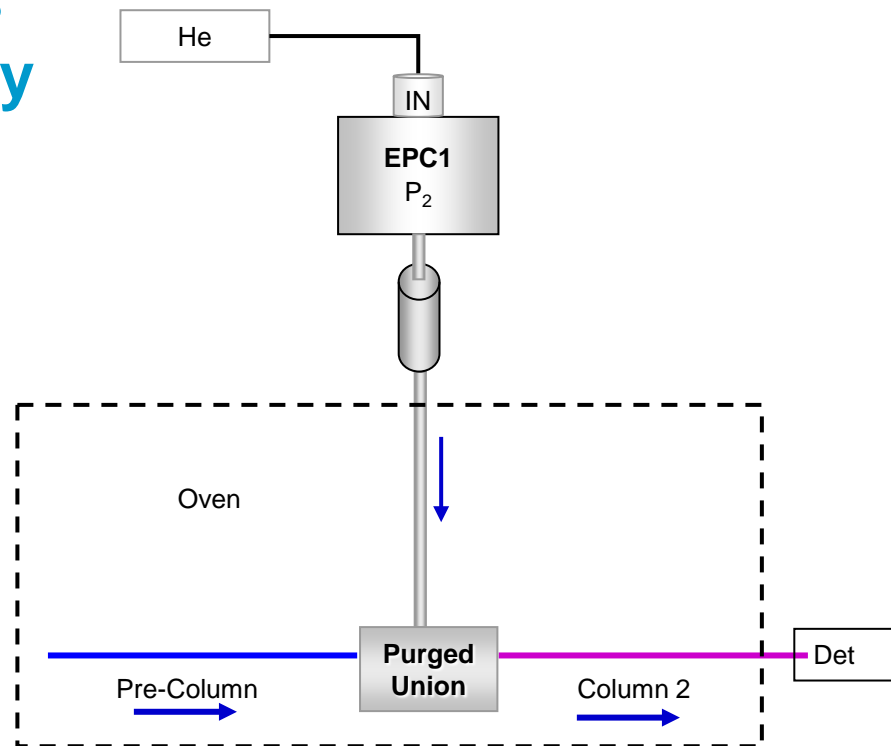


# Reliable Analysis of the Light Ends of Petroleum Fractions and Crude Oil Using Capillary Flow Backflush Technology

Roger L. Firor  
Agilent Technologies  
Chemical Analysis Group  
Wilmington, DE USA



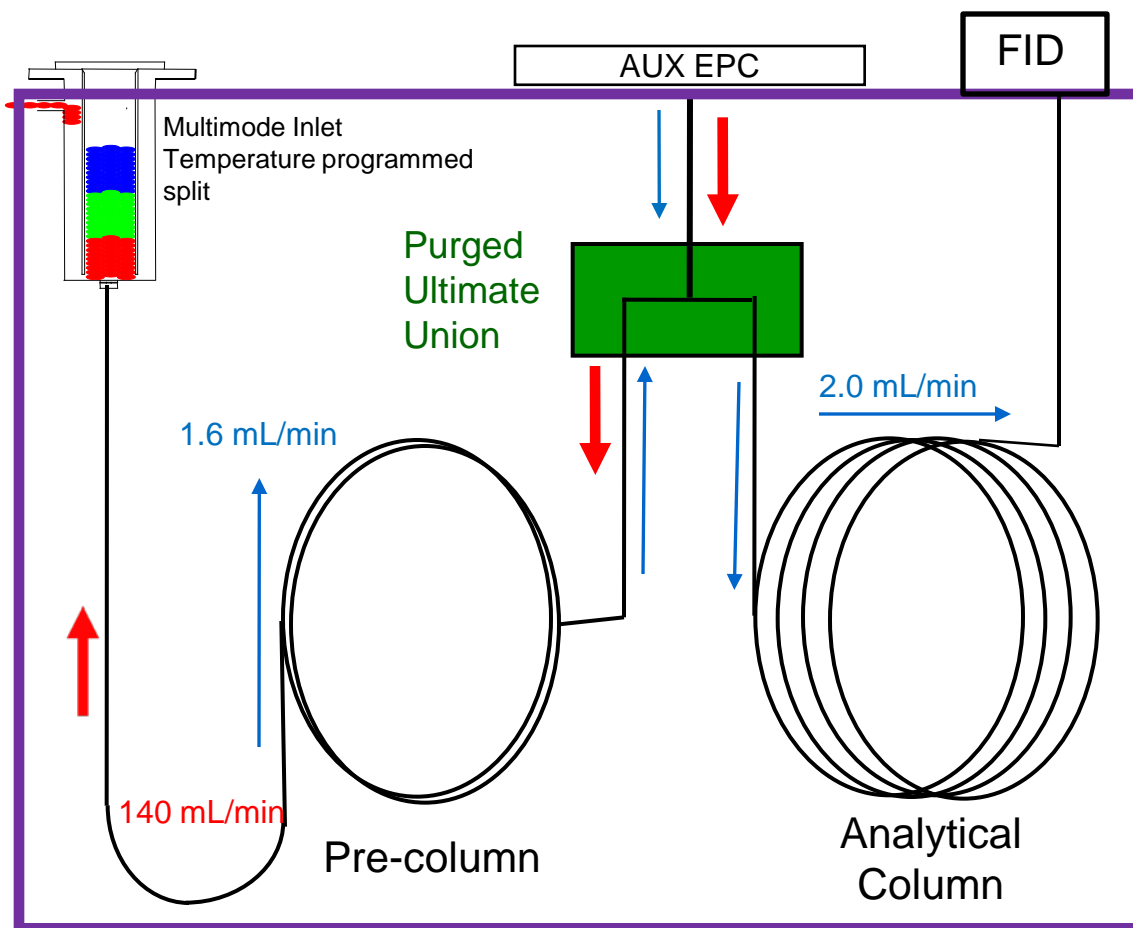
# Outline

- Why pre-column backflush
- System configuration with CFT (Capillary Flow Technology)
- Typical method setup
- Performance
- Application examples

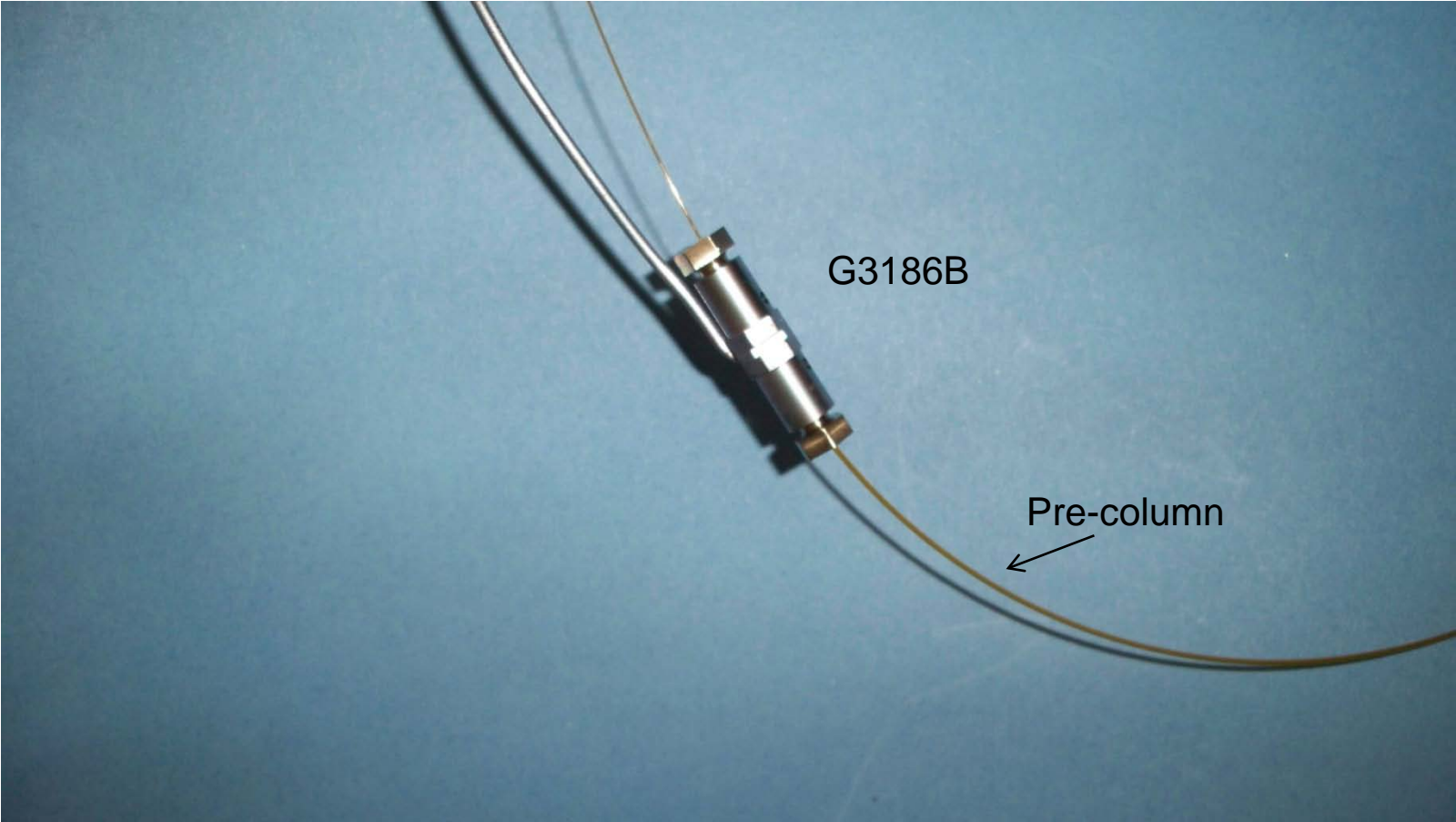
# Pre-column Backflush

- Eliminate bake out times at end of runs
- Dirty samples
- Late eluting compounds of no analytical interest
- Need to protect column from high boilers
- Interest in maximizing lab throughput and minimizing cost of operations
- Compatible with GC and GC/MSD systems

