

Practical Steps in GC Troubleshooting

Techniques, Tips, and Tricks

Simon Jones
Application Engineer
GC Columns & Supplies

What went wrong and how to fix it...

Common problems

Troubleshooting tools

Troubleshooting examples

**“Everything was just fine and then
this happened!”**

How do I go about
TROUBLESHOOTING?”

“Everything was just fine and then this happened!”

Logic = Something changed (slowly or sudden) =
Something is different

Track Events – log book

- Changed column, liner, septum, syringe, etc.
- Injected samples, other method, etc.
- Did maintenance, cut column, inlet flush, etc.

Logical Troubleshooting

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

Typical Problems of Optimized Methods becoming Unoptimized...and the Reason Why.

- Peak Tailing – Flow Path or Activity
- Bonus Peaks – In Sample or Back Flash (Carry Over)
- Split Peaks – Injector Problems, Mixed Solvent
- 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

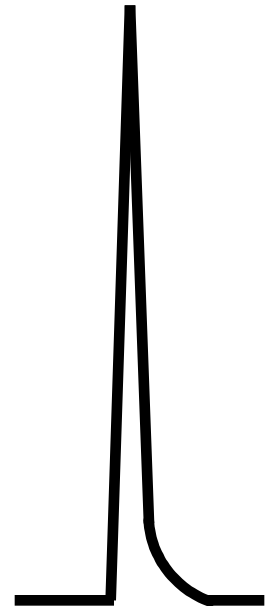
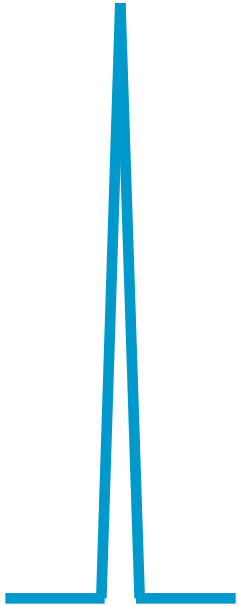
Peak Tailing

INJECTOR or COLUMN is Active

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

FLOW problem

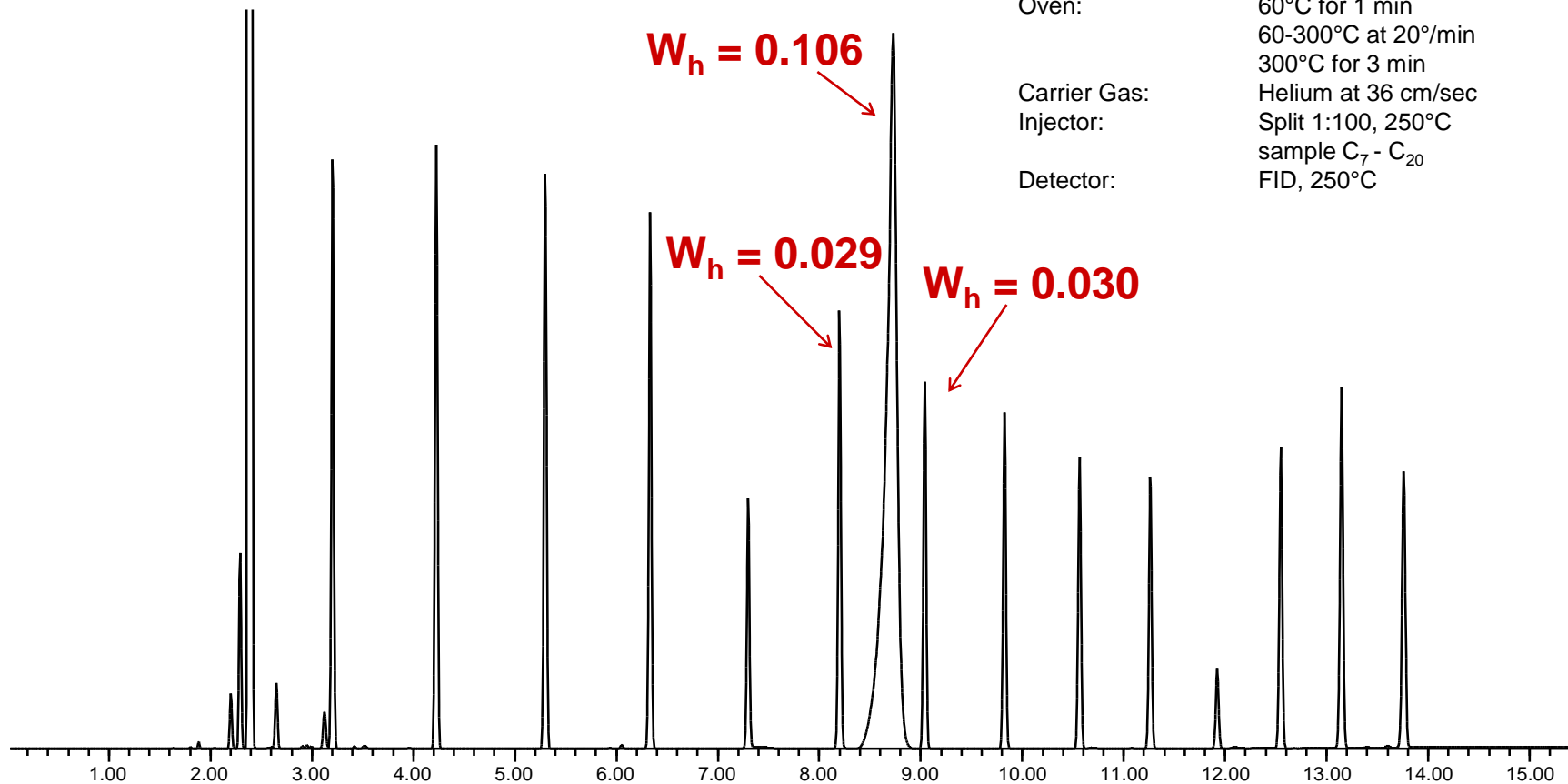
- dead volume, obstruction, poor installation, or
severe column contamination



Miscellaneous - 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.

Bonus Peaks



Column:

DB-5

J&W P/N:

**30 m x 0.53 mm I.D., 1.5 μ m
125-5032**

Oven:

60°C for 1 min
60-300°C at 20°/min
300°C for 3 min

Carrier Gas:

Helium at 36 cm/sec

Injector:

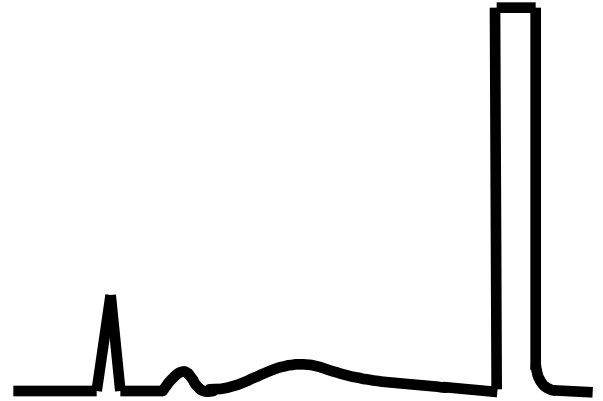
Split 1:100, 250°C

sample C₇ - C₂₀

Detector:

FID, 250°C

Bonus Peaks or Ghost Peaks

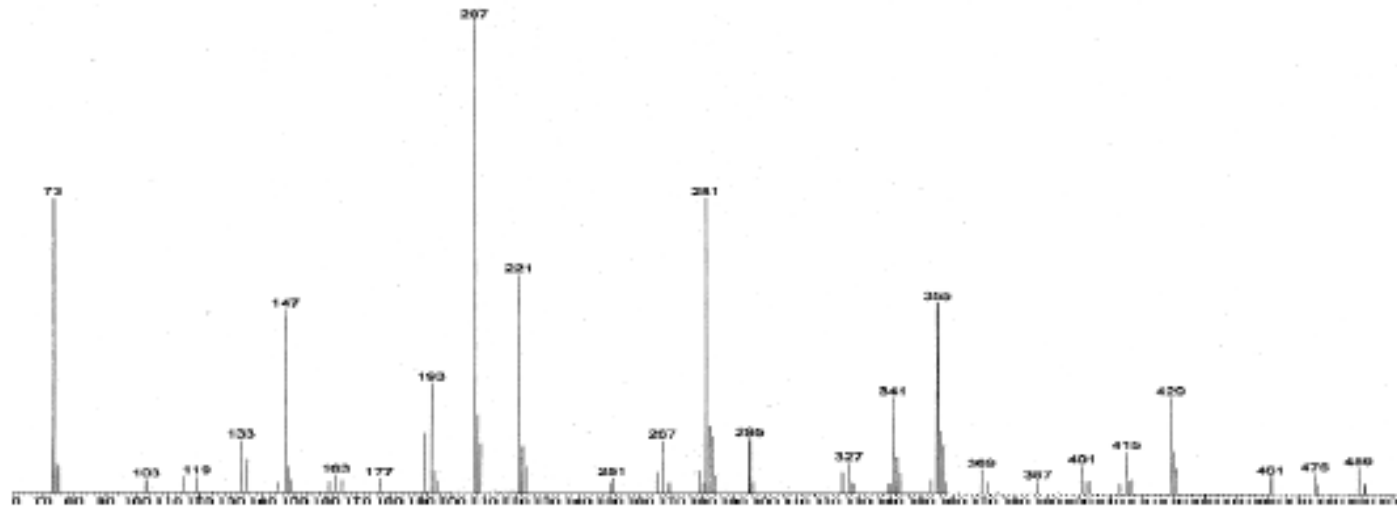
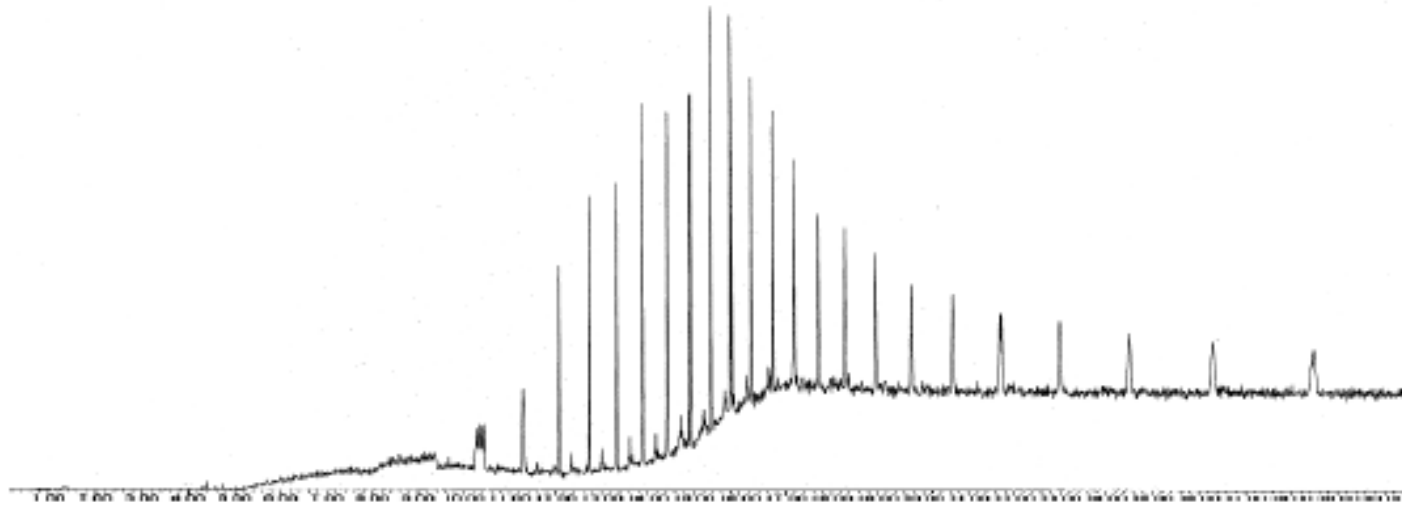


Contamination in INJECTOR, COLUMN or FLOW (carrier gas)

