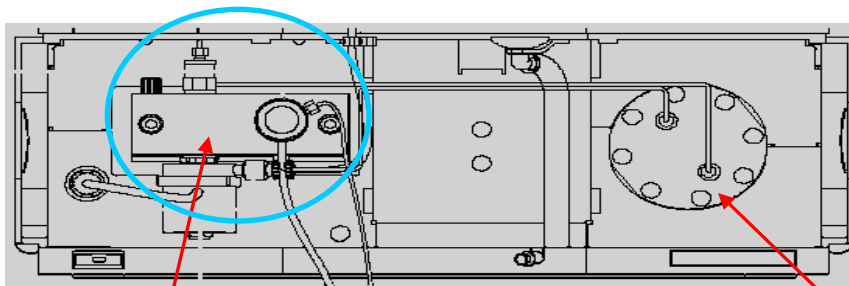
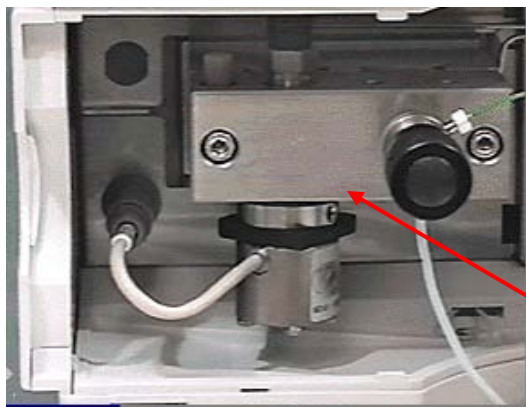
Operating Principle of the Dual Piston Pump



Pump - Main Components



Isocratic Pump

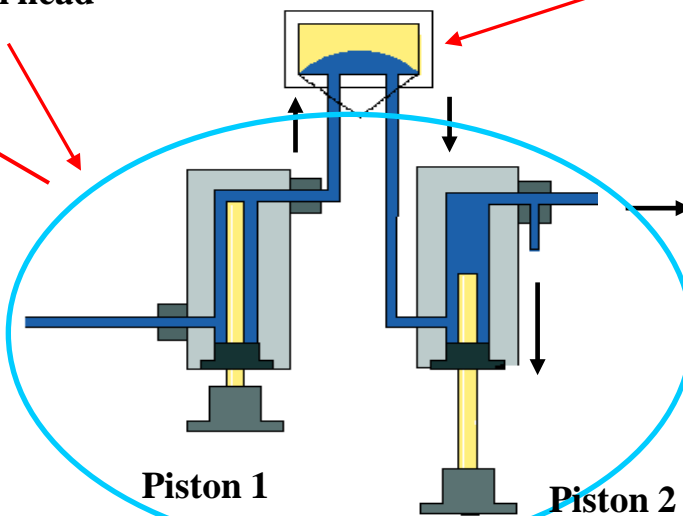


Piston head

Damper

(Pressure measurement)

Working Principle



Piston 1

Piston 2

Maintenance areas:

Seals,

AIV,

Outlet valve,

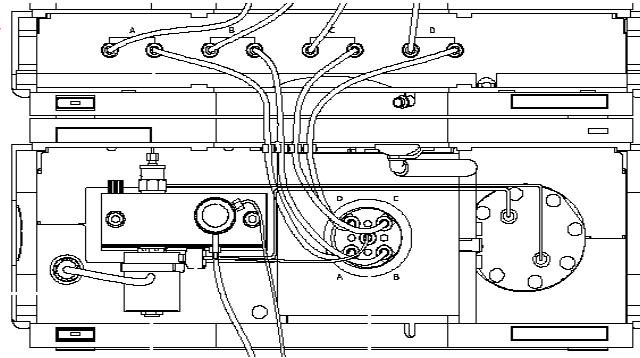
Purge valve,

Pistons.

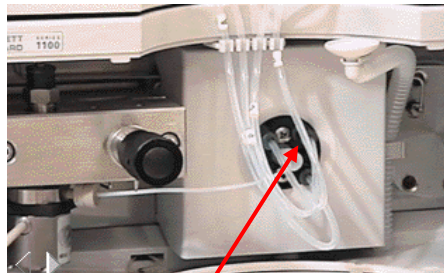
Quaternary Pump with Vacuum Degasser



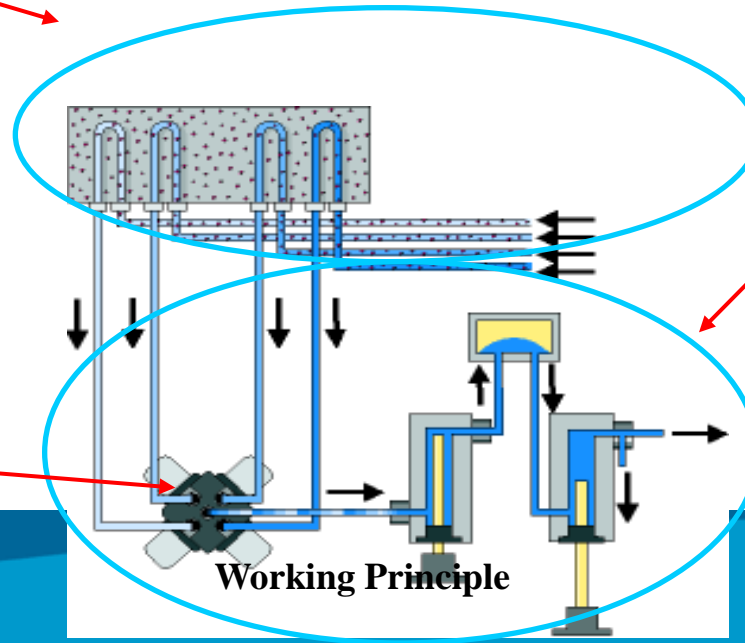
Vacuum degasser



Quaternary Pump



MCGV

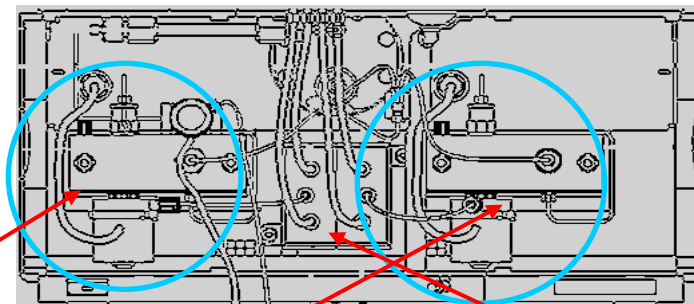


Working Principle

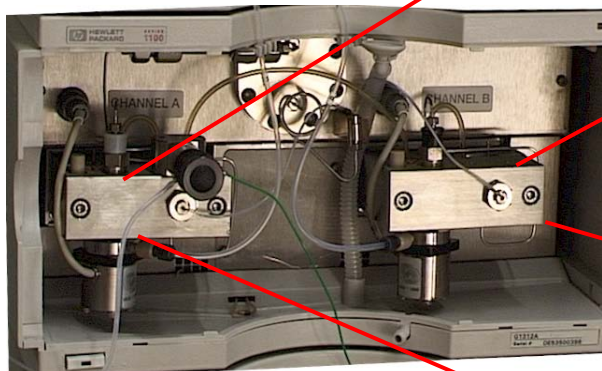
Maintenance areas:

- Seals,
- AIV,
- Outlet valve,
- Purge valve,
- Piston.

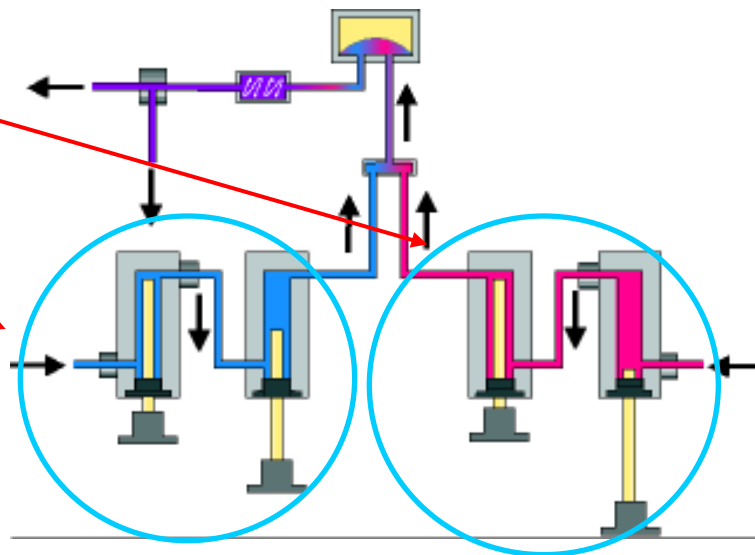
Binary Pump



Solvent Selection Valve (Option)



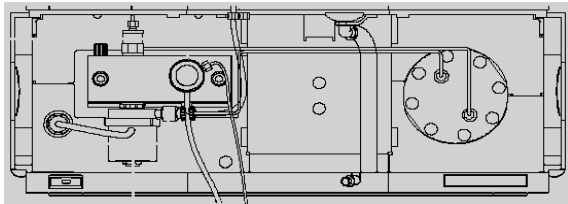
Pumpheads



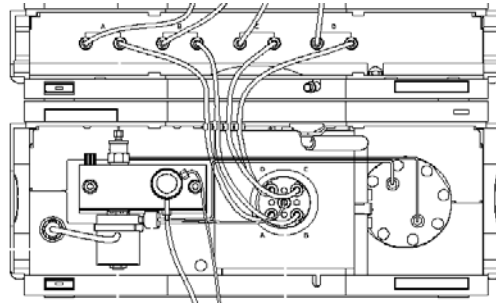
Maintenance areas:

- Seals,
- AIV,
- Outlet valve,
- Purge valve,
- Piston.

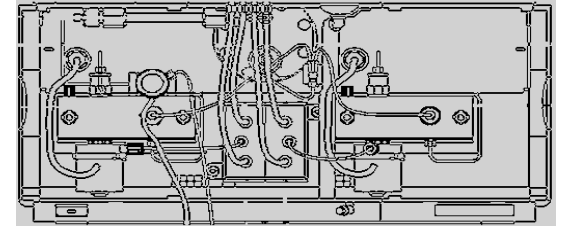
Pump Models for 1100 & 1200



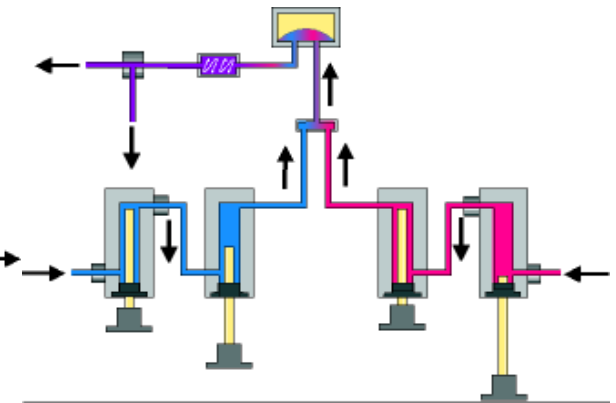
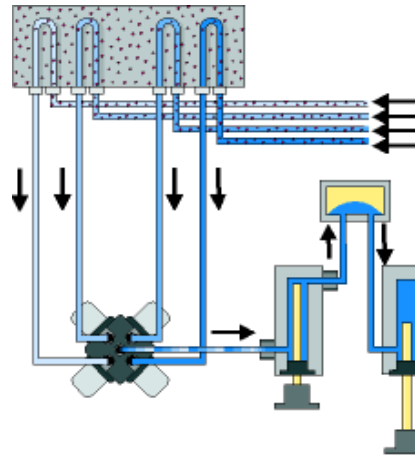
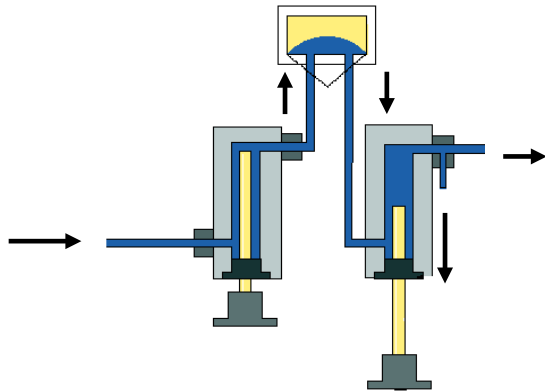
Isocratic Pump



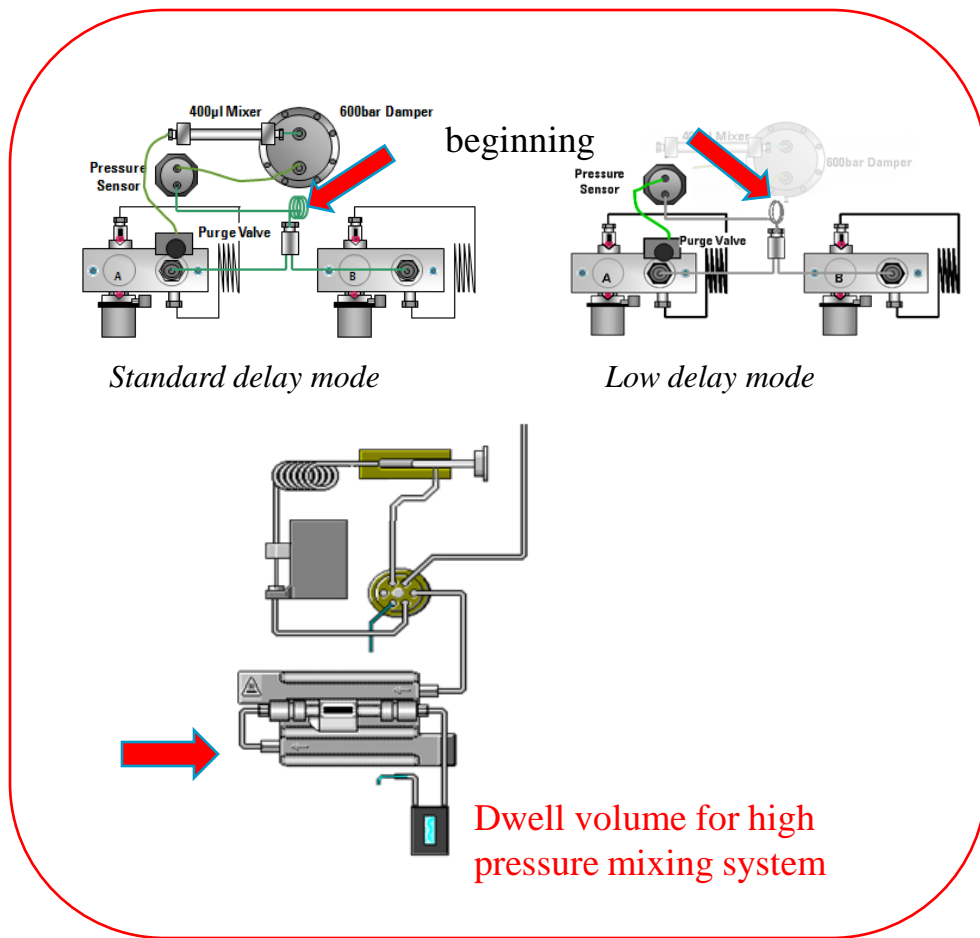
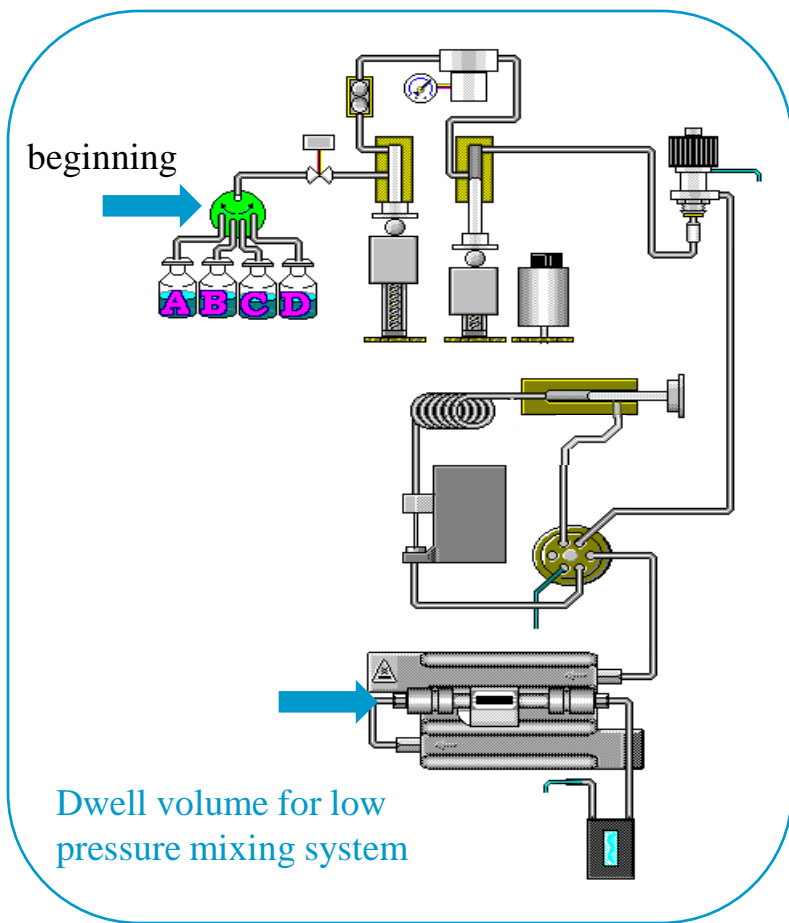
Quaternary Pump



Binary Pump



UHPLC Volume Effects



- Dwell volume = volume from formation of gradient to the column
- Behaves as isocratic hold at the beginning of gradient.

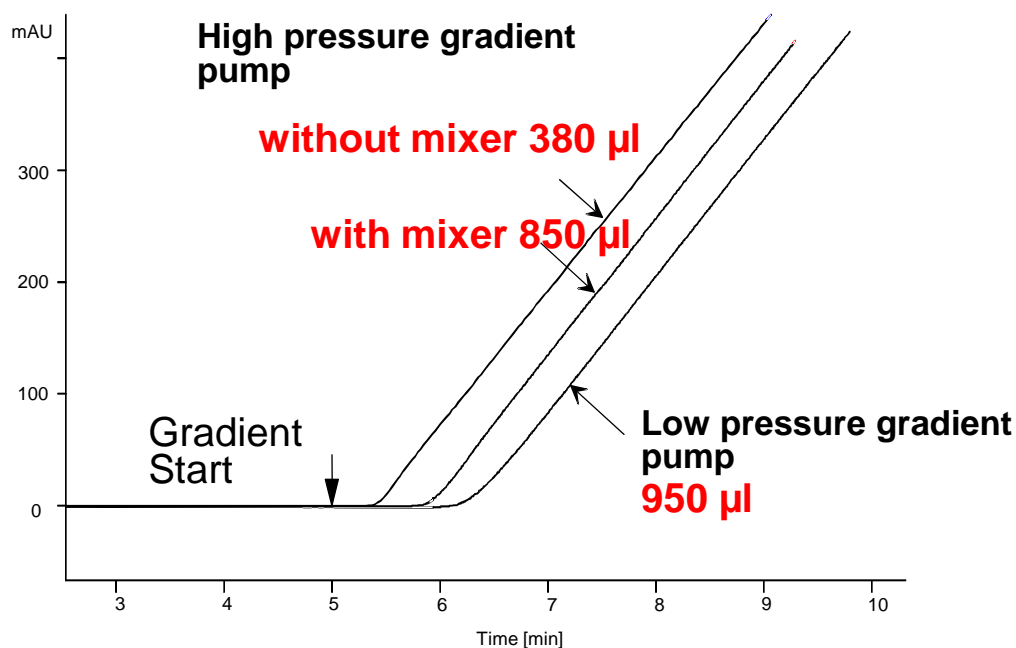
Comparison of System Delay Volumes

		1090	1050	1200 Quat.	1200 Bin.
Pump	w/o mixer	300-500	800-1200	800-1200	180-480
	w/ mixer	1050-1250	n/a	n/a	600-900
Mixer		750	n/a	n/a	420
		V (loop)	327 + V (inj)	300 + V (inj)	300 + V (inj)
Autosampler Standard		N/A	8	6.2	6.2
	Bypass	4.1 or 8.2 0	15 ul 0	3 or 6 0	3 or 6 0
Column compartment standard	Bypass				
min Range		304-504	1242-1442	1203-1406	189-489
Max Range		1058-1258			906-1206

