

Advanced Troubleshooting

Daron Decker / Eric Pavlich
Chromatography Technical Specialist



Troubleshooting Tips

1. Isolate the problem.

(Blank Runs, Inject Un-retained Compound, Know what it is not)

2. Change only one variable at a time.

3. Compare before/after chromatograms.

(Peak shape, response, retention, baseline rise, background, look for trends, etc.)



What Can Possibly Go Wrong?

INJECTOR – contamination, flow settings, flow path issues, overload, valve settings, faulty consumables

COLUMN – contamination, flow settings, flow path issues, damage (activity & bleed), breakage

DETECTOR – contamination, flow settings, flow path issues, electronics

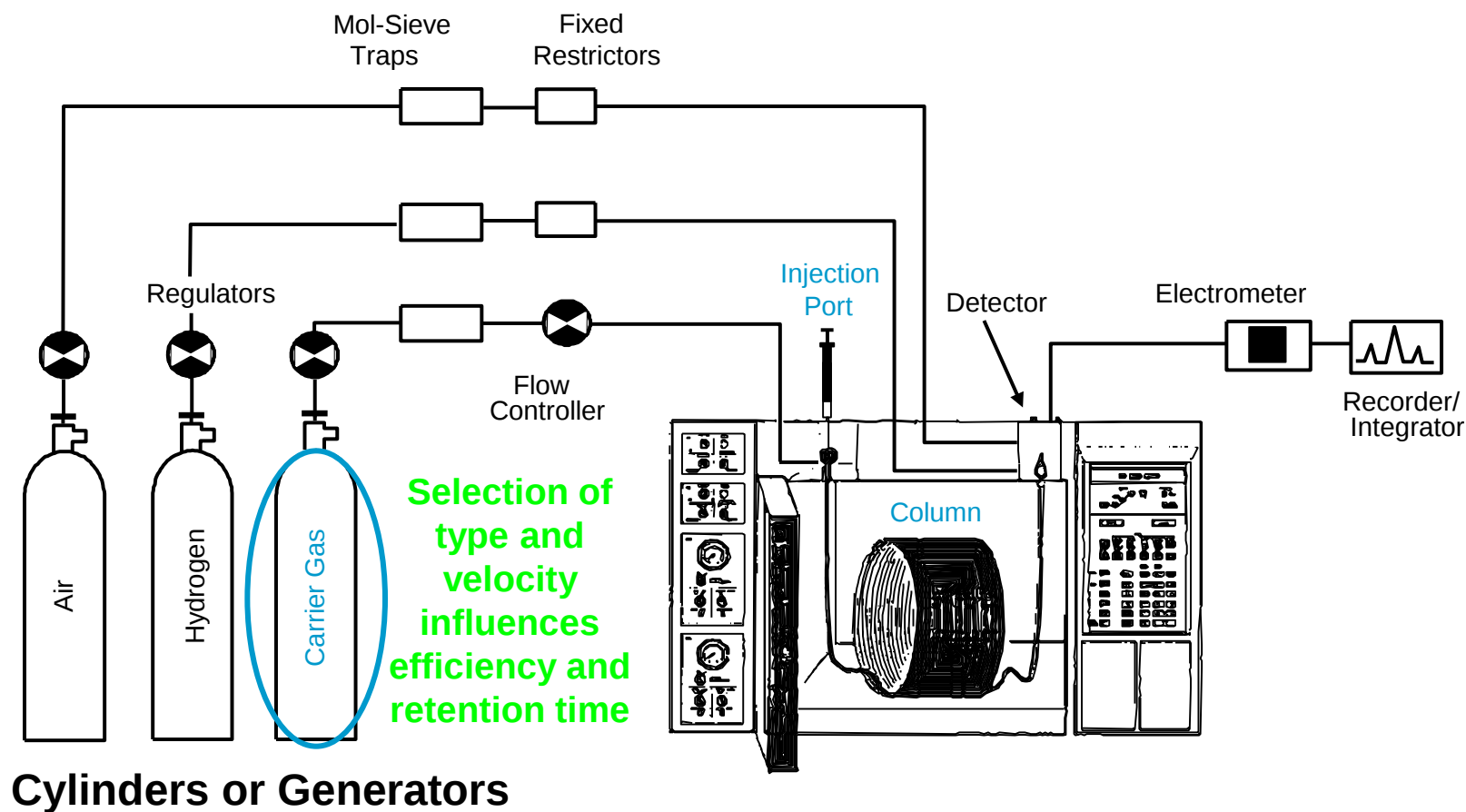


Logical Troubleshooting

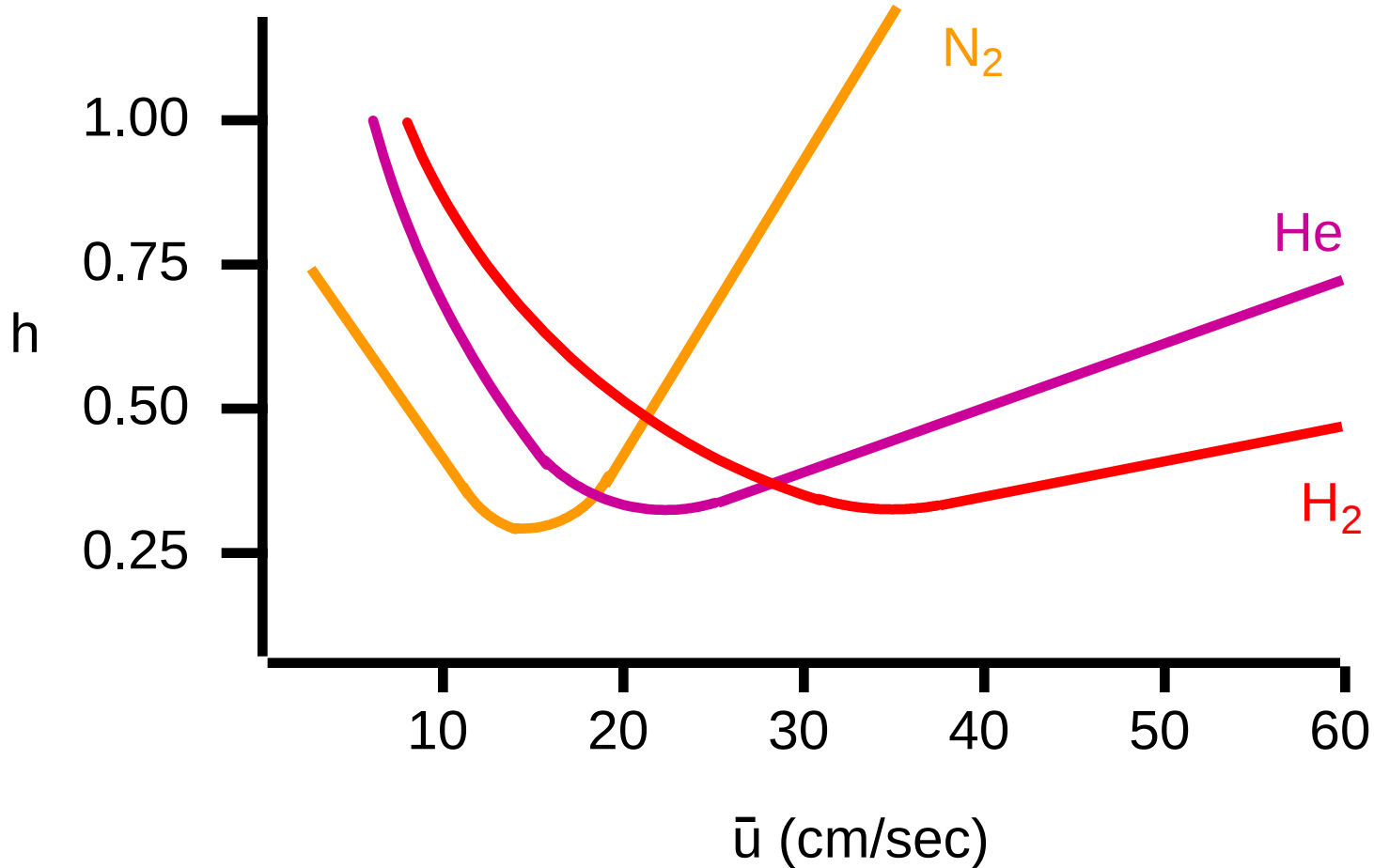
- **Troubleshooting Starts with Isolating the problem**
 - There are 5 basic areas from where the problem arises
 - FLOW
 - INJECTOR
 - COLUMN
 - DETECTOR
 - ELECTRONICS
(Temperature)
 - But of course it can always be some COMBINATION
- **Knowing what can & can't cause the symptom is the key**



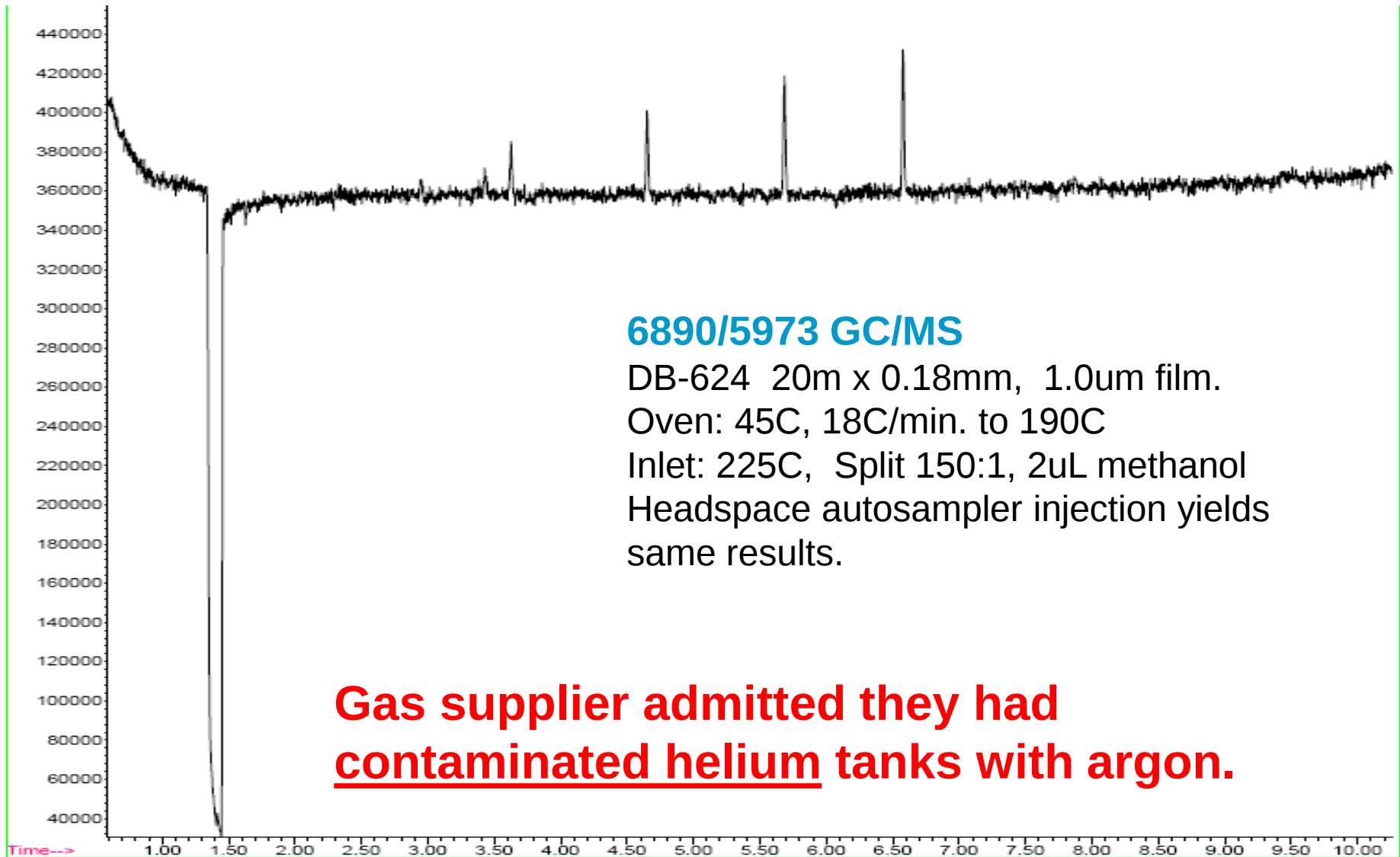
Typical Gas Chromatographic System



van Deemter Curves



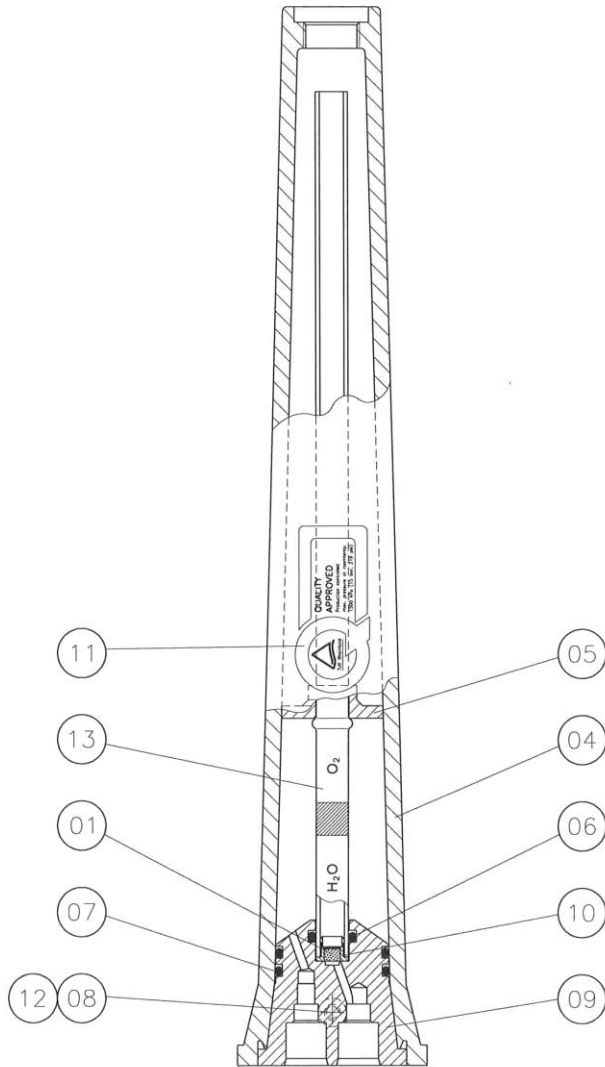
Noisy Baseline Even at Low Temperatures



Gas Clean Filters



Gas-Clean Filters Design



Sensitive indicators for moisture and oxygen

Short filter equilibration time

Fast, leak-free filter replacement

- without turning off the gas
- NO tools needed

Oxygen Damages GC Columns

Repair of the GC column is impossible

Replacements means re-calibration

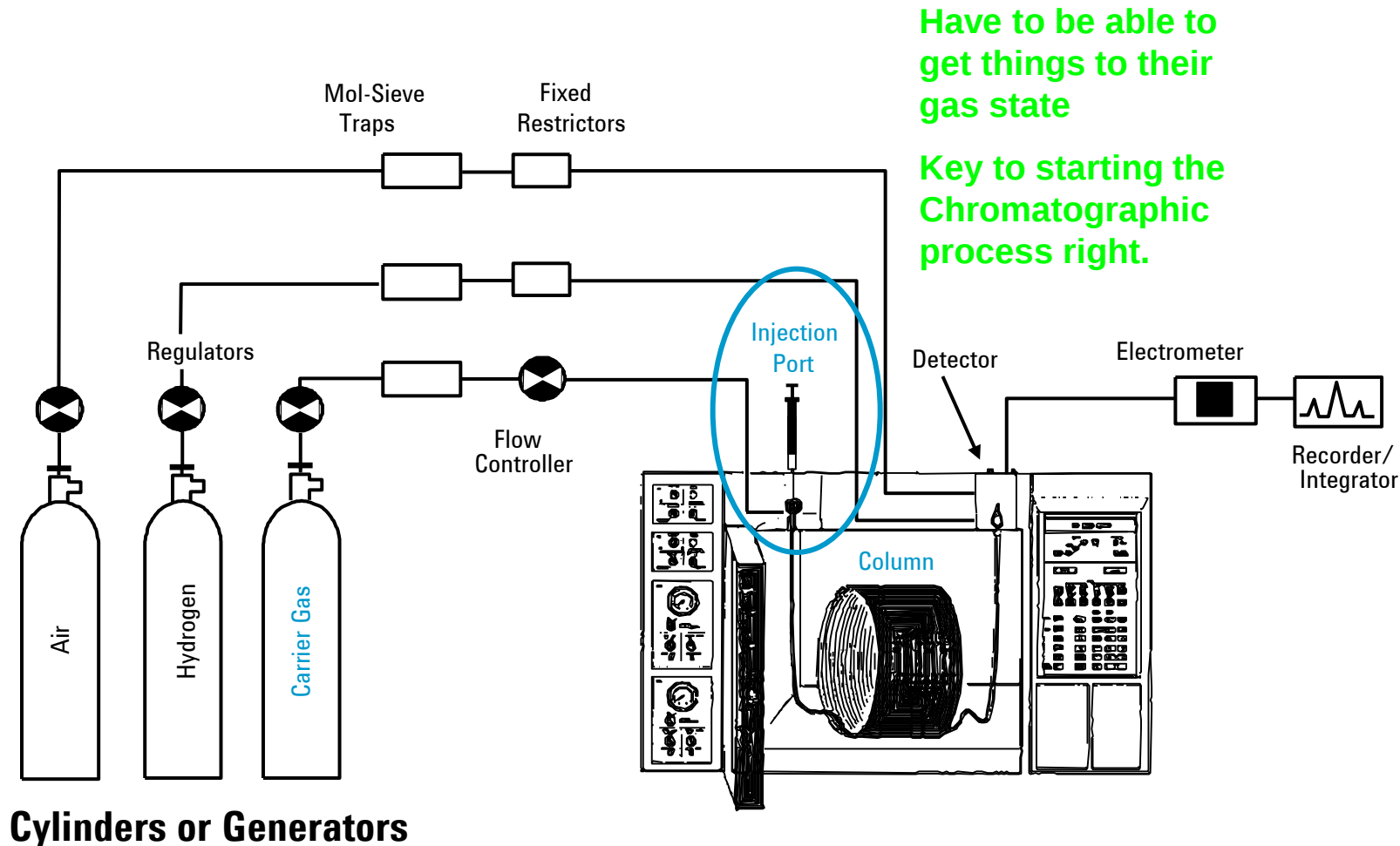
Replacement means delay

Replacement column costs money

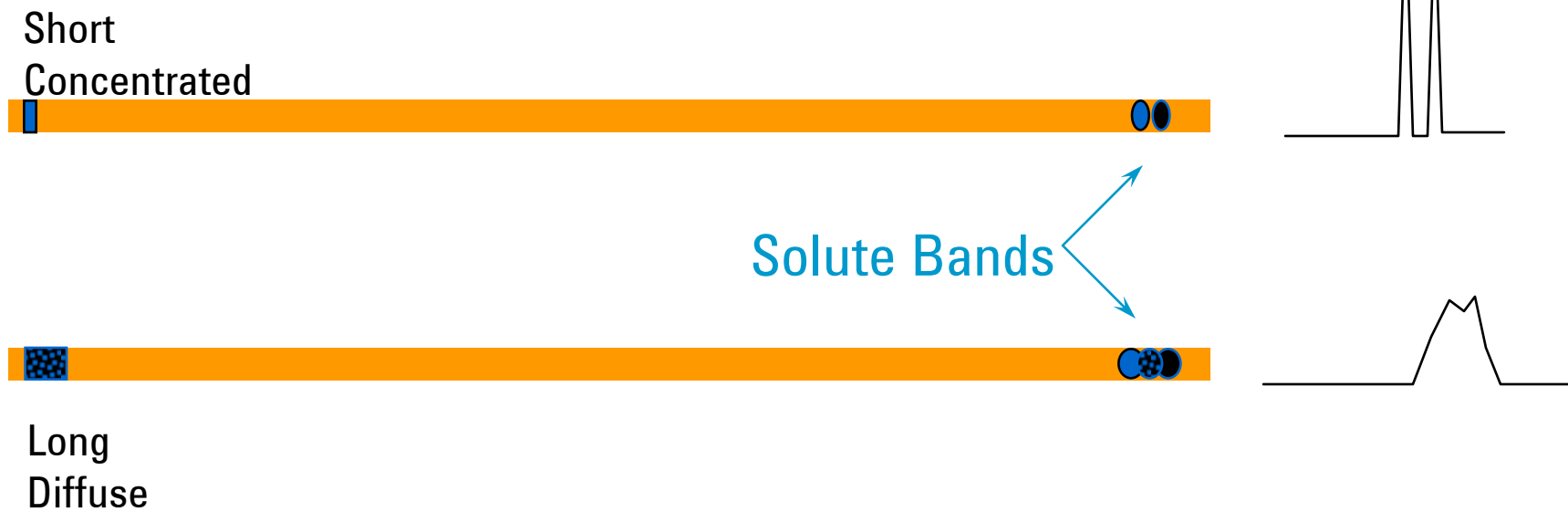
Solution = No Leaks!



Typical Gas Chromatographic System



Influence of Injection Efficiency



Same column, same chromatographic conditions

Solvent Vapor Volume Calculator

Vapor Volume Calculator

Solvent Properties

Acetonitrile

Boiling Point (°C) : 81.6

Density (g/cm³) : 0.782

Mol Wt. (amu) : 41.05

Injection Liner

5183-4647 single-tapered sy

Liner Volume (µL) : 850

Injection volume (µL)

1.00

Inlet Temperature (°C)

250

Inlet Pressure (gauge)

12.08

kPa ☒ psi ☐ bar

Estimated Volume

448 µL

% Capacity

52%

Solvents

Add Remove Defaults

Liners

Add Remove Defaults

Vapor Volume Calculator

Liner capacity exceeded! Choose a liner of greater volume or modify method parameters.

Solvent Properties

Water

Boiling Point (°C) : 100

Density (g/cm³) : 0.998

Mol Wt. (amu) : 18.02

Injection Liner

19251-60540 straight split

Liner Volume (µL) : 990

Injection volume (µL)

1.00

Inlet Temperature (°C)

280

Inlet Pressure (gauge)

14.993

kPa ☐ psi ☒ bar

Estimated Volume

1243 µL

% Capacity

125%

Solvents

Add Remove Defaults

Liners

Add Remove Defaults

The BIGGEST Problem in GC is...

There are more things that DON'T go through a GC than DO!

....therefore, don't inject anything and you'll never have problems.

OK, inject, but realize that everything just got dirty...deal with it!

Where Does it Get Dirty?

Here

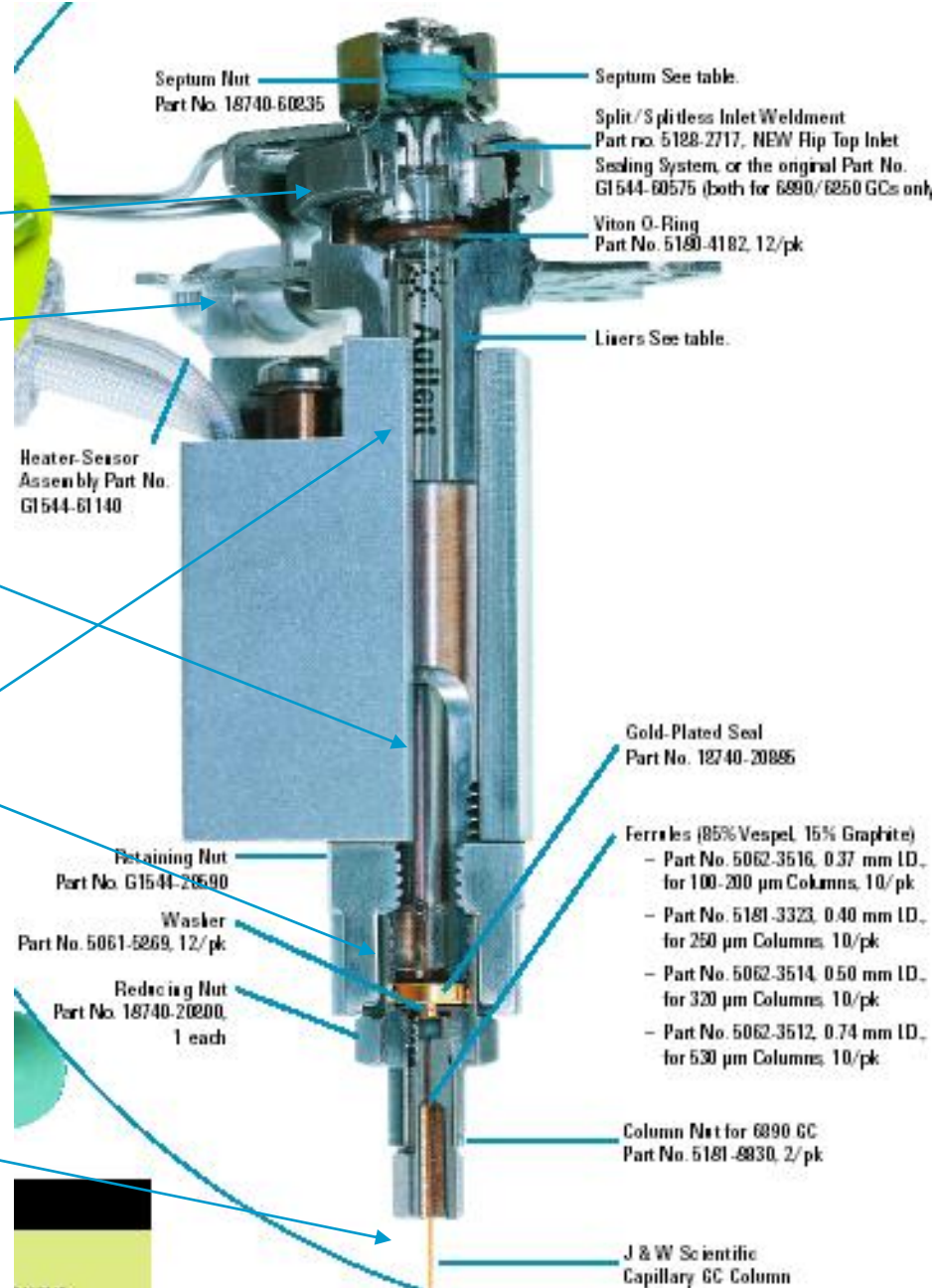
Here

Here

Here

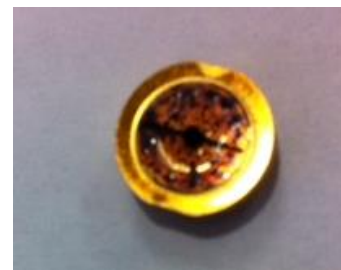
Here

Here



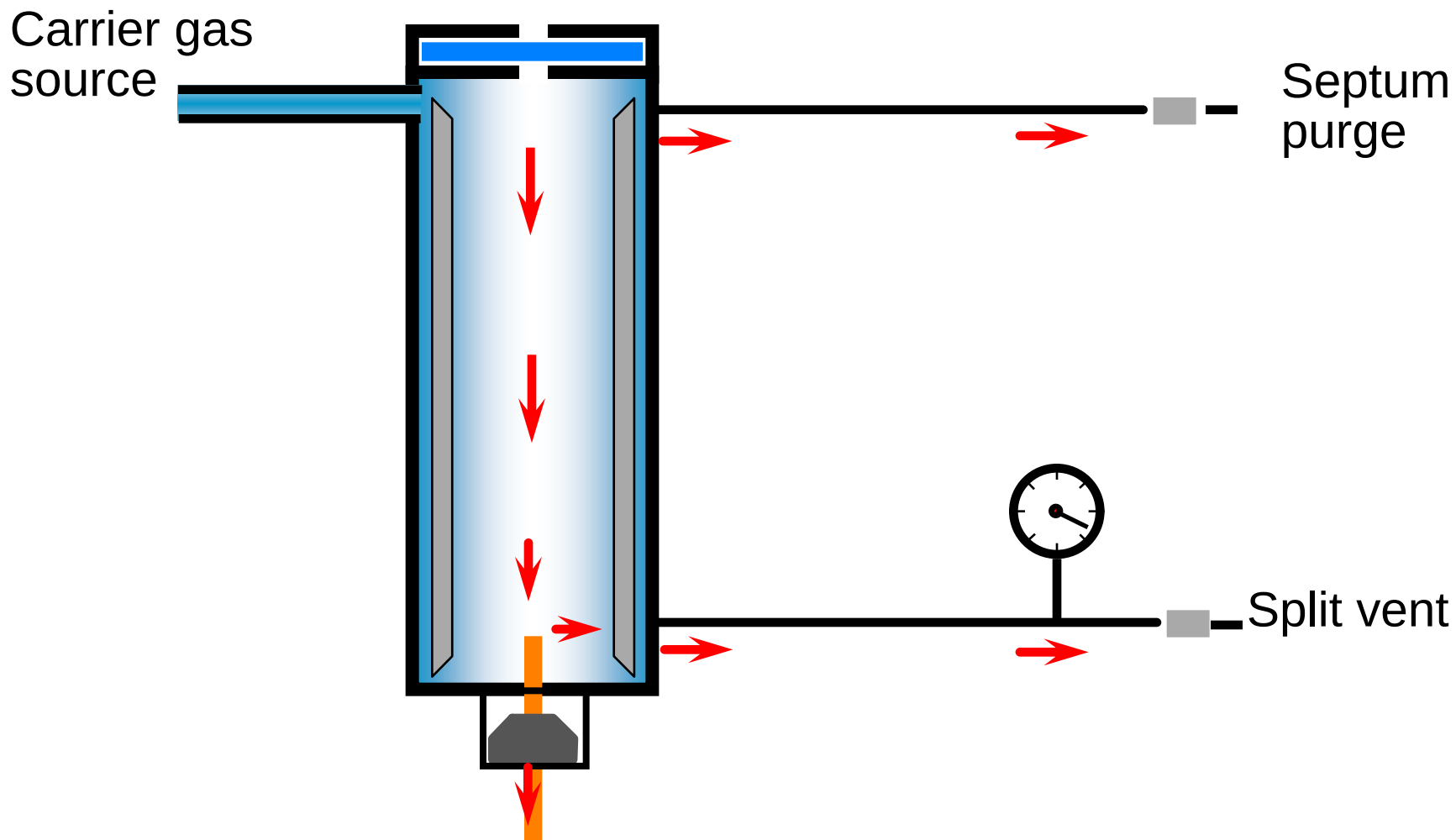
Agilent Technologies

What Are You Doing!?

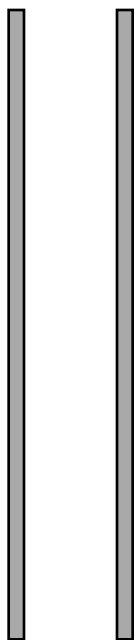


Split Injector

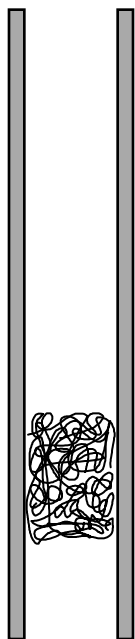
Flow Path



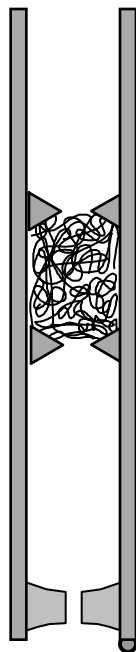
Split Liners – What's What?



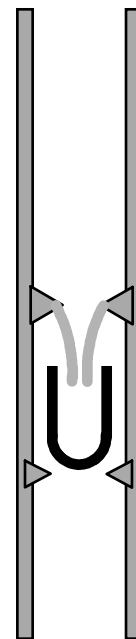
Straight
tube



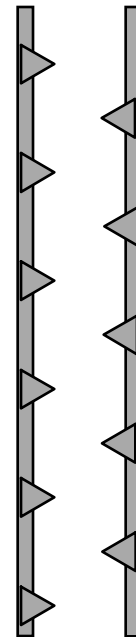
Straight
tube with
glass wool



Fixed glass
wool



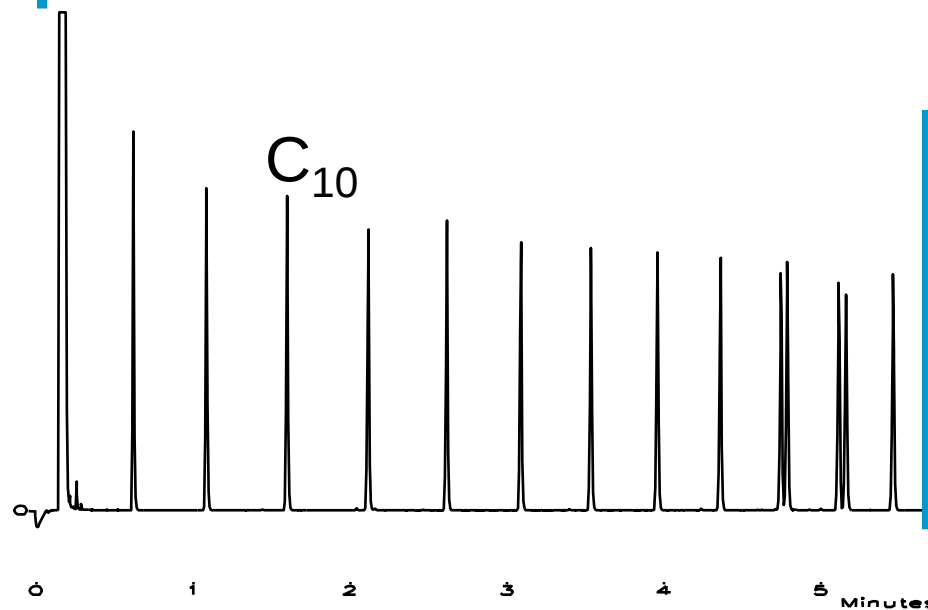
Inverted
cup



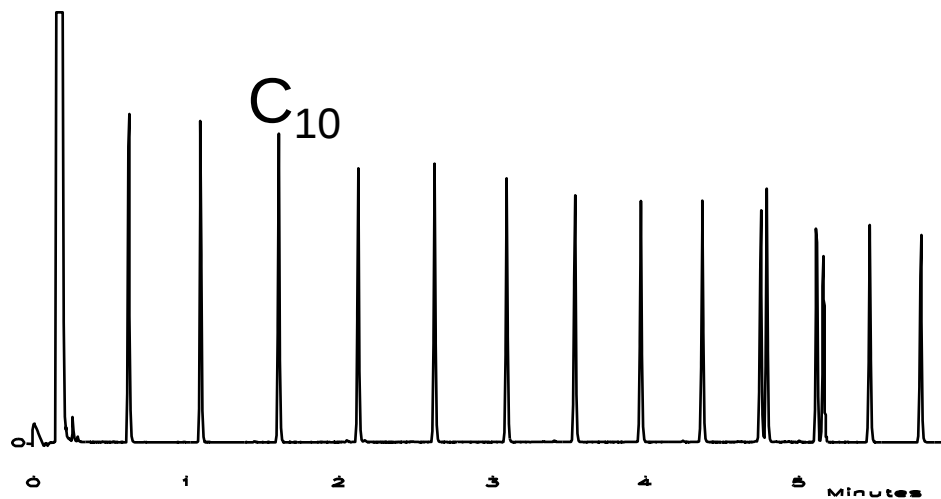
Baffle

Split Liner

Packed with Glass Wool



Without Glass Wool Packing

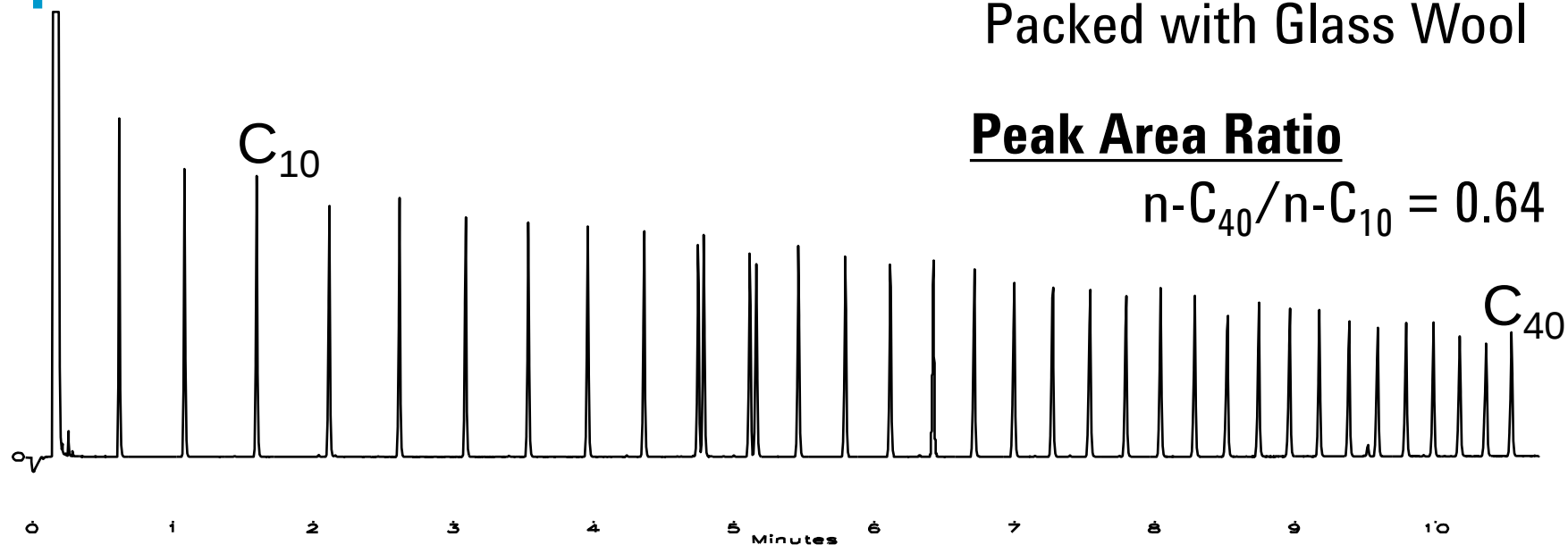


Split Liner

Packed with Glass Wool

Peak Area Ratio

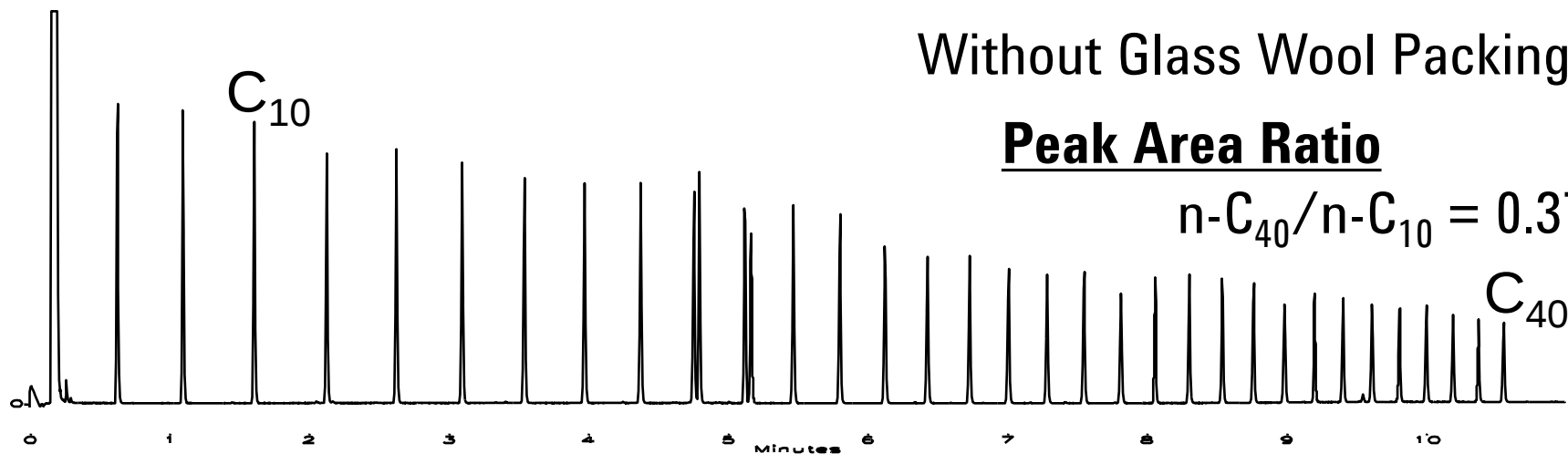
$$n\text{-C}_{40}/n\text{-C}_{10} = 0.64$$



