

# From Instrument to Column

## Tracking Down the Problem

Mark Powell

Applications Engineer

Columns and Supplies Technical Support



# Troubleshooting Topics

## System pressure

- Increased pressure
- Low pressure
- Pressure fluctuations

## Peak shape

- Tailing
- Broadening
- Fronting
- Peak splitting and doubling

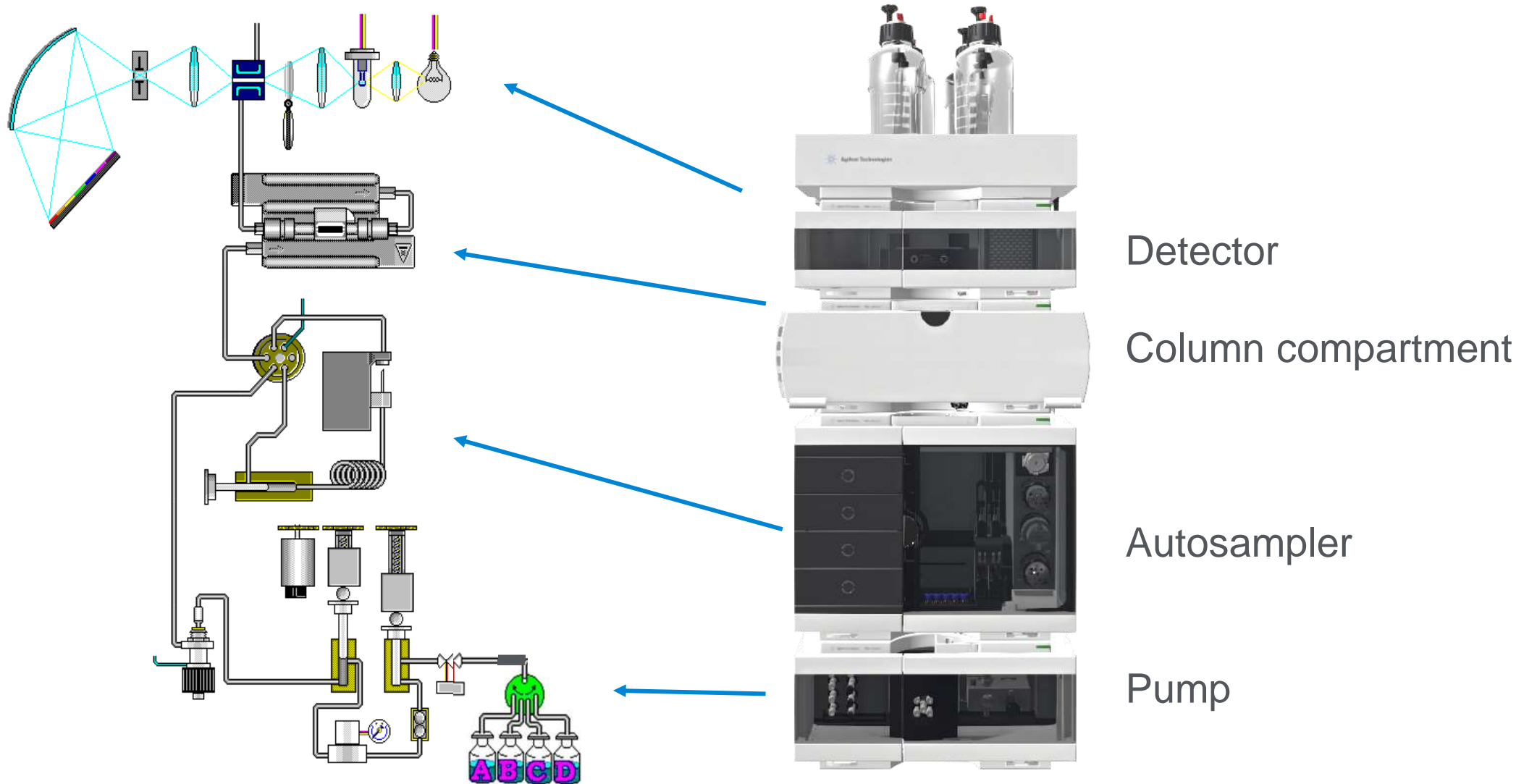
## Separation

- Changing retention time
- Loss of resolution

## Detection

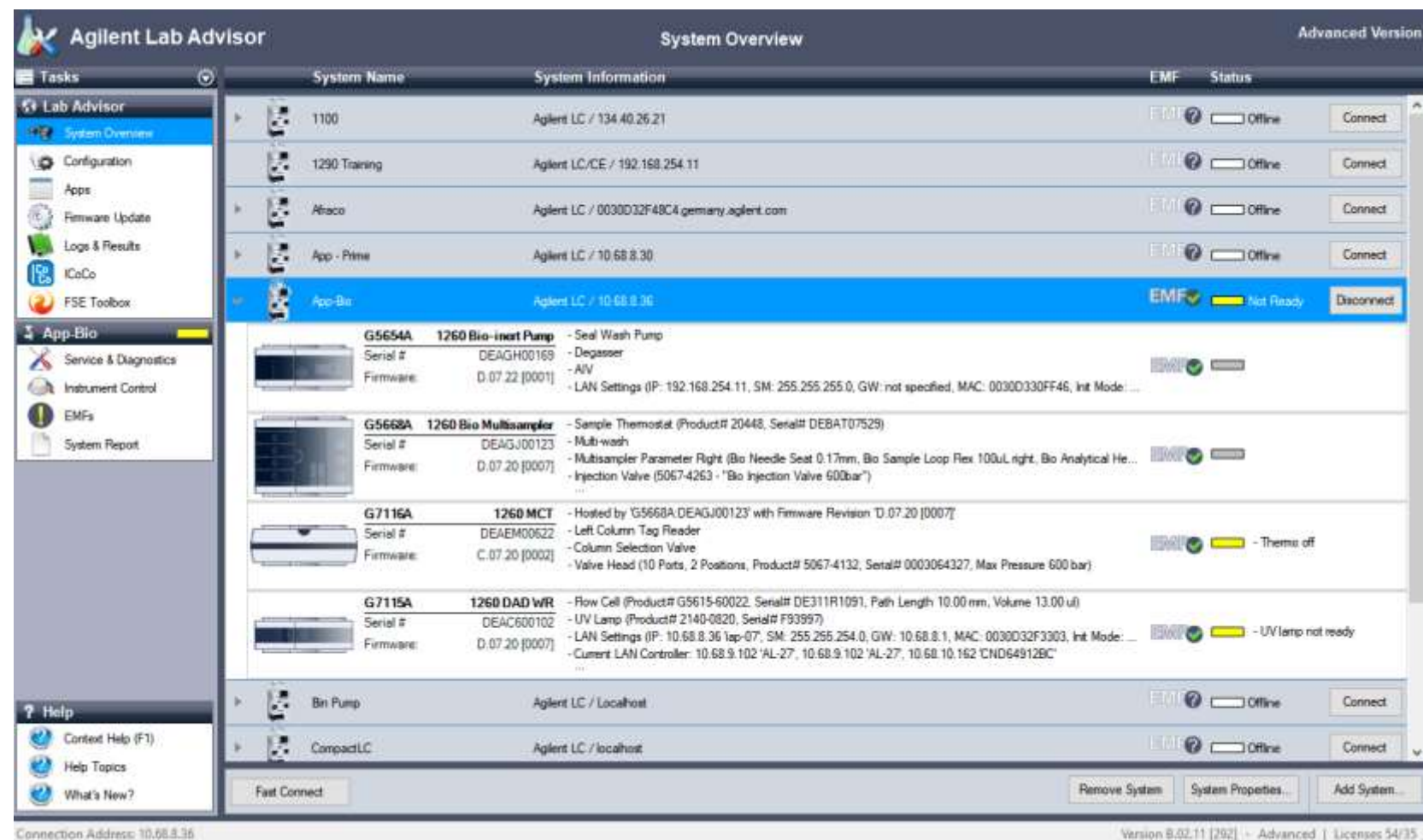
- Noisy baseline
- Reduced intensity or sensitivity
- Drifting baseline

# Understand Your LC System and Follow the Flow Path



# Agilent Lab Advisor

- Tools for calibration, diagnosis, and maintenance
- Daily instrument tests
- General calibration and maintenance procedures
- Advanced version also available for expert level troubleshooting
- EMF (Early Maintenance Feedback) shows the number of valve switches or pumped solvent, etc.

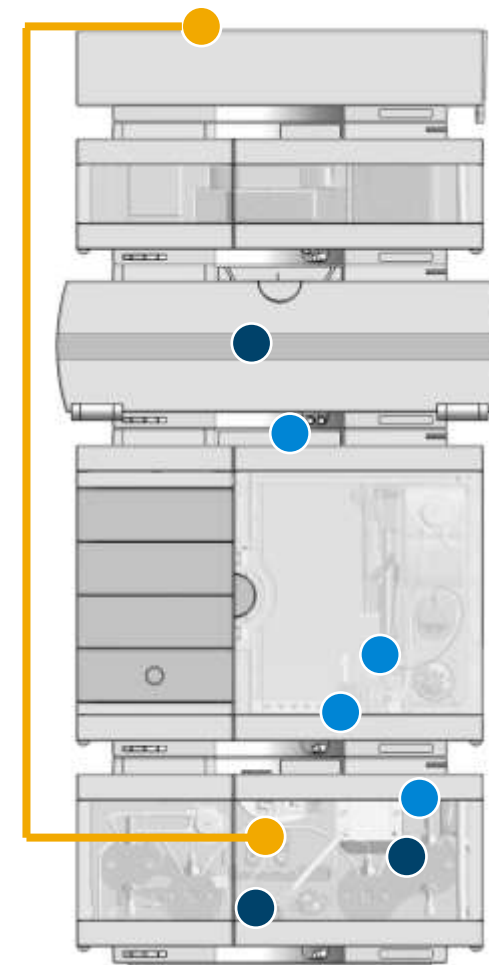


# System Pressure

# Changes in System Pressure

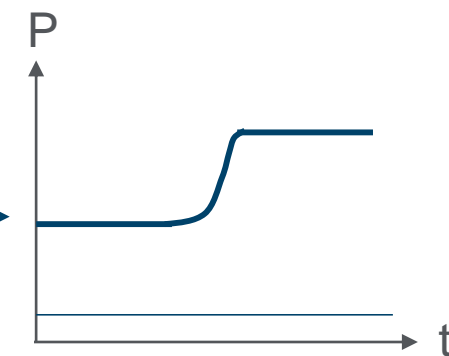
## Increased Pressure / Overpressure and Blockages

Potential cause		Recommended action
●	Clogging of filter frits in the high-pressure flow path	<ul style="list-style-type: none"><li>Identify the culprit by logical elimination process and replace affected part.</li><li>Use clean solvent (e.g. prefiltered solvent)</li><li>Prevent algae growth in water</li></ul>
●	Plugging of capillaries, needles and needle seats	
●	Wrong solvent	<ul style="list-style-type: none"><li>Check for correct mobile phase</li><li>Solvent reservoir and tube connections</li></ul>

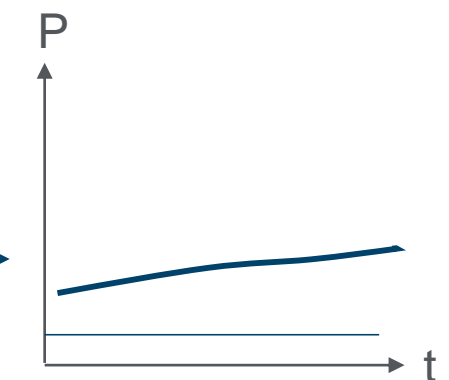


# Blockages and Clogging

Characteristics	
Parts affected	<p>Blockages:</p> <ul style="list-style-type: none"> <li>Capillaries, needle and needle seat</li> <li>Detector flow cells</li> </ul> <p>Clogging:</p> <ul style="list-style-type: none"> <li>Filter frits (inline filter, column filter)</li> </ul>
Characteristic	<div>●</div>
Identification	<ul style="list-style-type: none"> <li>Start by disconnecting the capillary at the column inlet</li> <li>Install test setup with restriction capillary</li> <li>Continue disconnecting capillaries, one-by-one, moving back toward the pump</li> </ul>
Possible Root Cause	<ul style="list-style-type: none"> <li>Debris from mechanically worn parts (needle seat material, rotor seal at injection valve)</li> <li>Coring of vial septa material</li> </ul>
Instant action / First aid	<ul style="list-style-type: none"> <li>Backflush affected part</li> <li>Replace part</li> </ul>
Preventive measures	<ul style="list-style-type: none"> <li>Replace wear parts in time; apply proper preventive maintenance schedules</li> <li>Use high quality septa</li> <li>Install inline filters</li> </ul>

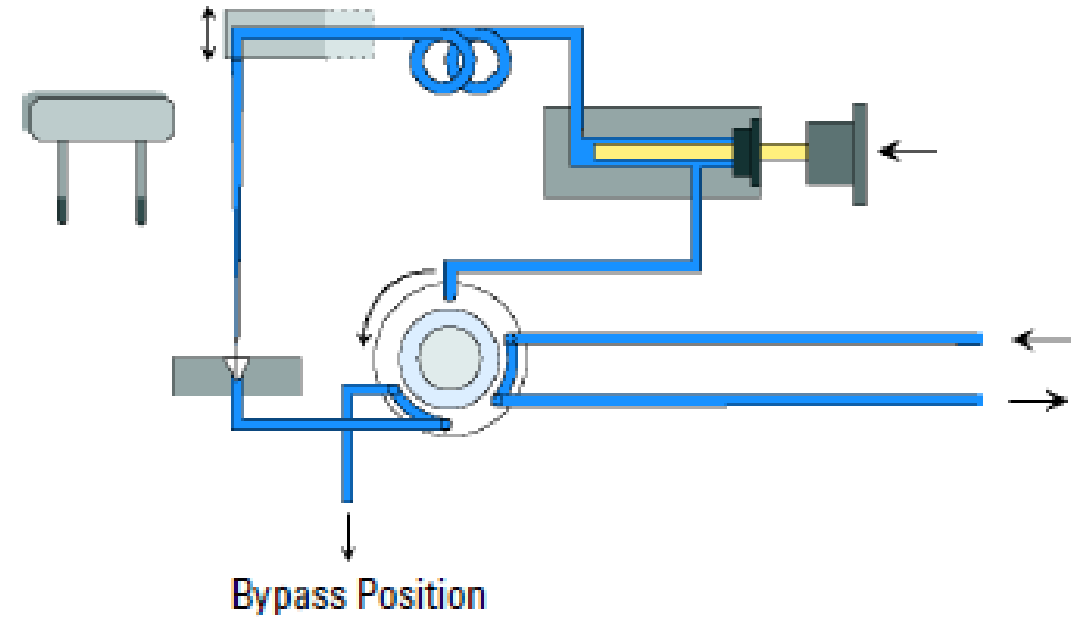
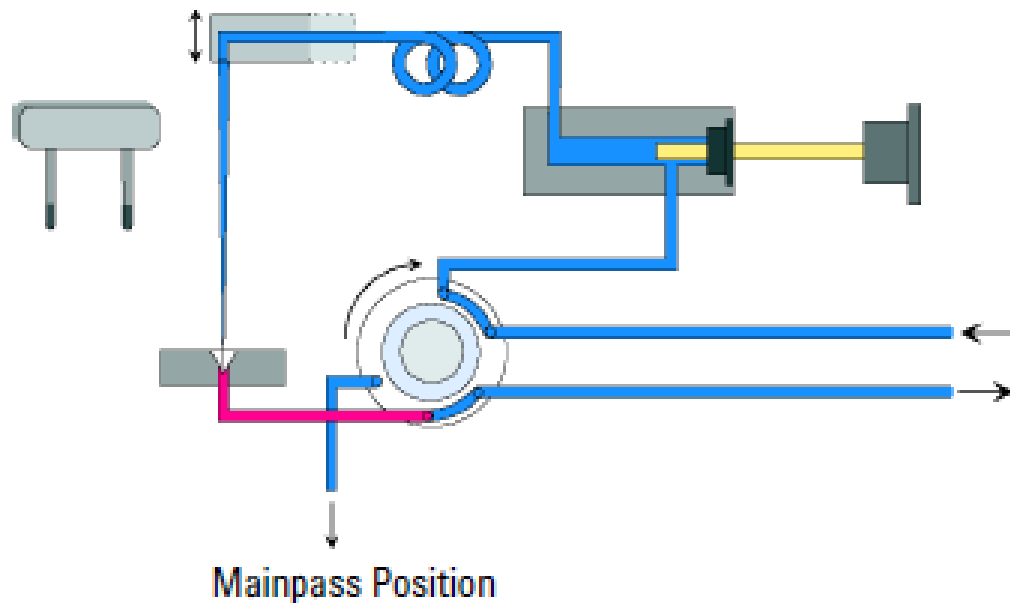