System Delay Volume with Low Pressure and High Pressure Gradient Pump

Agilent 1100 / 1200 Pumps

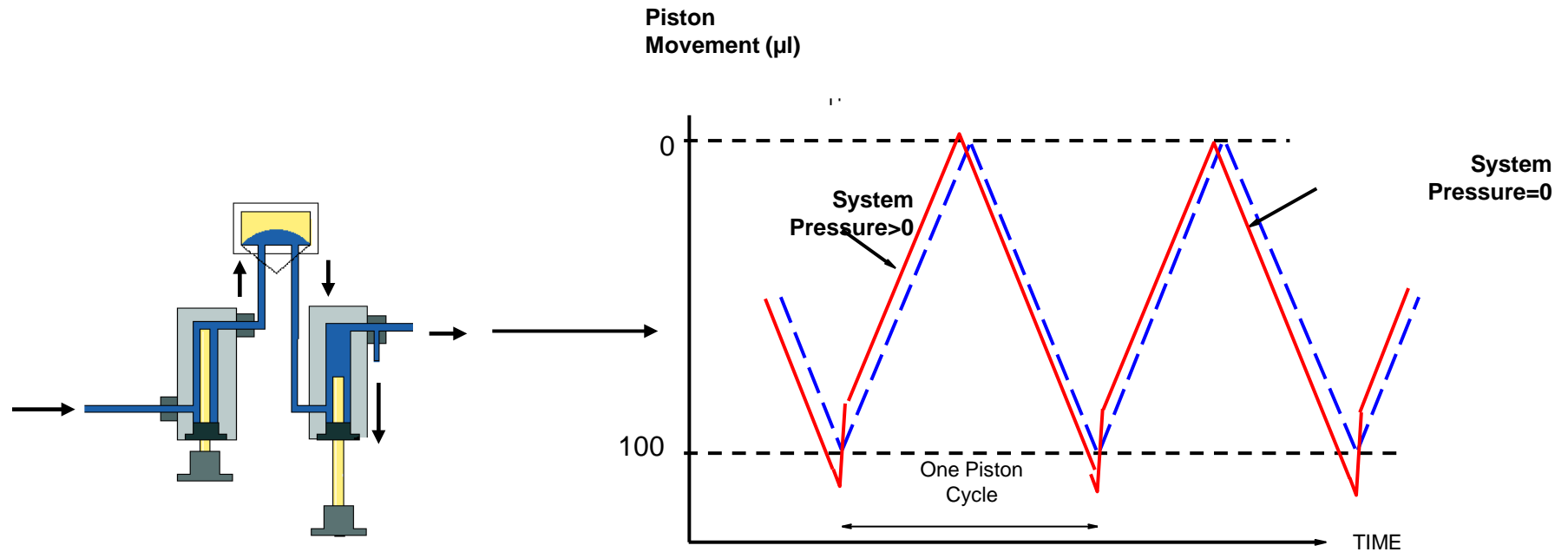
Performance



Solvent A: Water
Solvent B: Water+0.5%
Acetone

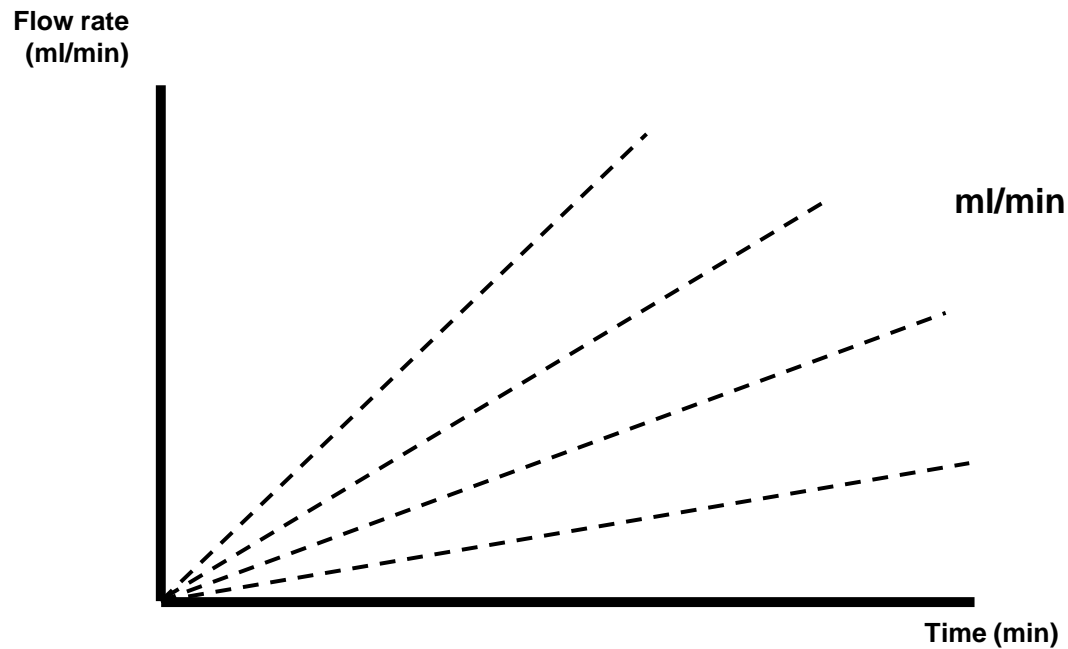
Flow rate: 1.0
ml/min
Pressure: 130 bar

Solvent Compressibility Corrections

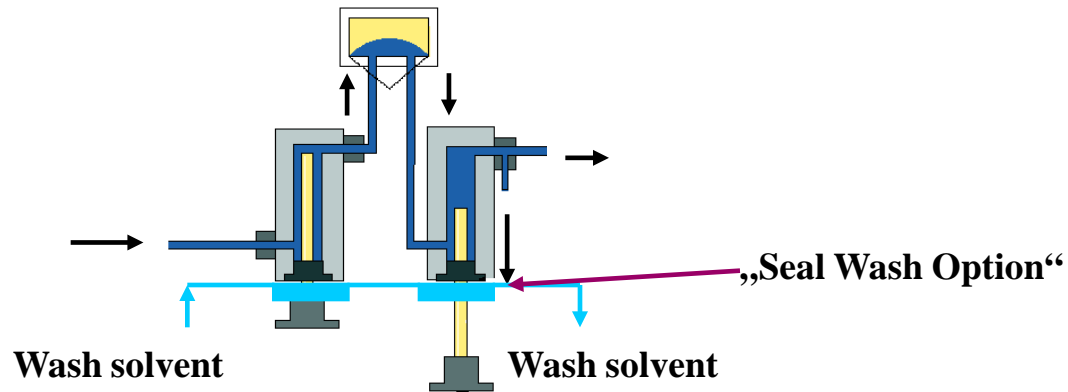


Column Protection: Soft Start Function

Flow Rate Ramp

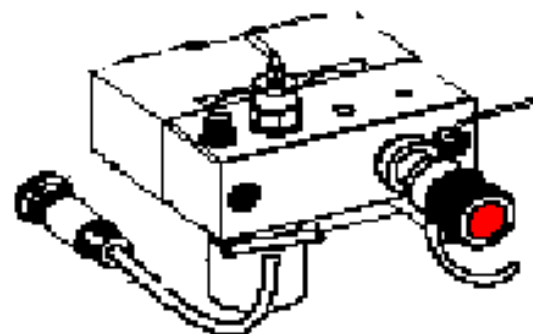
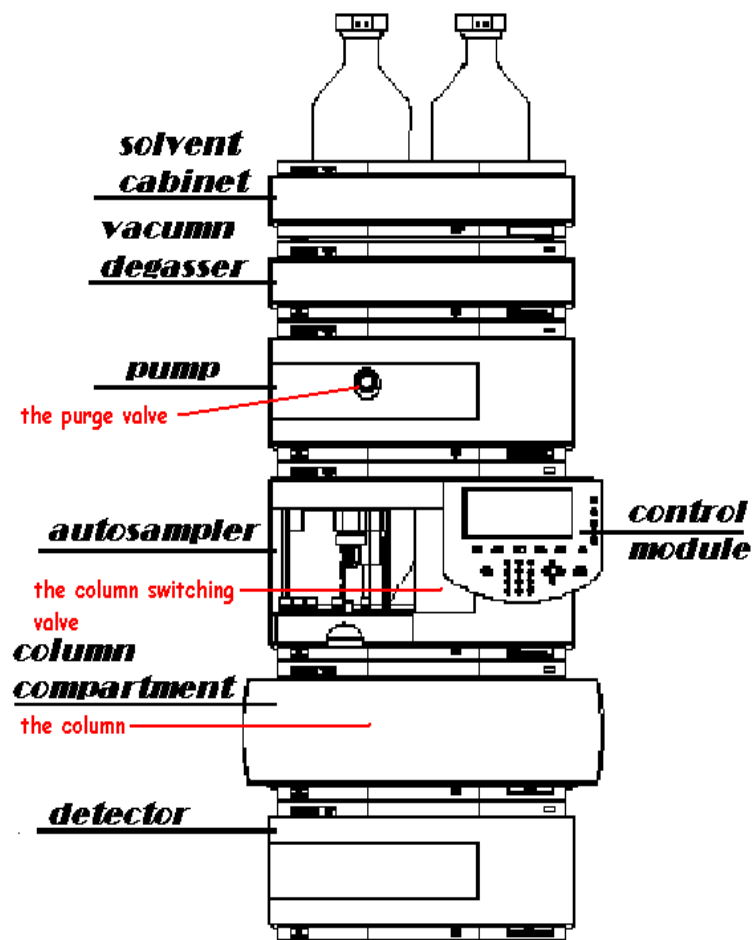


Seal Wash Option

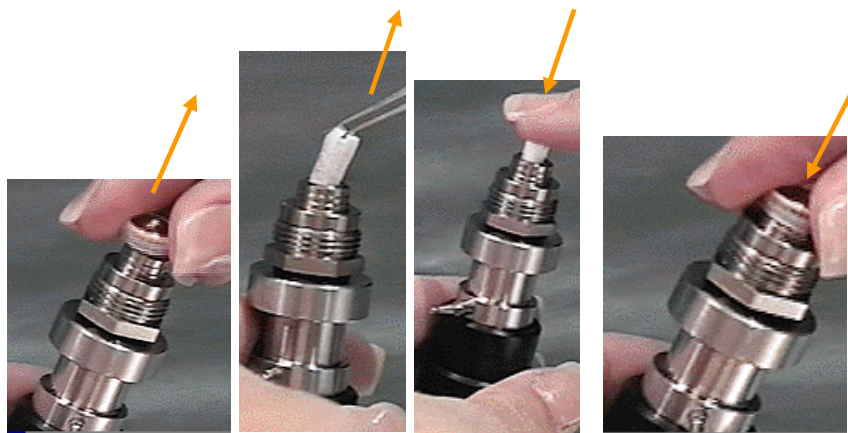


The *Seal Wash Option* can be installed to protect your pistons and seals from the routine use of highly concentrated aqueous buffer solutions (>0.01M)

The 3 point break

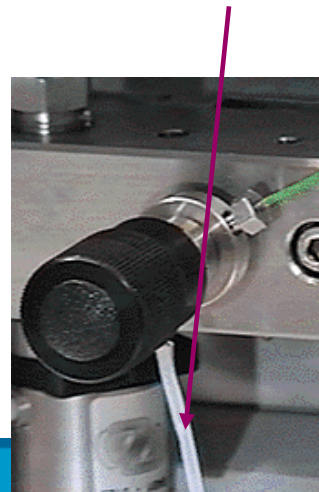
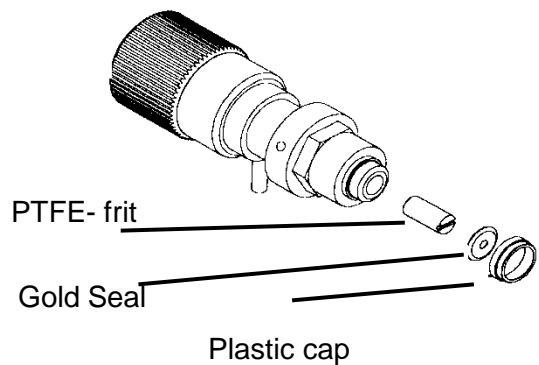


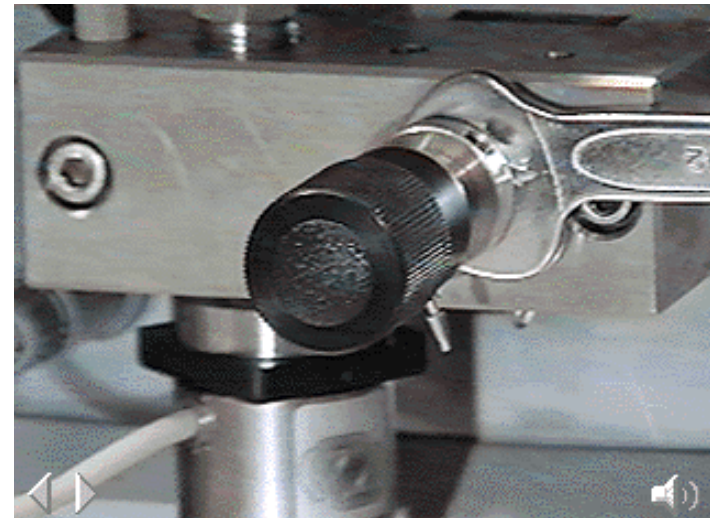
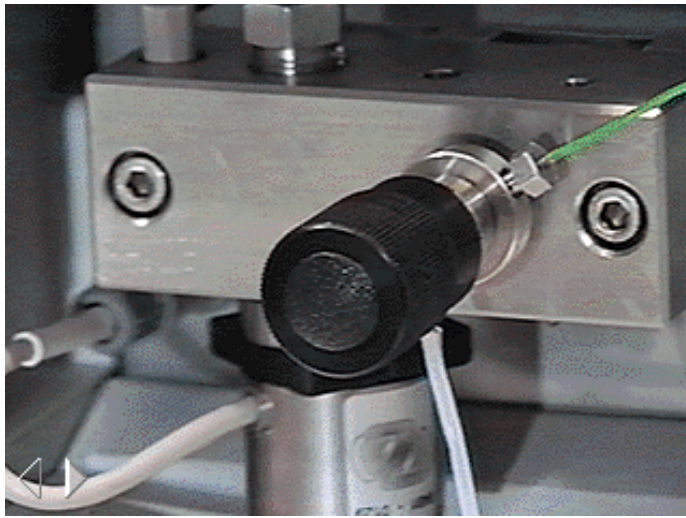
Purge Valve – Exchanging the PTFE Frit



1. Unscrew the valve using a 14 mm wrench
2. Remove the plastic cap and the gold seal
3. Take out the frit (tweezers)
4. Install a new frit
5. Replace the gold seal and the plastic cap
6. Install the valve

Note: Realign the waste tube in the correct orientation during installation.





The Outlet Ball Valve

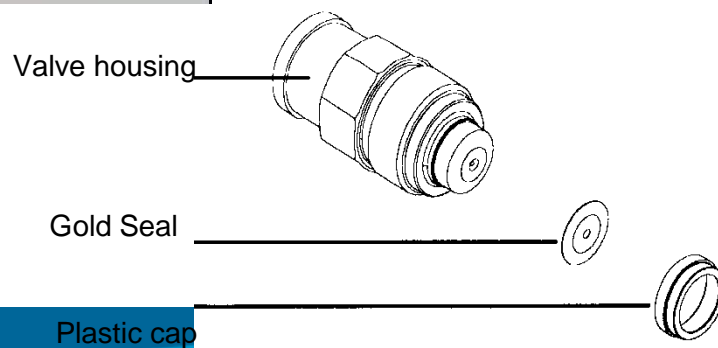


Outlet Valve – Cleaning or Exchanging the Outlet Ball Valve



1. Remove the capillary from the valve
2. Remove the valve using a wrench
3. Clean the ball valve in the ultrasonic bath or replace the ball valve
 - a. Take off the plastic cap and gold seal
 - b. Replace the ball valve
 - c. Replace the gold seal and cap
4. Reinstall

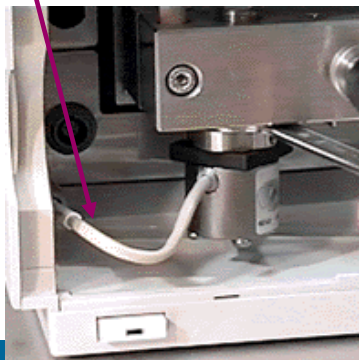
Note: The outlet ball valve of the binary pump has an additional sieve (5063-6505)



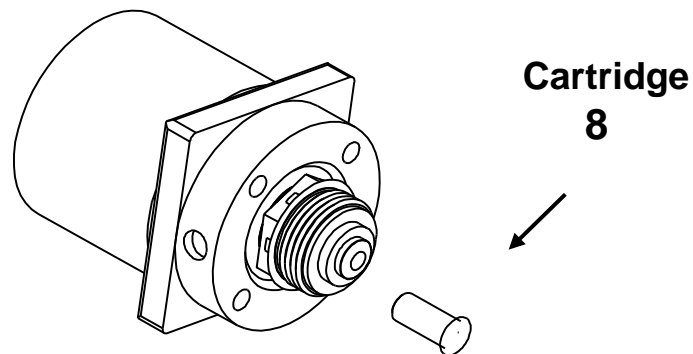
Active Inlet Valve (AIV) – Old/New

1. Remove the AIV using a 14 mm wrench
2. Change the cartridge (new design)
3. Change the gold seals when necessary
4. Reinstall the AIV

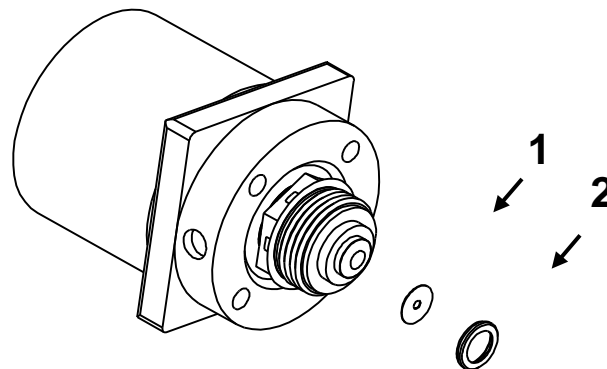
Note: Properly position the AIV cable when you reinstall the valve



New design



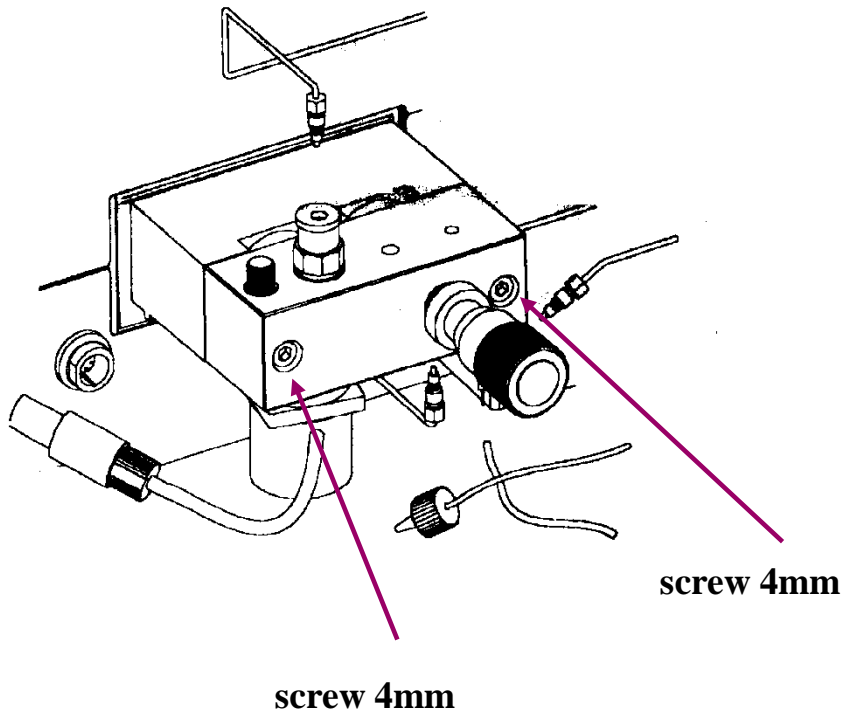
Old design



Active Inlet Valve (AIV)