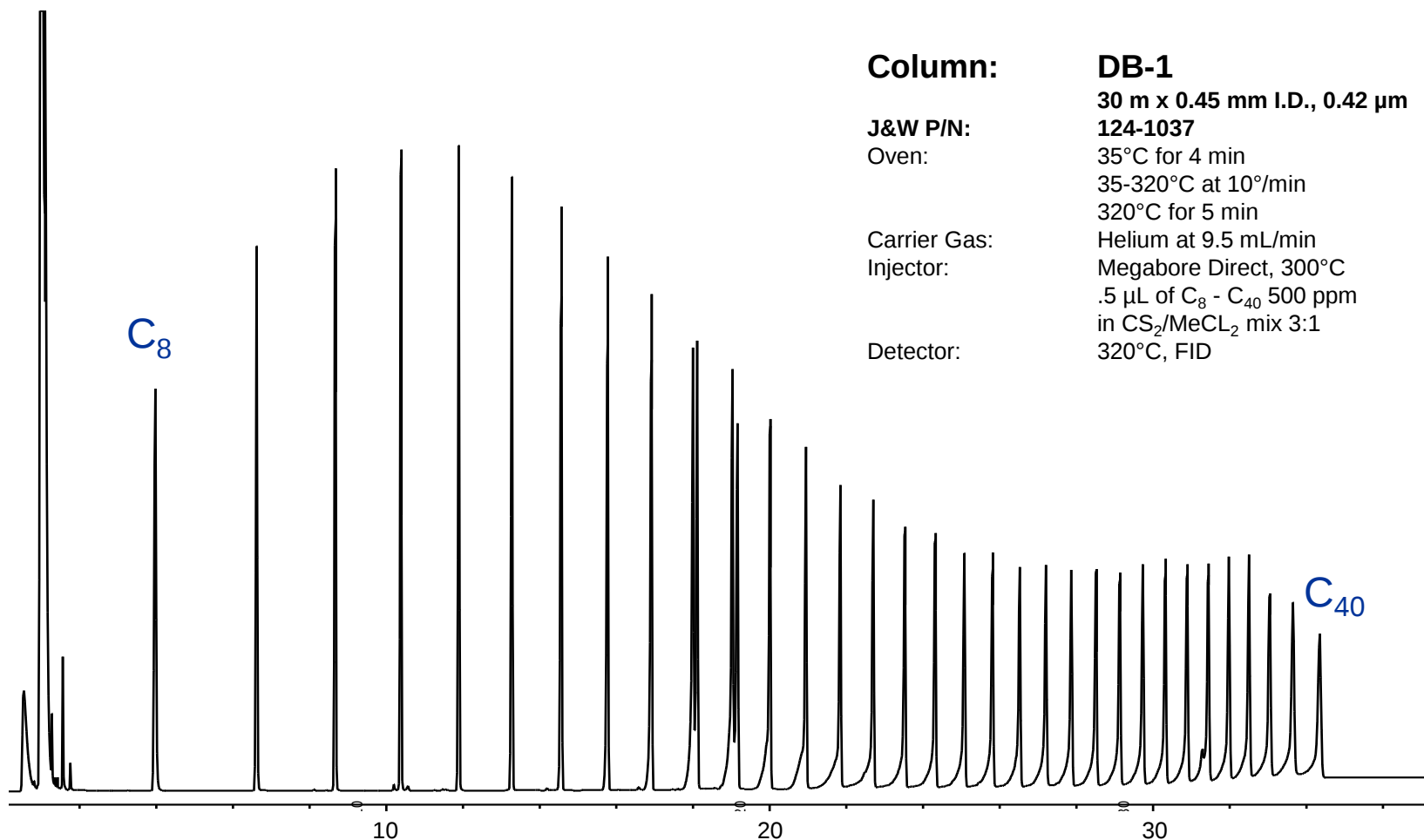
Without Glass Wool Packing

Peak Area Ratio

$$n\text{-C}_{40}/n\text{-C}_{10} = 0.37$$

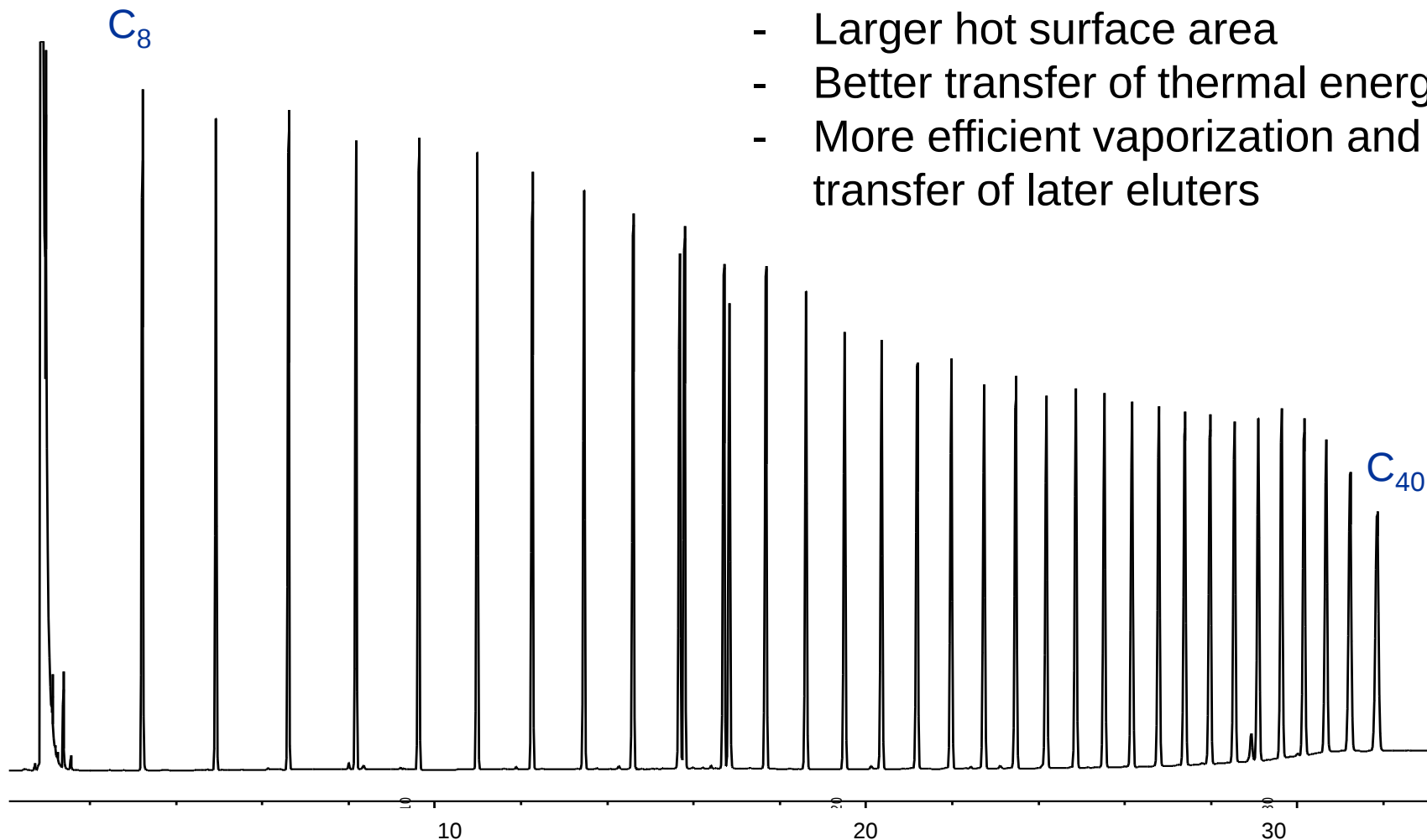


Fronting and Discrimination of Later Eluters



Large Glass Wool Plug Added to Liner

- Larger hot surface area
- Better transfer of thermal energy
- More efficient vaporization and transfer of later eluters



Discrimination

Considerations

Efficient heat transfer to injected sample

Efficient mixing of vaporized sample with carrier gas

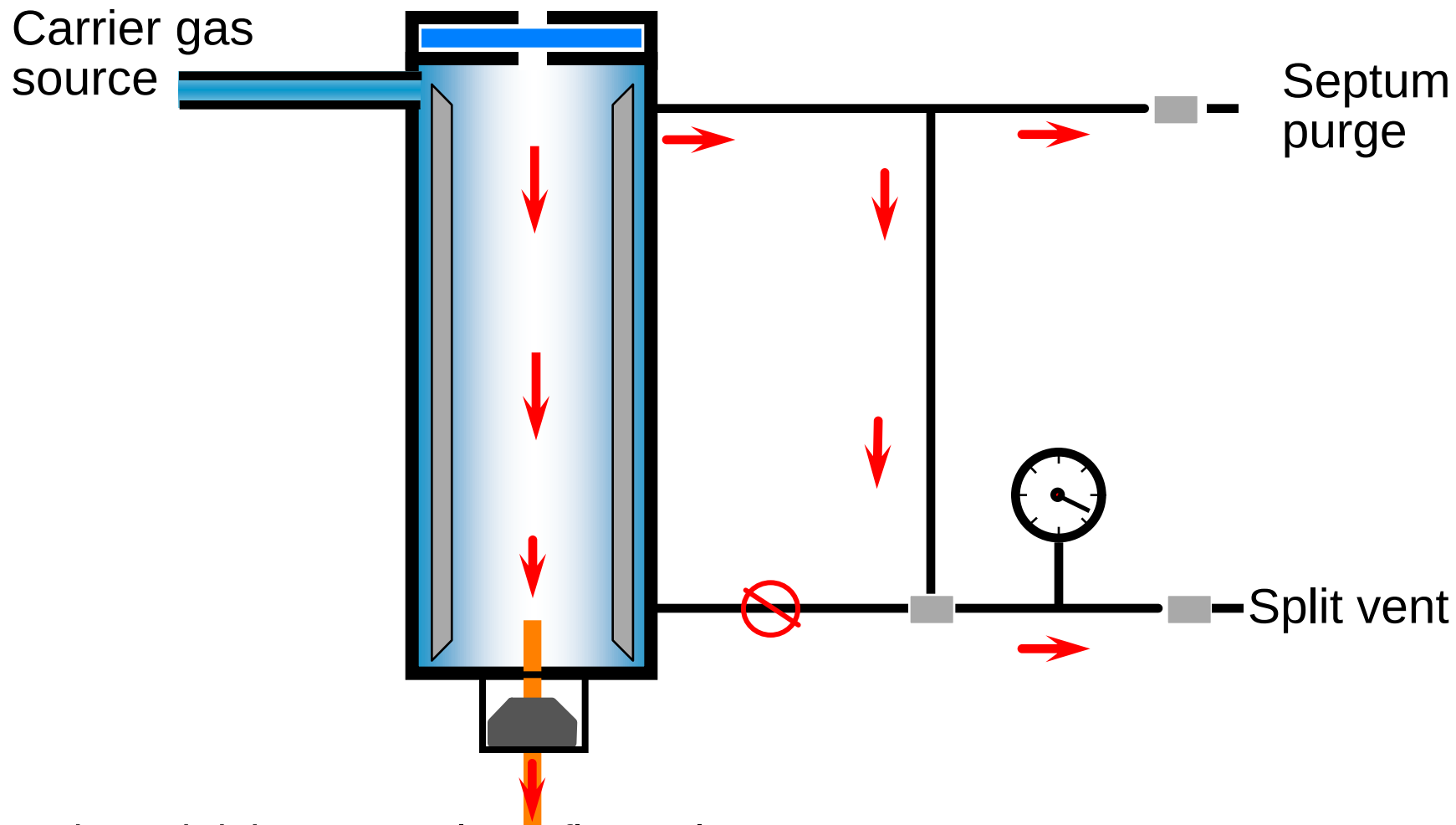
Column position in inlet

Consistent conditions are important



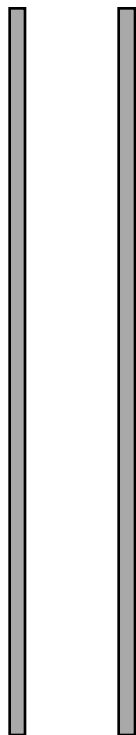
Splitless Injector

Purge Off At Injection

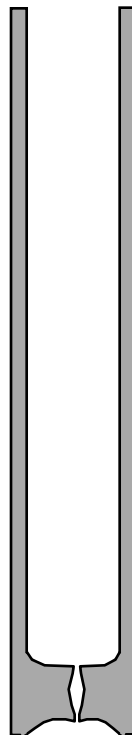


Splitless Injector

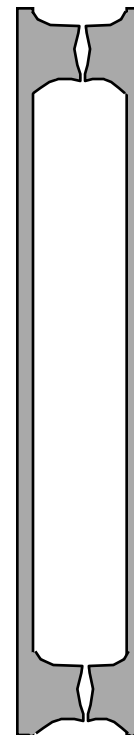
Liners



Straight
tube



Bottom
restriction



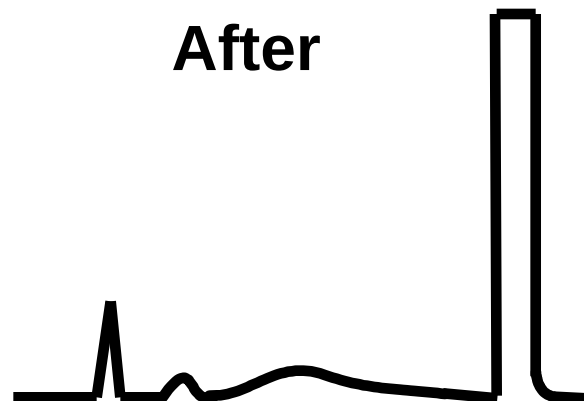
Dual
restriction

Bonus Peaks or Ghost Peaks

Before



After



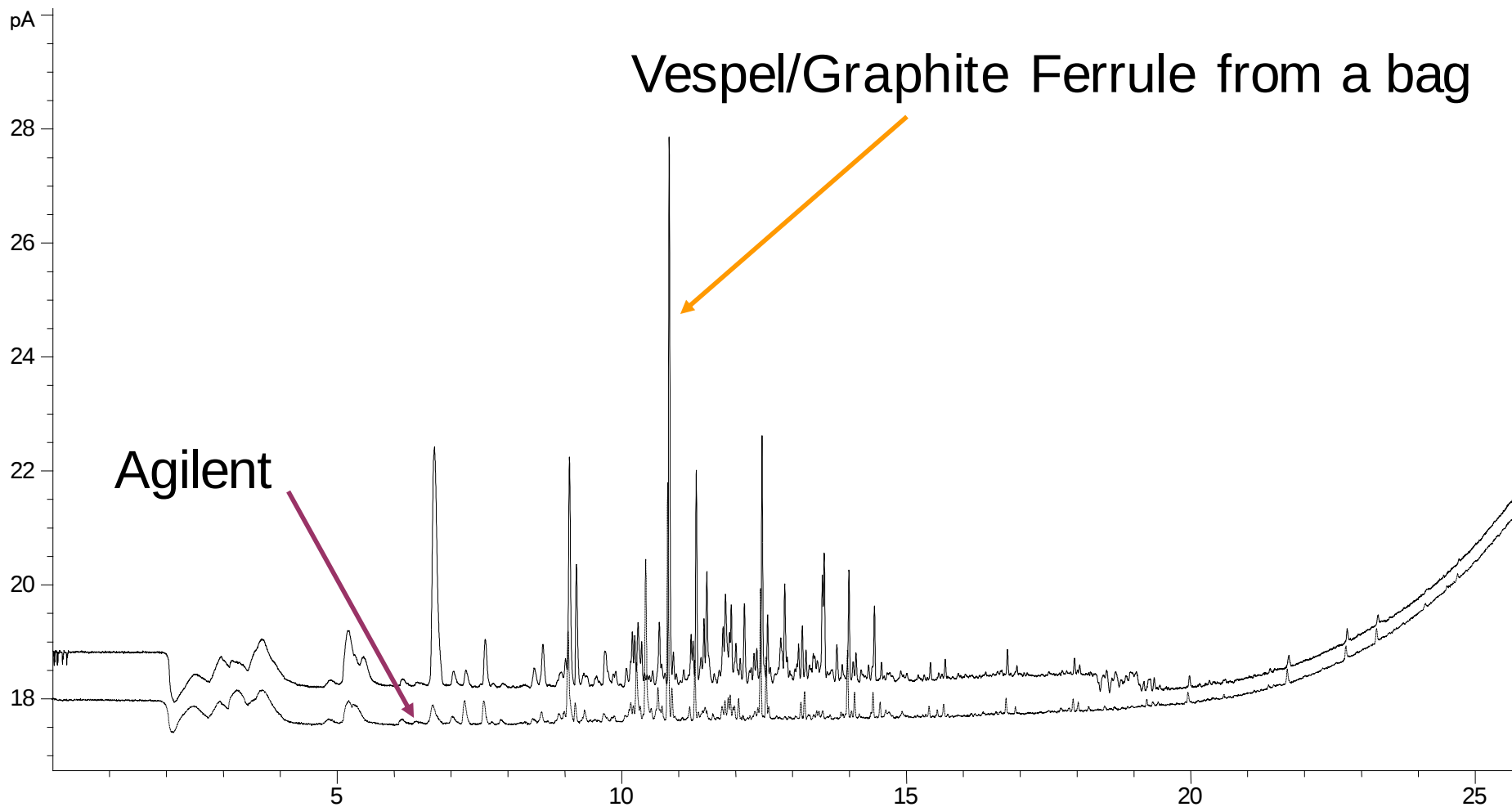
Contamination in INJECTOR or FLOW (carrier gas)

- Contaminated consumables
- Carryover from a backflash or previous sample
- Bad tank of gas or traps have expired
- Septum bleed

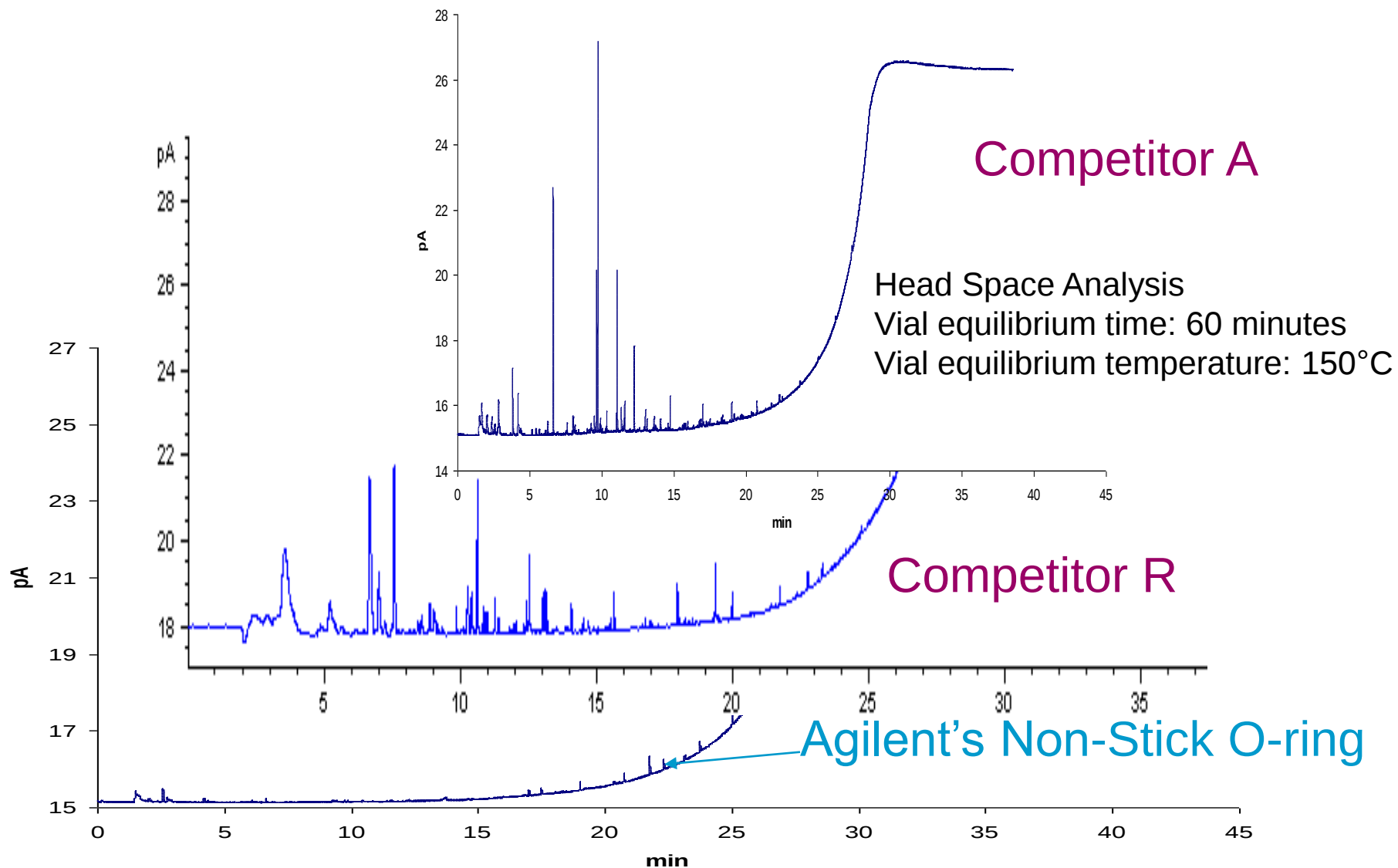
***TIP = Run a blank run...it should be blank!**



Bonus Peaks - Ferrule Contamination

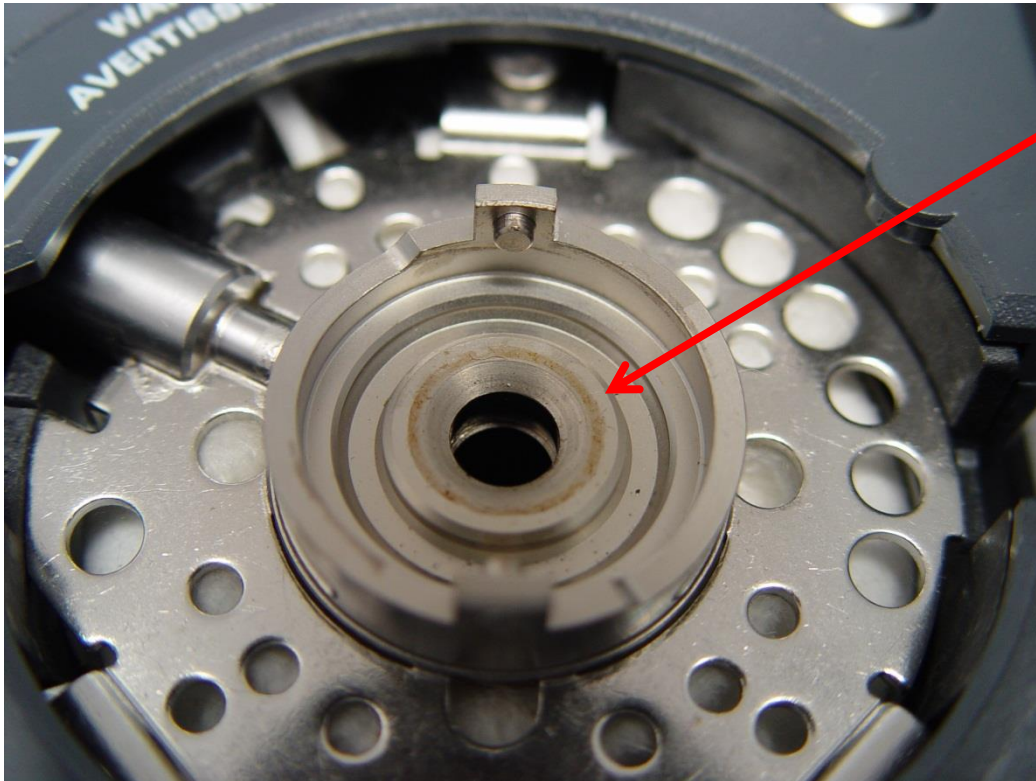


Bonus Peaks - O-Ring Contamination



More “Off-Brand” O-Ring Issues

Controlled Substances Analysis, H2 Carrier

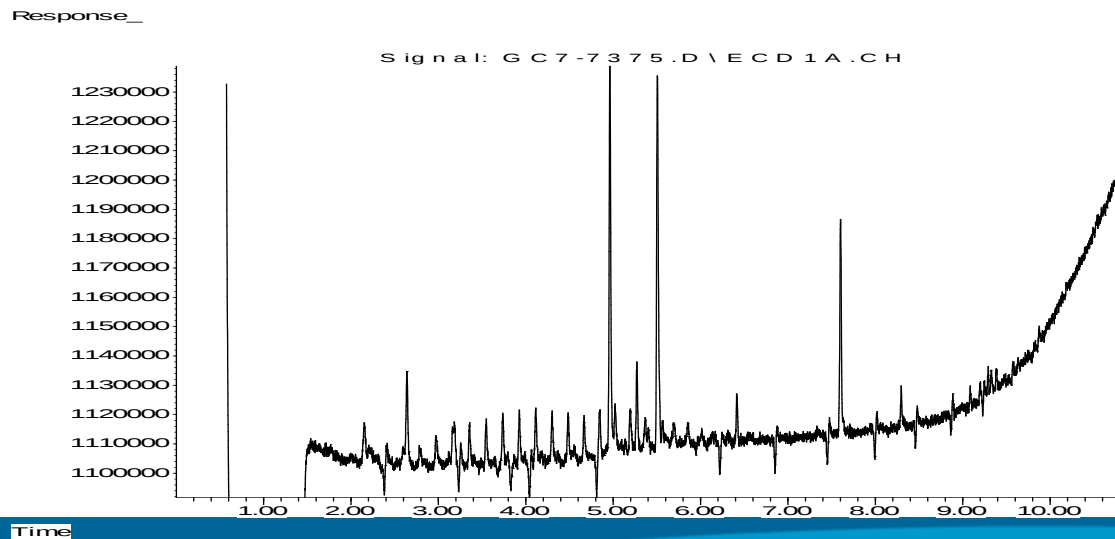
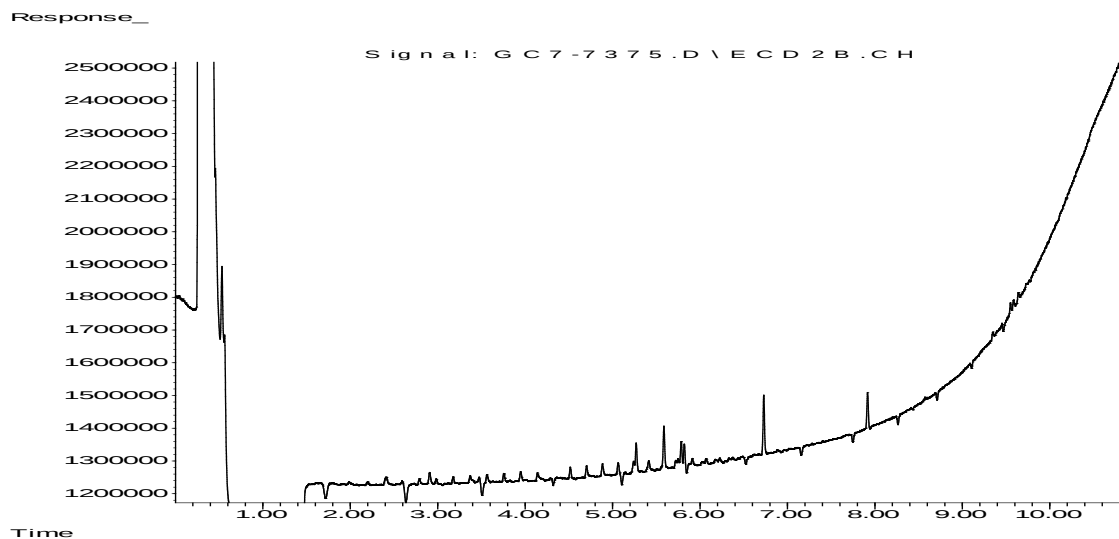


Residue on top of
inlet weldment

Problem Resolution:

Agilent Non-Stick Liner O-Ring
p/n 5188-5365, 10PK

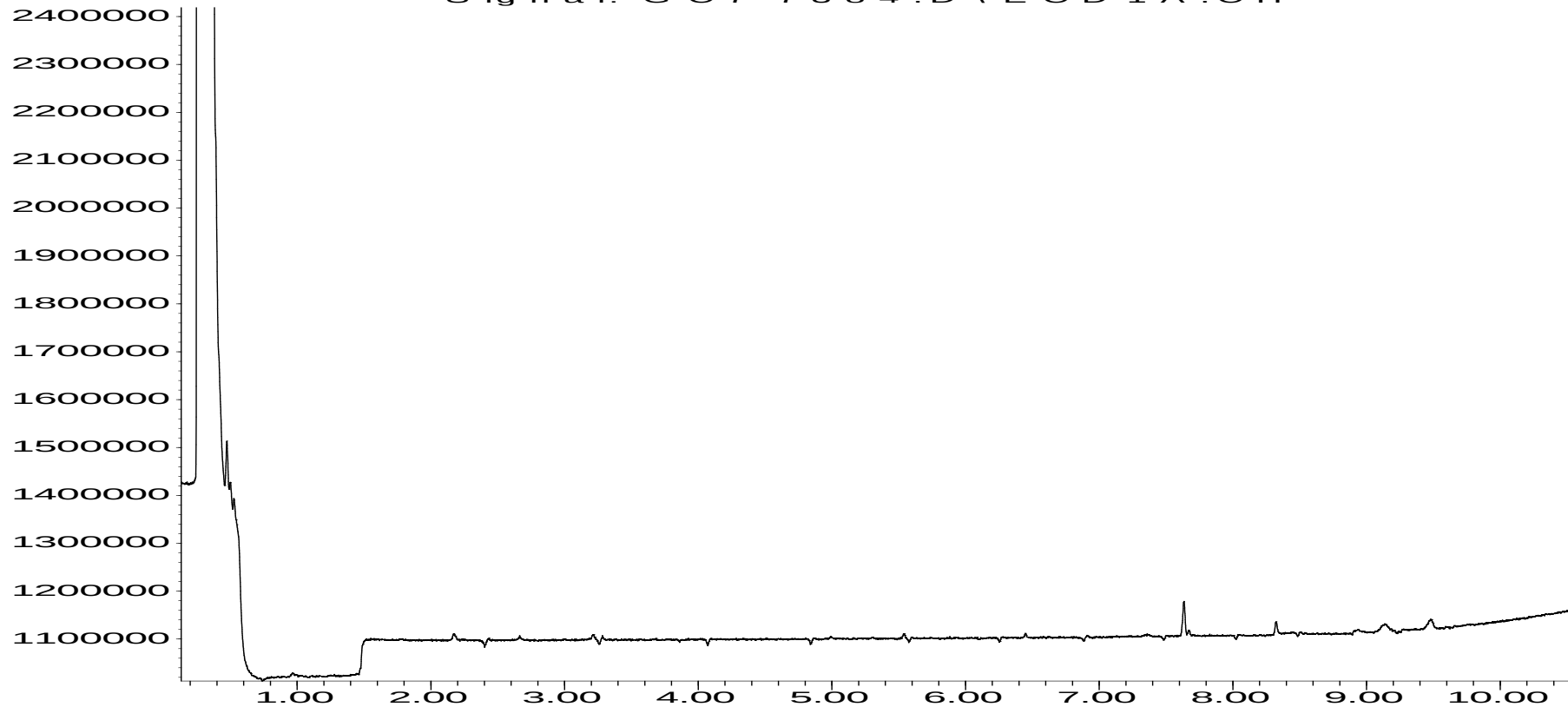
Contaminated Vial Inserts, New, Bag 1 and Bag 2



Hexane Rinsed Insert

Response_

Signal: GC7-7384.D \ ECD1A.CH



Time

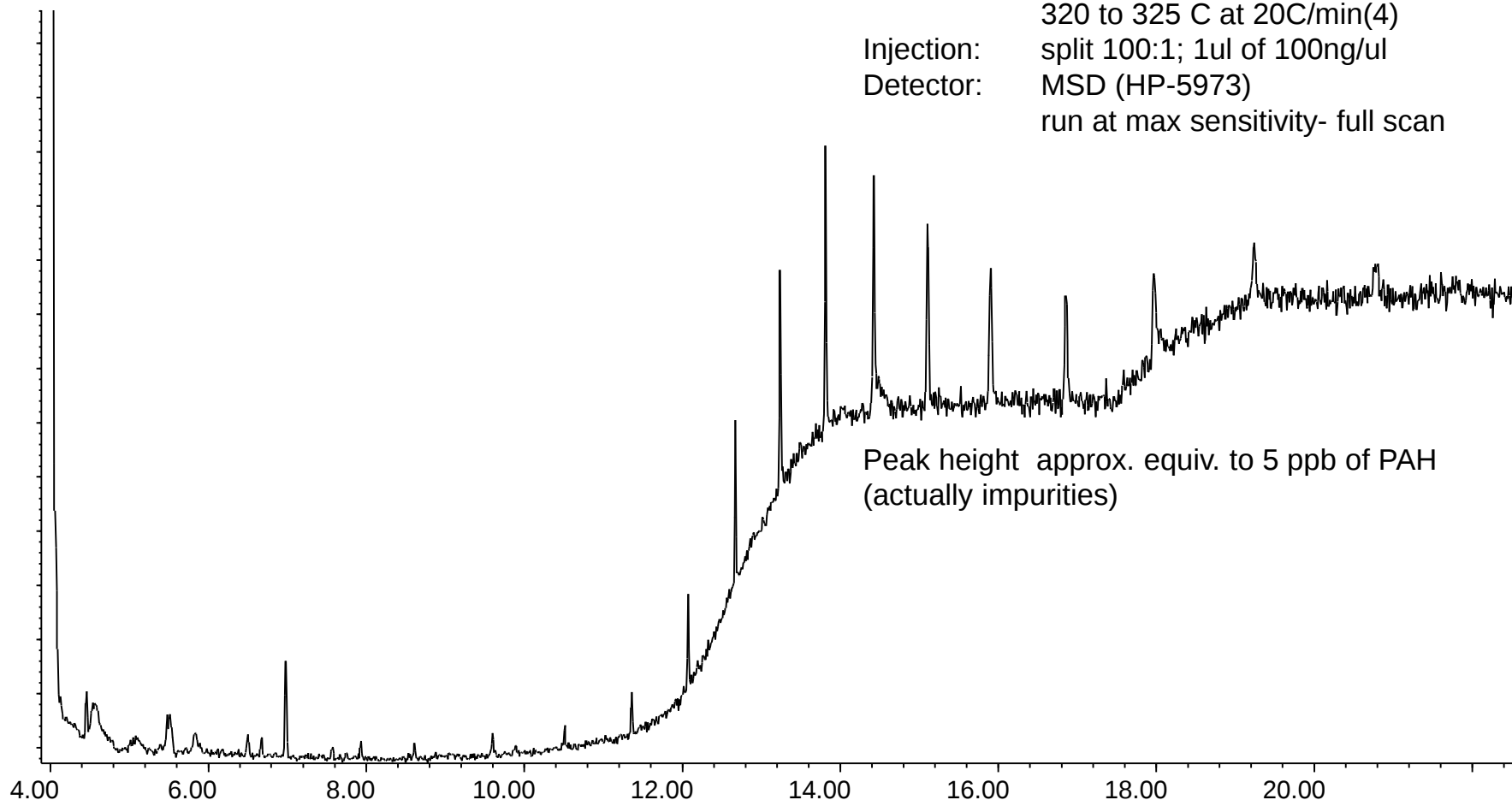


Agilent Technologies

Septa Bleed vs Column Bleed (MSD)