**Blockages:** instant pressure increase step



**Clogging:** constant pressure increase over time

# Checking for Blocked Needle Seat



# Checking for Blocked Needle Seat

- Track the typical operating pressure for a given application
- To troubleshoot, create high pressure on the system by turning on flow
- While the pressure is climbing, move the sampler to the “Bypass” position
- Watch the pressure when the valve switches to “Bypass”
- If the pressure drops immediately, then the source of the high pressure is in the portion of the flow path specific to the bypass position.
- Needle Seat
  - The most commonly clogged piece of tubing in the LC
  - Where sample first meets the separation solvent
- Needle
  - Less commonly clogged
  - Watch for issues with septa

The screenshot displays the 'Control' panel of an Agilent instrument. Under the 'Method Parameters' section, the 'Injection valve (Single Needle)' is set to 'Mainpass' (indicated by a selected radio button). Below this, the 'Needle Wash' section is expanded, revealing 'Set Needle Wash Multi Mode Parameters'. This section includes a dropdown for 'Solvent 1', a 'Channel' dropdown set to 'A', a 'Time [s]' input field set to '30', a 'Seat back flush' dropdown set to 'Off', a 'Needle wash at flush port' dropdown set to 'On', and a 'Solvent Name' input field.

# Checking for Blocked Needle Seat

## Loop

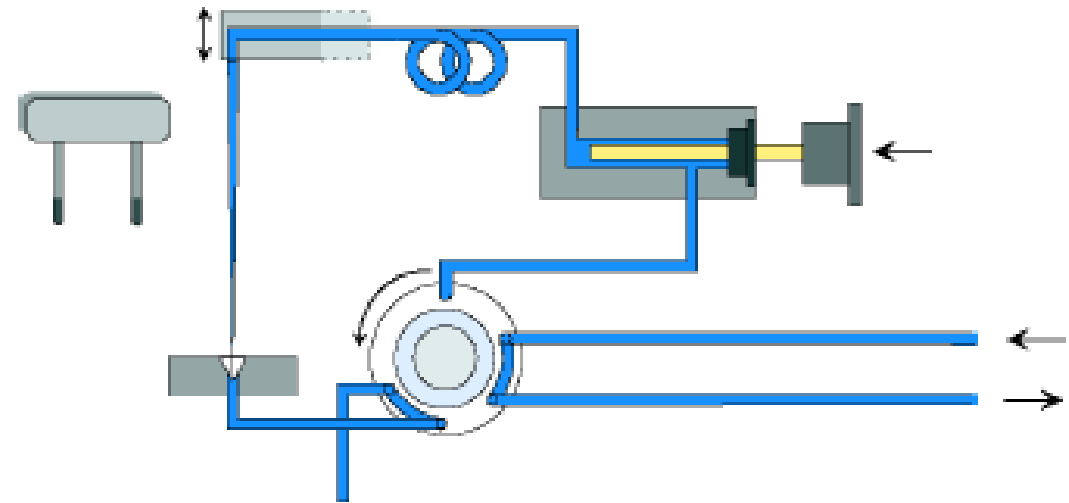
- not commonly clogged
- watch for issues with sample

## Metering Head

- never exposed to sample
- consider solvent issues

## Injection Valve

- most commonly the rotor seal
- look for scratches on stator face



# Checking for Blocked Needle Seat

- In well plate samplers, bottom sensing gives the most consistent position
- But this isn't recommended if there's debris in the bottom of the sample vial
- Driving the needle into debris may result in a clogged needle or seat
- For vialsamplers, a zero offset is approximately 2 mm from the bottom of a 2 ml vial

Needle height position without bottom sensing:

G1367E/G4226A 54 vial tray = 4 mm


G1367E/G4226A 100 vial tray = 2.5 mm

G7167X 54 vial tray = 5 mm

# Locating a Clog

- If the pressure doesn't drop when the valve switches to "Bypass" the issue is likely outside the sampler
- Other easily accessible points in the flow path to check are:
  - With 1260 model pumps open the manual purge valve, the pressure should drop to between 0 and 5 bar. If the pressure is higher the PTFE filter may be clogged.
  - With 1290 model pumps purging is done through an automated valve activated through software. 1290 Binary pumps have the same PTFE filter, 1290 quaternary pumps have a 5 µm filter frit.



**G7120A 1290 High Speed Pump**

Serial #	DEA0000000
Firmware:	B.07.20 [0007]

**Controls**

**Control**

Pump: ☐ On ☒ Off ☐ Standby ☐ Initializing

Purge + Prime

Purge Process: ☐ On ☒ Off

Prime: ☐ On ☒ Off

# Locating a Clog



PTFE replacement on a 1260 pump:

1. Remove pump outlet and purge waste tubing
2. Unscrew the purge valve using a 14mm wrench
3. Remove the gold seal cap
4. Remove the frit
5. Install the new frit, slot side up
6. Replace the gold seal cap
7. Reinstall the valve

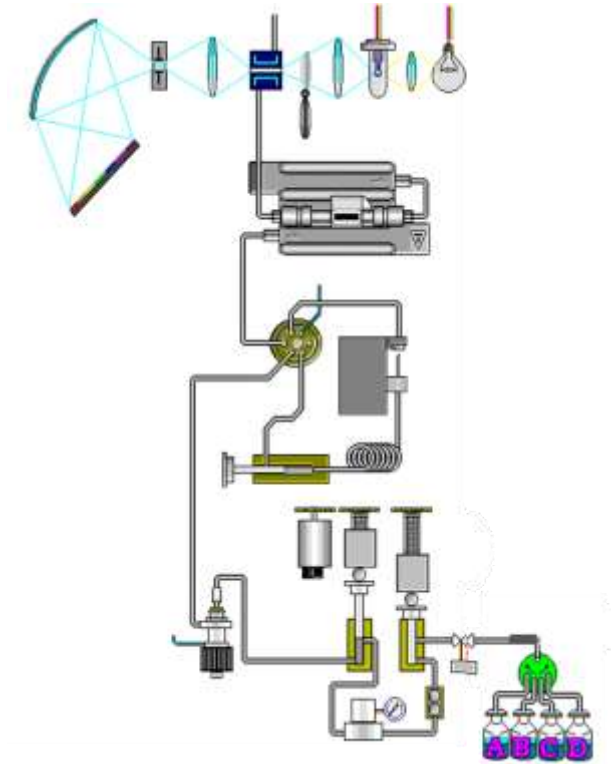
Realign the waste tubing in the correct orientation during installation.

# Locating a Clog

- If the pressure doesn't drop when the valve switches to "Bypass" the issue is likely outside the sampler

Other easily accessible points in the flow path to check are:

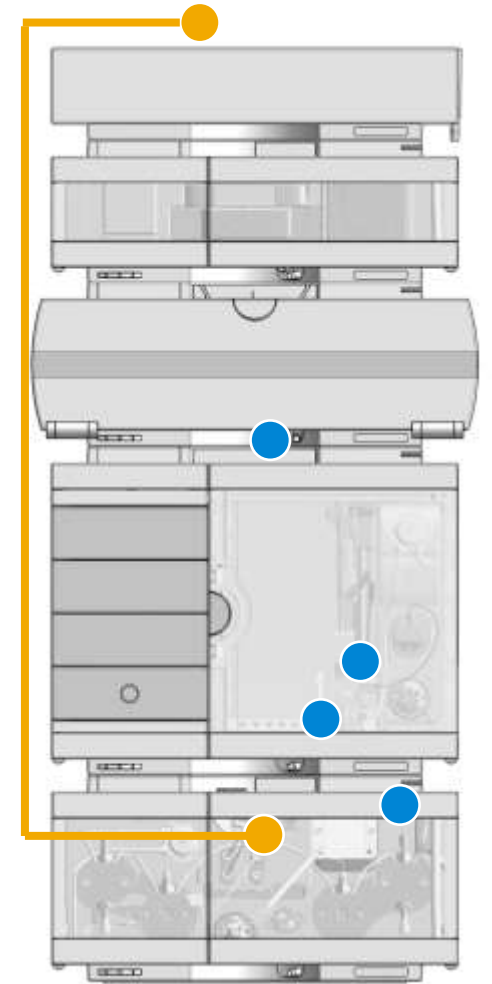
- Open the fitting at the inlet of the column. Pumping 1 ml/min of water through an Agilent LC with 0.17mm id tubing typically shows a pressure of 40 bar. If the pressure is much higher than this a capillary may be clogged. If the pressure appears "normal" the issue may be with the column.
- Clogs are located by opening a fitting, typically at most a half turn. If the pressure drops the clog is downstream from the fitting or towards the detector. If pressure remains high the clog is upstream or towards the pump.



# Changes in System Pressure

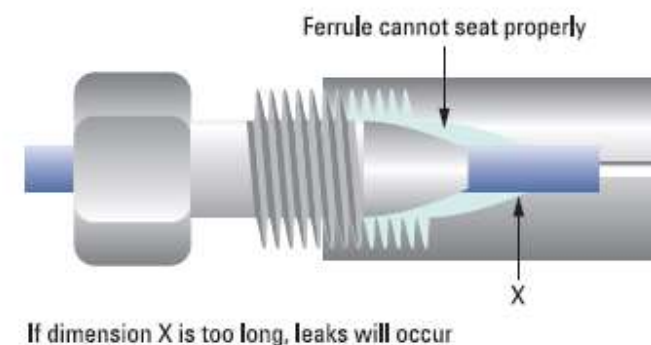
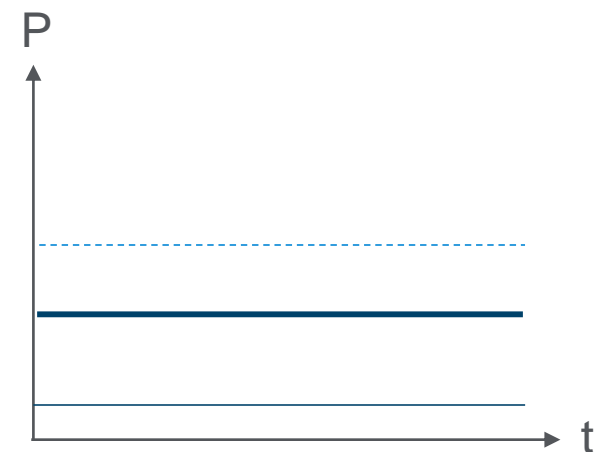
## Low pressure

	Potential cause	Recommended action
●	Leak in high-pressure flow path	<ul style="list-style-type: none"><li>• Visual inspection of flow path</li><li>• Instrument diagnostic tests</li></ul>
●	Wrong mobile phase	<ul style="list-style-type: none"><li>• Check for correct mobile phase</li><li>• Solvent reservoir and tube connections</li></ul>



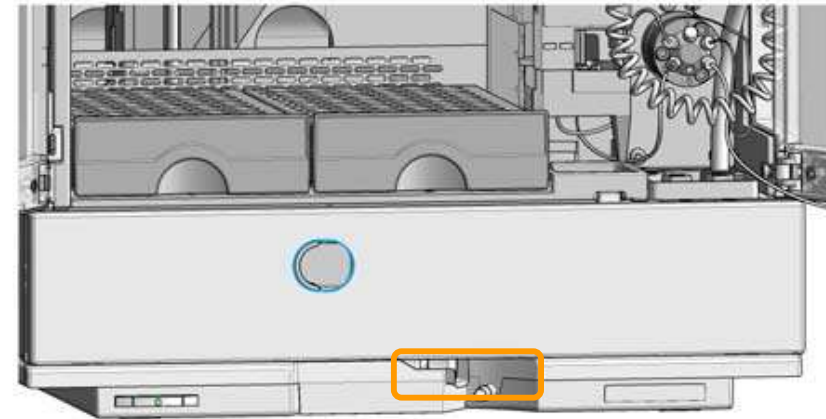
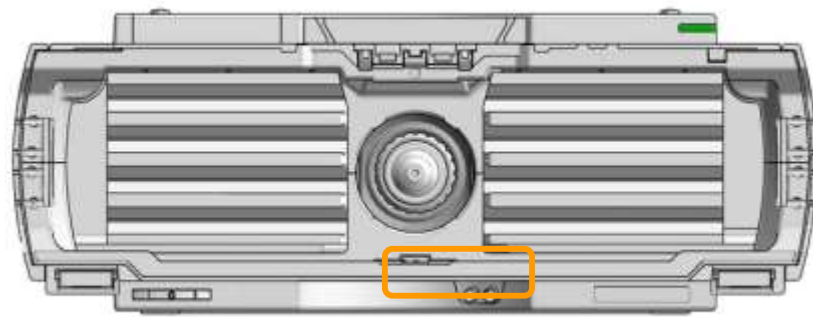
# Leaks

Characteristics	
Parts affected	<ul style="list-style-type: none"> <li>Potentially all parts in the flow path</li> <li>High potential at frequently operated fitting connections (e.g. column inlet) and parts with high mechanical stress (rotor seal, needle, and needle seat)</li> </ul>
Characteristic	<ul style="list-style-type: none"> <li>Lower pressure</li> <li>Potentially impacting retention times and peak shape</li> </ul>
Identification	<ul style="list-style-type: none"> <li>Drops of solvent or residues of salt</li> <li>System diagnostic tests</li> </ul>
Possible root cause	<ul style="list-style-type: none"> <li>Loose or bad fitting connections</li> <li>Cracked capillaries</li> <li>Worn needle and needle seat</li> </ul>
Instant action/first aid	<ul style="list-style-type: none"> <li>Replace affected parts</li> <li>Renew or redo fitting connection</li> </ul>
Preventive measures	<ul style="list-style-type: none"> <li>Use proper fitting connections</li> <li>Replace fittings and wear parts in time</li> </ul>