# Basic System Configuration



# Purged Union with Columns Installed



# Advantages of Pre-column Backflush

- Backflush is carried out while analysis is occurring on the analytical column
- Midpoint pressure control allows analytical column to run at desired flow while pre-column is backflushed
- Use of an uncoated pre-column allows transfer of components (in this application up to C12 – C13) at a relatively low temperature (i.e. 35 C) and faster backflushing of the heavier components
- Sample matrix has little influence
- Wide choice of pre-column and analytical column dimensions and phase ratios possible to meet specific application needs

# System Setup and Operation

- A pre-column, typically a deactivated retention gap of 2 meter length is used.
- The inlet of the pre-column is defined as the MMI (multimode inlet) and the outlet as a Aux module channel
- The analytical column (100m x 0.20mm x 0.5um HP1 for crude oil analysis) uses an Aux channel (no restrictor) as inlet and FID as outlet
- At the specified runtime the inlet pressure is dropped to reverse flow in the pre-column while the Aux maintains forward flow in the analytical column
- At backflush runtime, the MMI's temperature is ramped to 425 C to clear out the inlet. It is held at this temp to end of run
- Single taper liner with glass wool is used
- Pressure zeros should be set on the GC for the inlet and Aux channel before use

# Backflush: Column Settings for Polywax Carbon Number Specific Backflushing

PW 500

Pre-column →

Analytical column →

Description	
1	agilent RG-1: 450 °C: 1 m x 530 μm x 0 μm In: Back PTV Inlet He Out: Aux EPC 1 He
2	J&W 145-1001: 400 °C: 5 m x 530 μm x 0.15 μm In: Aux EPC 1 He Out: Front Detector FID
	Front SS Inlet He
	Aux EPC 2 He
	Aux EPC 3 He

Control Mode  On

Flow      12 mL/min  
 Pressure      1.5871 psi  
 Average Velocity      93.246 cm/sec  
 Holdup Time      0.089369 min

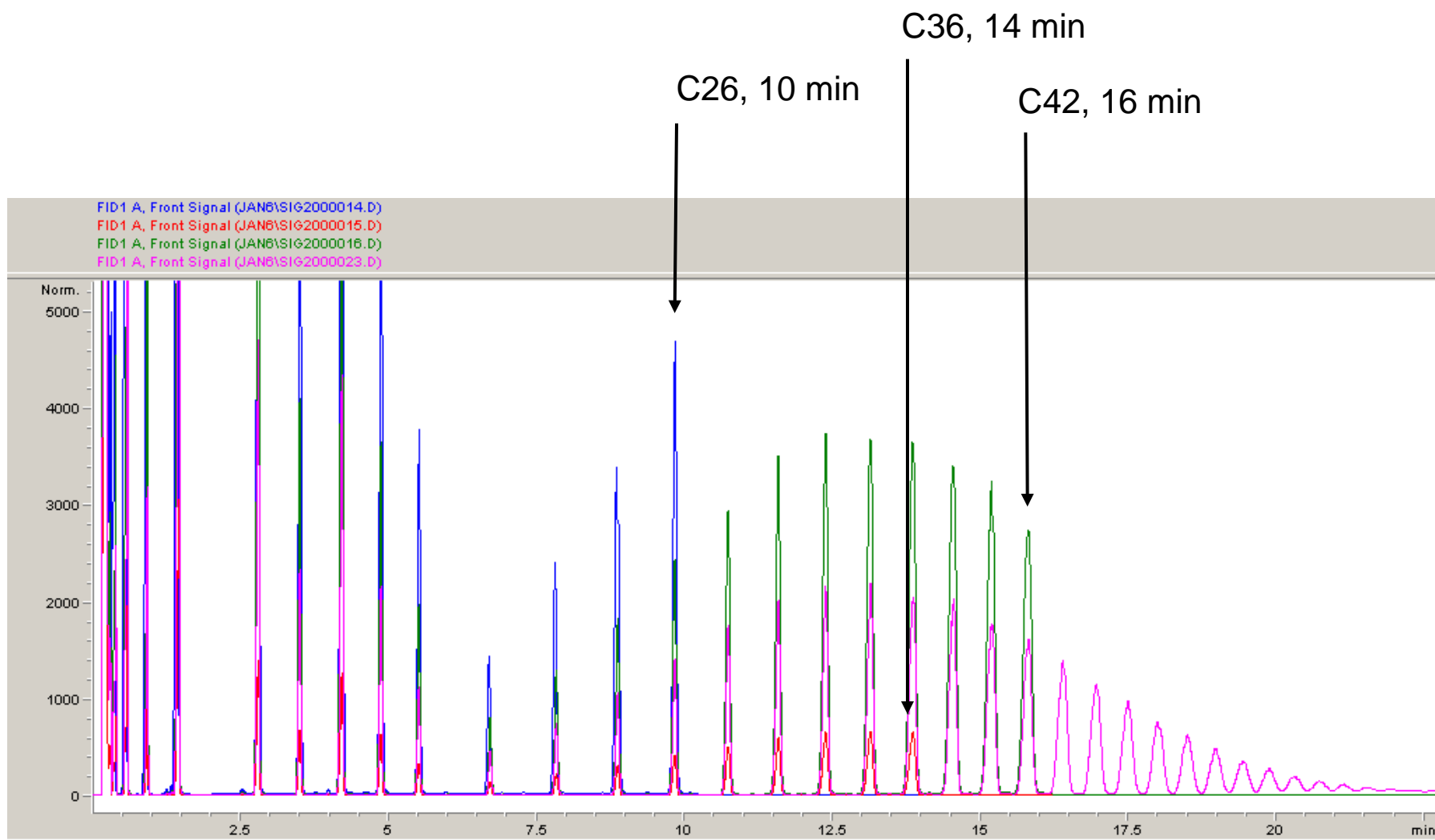
(Initial): 0 min  
He @ 50 °C Oven  
Out: Ambient Pressure  
5 m x 530 μm x 0.15 μm

	Rate mL/min per min	Value mL/min	Hold Time min	Run Time min
▶ (Initial)		12	0	28.333
*				

