



# NEW 7890B Gas Chromatograph + 5977 Single Quadrupole

## January 2013

# Today's Challenges

## Economic - Routine -

- Productivity
- Cost of Ownership
- Investment Protection
- Ease-of-Use

## Economic - Method - Development

- Time = Money
- Skilled Resources  
... Harder to Find

## Sensitivity

- Trend – Lower LOD's
- "Imperfect" Samples



# Agilent's Response - 7890B GC & 5977A MSD - OpenLAB & MassHunter

## Economic - Routine -

- Productivity
- Investment Protection
- Cost of Ownership
- Ease-of-Use

**Integrated GC/MSD Communications**  
OpenLAB, MassHunter  
**Integrated GC/MSD Communications**  
Gas/Power Resource Management  
OpenLAB, MassHunter

## Economic - Method - Development

- Time = Money
- Skilled Resources  
... Harder to Find

**Integrated Intelligence**

## Sensitivity

- Trend – Lower LOD's
- "Imperfect" Samples

**5977A MSD**  
**GC FPD "Plus"**  
**Inert Flow Path**

# 7890B ↔ 5977A Direct, 2-Way (LAN) Communication ... Enhances and Protects Your GC/MSD Investment

## 5977A → 7890B

### If the MSD vents

- Increases GC gas flows
- Cools MSD faster
- **40% Faster Vent Times**

### If a Pump Event/Failure

- Stops GC Flows
  - Don't waste expensive He
  - Don't build-up H<sub>2</sub>



## 5977A ← 7890B

**GC monitors  
GC/MSD connection**

**If communication is  
lost...**

- GC shuts down various (GC) thermal zones

# 7890B – Better GC Selective Detector MDL Specs ... uECD, NPD

- No Changes in Designs
- Improved Manufacturing Processes/Controls

→ Less Variability

→ Lower MDL Specifications Guaranteed

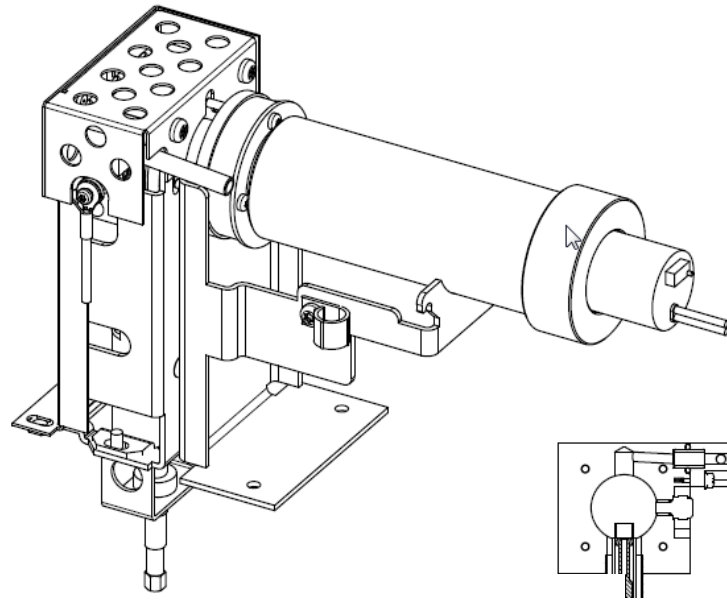
**uECD**    fg/sec                      5.5 → 4.5