Source of peaks from outside of the column

Columns: HP 5MS
30mx0.25mmx0.25um
Oven: 80 to 160C at 25 C/min,
160 to 320 C at 3 C/min(4),
320 to 325 C at 20C/min(4)
Injection: split 100:1; 1ul of 100ng/ul
Detector: MSD (HP-5973)
run at max sensitivity- full scan



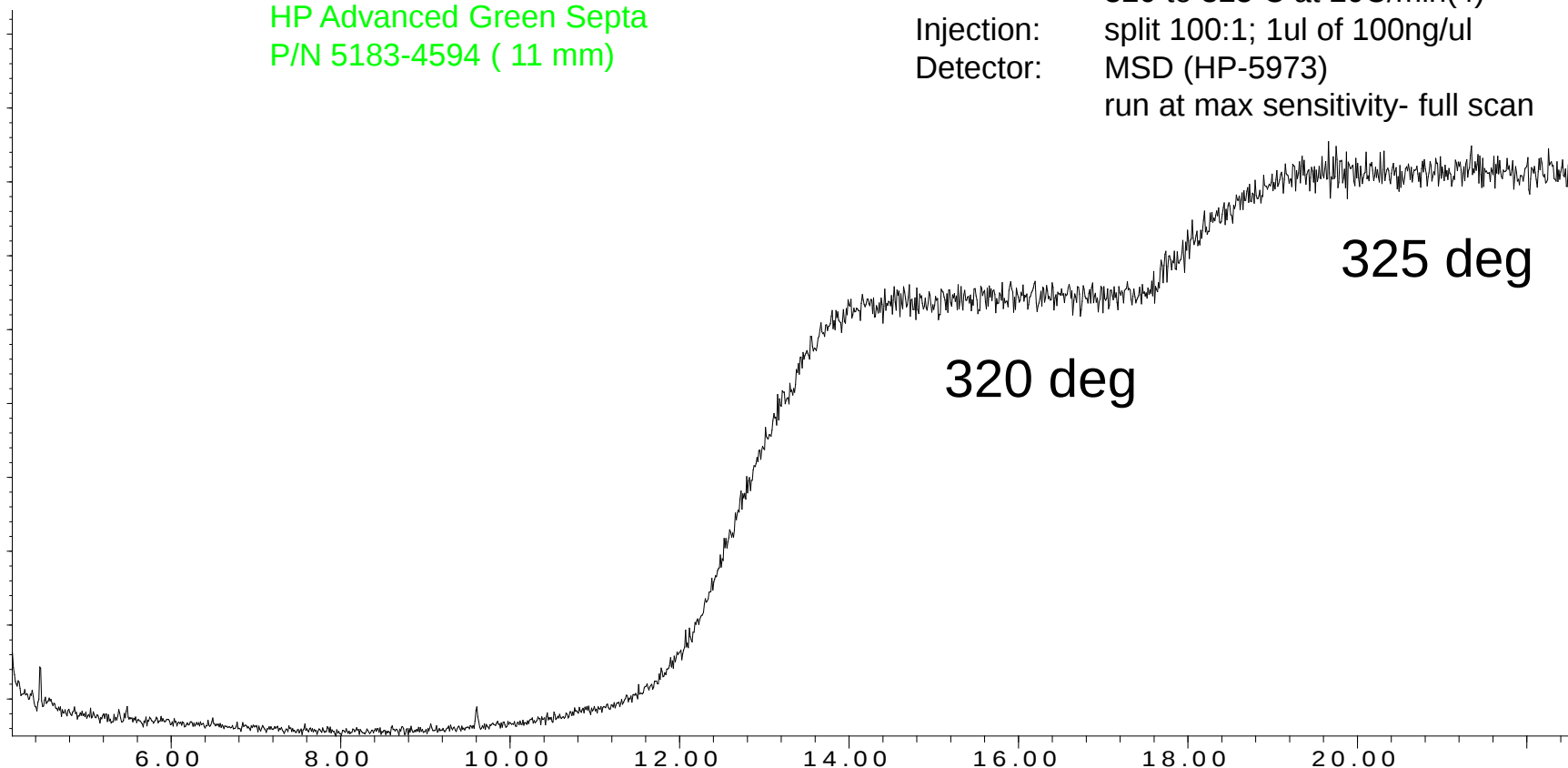
Peak height approx. equiv. to 5 ppb of PAH
(actually impurities)

Septa Bleed vs Column Bleed (MSD)

Source of peaks from outside of the column *ELIMINATED*

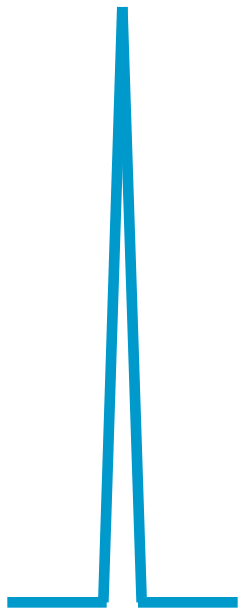
New Septa Installed
HP Advanced Green Septa
P/N 5183-4594 (11 mm)

Columns: HP 5 MS
30mx0.25mmx0.25um
Oven: 80 to160 C at 25 C/min,
160 to 320 C at 3 C/min(4)
320 to 325 C at 20C/min(4)
Injection: split 100:1; 1ul of 100ng/ul
Detector: MSD (HP-5973)
run at max sensitivity- full scan



Peak Tailing

Before



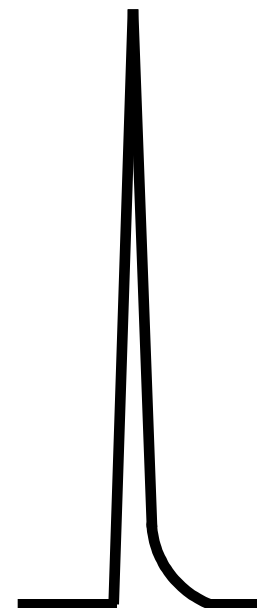
INJECTOR or COLUMN is Active

- Reversible adsorption of active compounds
(-OH, -NH, -SH)

FLOW problem

- dead volume, obstruction, poor installation, or
severe column contamination

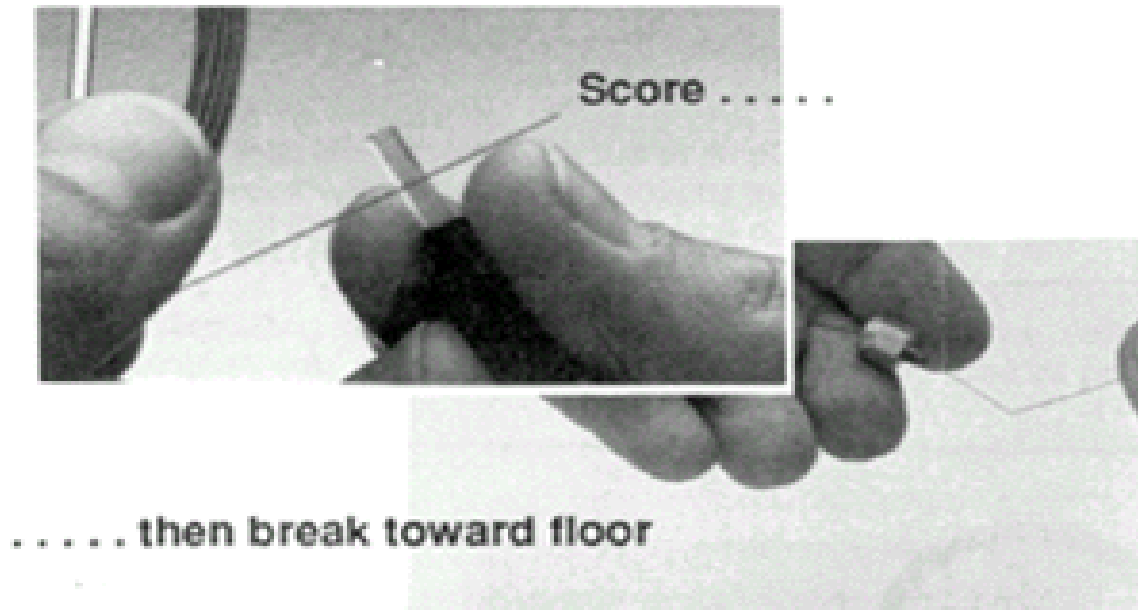
After



Miscellaneous – temperature issues for late eluters, overloading of PLOT columns, co-elution, polarity mismatch between phase, solute or solvent, and some compounds always tail

***Tip = Inject a light hydrocarbon, should not tail unless flow path problem.**

Flow Path Matters

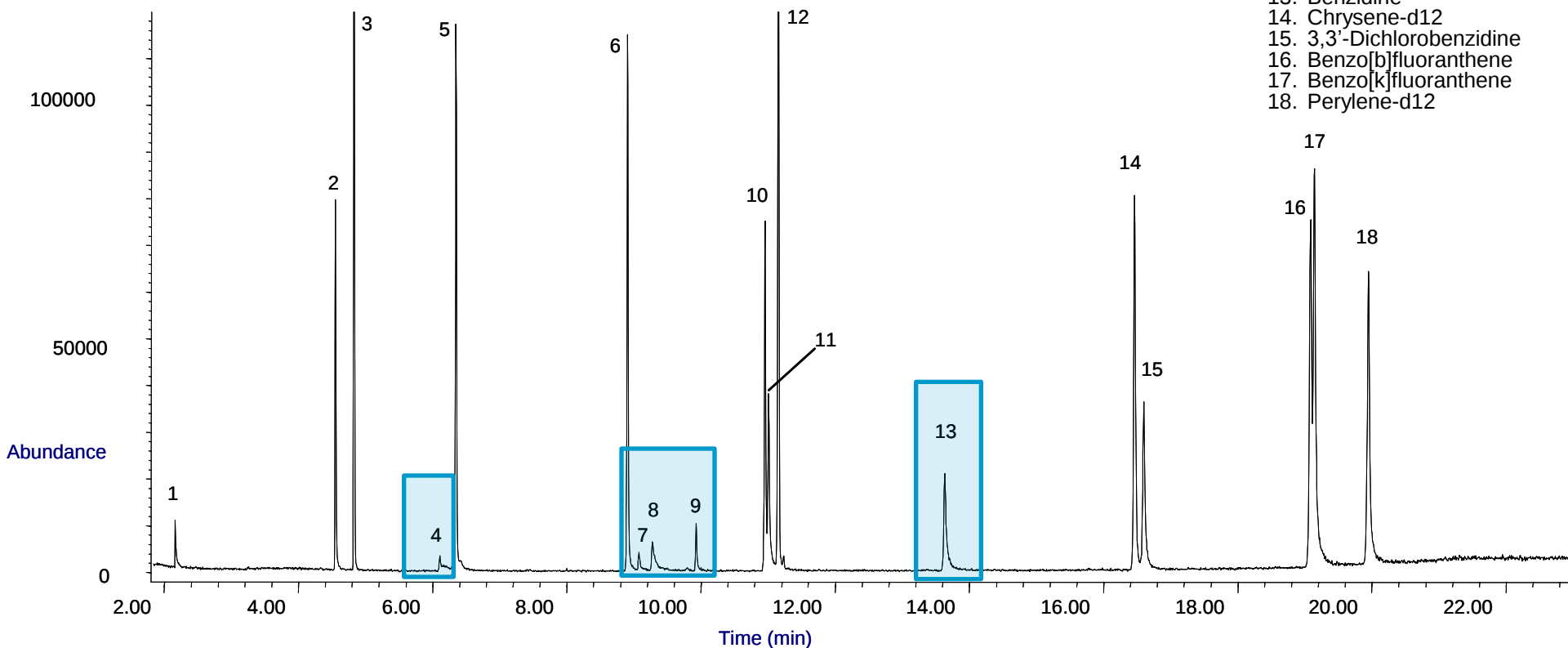


A flat, 90° square cut will be optimal for all connections

Symptom – Tailing of Active Compounds

Sample: 0.5 ng on column loading with ISTD
Column: 20m 0.18mm 0.18µm
Carrier: Helium 37cm/sec, Ramped flow; 0.7ml/min (0.1min) to 1.3ml/min (15ml/min²)
Oven: 35°C (2.5 min) to 80°C (40°C/min), 15°C/min to 200°C, 8°C/min to 275°C (2 min)
Injection: 0.5µl, Splitless, 280°C, purge flow 30ml/min at 0.75 min
MSD: Transfer Line 290°C, Source 300°C, Quad 180°C

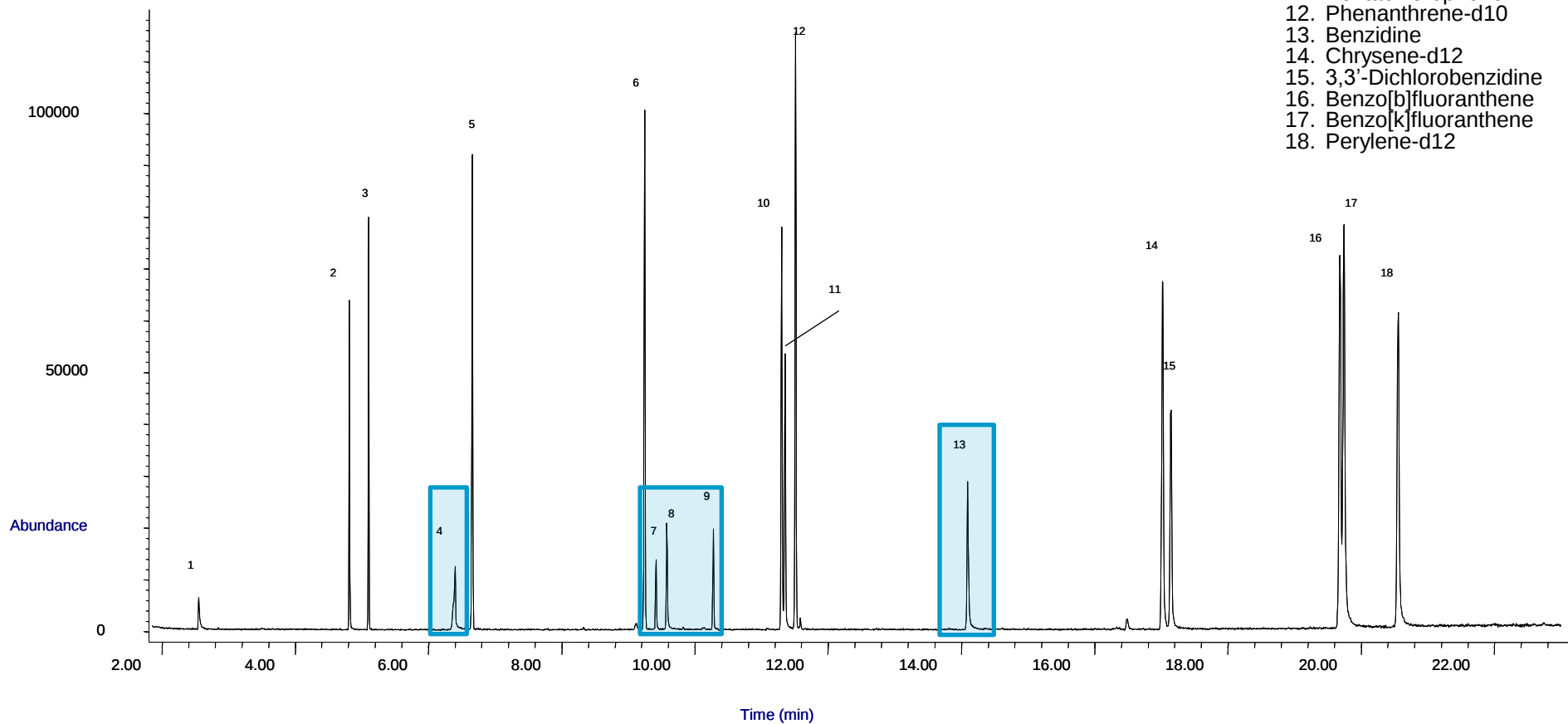
1. n-Nitrosodimethylamine
2. Aniline
3. 1,4-Dichlorobenzene-d4
4. Benzoic Acid
5. Naphthalene-d8
6. Acenaphthene-d10
7. 2,4-Dinitrophenol
8. 4-Nitrophenol
9. 2-Me-4,6-dinitrophenol
10. 4-Aminobiphenyl
11. Pentachlorophenol
12. Phenanthrene-d10
13. Benzidine
14. Chrysene-d12
15. 3,3'-Dichlorobenzidine
16. Benzo[b]fluoranthene
17. Benzo[k]fluoranthene
18. Perylene-d12



Solution – Ultra Inert GC Column

Sample: 0.5 ng on column loading with ISTD
Column: 20m 0.18mm 0.18µm
Carrier: Helium 37cm/sec, Ramped flow; 0.7ml/min (0.1min) to 1.3ml/min (15ml/min²)
Oven: 35°C (2.5 min) to 80°C (40°C/min), 15°C/min to 200°C, 8°C/min to 275°C (2 min)
Injection: 0.5µl, Splitless, 280°C, purge flow 30ml/min at 0.75 min
MSD: Transfer Line 290°C, Source 300°C, Quad 180°C

1. n-Nitrosodimethylamine
2. Aniline
3. 1,4-Dichlorobenzene-d4
4. Benzoic Acid
5. Naphthalene-d8
6. Acenaphthene-d10
7. 2,4-Dinitrophenol
8. 4-Nitrophenol
9. 2-Me-4,6-dinitrophenol
10. 4-Aminobiphenyl
11. Pentachlorophenol
12. Phenanthrene-d10
13. Benzidine
14. Chrysene-d12
15. 3,3'-Dichlorobenzidine
16. Benzo[b]fluoranthene
17. Benzo[k]fluoranthene
18. Perylene-d12



Agilent Ultra Inert Liners

Touchless packaging

Easy installation of new, clean liner

without risk of contamination from touching

Includes non-stick plasma treated O-ring



Instructions for Use



1 Squeeze cap sides tightly to hold liner as you remove plastic tube.



2 Align liner with inlet and gently release.

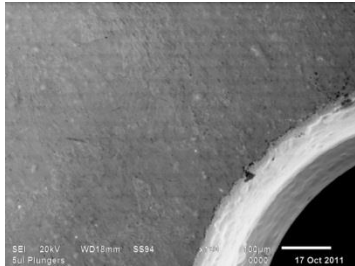


3 Use cap edge to press liner all the way down.

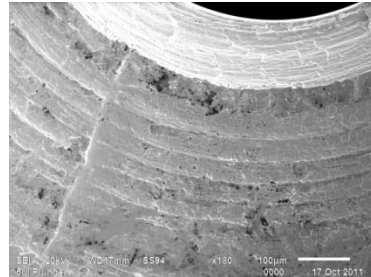


Agilent UI Gold Seal: Deactivated Gold Surface

- Soft gold plating is essential for proper sealing
- Ultra Inert chemistry blocks active sites (gold is NOT inert)
- Smooth surface doesn't leak



Agilent MIM seal



Competitor's
machined seal

*Reliable ppb and ppt
measurements require
attention to the little things!*

New Product: Agilent Inert Inlet

UltiMetal Plus treatment
creates inert surface 7890 inlet weldment & shell

- Limits adsorption/degradation of active analytes
- Optimum for Trace GC/MS and GC-ECD pesticide analyses



***Agilent's proven
proprietary UltiMetal
Plus surface
treatment***

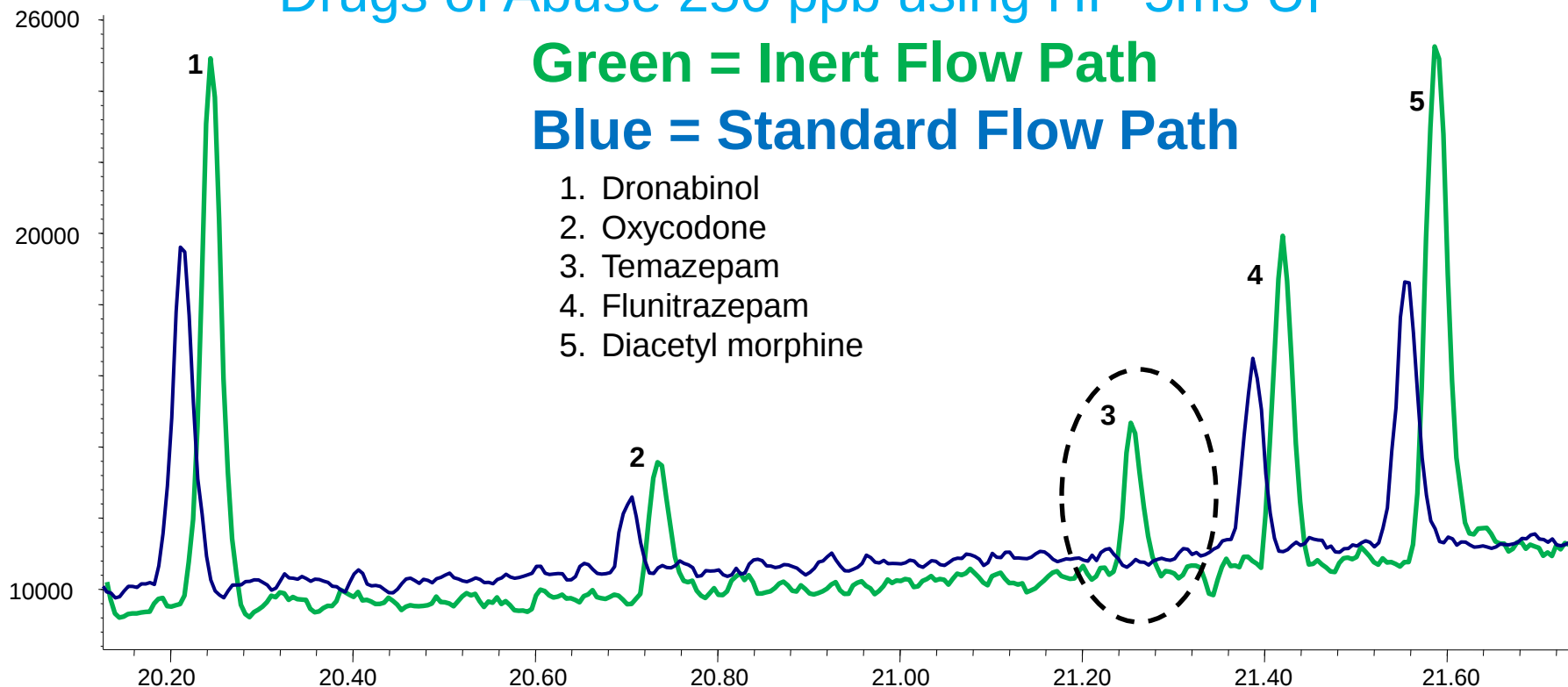
Ultra UltiMetal Plus Inert Inlet

Drugs of Abuse 250 ppb using HP-5ms UI

Green = Inert Flow Path

Blue = Standard Flow Path

1. Dronabinol
2. Oxycodone
3. Temazepam
4. Flunitrazepam
5. Diacetyl morphine

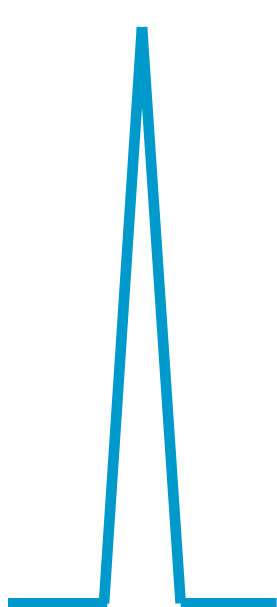


Column: Agilent J&W HP-5ms UI 30 m x 0.25 mm x 0.25 μ m
Oven: 100°C 4 min hold, 10°/min to 280°C, 6 °/min to 300°C (4.67 min hold),
Carrier : Helium 52.7 cm/s (2 mL/min) set at 100°C, EPC-Constant Flow
Inlet: Pulsed Splitless 35 PSI pulse until 0.73 min, 0.75 min purge 50 ml/min, gas saver 20, ml/min at 2 min
Inlet liner: Ultra Inert with wool / Standard single taper liner with wool (p/n 5190-3165)
Gold Seal: UI Gold Seal / Standard gold seal
Detector: MSD Scan mode 40 to 450 m/z, 230 °C source temp, 150 °C Quad temp, 310 °C transfer line



Agilent Technologies

Split Peaks

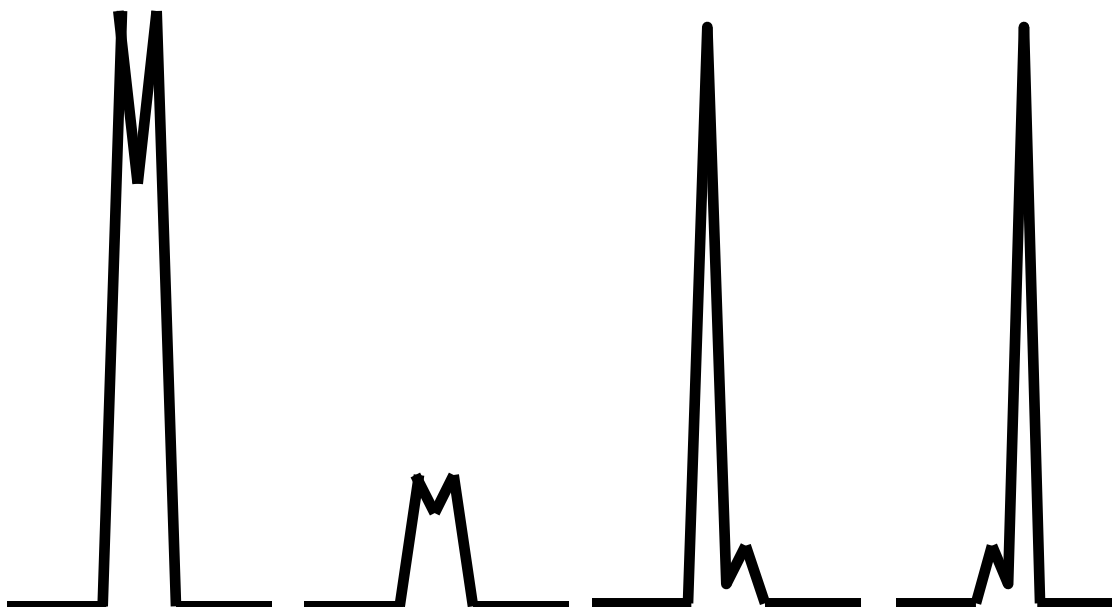


INJECTOR (poor sample introduction)

- Injecting the sample twice (some how?)
- Mixed sample solvent (polarity difference)
- Sample in syringe needle (manual inject)

INJECTOR (activity)

- Breakdown (not really a split peak, 2 peaks)
- Sample degradation in injector



VOLATILITY

High boilers dropping out
on Cold Spots

- Transfer line temps
- Unions or fittings not
tracking column temp



PAH's in Toluene

Splitting Peaks Immediately After Installation

Column: Agilent J&W DB-1ms

30 m x 0.25 mm I.D., 0.25 µm

P/N: 123-1522

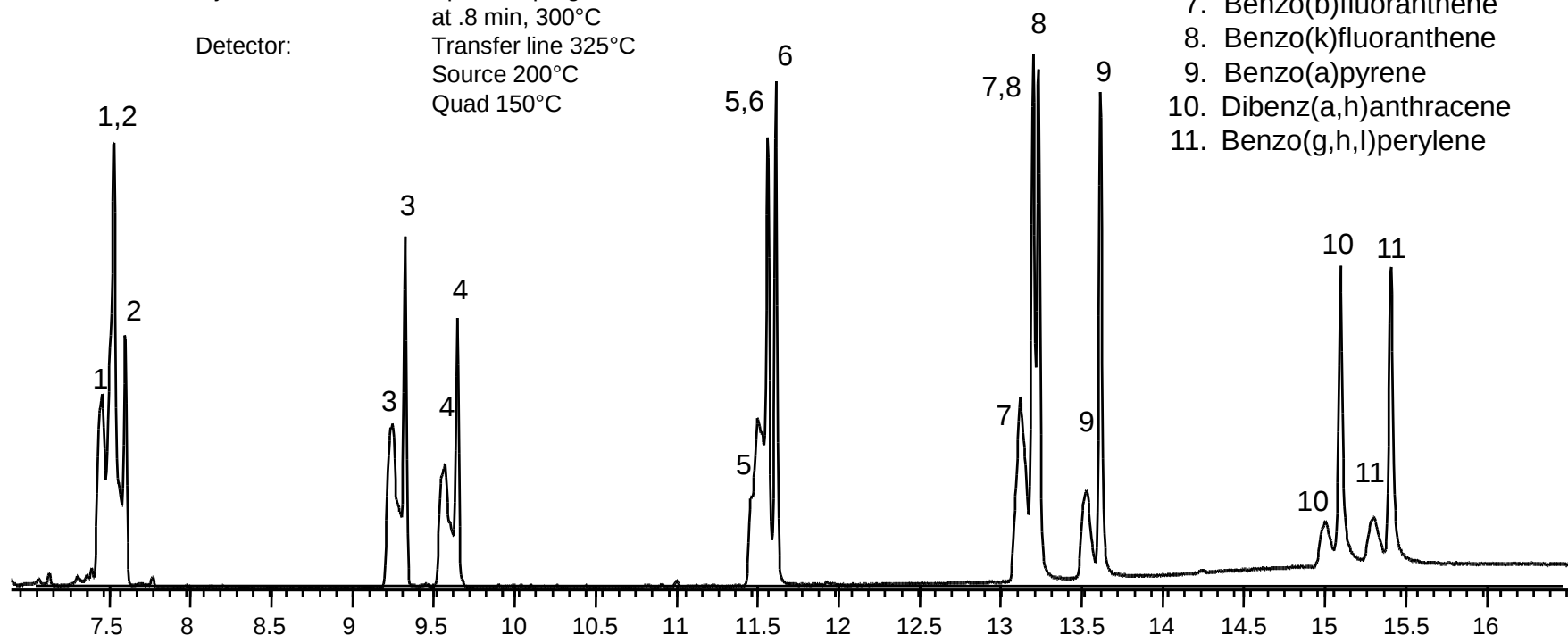
Test Temperature: 80°C for 0.8 min
80-325°C at 15°/min
325°C for 2 min

Carrier Gas: Helium at 40 cm/sec

Injector: Splitless, purge on
at .8 min, 300°C

Detector: Transfer line 325°C
Source 200°C
Quad 150°C

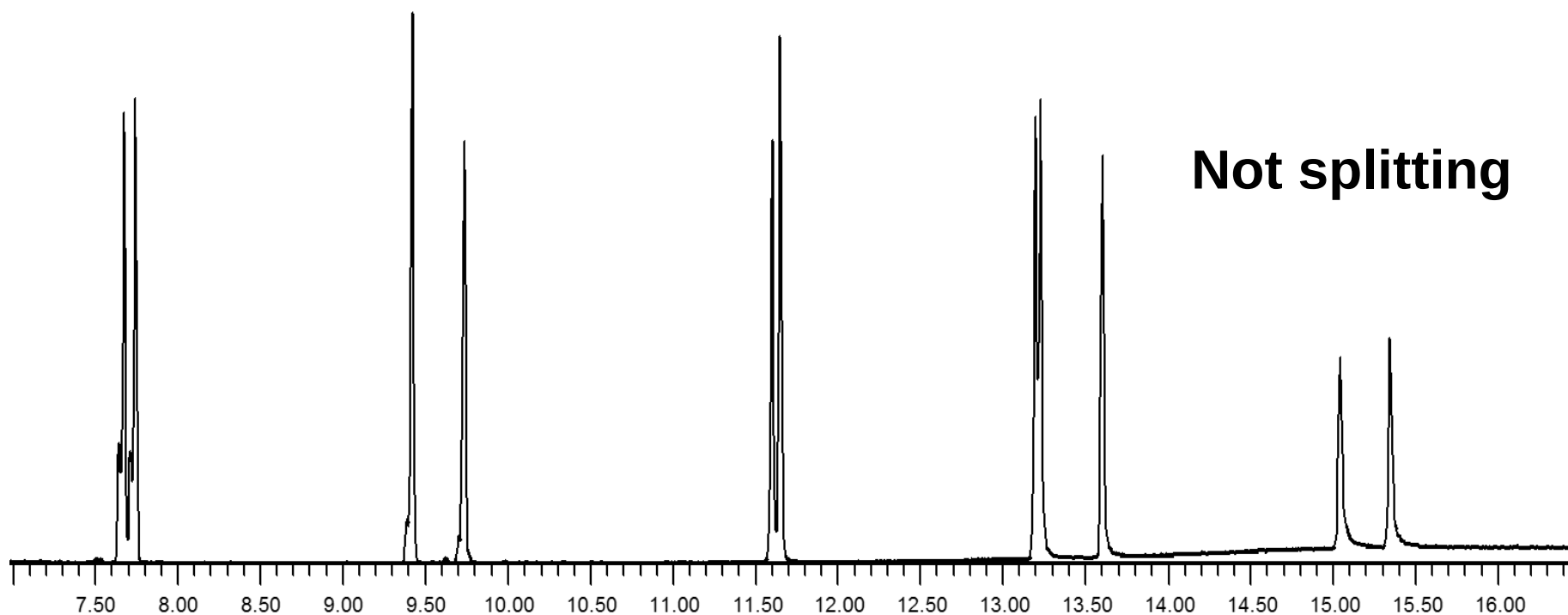
1. Phenanthrene
2. Anthracene
3. Fluoranthene
4. Pyrene
5. Benzo(a)anthracene
6. Chrysene
7. Benzo(b)fluoranthene
8. Benzo(k)fluoranthene
9. Benzo(a)pyrene
10. Dibenz(a,h)anthracene
11. Benzo(g,h,i)perylene



Reinstalled the Column

Distance was off - previously installed too far

Improved but still splitting



Broad Peaks

INJECTOR

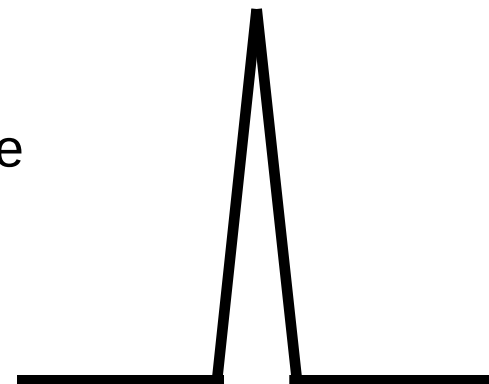
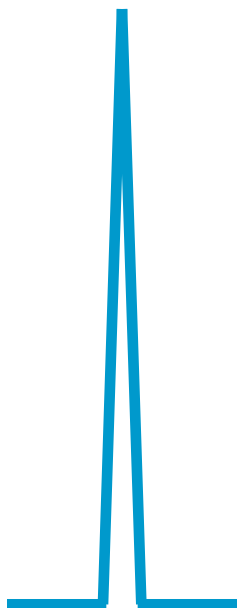
- Poor Installation
- Change in settings (temps/flows)
- Poor sample focusing
- Large change in sample concentration