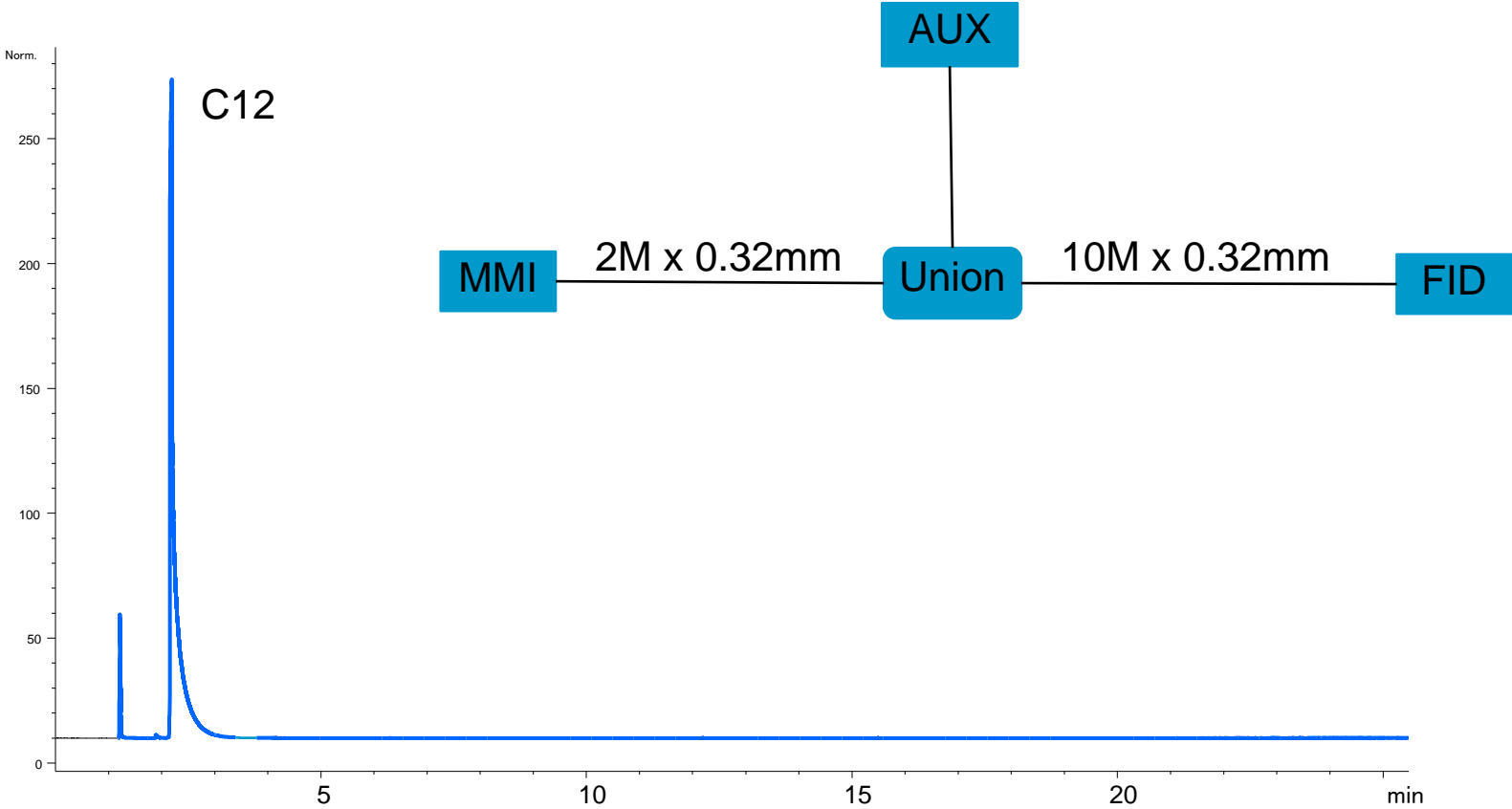


# Backflush with Purged Tee: Polywax 500

Sample: C5-C18 plus PW500

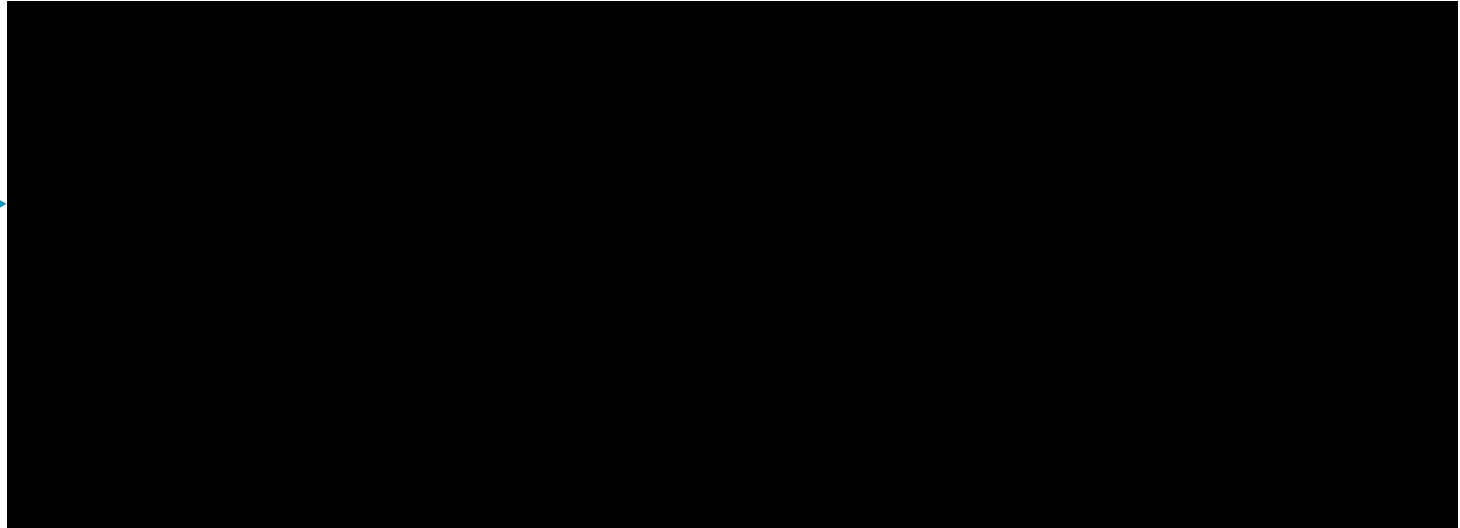


# Peak Shapes on Retention Gaps: C12 Only



# Setup for Pre-column and Analytical Column

Pre-column →

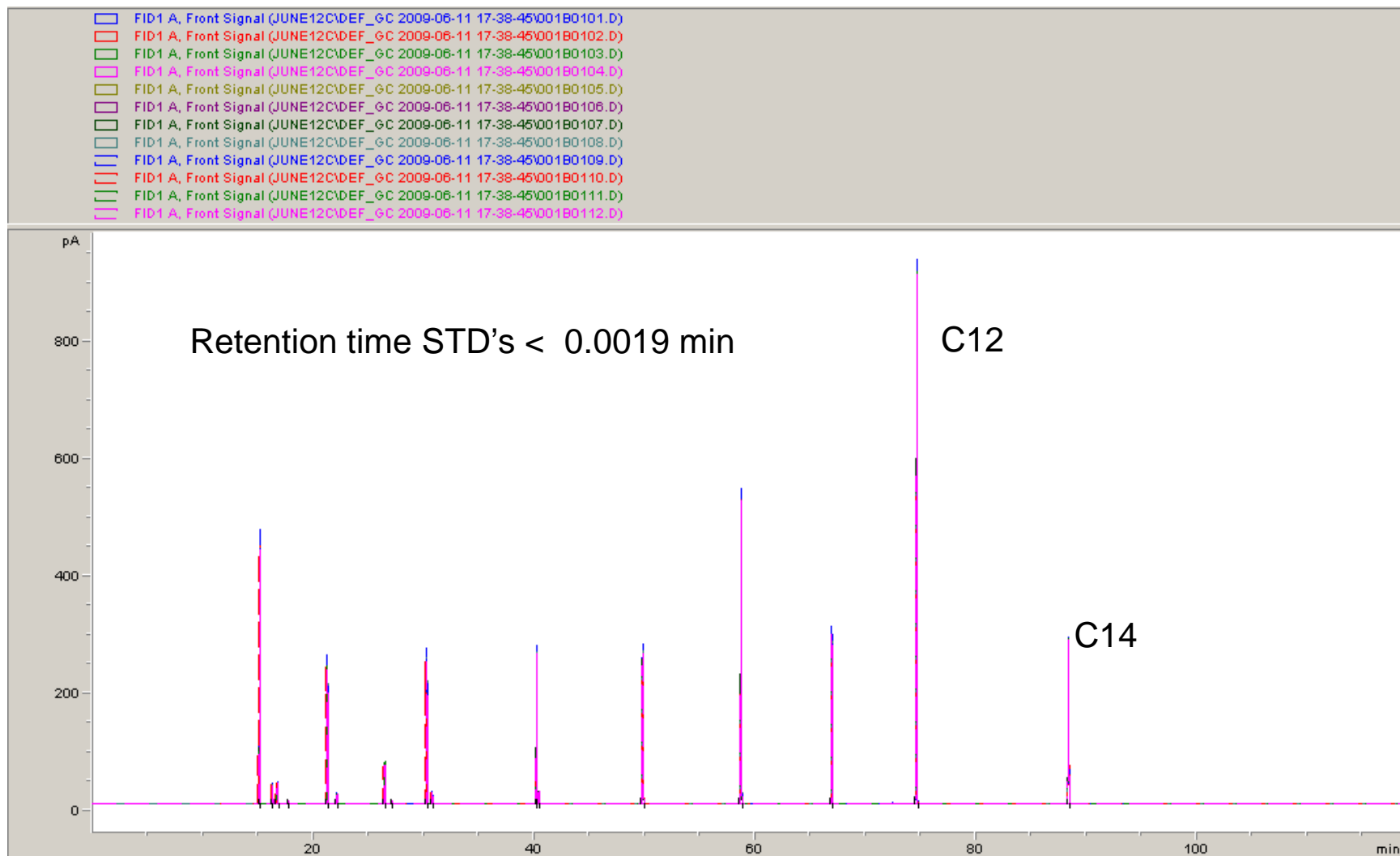


Analytical →

	Rate mL/min per min	Value mL/min	Hold Time min	Run Time min
▶ (Initial)		2	0	42.25
*				