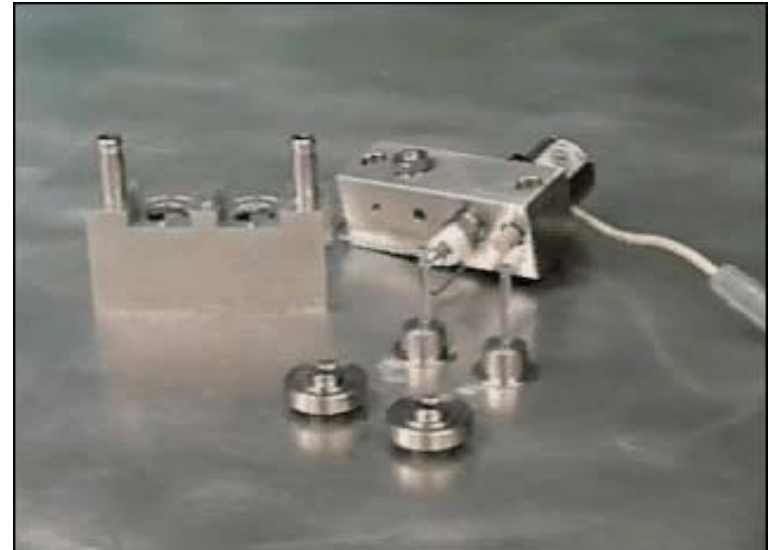
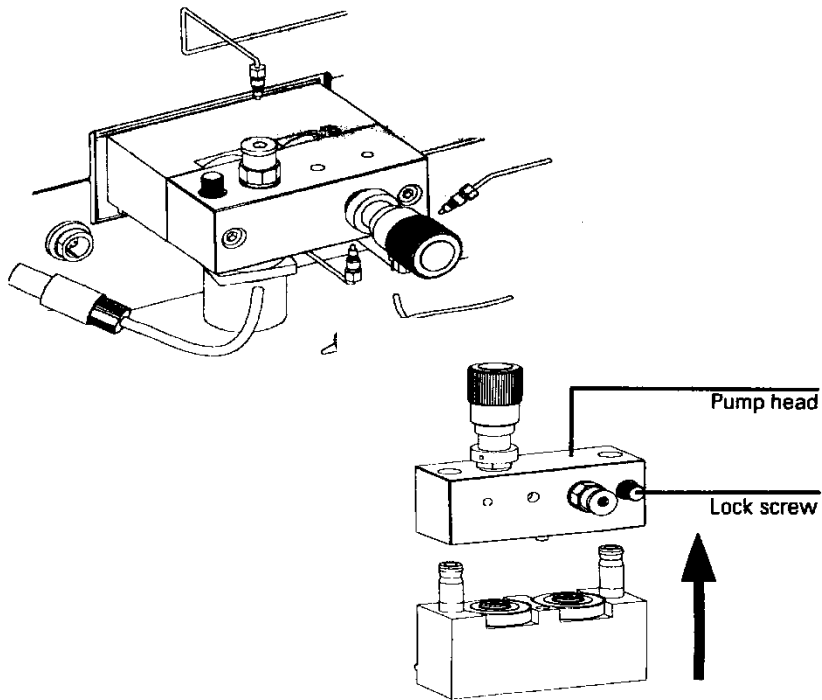


Removing the Pump Head



- Remove all capillaries
- Disconnect the AIV supply cable
- Remove both pump head hexagonal screws
- Remove the pump

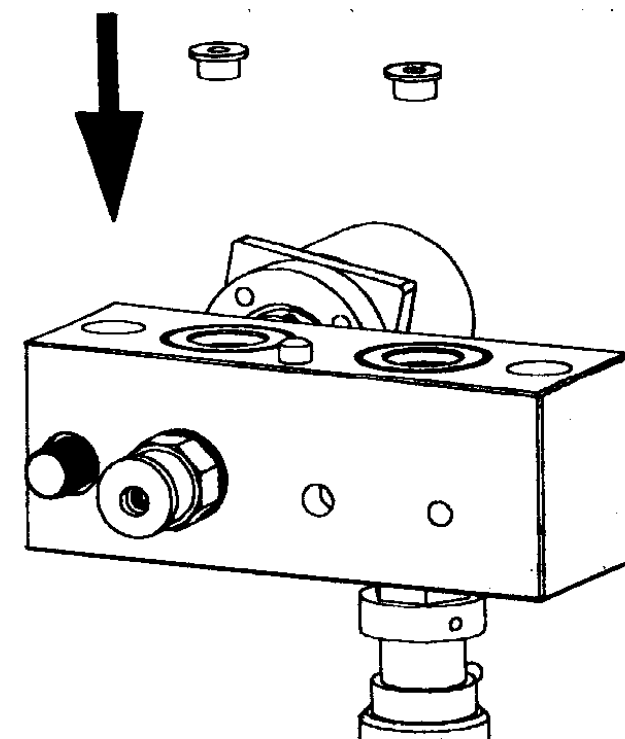
Disassembling the Pump Head



The two parts of the pump head are disconnected by releasing the lock screw.

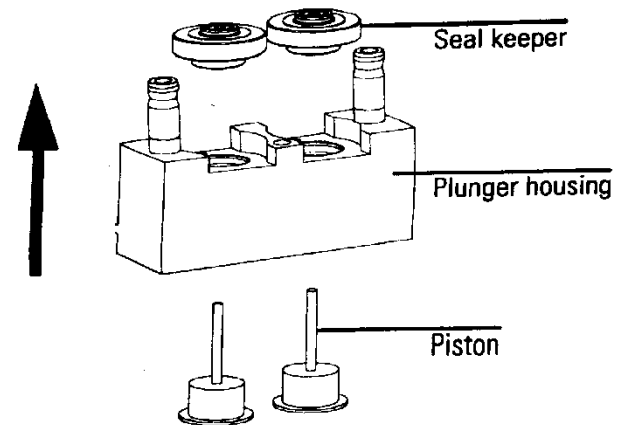
Changing the Pump Seals

1. Remove the old seals.
2. Remove the wear retainers, if present.
3. Clean the pump chambers.
4. Insert new seals.
5. Reassemble the pump head.
6. Perform seal wear-in procedure for standard seals

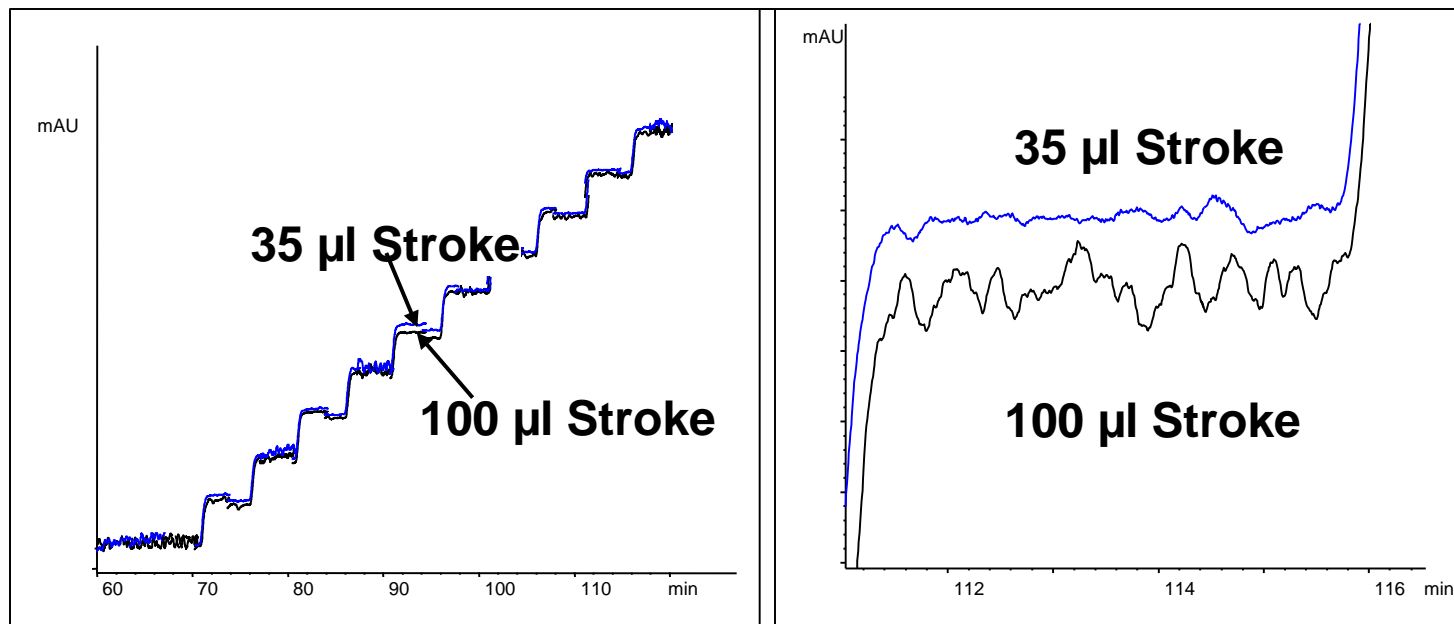


Changing and Inspecting the Sapphire Pistons

1. **Disassemble the pump head assembly**
2. **Check the plunger surface and remove any deposits or layers with alcohol or tooth paste**
3. **Replace the pistons if scratched**
4. **Reassemble the pump head**
5. **Check to make certain there are not any fractures in the springs**
6. **Put in the pistons**



Agilent 1200 Series Quaternary Pump: Variable Stroke Influence on Composition Ripple

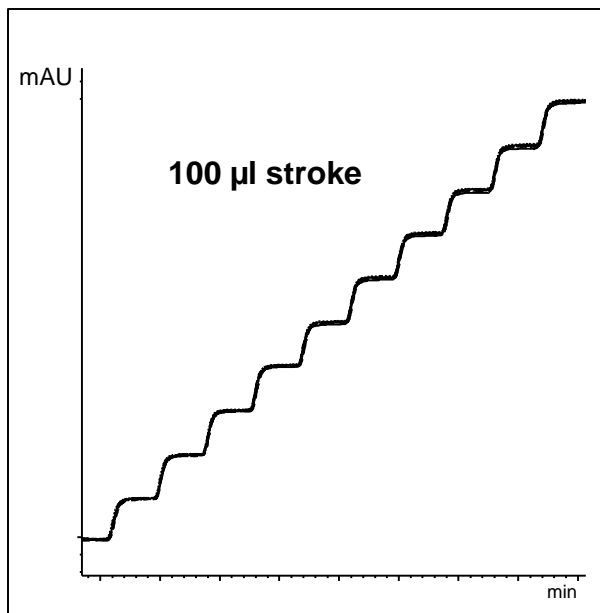


Tracer	Water/Methanol +0.5% Acetone
Flow rate	1 ml/min
Gradient	55-65 %

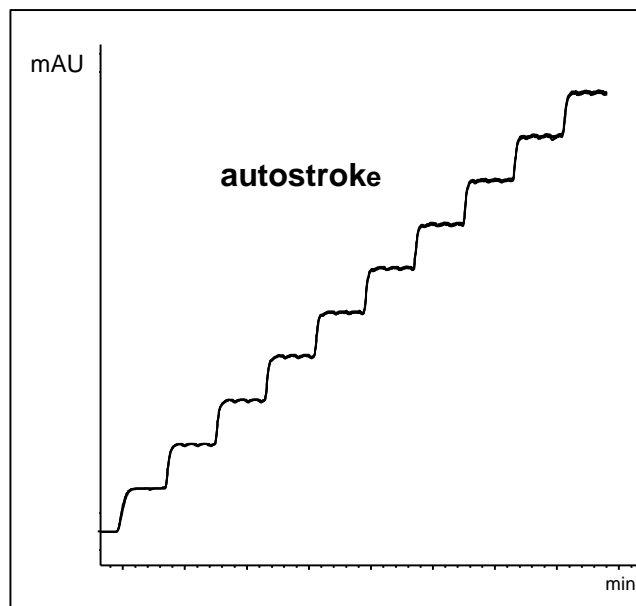
Agilent 1200 Series Pumps:

Gradient Performance 0-10%B

Quaternary Pump



Binary Pump



Tracer

**Water/Water
+0.5% Acetone**

Flow rate

1 ml/min

Gradient

0-10 %

Injector

- **Important performance characteristics**

- Injection volume precision
- Wide linearity
- Minimum carry over
- Wide dynamic injection volume

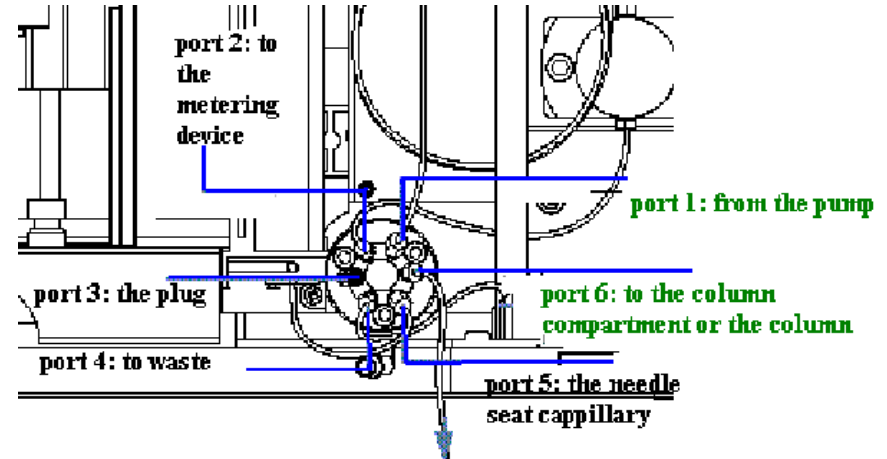
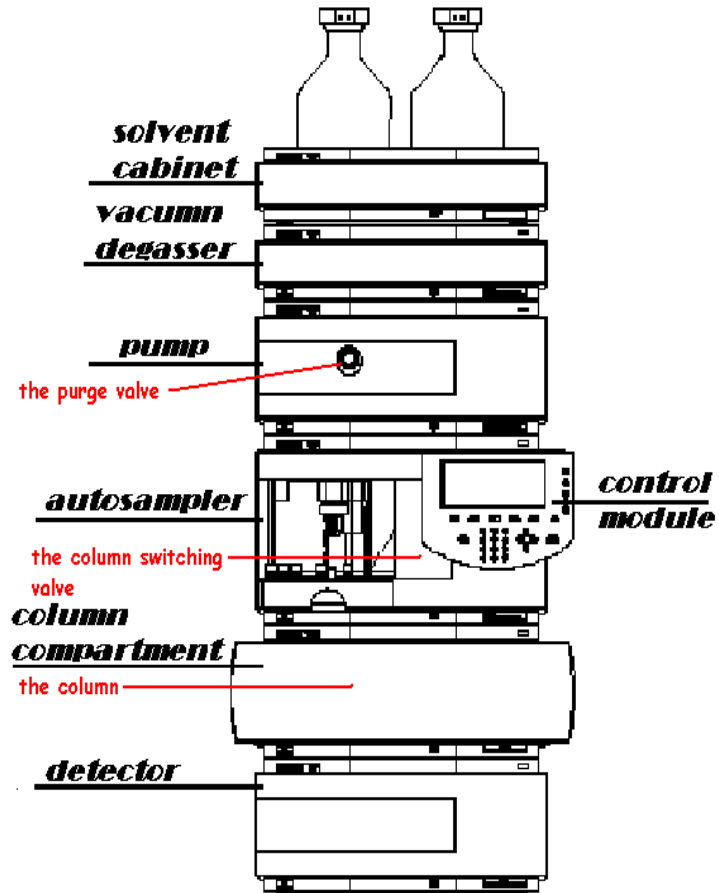


- **Influence on...**

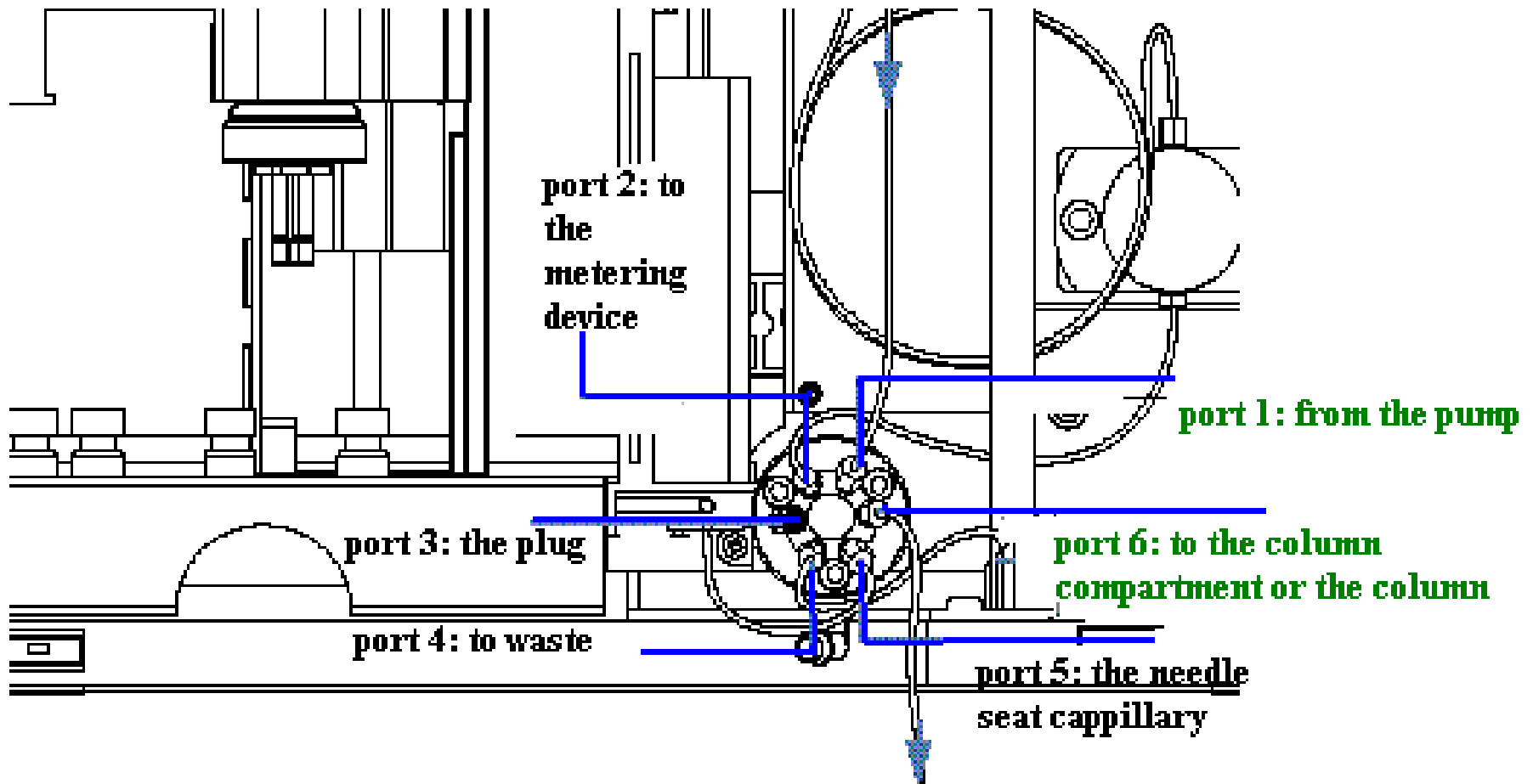
- Precision of peak area/height
- Accuracy of peak area/height (when using different injection volumes)
- Precision of peak area/height
- Versatility, application range