FLOW

- Change in gas velocity
- Constant Flow vs Constant Pressure

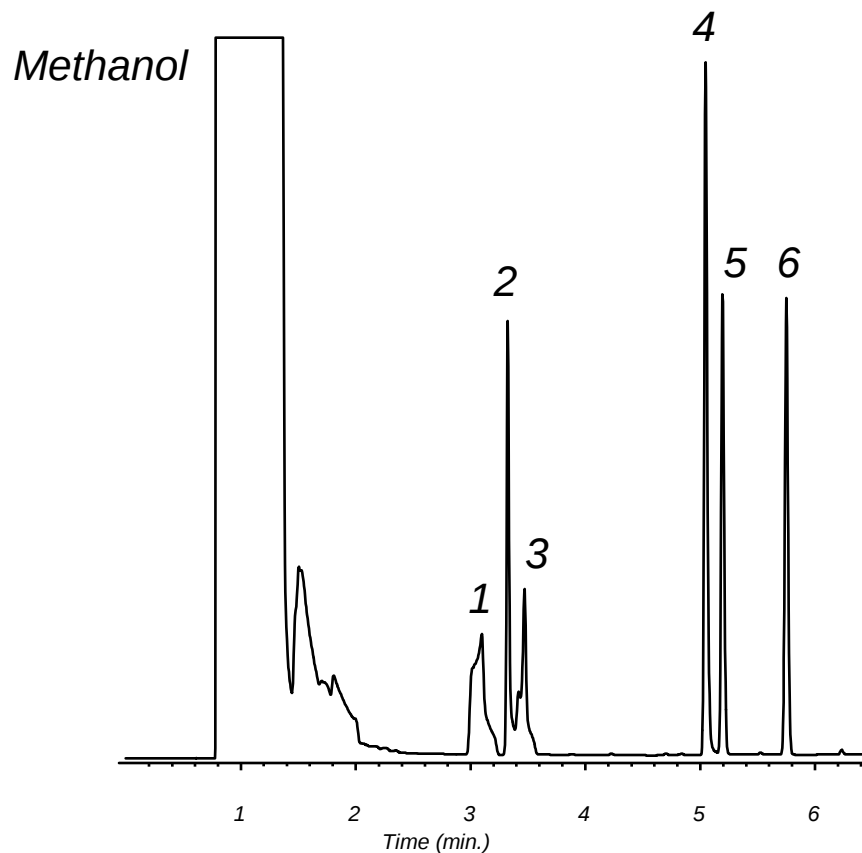
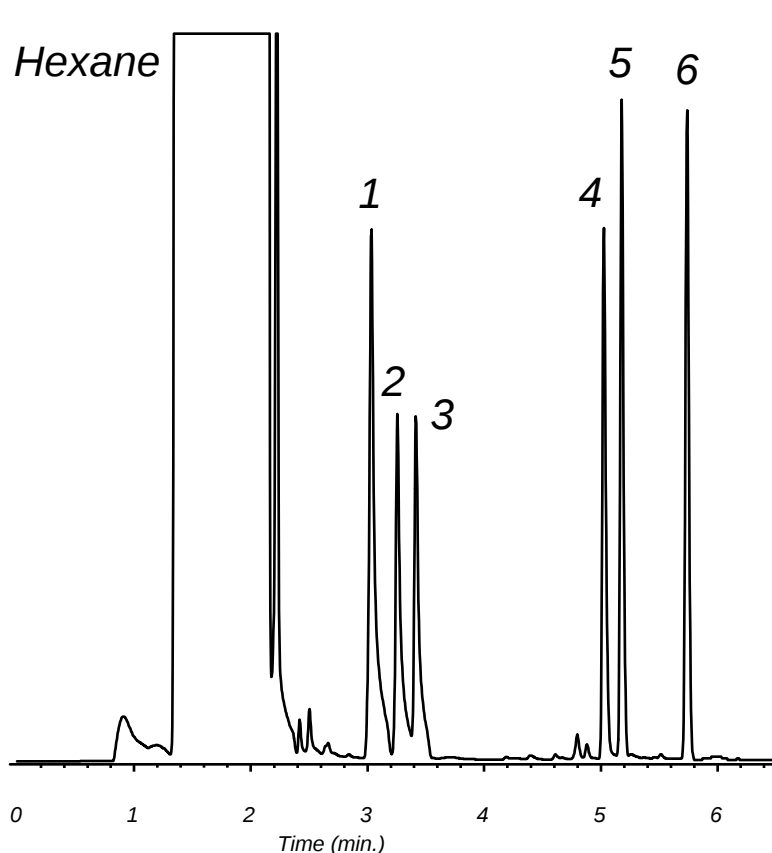
COLUMN

- Contamination
- Damaged/old stationary phase
- Reverse Solvent Effect



Splitless Injection

Reverse Solvent Effect



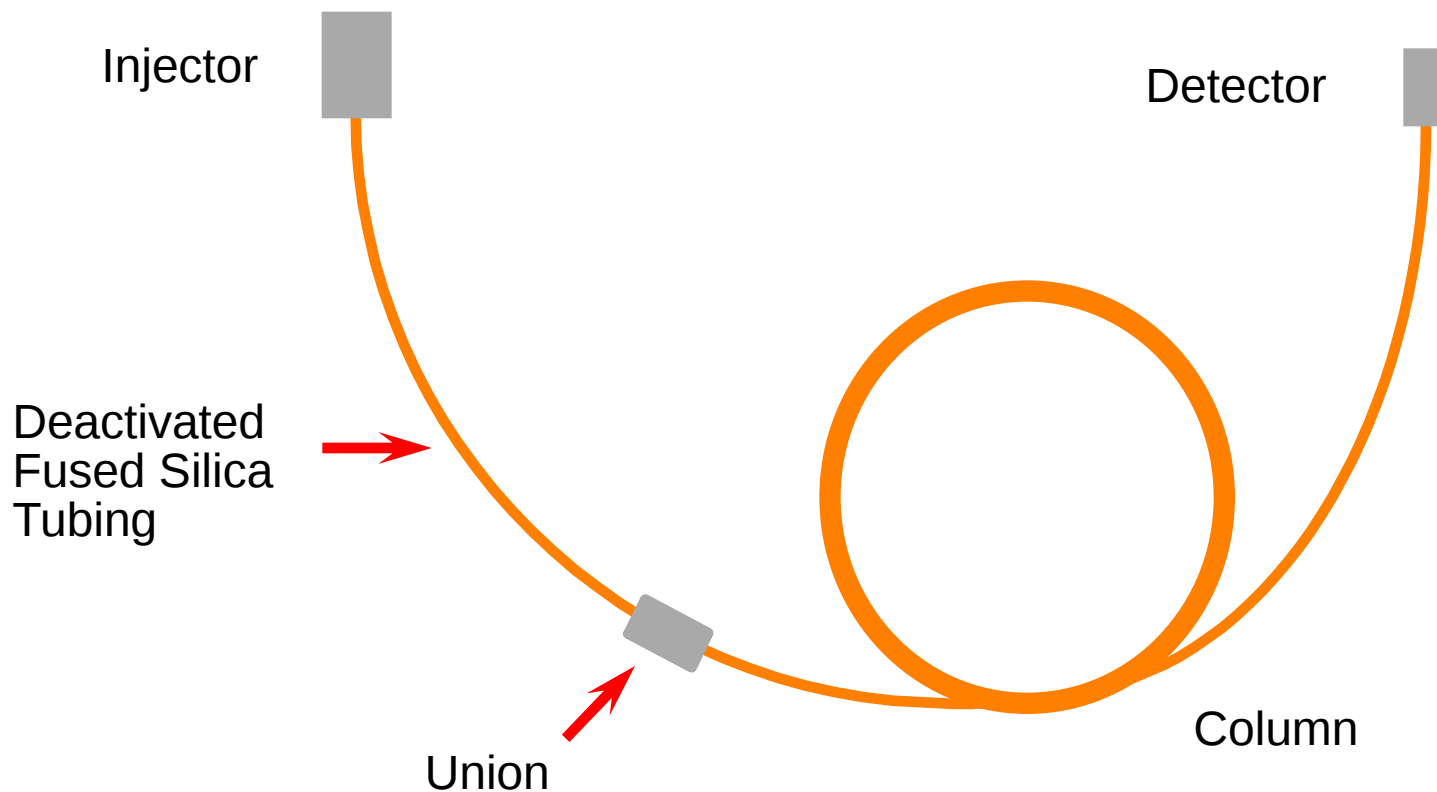
DB-1, 15 m x 0.25 mm I.D., 0.25 μ m

50°C for 1 min, 50-210°C at 20°/min; Helium at 30 cm/sec

1. 1,3-DCP 2. 3-hexanol 3. butyl acetate 4. 1-heptanol 5. 3-octanone 6. 1,2-dichlorobenzene

Retention Gap

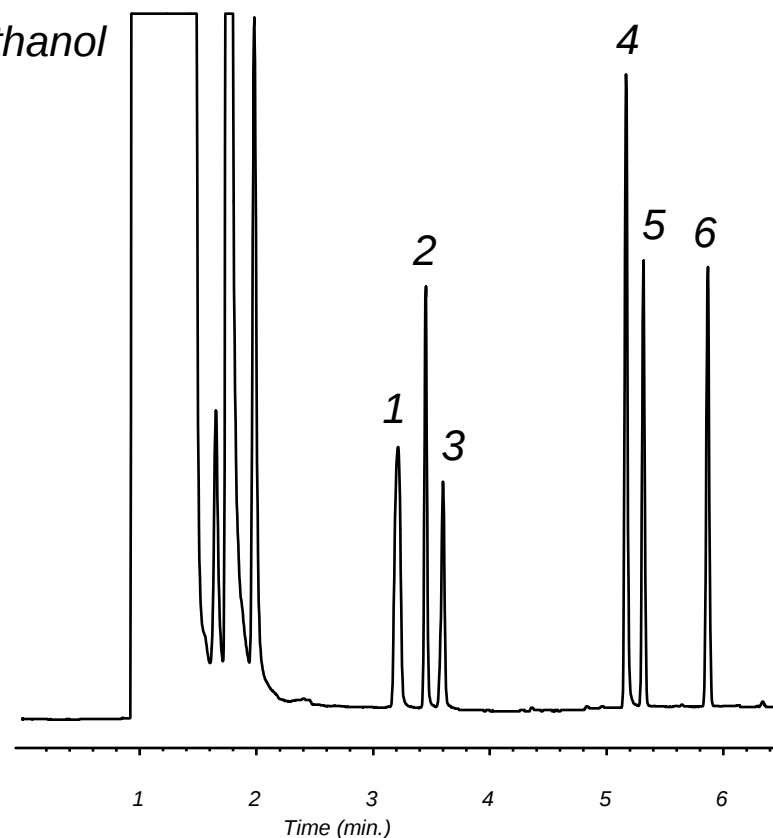
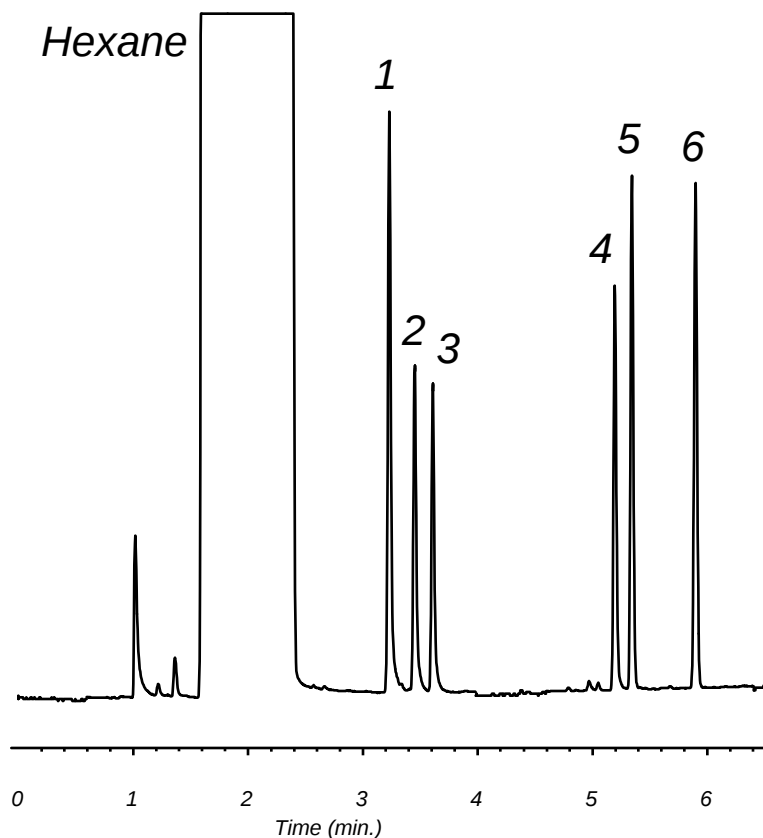
Also Called A Guard Column



Usually 2-10 meters long and same diameter as the column
(or larger if needed)

Fixing the Reverse Solvent Effect

3 m x 0.25 mm I.D. Retention Gap



DB-1, 15 m x 0.25 mm I.D., 0.25 μ m

50°C for 1 min, 50-210°C at 20°/min; Helium at 30 cm/sec

1. 1,3-DCP 2. 3-hexanol 3. butyl acetate 4. 1-heptanol 5. 3-octanone 6. 1,2-dichlorobenzene

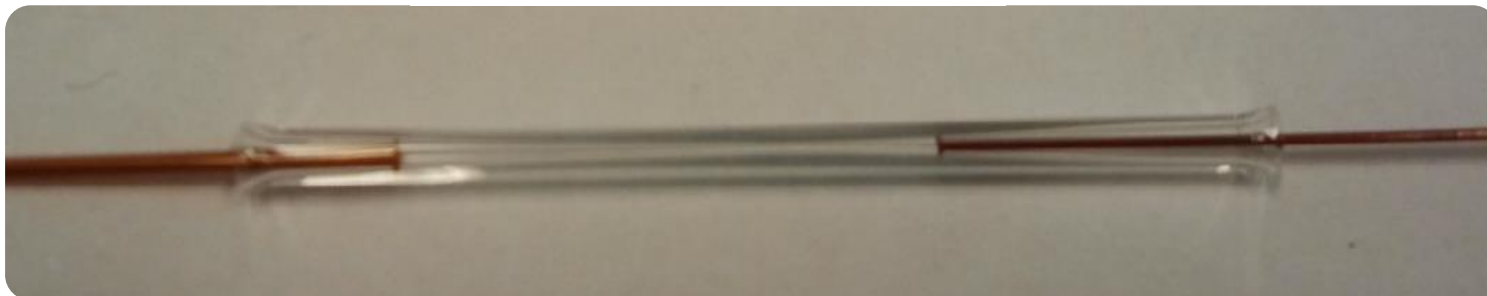


Press Fit Connectors

Competitive Press-Tight

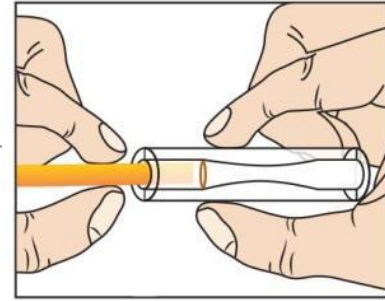


Agilent Ultra Inert Press Fit



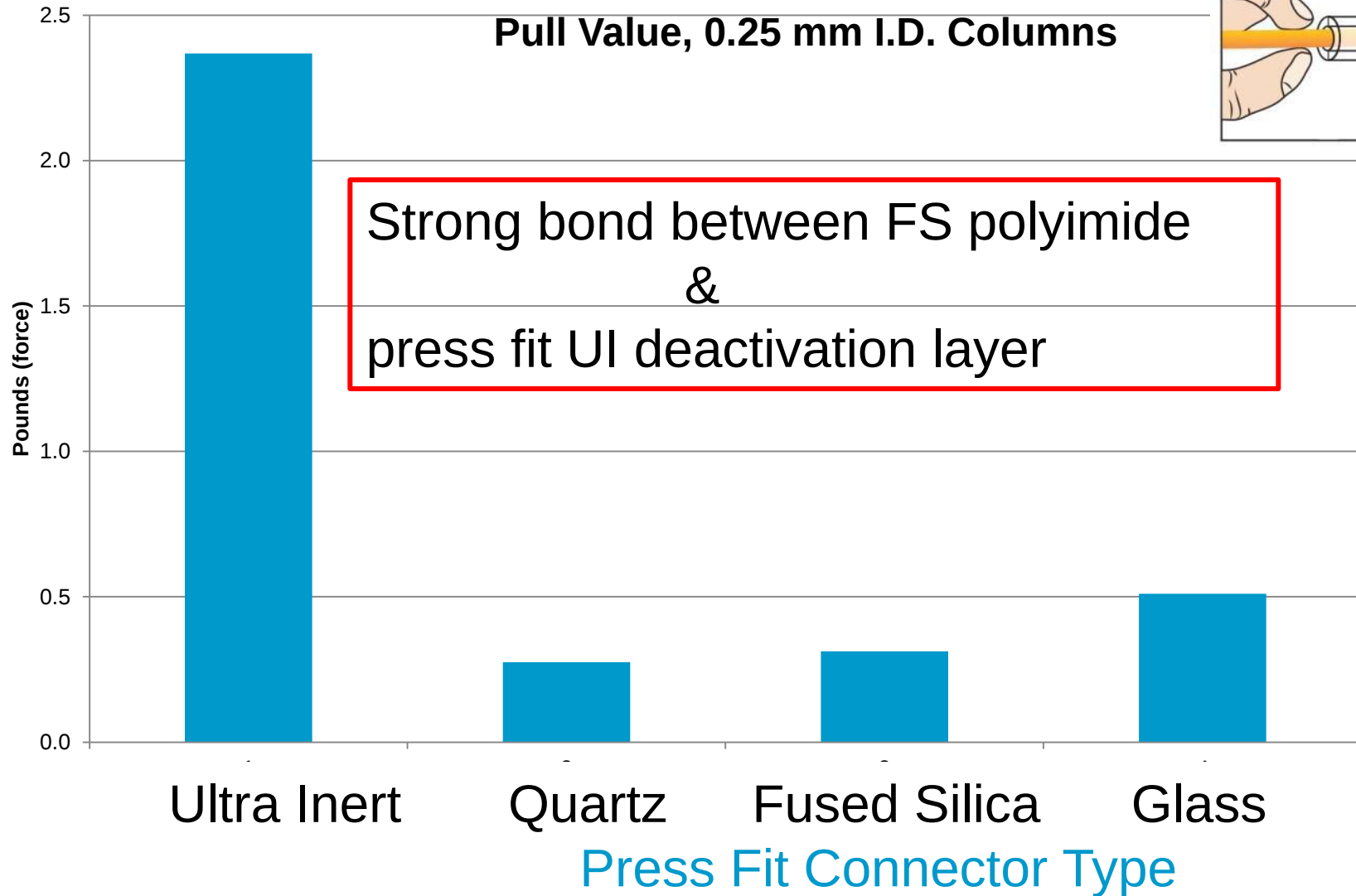
Inertness, seal and transparency are key parameters

Maximum Force Required to “pull apart”

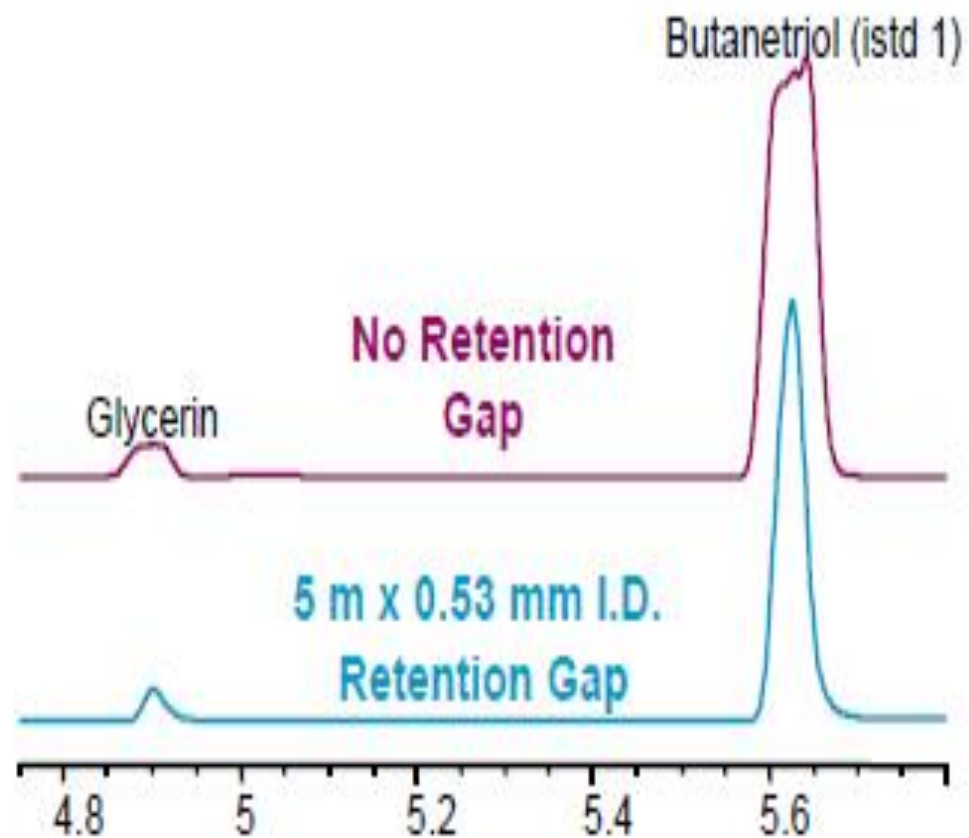


Pull Value, 0.25 mm I.D. Columns

Strong bond between FS polyimide
&
press fit UI deactivation layer



Ultimate Union and Retention Gap



Two Connection Options from Agilent

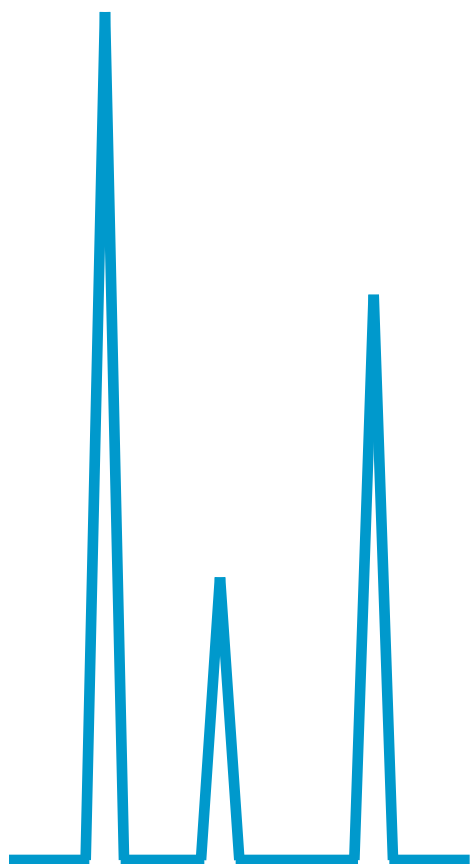


Ultra Inert Press Fit Connectors



Capillary Flow Technology

No Peaks



DETECTOR (not on or not operational)

INJECTOR (not working)

- Plugged syringe/plunger not moving

- Wrong injector (or detector)

- Huge leak/no carrier gas flow (older systems)

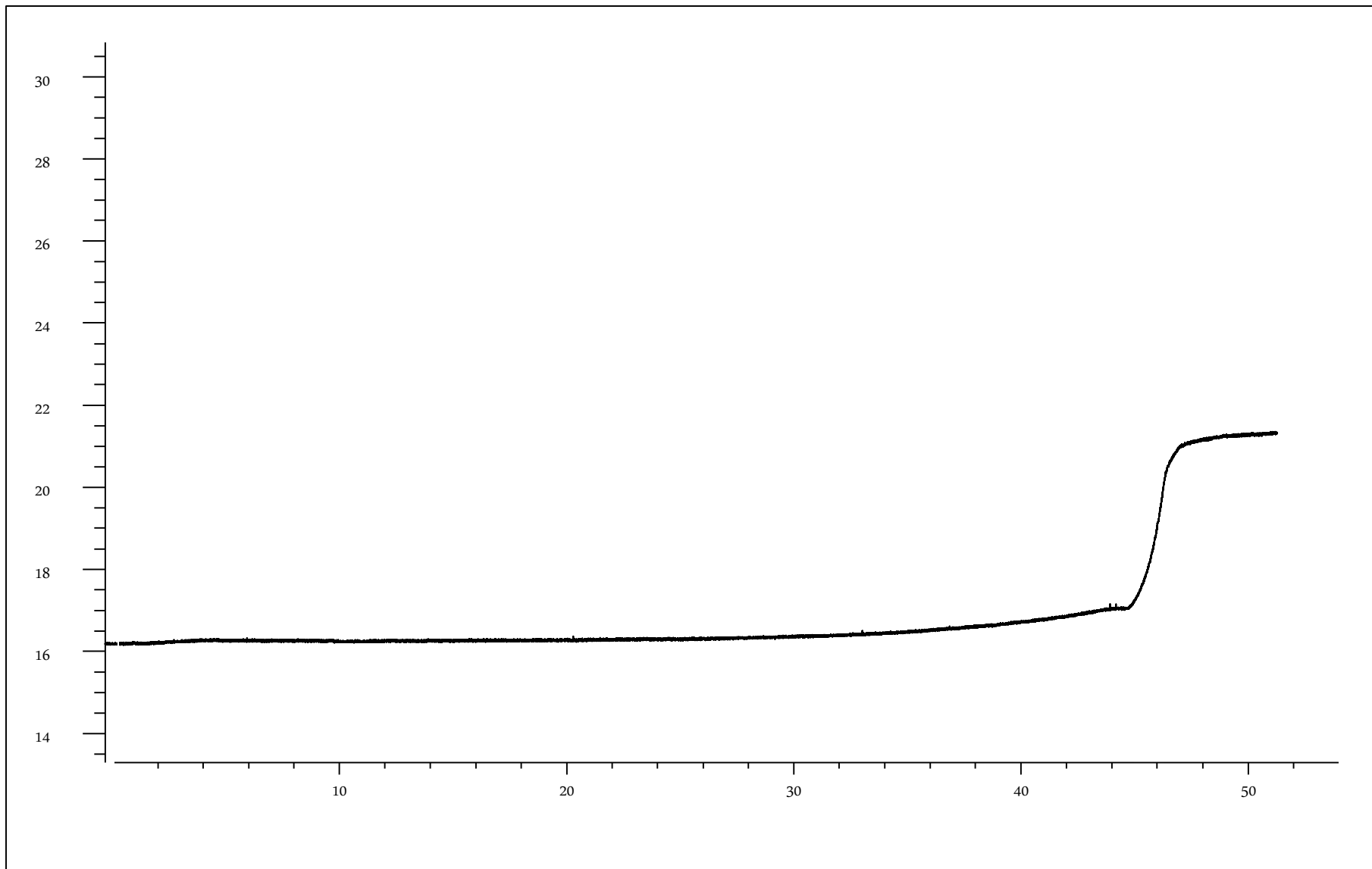
NOT the COLUMN Unless...

- Broken column

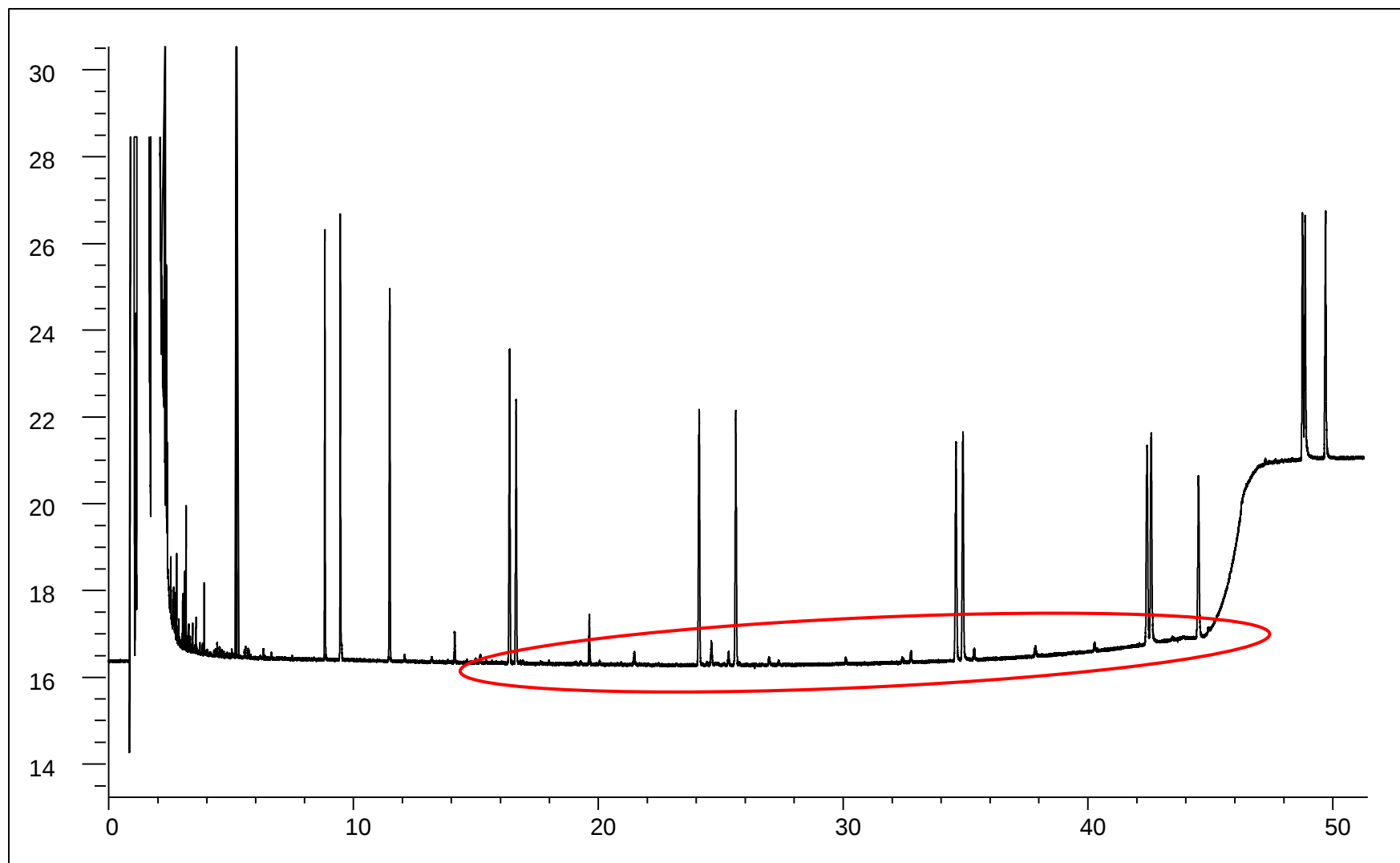
- No column



Symptom – No Peaks



Solution - Unplugged Syringe



Peak Response

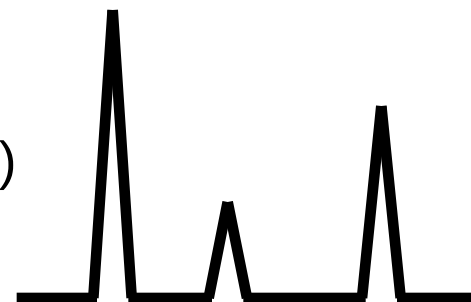
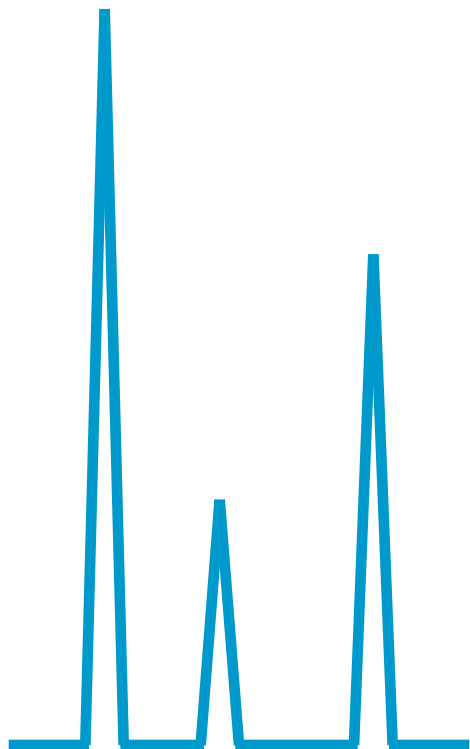
All Change in Size

INJECTOR

- Leaky syringe
- Split ratio set incorrectly
- Wrong purge activation time
- Septum purge flow too high
- Injector temperature too low*

DETECTOR (response problem)

- Settings or flows changed
- Electronics failing

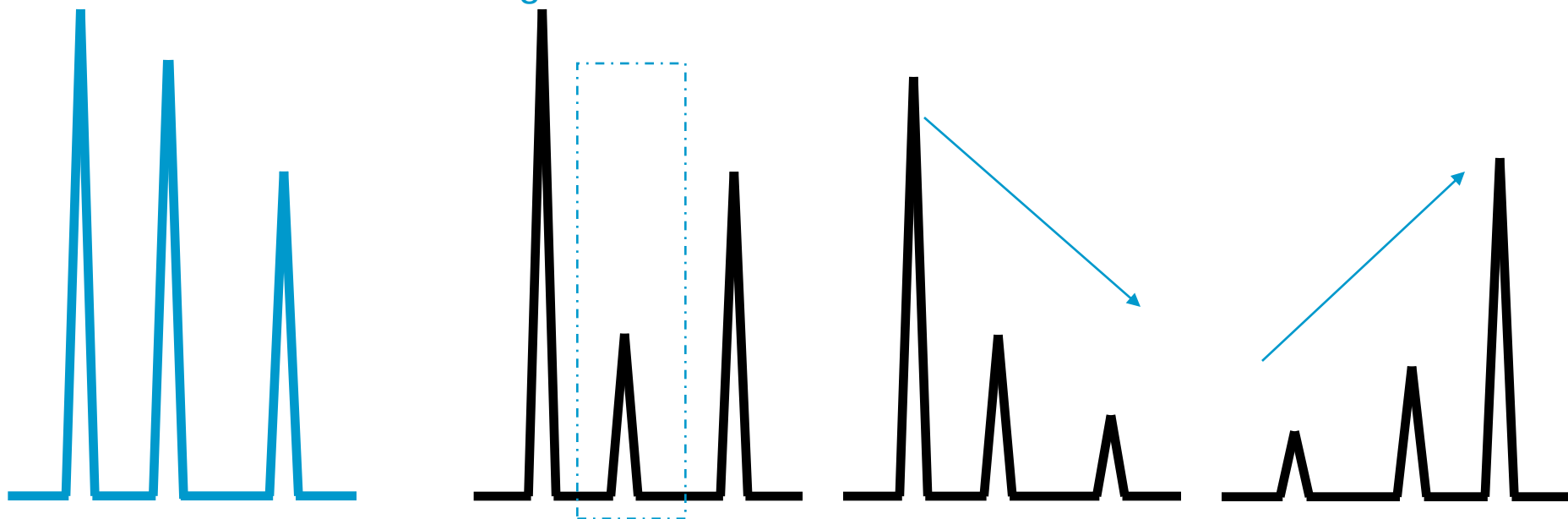


*Tip = Ask is it all of them or some of them, if all then injector or detector



Peak Response

Some Change in Size

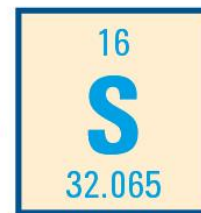
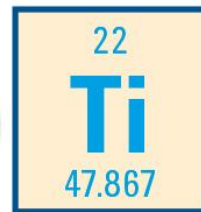


INJECTOR or COLUMN is active/contaminated

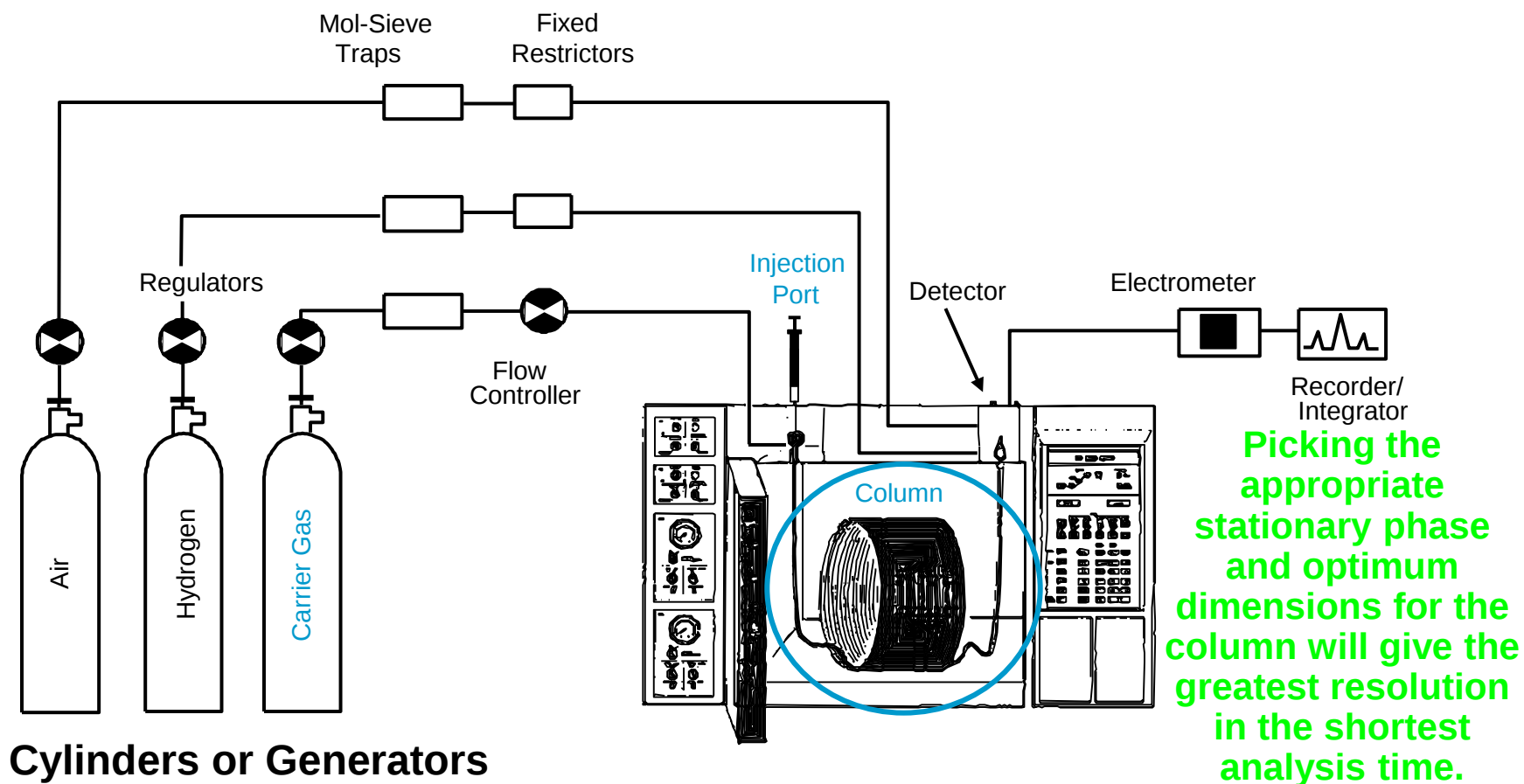
- Irreversible adsorption of active compounds (-OH, -NH, -SH)
- Decomposition of sample
- Temperature Change – Discrimination
- Evaporation from sample

*Tip = If only some change, then ask which ones? If active compounds then activity. If tracks volatility then cold spots or inlet discrimination.