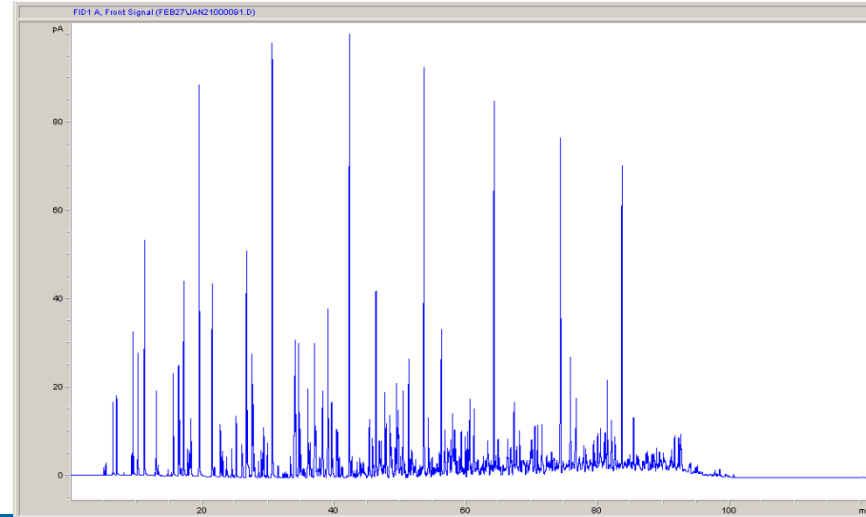
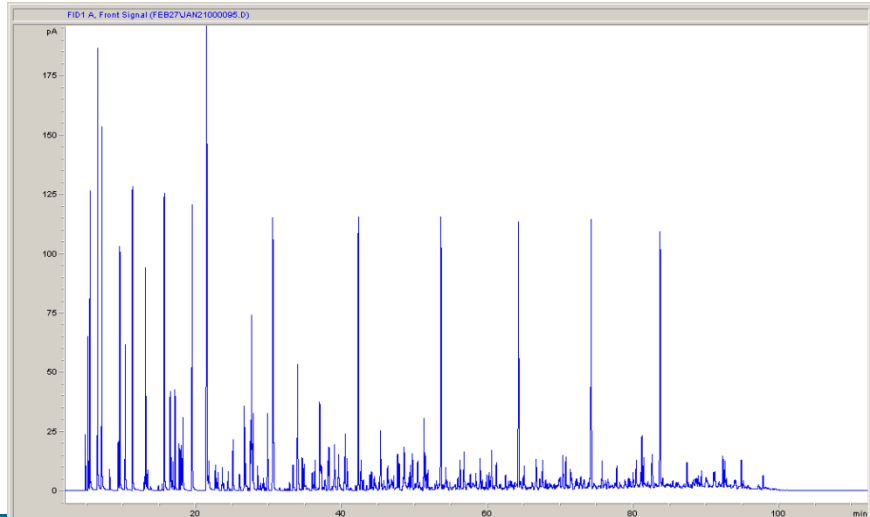
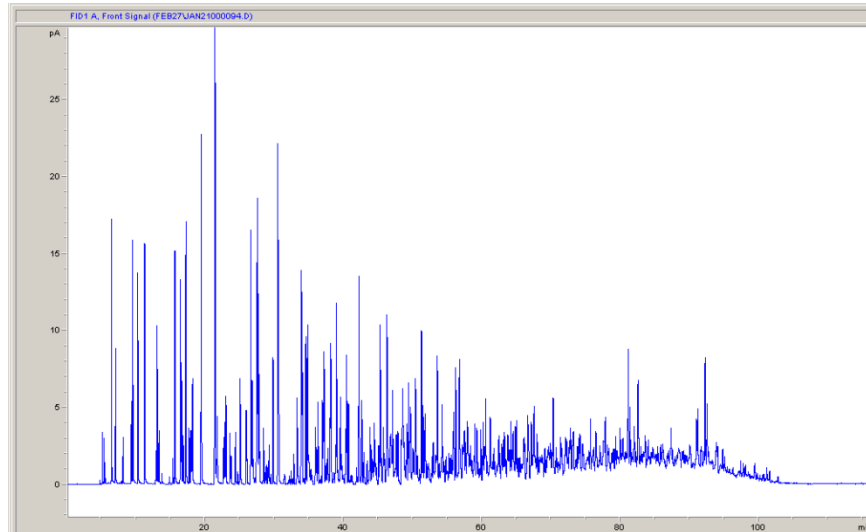
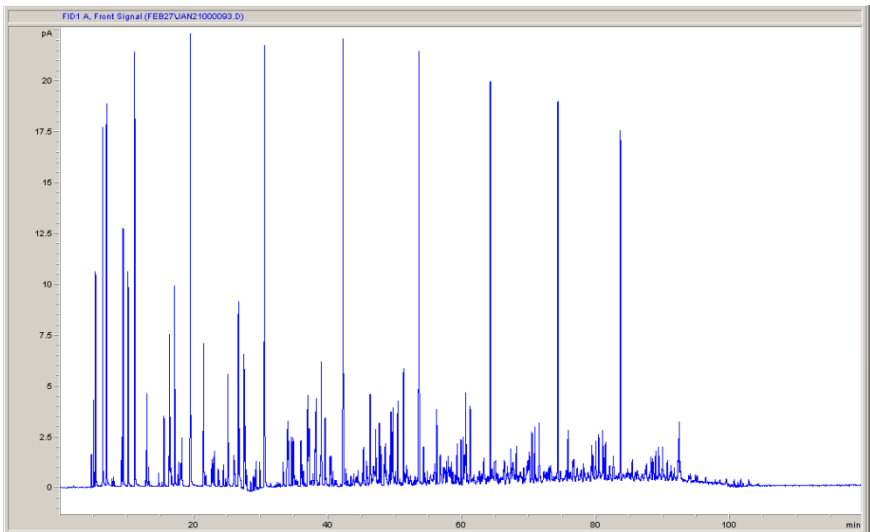
# Pre-column Backflush of C5-C18 Mix After C14

1 m x 0.32 mm retention gap, 100 m x 0.25 mm x 0.50 um DB-PETRO



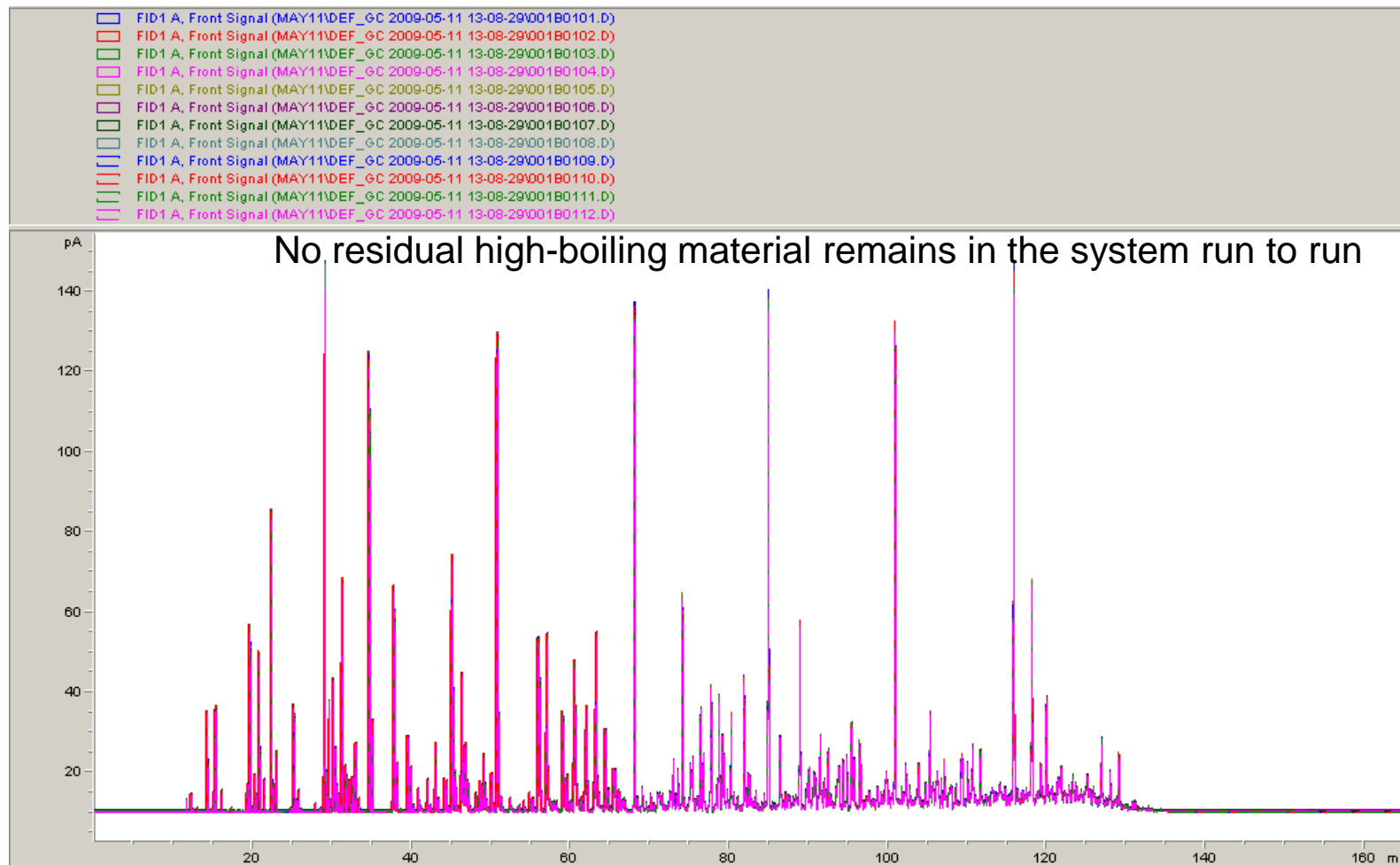
# Four Crudes From Different Regions: Backflush after C12