The 3 point break

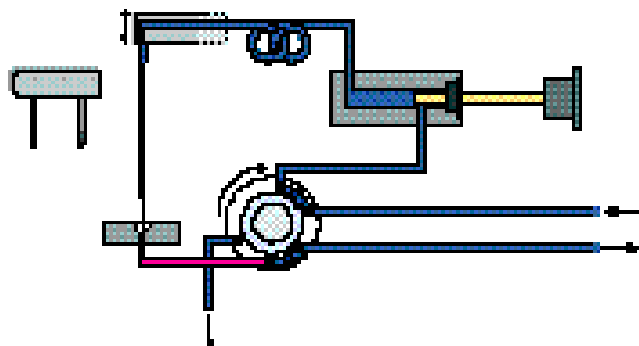


The autosampler connections

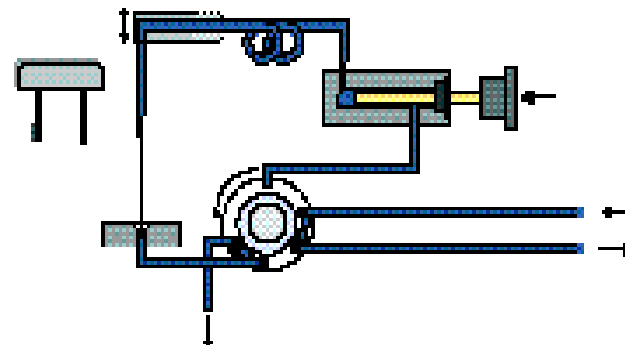


The injection sequence

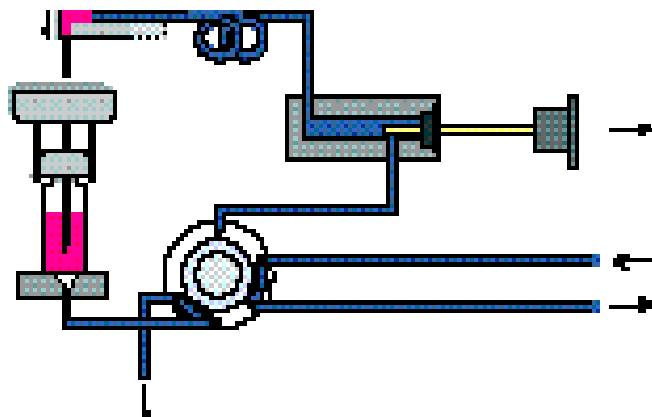
Mainpass Position



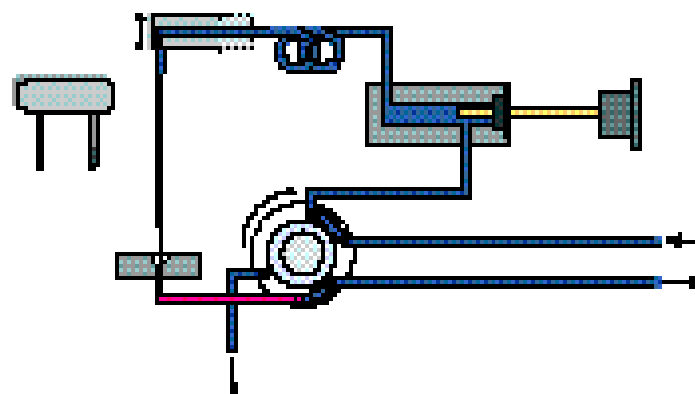
Bypass Position

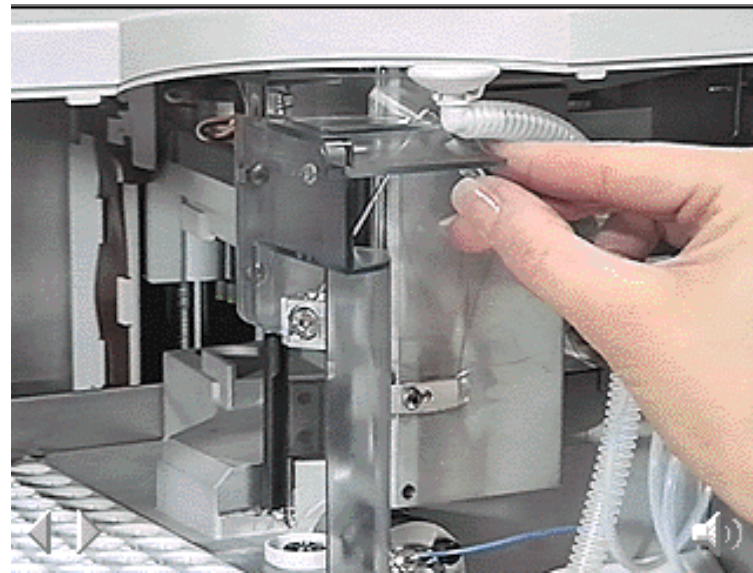
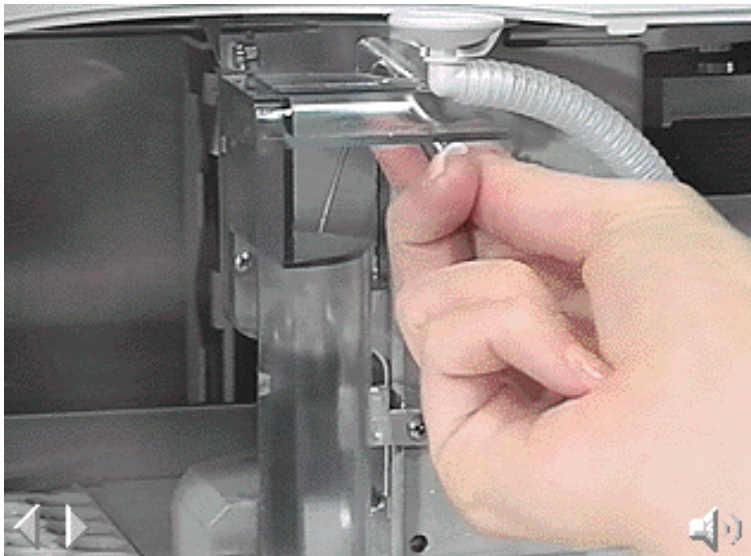
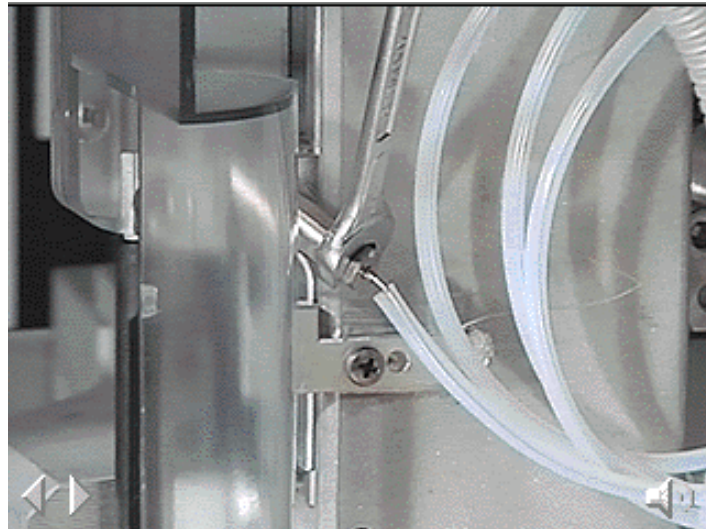
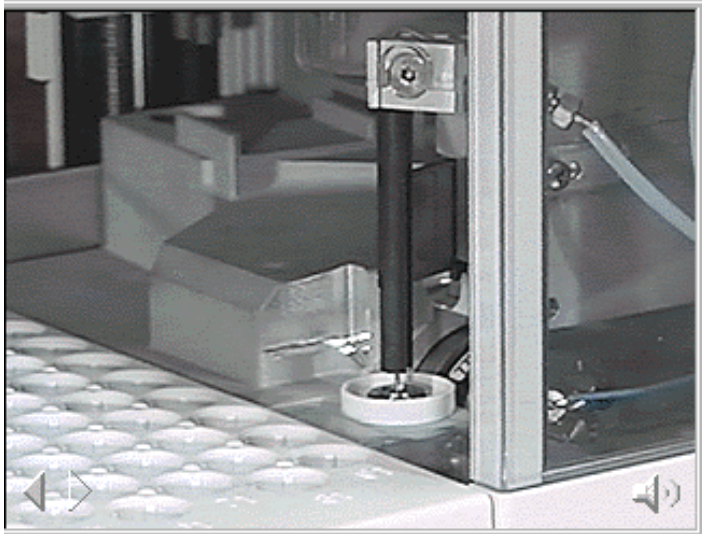


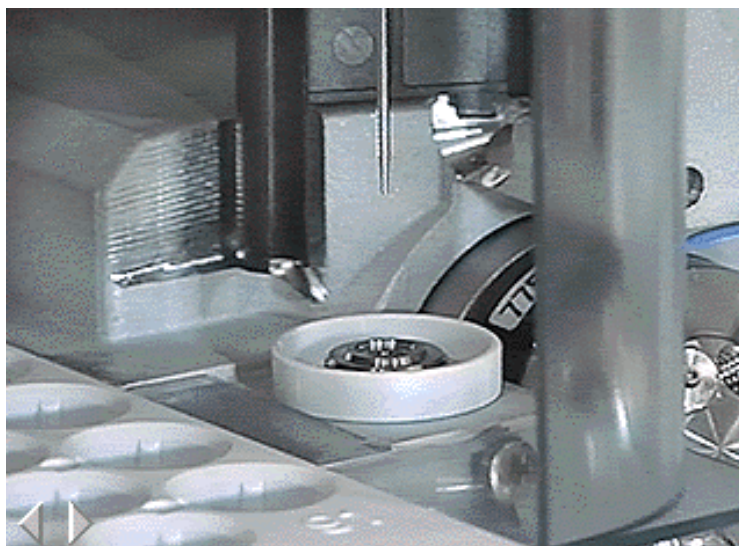
Drawing the Sample



Mainpass Position (Sample Injection)







Thermostatted Column Compartment

- **Important performance characteristics**

- Excellent temperature accuracy



- Excellent temperature precision



- **Influence on...**

- Elution order

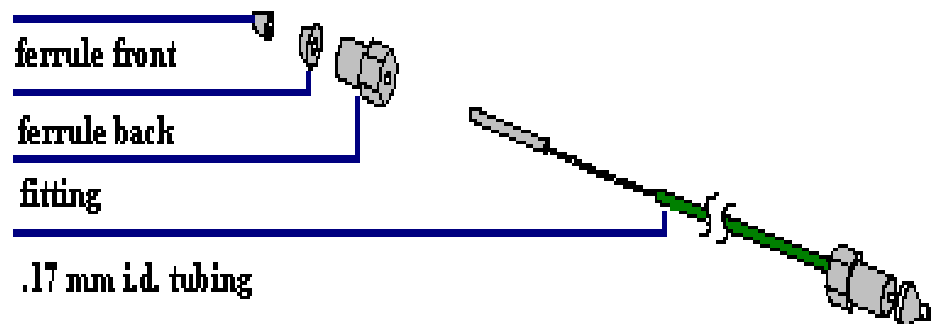
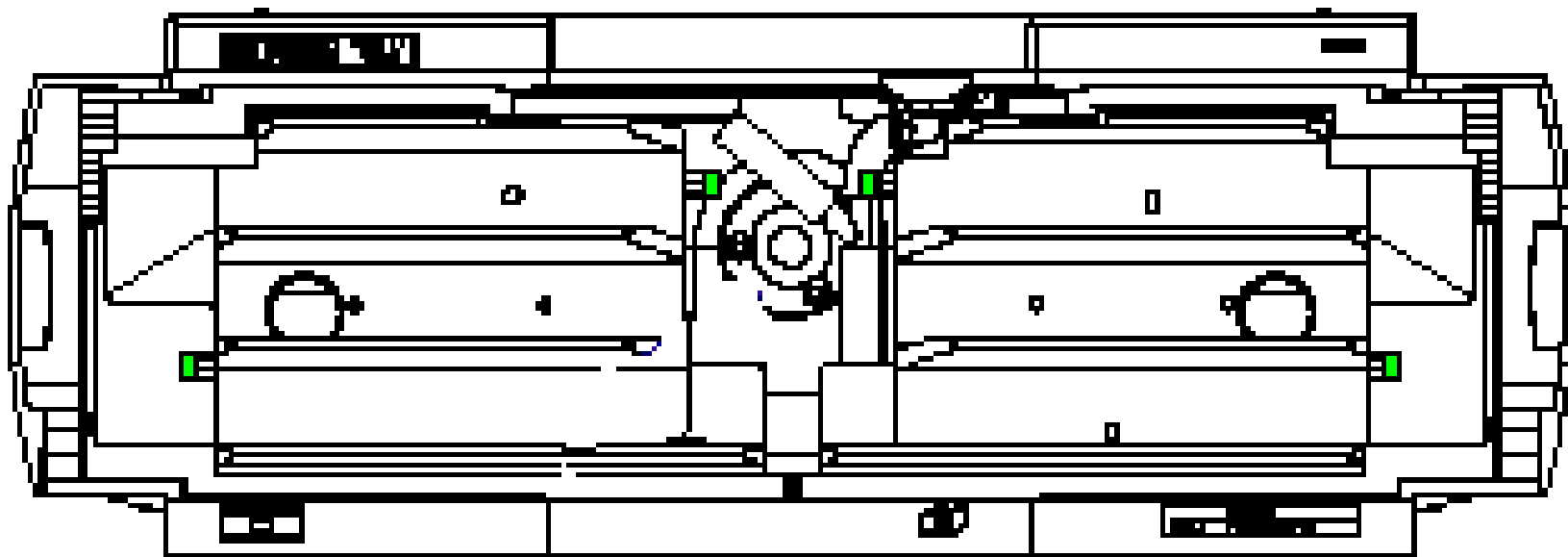
- Peak identification

- Elution order

- Retention time precision

- Peak identification





The column compartment is an optional module the output of the autosampler (port 6) is attached to the right side, the left side or the column switching valve.

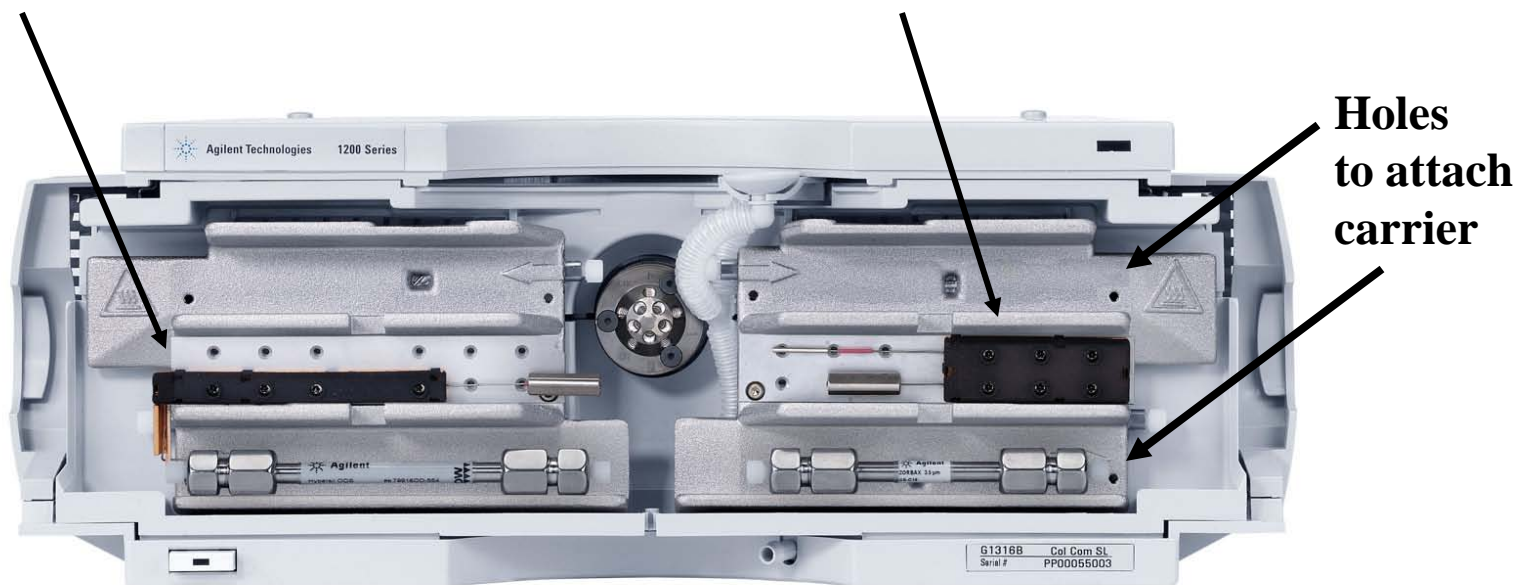


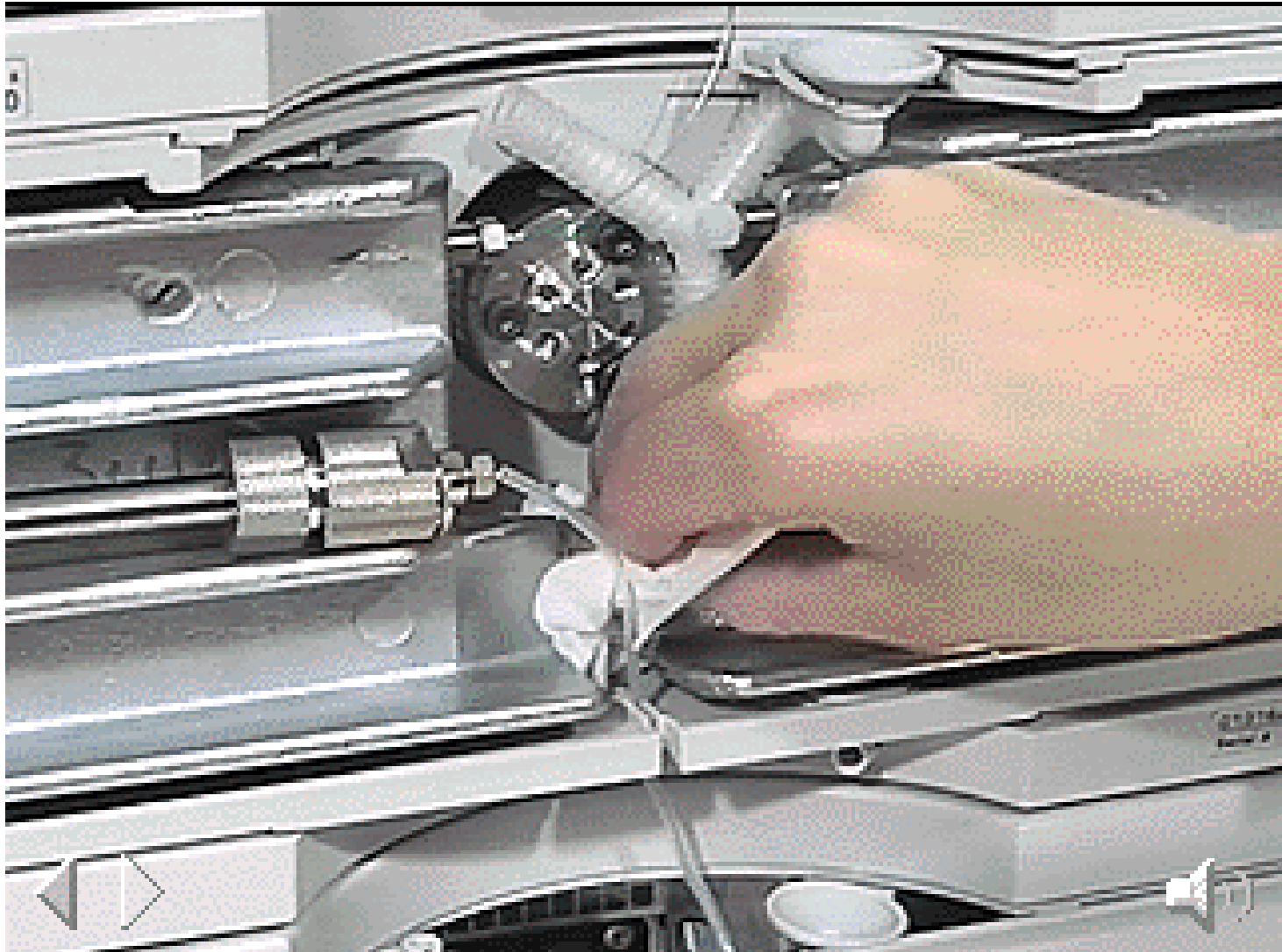
1200 Series TCC SL

Pre-column Heater and Post-column Cooler

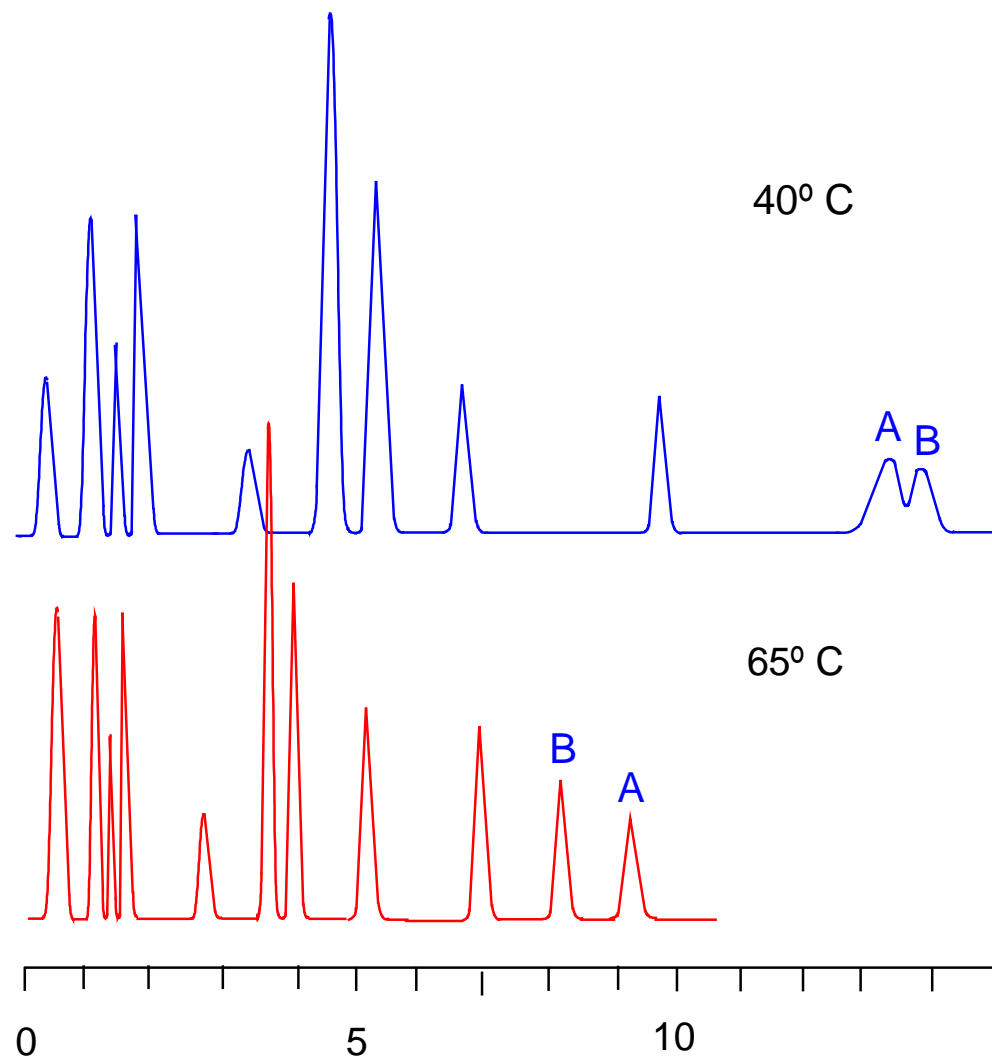
- L-shaped pre-column heater
- Volume: 1.6 μ l
- Mounted on carrier