Advanced
Troubleshooting
Part 2 The Rest of the Story



Typical Gas Chromatographic System



Peak Fronting

Shark Fin Shaped or Just Slight

COLUMN (contaminated)

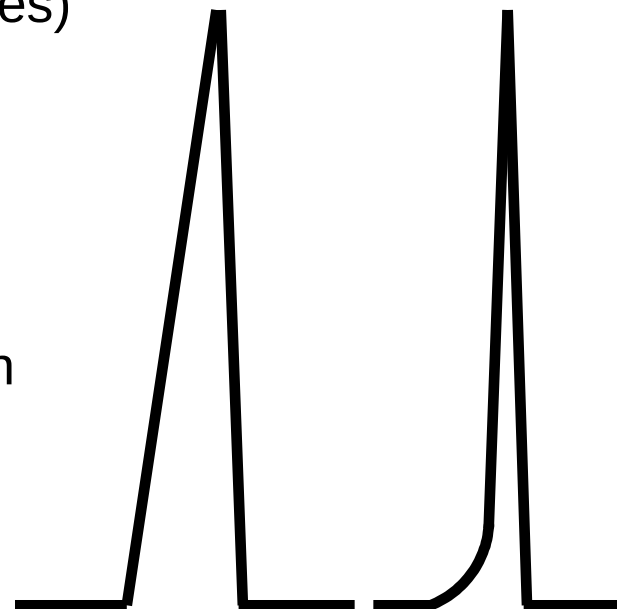
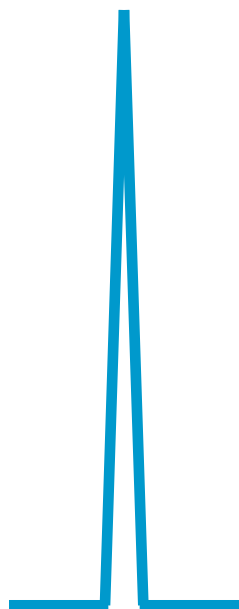
- Overload (More pronounced with large solute and phase polarity differences)

INJECTOR

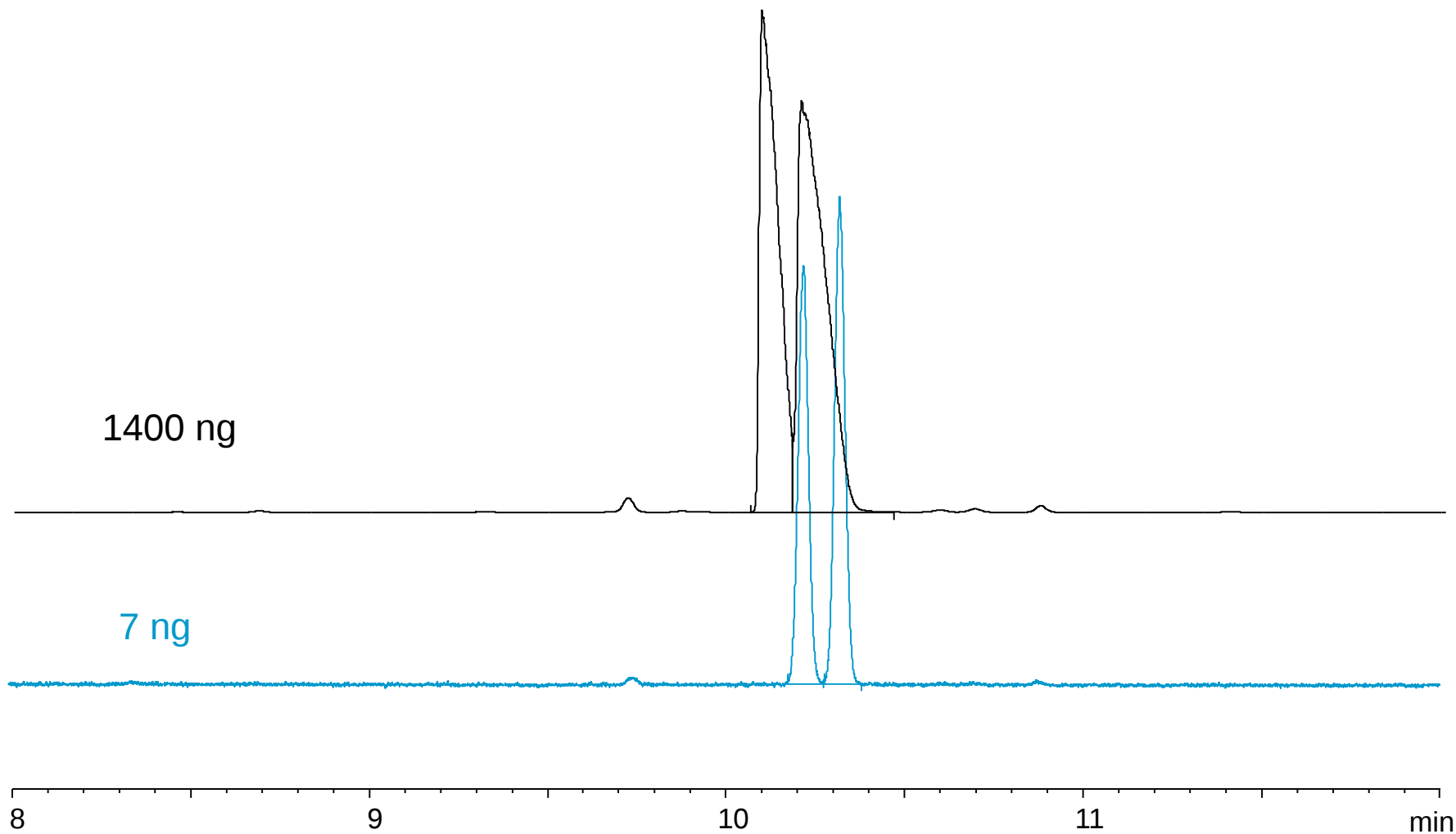
- Poor efficiency (flow/temp)
- Column installation
- Compound very soluble in injection solvent (need retention gap)
- Mixed sample solvent

OTHER

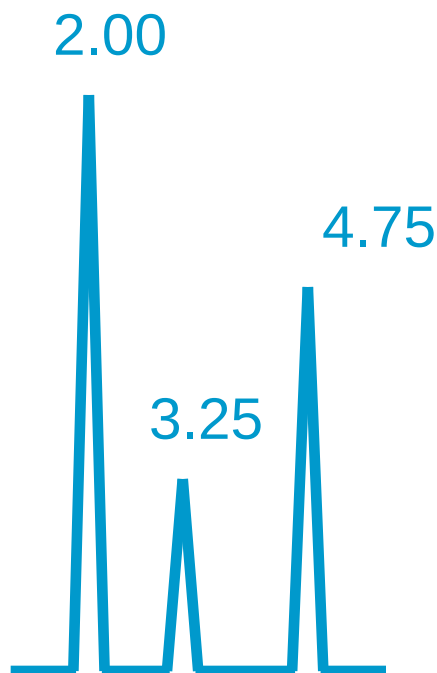
- Co-elution
- Breakdown



Effect of Sample Overload on Retention Time and Peak Shape



Retention Time Shift



INJECTOR

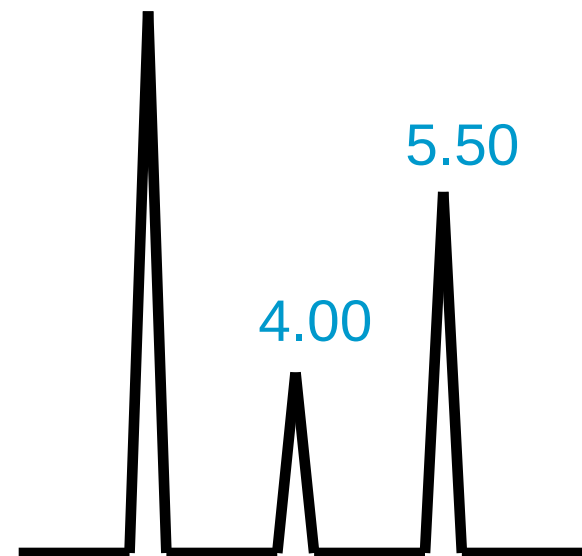
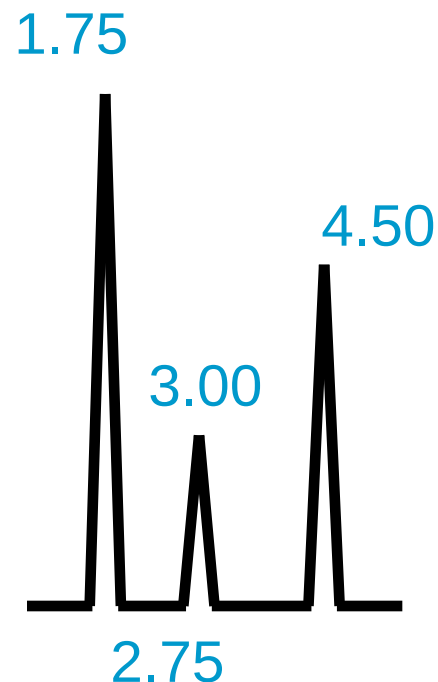
- Change in injection solvent
- Large change in sample concentration

FLOW

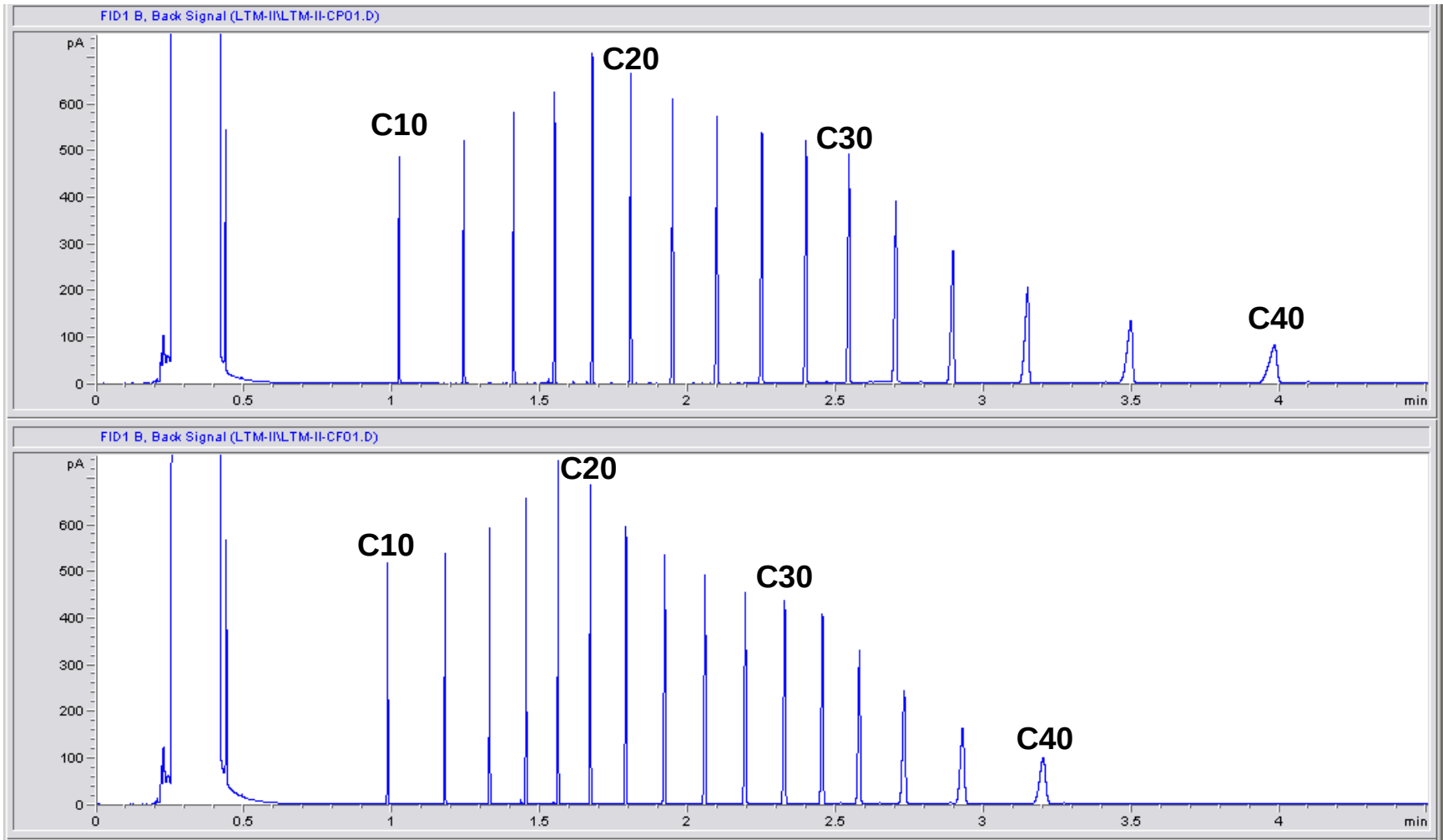
- Leak in the septum
- Change in gas velocity

COLUMN

- Contamination
- Damaged stationary phase
- Loss of stationary phase
- Change in temperature

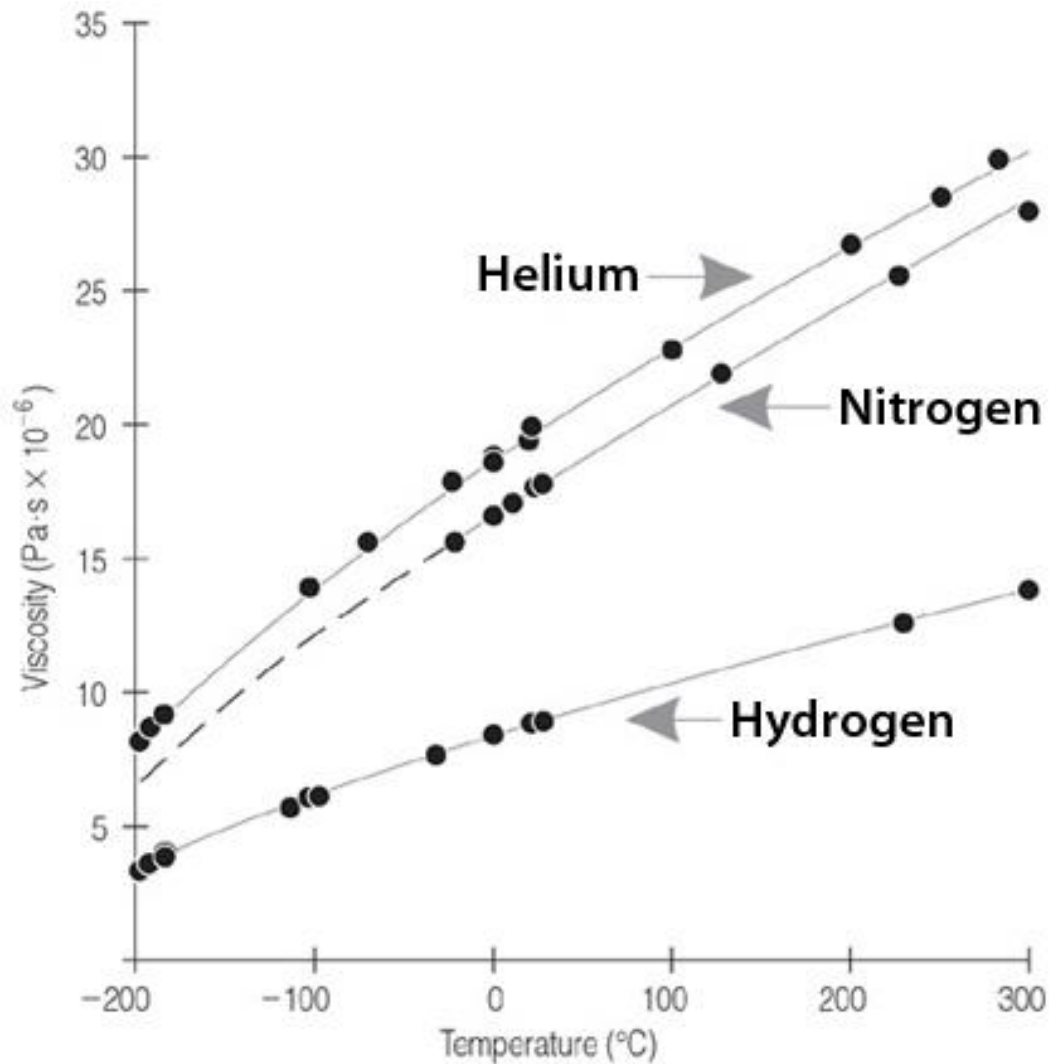


Constant Pressure vs Constant Flow Retention



Under constant pressure conditions, flow decreases as temperature increases.
(viscosity of a gas increases as temperature increases)

Gas Viscosity vs Temperature



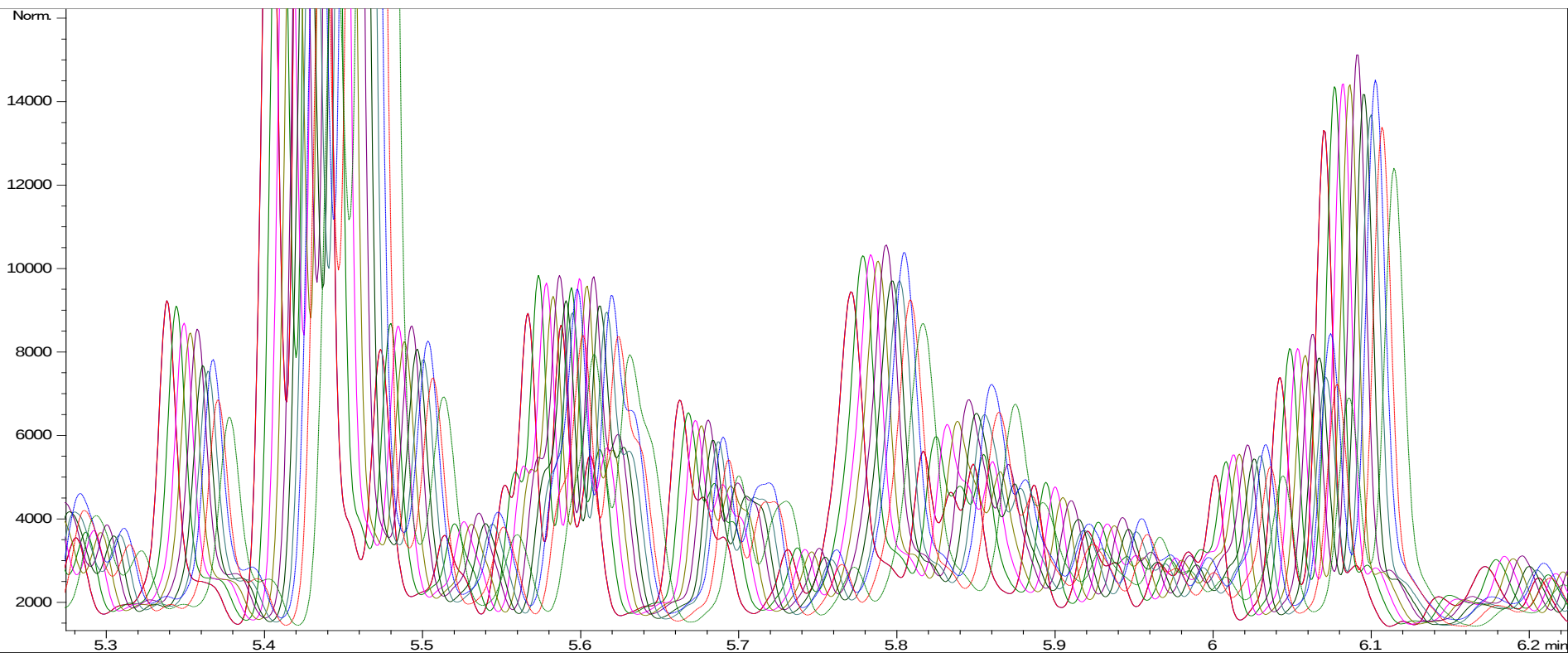
J.V. Hinshaw, Column Connections, LCGC Asia Pacific, 12(2), 1100 (2009).



Agilent Technologies

Retention Times Shift 4-5 sec During the First 10 Runs

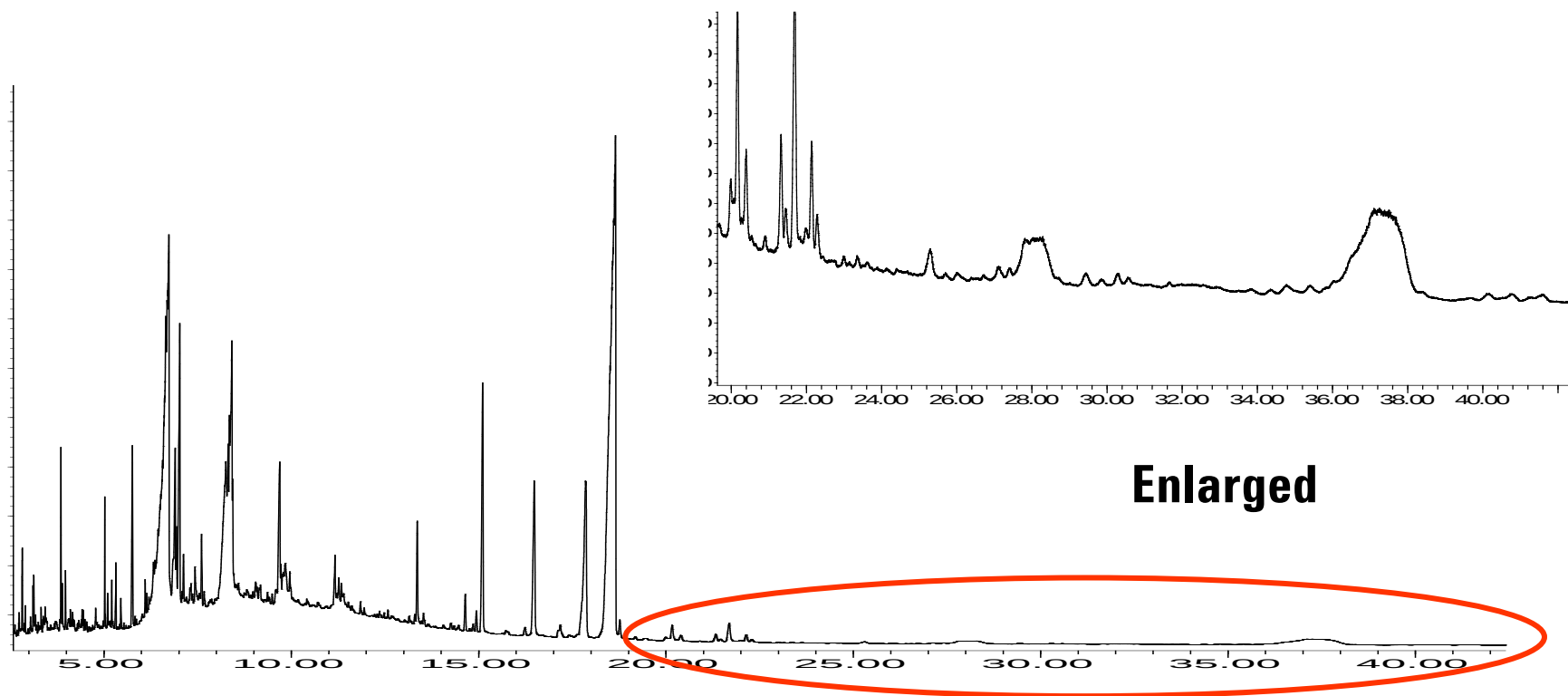
Overlay of first 10 runs



GC-FID Chromatogram of Fish Oil

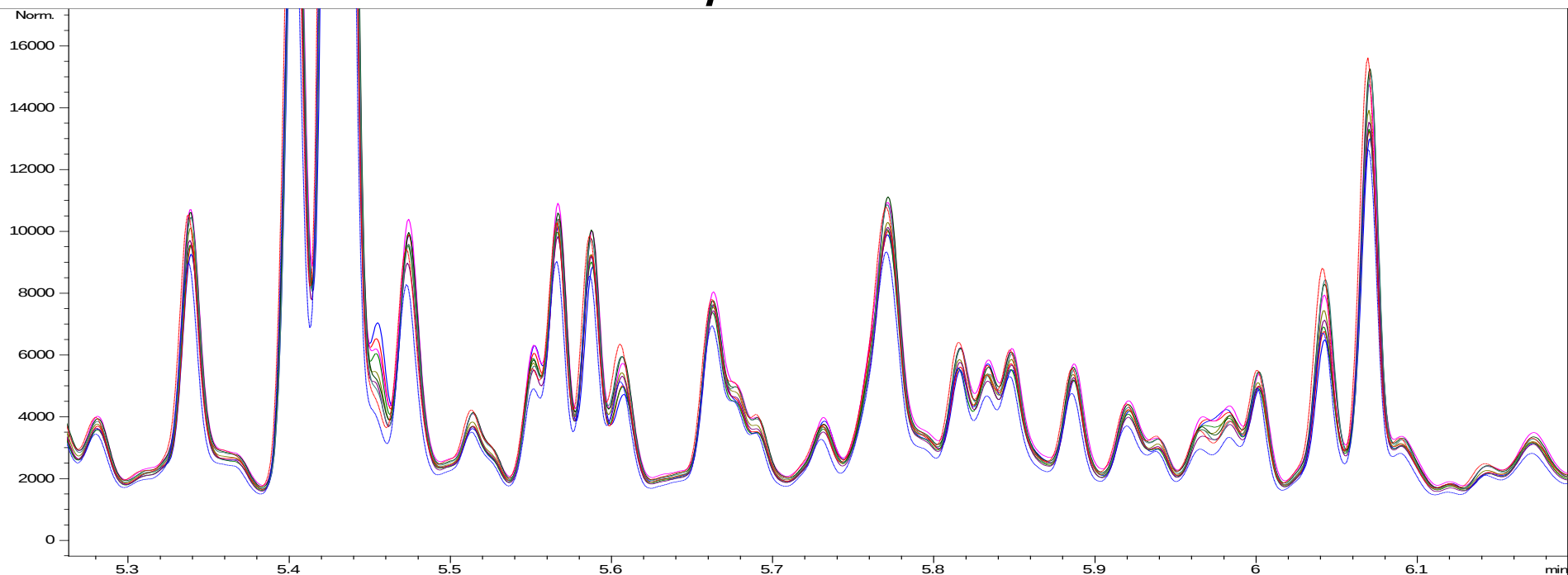
1- μ L splitless injection of 10% fish oil

GC oven was held at 290 °C for 30 min at the end of the run.



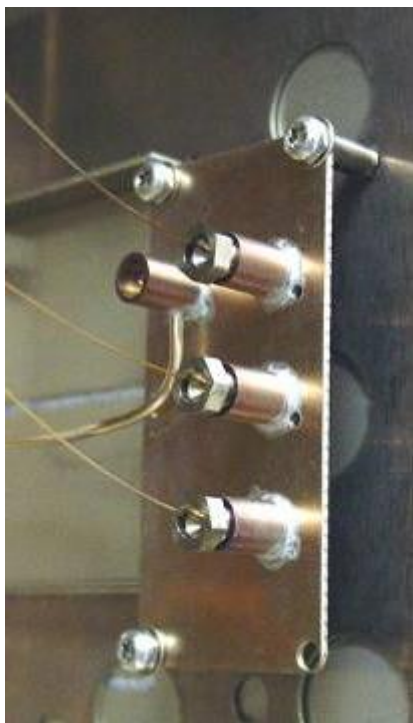
Get High M.W. Fish Oil Off the Column After Every Run = No More Shifting Retention

Retention Time Precision with Backflushing Overlay of 10 Runs

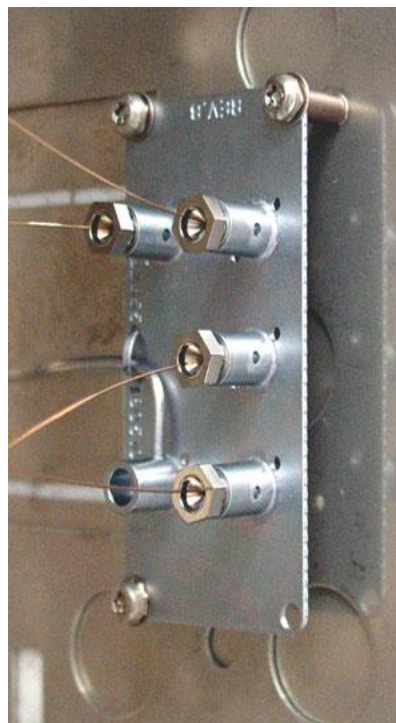


Purged Capillary Flow Devices

**2-Way Splitter
with Makeup**



**3-Way Splitter
with Makeup**



Deans Switch

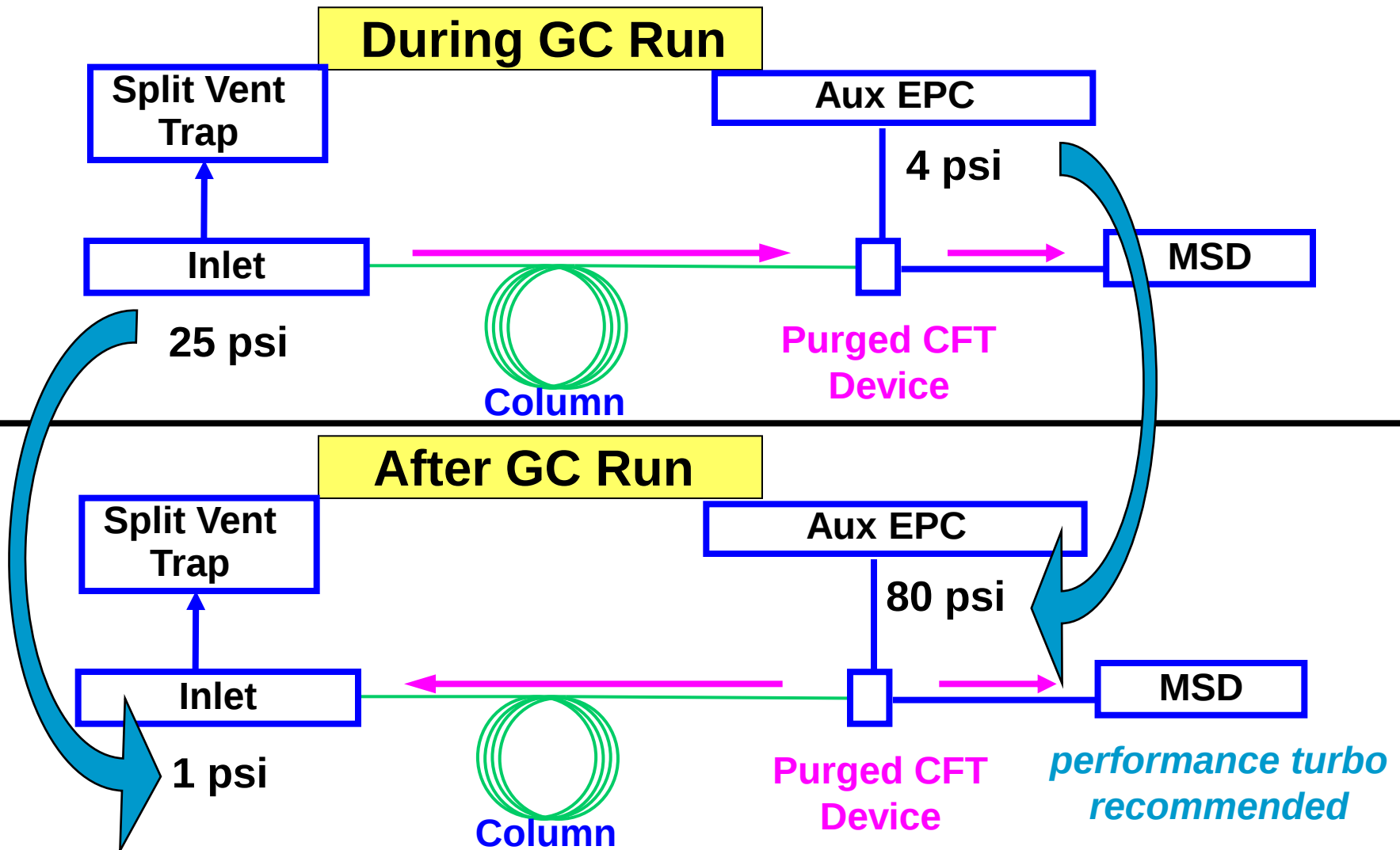


**Purged Union
(most recent)**



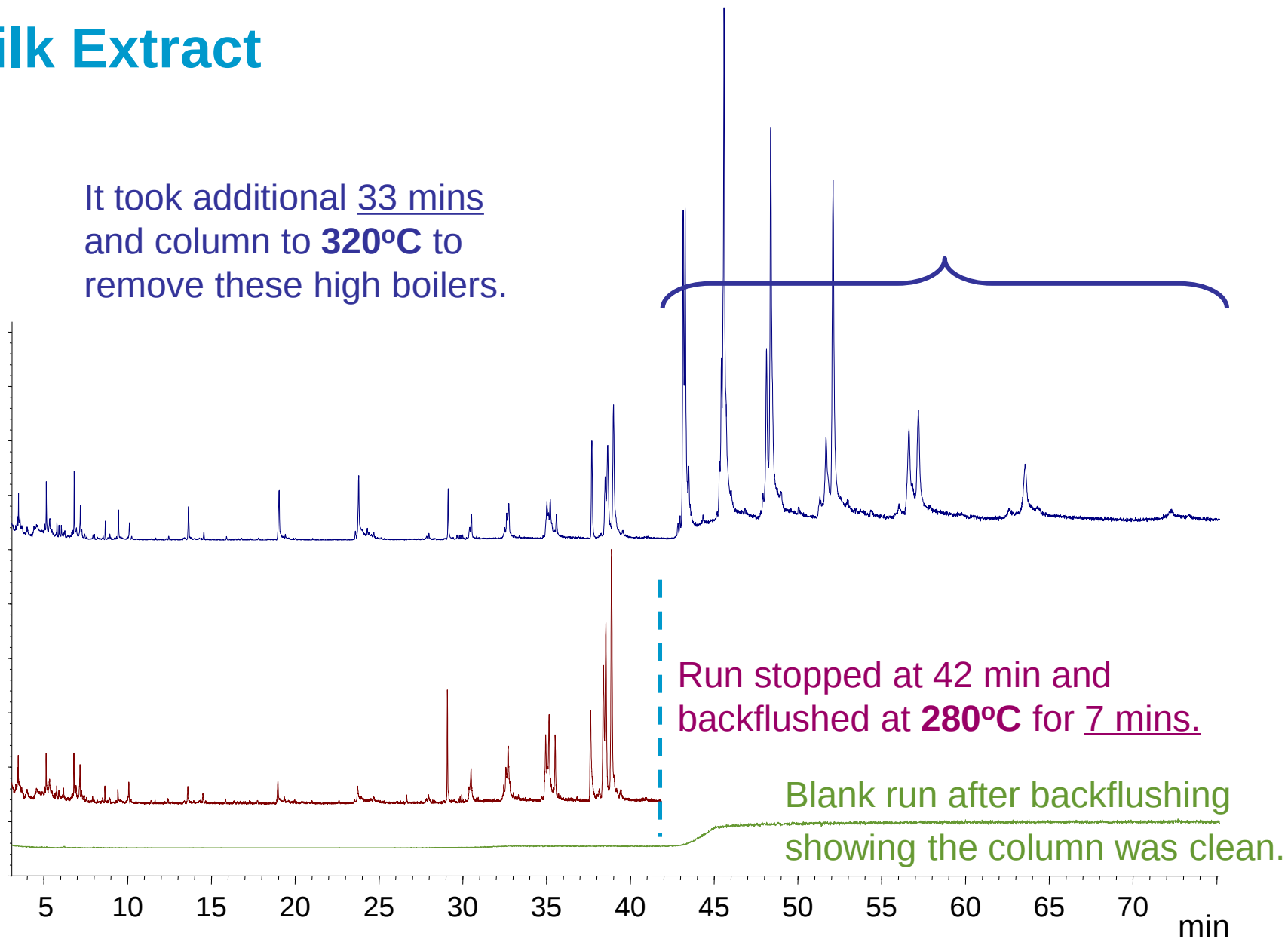
All Purged Capillary Flow Devices are Capable of Backflushing