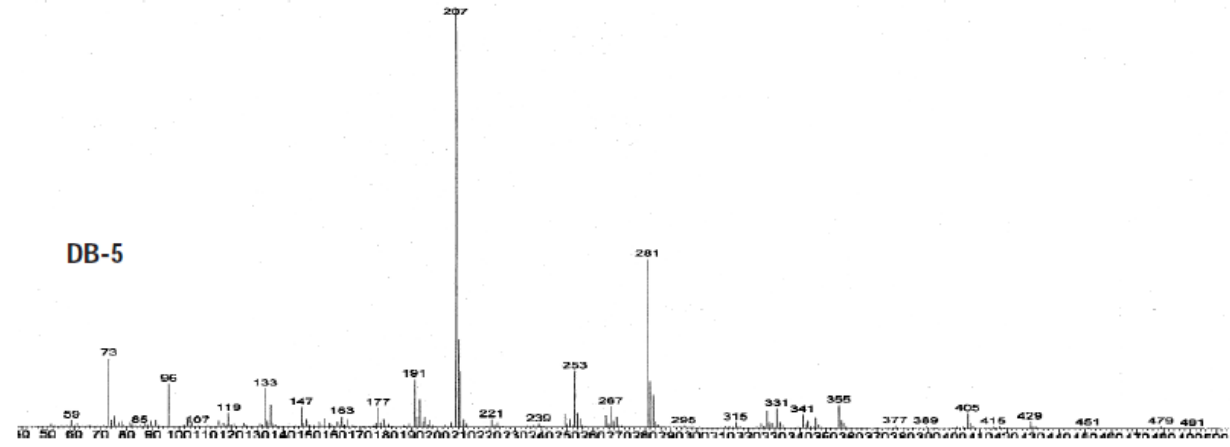
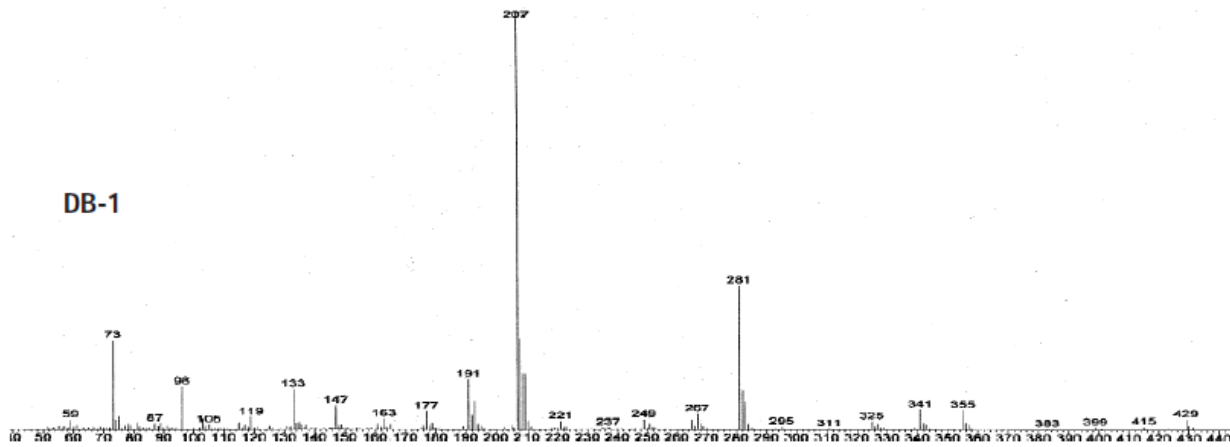
- Carry-over from a backflash or previous sample
- Bad tank of gas or traps have expired
- Septum bleed

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

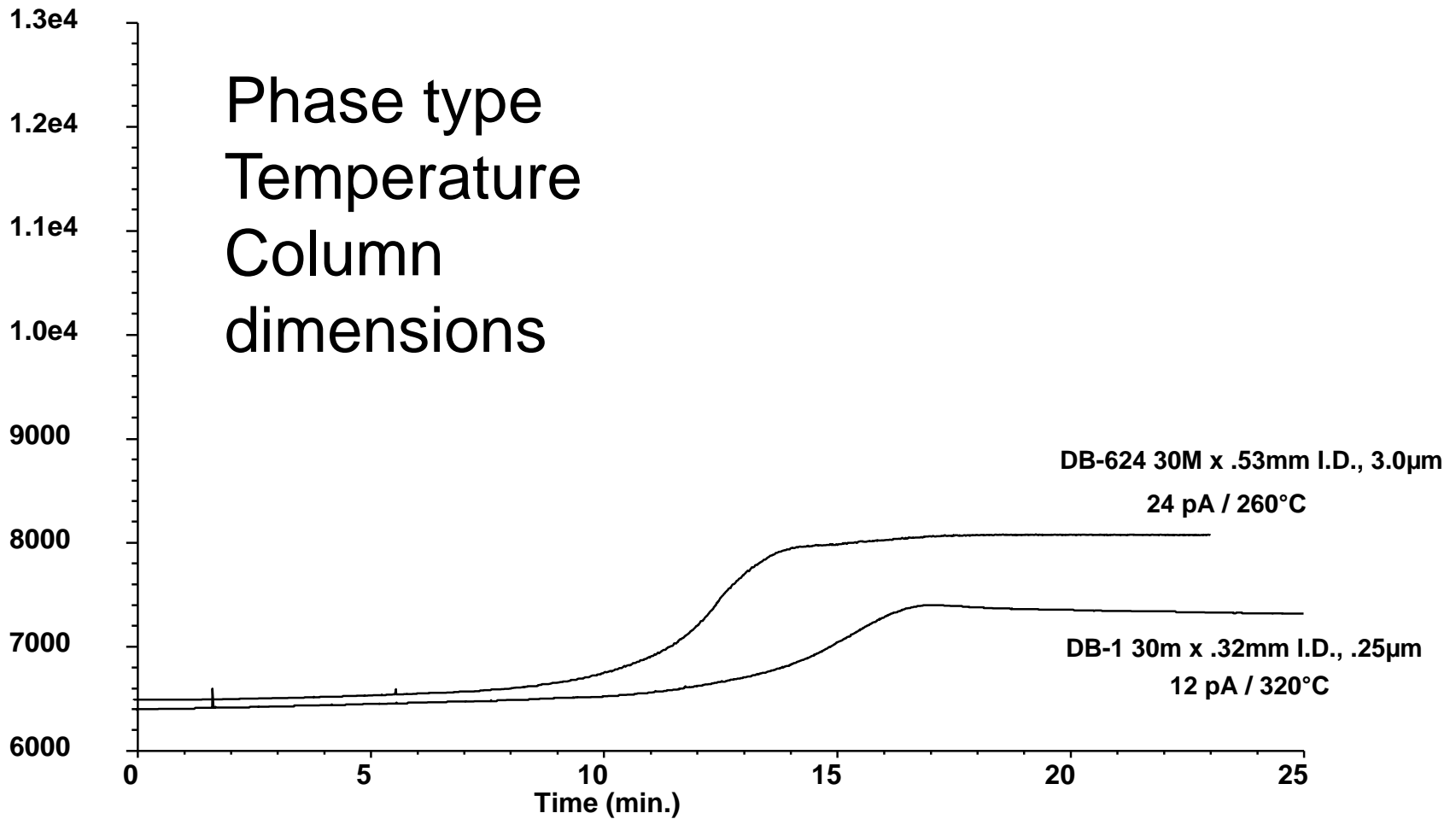
Bonus 'Siloxane' Peaks



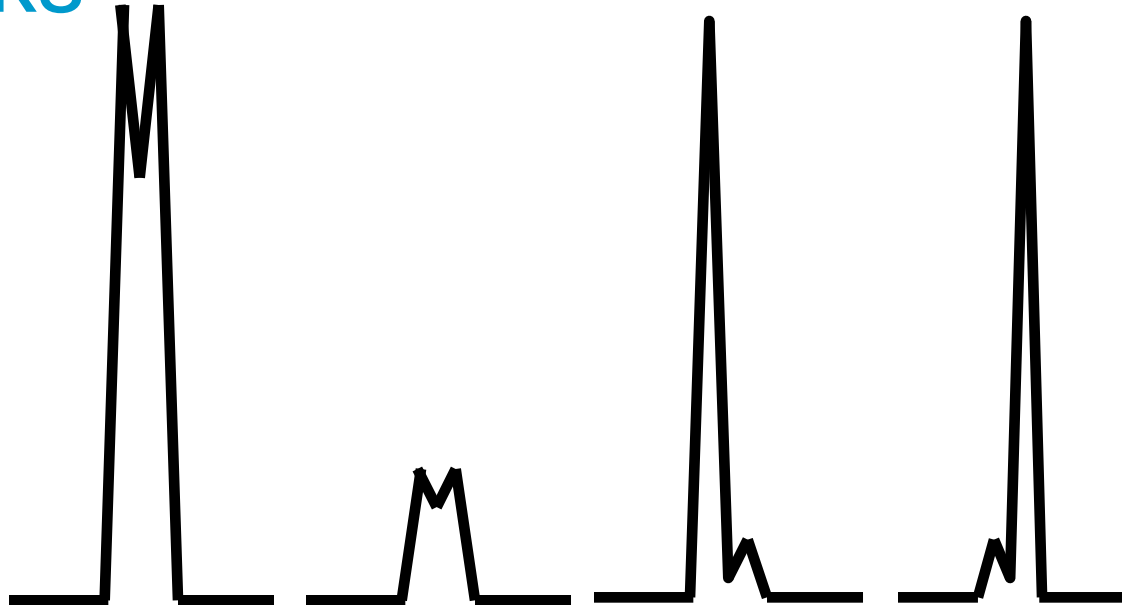
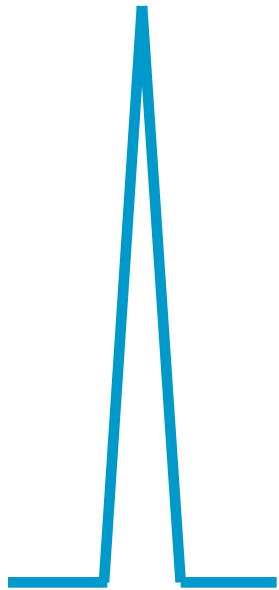
GC Column Bleed Ions



Column Bleed is Influenced by:



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

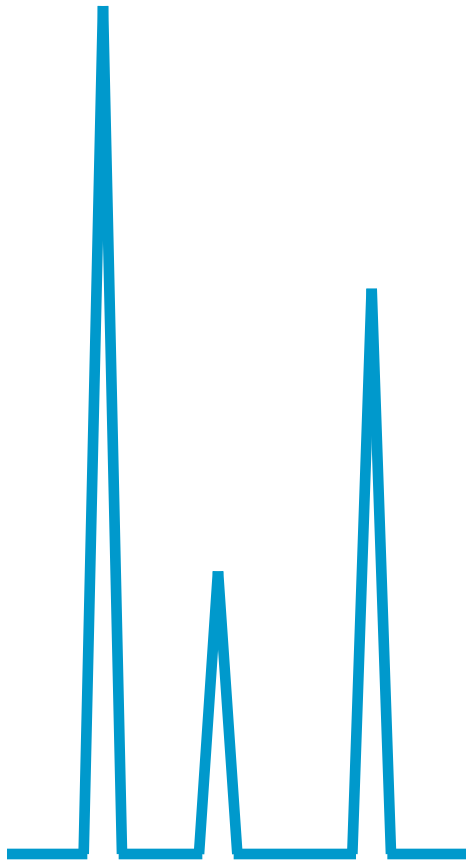
VOLATILITY

High boilers dropping out on Cold Spots

-Transfer line temps

-Unions or fittings not tracking column temp

No Peaks



DETECTOR (not on or not operational)

INJECTOR (not working)

-Plugged syringe/plunger not moving

-Wrong injector (or detector)

-Huge leak (older systems)

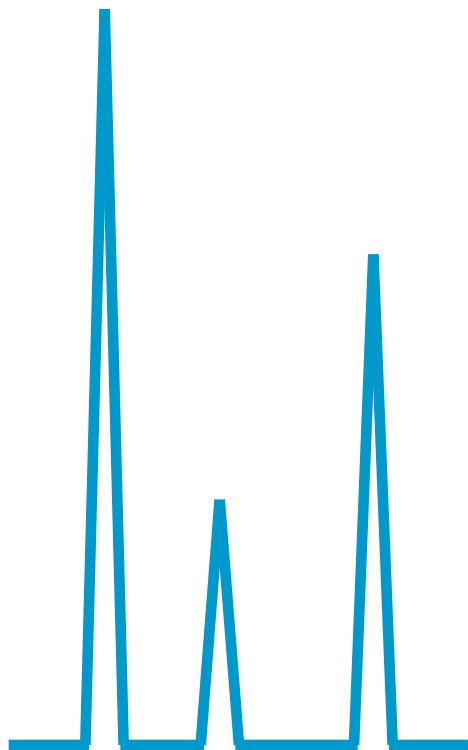
-No carrier gas flow

NOT the COLUMN Unless...