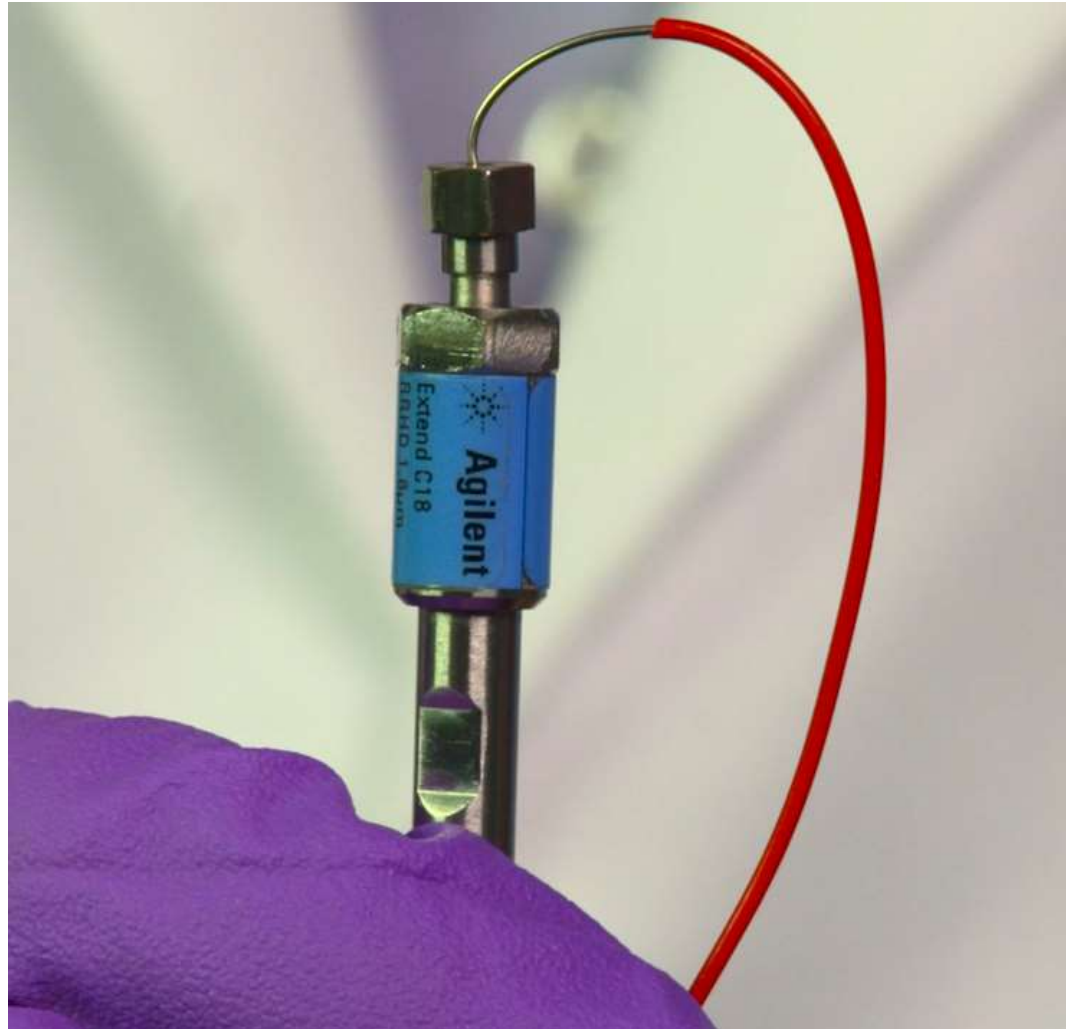


# How do I Locate a Leak?

- Each Agilent LC module is equipped with a leak sensor
- If liquid is detected the entire LC stack will shut down
- The LC will not start up again until the sensor has been dried and returned to temperature



# Overtightened Fittings



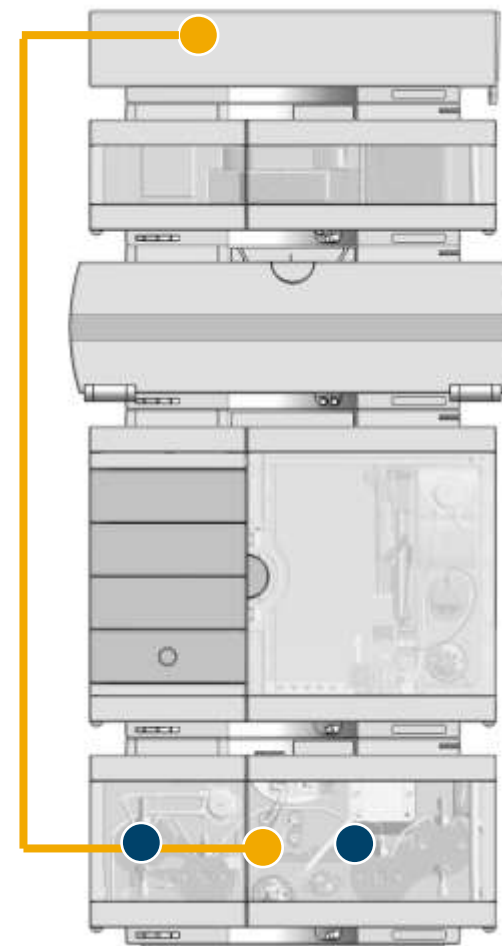
# Changes in System Pressure

## Pressure fluctuations

Potential Cause		Recommended Action
●	Air in the system	<ul style="list-style-type: none"><li>• Prime and flush instrument</li><li>• Check for sufficient solvent supply</li><li>• Check for correct plumbing (SSV/MCGV)</li><li>• Check for correct degassing</li></ul>
●	Malfunctions at pump head	<ul style="list-style-type: none"><li>• Perform pump head diagnostic tests LA</li><li>• Replace defective parts</li><li>• Implement proper maintenance schedule</li></ul>
●	Cavitation effects	<ul style="list-style-type: none"><li>• Check for flow restrictions (solvent bottle to pump head inlet)</li><li>• Clean or replace parts</li><li>• Verify that solvent supply is positioned above pump inlet</li></ul>

### Important to know

Pressure fluctuations typically also will impact the UV-signal due to refractive index effects.



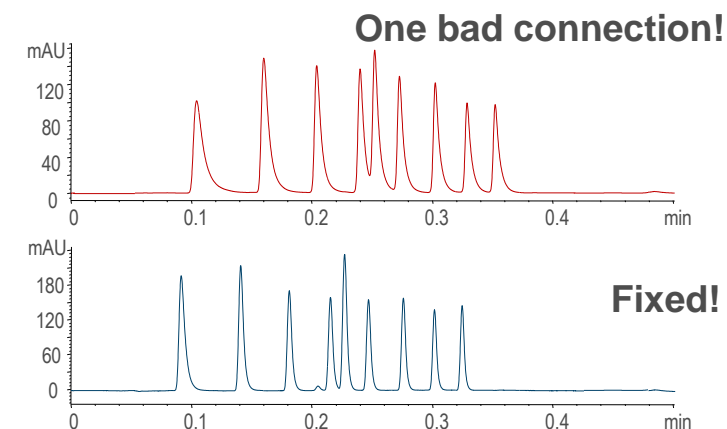
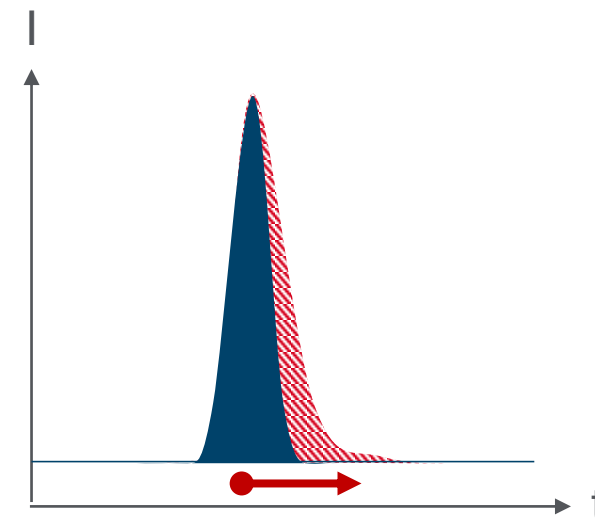
# Peak Shape

# Changes in Peak Shape

## Peak Tailing

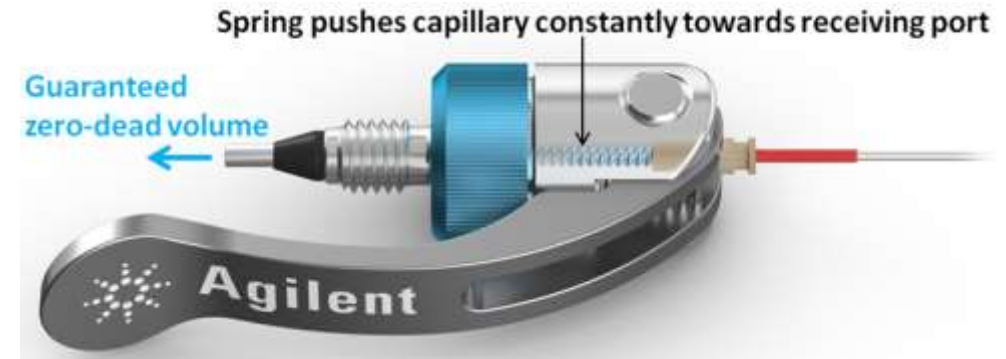
If applicable to some peaks	Recommended Action
Secondary interactions	<ul style="list-style-type: none"> <li>• Change pH</li> <li>• Change stationary phase</li> </ul>
Small peak eluting on tail of larger peak	<ul style="list-style-type: none"> <li>• Change selectivity (column, mobile phase)</li> <li>• Switch to methods with higher resolution (UHPLC, 2D-LC)</li> </ul>

If applicable to all peaks	Recommended Action
Poor tubing connections; high dispersion volume	<ul style="list-style-type: none"> <li>• Minimize number of connections</li> <li>• check connections / fitting condition and proper seat of fittings</li> <li>• use fittings with spring-load function</li> </ul>
Column damage	<ul style="list-style-type: none"> <li>• Use specialty, polymeric or sterically protected column</li> <li>• Column cleaning</li> </ul>



# InfinityLab Quick Connect and Quick Turn Fittings

- Spring loaded design
- Easy! **No tools needed**
- Works for all column types
- Reusable
- Consistent ZDV connection



## Quick Connect Fitting

- Finger tight up to 1300 bar
- Hand tighten the nut, then depress the lever

## Quick Turn Fitting

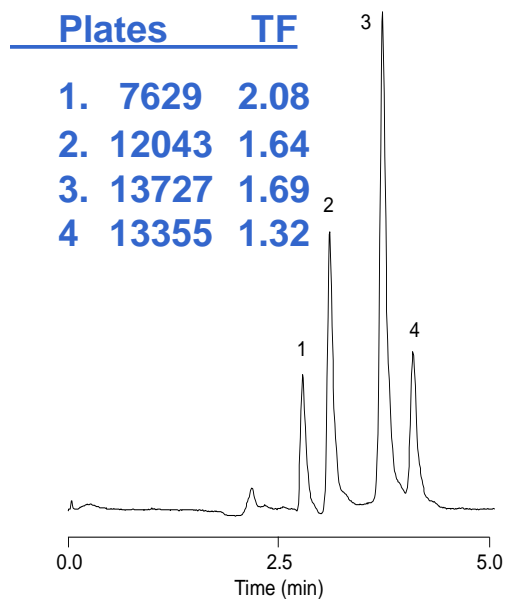
- Finger tight up to 400 bar
- Up to 1300 bar with a wrench
- Compact design



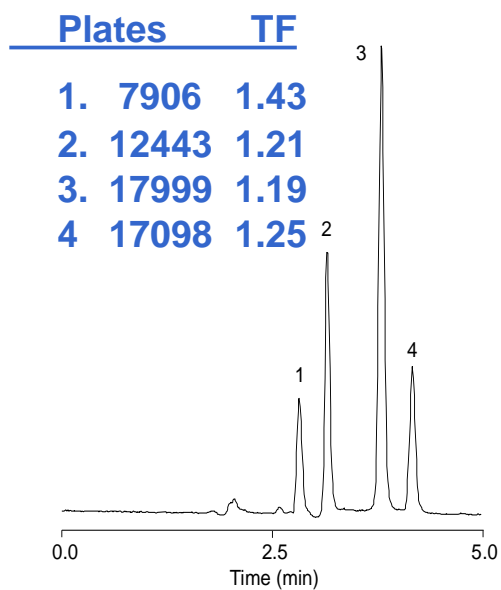
# Peak Tailing Column Contamination

Column: StableBond SB-C8, 4.6 x 250 mm, 5 mm      Mobile phase: 20% H<sub>2</sub>O : 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

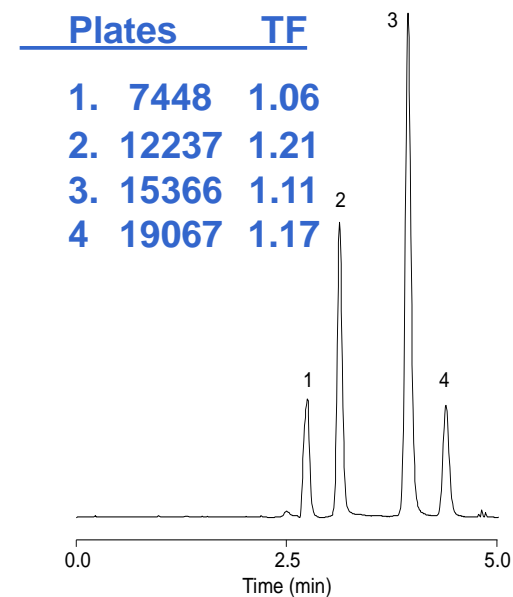
## QC test forward direction



## QC test reverse direction



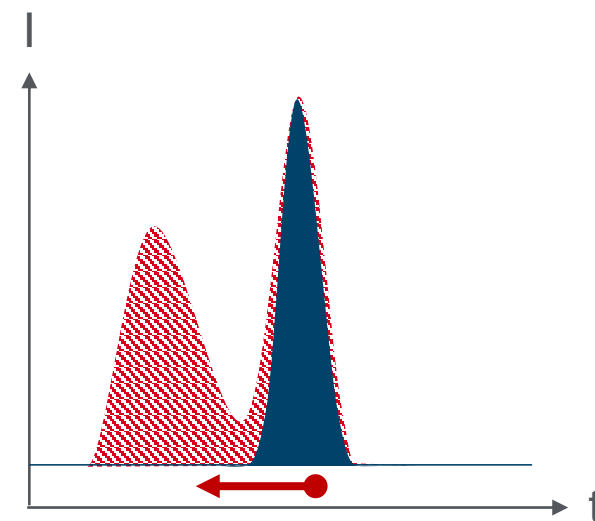
## QC test after cleaning 100% IPA, 35°C