Backflush with Purged Union or Any Purged Capillary Flow Technology Device

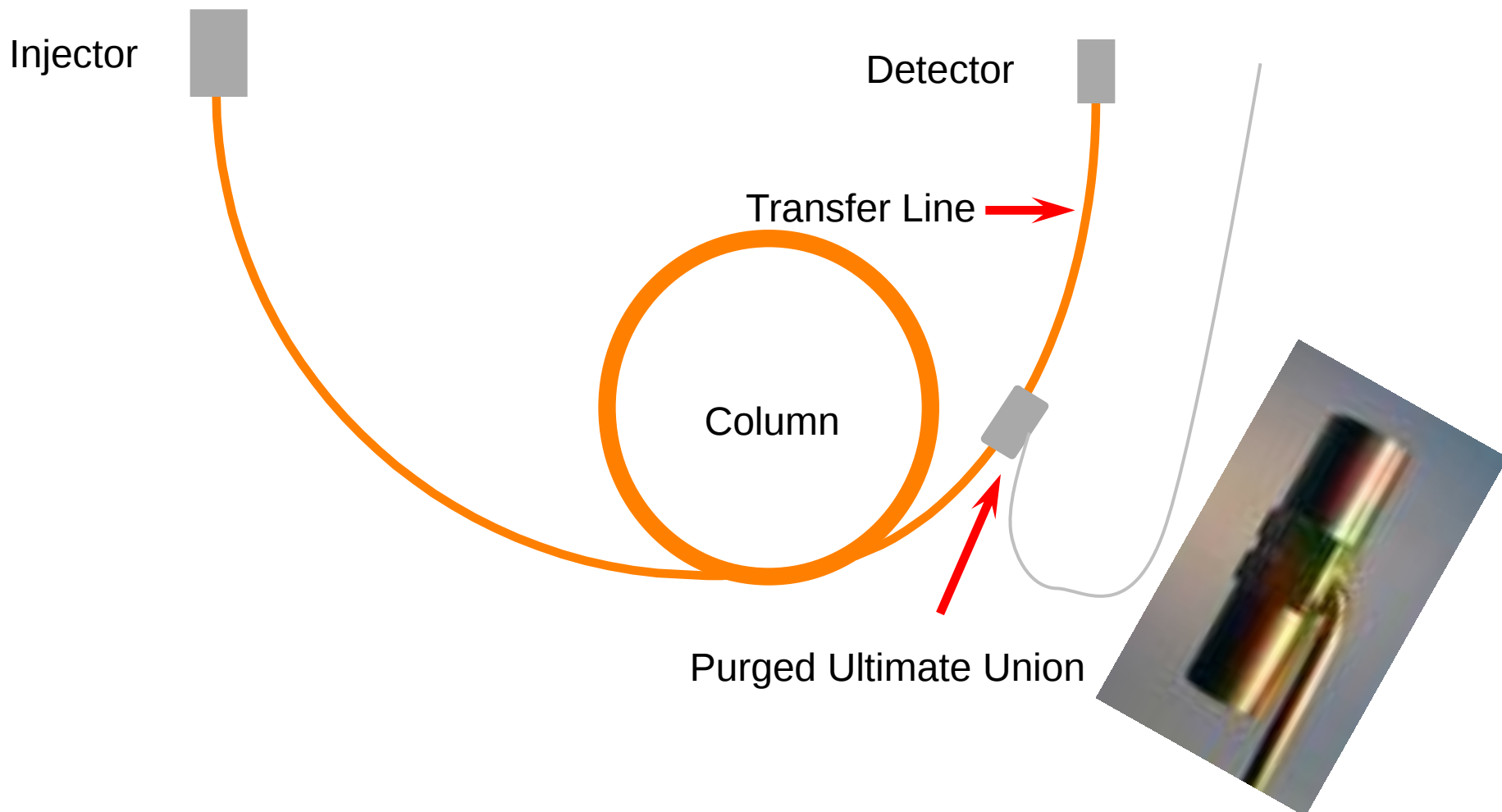


Milk Extract

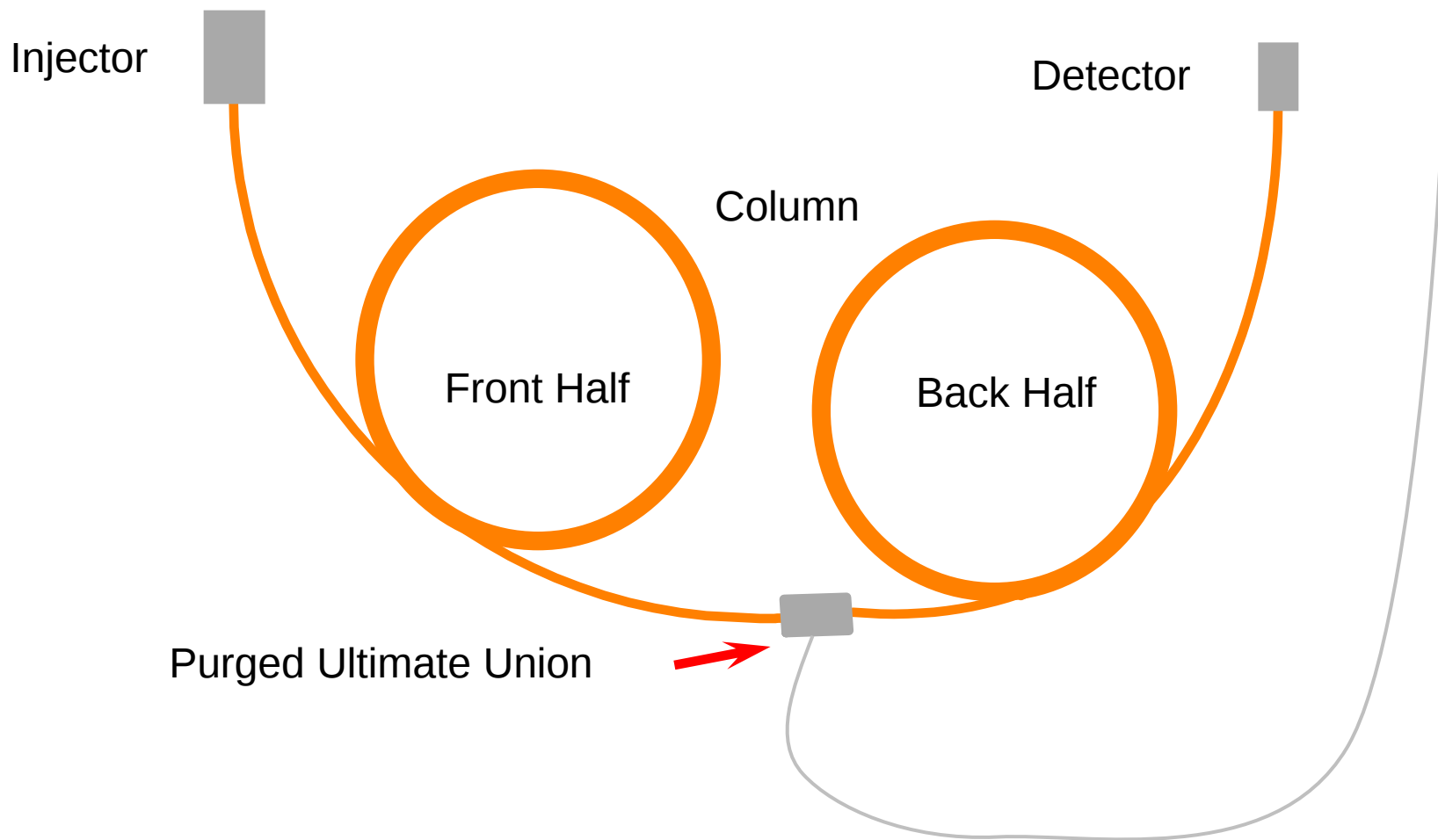
It took additional 33 mins
and column to **320°C** to
remove these high boilers.



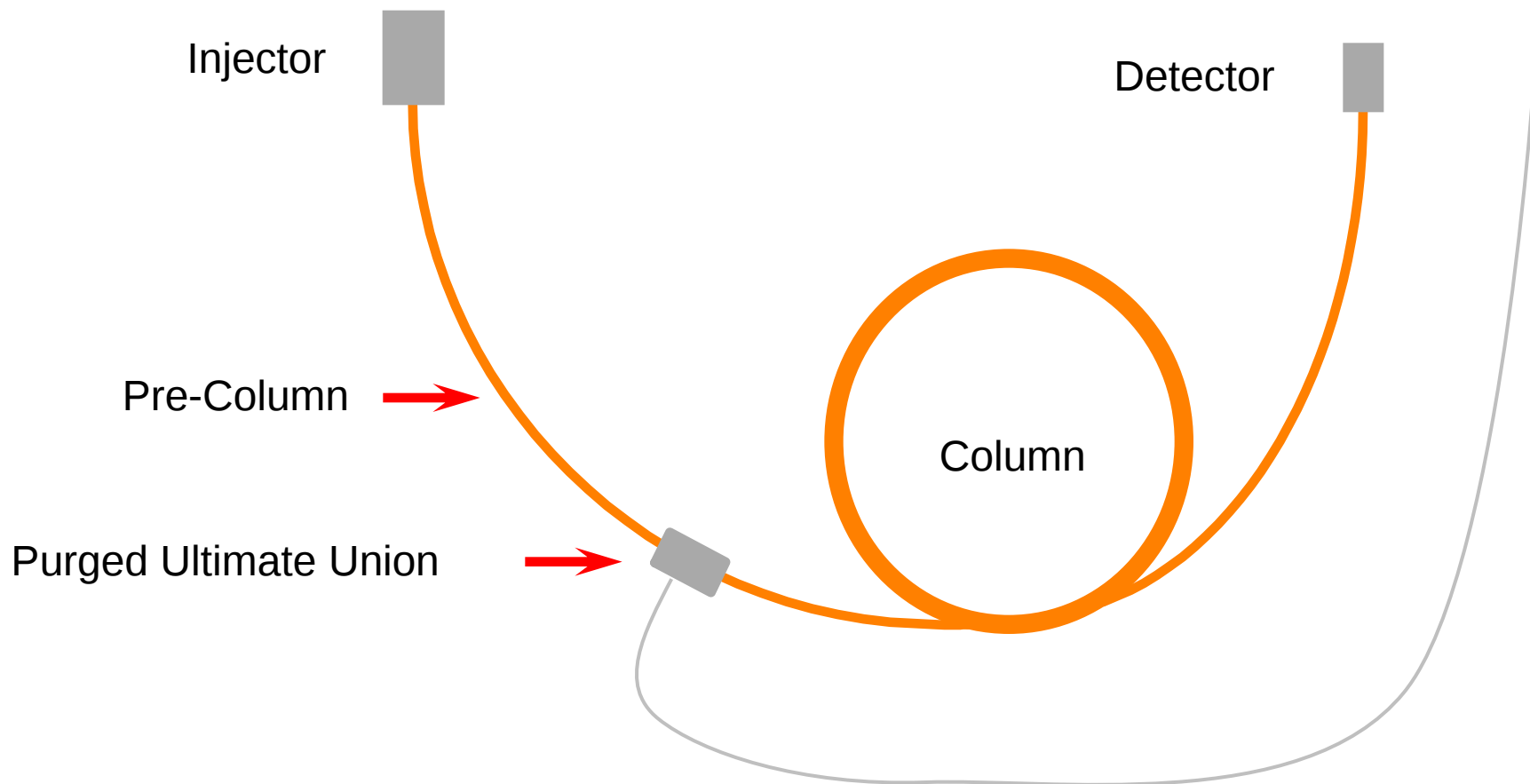
Post-Column Backflush



Mid-Column Backflush



Pre-Column Backflush



Without Backflush: A Serious Problem

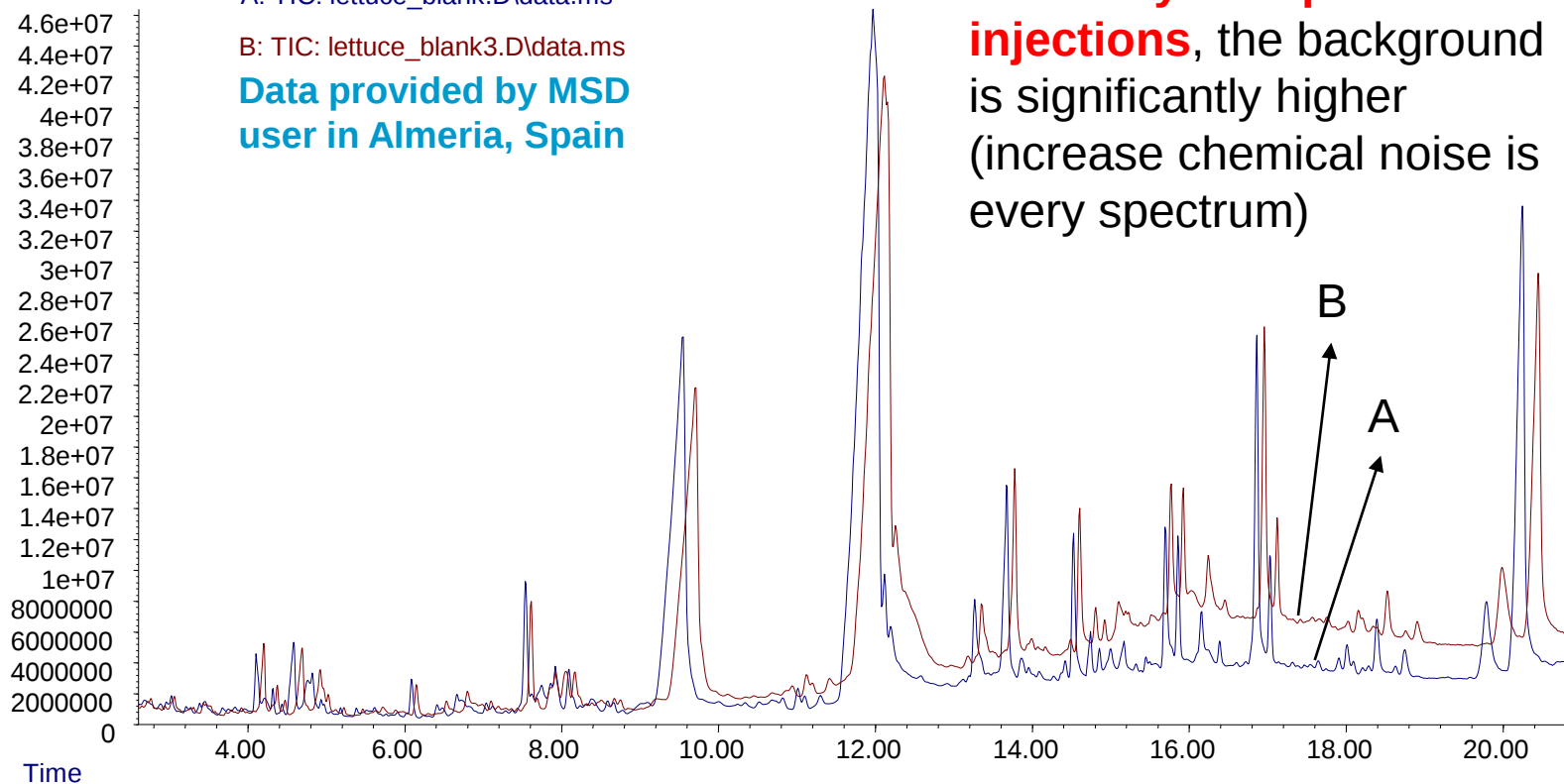
Abundance

A: TIC: lettuce_blank.D\data.ms

B: TIC: lettuce_blank3.D\data.ms

Data provided by MSD
user in Almeria, Spain

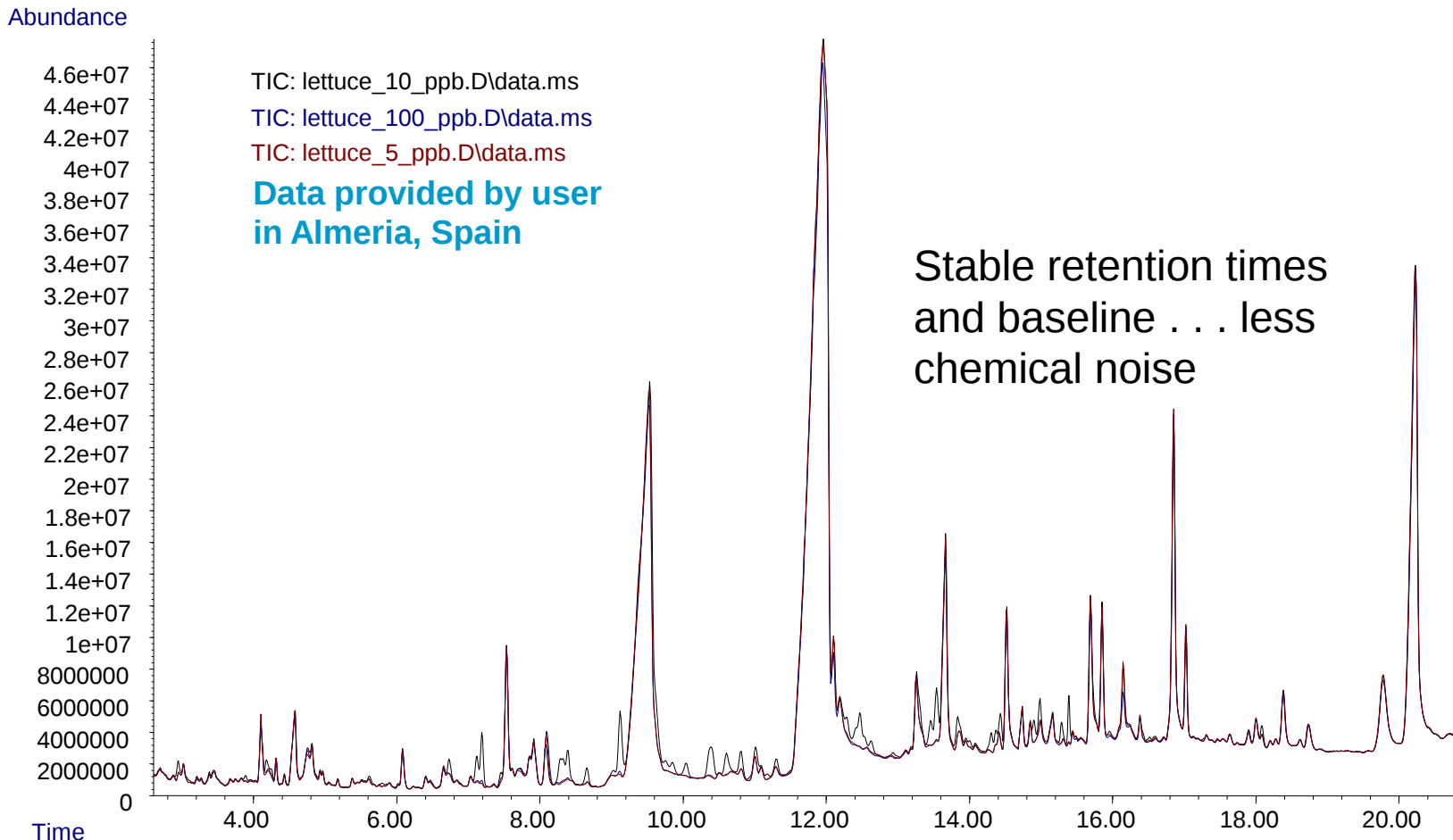
After **only 3 10- μ L injections**, the background is significantly higher (increase chemical noise is every spectrum)



Overlay of two chromatograms of a blank extract injected BEFORE (A) and AFTER (B) three injections without backflush

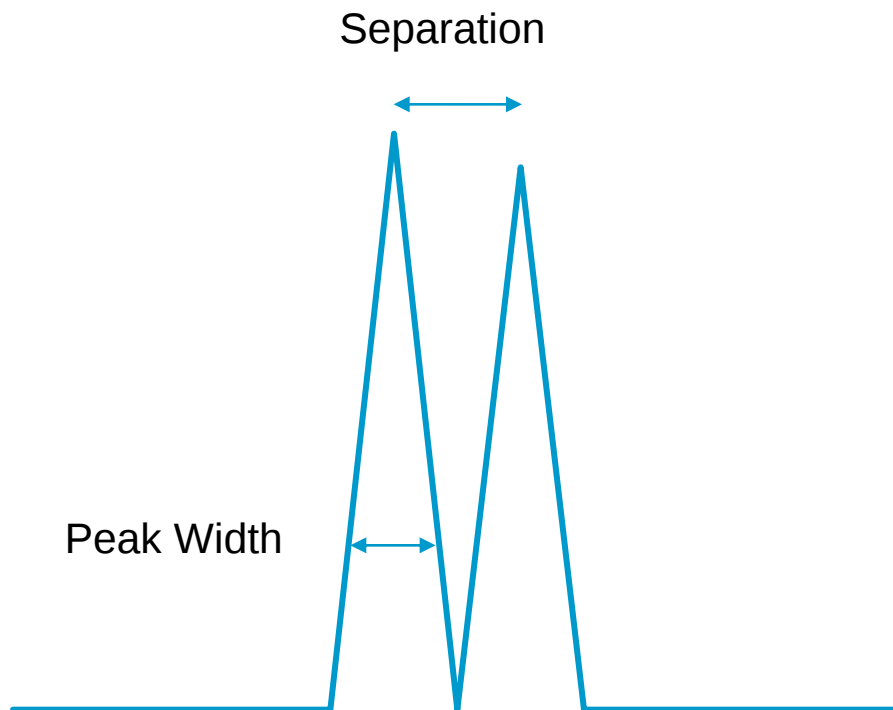


With Backflush: No Increased Background (Less Spectral Noise) and Consistent Retention Times



Overlay of three chromatograms of lettuce extract run with 2 min of back flush

Loss of Resolution

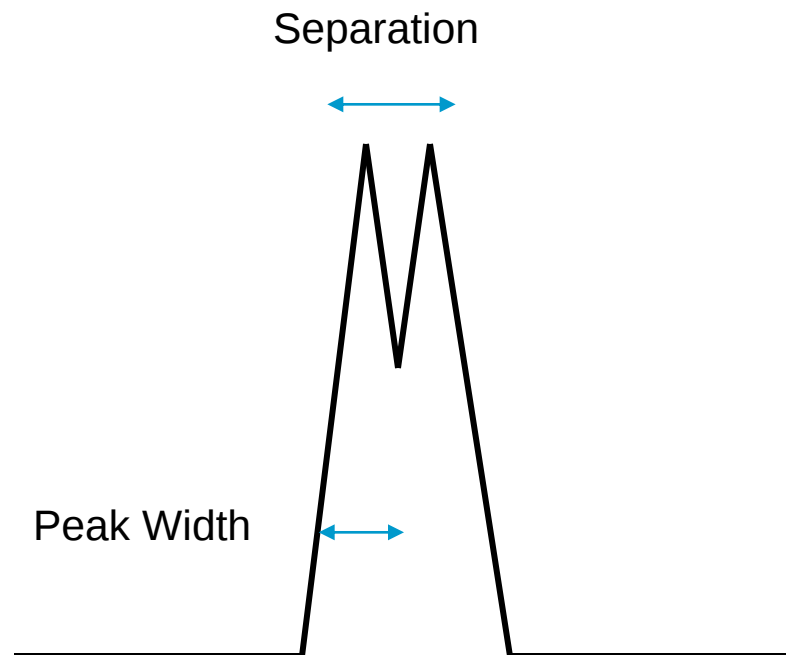


Resolution is a function of separation and peak width

Loss of Resolution - Separation Decrease

COLUMN

- Different column temperature
- Contamination (more phase?)
- Matrix components co-eluting
- Different column phase?



Loss of Resolution - Peak Broadening

FLOW

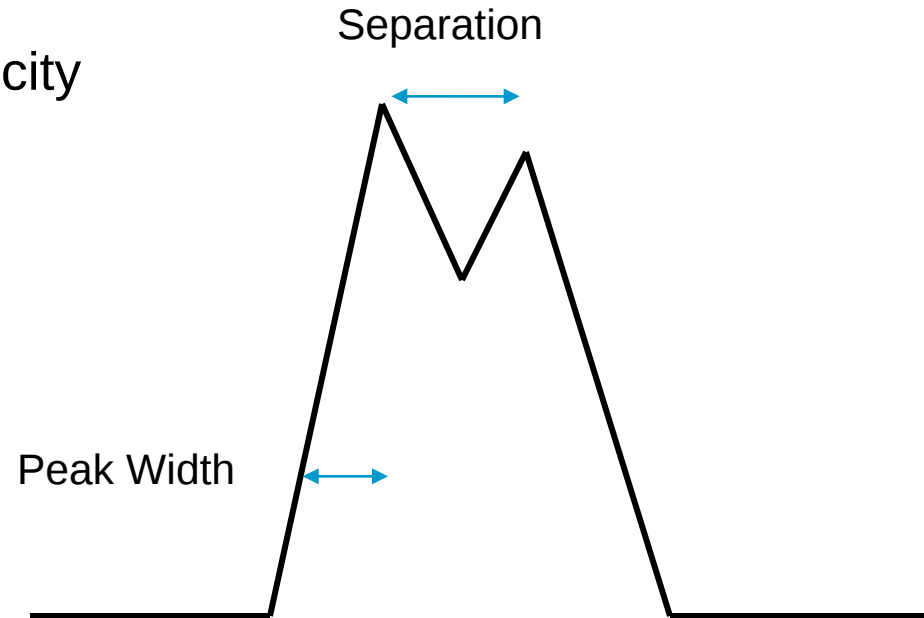
- Change in carrier gas velocity
- Make-up gas

COLUMN

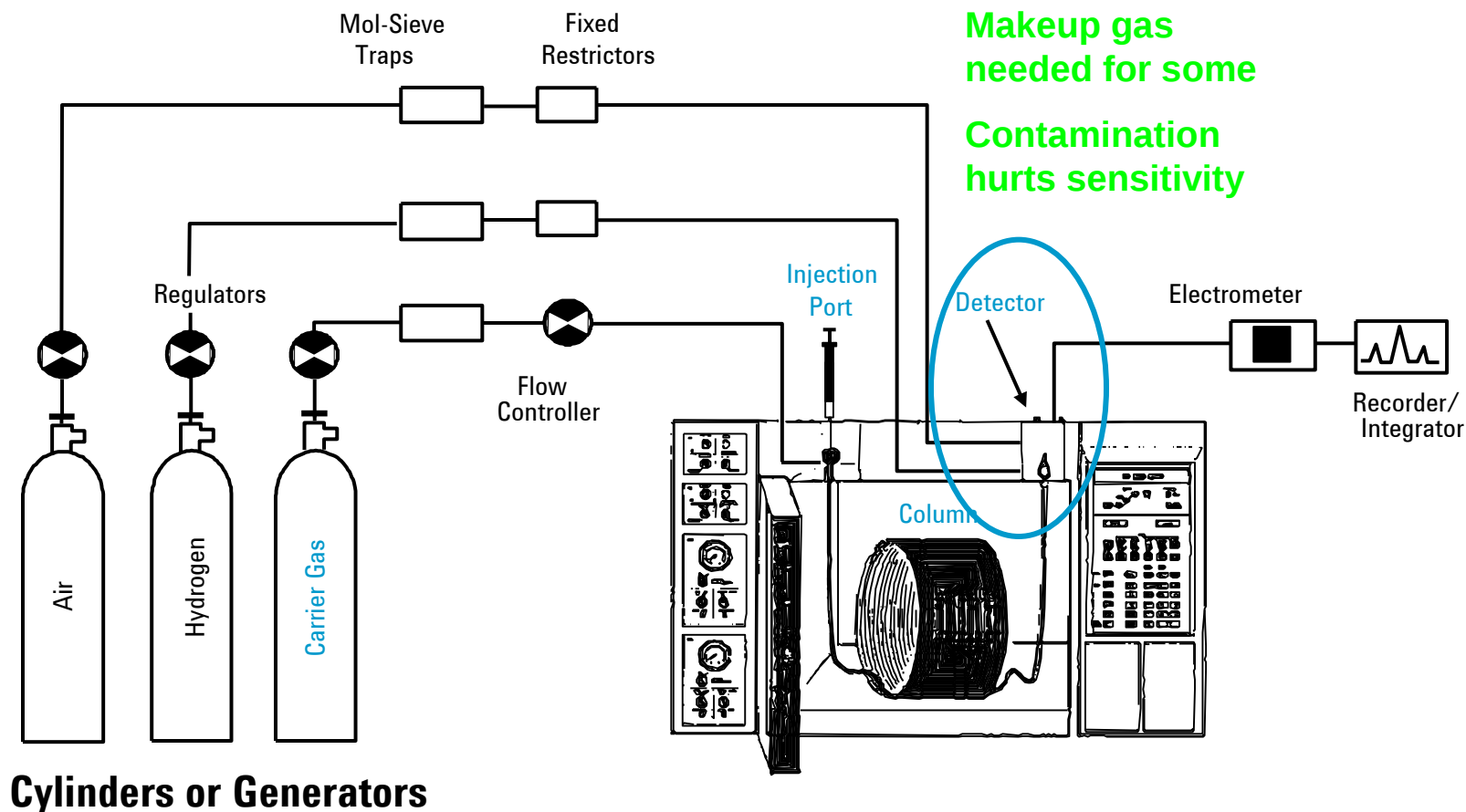
- Contamination
- Phase degradation

INJECTOR (efficiency)

- Settings, Liner, Installation, etc.



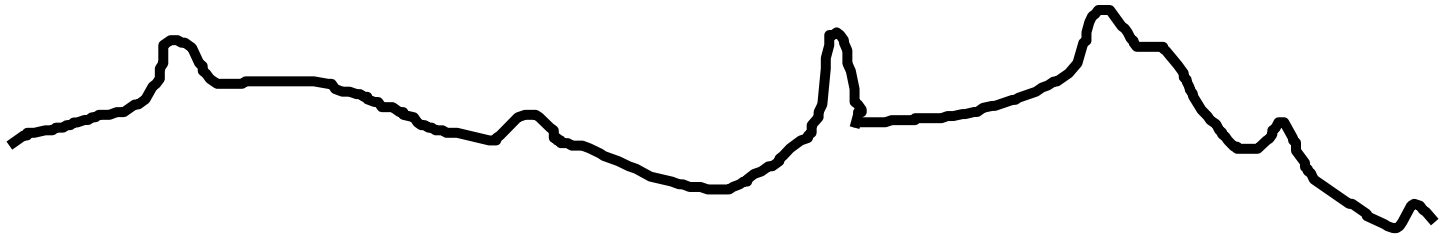
Typical Gas Chromatographic System



Baseline Disturbances

Sudden Changes, Wandering, or Drifting

WANDER



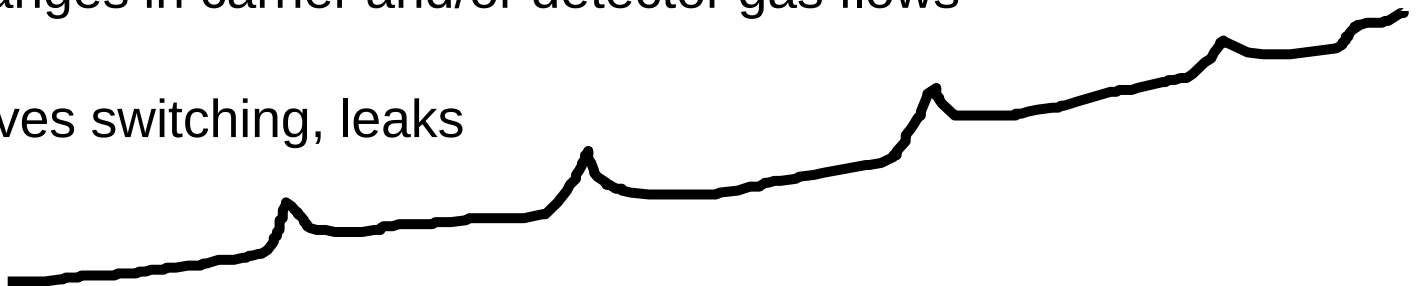
COLUMN or DETECTOR

- Not fully conditioned or stabilized (electronics)
- Contamination

FLOW

- Changes in carrier and/or detector gas flows
- Valves switching, leaks

DRIFT



Noisy Baseline

MILD



SEVERE



FLOW

- Contaminated gas
- Incorrect detector settings

COLUMN

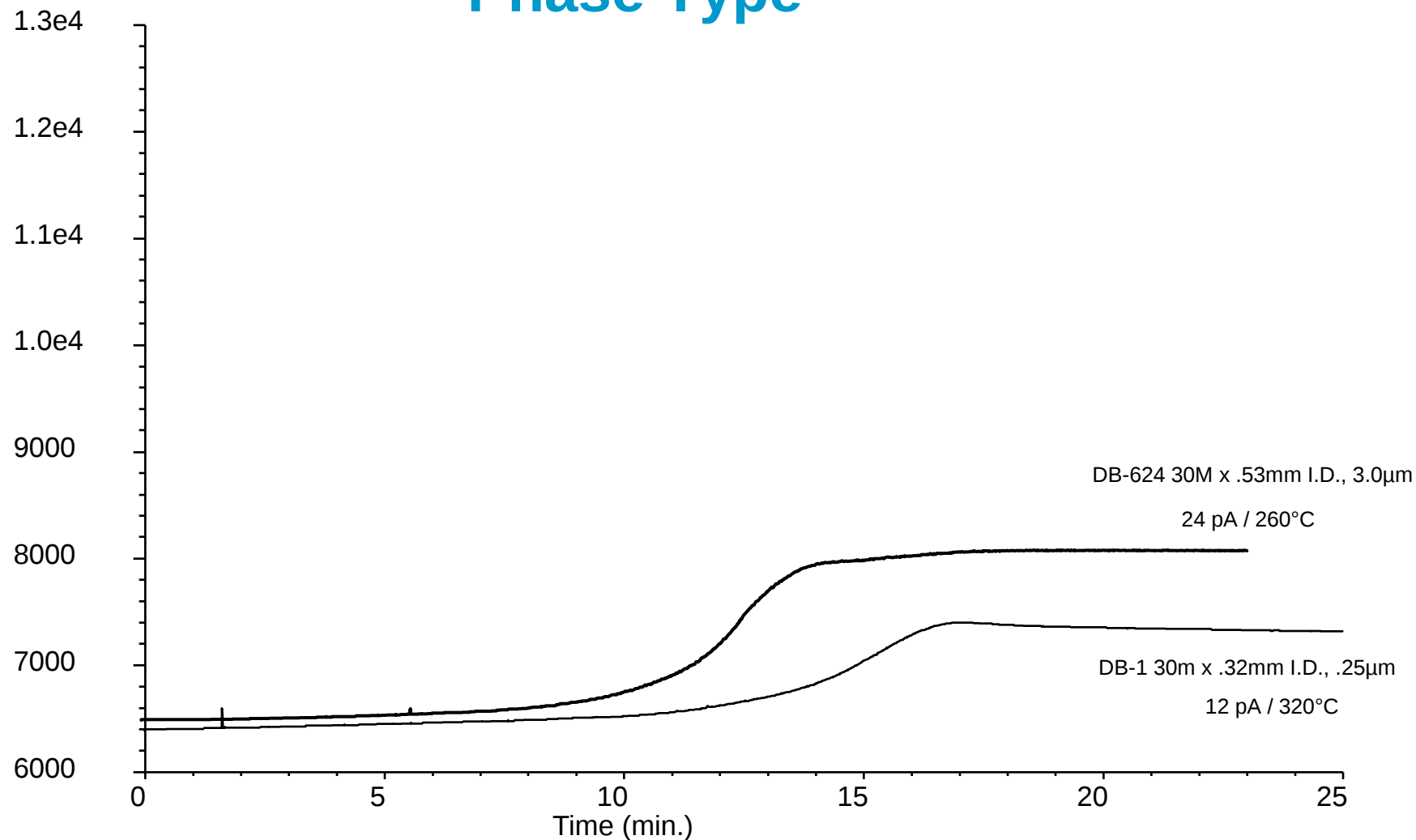
- Bleed if at high temperature
- In detector flame (poor installation)