-Broken column or No column

Peak Response

All Change in Size

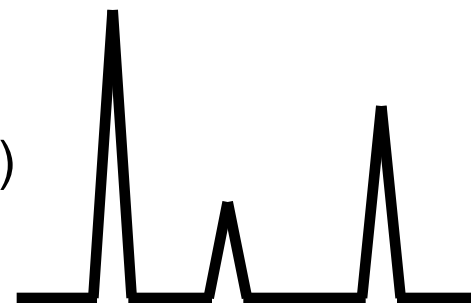


INJECTOR

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

DETECTOR (response problem)

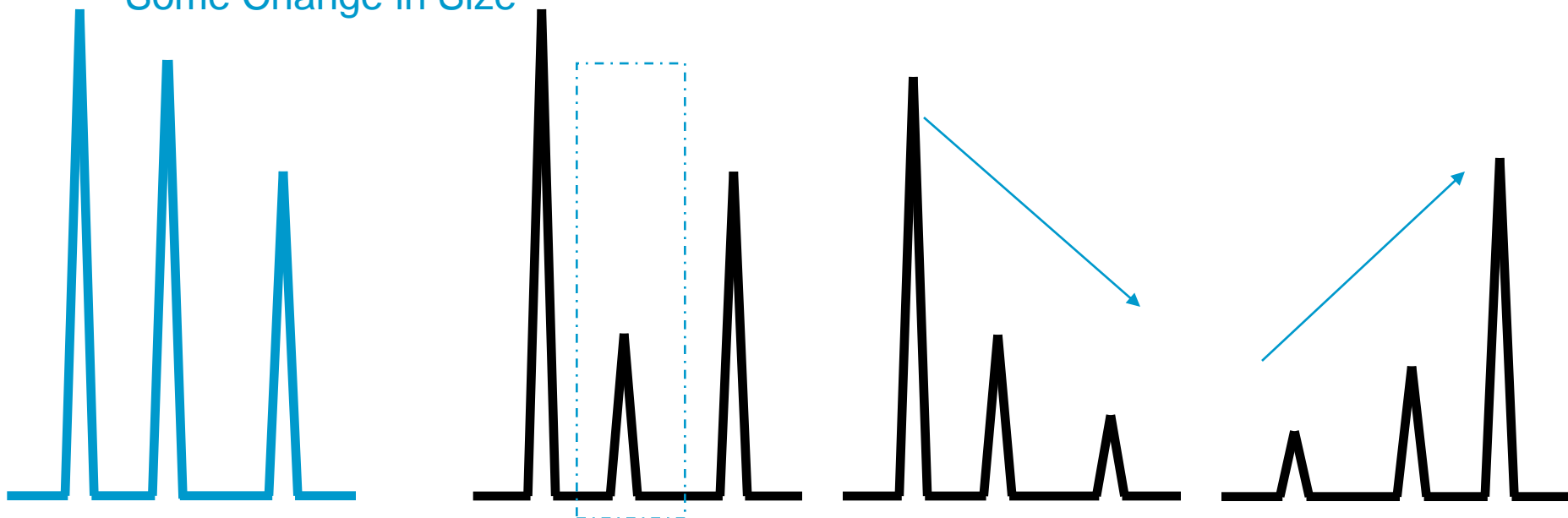
- Settings or flows changed
- Electronics failing



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

Peak Response

Some Change in Size



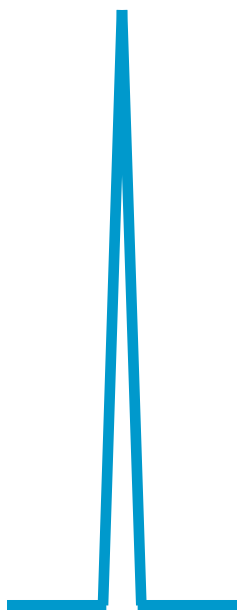
INJECTOR or COLUMN is active/contaminated

- 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.

Peak Fronting

Shark Fin Shaped or Just Slight



COLUMN (contaminated)

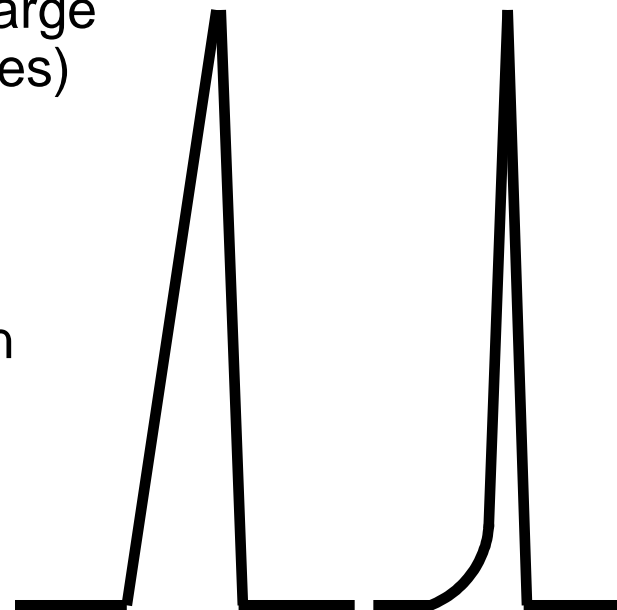
- Overload (More pronounced with large solute and phase polarity differences)

INJECTOR

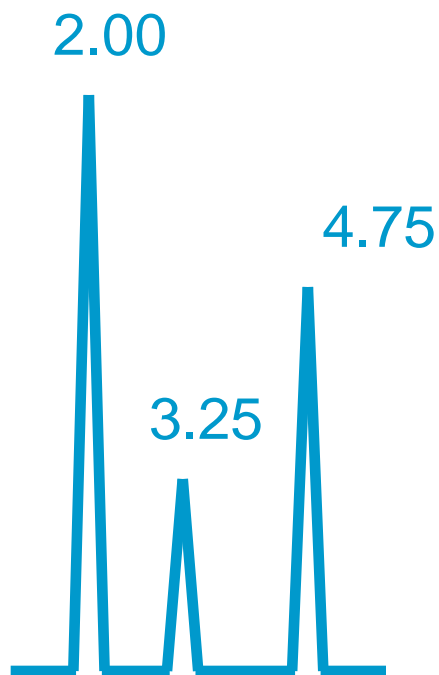
- Column installation
- Compound very soluble in injection solvent (need retention gap)
- Mixed sample solvent

OTHER

- Co-elution
- Breakdown



Retention Time Shift



INJECTOR

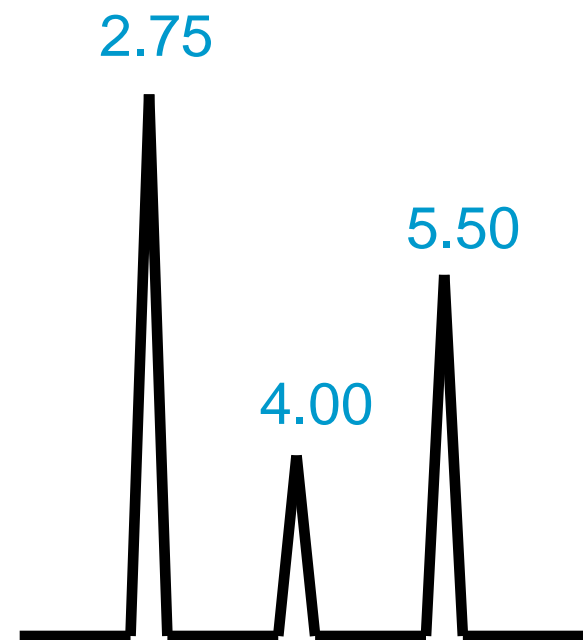
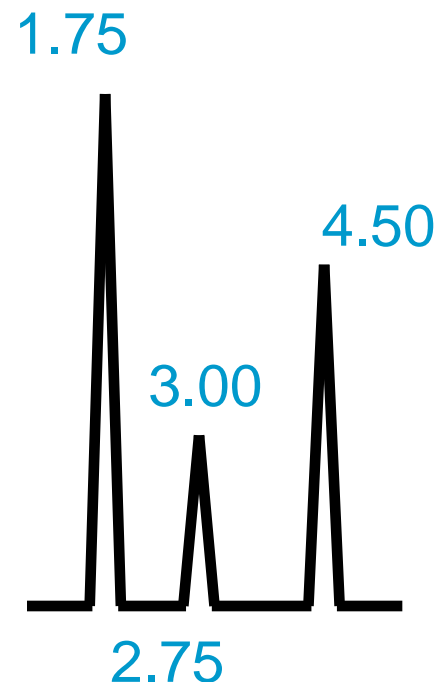
- Leak in the septum
- Change in injection solvent
- Large change in sample concentration