# Changes in Peak Shape

## Peak splitting / doubling

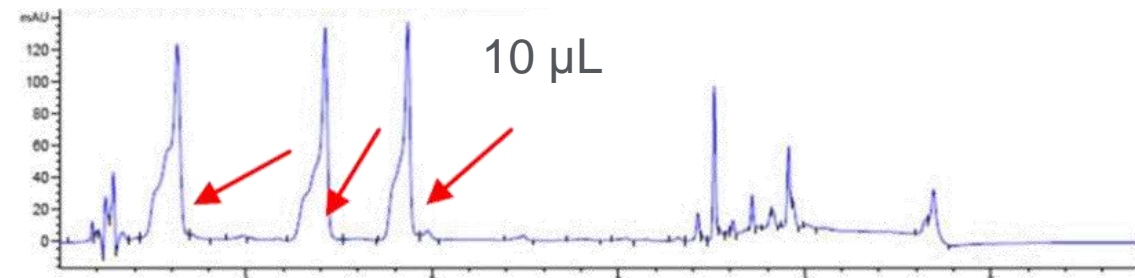
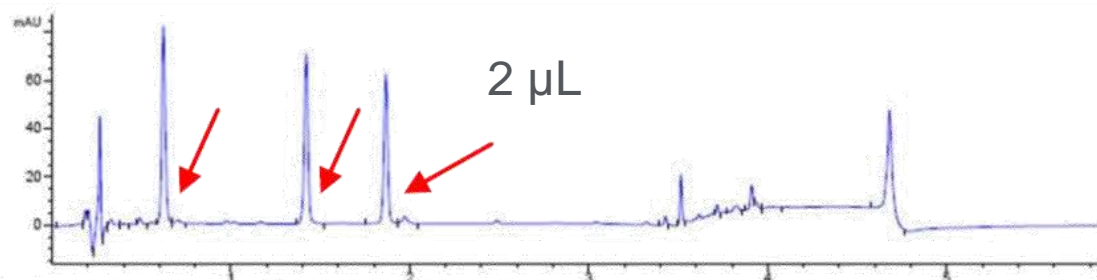
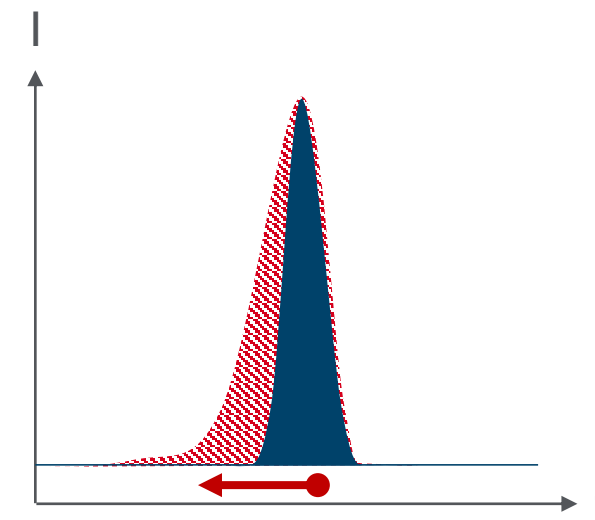
Potential Cause	Recommended Action
Partially plugged column frit	<ul style="list-style-type: none"><li>• Backflush column (if applicable)</li><li>• use inline filter</li><li>• use guard column</li></ul>
Column void	<ul style="list-style-type: none"><li>• Replace column</li><li>• use guard column</li><li>• use less aggressive mobile phase conditions</li></ul>
Sample volume overload	<ul style="list-style-type: none"><li>• Use smaller injection volume</li></ul>
Sample solvent incompatibility with mobile phase	<ul style="list-style-type: none"><li>• Use mobile phase or weaker miscible solvent as injection solvent</li></ul>
Issues with injection valve	<ul style="list-style-type: none"><li>• Check injector valve parts</li><li>• replace worn parts</li></ul>



# Changes in Peak Shape

## Fronting

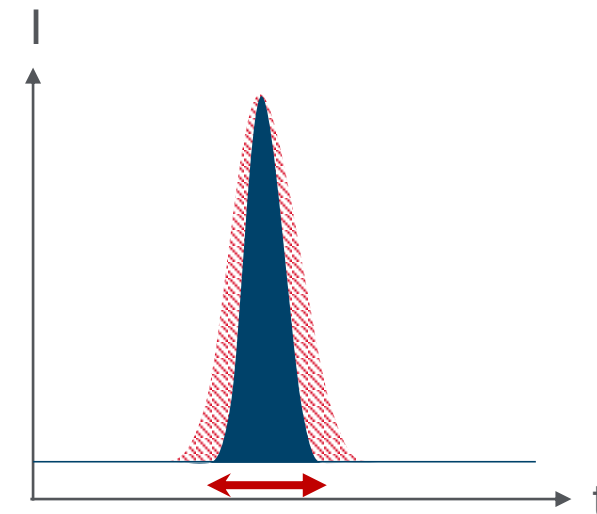
Potential Cause	Recommended Action
Channeling in column	<ul style="list-style-type: none"> <li>• Replace column</li> <li>• use guard columns</li> </ul>
Column overload	<ul style="list-style-type: none"> <li>• Use higher capacity column (increase length, diameter or change to high-capacity material)</li> <li>• decrease sample amount</li> </ul>



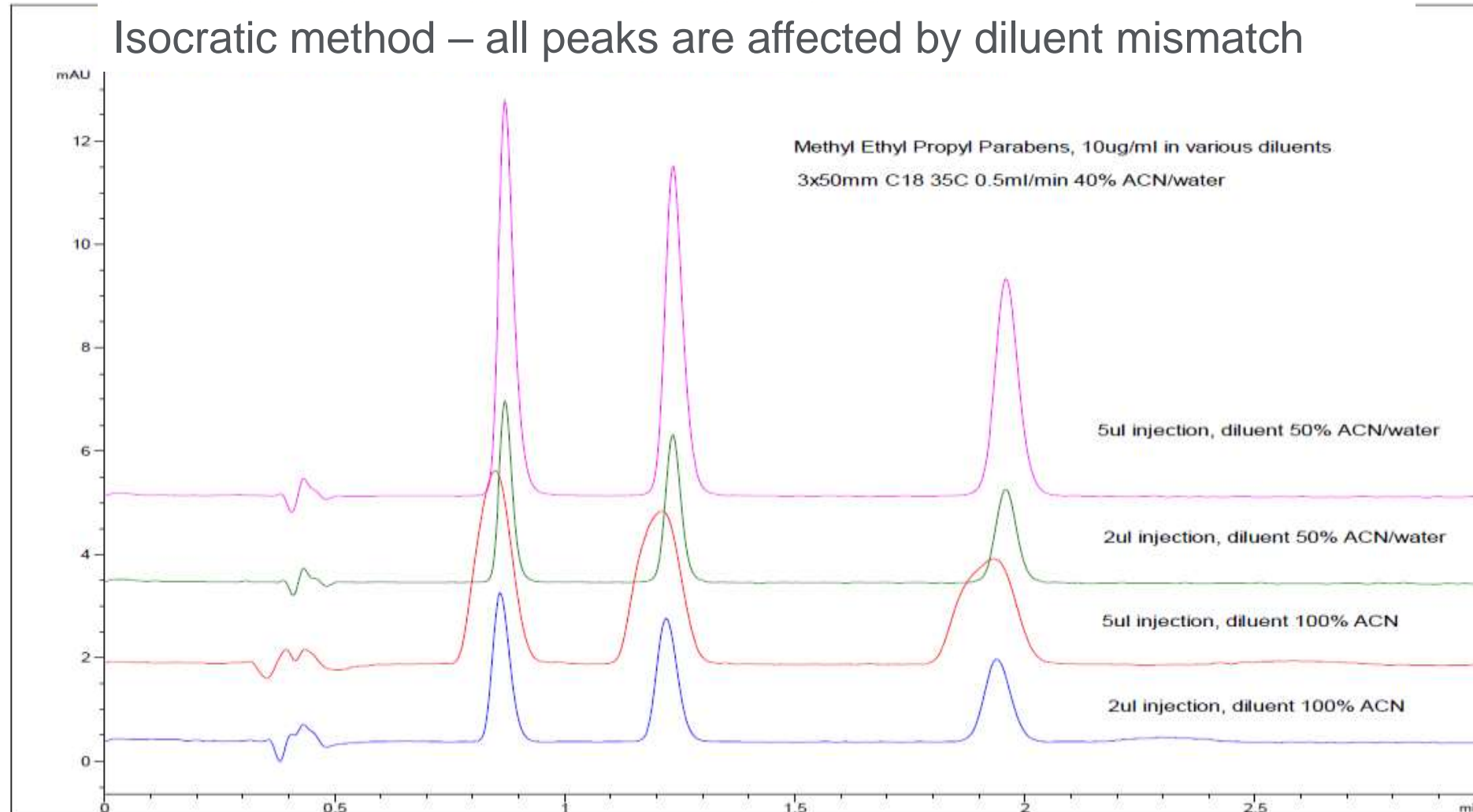
# Changes in Peak Shape

## Peak broadening

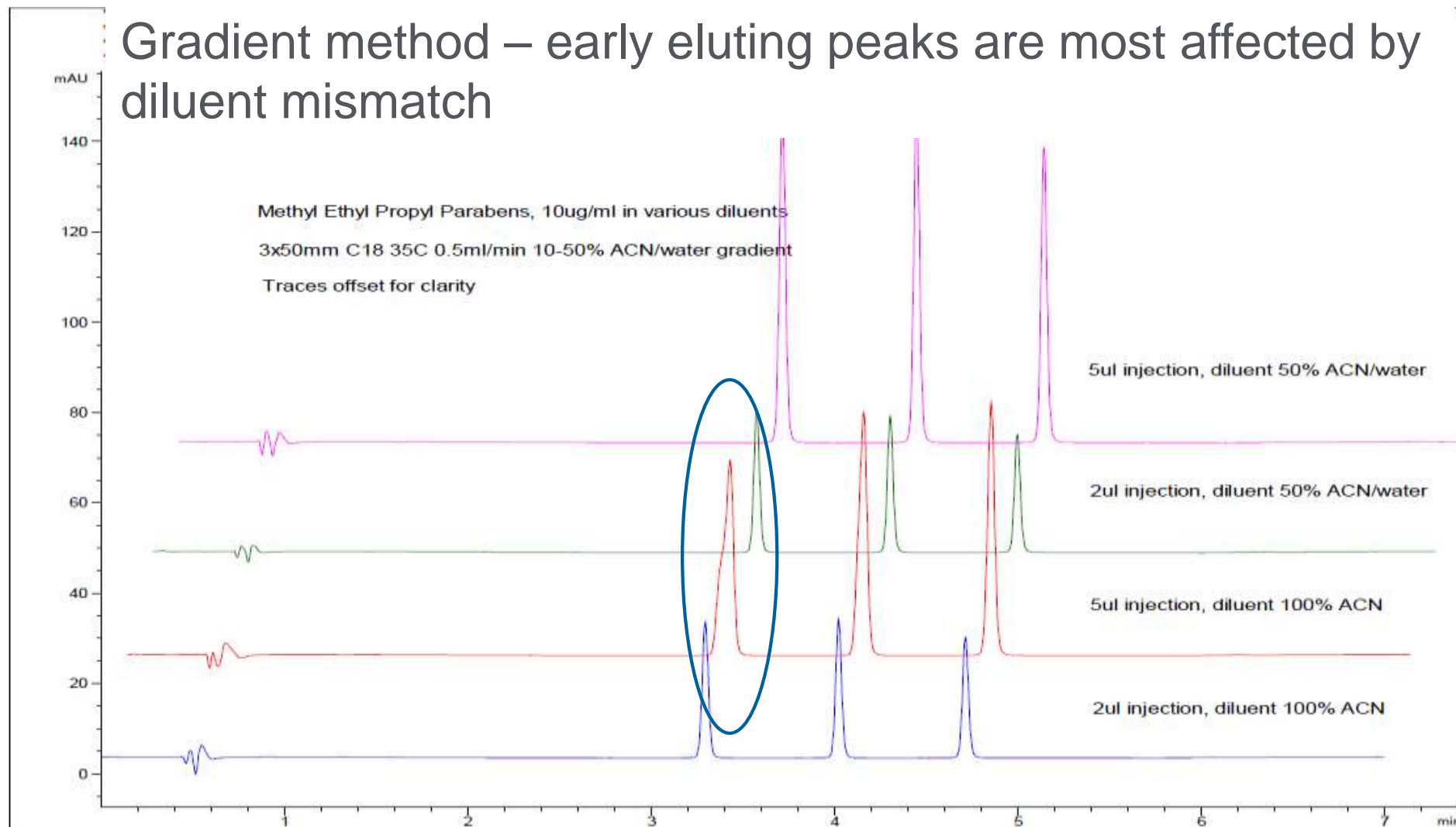
Potential Cause	Recommended Action
Injection volume too large	<ul style="list-style-type: none"> <li>Decrease injection volume</li> </ul>
Long retention times	<ul style="list-style-type: none"> <li>Use gradient elution or stronger mobile phase</li> </ul>
System settings	<ul style="list-style-type: none"> <li>Check data collection rate:</li> <li>Adjust the detector setting and / or time constant to the fastest possible value without compromising signal-to-noise.</li> </ul>
Viscosity of mobile phase too high	<ul style="list-style-type: none"> <li>Increase column temperature</li> </ul>
Detector cell volume too large	<ul style="list-style-type: none"> <li>Use smallest possible cell volume</li> </ul>
Improper fittings / connections	<ul style="list-style-type: none"> <li>Ensure that your fitting connections are made correct</li> </ul>
Extra tubing volume on system	<ul style="list-style-type: none"> <li>Ensure that the tubing is narrow and as short as possible to avoid extra volume.</li> </ul>
Sample diluent too strong	<ul style="list-style-type: none"> <li>Reduce diluent strength</li> </ul>