Pre-column: 1.5 m x 0.32 mm deactivated RG, Analytical column: 100 m x 0.25 mm x 0.50u DB-PETRO



# 12 runs of Crude Oil Overlay: Pre-column Backflush After C12

2 m x 0.32 mm RG

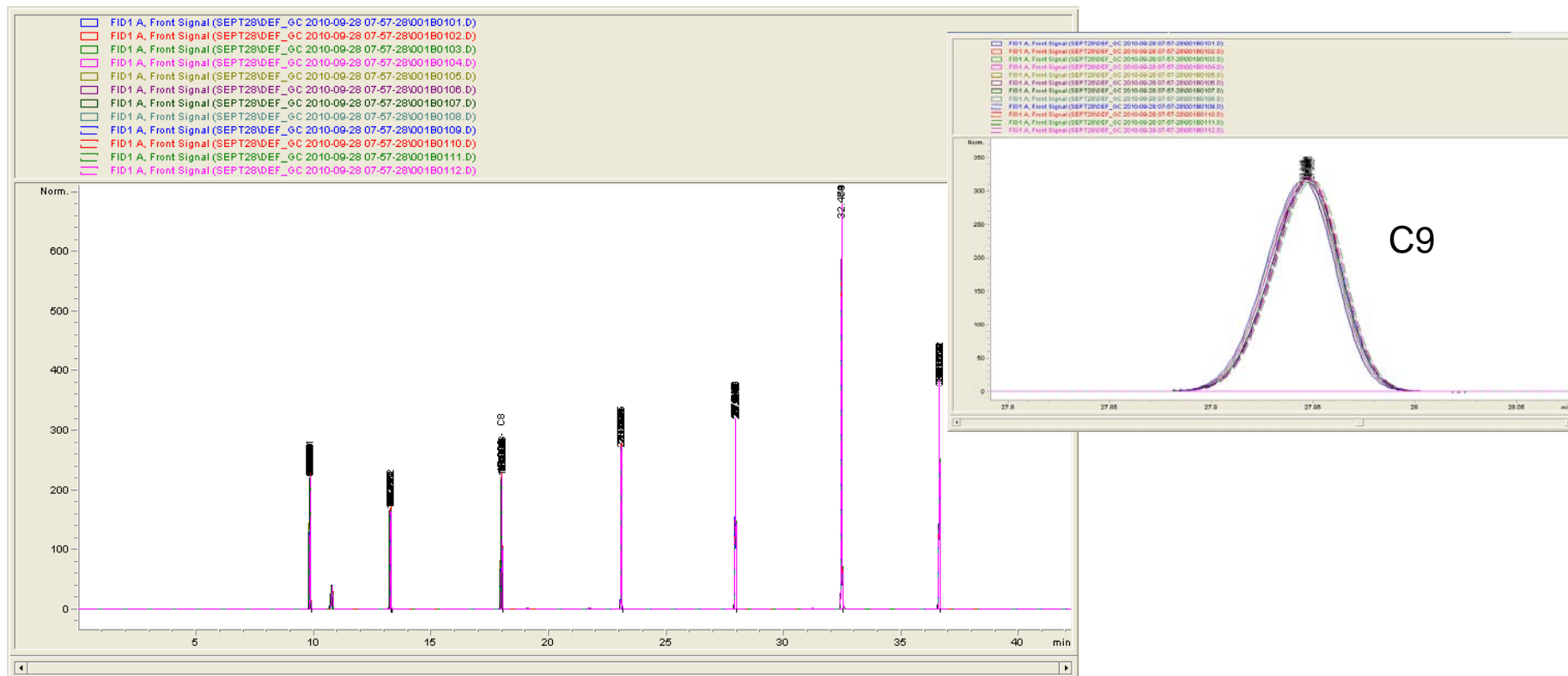


# Backflush with Coated Pre-columns

- Pre-columns with thin coating (i.e. 0.10  $\mu\text{m}$  films) can be used
- Provides finer control of compounds that are backflushed
- Backflush times will be much longer typically
- Backflush of heavy compounds remaining on the pre-column will take longer
- Better inertness

# Backflush of C5-C40 Mix After C11

Overlay of 12 runs, 2 m x 0.32 mm x 0.10  $\mu$ m pre-column





# Overlay of Souide Crude Oil

C10 Retention time STDEV = 0.0010, % STDEV = 0.315

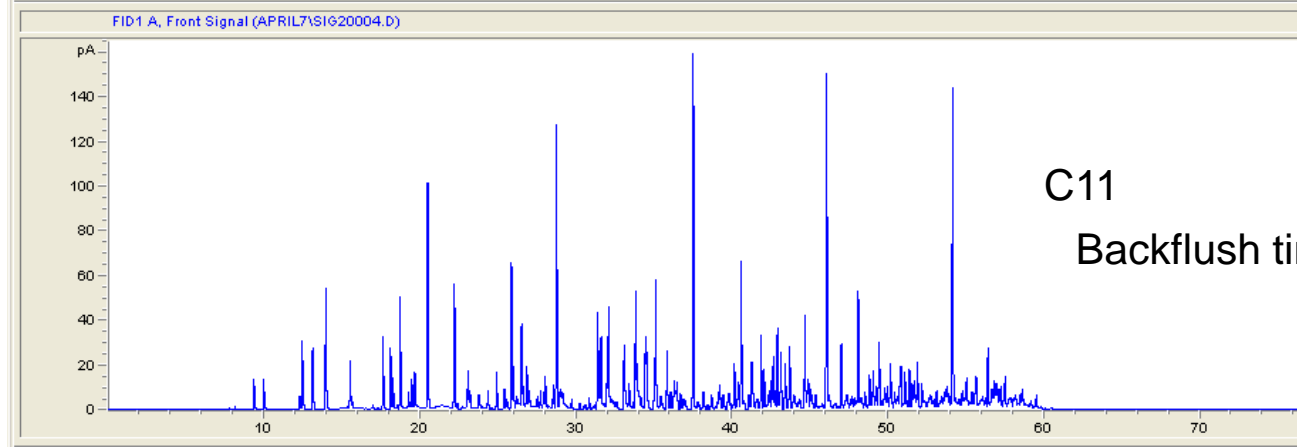
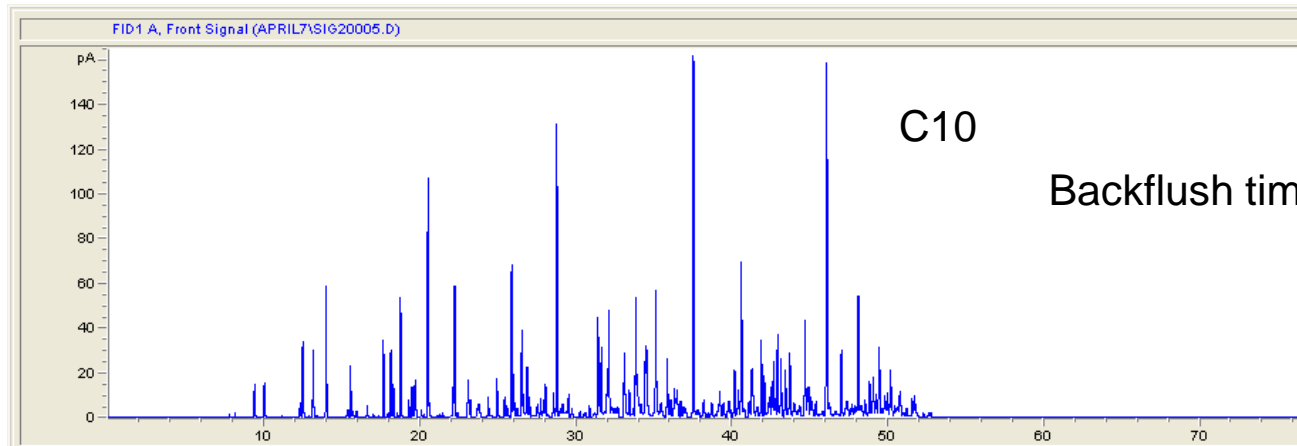