FLOW

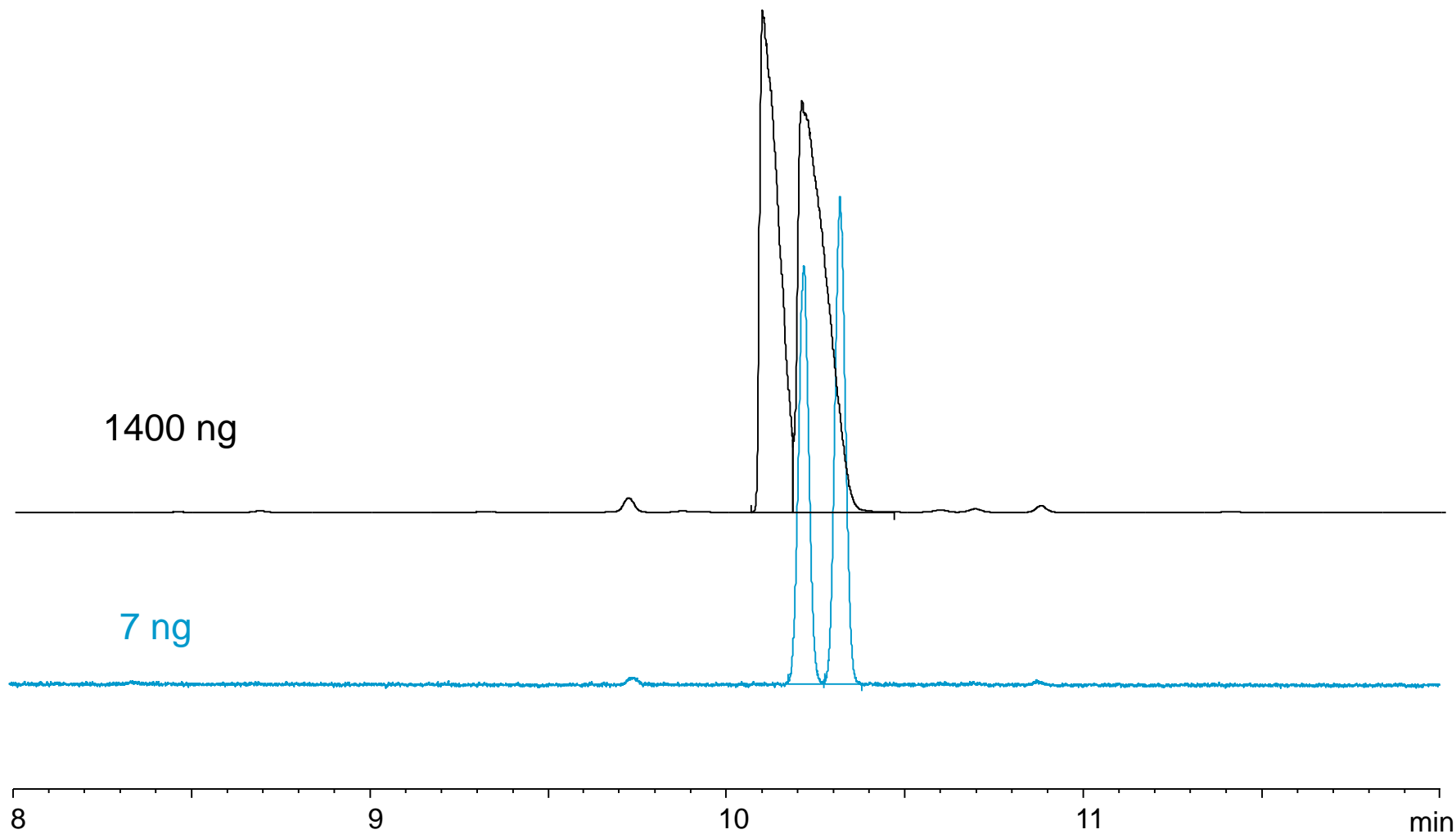
- Change in gas velocity

COLUMN

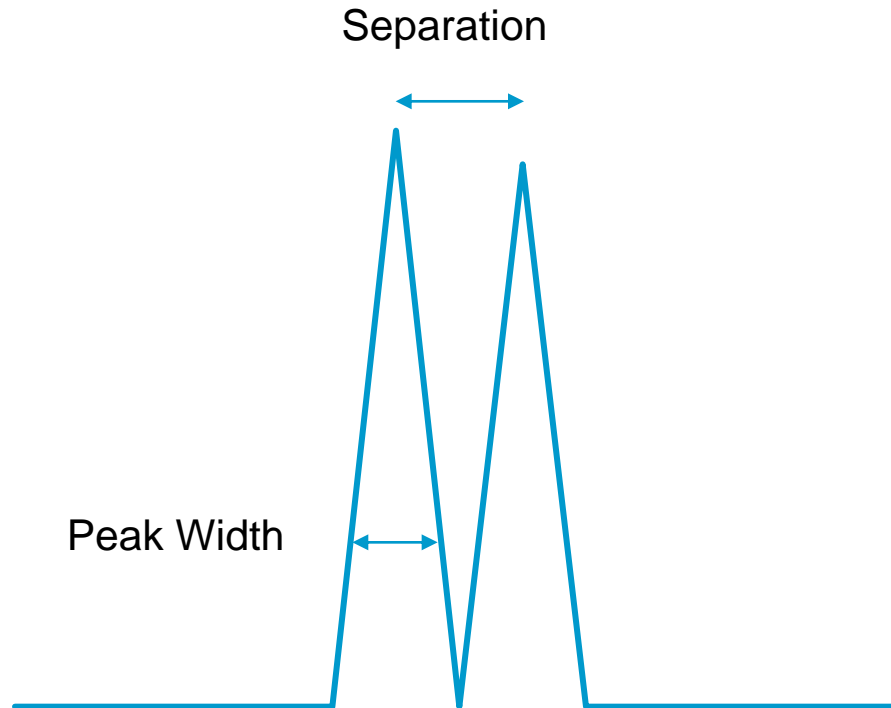
- Contamination
- Damaged stationary phase
- Loss of stationary phase
- Change in temperature



Effect of Sample Overload on Retention Time and Peak Shape



Loss of Resolution

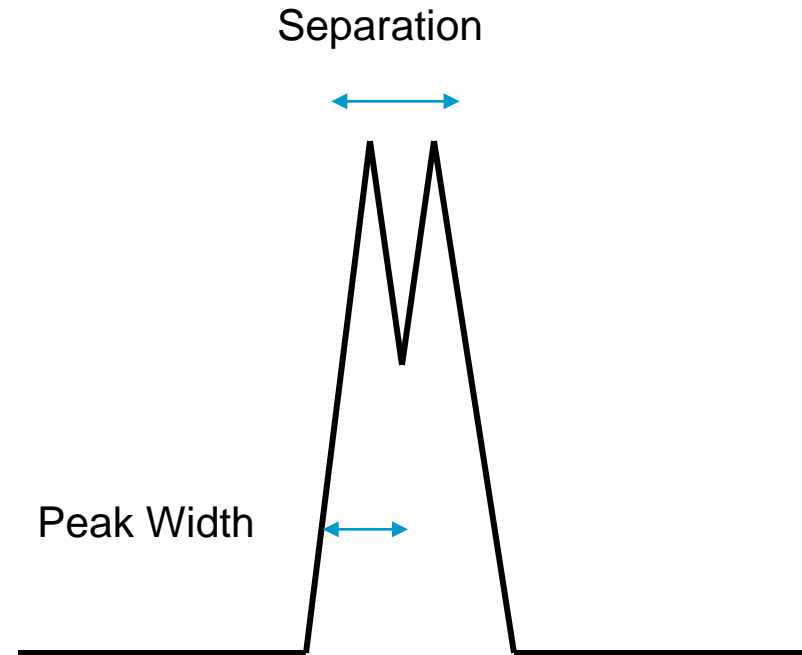


Resolution is a function of separation and peak width

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

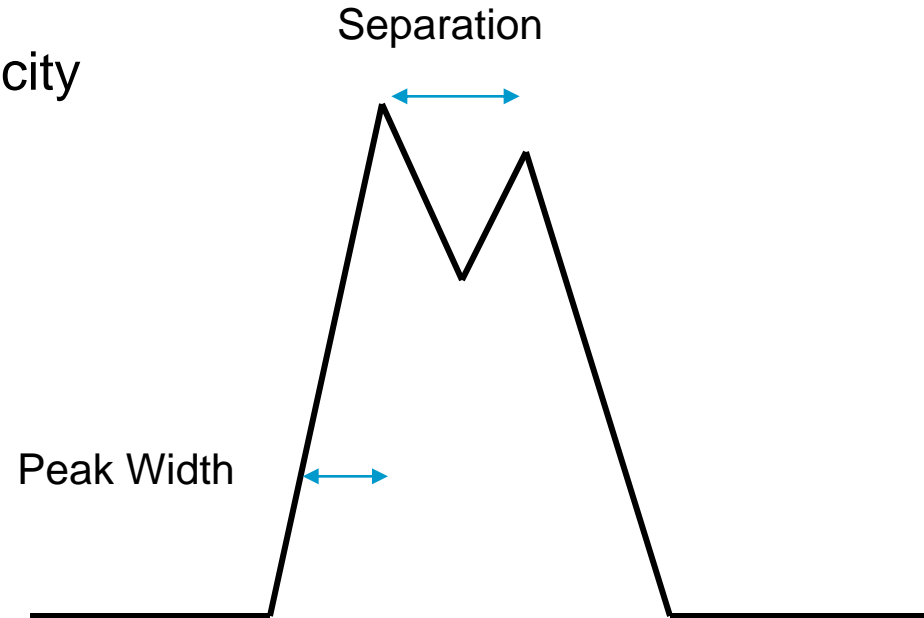
- Change in carrier gas velocity
- Make-up gas

COLUMN

- Contamination
- Phase degradation

INJECTOR (efficiency)

- Settings, Liner, Installation, etc.



Baseline Disturbances

Sudden Changes, Wandering, or Drifting

WANDER



COLUMN or DETECTOR

-Not fully conditioned or stabilized (electronics)

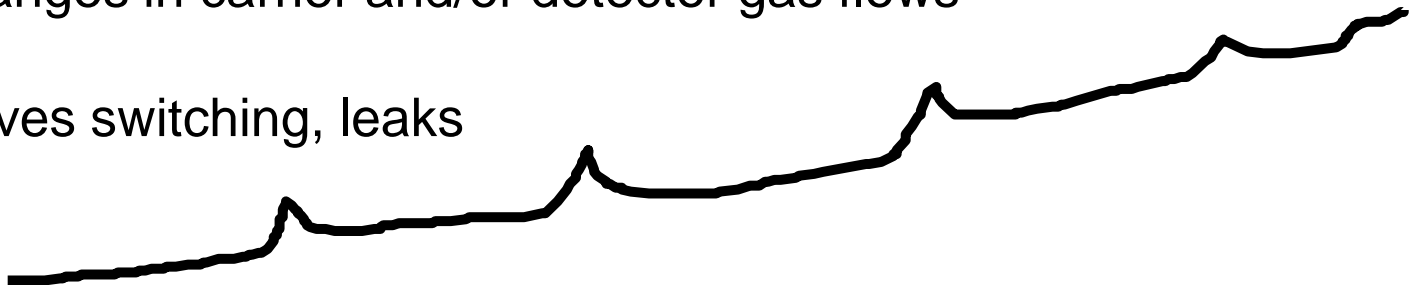
-Contamination

FLOW

-Changes in carrier and/or detector gas flows

-Valves switching, leaks

DRIFT



Noisy Baseline

MILD



SEVERE



FLOW

- Contaminated gas
- Incorrect detector settings

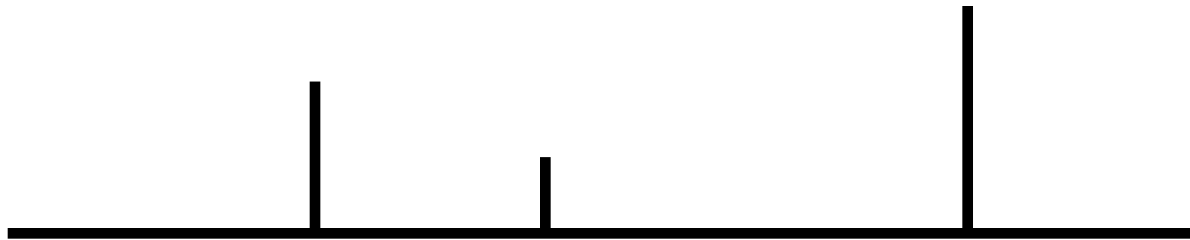
COLUMN

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

DETECTOR

- Air leak - ECD, TCD
- Electronics malfunction

Spiking Baseline



DETECTOR

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

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

Troubleshooting “Tools”