# Strong Diluents can Disrupt Equilibration – Isocratic Method

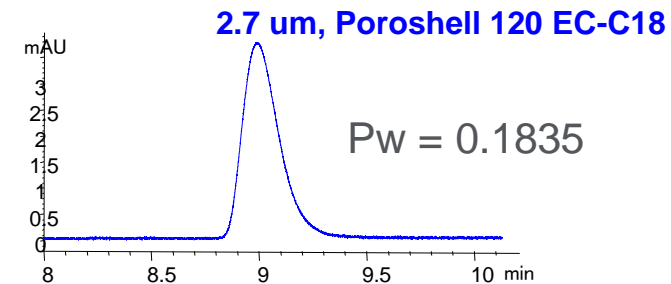
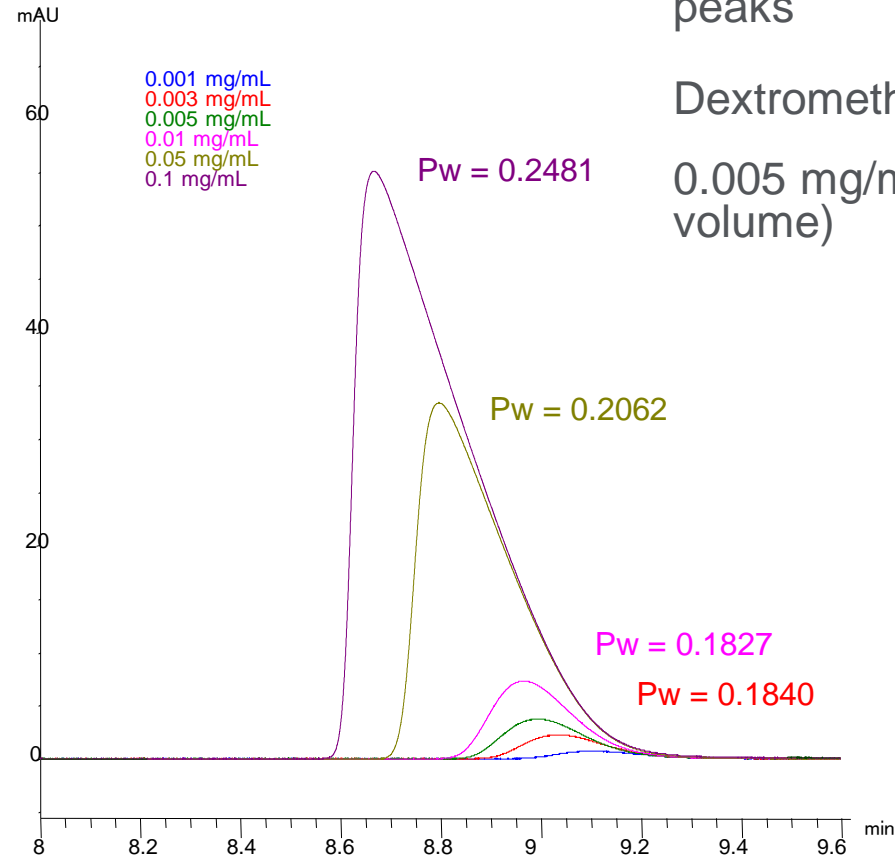


# Strong Diluents can Disrupt Equilibration – Gradient Analysis



# Comparison of Peak Shape at Low and High Loads

## Broadening and Tailing

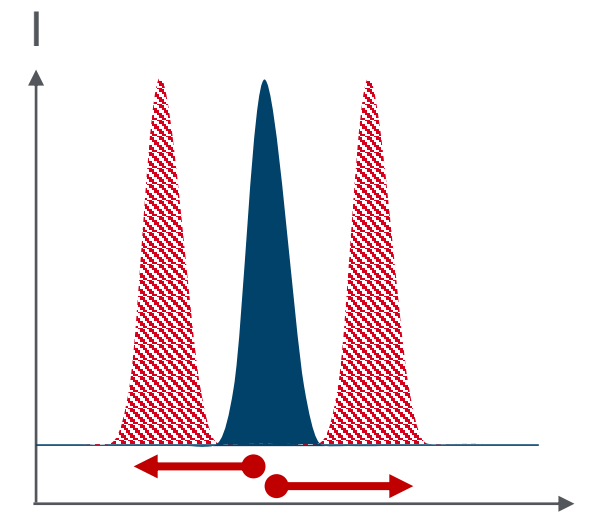


# Changes in Separation

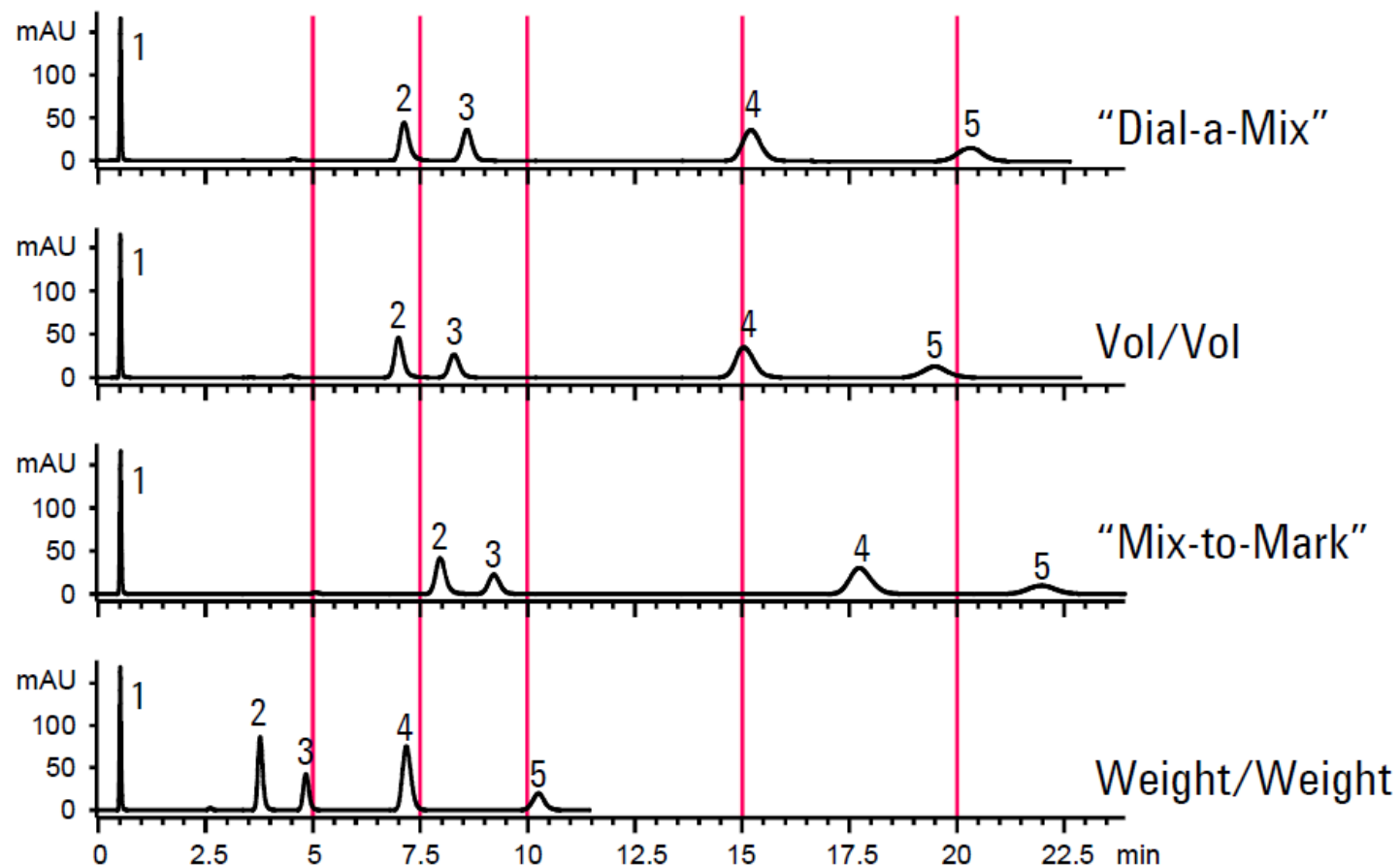
# Changes in Separation

## Retention time changing

Potential Cause	Recommended Action
Inconsistent on-line mobile phase mixing	Ensure gradient system delivering constant composition check vs. manual prep of mobile phase
Flow rate changing	Check 'Pressure fluctuation'
Column temperature varying	Thermostat column and ensure constant lab temperature
Equilibration time insufficient with gradient run or change in isocratic mobile phase	Flush with at least 10 column volumes after solvent change or gradient conclusion
Selective evaporation of mobile phase component	keep solvent reservoirs covered prepare fresh mobile phase
Buffer capacity insufficient	Use > 20 mM concentration of buffer
Contamination buildup	Occasionally flush column with strong solvent to remove contaminants
First few injections – adsorption on active sites	Condition column by initial injection of concentrated sample
Column overloaded with sample	Decrease injection volume or concentration
Mobile phase composition changing	Follow 'best practices'



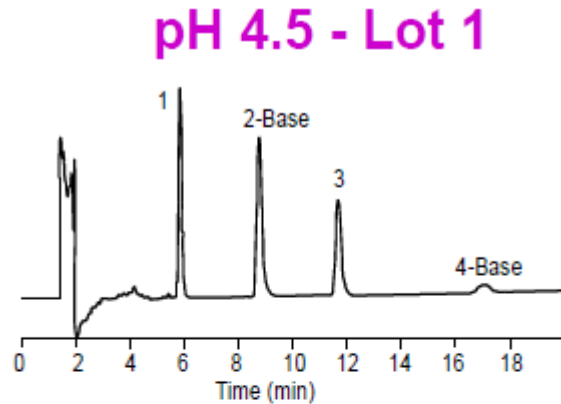
# Mobile Phase Preparation



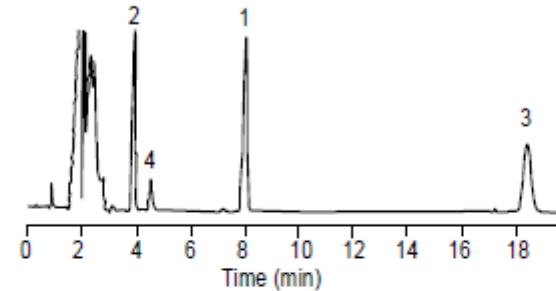
Agilent 1100 with quaternary pump  
ZORBAX Eclipse XDB-C8 Rapid-Resolution (3.5 $\mu$ m), 4.6 x 50 mm  
Agilent Part No. 935967-906  
Dial-a-Mix= A: water B: MeOH, pump 50% B  
Vol/Vol=250 mL water + 250 mL MeOH, pump 100%  
Mix-to-Mark = 250 mL MeOH, fill to 500 mL with water, pump 100%  
Premixed (w/w) = 200 g MeOH + 200 g water, pump 100%  
UV 254 nm  
1 mL/ min.

# Retention Time Shift – Selectivity differences due to incorrect pH

pH 4.5 shows selectivity change from lot-to-lot for basic compounds

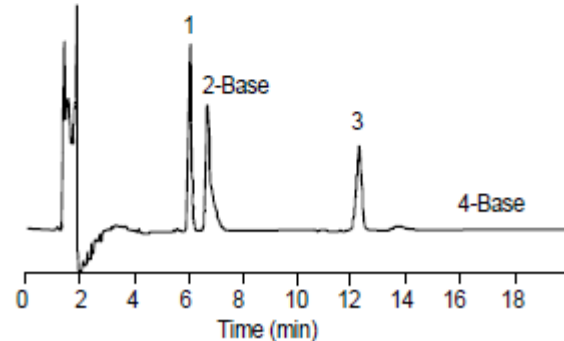


pH 3.0 - Lot 1

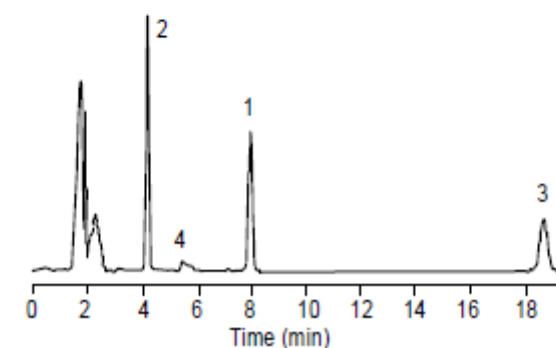


pH 3.0 shows no selectivity change from lot-to-lot

pH 4.5 - Lot 2



pH 3.0 - Lot 2



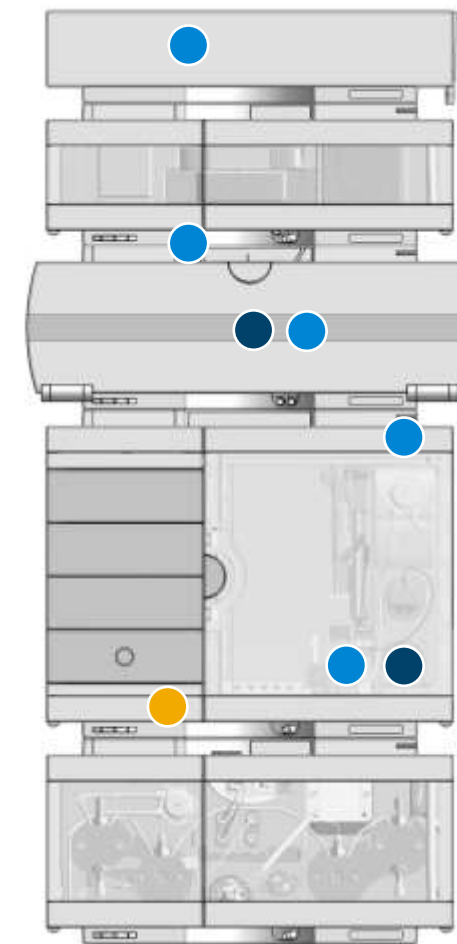
- For Method Ruggedness
  - Test 3 different column lots
  - Compare  $R_s$  for the 3 lots
    - If  $\Delta R_s$  is too large, modify method

# Changes in Separation

## Ghost peaks, carry over

	Potential Cause	Recommended Action
●	Peaks from previous injection	<ul style="list-style-type: none"><li>• Flush column to remove contaminants</li><li>• Check with blank injection</li></ul>
●	Specific interaction with metal surfaces	<ul style="list-style-type: none"><li>• Passivate instrument</li><li>• Use InfinityLab Deactivator Additive</li><li>• Use bio-inert LC equipment</li></ul>
●	Contamination or unknown interferences in samples	<ul style="list-style-type: none"><li>• Proper sample clean-up</li></ul>

**BIO  
INERT**



# Mobile Phase Hygiene

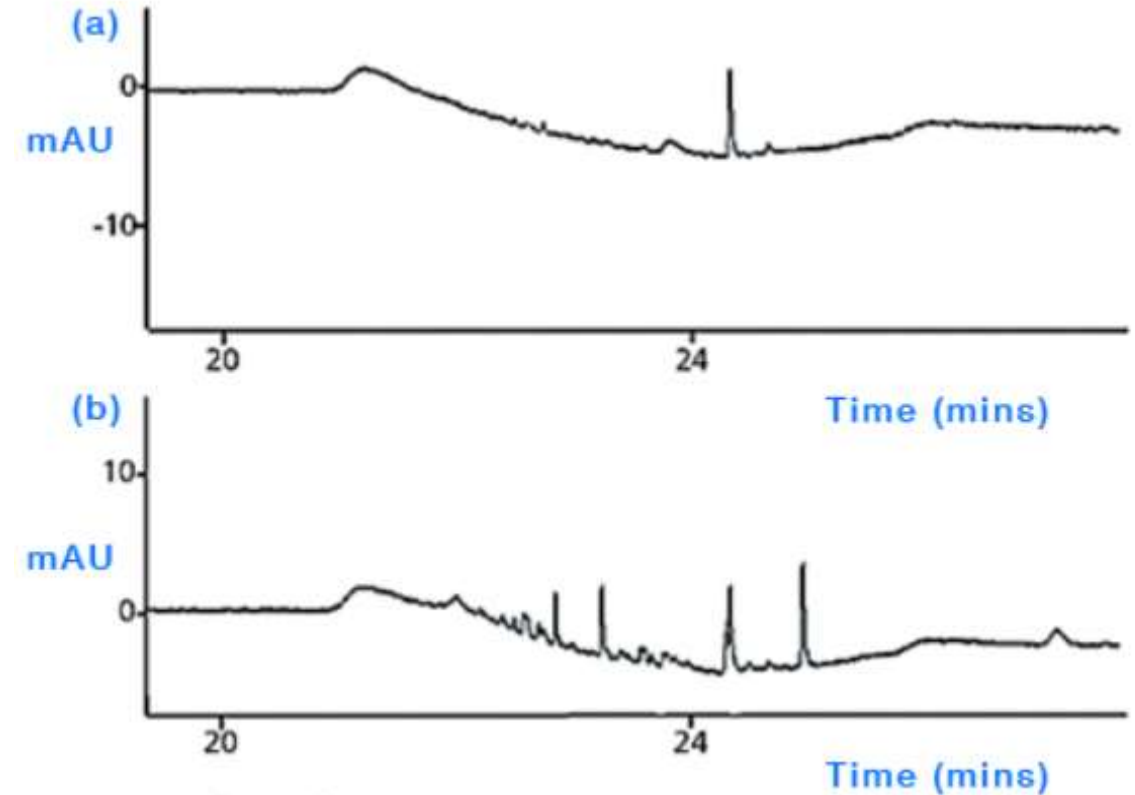
Contaminated mobile phases can cause:

- Lower sensitivity
- Rising/drifted baselines
- Higher noise
- Ghost peaks with gradient separations

Often the issue is confused with Autosampler carryover.