DETECTOR

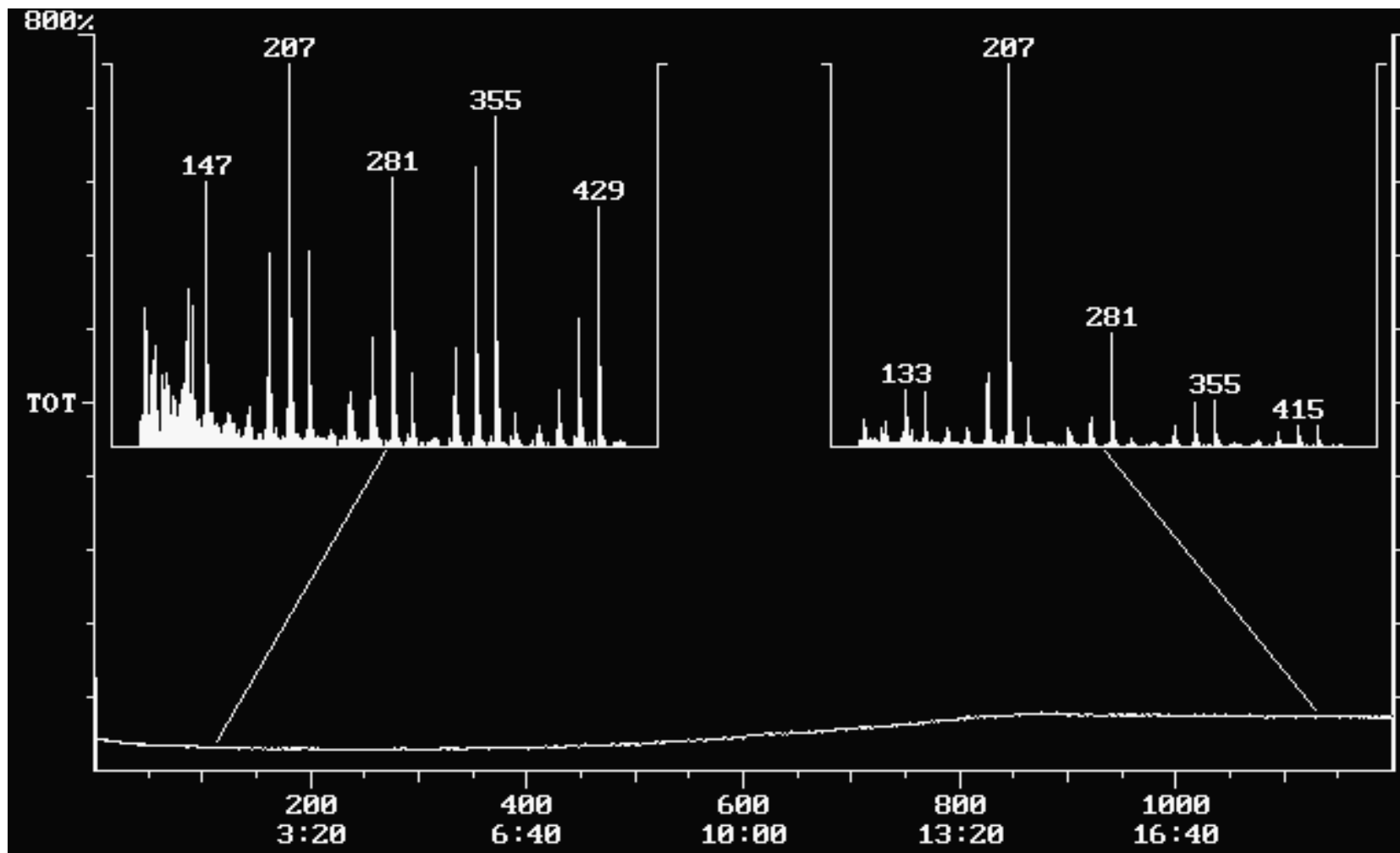
- Air leak - ECD, TCD
- Electronics malfunction



Column Bleed – Affected by Dimensions and Phase Type



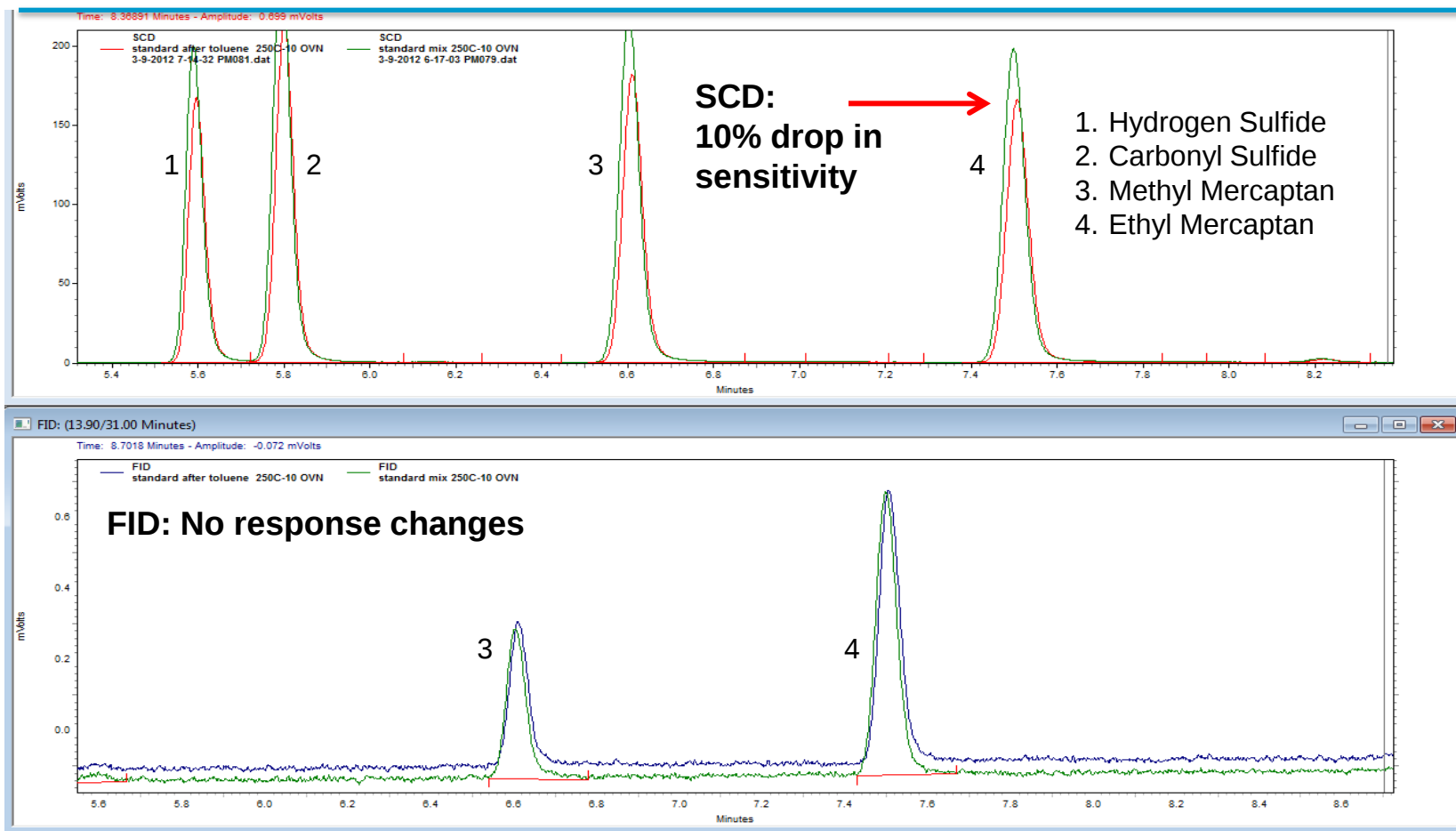
Mass Spectrum of Phenylmethylpolysiloxane Column Bleed (Normal Background)



Mass spectral library search is not always accurate

Traditional PDMS column- Coking (desensitization) of Reactor Tubes

Overlay of before (green) and after 2 x 2 uL neat toluene Injection (red)



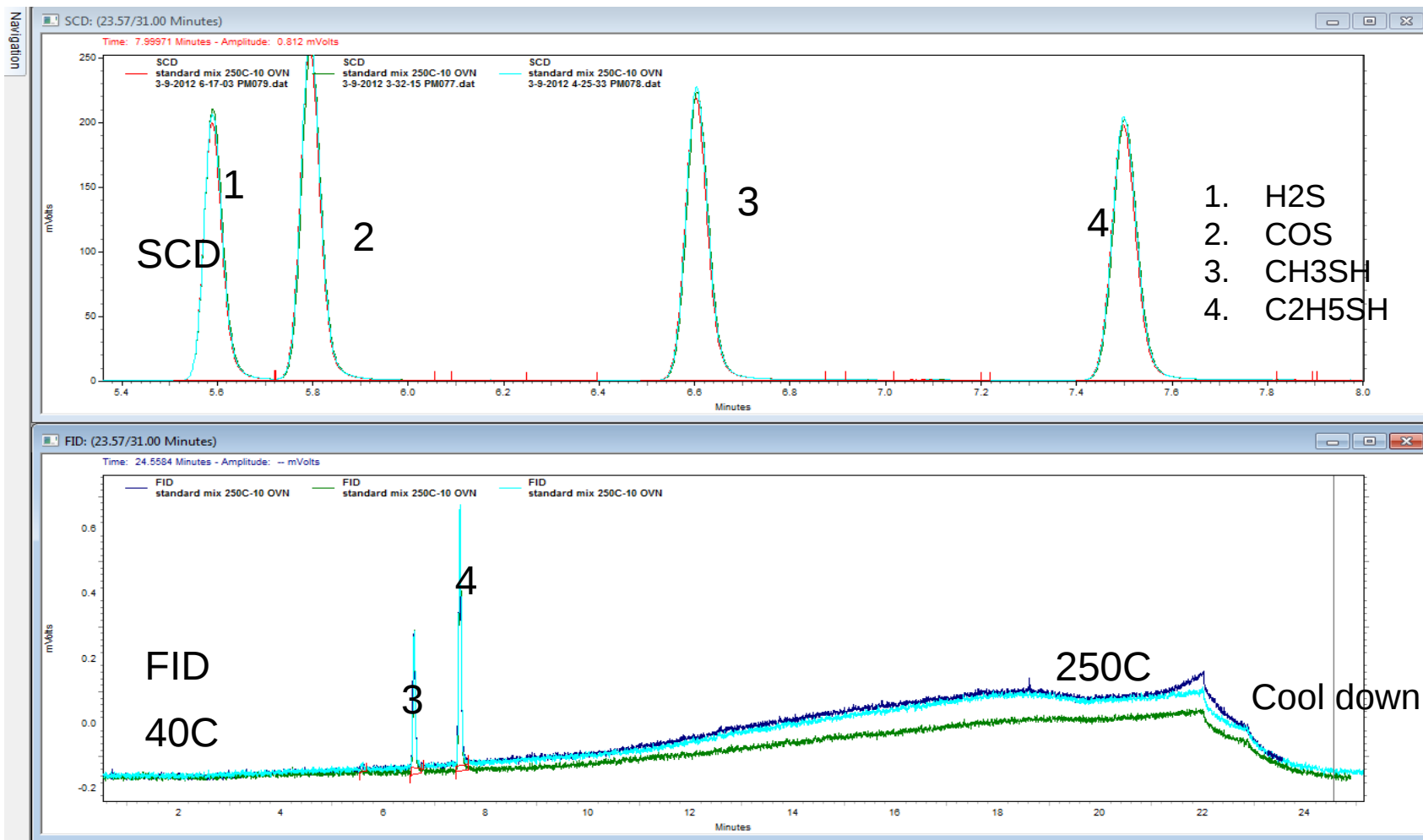
Data courtesy of Jim Luong, Ronda Gras, Myron Hawryluk of Dow Chemical Canada



Agilent Technologies

New DB-Sulfur SCD column

Last Three Runs of the day (n=20, 100 ppmv std)

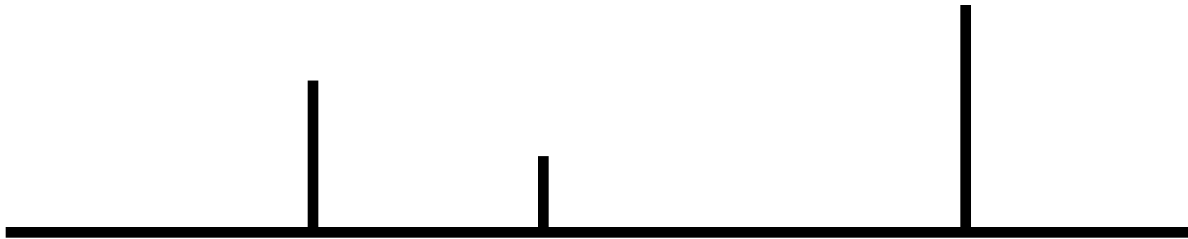


Data courtesy of Jim Luong, Ronda Gras, Myron Hawryluk of Dow Chemical Canada



Agilent Technologies

Spiking Baseline



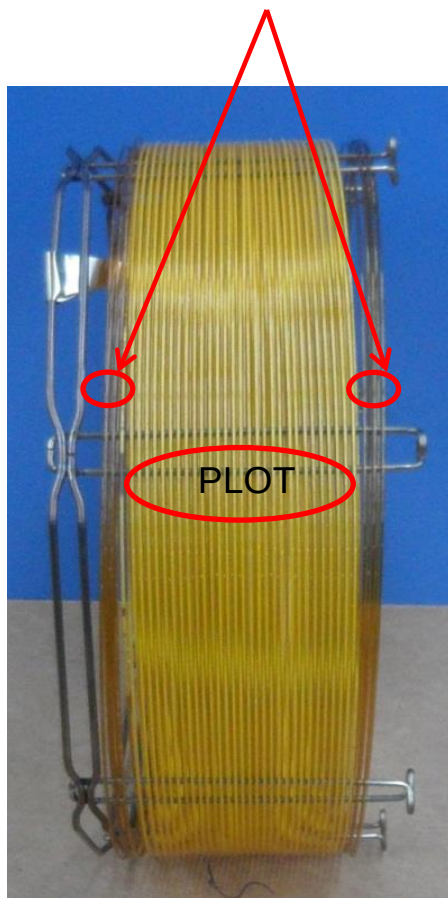
DETECTOR

- Particles entering the detector
- Random: poor connection
- Regular: nearby "cycling" equipment (electronics)



NEW Integrated Particle Trap PLOT Column

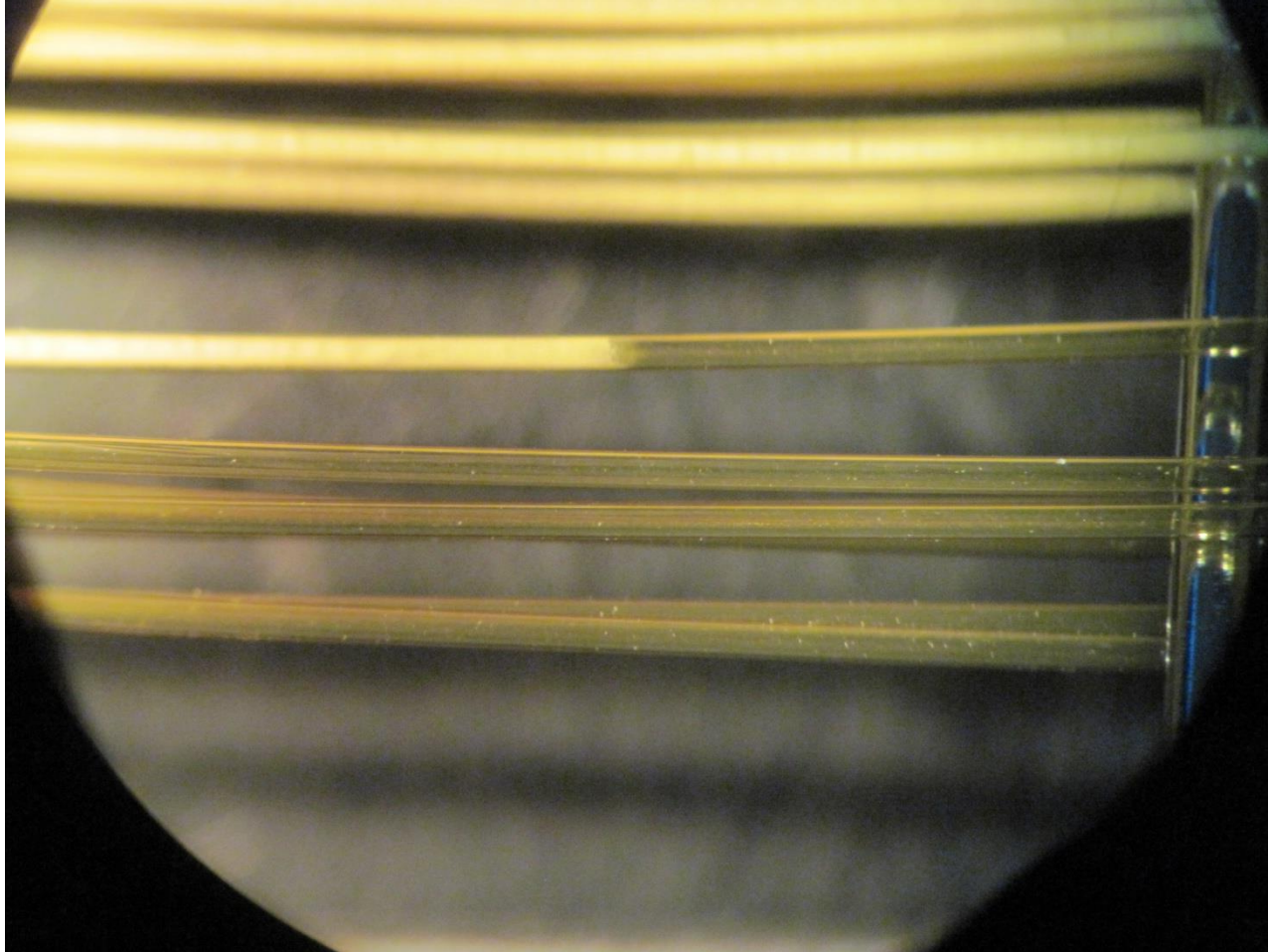
2.5 meter integrated particle traps on both ends



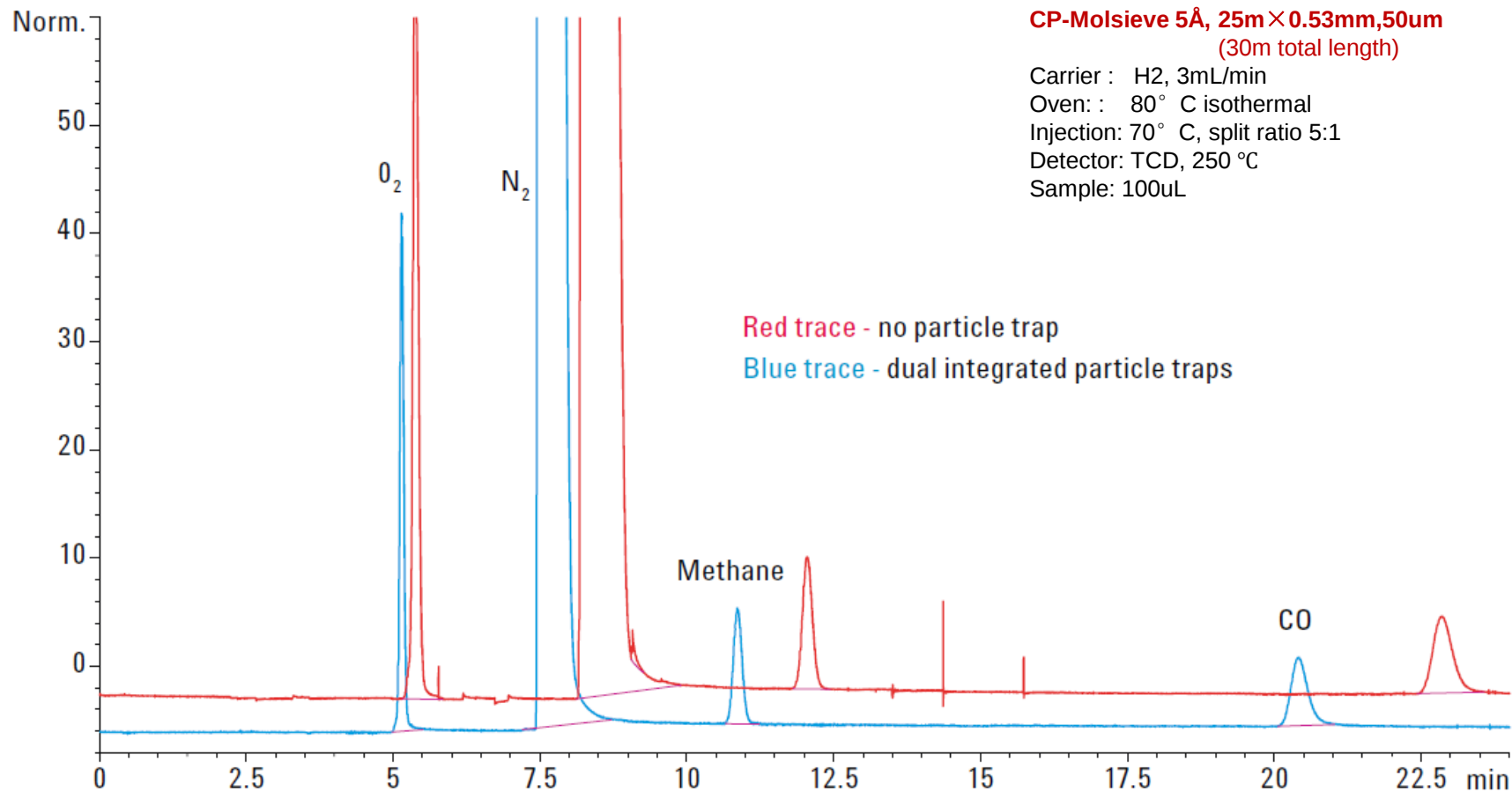
Compatible with GC/MS and valve switching systems,
and Capillary Flow Technology

Very similar selectivity, plates and peak shape to
existing Agilent PLOT columns

The PLOT Column and Integrated Particle Trap



No spikes at Fixed Gases Analysis on CP-Molsieve 5Å PLOT PT column



Quantitation Problems

DETECTOR

- Poor stability (electronics) or Baseline disturbances (contamination)
- Outside detector's linear range or wrong settings

Activity (adsorption) in INJECTOR or COLUMN

INJECTOR

- Technique, settings, conditions
- Syringe worn

OTHER

- Co-elution
- Matrix effects
- Sample evaporation – leaky vials
- Sample decomposition



Agilent 5975C GC/MS

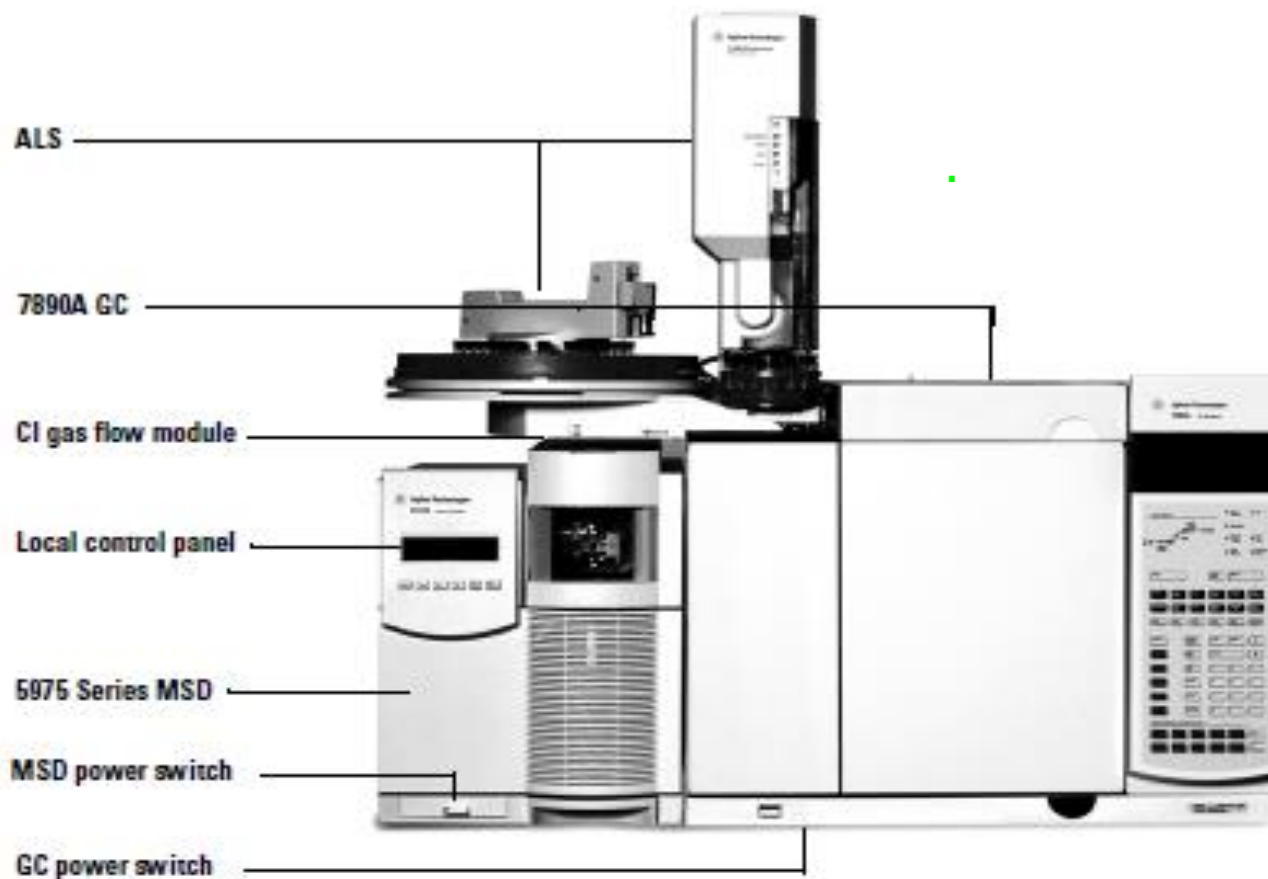
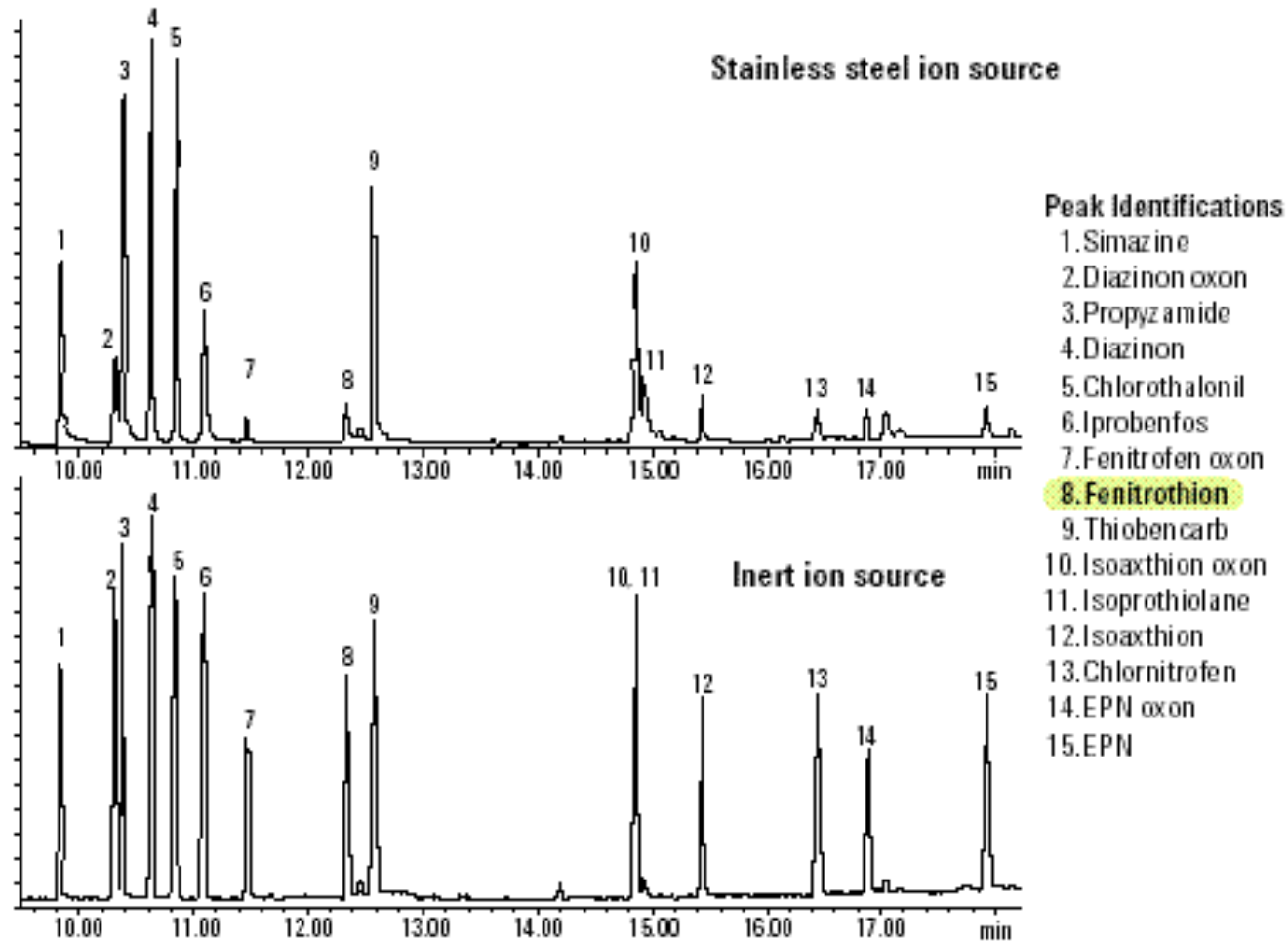


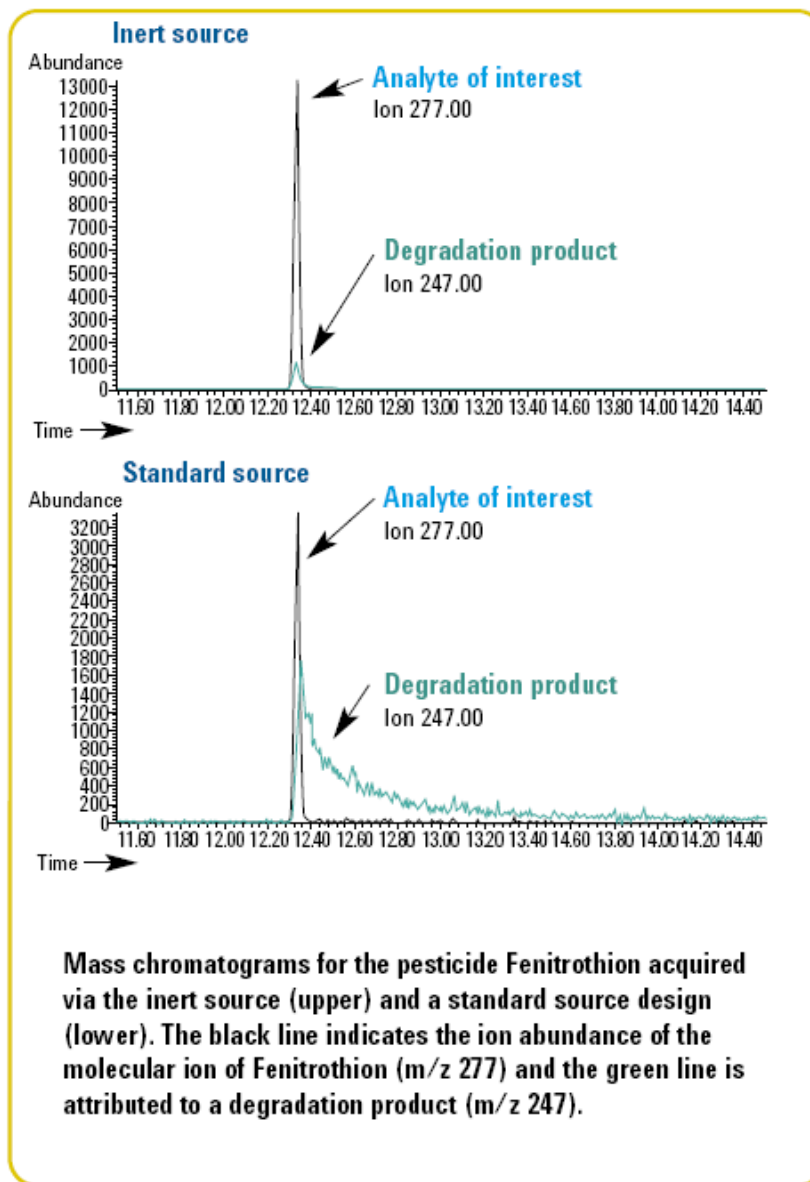
Figure 1 5975 Series GC/MSD system

MSD Source Inertness – an old problem revisited

Inert MSD source increases response of chemically active analytes



Inert MSD Source



Mass Spectral Problems

High background

- Excessive pressure in the analyzer chamber
- Air leak
- Contamination

Mass Spectral Problems

High abundances at m/z 18,28,32 and 44

- System recently vented (pumping down)
- Masses at 14,16 and 19 are indicative of a massive leak



Common Contaminant Ions

Ions (m/z)	Compound	Possible source
18, 28, 32, 44 or 14, 16	H ₂ O, N ₂ , O ₂ , CO ₂ or N, O	Residual air and water, air leaks, outgassing from Vespel ferrules
31, 51, 69, 100, 119, 131, 169, 181, 214, 219, 264, 376, 414, 426, 464, 502, 576, 614	PFTBA and related ions	PFTBA (tuning compound)
31	Methanol	Cleaning solvent
43, 58	Acetone	Cleaning solvent
78	Benzene	Cleaning solvent
91, 92	Toluene or xylene	Cleaning solvent
105, 106	Xylene	Cleaning solvent
151, 153	Trichloroethane	Cleaning solvent
69	Foreline pump oil or PFTBA	Foreline pump oil vapor or calibration valve leak
73, 147, 207, 221, 281, 295, 355, 429	Dimethylpolysiloxane	Septum bleed or methyl silicone column bleed
77, 94, 115, 141, 168, 170, 262, 354, 446	Diffusion pump fluid and related ions	Diffusion pump fluid
149	Plasticizer (phthalates)	Vacuum seals (O-rings) damaged by high temperatures, vinyl gloves
Peaks spaced 14 m/z apart	Hydrocarbons	Fingerprints, foreline pump oil



But at the End of the Day...It's in the Inlet!