Bleed Profile: *baseline problems*

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

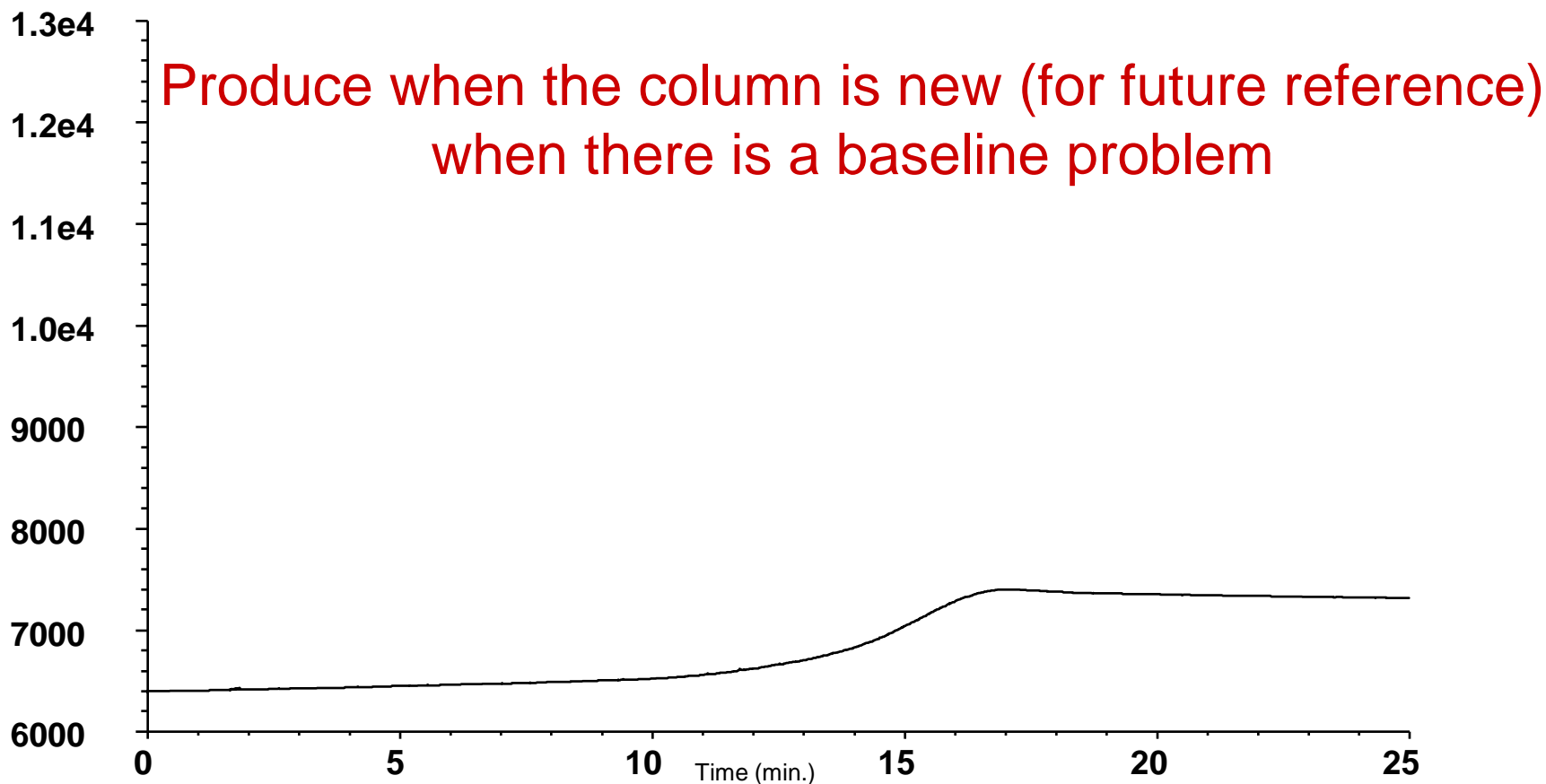
Test mix: *all problems*

Isolate the components: *all problems*

Condensation Test: *baseline problems*

Jumper Tube Test: *baseline problems*

Generating a Bleed Profile

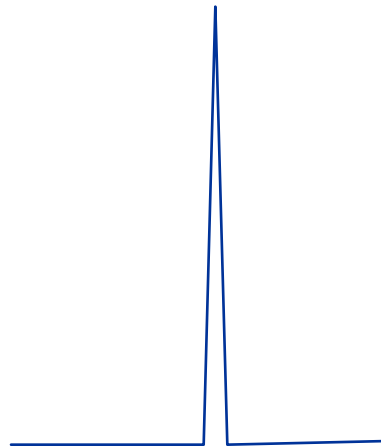


*DB-1 30m x .32mm I.D., .25 μ m

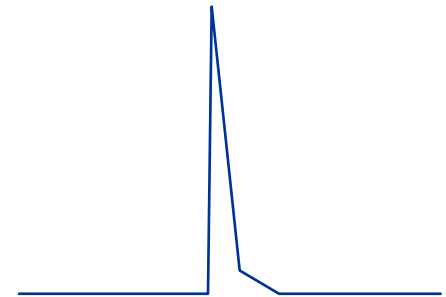
Temperature program // 40°C, hold 1 min // 20°/min to 320°C, hold 10 min.

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

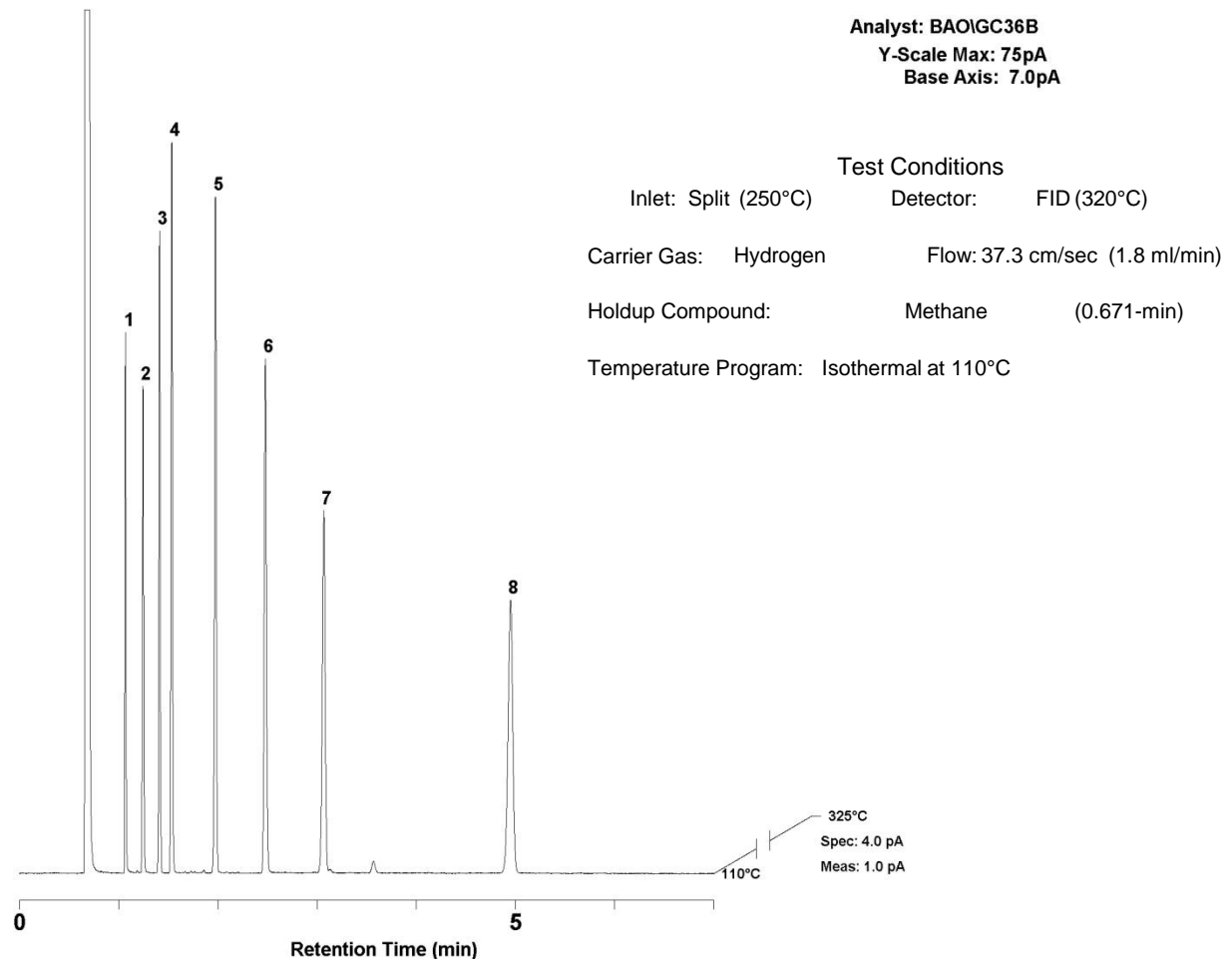
Liner problem

(broken, leaking, misplaced)

Column position in injector and detector

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 to your application
- Selective detectors
- Concentrations specific to your application
- Use same instrument conditions
- Easiest to simply inject a calibration standard
- Store for future measure of column performance

Isolate the Components

Simplify the system:

- example - Direct injection instead of P&T sample introduction

Put in a known good column

Move column to a different GC, inlet or detector

Condensation Test



Used* to isolate the cause of:

- Erratic baselines
- Ghost peaks or carryover

***Use when problems are worse after periods of GC non-use**

Condensation Test

Procedure

Leave GC at 40-50°C for > 8 hours

Blank run

Repeat a blank run immediately after the first blank run is complete

Compare the two blank runs

Condensation Test

Results

First blank run is worse:

- Contaminants (from injector, lines, traps or carrier gas) carried into the column

Blank runs the same: *contaminants are not strongly focused on the front of the column*

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

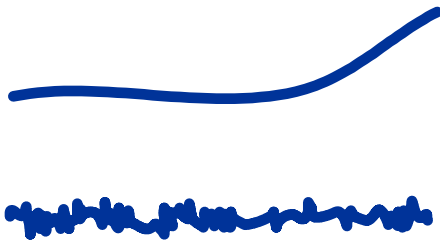
Blank run

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

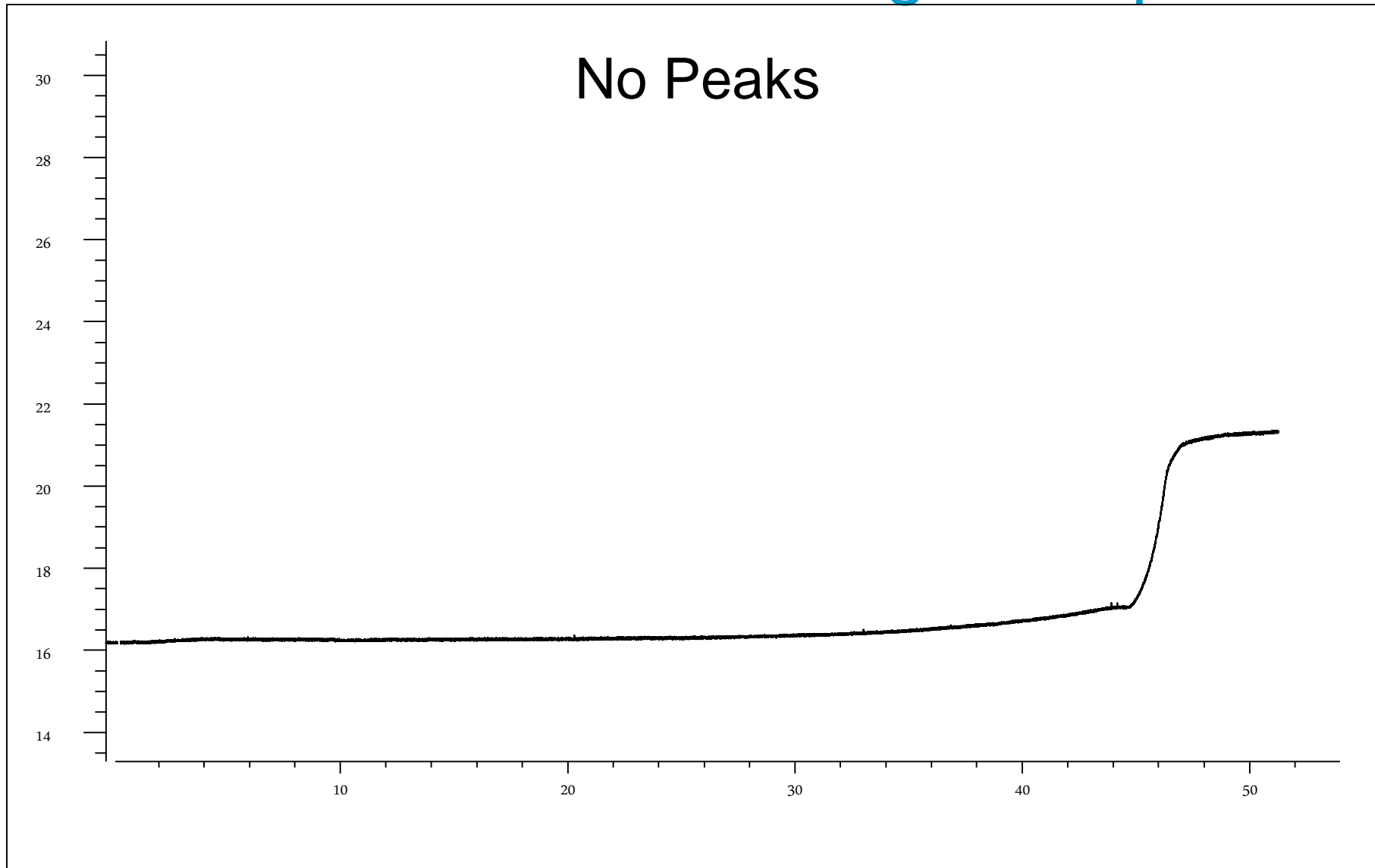
Problem gone: Previous leak, solid debris, or installation problem

And Now Let's do Some

TROUBLESHOOTING

Troubleshooting-Example #1

A Real Troubleshooting Example



Logical Steps Taken to Find Peaks

(most of our problems are leaks and plugs)

Is the flame Lit?

- put glass piece over FID outlet----*Answer in this case, Water condenses*
- look at output in instrument guage-- is the digital value greater than 0.0?
Answer in this case is approximately 16.2 pico amps

Is there flow through the column?