# Crude Oil Backflush with Coated 2 m x 0.32 mm x 0.10 um DB-1 Pre-column

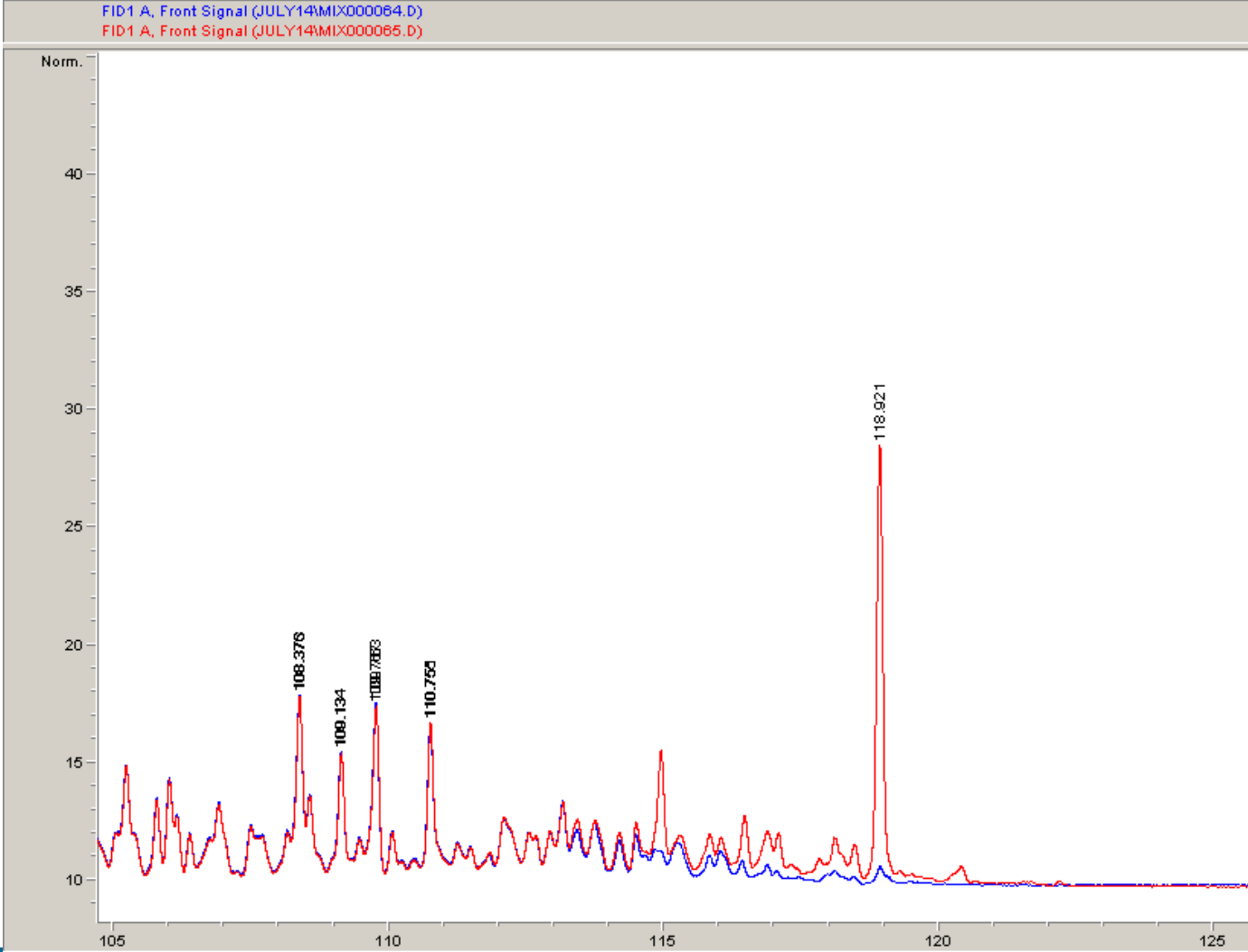
Conditions: Pre-column flow 1.6 ml/min, Analytical column flow 2.0 ml/min