- U-shaped Post-column cooler
- Volume: 1.5 ul
- Mounted on Carrier





Effect of Temperature on Separation



Time in
Minutes

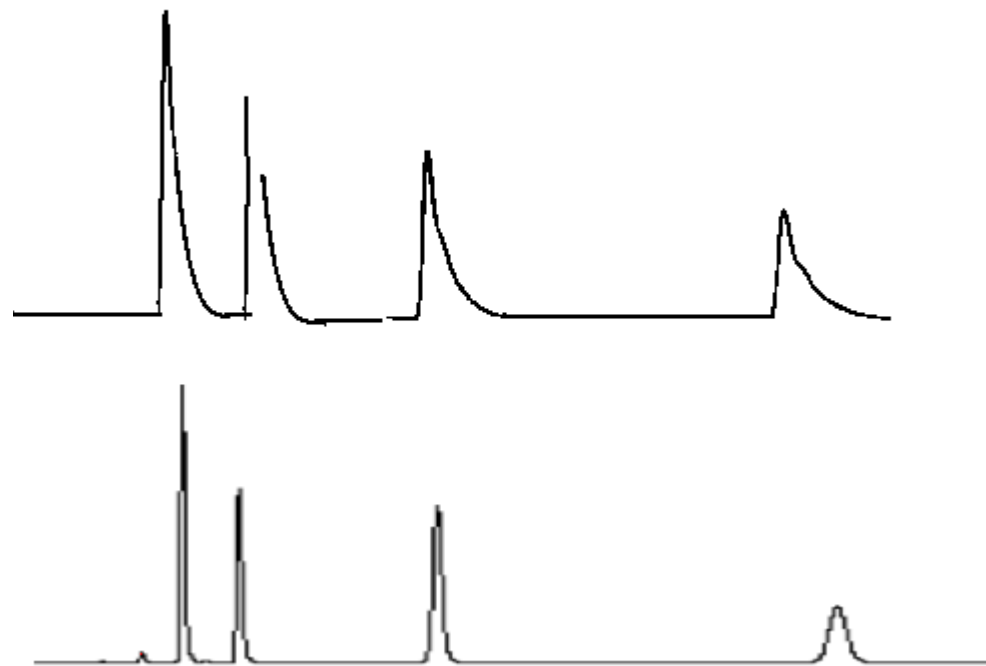


Agilent Technologies

Split Peaks

Can be caused by:

- Column contamination
- Partially plugged frit
- Column void
- Injection solvent effects



Peak Tailing, Broadening and Loss of Efficiency

May be caused by:

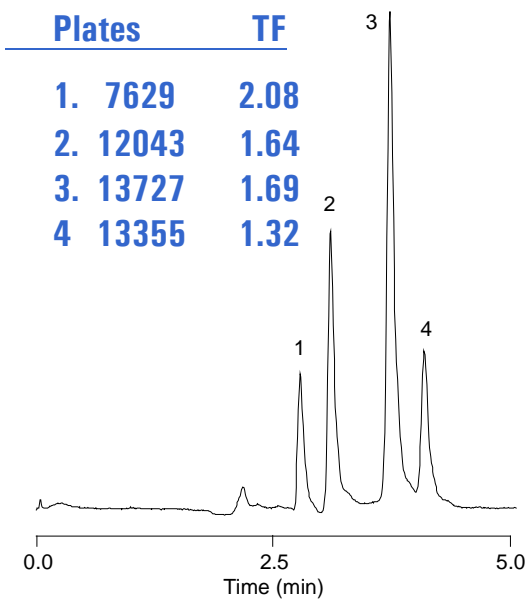
- Column “secondary interactions”
- Column void
- Column contamination
- Column aging
- Column loading
- Extra-column effects

Peak Tailing

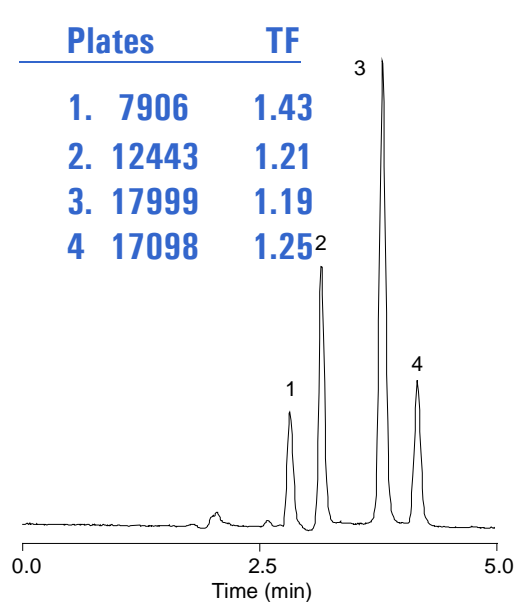
Column Contamination

Column: StableBond SB-C8, 4.6 x 250 mm, 5 μ m Mobile Phase: 20% H₂O : 80% MeOH Flow Rate: 1.0 mL/min
 Temperature: R.T. Detection: UV 254 nm Sample: 1. Uracil 2. Phenol 3. 4-Chloronitrobenzene 4. Toluene

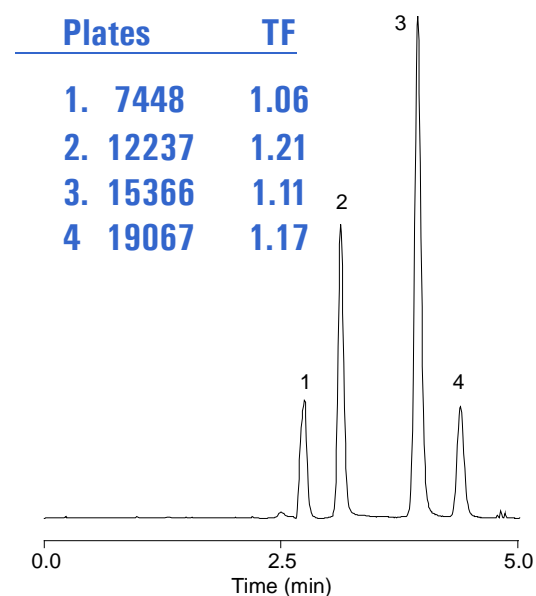
QC test forward direction



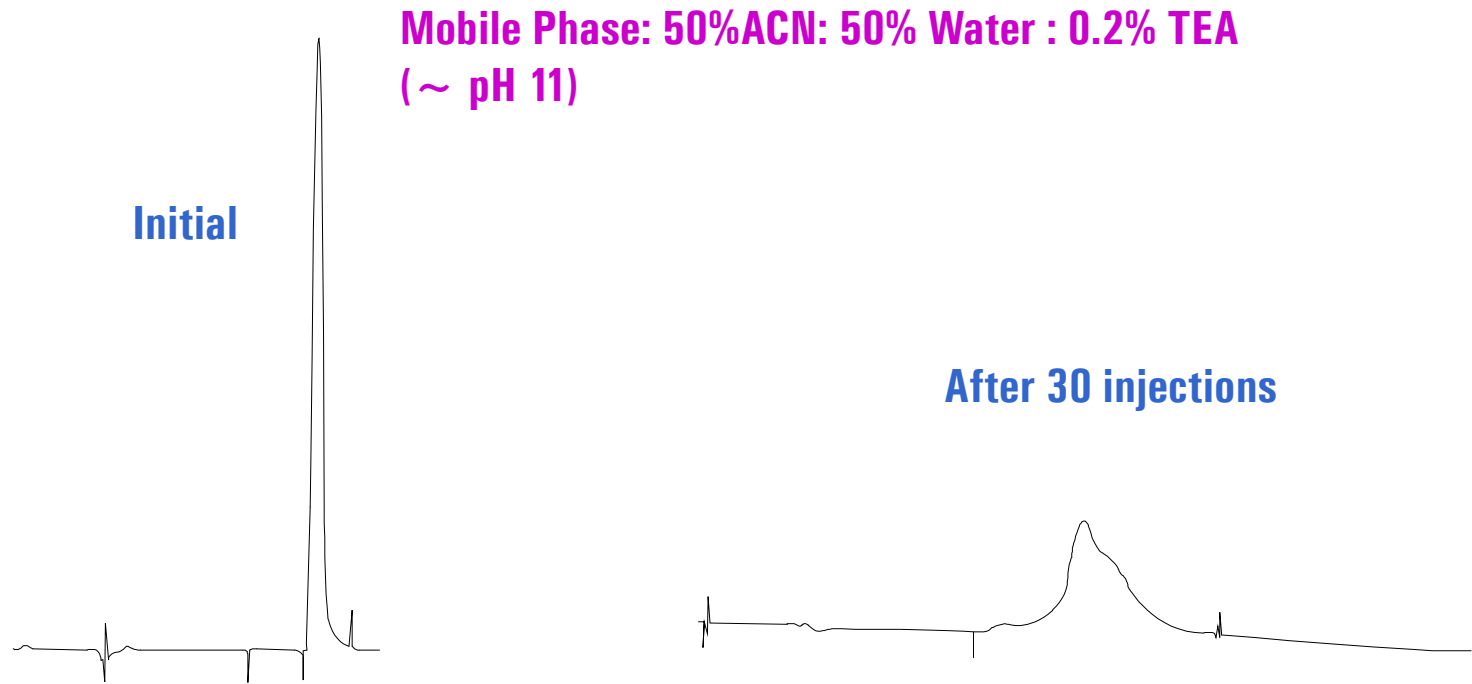
QC test reverse direction



QC test after cleaning 100% IPA, 35°C



Peak Broadening & Splitting



- Multiple peak shape changes can be caused by the same column problem. In this case a void resulted from silica dissolved at high pH.

HPLC UV/Vis Detectors

- **Important performance characteristics**

- Variable Wavelength and Diode Array Detector

- Low noise, wander and drift
- Wide linear range

- Very good wavel. accuracy
- Excellent wavel. precision

- Diode Array Detector only

- High spectral resolution

- Excellent spectral sensitivity

- **Influence on...**

- Variable Wavelength and Diode Array Detector

- Detection limit, quantitation limit
- Confidence in quantitation at high and low concentrations
- Accuracy of peak areas/heights
- Precisions of peak areas/heights

- Diode Array Detector only

- Accuracy of spectra, peak identification by spectra
- Accuracy of spectra, peak identification by spectra at low concentrations

Flow Cells with longer path lengths yield higher signals

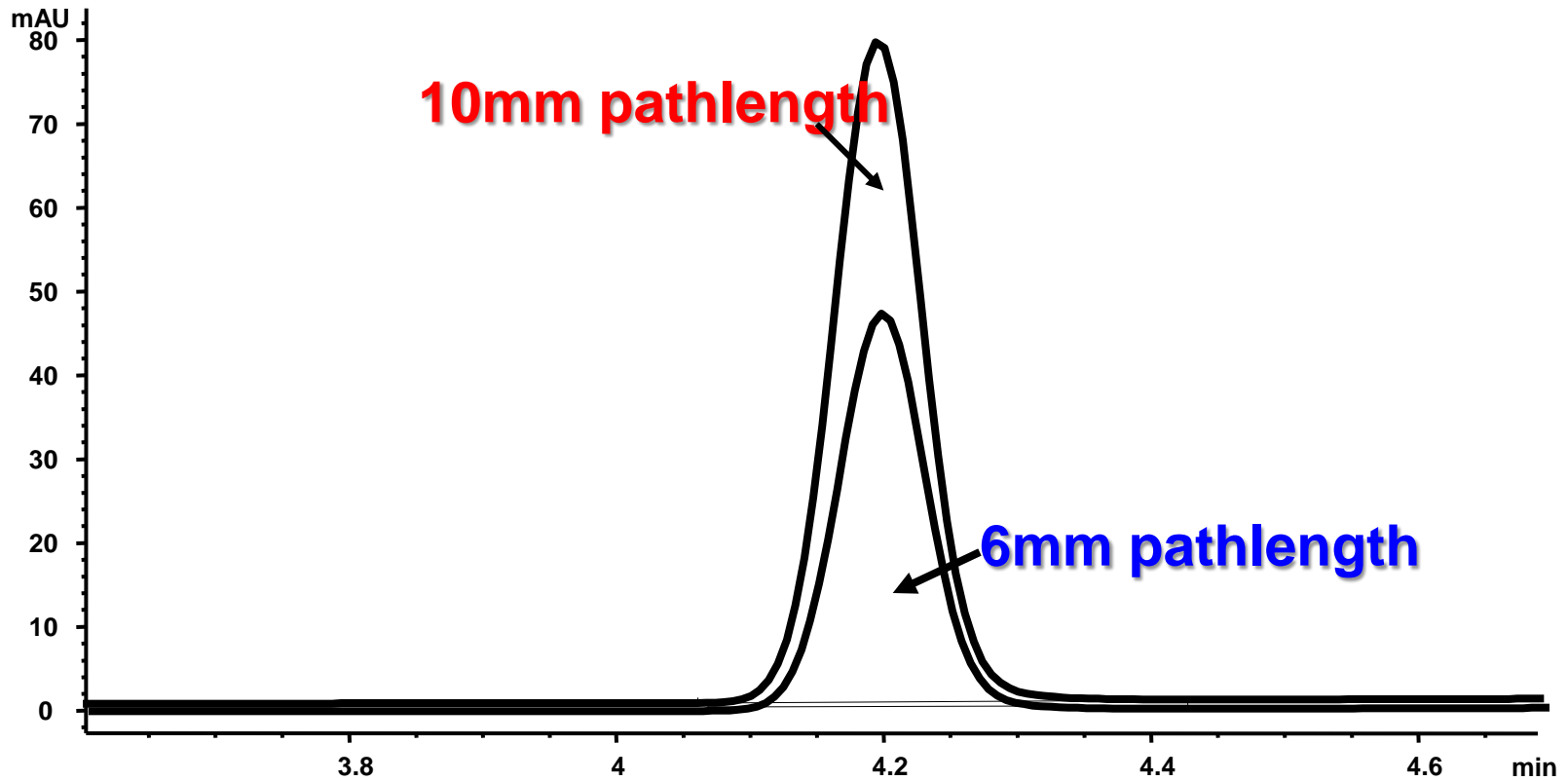
VWD STANDARD CELL

Standard: 14- μ l volume, 10-mm cell path length and 40 bar

DAD STANDARD CELL

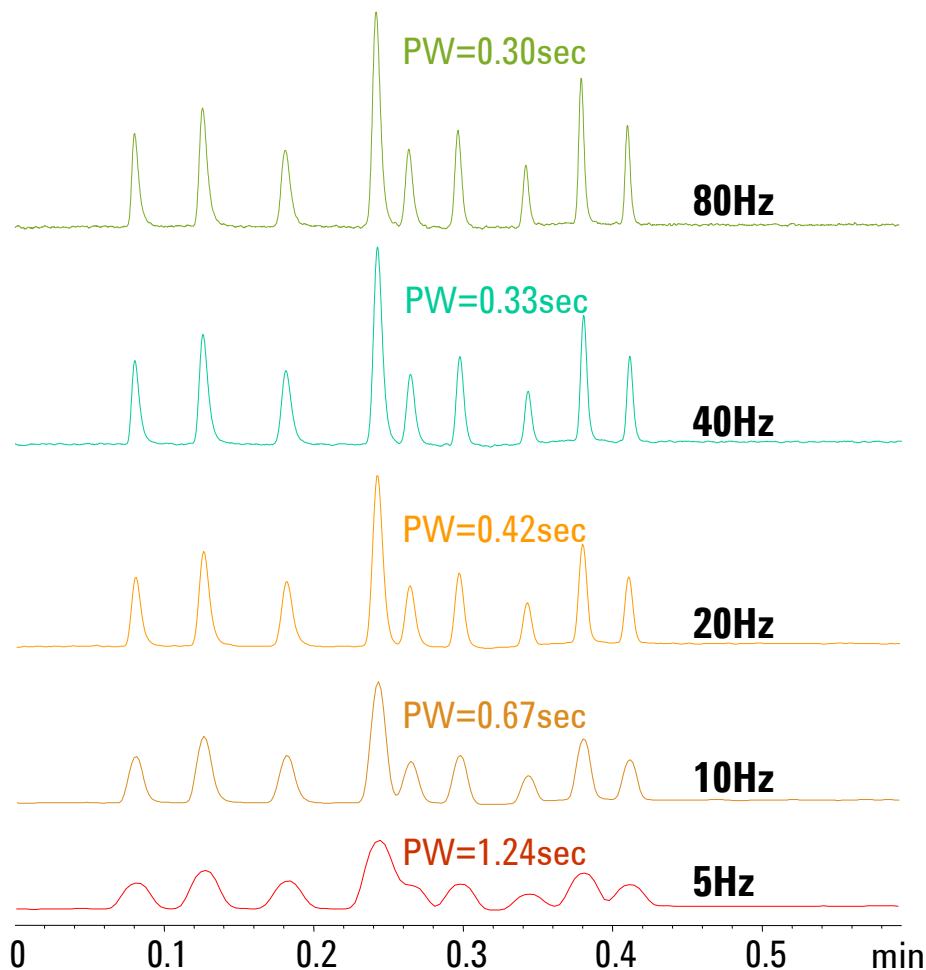
Standard: 13 μ l volume, 10 mm cell path length and 120 bar

Influence of Pathlength on Signal Sensitivity



Benefit of 80Hz Data Acquisition Rate

Peak Width, Resolution and Peak Capacity in Ultra-Fast LC



20Hz versus 80Hz

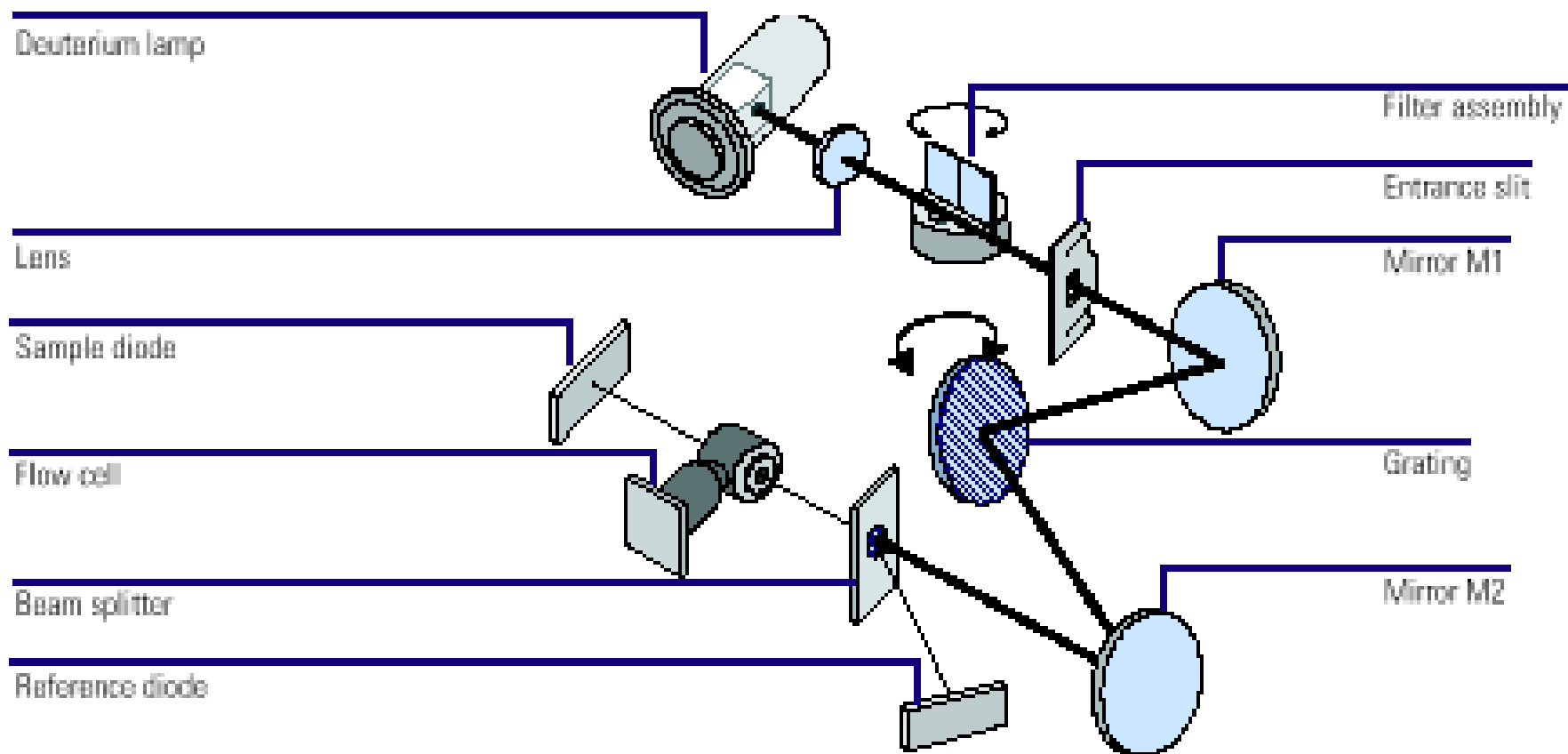
- + 40% Peak Width
- 40% Peak Capacity
- 30% Resolution
- 70% Apparent Column Efficiency