- disconnect column from detector and measure flow with bubble solution or meter
Answer in this case was YES THERE IS FLOW

Assess the observations

- *Flame is lit and we have flow from end of column*
- *Hypothesis: Sample not getting on column-syringe plugged?*

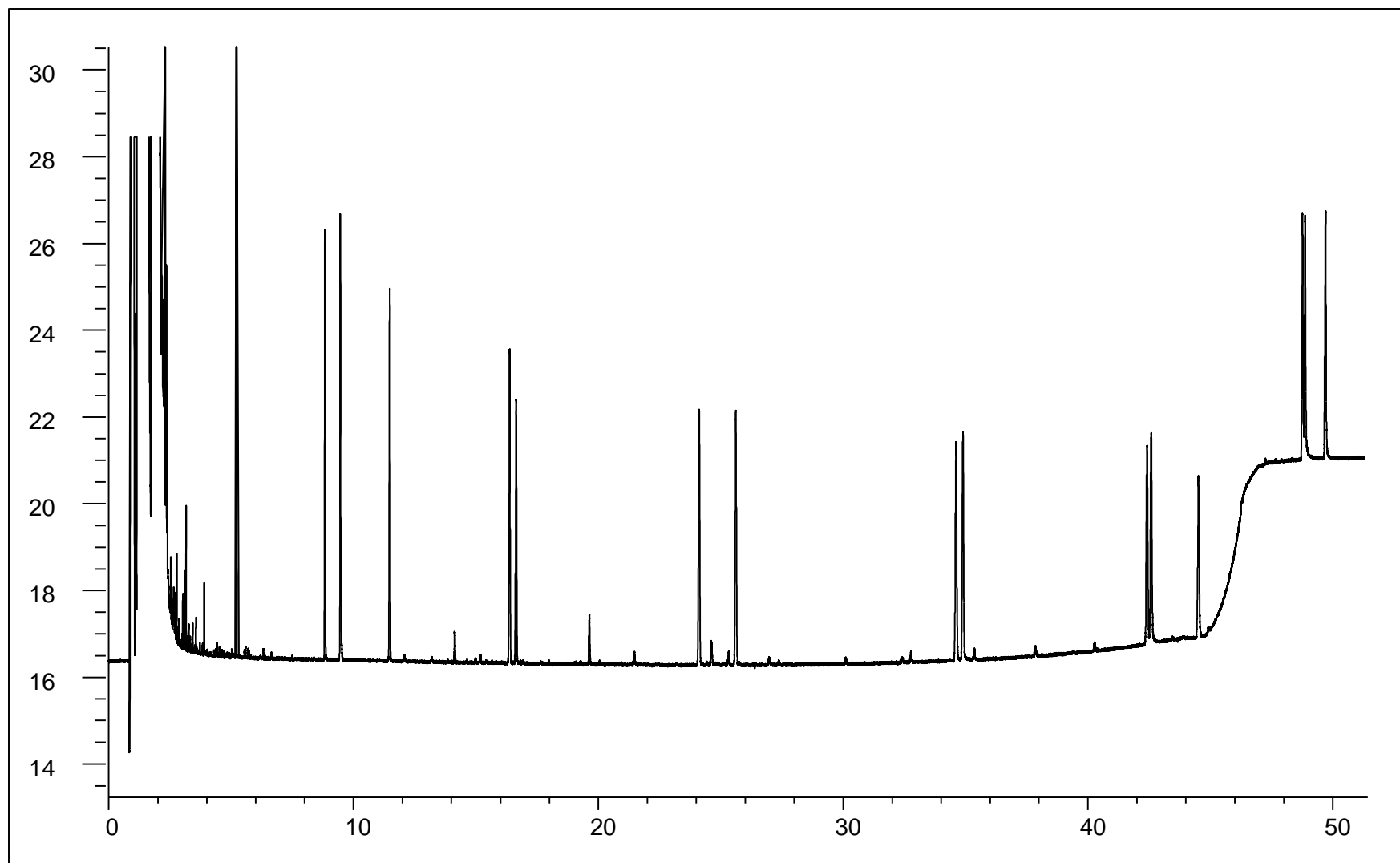
Take syringe out and make injection manually on a dry paper towel

Answer – towel stays dry (Syringe was clogged with septum)

Pull plunger out top, add solvent and replace plunger will usually dislodge septum particle (should hear a little pop) If you can't dislodge plug, Replace syringe

Reassemble the Injector & Re-inject

Peaks !!