MMI: 250 C(0.5 min) to 425 C @ 200 C/min

Oven Program: 35 C (4 min) to 200 C (1 min) @ 2.0 C/min

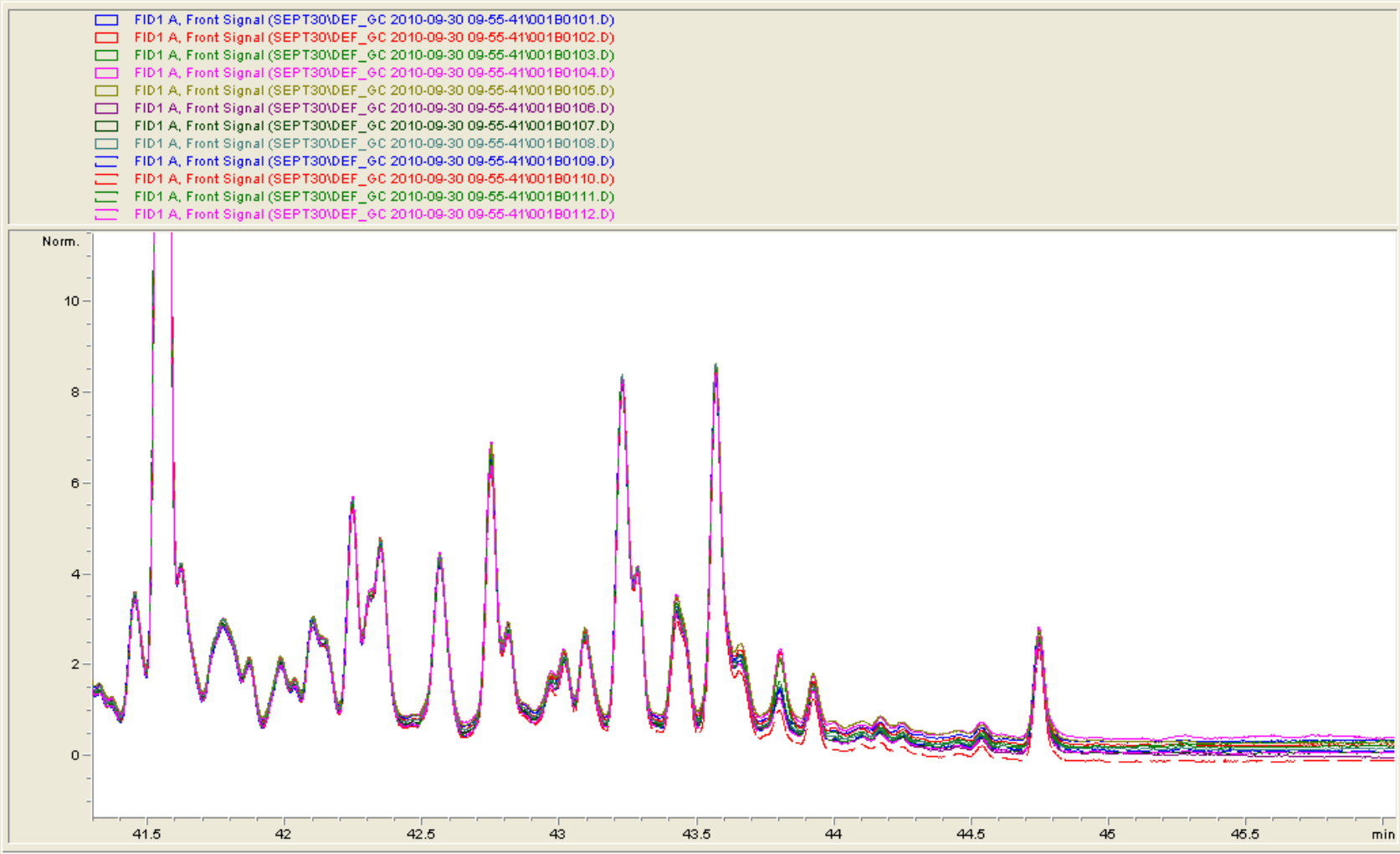


# 2 m x 0.32 mm x 0.10 um Pre-column: Blue Backflush at 14 min, Red – Backflush at 16 min



# Zoom on C11 Plus Hydrocarbons

Stable baseline shows that column is clean after each run



# Combining Pre-column Backflush with GCxGC

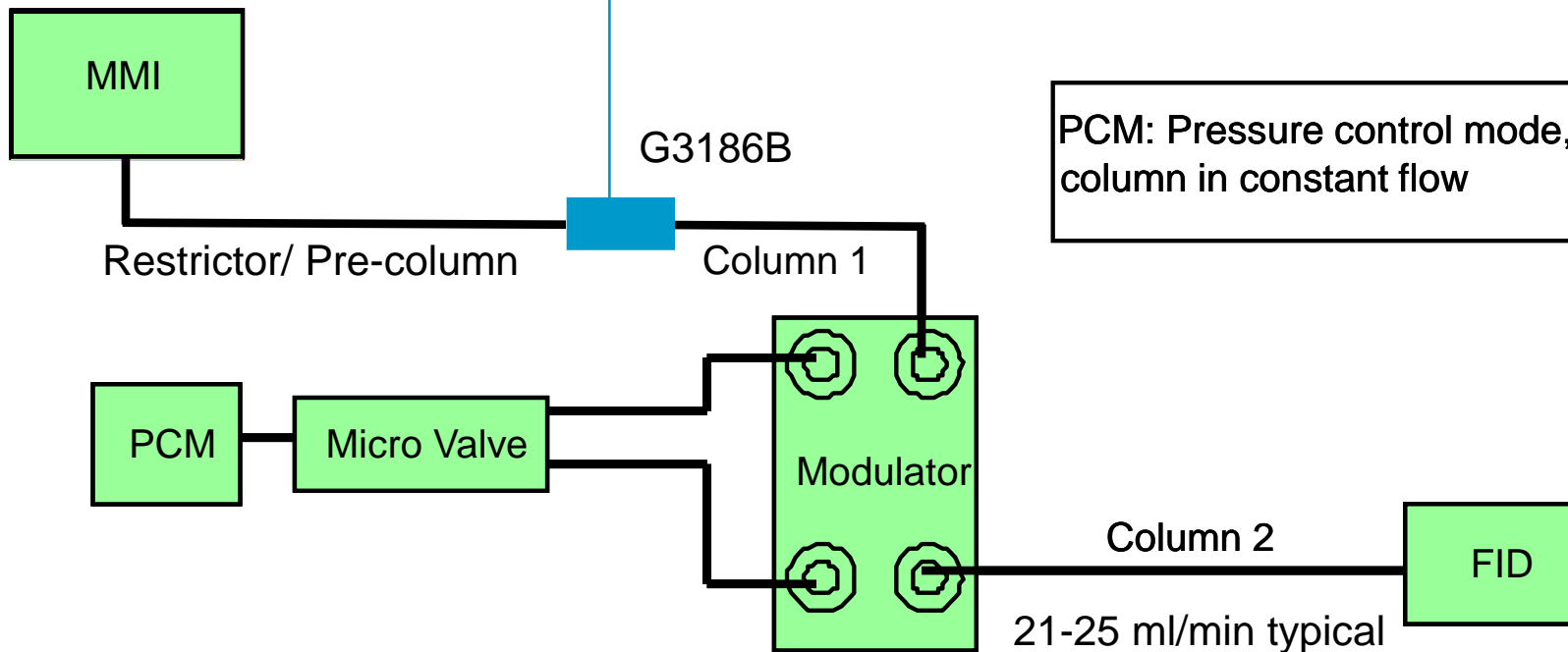
- Investigate possibility of analyzing heavy crudes with comprehensive gas chromatography
- Apply chemometrics to identify source or contamination
- Process 50 Hz raw data files
- Analyze four different crudes
- Column set
  - First Dimension: 30 m x 0.25 mm x 0.10  $\mu$ m DB-5HT
  - Second Dimension : 5 m x 0.25 mm x 0.15  $\mu$ m DB-17HT

# Flow Modulator Diagram: Backflush

MMI: ramped flow,  
negative ramp at backflush time

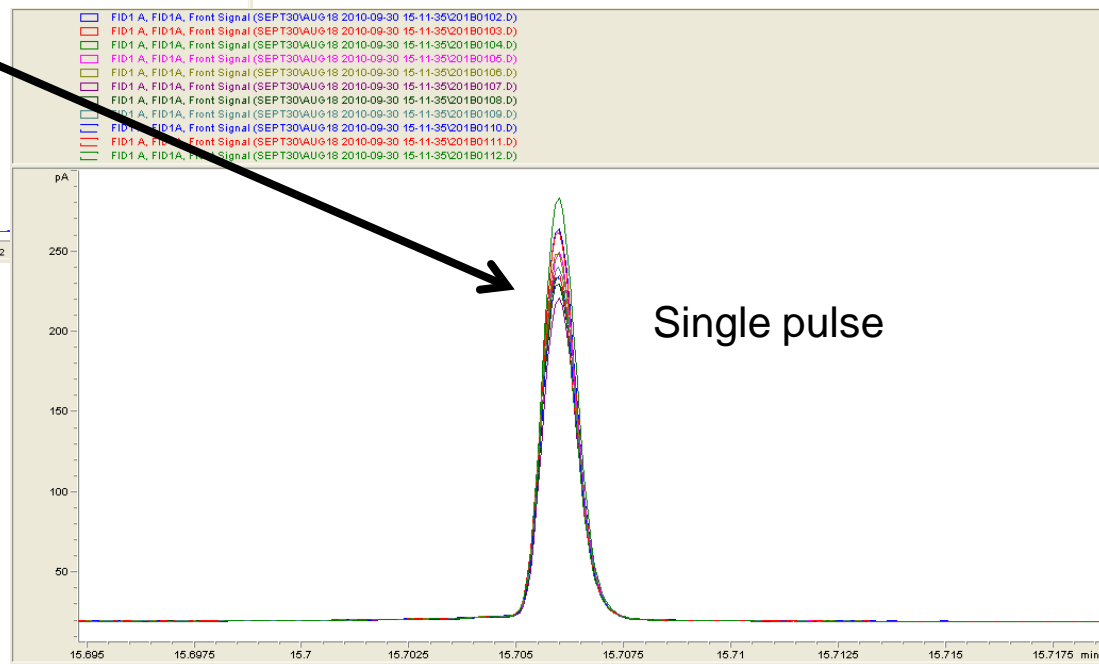
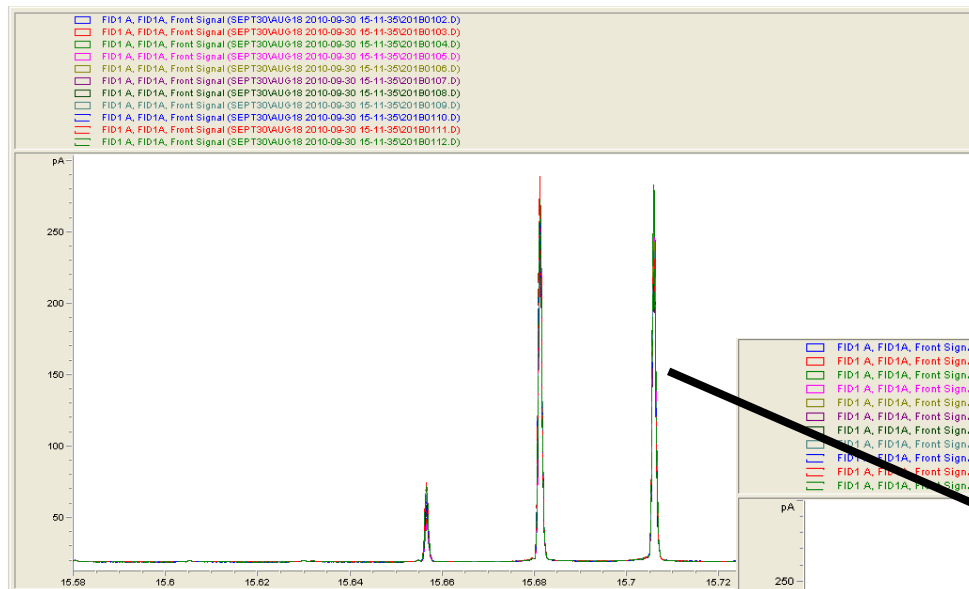
AUX

Aux: Controls Column 1 in  
constant flow



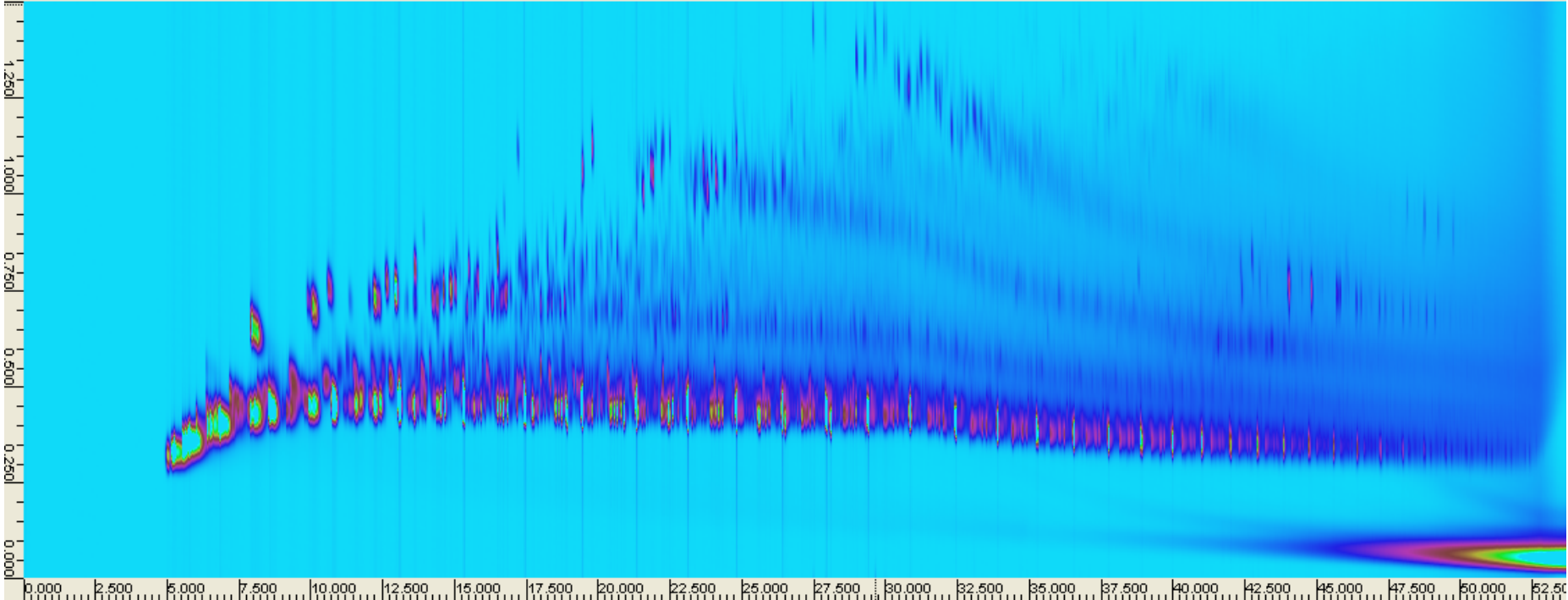
# RT Stability and Effect on Modulated Area: C13