**NPD**    pg N/sec                      0.1 → 0.08

            pg P/sec                      0.03 → 0.01

# New FPD "Plus"

- Advanced Photomultipliers
- Deactivation Treatments
- Design Improvements

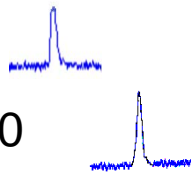


7890A  
FPD

Max Temp: 250C

Sulfur MDL (pg S/sec) : 3.6

Phosphorus MDL (fg P/sec): 60



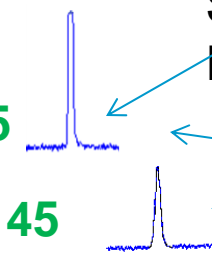
7890B  
FPD+  
400C

2.5

45

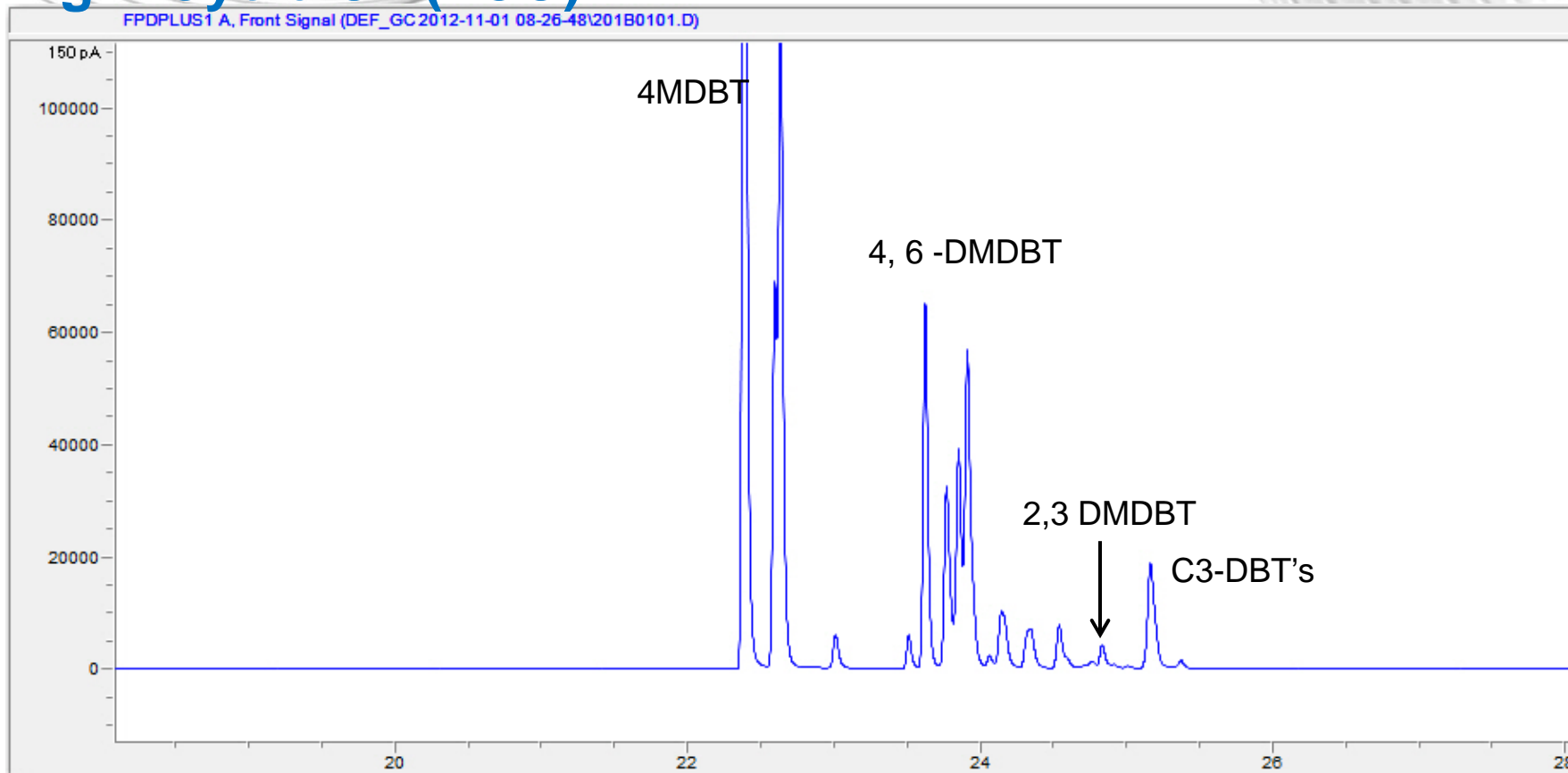
Higher Sulfur Response

Lower Noise



[Agilent SCD = 0.5 pg S/sec]

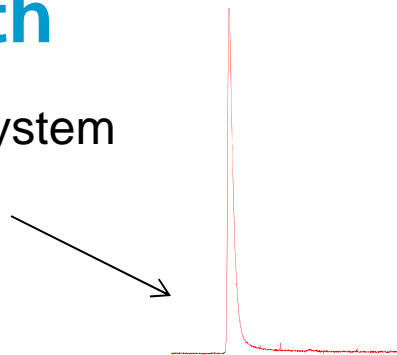
# Analysis of Substituted Dibenzothiophenes in Light Cycle Oil (LCO)