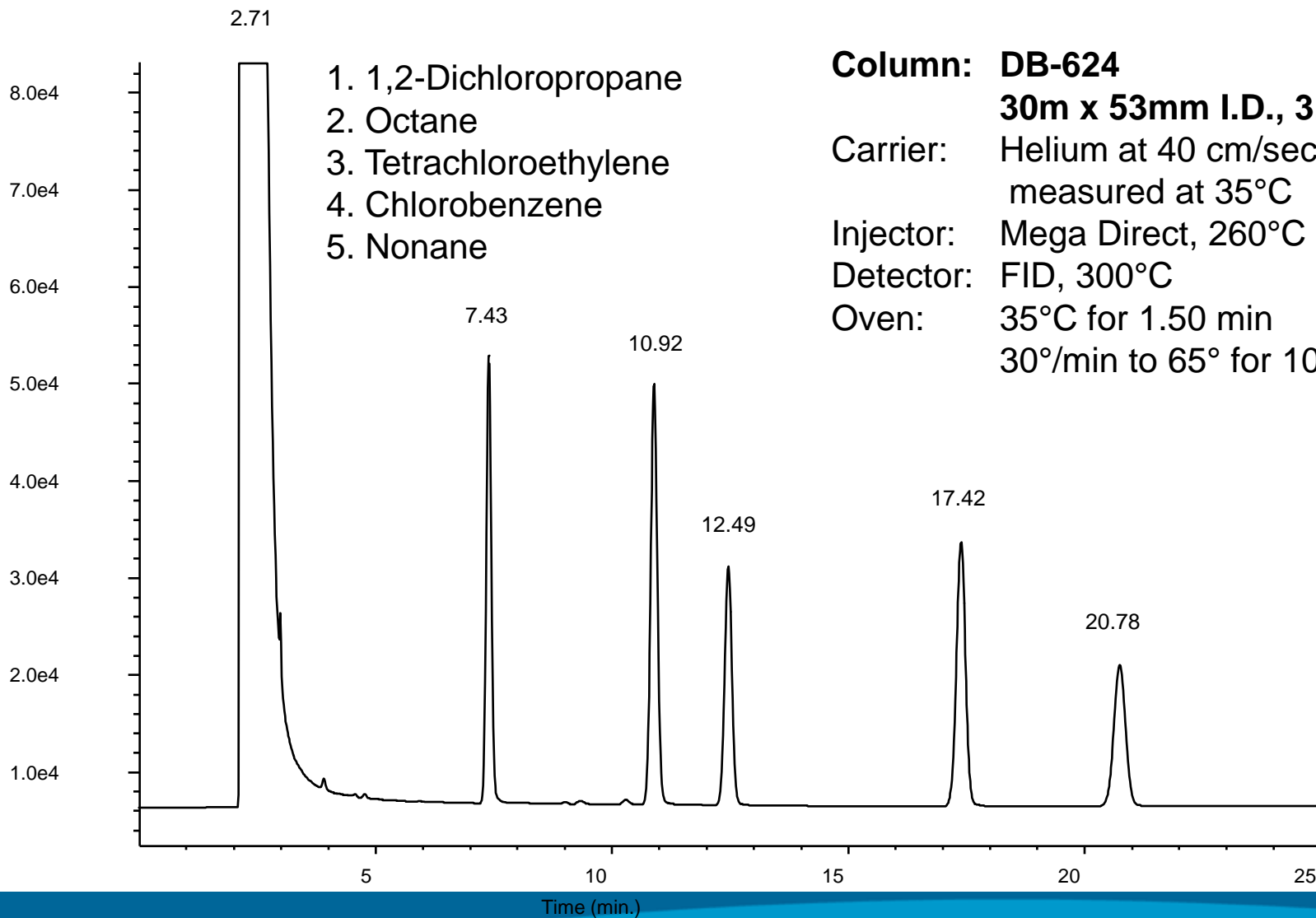


Troubleshooting-Example #2

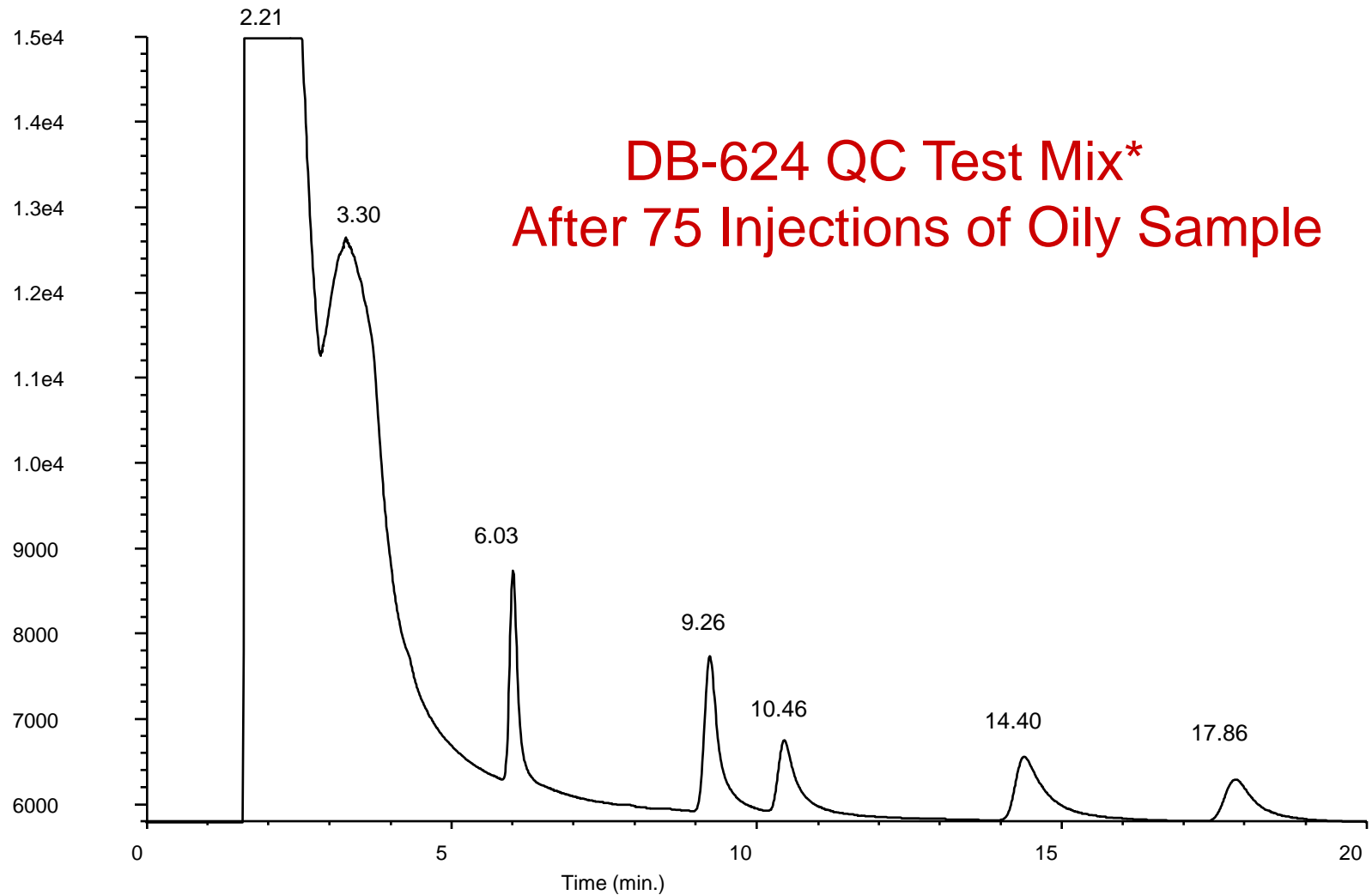
DB-624 COLUMN

QC Test Mix

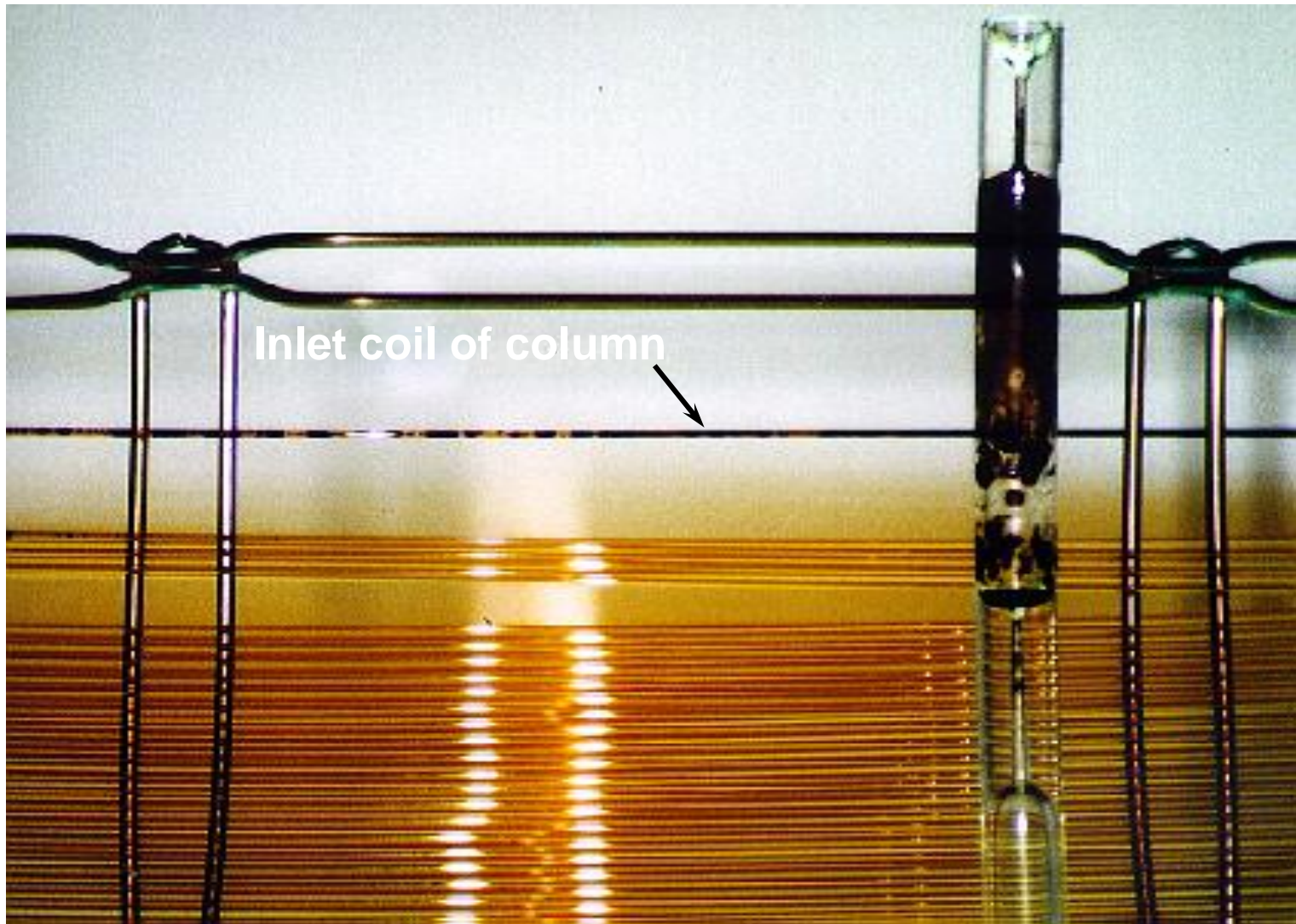
Column: DB-624
30m x 53mm I.D., 3.0µm
Carrier: Helium at 40 cm/sec
measured at 35°C
Injector: Mega Direct, 260°C
Detector: FID, 300°C
Oven: 35°C for 1.50 min
30°/min to 65° for 10 min



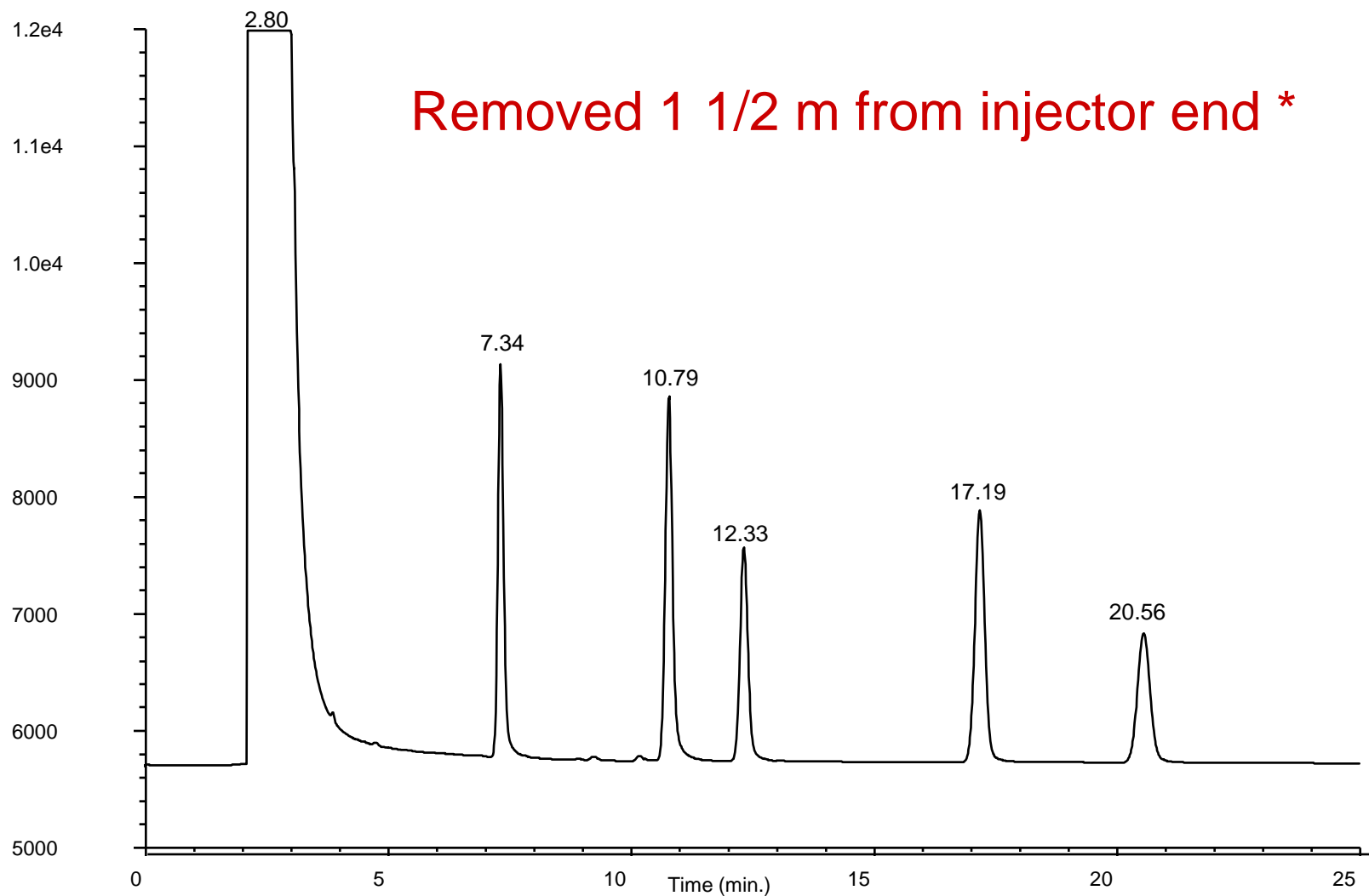
Example of Column Contamination



Column and Liner Contamination



Example of Column Contamination



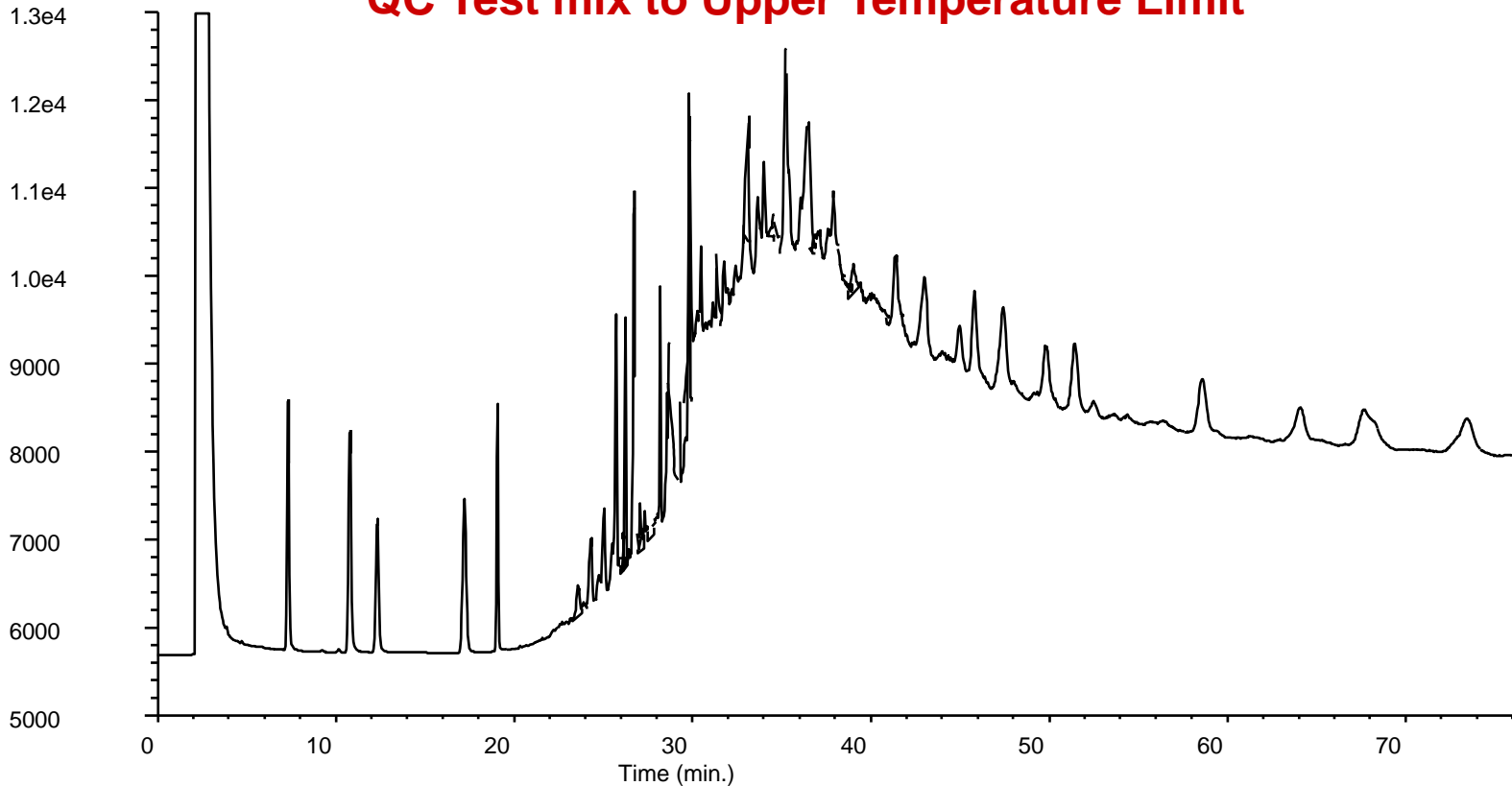
*Before Column rinse and bake

Temperature program // 35°C hold 1.50 min // 30°/min to 65°C, hold 10 min

Looks Fixed Doesn't it?