It can be identified by repeating the gradient run without sample injection and/or increasing the pre-run equilibration.

Always run multiple blanks before standards or samples to distinguish gradient artifacts from possible carryover.



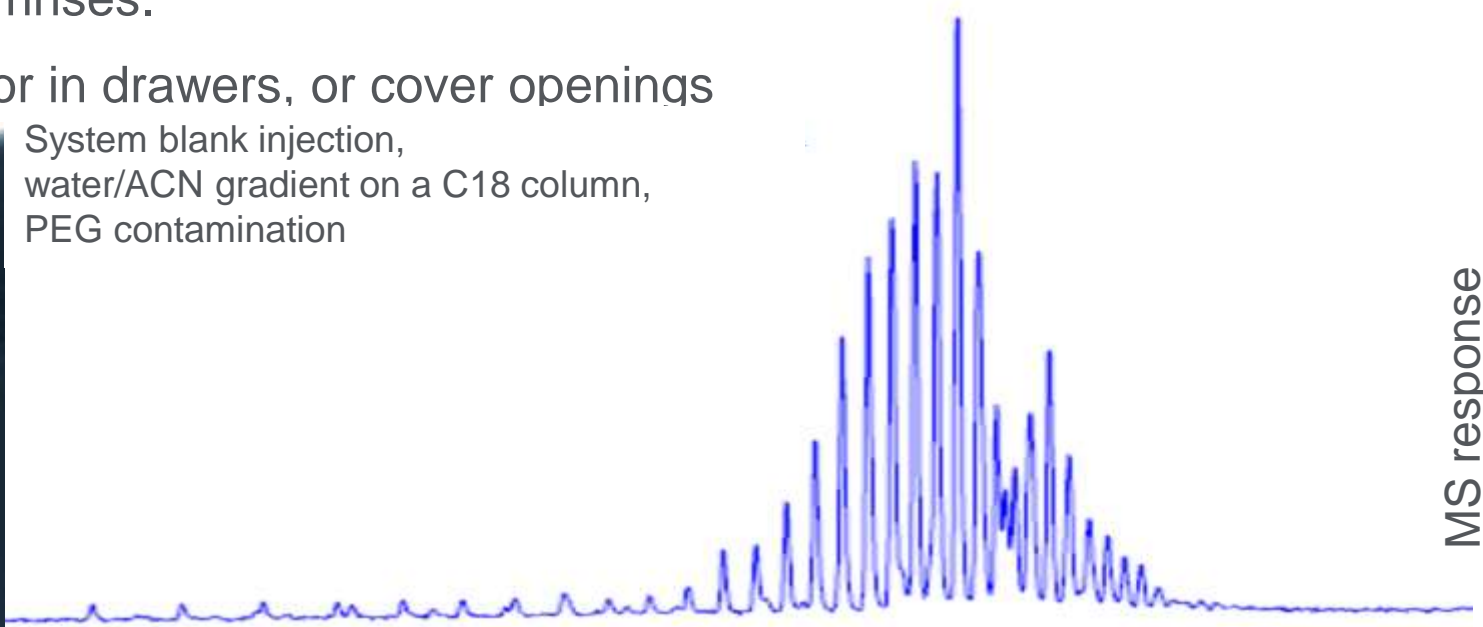
# Mobile Phase Hygiene: Glassware

## Improper cleaning of solvent bottles can cause contamination of mobile phases and result in gradient artifacts

- Wash solvent bottles with hot water, deionized water, and organic solvent (IPA or acetonitrile).
- Leave glassware inverted on paper towels on bench or on clean pegboard dowels to dry.
- **Avoid using detergents!** If it is necessary to use detergents to get glassware clean, re-wash with plenty of hot water and cold water so that all detergent residues are removed. Follow with deionized water and organic (IPA or acetonitrile) rinses.
- Store glassware inverted on shelves or in drawers, or cover openings



System blank injection,  
water/ACN gradient on a C18 column,  
PEG contamination



# Mobile Phase Hygiene: Solvent Purity and Buffer Preparation

- Use HPLC grade Organic mobile phases
- Use HPLC grade water or Milli Q DI water
- Use HPLC grade reagents including salts, ion pair reagents, and base and acid modifiers
- Always rinse pH electrode thoroughly when measuring/adjusting pH of mobile phase
- Prepare fresh buffers to avoid contaminants from the growth of bacteria or algae
- Filter your mobile phase buffer with 0.45  $\mu\text{m}$  filter before use
- Solvent filters installed at the end of solvent lines should be replaced periodically



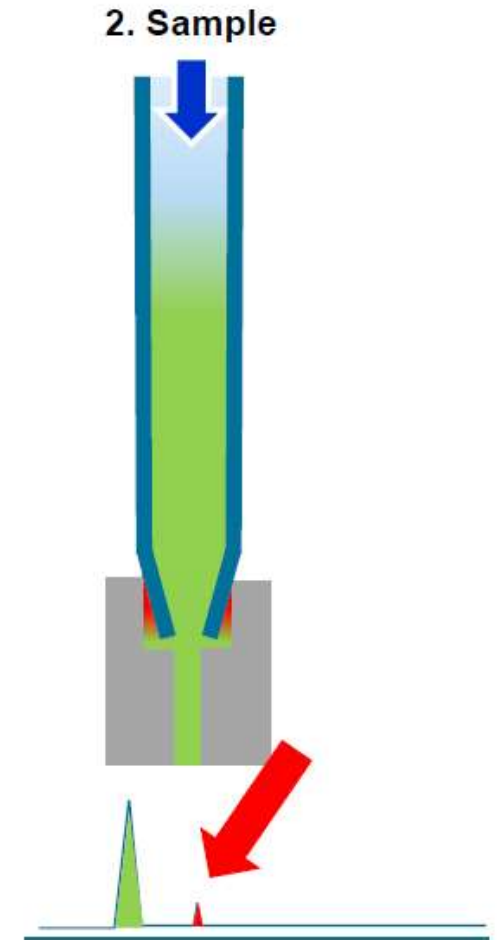
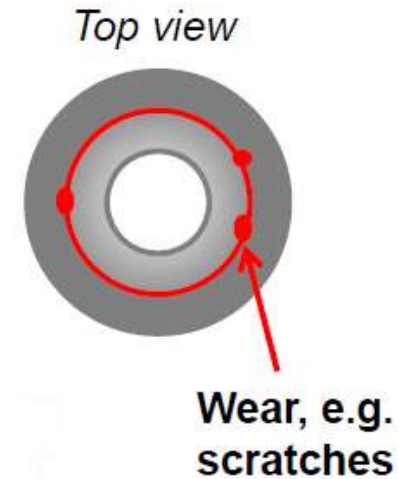
100% ACN

90%ACN+10% buffer  
(10mM phosphate)

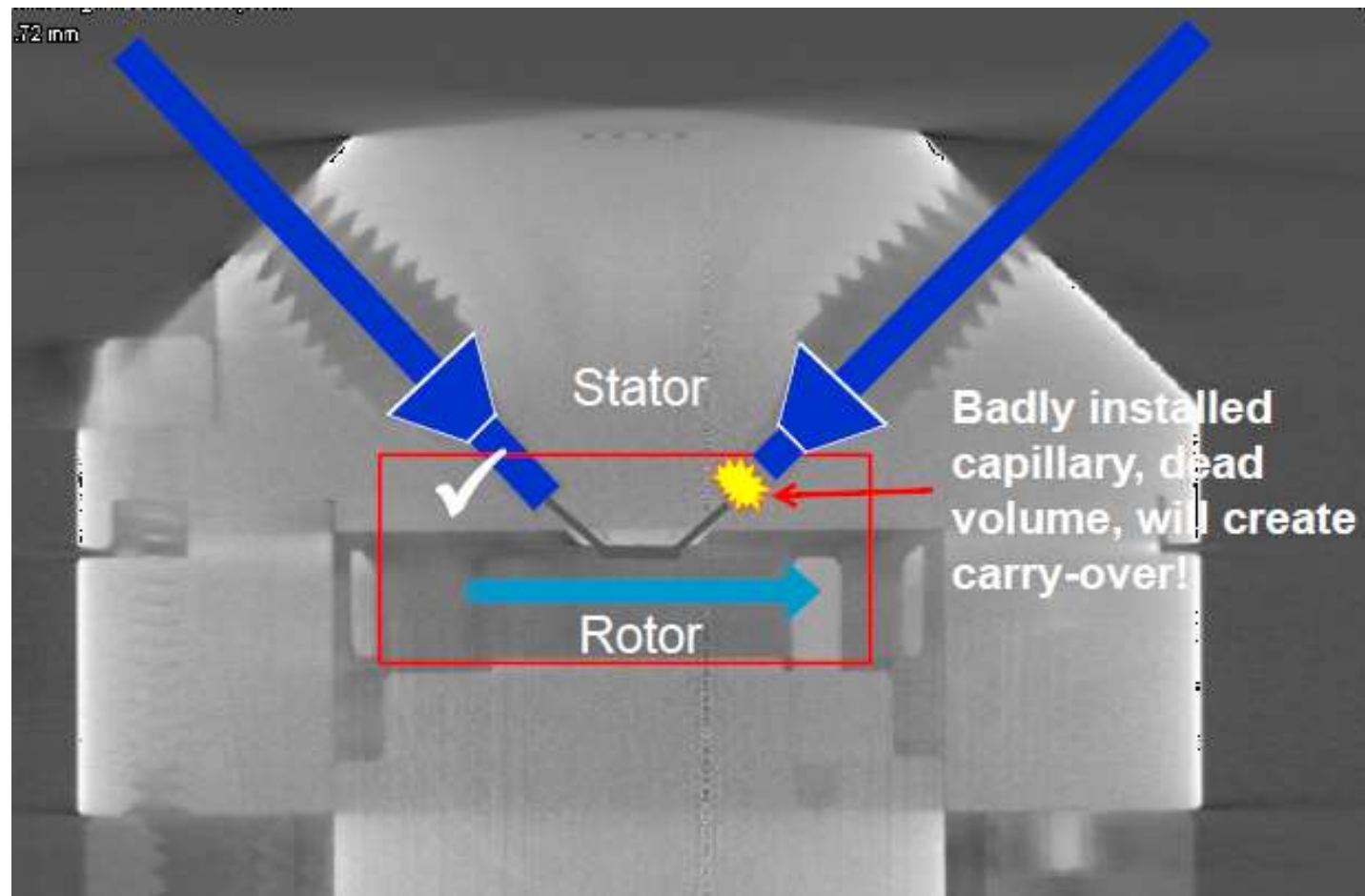
# Autosampler Carryover

## Common sources:

- Exterior of needle (use needle wash)
- Worn needle seat
- Worn rotor seal
- Poorly made fitting



# Autosampler Carryover

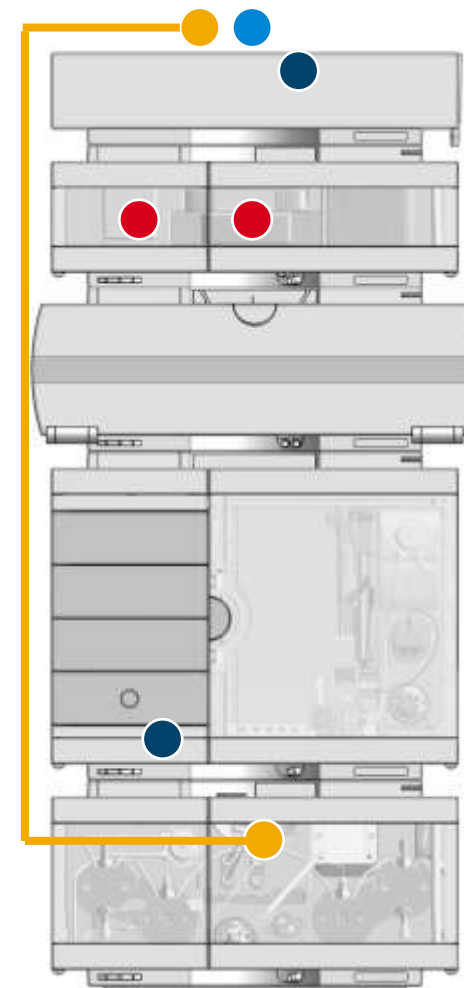


# Changes in Detection

# Changes in Detection

## Noisy baseline

	Potential Cause	Recommended Action
●	Gas bubbles in mobile phase	<ul style="list-style-type: none"><li>• Apply degassing</li><li>• check degasser performance</li></ul>
●	Low difference between sample and mobile phase absorbance	<ul style="list-style-type: none"><li>• Check absorbance values of sample vs. mobile phase</li></ul>
●	Contamination	<ul style="list-style-type: none"><li>• Use degassed HPLC-grade solvents</li><li>• flush system</li><li>• Clean up the sample</li></ul>
●	Detector optics	<ul style="list-style-type: none"><li>• Perform intensity test</li><li>• Check signal with flow cell removed if possible</li><li>• Replace lamp</li></ul>
	Pressure instability	<ul style="list-style-type: none"><li>• Check 'Pressure fluctuation'</li></ul>







# UV Lamp Tests

How do I know if my UV lamp is good?