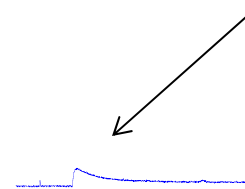
- The detector must be operated at temperatures above 300 C
- Agilent FPD Plus' higher temperature /improved sensitivity expands S-applications addressable by the easy-to-use FPD

# Improving Performance: Focus on sample flowpath

Inert system



Active system



Active or contaminated components in the flow path, before the detector, can reduce the overall sensitivity of the system



## Consider the following:

- Inlet liner
- Gold seal
- SSL inlet (weldment)
- Column
- Capillary Flow Device
- Detector

# Technologies that matter: Only from Agilent

Agilent has developed core competency in chemical vapor deposition surface deactivation resulting in proprietary processes to dramatically improve GC inertness.

## Deactivation processes for:

- Glass
- Steel
- Gold
- Glass wool



# Agilent Inert Flow Path ... From Injection to Detection

Ensures superior performance for trace level analysis

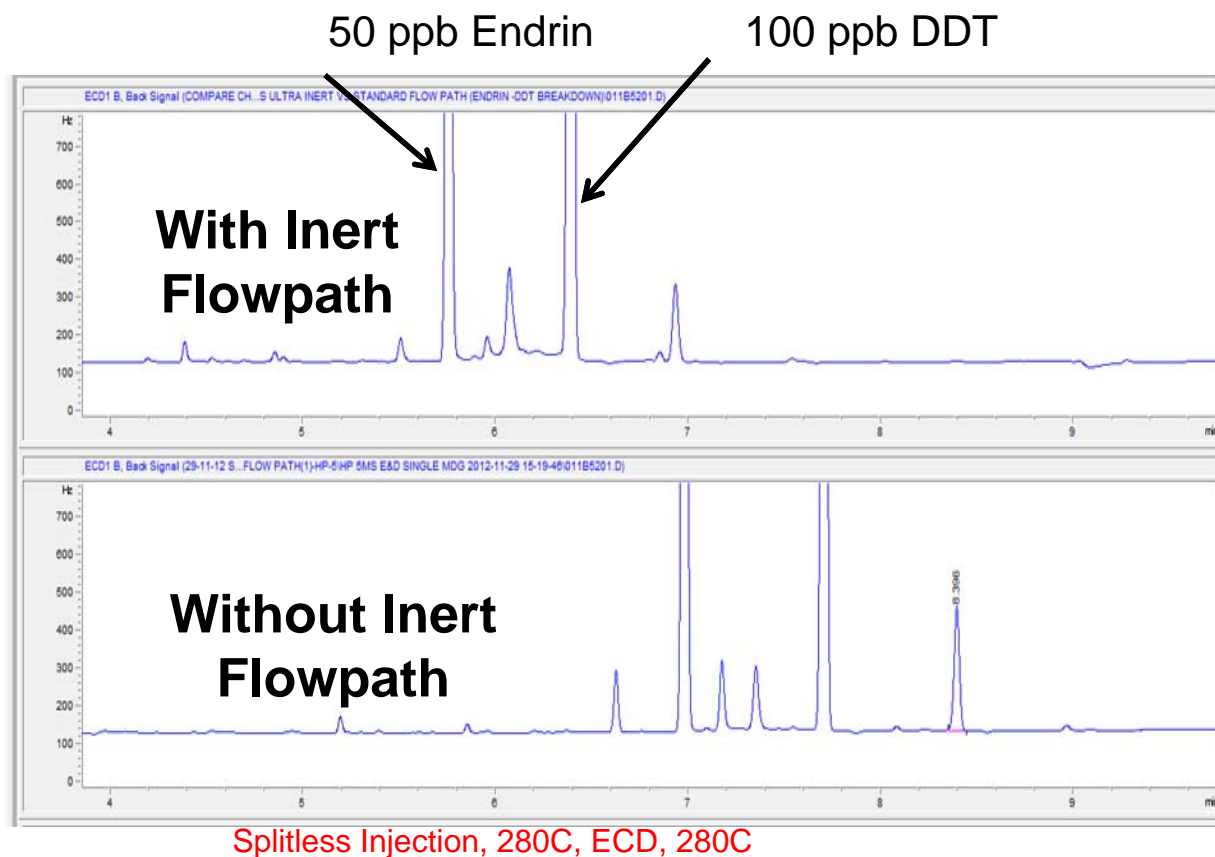


# Agilent's Inert Flowpath..From Injection to Detection

*Clearly less breakdown of a ppb level standard after 100 injections!*

## Consider the following:

- Inlet liner
- Gold seal
- Inlet weldment
- Column
- Capillary Flow Device
- Detector




# Integrated Intelligence ... A Smarter GC System

Integrated Intelligence	Benefit
Integrated GC calculators	Develop methods easier
Barcode Scanning with Auto-Input	Easier configuration
Integrated early-maintenance-feedback	Manage assets more efficiently
GC:MSD communication	Faster MSD vent Protect system better
3D Interactive Consumable and Parts Finder	Find parts faster & order easier
Sleep-Wake Modes ... Also coordinated with 5977A MSD	Lower cost of ownership

# 7890B Sleep/Wake Modes → Reduce Power/Gas Costs

Agilent 7890B Configuration: Instrument 1

Connection | Configuration | **Resource Conservation**

 Reduce gas and power consumption by setting gas saver and instrument schedule options

### Instrument Schedule

Select a schedule that best matches how you use this instrument:

Custom