Example of Column Contamination

1 1/2 mtrs removed*
QC Test mix to Upper Temperature Limit

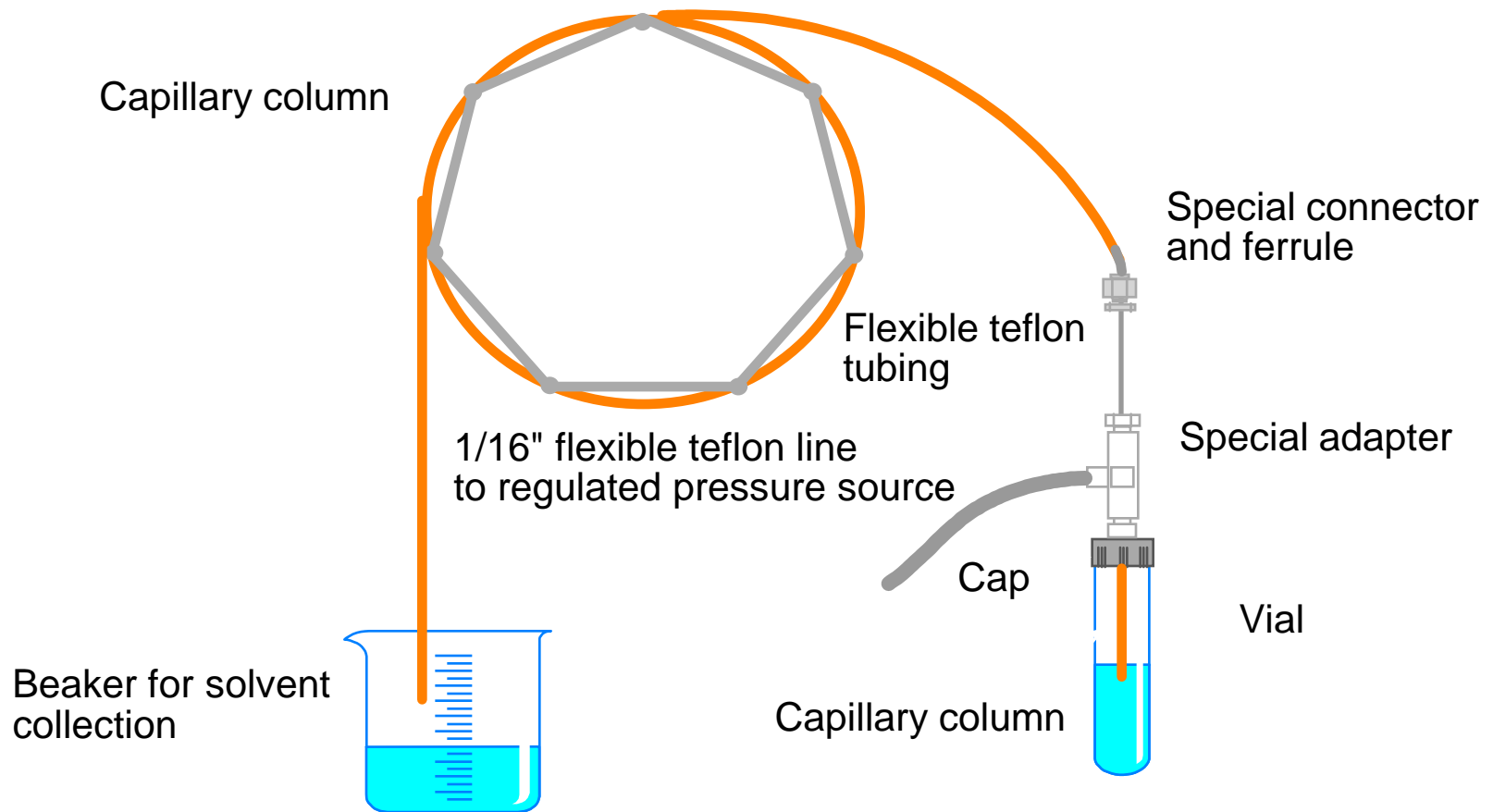


***Before Column rinse and bake.**

**Temperature program // 35°C, hold 1.50 min // 30°/min to 65°C,
hold 15 min // 20°/min to 260°, hold 50 min**

Backflush Column

Rinse with 10ml each:
Methanol, Methylene Chloride, Hexane

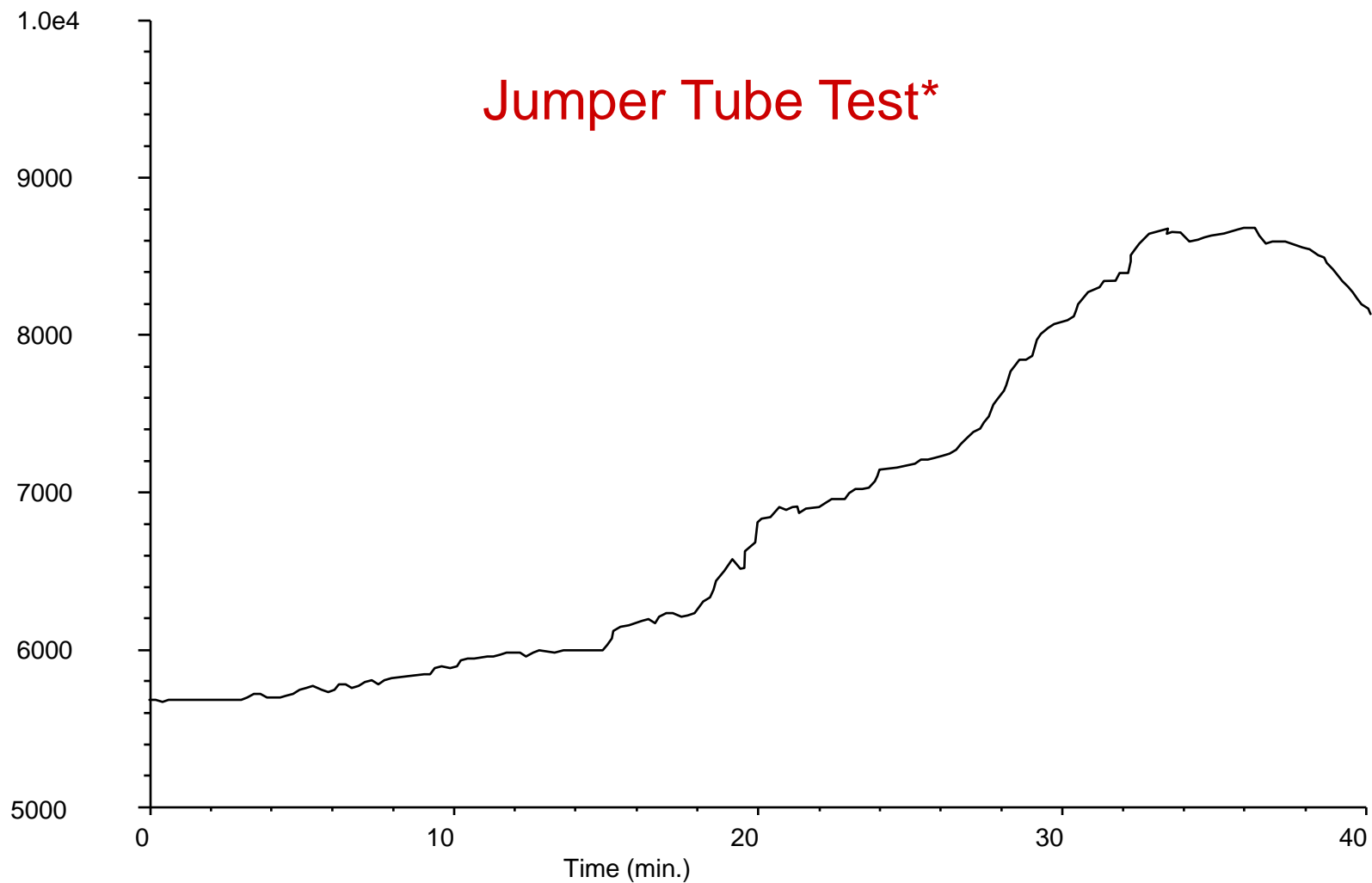


Jumper Tube Test

Used to Isolate Source of Contamination

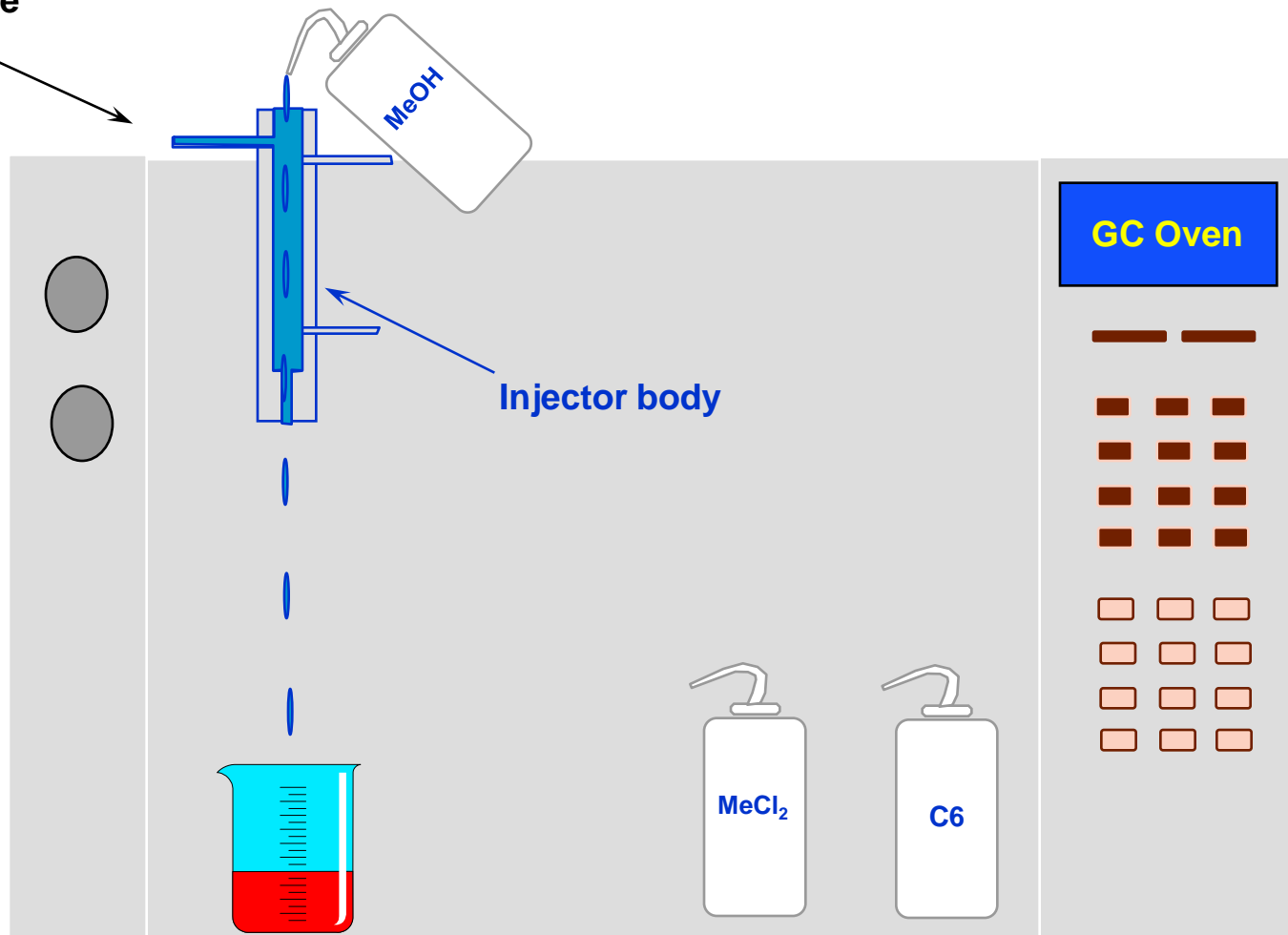
- . Cap off the detector and establish normal gas flows and temperature.
- . Plot the baseline using a temperature program. If flat.....
- . Connect 1 meter of deactivated tubing between the injector and detector
- . Plot the baseline using a temperature program. If flat.....
- . Install the column.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

Contaminated Inlet



Rinsing Injector

Carrier gas line



Troubleshooting Tips

1. Isolate the problem.

(Blank Run, Inject Un-retained Compound, Jumper Tube Test)

2. Change only one variable at a time.

3. Compare before/after chromatograms.

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

4. Utilize Technical Support.

Remember

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

Multiple cause and effect

Do not change too many variables at once

TECHNICAL SUPPORT

1-800-227-9770, #3, #3, #1

E-mail: gc-column-support@agilent.com

866-422-5571 (FAX)