Do Inlet Maintenance

- Trim the Column
- Replace the Liner
- Replace the Gold Seal
- Replace the Syringe
- Replace the Split Vent Trap, Split Vent Line
- Rinse the Inlet



Common Care and Maintenance Scheme for GC Columns

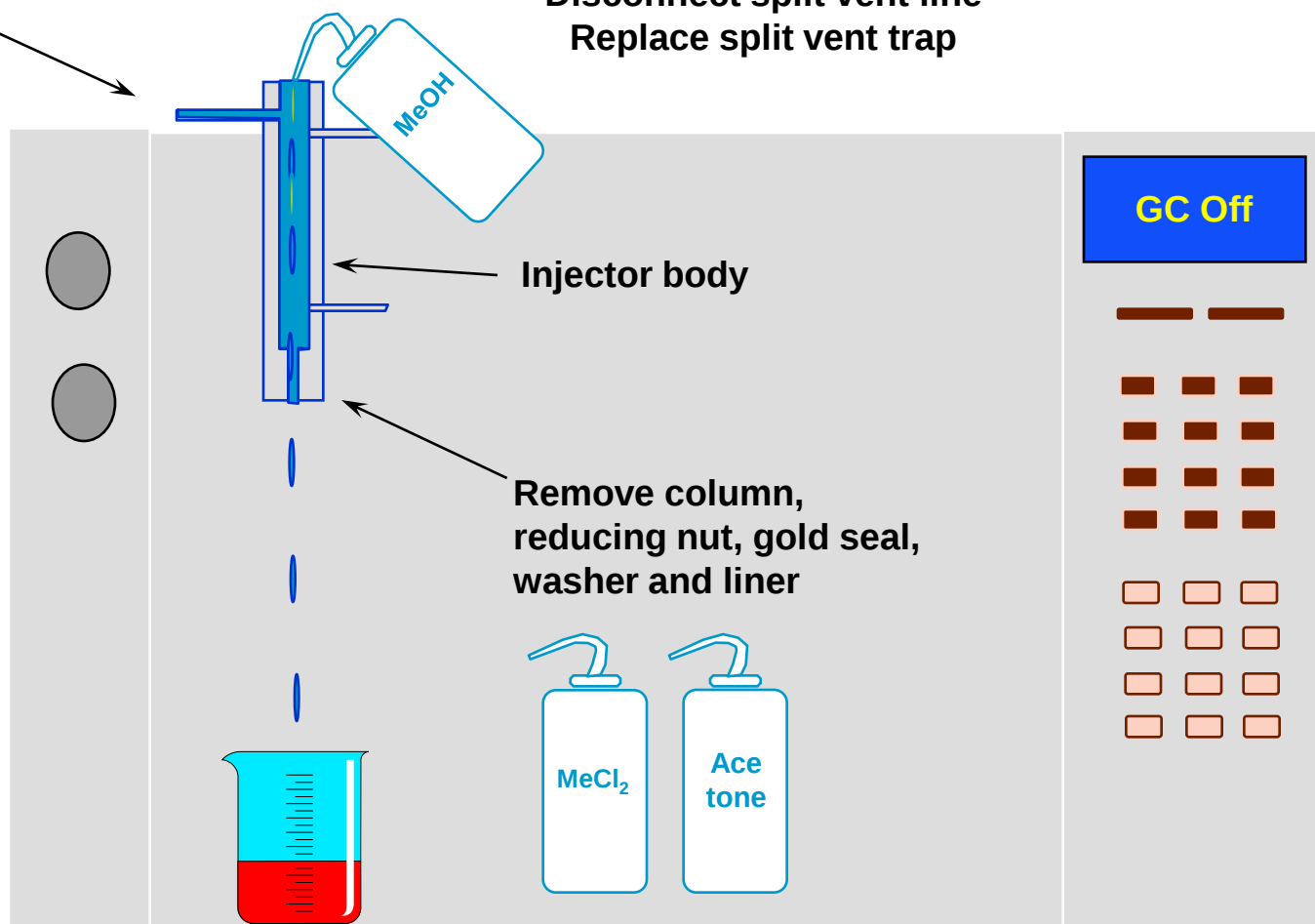
1. Bake out the column for no more than 2 hours.
2. Cut off 6"-1ft of the inlet end of the column.
3. Cut off more column. (repeat as necessary)



Cleaning the Split/Splitless Injector

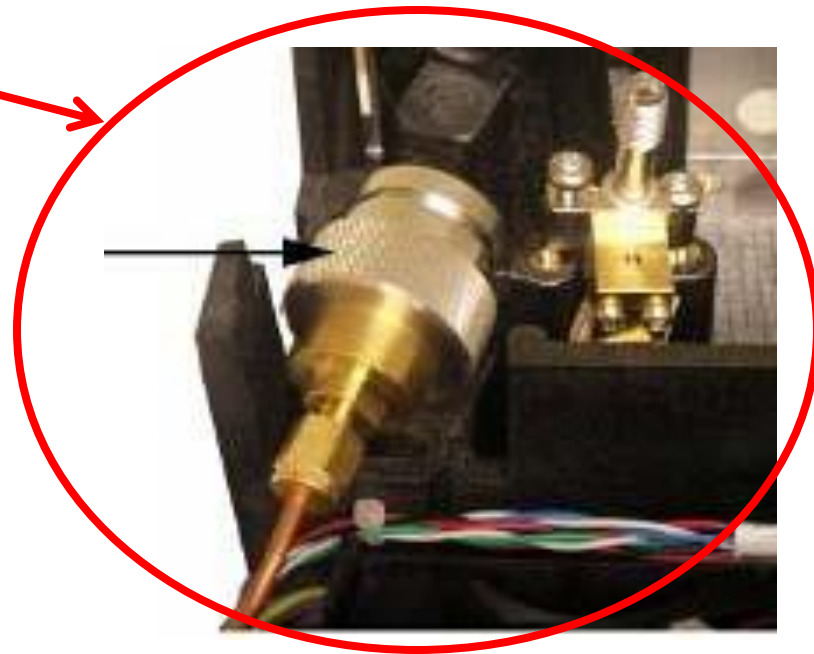
Carrier gas flow off

Disconnect split vent line
Replace split vent trap



Finding the Split Vent Trap

Remove cover at Split Vent

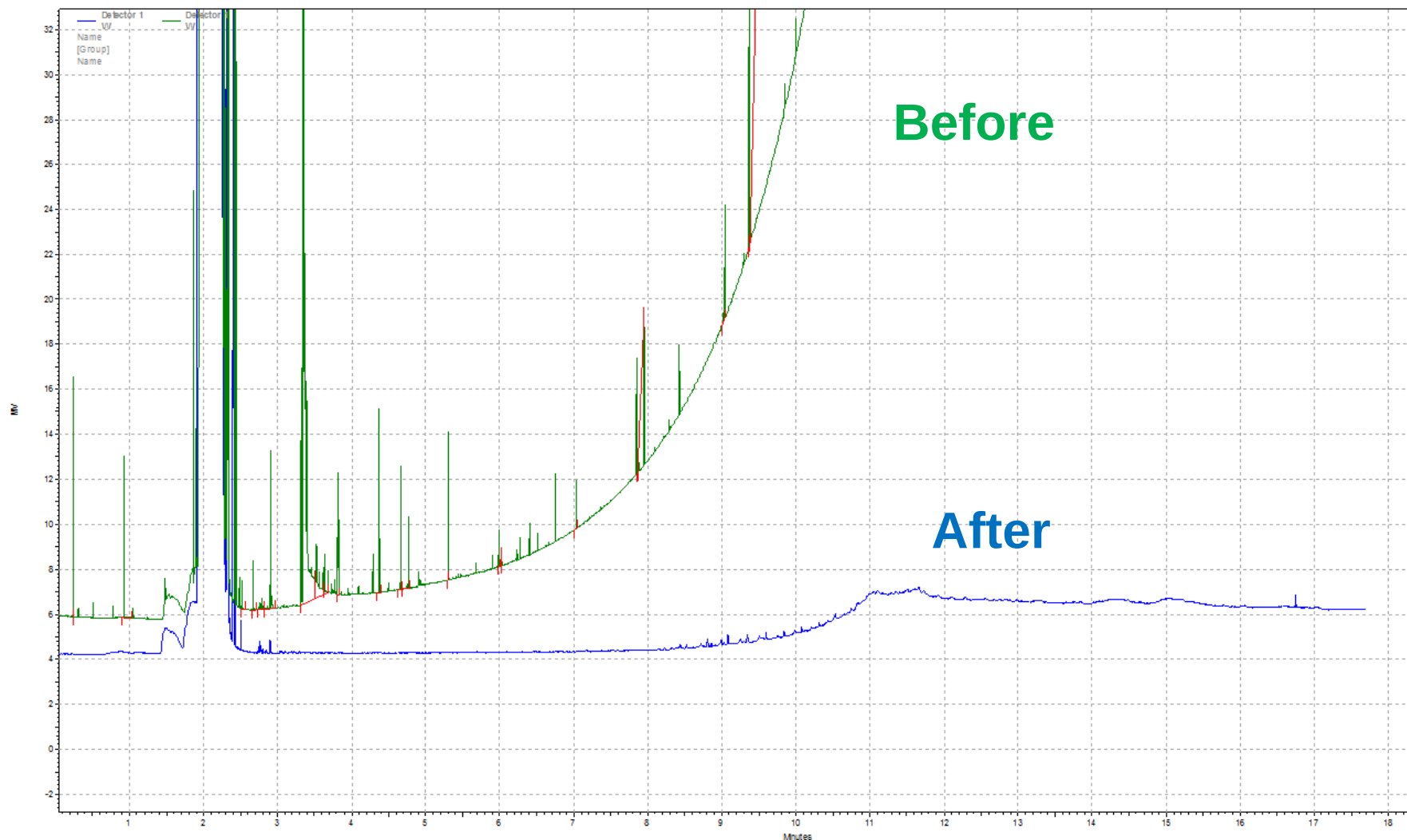


Replacing the Split Vent Trap

Finger Tight Knurled Nut



Split Vent Trap Changed (Column Bleed?!?)

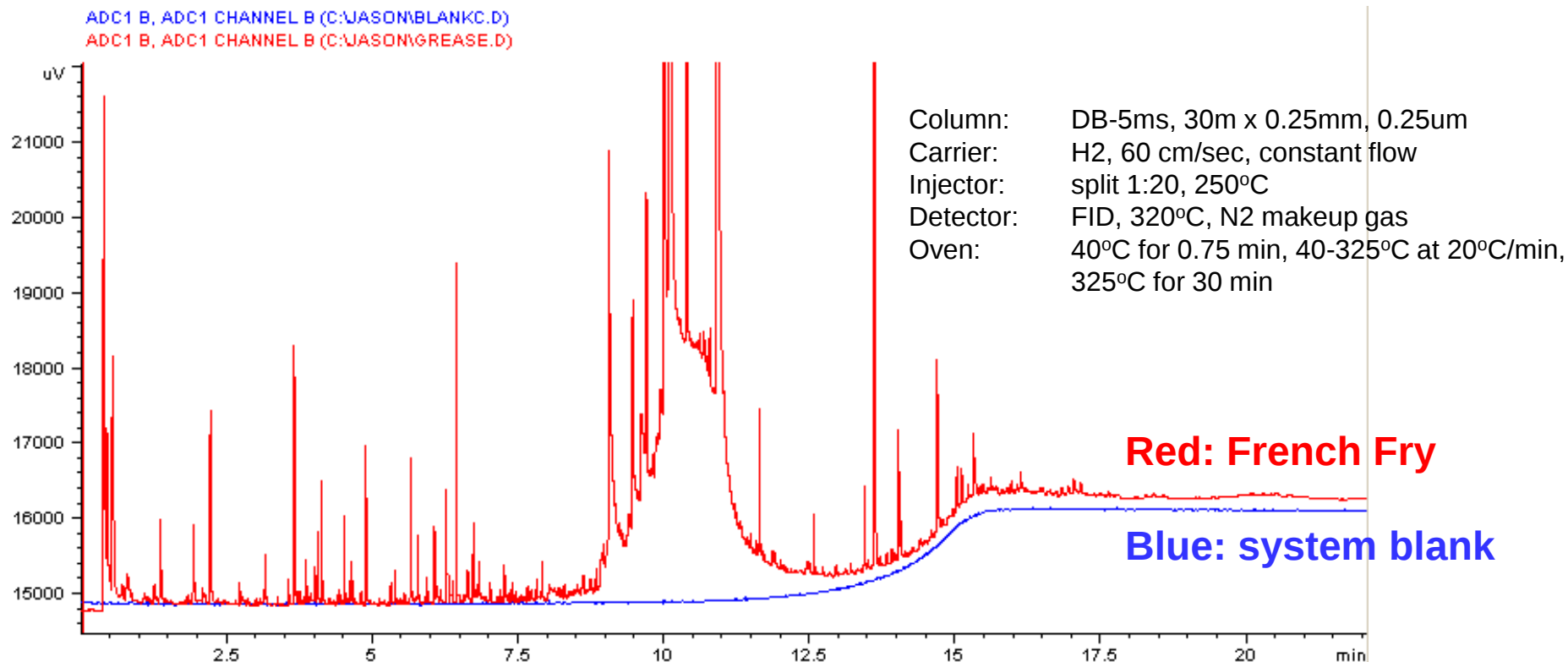


Be Careful When Doing Maintenance...

You may be the CONTAMINATOR!



Contamination of system by residue on fingers during column installation

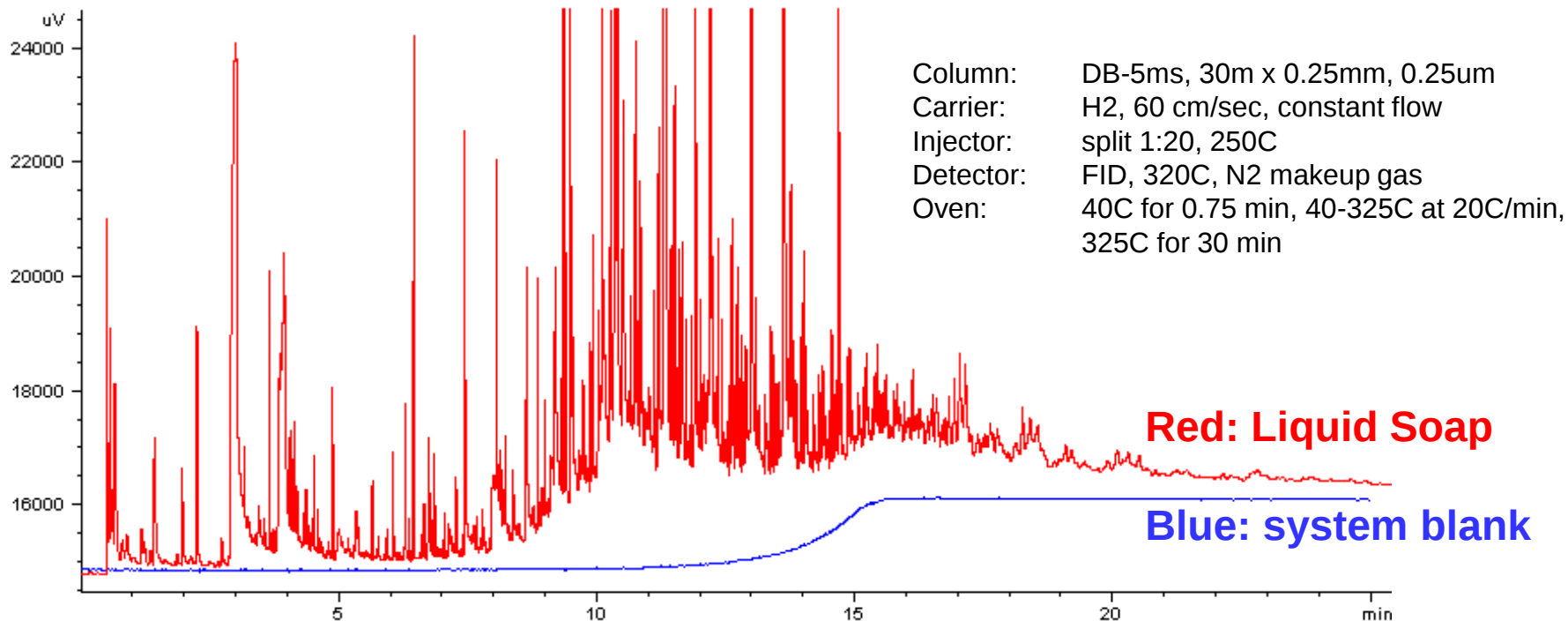


Procedure:

- (1) Held French fry for 5 seconds.
- (2) Fingertip was wiped with paper towel to remove as much of the offending material as possible.
- (3) Lightly touched the part of the column sticking up above the ferrule.
- (4) Installed column into injector.
- (5) Set oven temperature to 40°C.
- (6) Started oven temperature program as soon as oven reached 40°C.



Contamination from Liquid Soap

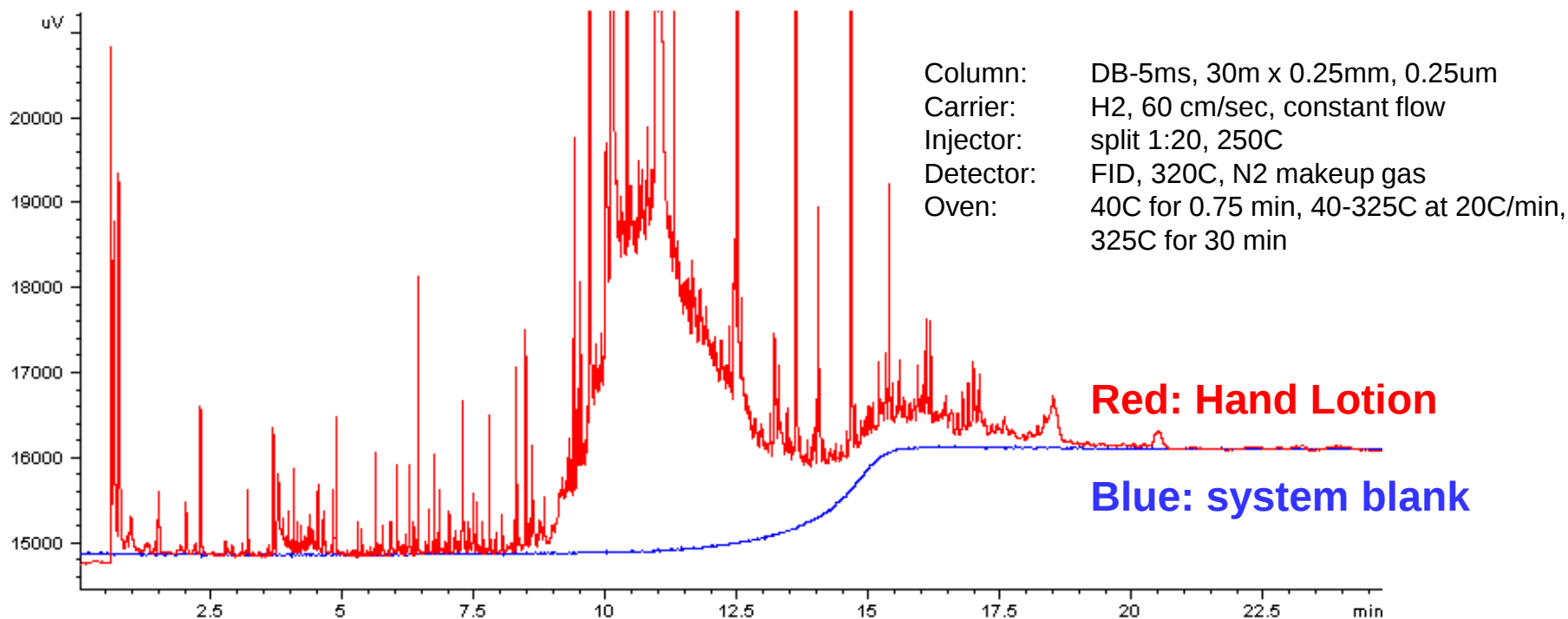


Procedure:

- (1) One very small drop of liquid soap placed on one fingertip.
- (2) Fingertip was wiped with paper towel to remove as much of the offending material as possible.
- (3) Lightly touched the part of the column sticking up above the ferrule.
- (4) Installed column into injector.
- (5) Set oven temperature to 40C.
- (6) Started oven temperature program as soon as oven reached 40C.



Contamination from Hand Lotion

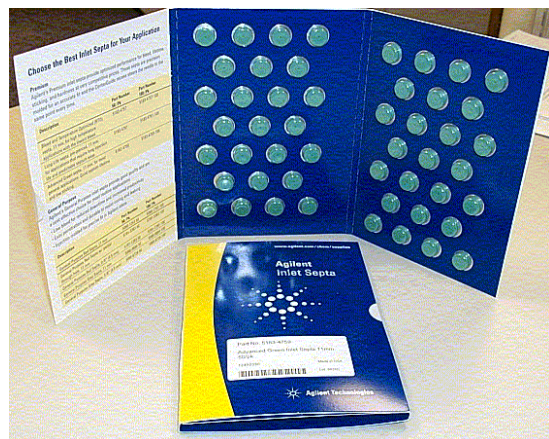


Procedure:

- (1) One very small drop of hand lotion placed on one fingertip.
- (2) Fingertip was wiped with paper towel to remove as much of the offending material as possible.
- (3) Lightly touched the part of the column sticking up above the ferrule.
- (4) Installed column into injector.
- (5) Set oven temperature to 40C.
- (6) Started oven temperature program as soon as oven reached 40C.



More and More “Touchless” Packaging



Troubleshooting “Tools”

Bleed Profile: *baseline problems*

Inject a non-retained peak: *peak shape problems*

Leak Test: *all problems*

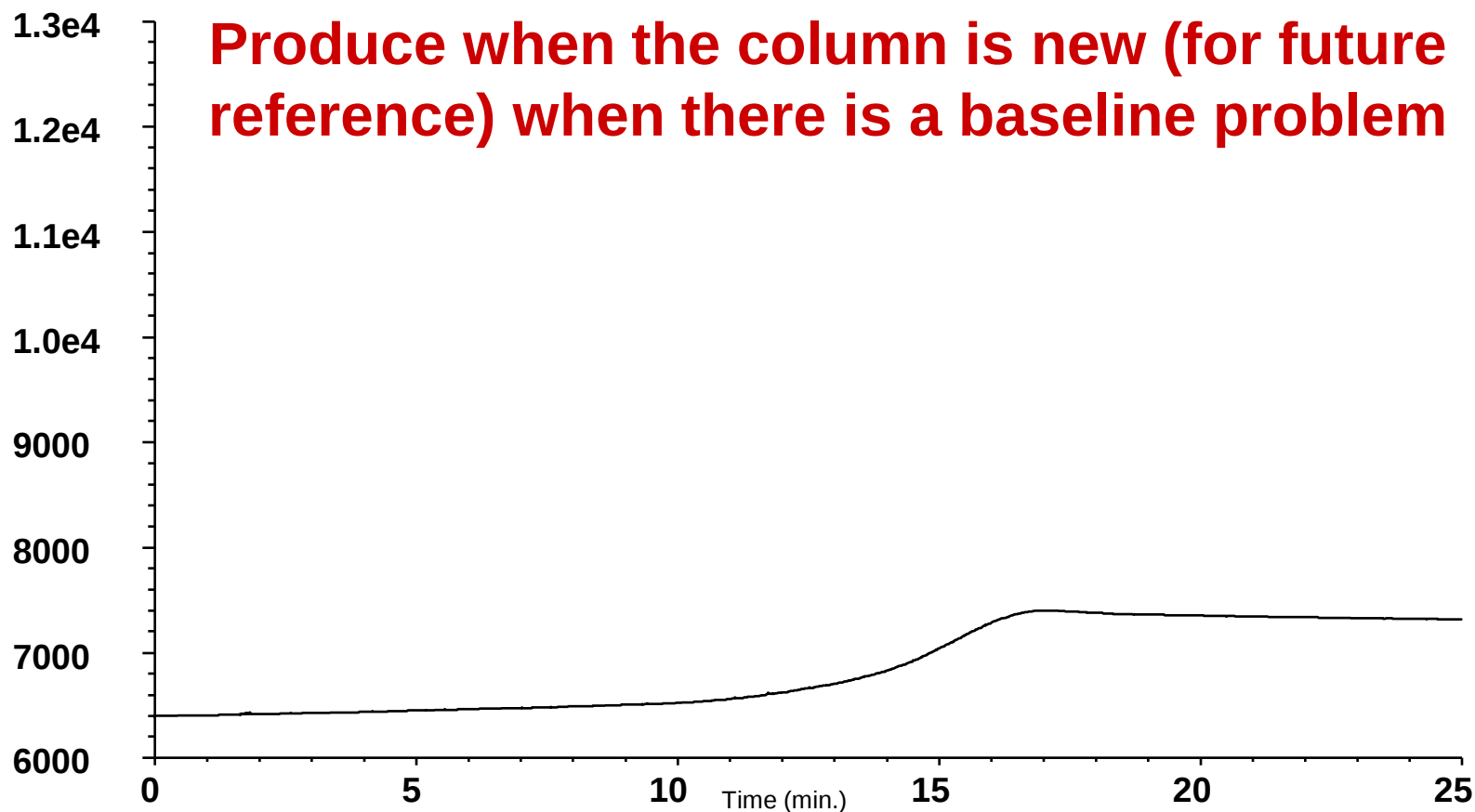
Test mix: *all problems*

Isolate the components: *all problems*

Jumper Tube Test: *baseline problems*



Generating a Bleed Profile

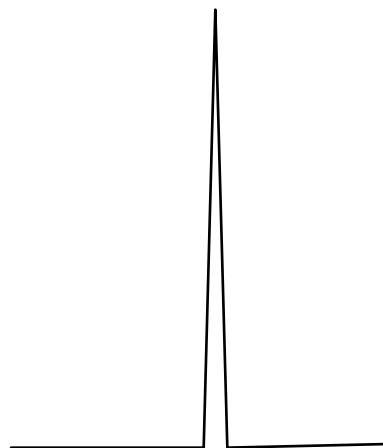


***DB-1 30m x .32mm I.D., .25 μ m**

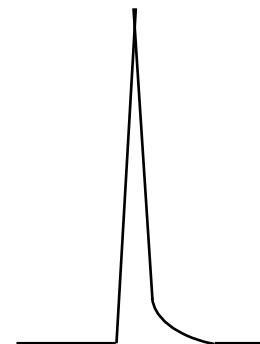
Temperature program // 40°C, hold 1 min // 20°/min to 320°C, hold 10 min.

Non-Retained Peak Shapes

**Used to Check
Flowpath**



Good Installation



Improper Installation or
Injector Leak

Potential problems:

- Injector or septum leak

- Too low of a split ratio

- Liner problem

 - (broken, leaking, misplaced)

- Column position in injector and detector



Leak Test

- MS – Air/Water
- Leak Detector (G3188B)
 - Portable, handheld unit – only 105g
 - USB cable for recharging
 - Audible and visual alerts for 12 gases
 - Minimum detection limit of 0.0005 mL/minute for hydrogen and helium
 - Rechargeable lithium ion battery with over 5 hours of life



Ferrule Pre-swaging & MS Interface Installation Tools

- Ensures proper length of column into the fittings, every time
- For graphite or metal ferrules

Metal ferrule
G3440-80218

Graphite ferrule
G3440-80217



Self-Tightening Column Nuts

Innovative spring-driven piston continuously presses against ferrule

- **Less wasted time:** No retightening needed after repeated thermal cycles
- **Ease of use:** Finger-tight, consistent connections *without tools*
- **Leak Free = Lower column bleed:** Longer column life

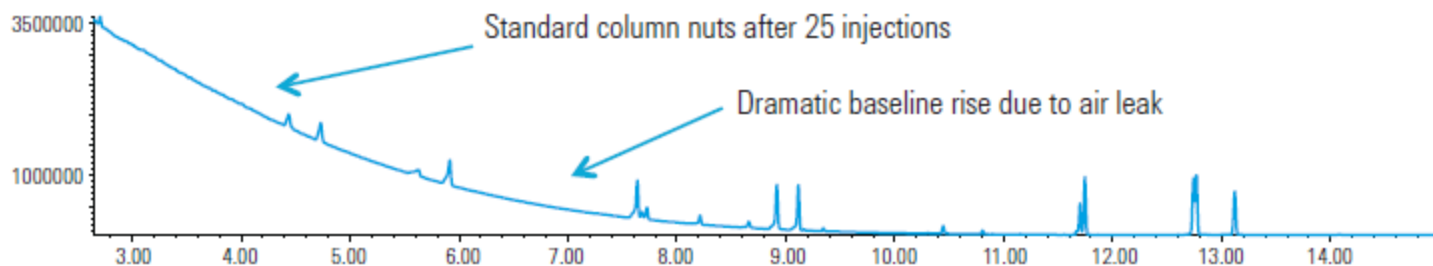
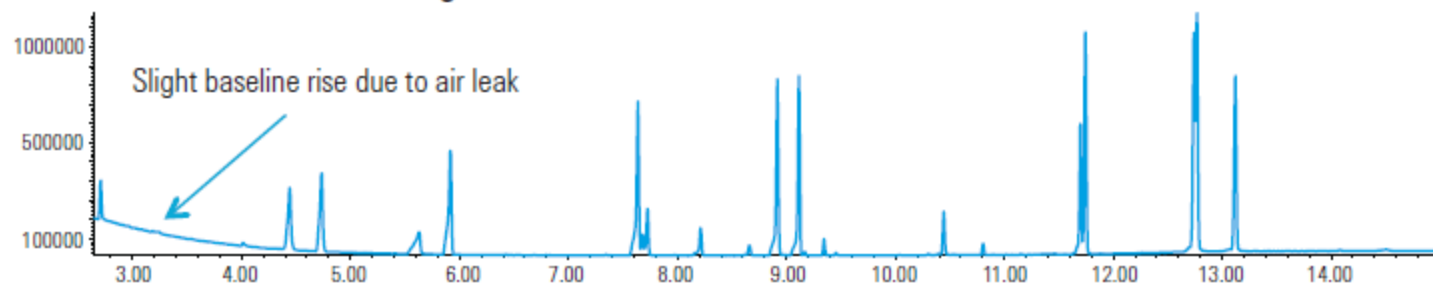


Video at [agilent.com/chem/STnutvideo](https://www.agilent.com/chem/STnutvideo)

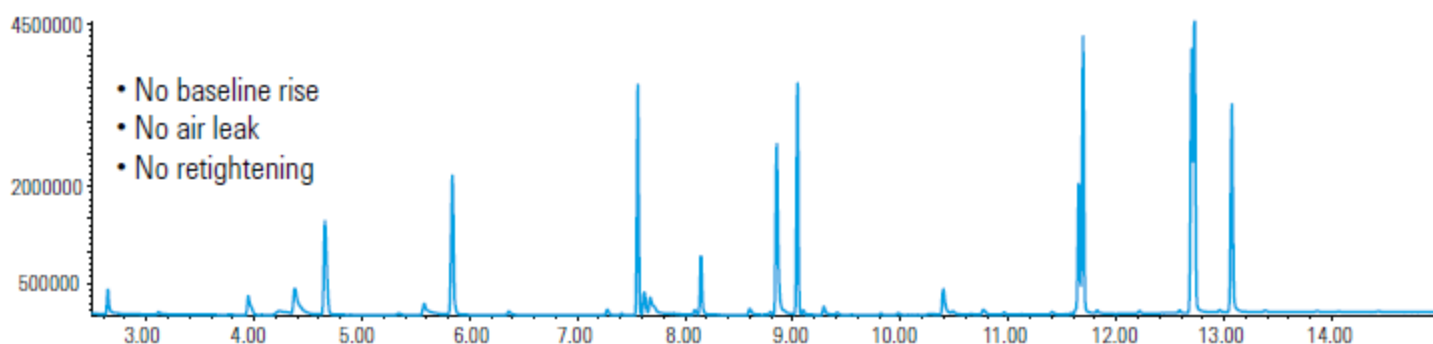


Self-Tightening Column Nuts

Standard column nuts new fitting

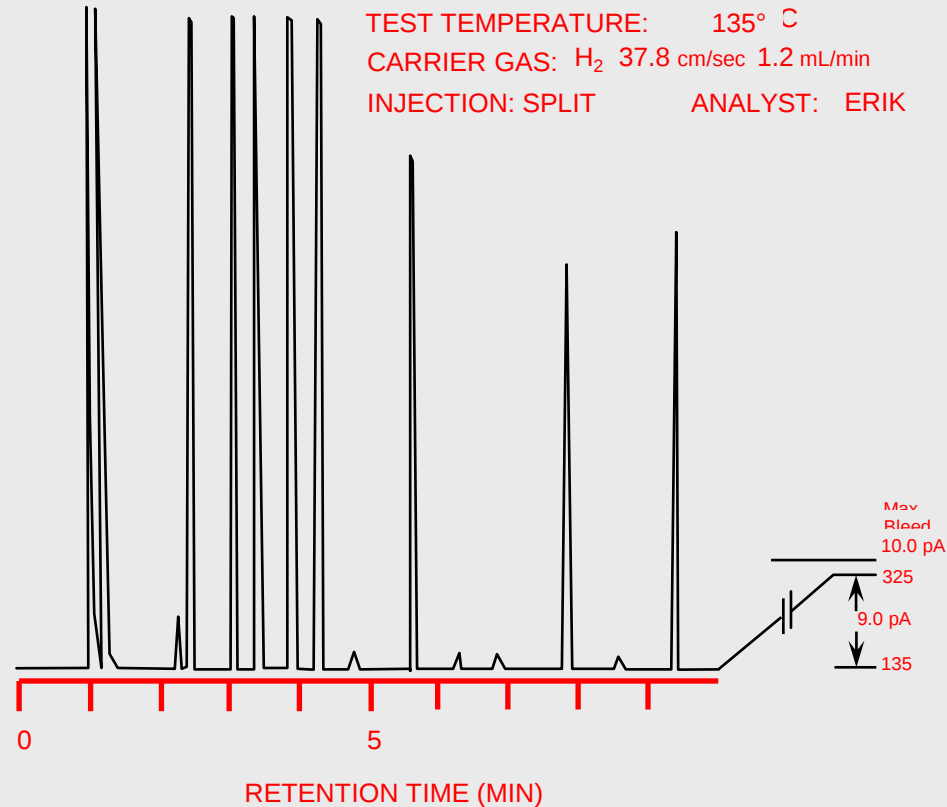


Agilent Self Tightening Column Nuts after 400 injections



Test Mix

Used to determine how “good” the column is or if the problem is related to the chemical properties of the analytes.



Test Mixture Components

Compounds

Hydrocarbons

Alcohols

FAME's, PAH's

Acids

Bases

Purpose

Efficiency

Retention

Activity

Retention

Acidic Character

Basic Character



Own Test Mixture

More specific

Selective detectors

Actual concentrations

No conditions or instrument changes



Jumper Tube Test

Purpose

Helps to locate the source of contamination or noise

Isolates GC components



Jumper Tube Test

Isolate the Detector

Remove column from the detector

Cap detector and turn on

Blank run



Jumper Tube Test

Isolation of Detector - Results



Detector OK



Detector is the problem



Jumper Tube Test

Isolate the Injector

Connect the injector and detector

- 1-2 meters deactivated fused silica tubing

Turn on carrier gas

Blank run



Jumper Tube Test

Isolate the Injector - Results



Injector OK



**Injector, lines or carrier
gas contaminated**

Jumper Tube Test

Isolate the Column

Reinstall the column

Setup as before

Blank Run



Jumper Tube Test

Isolate the Column - Results

Problem returns: It's the column

**Problem gone: Previous leak, solid debris,
or installation problem**



Jumper Tube Test

Used to Isolate Source of Contamination

- . Cap off the detector and establish normal gas flows and temperature.
- . Plot the baseline using a temperature program. If flat.....
- . Connect 1 meter of deactivated tubing between the injector and detector
- . Plot the baseline using a temperature program. If flat.....
- . Install the column.
- . Plot the baseline using a temperature program.



Remember

**Complete system = Carrier Gas + Injector +
Column + Detector + Data System**

Multiple cause and effect

**Do not change too many variables at
once**



Symptoms and Reasons (Cheat Sheet)

**Solution = Get Rid of the Reason*

- Peak Tailing – Flow Path or Activity
- Bonus Peaks – In Sample or Back Flash (Carry Over)
- Split Peaks – Injector Problems, Mixed Solvent, Cold Spots
- Broad Peaks – Injection Problems, Temperature/Flow, Solubility Problems
- No Peaks – Wasn't Introduced, Wasn't Detected
- Response Changes – Activity, Injector Discrimination, Detector Problem
- Peak Fronting – Overload or Solubility Mismatch, Injector Problems
- Shifting Retention – Leaks, Column Aging, Contamination or Damage
- Loss of Resolution – Separation Decreasing, Peak Broadening
- Baseline Disturbances – Column Bleed, Contamination, Electronics
- Noisy or Spiking Baseline – Electronics or Contaminated Detector
- Quantitation Problems – Activity, Injector or Detector Problems



Troubleshooting Resources

Online Troubleshooting and Maintenance Videos

<http://www.chem.agilent.com/en-US/Technical-Support/Instruments-Systems/Gas-Chromatography/Pages/troubleshootingvideos.aspx>

GC Troubleshooting Guide

<http://www.chem.agilent.com/en-US/Products-Services/Instruments-Systems/Gas-Chromatography/pages/gp6770.aspx>

Method Translation Software

<http://www.chem.agilent.com/en-US/Technical-Support/Instruments-Systems/Gas-Chromatography/utilities/Pages/gcmethodtranslation.aspx>



Agilent Better GC Connections

www.agilent.com/chem/betterGCconnections

Order the poster...

SIX TIPS FOR TIGHTER GC CONNECTIONS AND BETTER RESULTS

Inspecting your GC column connections is a key part of good preventive maintenance – and a laboratory practice that you simply cannot afford to overlook. That's because poor, leaky connections can cause:

- Empty baselines
- Loss of expensive, high-quality gas
- Shorter column and detector life
- Decreased system sensitivity
- Reduced system productivity

This poster highlights the critical GC connection 'hotspots' to help you fix problems before they compromise your results.

The Mechanics of Confidence

There is a lot that often goes wrong when the quality and consistency of your data.

To learn more about creating and maintaining leak-free GC connections, go to www.agilent.com/chem/betterGCconnections

Agilent Technologies

The poster features a diagram of a GC system with numbered hotspots (1-6) and six numbered tips on the right side. The tips are: 1. Inspect the ferrule and nut, and use the appropriate tool for your application. 2. Inspect the ferrule and nut, and use the appropriate tool for your application. 3. Inspect the ferrule and nut, and use the appropriate tool for your application. 4. Inspect the ferrule and nut, and use the appropriate tool for your application. 5. Inspect the ferrule and nut, and use the appropriate tool for your application. 6. Inspect the ferrule and nut, and use the appropriate tool for your application.

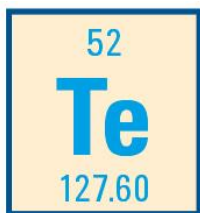
View the video...

Better_GC_Connections_08c-FINAL_3280x720

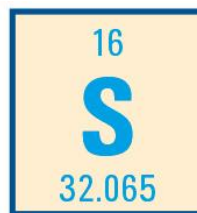
Agilent Self Tightening Column Nut with short graphite polyimide blend ferrule

00:52

The video shows a close-up of a GC column connection. A hand is shown using a tool to tighten a column nut. The video is titled 'Better GC Connections' and shows a close-up of a GC column connection. A hand is shown using a tool to tighten a column nut. The video is titled 'Better GC Connections' and shows a close-up of a GC column connection. A hand is shown using a tool to tighten a column nut.



chnical



upport

1-800-227-9770, #3

1 (972) 699-6423 (Daron)

1 (214) 883-2260 (Eric)



E-mail:

gc-column-support@agilent.com

Eric.pavlich@agilent.com

Daron_Decker@agilent.com



Agilent Technologies