## Analysis of modulated C13



# Crude Oil GCxGC: Pre- column Backflush

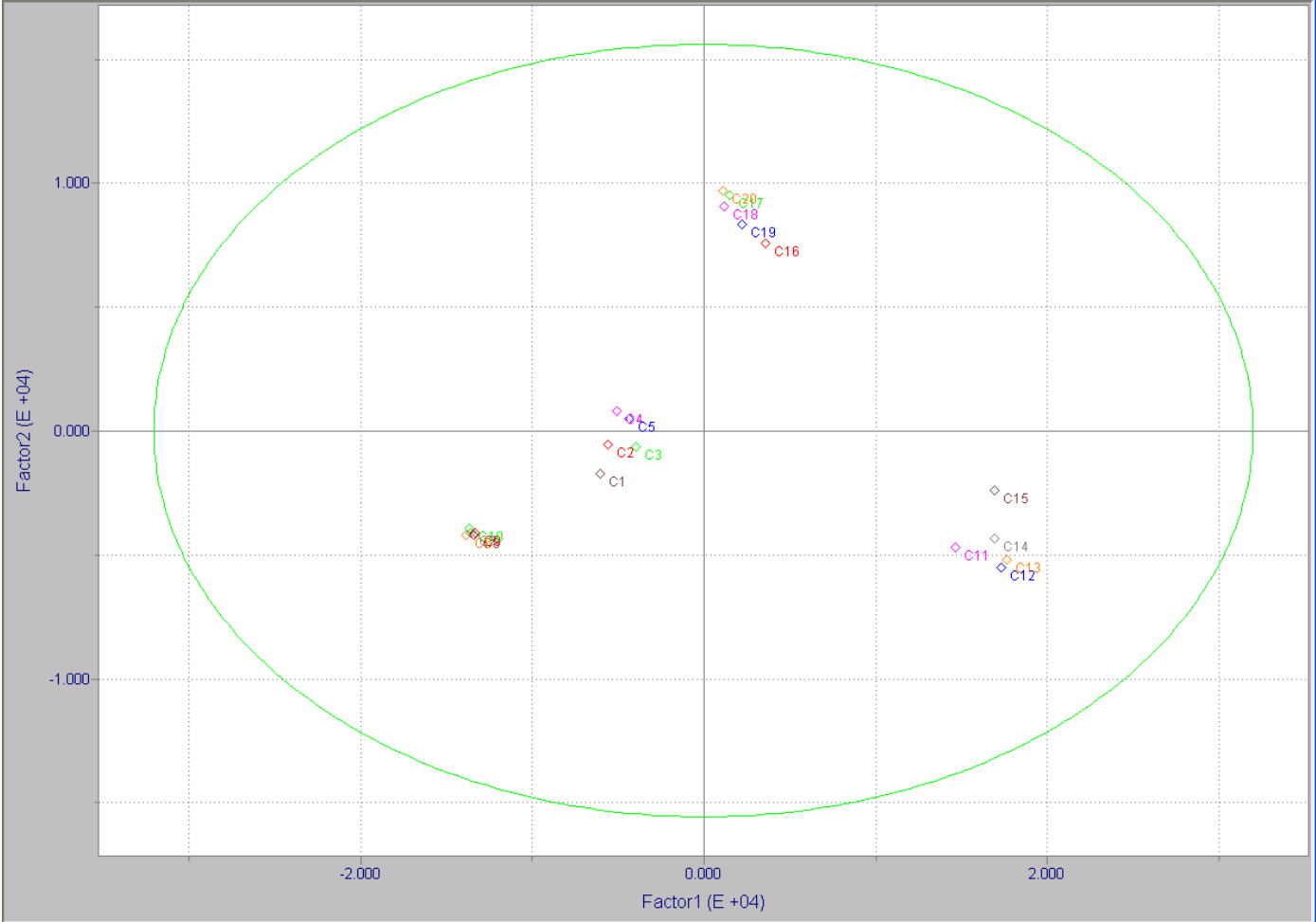
BF at 30 min, Oven programmed to 390 C



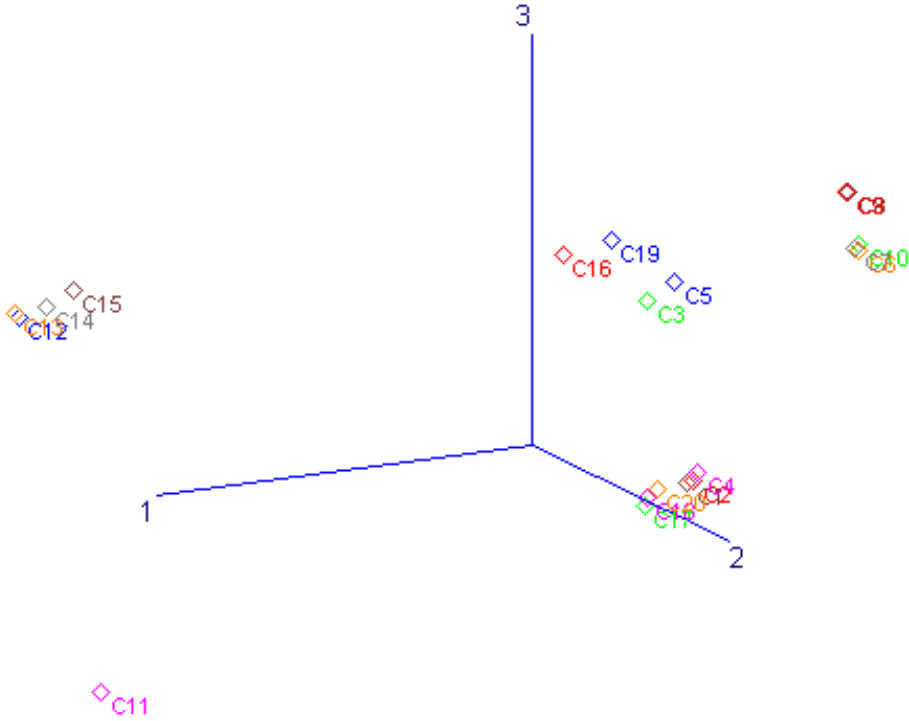


# Principal Component Analysis: 2D view

Four crudes



# Principal Component Analysis: 3D view



# Summary

- The CFT purged union offers a high performance alternative to traditional valved pre-column backflush systems
- Relies on precise EPC midpoint pressure control
- Backflush is concurrent while analytical column runs at constant flow
- Excellent thermal control since all parts are in the oven
- Heavy petroleum including crude oils can be routinely analyzed with complete protection of the high resolution capillary column
- The specific carbon number where backflush beginning can be easily controlled and fine tuned
- Pre-column backflush is compatible with GCxGC
- Coated pre-columns can be used for fine tuning
- Chemometrics can be applied to identify oils

# Acknowledgement

Pirouette® Chemometric Software provided by Infometrix, Inc.

# Capillary Flow Technology (CFT) Based Pre-fractionator: Pre-column Backflush System

- In oven CFT device is used in a relatively simple hardware configuration
  - CFT “purged union”
  - The multimode inlet is used to assist with backflushing heavy material from the inlet. Inlet is programmed to 425 C after backflush is started
  - Run time backflush is easily programmed from the ChemStation
  - Analytical column flow controlled from an Aux module
- Key Application: Front-end detailed analysis (DHA) of crude oils

# Flow Calculator Parameters

## 100M Column

**Column Pressure/Flow Calculator**

**Column Parameters**

Length (m)

i.d. (mm)

Temp (C)

**Carrier Gas Parameters**

Inlet Pressure (gauge)

Outlet Flow (mL/min)

Average Velocity (cm/s)

Outlet Pressure (Absolute)

## 2.0M RG

**Column Pressure/Flow Calculator**

**Column Parameters**

Length (m)

i.d. (mm)

Temp (C)

**Carrier Gas Parameters**

Inlet Pressure (gauge)

Outlet Flow (mL/min)

Average Velocity (cm/s)

Outlet Pressure (Absolute)

1 Atm  Vacuum  Other

## 1.5M RG

**Column Pressure/Flow Calculator**

**Column Parameters**

Length (m)

i.d. (mm)

Temp (C)

**Carrier Gas Parameters**

Inlet Pressure (gauge)

Outlet Flow (mL/min)

Average Velocity (cm/s)

Outlet Pressure (Absolute)

# 2 m x 0.32 mm x 0.10 um Pre-column: 12 Run Overlay of Crude Oil