- Visual inspection of an equilibrated baseline
- Accumulated UV lamp on-time from RFI tag or Lab Advisor
- Lab Advisor intensity test
- Lab Advisor ASTM drift and noise test
- Lab Advisor cell test

Filter

☒ All Counters ☐ Counters with Limit

	Title	Value	Unit	Limit	Progress		
 <b>G4220A</b> Serial # DEBAA00157	<b>1290 Bin Pump</b>	0	Hour	3000	0%	★ ↺	
		3.45	Liter	50	6%	★ ↺	
		8.64	Liter	50	17%	★ ↺	
		631	Count	15000	4%	★ ↺	
	Liquimeter (A+B)	12.09	Liter	0	0%	★ ↺	
 <b>G4226A</b> Serial # DE93000560	<b>1290 ALS</b>	0	Count	1000	0%	★ ↺	
		Needle into seat counter	1191	Count	1500	79%	★ ↺
			1.53	Hour	3000	0%	★ ↺
	Valve switching counter	2418	Count	60000	4%	★ ↺	
 <b>G4212A</b> Serial # DEBAF00163	<b>1290 DAD</b>	Accumulated UV lamp on-time	2519.65	Hour	2000	100%	★ ↺
		UV lamp ignition counter	28	Count	1500	1%	★ ↺
		UV lamp on-time	360.65	Hour	0	0%	★ ↺
 <b>G4208A</b> Serial # PP55055002	<b>1200 Instant Pilot</b>						

# UV Lamp Tests

Diode array and  
multiple wavelength

Counters and hours

		Title	Value	Unit	Limit	Progress	
		G7117B					
	<b>G7117B</b>	<b>1290 DAD</b>	Accumulated UV Lamp On-Time	3.17	h	<input type="text" value="0"/>	0%
	Serial #	PPBAW00058	Number of UV Lamp Ignitions	2	Count	<input type="text" value="0"/>	0%

The useable lifetime of a deuterium lamp will depend on it's use:

- How many hours has it been on?
- How many times has it been ignited?
- What wavelength is being used?

# UV Lamp Tests

Diode array and  
multiple wavelength

Intensity test

Intensity Test

1260 HPLC » 1260 HPLC » G7115A:DEAC600377

GeneralLimitsSignals

Test Name

Intensity Test

Description

The test scans the Intensity spectrum generated by the UV and VIS Lamp.

Module

G7115A:DEAC600377 (1260 DAD WR)

Status

Passed

Start Time

2/21/2019 3:21:31 PM

Stop Time

2/21/2019 3:22:15 PM

Test Procedure

✓ 1. Check Prerequisites...

✓ 2. Remove Flow Cell.

✓ 3. Scan Intensity Spectrum...

✓ 4. Evaluate Data...

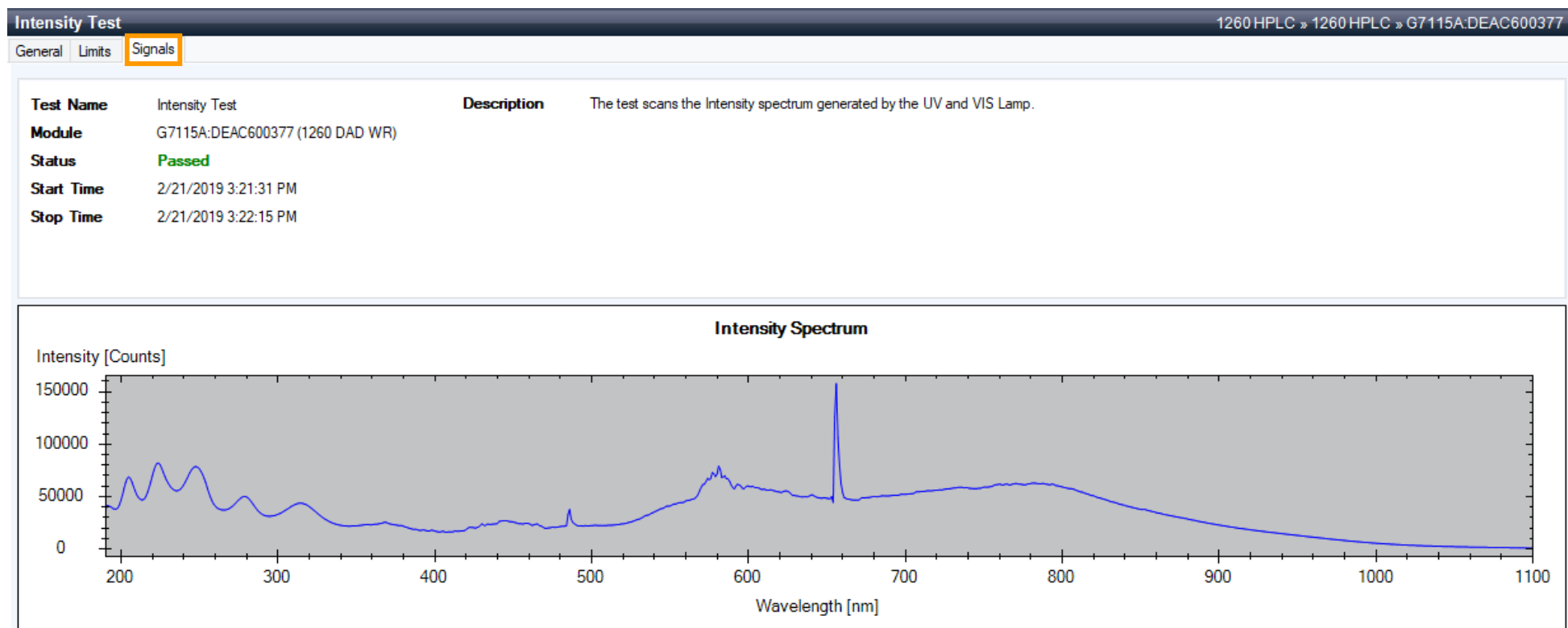
Result

Name	Value
Accumulated UV Lamp Burn Time	126.35 h
UV Lamp On-Time	0.02 h
Accumulated Vis Lamp Burn Time	262.45 h
Vis Lamp On-Time	0.02 h
Lowest Intensity in Range 190 - 220 nm	36289 Counts
Lowest Intensity in Range 190 - 220 nm	2000 Counts
Lowest Intensity in Range 221 - 350 nm	21963 Counts
Lowest Intensity in Range 221 - 350 nm	5000 Counts
Lowest Intensity in Range 351 - 500 nm	16150 Counts
Lowest Intensity in Range 351 - 500 nm	2000 Counts
Lowest Intensity in Range 501 - 950 nm	13102 Counts
Lowest Intensity in Range 501 - 950 nm	2000 Counts
Highest Intensity in Range 190 - 350 nm	81934 Counts
Highest Intensity in Range 190 - 350 nm	2000 Counts
Highest Intensity in Range 700 - 950 nm	62919 Counts
Highest Intensity in Range 700 - 950 nm	2000 Counts
Highest Intensity for D2 Alpha Line (600 - 700 nm)	157676 Counts
Highest Intensity for D2 Alpha Line	2000 Counts
Spectrum Integral	31863112
UV Integral (190 - 349 nm)	7384053

# UV Lamp Tests

Diode array and  
multiple wavelength

Intensity test



The profile of the intensity scan changes as a lamp ages

# UV Lamp Tests

Diode array and  
multiple wavelength

ASTM drift and noise

ASTM Drift and Noise Test

1260 HPLC » 1260 HPLC » G7115A:DEAC600377

GeneralLimitsSignals

Test Name

ASTM Drift and Noise Test

Description

The test performs ASTM drift and noise evaluation without reference.

Module

G7115A:DEAC600377 (1260 DAD WR)

Status

Passed

Start Time

2/21/2019 4:56:26 PM

Stop Time

2/21/2019 5:16:47 PM

Test Procedure

✓ 1. Check Prerequisites...

✓ 2. Remove Flow Cell.


✓ 3. Measure Noise...

✓ 4. Evaluate Data...

Result

Name	Value
Accumulated UV Lamp Burn Time	127.94 h
UV Lamp On-Time	1.60 h
Minimum Lamp On-Time	1 h
Accumulated Vis Lamp Burn Time	264.03 h
Vis Lamp On-Time	1.61 h
Minimum Lamp On-Time	1 h
Signal Drift value at 254 nm (UV)	-0.138 mAU/h
Maximum Allowed Drift	-1 ... 1 mAU/h
Signal Noise value at 254 nm (UV)	0.008 mAU
Maximum Allowed Noise	0 ... 0.02 mAU
Signal Drift value at 750 nm (Vis)	-0.415 mAU/h
Maximum Allowed Drift	-1 ... 1 mAU/h
Signal Noise value at 750 nm (Vis)	0.013 mAU
Maximum Allowed Noise	0 ... 0.02 mAU

ASTM Drift and Noise Test



Remove the flow cell from the lightpath of the detector.

OK

Cancel

# UV Lamp Tests

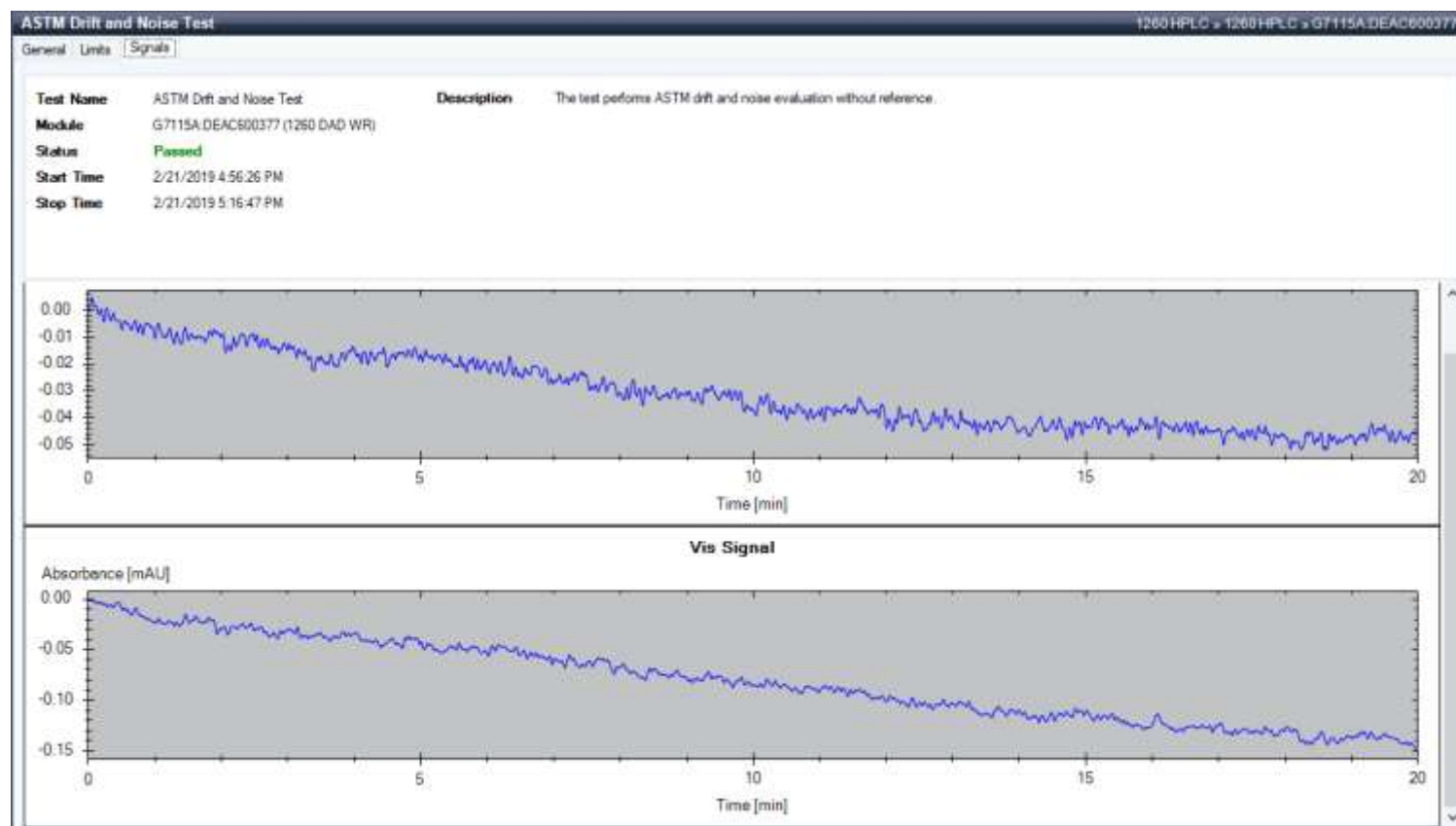
Diode array and  
multiple wavelength

ASTM drift and noise

Run on a monthly basis this test  
can help track the natural  
decline of the lamp and perhaps  
raise awareness of a dirty cell.

Name	Value	Description
Minimum Lamp On-Time	1 h	The minimum lamp on-time to perform a noise check.

Name	Lower limit	Upper limit	Description
Maximum Allowed Noise	0 mAU	0.02 mAU	The maximum allowed Signal noise in mAU.
Maximum Allowed Drift	-1 mAU/h	1 mAU/h	The maximum allowed Signal drift in mAU.



# UV Lamp Tests

Diode array and  
multiple wavelength

Cell test

Cell Test1260 HPLC » 1260 HPLC » G7115A:DEAC600377

GeneralLimitsSignals

Test Name

Cell Test

Description

The test compares the lamp intensity with and without the flow cell installed. The intensity ratio is an indicator of the amount of light absorbed by the flow cell.

Module

G7115A:DEAC600377 (1260 DAD WR)

Status

Passed

Start Time

2/21/2019 3:57:24 PM

Stop Time

2/21/2019 3:59:37 PM

Test Procedure

✓ 1. Check Prerequisites...

✓ 2. Remove Flow Cell.

✓ 3. Scan Intensity Spectrum...

✓ 4. Insert Flow Cell.

✓ 5. Scan Intensity Spectrum...

✓ 6. Evaluate Data...