10Hz versus 80Hz

- + 120% Peak Width
- 120% Peak Capacity
- 90% Resolution
- 260% Apparent Column Efficiency

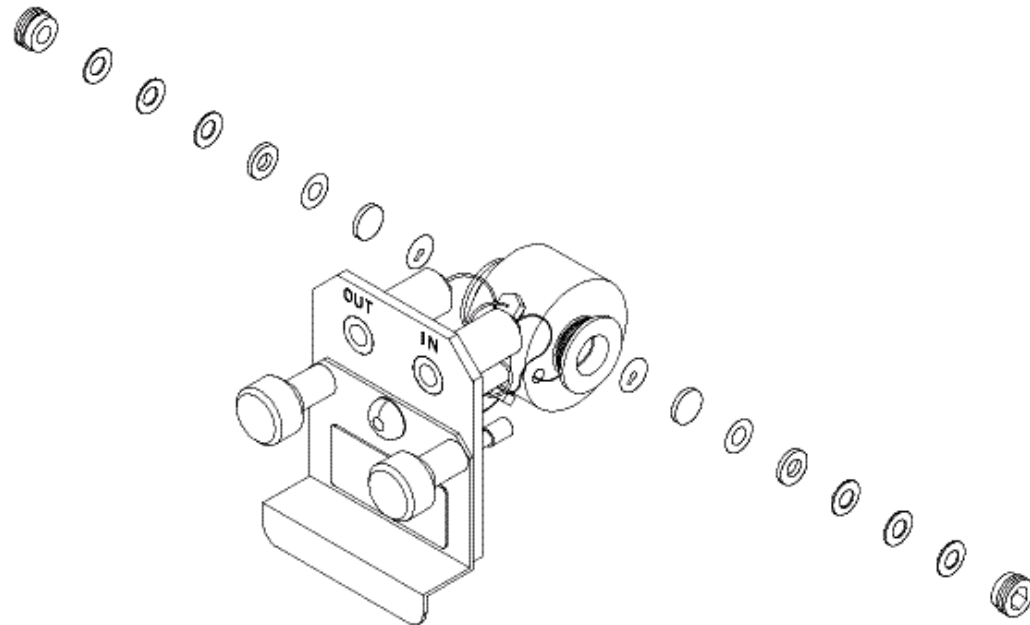
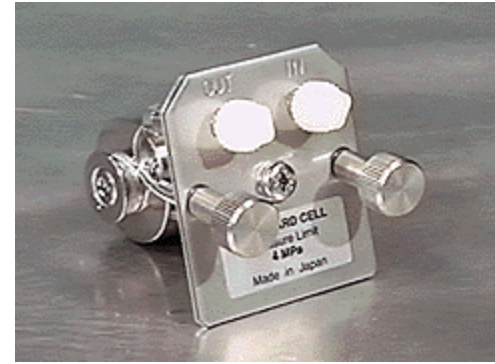
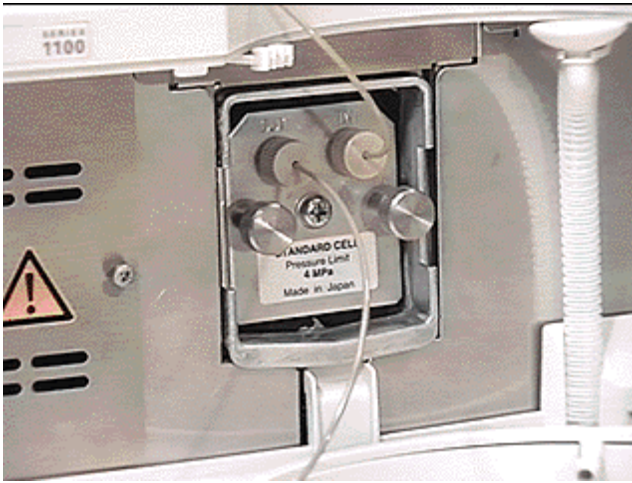
- Sample: Phenone Test Mix
- Column: Zorbax SB-C18, 4.6x30, 1.8 μ m
- Gradient: 50-100% ACN in 0.3min
- Flow cell: 5 μ l

Figure 6 **Optical Path of the Variable Wavelength Detector**

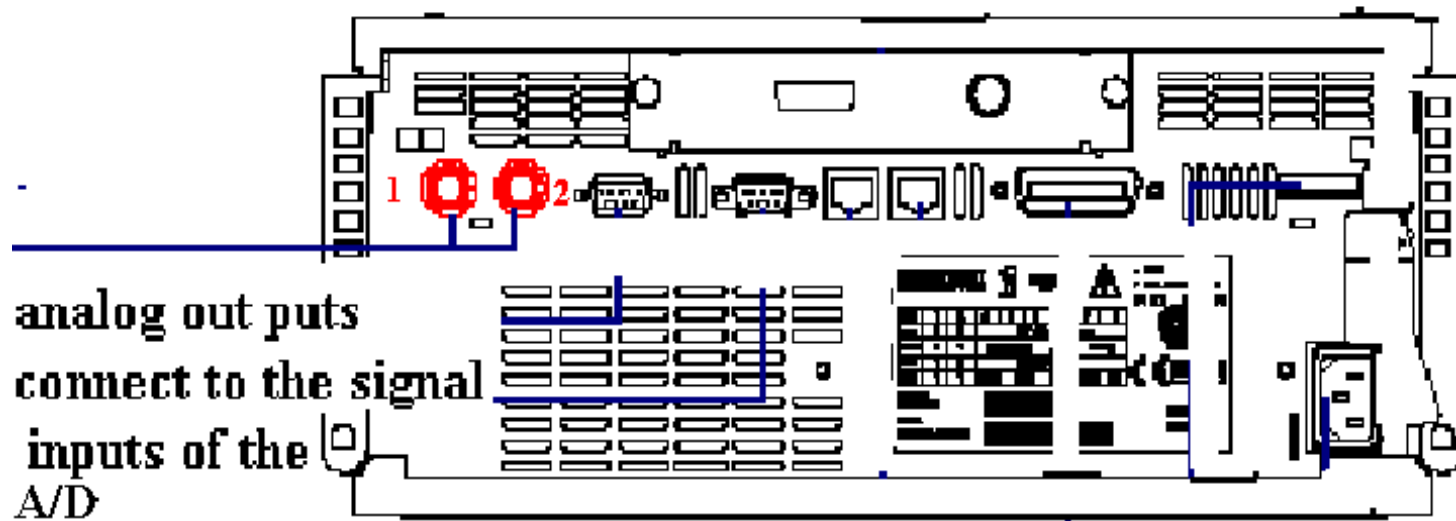


Changing the VWD Lamp





Connections to an A/D box



Analog connections

General purpose (spade lugs)

01046-60105

Please note the analog cable and connections are the same for all 1200 detector outputs to a non HP data system

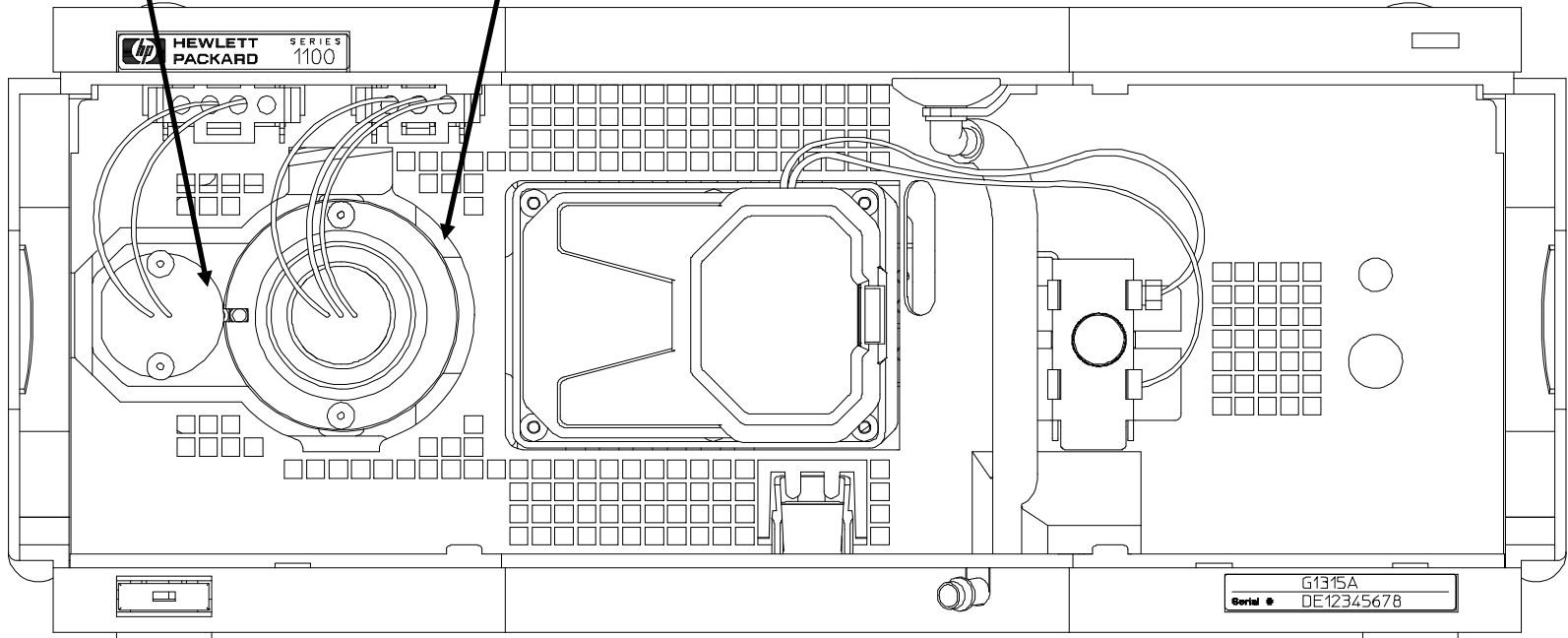
HP 1200 Diode Array

- **Wavelength Range:** 190 - 950 nm
- **Lamps:** Shine-through deuterium lamp (uv-range)
Tungsten lamp (vis-range)
- **Slits:** Programmable electromechanical; 1, 2, 4, 8 and 16 nm
- **Noise:** 2×10^{-5} AU at 254 nm and 750 nm
- **Flow Cells:** STD 10 mm 13 ul, Semi-micro 6 mm 5 ul, Micro/HP 10 mm 1.7 ul, 500 nl flow cell

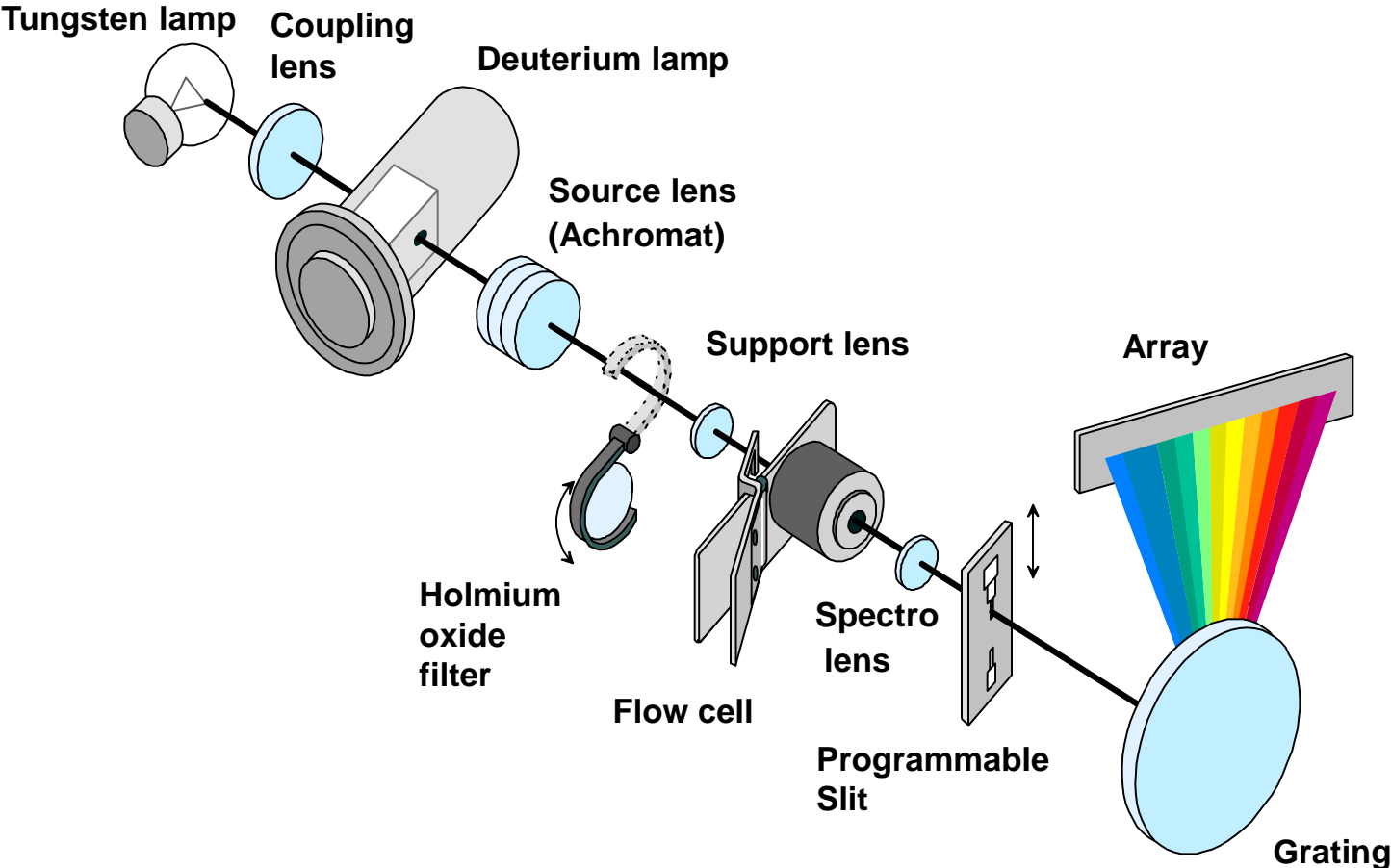
Frontal View – DAD/MWD

Tungsten Lamp

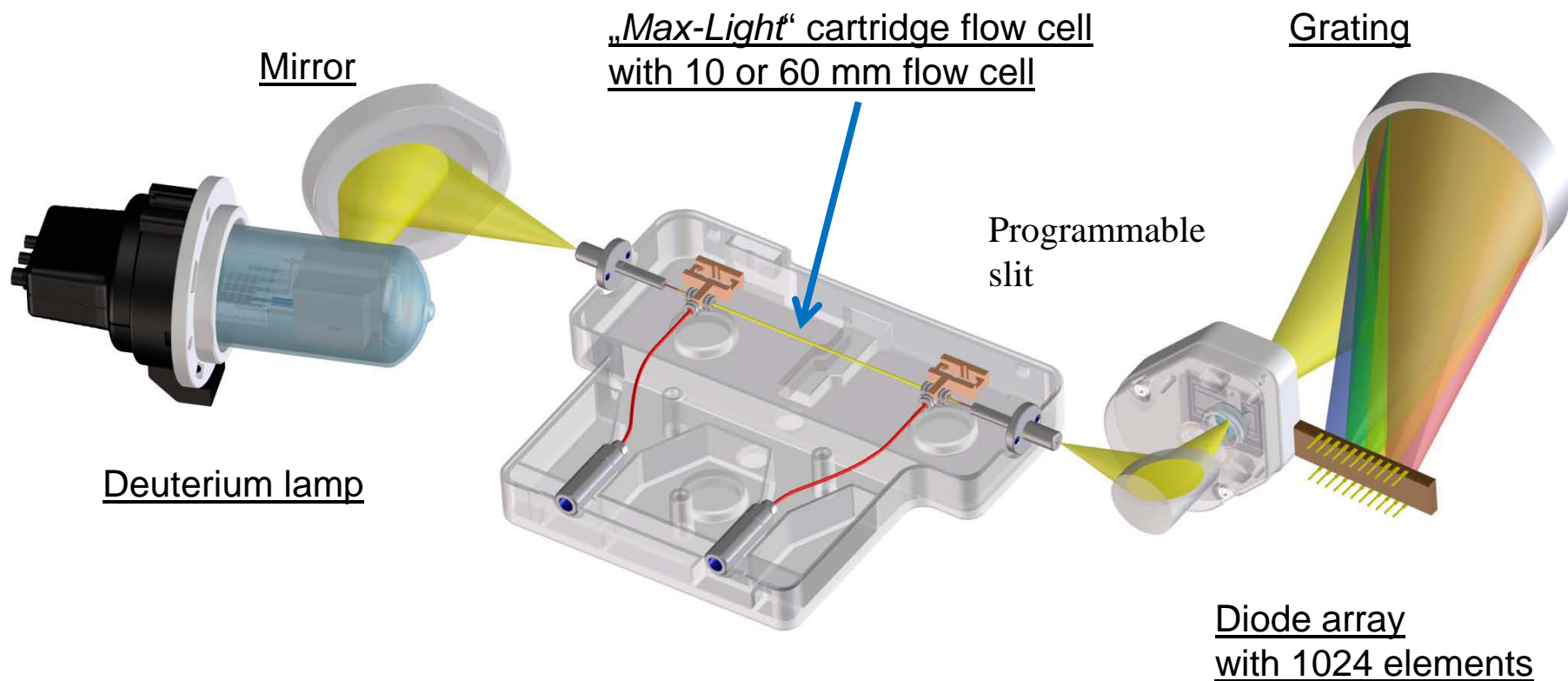
Deuterium Lamp



Optical Path - HP 1200



1260 & 1290 Infinity Diode Array Detectors



Optofluidic Waveguides: Max-Light Flow Cells

Total-internal reflection in a non-coated fused silica fiber

Diode Array Parameters - Review

Signal Storage

Spectral Storage Parameters

Response Time

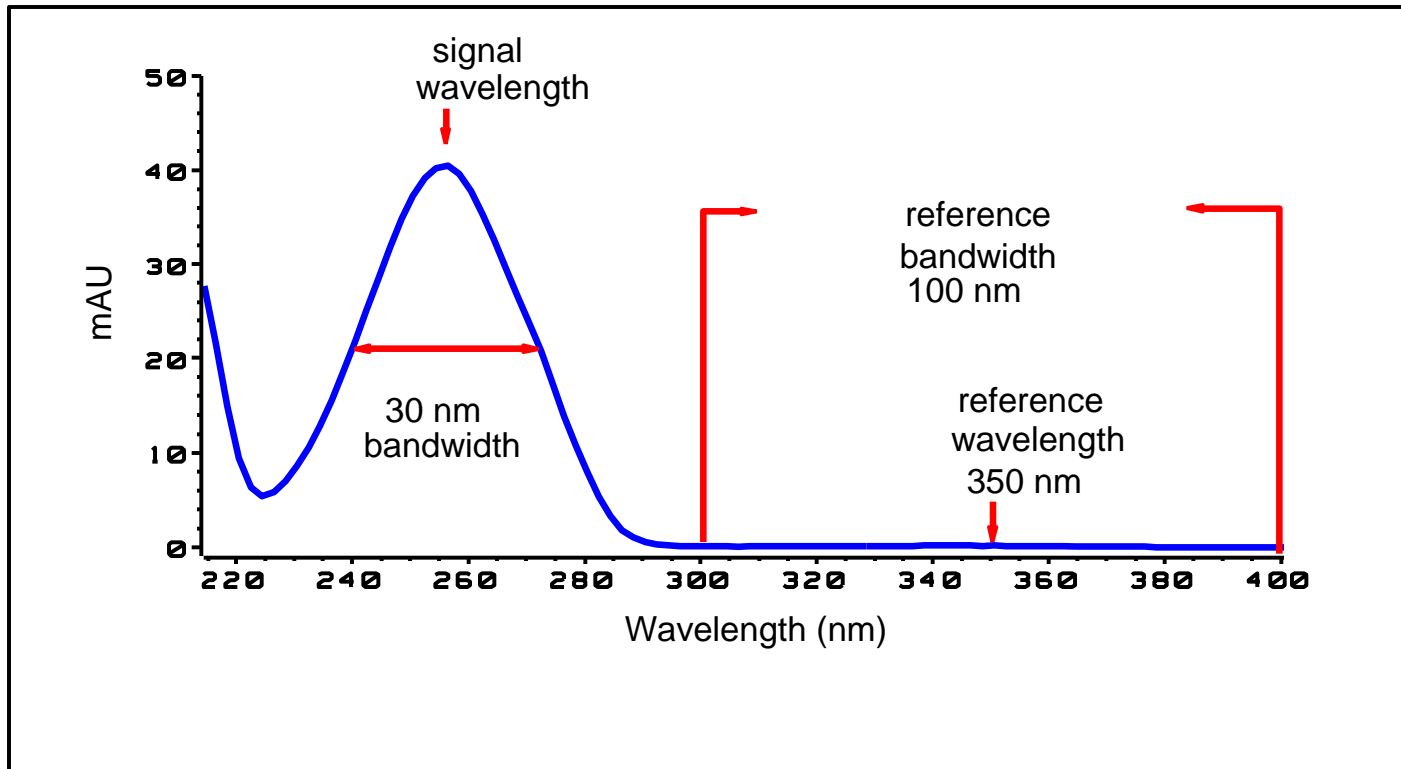
The screenshot shows the 'DAD Signals : Instrument 2' dialog box with the following settings:

- Signals:**

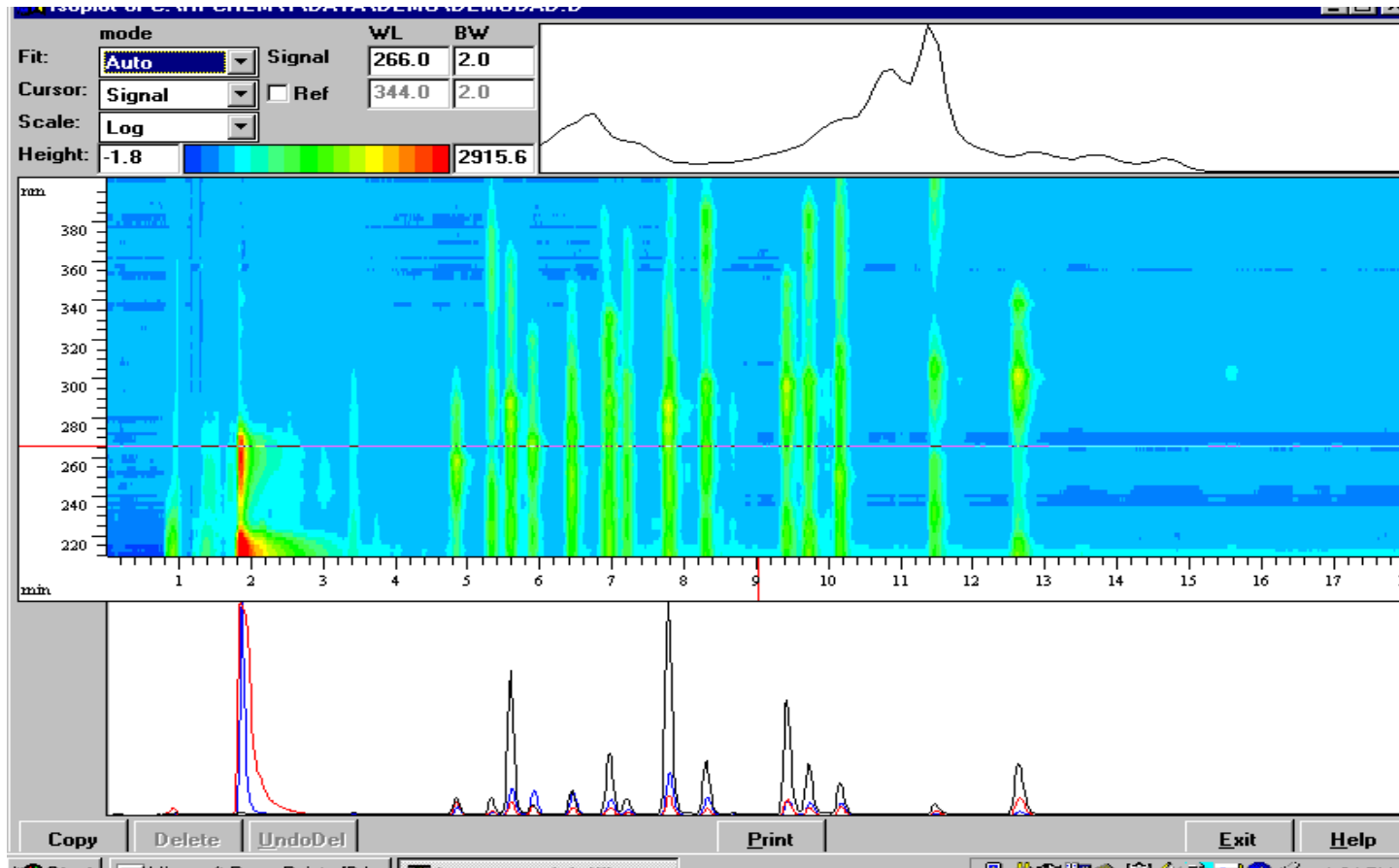
Store	Sample_Bw	Reference_Bw	nm
A: <input checked="" type="checkbox"/>	254	20	450 100
B: <input checked="" type="checkbox"/>	230	20	360 100
C: <input type="checkbox"/>	210	8	360 100
D: <input type="checkbox"/>	230	16	360 100
E: <input type="checkbox"/>	280	16	360 100
- Time:**
 - Stoptime: as Pump no Limit min
 - Posttime: Off min
- Required Lamps:**
 - UV
 - Vis
- Spectrum:**
 - Store: All
 - Range: 190 to 400 nm
 - Step: 2.0 nm
 - Threshold: 1.000 mAU
- Peakwidth (Responsetime):** > 0.1 min (2 s)
- Autobalance:**
 - Prerun
 - Postrun
- Slit:** 4 nm
- Margin for negative Absorbance:** 100 mAU

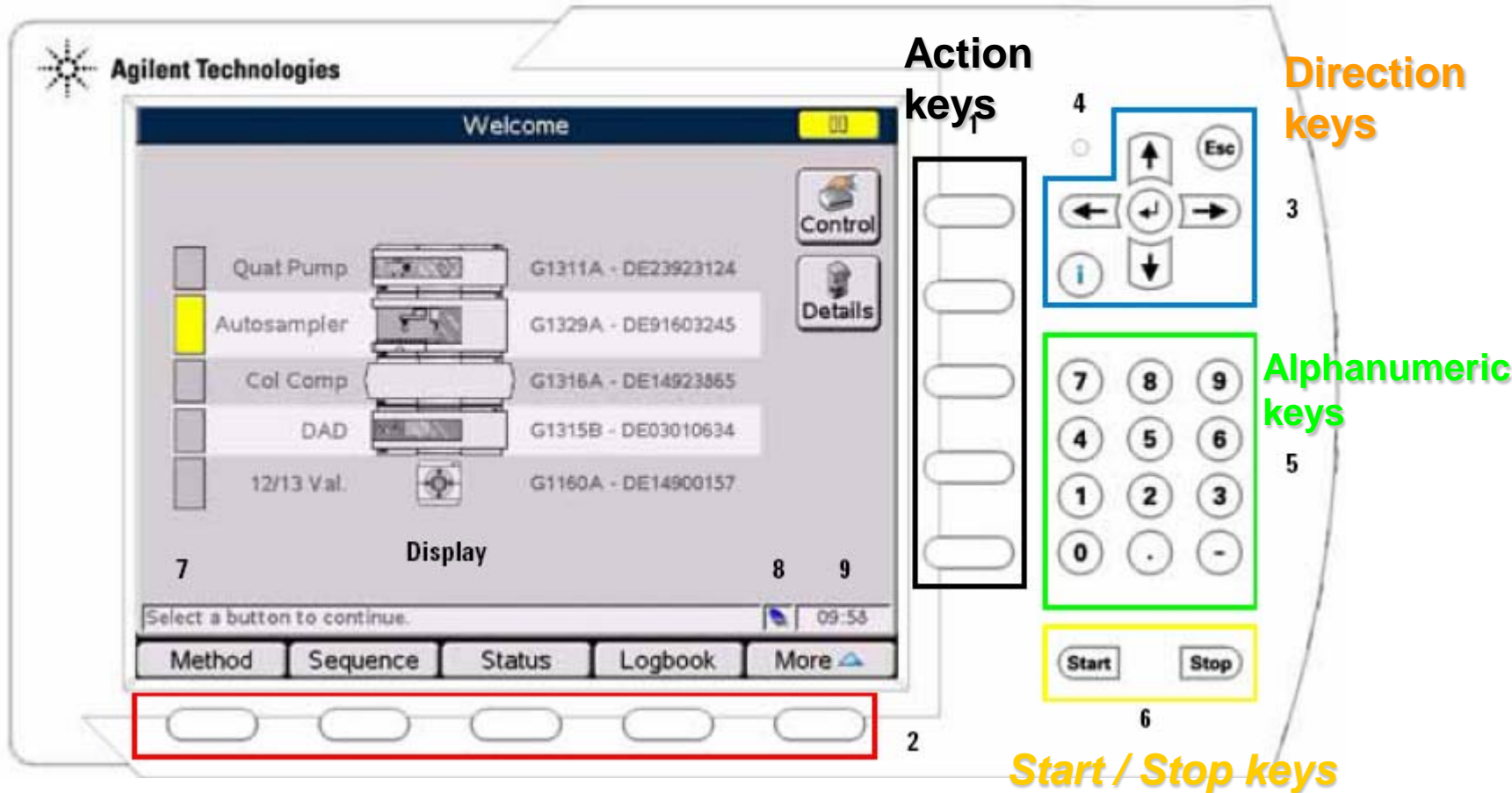
Buttons: Timetable ..., Total Lines: 1, OK, Cancel, Help

Sample Signal and Reference Optimization



Diode Array: Signal Extraction with Isoplot

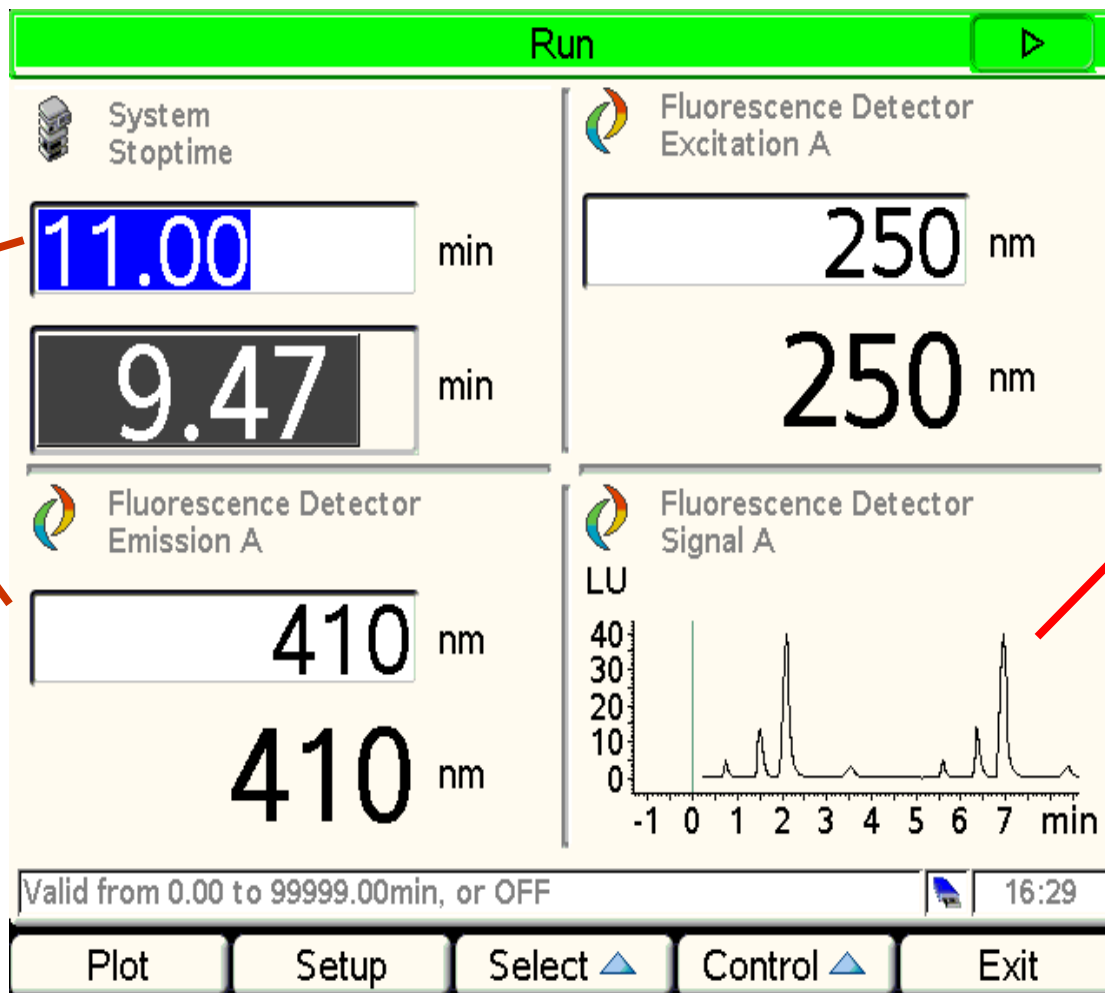




Navigation keys

Customizable Status Display

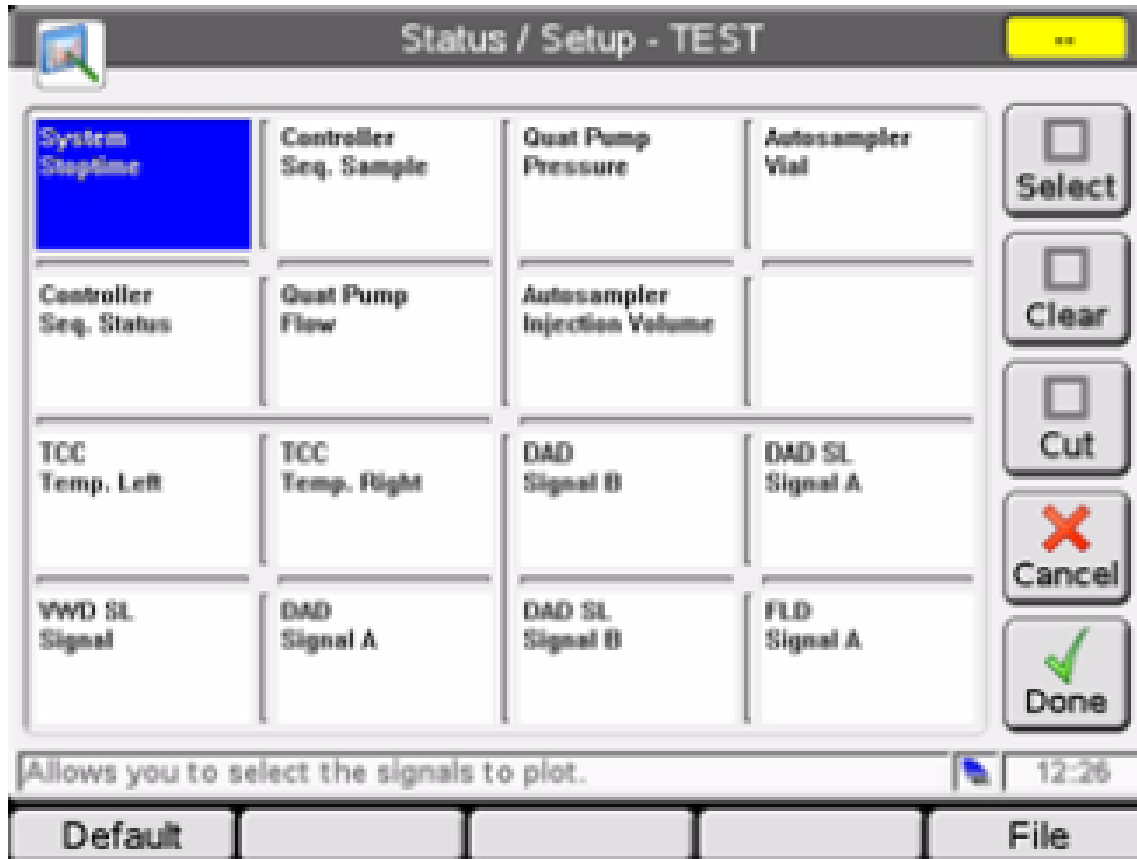
Direct
parameter
change



Signal
Plot

Tiles customizable by user

Press the **Setup** button.



allows the selection of a signal/parameter.

clears a selected field.

cuts a selected field to be pasted to another position.

leaves this screen without changes.

leaves this screen with all changes.

Default: takes signals from all modules

File: load/save a setup.

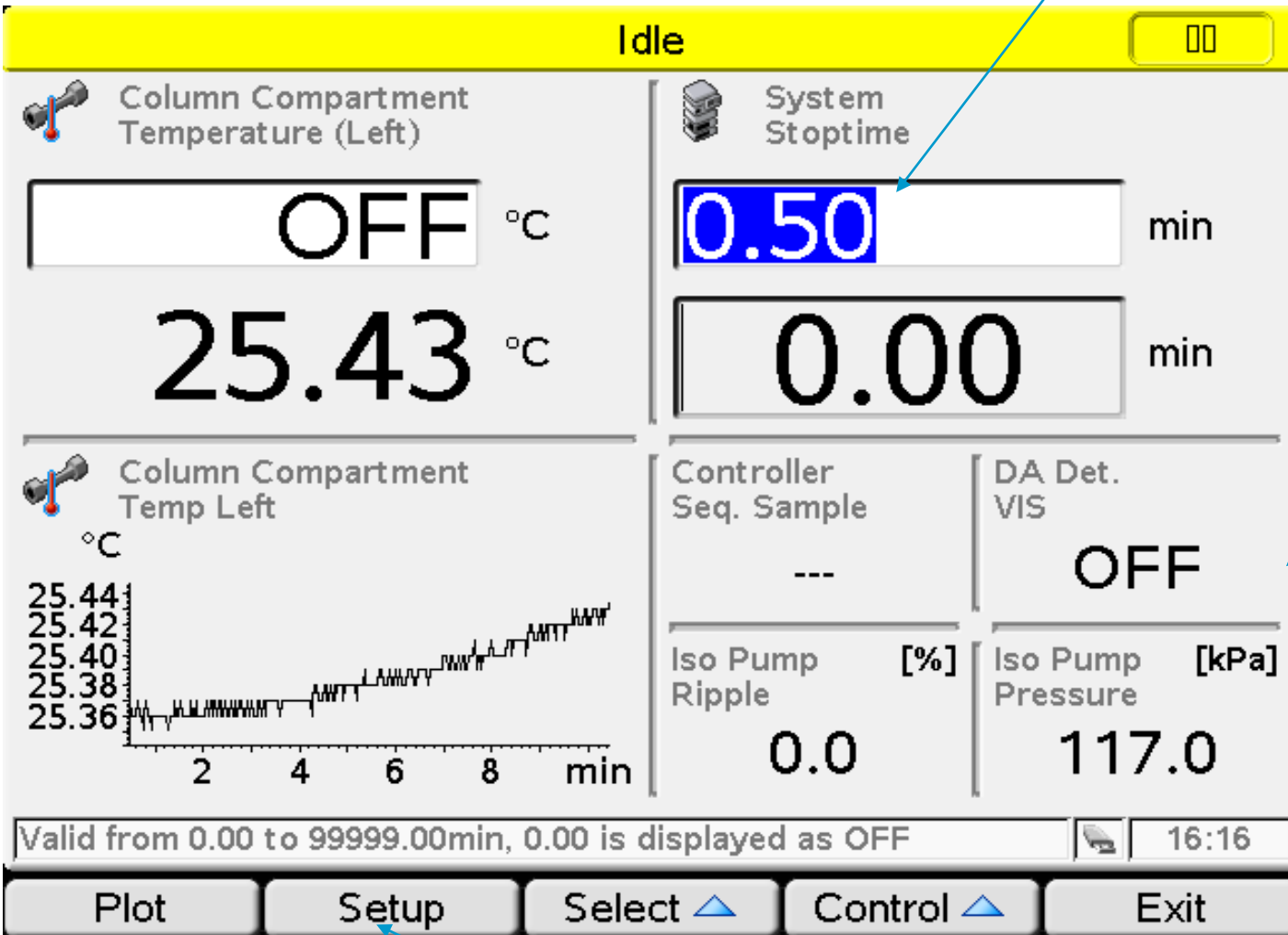
Setup of a Status Information Screen

Idle			
System Stoptime [min] 0.00	Controller Seq. Sample ---	Quat Pump Pressure [bar] 0.1	Autosampler Vial 0
Controller Seq. Status Idle	Quat Pump Flow [...] 0.000	Autosampler Injection Volume [µl] 0.00	
TCC Temp. Left [°C] 27.17	TCC Temp. Right [°C] 26.62	DAD Signal B [mAU] -1.702	DAD SL Signal A [mAU] 1.289
VWD SL Signal [mAU] 0.000	DAD Signal A [mAU] -35.75	DAD SL Signal B [mAU] -3.819	FLD Signal A [LU] 0.000
Allows you to modify the network configuration.			12:25
Plot	Setup	Select ▲	Control ▲
		Exit	

When the Status Information screen has not been setup before, it will show from each module in the system one or more signals/parameters

Customizable Status Display

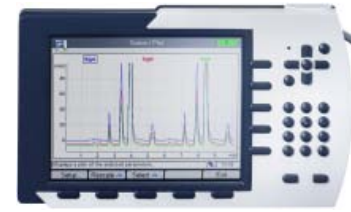
Direct parameter change



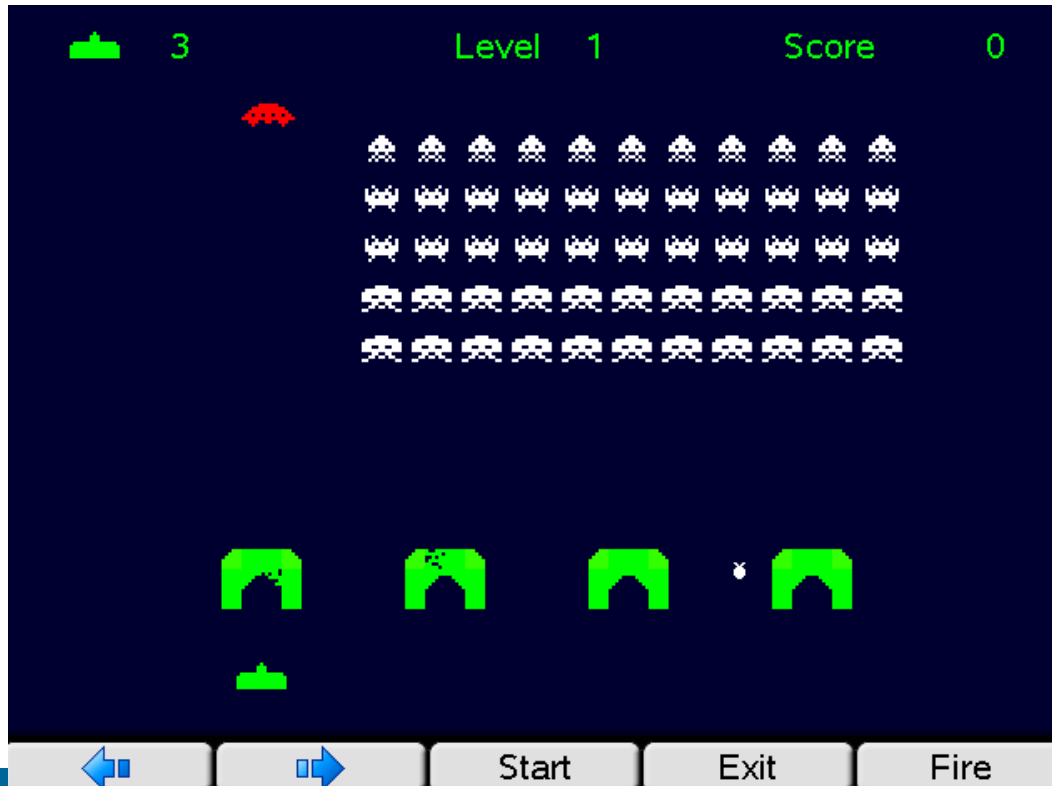
Tile with 4 subtiles

Screen customizable

Agilent 1200 Series Instant Pilot



Unsupported Feature – Space Invaders



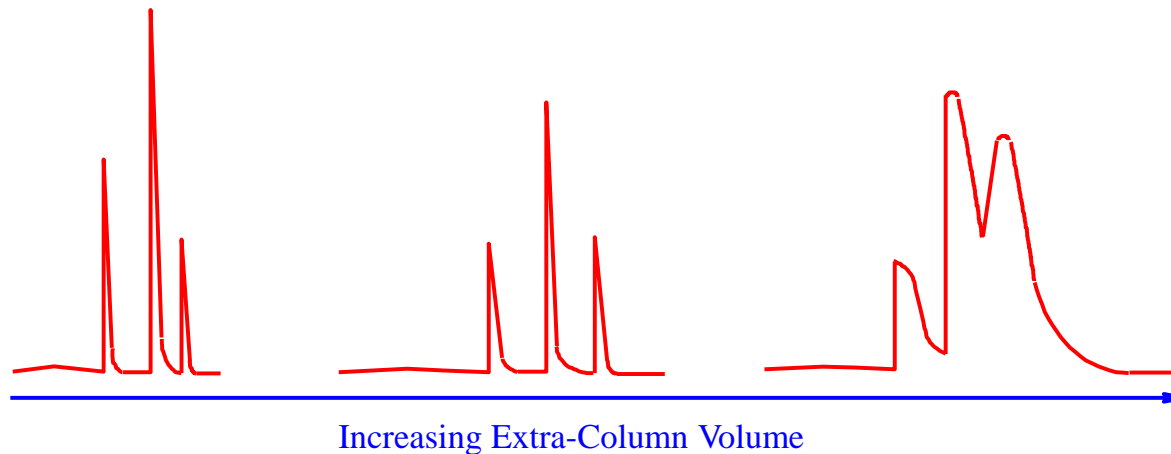
Welcome screen

Enter

3 5 1 1 3 7

H O B B I T

Extra-Column Dispersion



- Use short, small internal diameter tubing between the injector and the column and between the column and the detector.
- Make certain all tubing connections are made with matched fittings.
- Use a low-volume detector cell.
- Inject small sample volumes.

Dispersion is effected by

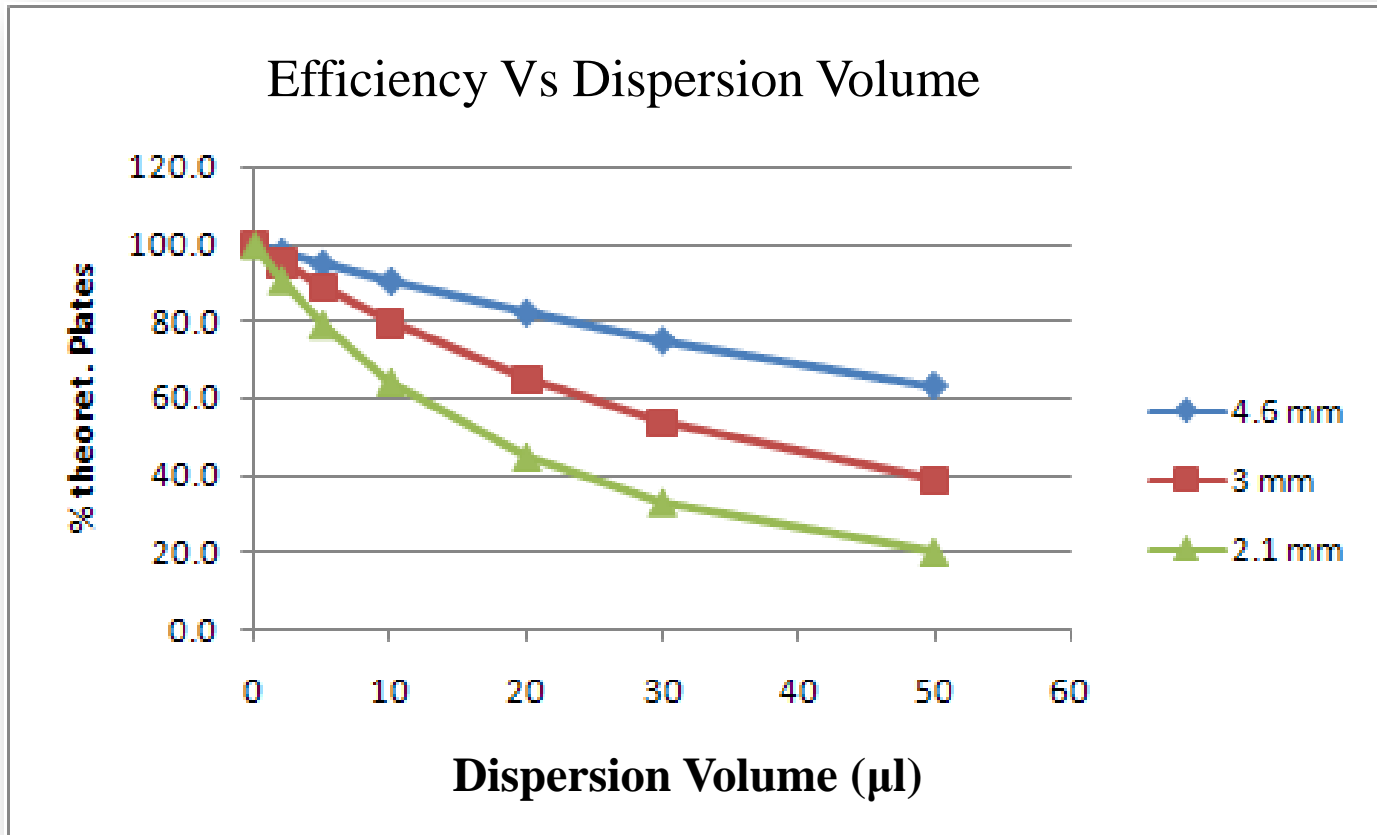
Tubing (amount and id)

Connections (number of)

Flow cell volume (larger >)

The goal is to have the smallest id tubing, least number of connections, and the lowest flow cell volume. This will give the best number for dispersion.

Dispersion



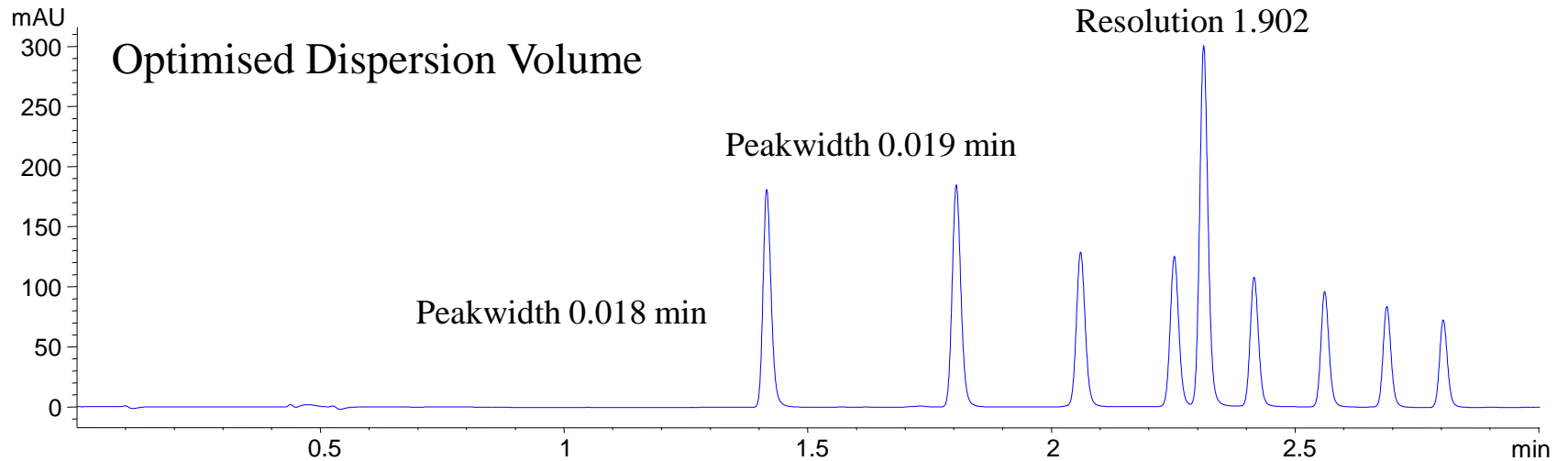
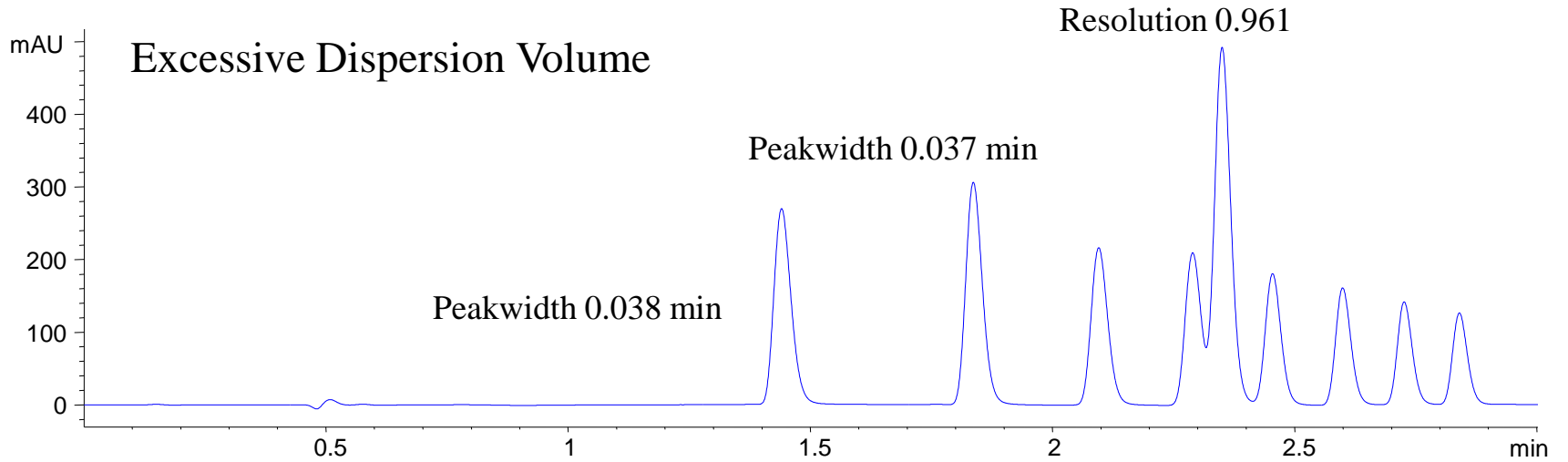
Flow rates converted for Column i.d.

4.6 mm Column	1.00 mL/min
3.0 mm Column	0.43 mL/min
2.1 mm Column	0.21 mL/min

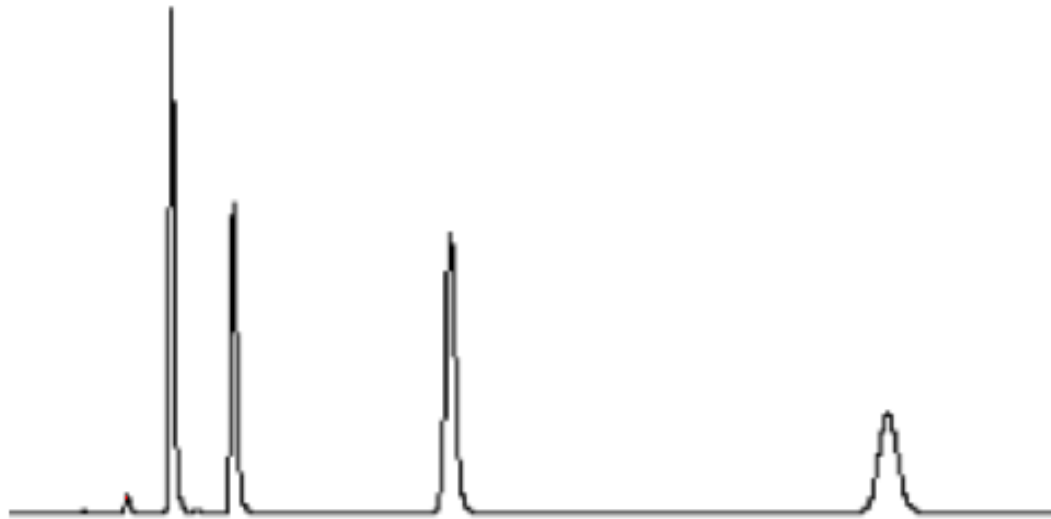
Column Dimension

Length = 50 mm
Particle size = 1.8 µm

Dispersion

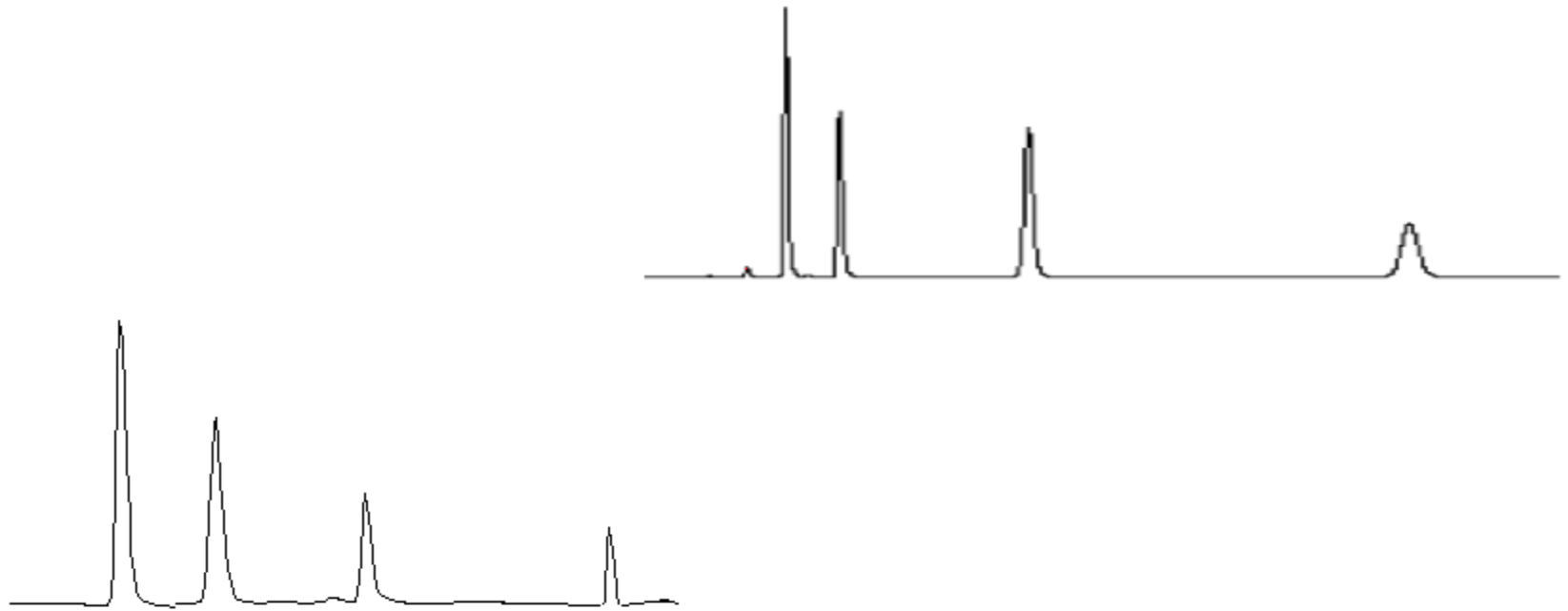


A typical chromatogram



Peak Tailing

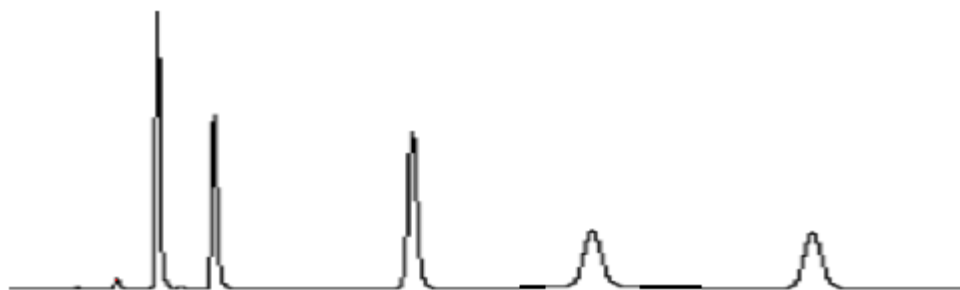
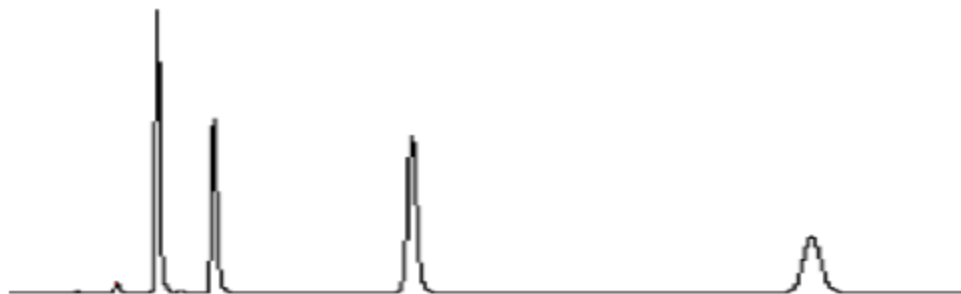
Injector Seal Failure



- Overdue instrument maintenance can cause peak shape problems.

Broad Peaks

Unknown “Phantom” Peaks



- The extremely low plates are an indication of a very late eluting peak from the preceding run.

Determining the Cause of Peak Tailing

- Evaluate mobile phase effects - alter mobile phase pH and additives to eliminate secondary interactions
- Evaluate column choice - try column with high purity silica or different bonding technology
- Reduce sample load
- Eliminate extra-column effects
- Flush column and check for aging/void

Questions?