Result

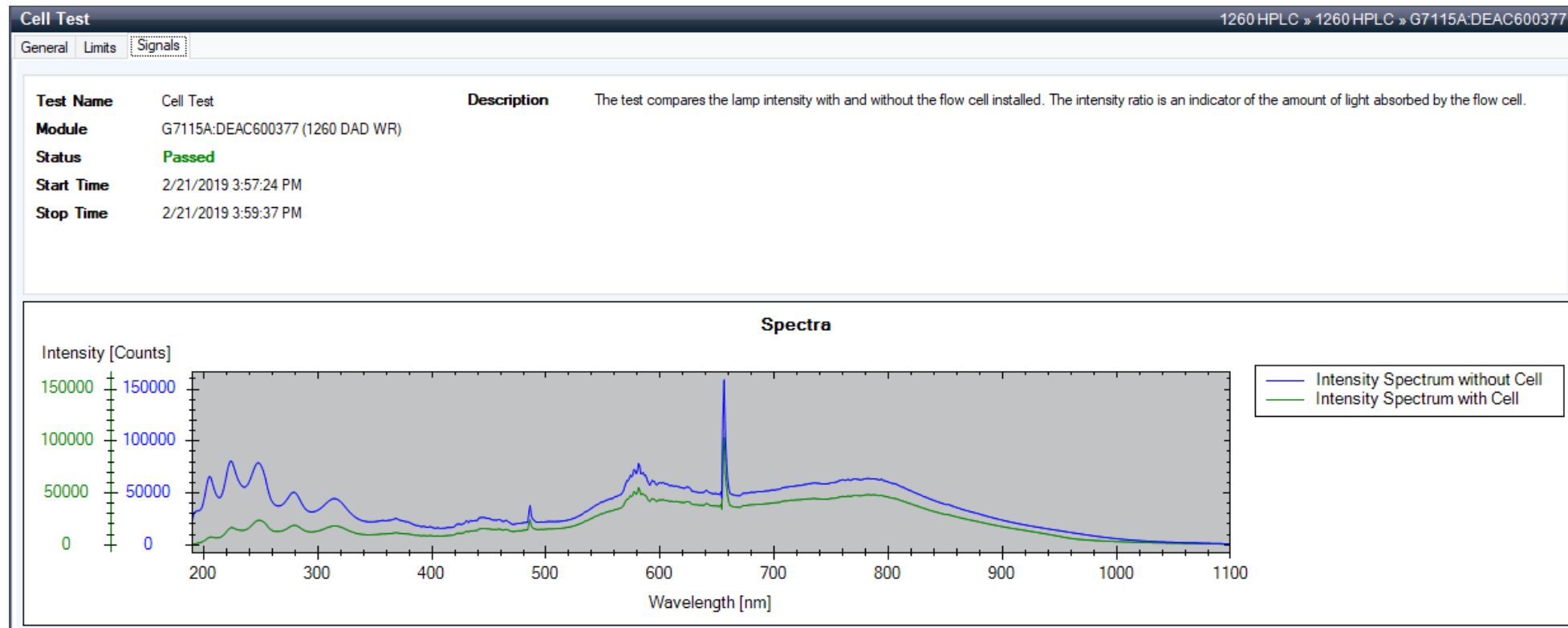
Name	Value
Accumulated UV Lamp Burn Time	126.95 h
UV Lamp On-Time	0.62 h
Accumulated Vis Lamp Burn Time	263.05 h
Vis Lamp On-Time	0.62 h
Intensity Integral without Flow Cell	32,088,720
Intensity Integral with Flow Cell	19,830,098
Intensity Ratio	0.62
Minimum Intensity Ratio	0.3

Diode array detectors with the fiberoptic style flow cell require a Max Light test cell for this test - part number G4212-60011.

# UV Lamp Tests

Diode array and  
multiple wavelength

Cell test

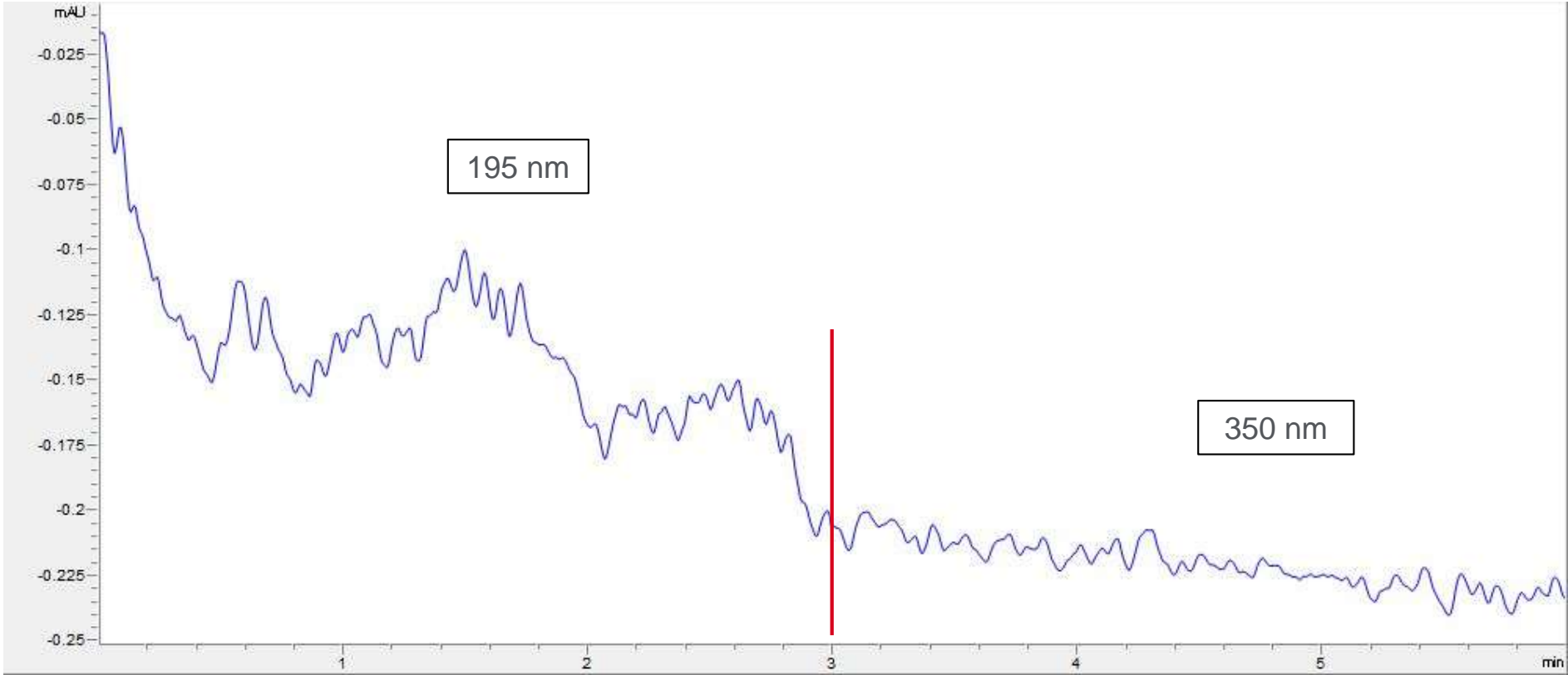


Example of scans with and without cell installed.

# How do I know if my UV lamp is good?

Diode array and  
multiple wavelength

Baseline inspection



A number of factors contribute to the specific amplitude and pattern of baseline noise and drift, including the specific wavelength, mobile phase, room temperature and data rate.

# How do I know if my UV lamp is good?

Diode array and  
multiple wavelength

## Long cycle wave

This is a rhythmic change in the baseline where the periodicity may be hours.

- environmental influences

## Baseline inspection

## Short cycle wave

This is a rhythmic change in the baseline where the periodicity may be seconds or minutes.

- solvent mixing noise
- mechanical issue in pump

If the cycle of the wave does not appear to be mixing noise, evaluate the health of the lamps through Lab Advisor intensity and noise and drift tests. Also, the cleanliness of the flow cell through the cell test.

## Excessive drift

In a UV baseline light scattering shows up as drift. If the baseline is drifting more than expected, empty and rinse the solvent bottles, refilling with fresh solvent. Perform a cell test to check the cleanliness of the flow cell.

# Useful parts

Parts that address potential issues and help to ease your daily tasks

Part Description	Information	Part number
InfinityLab Stay Safe caps	Prevents solvent evaporation; changes in mobile phase concentration	Various <a href="http://www.agilent.com/chem/staysafecaps">www.agilent.com/chem/staysafecaps</a>
InfinityLab Quick Connect and Quick Turn fittings	With spring-load function for optimized dead volume reduction	Various <a href="http://www.agilent.com/chem/infinitylabfittings">www.agilent.com/chem/infinitylabfittings</a>
Blank nut, long, 10-32	Blank nut, PEEK with steel core; for system diagnostic tests; finger tight up to 1300 bar, easy to use and gentle to receiving port	5043-0277
Agilent Captiva syringe filters	Solve issues like inlet clogging, increased backpressure, and retention time shift by filtering your samples	Various <a href="http://www.agilent.com/chem/filtration">www.agilent.com/chem/filtration</a>
InfinityLab Poroshell 120 columns	High efficiency and high resolution; available in 18 chemistries	Various <a href="http://www.agilent.com/chem/discoverporoshell">www.agilent.com/chem/discoverporoshell</a>



InfinityLab Poroshell 120 columns



InfinityLab Stay Safe cap on solvent bottle



InfinityLab Quick Connect fitting



InfinityLab Quick Turn fitting



Blank nut, 5043-0277

# LC Troubleshooting Poster Available

## LC Troubleshooting Guide

Your guide to solving common problems and staying productive

Agilent  
InfinityLab

### Places to Start

#### Solvents

- Use brown borosilicate bottles to avoid algae growth
- Prepare solvent volume to be used up within 1 to 2 days
- Use only HPLC-grade solvents filtered through 0.2 µm filters

#### Preparing and powering up the pump

- Inspect solvent bottles and inlet filters for damage or clogging
- Always use seal wash when installed and purge the pump
- Use the appropriate system conditioning method

#### Daily tasks

- Replace aqueous and organic mobile phases every second day
- Check seal wash solvent
- Flush the system with the composition of your application

#### Weekly tasks

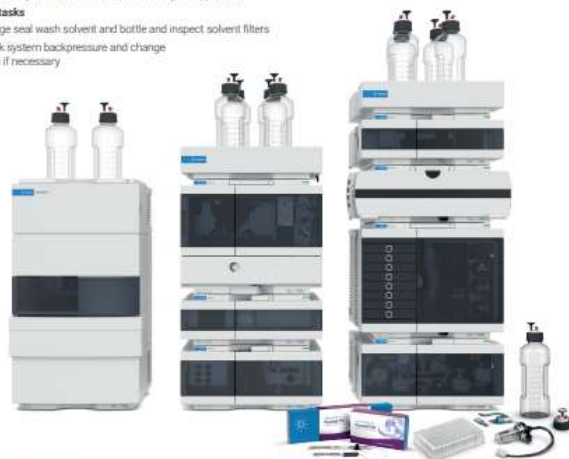
- Change seal wash solvent and bottle and inspect solvent filters
- Check system backpressure and change filters if necessary

#### Pump shutdown

- Flush all channels to remove salt deposits and particulate matter
- Flush the system with appropriate storage solvent and power down the system

#### Handling of acetonitrile

- If possible, use 5 to 10% of water in your mobile phase
- Be sure to avoid ACN evaporation
- Don't leave ACN on the system for more than 2 to 3 days
- Perform a periodic warm water wash (50 to 70 °C) if you face problems



### Maintenance

Agilent Lab Advisor software helps you manage your Agilent LC instruments to achieve high-quality chromatographic results in the most efficient way by ensuring high instrument performance, productivity, and reliability. It is available free-of-charge.

- Diagnostic tests to evaluate performance
- Easier maintenance of all Agilent LC modules
- Comprehensive reports generated to ease communication with Agilent service

#### Retention Time Drift



Possible Cause	Solution
Inconsistent online mobile phase mixing	Ensure gradient system delivers constant composition; compare with manual preparation of mobile phase
Variation in column temperature	Thermostat or insulate column; ensure constant lab temperature
Insufficient equilibration time with gradient run or change in isocratic mobile phase	Make sure at least 10 column volumes pass through column after sample run
Selective evaporation of mobile phase component	Less volatile helium sparging; keep solvent reservoirs covered; prepare fresh mobile phase
Contamination buildup	Occasionally flush column with strong solvent
Column overloaded with sample	Decrease injection volume or concentration

#### Pressure Fluctuation



Possible Cause	Solution
Leak in the system	Identify the channel and clean or replace check valves; replace pump seals
Buildup of particulates	Filter sample and mobile phase
Bubbles in pump	Perform solvent degassing; sparge solvent with helium

#### Pressure Increase



Possible Cause	Solution
System blockage	Check flowpath (needle seat, capillaries, filter and filter)
Water/organic systems: buffer precipitation	Test buffer/organic mixtures to ensure compatibility

#### High Column Backpressure



Possible Cause	Solution
Column blockage	Bottle sample cleanup; use guard column
Mobile phase viscosity too high	Use lower viscosity solvents or higher temperature
Particle size too small	Use larger d <sub>p</sub> packing
Plugged inlet frit	Replace column

#### Drifting Baseline



Possible Cause	Solution
Positive/negative direction: contaminant buildup/cleanup	Flush column; clean up sample; use pure solvents
Positive/negative: difference in refractive index of injection solvent	Use mobile phase for sample solvent
Temperature changes	Insulate and thermostat column and tubing

#### Noisy Baseline



Possible Cause	Solution
Contamination	Use degassed HPLC-grade solvents; flush system; clean up sample
Detector problems	Check number of hours of UV lamp; replace UV lamp or flow cell

#### Ghost Peaks



Possible Cause	Solution
Peaks from previous injection	Flush column to remove contaminants; check with blank injection
Contaminant; unknown interferences in samples	Proper sample cleanup
Ion pair: diazepam	Prepare sample in actual mobile phase to minimize disturbance
Contaminated mobile phase	Check your mobile phase
Bubbles in solvent	Check and degas your solvents

#### Peak Tailing



Possible Cause	Solution
Unwashed dead volumes	Minimize number of connections; ensure injector seal is tight; ensure fittings are properly seated
Column performance	Change mobile phase; replace column
Silica-based: column degradation	Use specialty, polymeric, or sterically protected column
Silica-based: basic interactions with stationary phase	Use stronger mobile phase or add appropriate base (e.g., TEA)

#### Peak Broadening



Possible Cause	Solution
Injection volume too large	Decrease injection volume or solvent strength of injection solvent; use gradient methods
Low sampling rate of data system	Increase data rate
Detector cell volume too large	Use smallest possible cell volume
Injection volume too large	Decrease injection volume

#### Sensitivity Problems



Possible Cause	Solution
Peaks are outside of sensitivity range of detector	Dilute/concentrate sample to bring into linear region
Sample-related losses during preparation	Use internal standard during sample preparation; optimize sample preparation method

#### Leaks



Possible Cause	Solution
White powder at fittings; loose fitting	Tighten fittings; replace capillaries
System leak	Identify location checking leak sensors/sensors; check flow cell

Discover more best practices for using an Agilent LC system:  
<https://www.agilent.com/chem/lc-best-practices>



Training courses are available at:  
<https://www.agilent.com/crosslab/university>



Get answers. Share insights. Join the Agilent Community at:  
<https://community.agilent.com>



For Lab Advisor software, please visit:  
<https://www.agilent.com/chem/lab-advisor>



Request yours today at  
[www.agilent.com/chem/troubleshootLC](https://www.agilent.com/chem/troubleshootLC)

# Resources for Support

- *New!* LC Troubleshooting poster: 5994-0709EN
- Resource page: <http://www.agilent.com/chem/agilentresources>
  - Quick reference guides
  - Catalogs, column user guides
  - Online selection tools, how-to videos
- InfinityLab Supplies catalog: [5991-8031EN](#)
- LC handbook: [5990-7595EN](#)
- YouTube – [Agilent channel](#) (maintenance videos)
- Agilent service contracts



# Contact Agilent Chemistries and Supplies Technical Support



1-800-227-9770 Option 3, Option 3:

Option 1 for GC and GC/MS columns and supplies

Option 2 for LC and LC/MS columns and supplies

Option 3 for sample preparation, filtration and QuEChERS

Option 4 for spectroscopy supplies

Option 5 for chemical standards

**Available in the USA & Canada 8-5 all time zones**

[gc-column-support@agilent.com](mailto:gc-column-support@agilent.com)

[lc-column-support@agilent.com](mailto:lc-column-support@agilent.com)

[spp-support@agilent.com](mailto:spp-support@agilent.com)

[spectro-supplies-support@agilent.com](mailto:spectro-supplies-support@agilent.com)

[chem-standards-support@agilent.com](mailto:chem-standards-support@agilent.com)