#### Schedule

Day	Set Wake Method	Wake Time	Set Sleep Method	Sleep Time
Sunday	<input type="checkbox"/>		<input type="checkbox"/>	
Monday	<input type="checkbox"/>		<input type="checkbox"/>	
Tuesday	<input type="checkbox"/>		<input type="checkbox"/>	
Wednesday	<input type="checkbox"/>		<input type="checkbox"/>	
Thursday	<input type="checkbox"/>		<input type="checkbox"/>	
Friday	<input type="checkbox"/>		<input type="checkbox"/>	
Saturday	<input type="checkbox"/>		<input type="checkbox"/>	

Wake Method:       Sleep Method:

Wake to last active method before sleep

Perform a conditioning run on Wake

# Early Maintenance Feedback

## Improving Operational Efficiency

Maintenance

Early Maintenance Feedback Counters | Maintenance Log | Maintenance Actions

Front Inlet | Back Inlet | Front Detector | Back Detector | Third Detector | Front Injector | Column 1 | Column 2

NPD

Collector assembly injections	0	Injections	<input type="text"/>	+
Ceramics injections	2	Injections	<input type="text"/>	+
Bead baseline offset	0	$\mu A$	<input type="text"/>	+
Bead baseline voltage	0	Volts	<input type="text"/>	+
Bead integral of current	0	Unspecified	<input type="text"/>	+
Bead on time	:	0	<input type="text"/>	+
Bead injections	2	Injections	<input type="text"/>	+

[Add User Defined EMF](#) | [Hide selected EMFs on this page](#) | [Show all hidden EMFs](#) | [Print](#) | [Options](#)

# Method Translation Tools

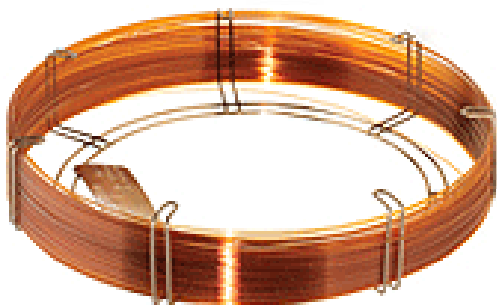
Completely Updated with AutoImport to Method

**GC Method Translation**

Criterion:  Translate Only  Best Efficiency  Fast Analysis  None **Speed gain: 2.50000**

	Original Method	Translated Method																														
<b>Column</b>																																
Length, m	30	<input type="checkbox"/> 20																														
Internal Diameter, $\mu\text{m}$	250	<input type="checkbox"/> 180																														
<b>Film</b>		<input type="radio"/> Unlock																														
Thickness, $\mu\text{m}$	0.25	<input type="radio"/> 0.180																														
Phase Ratio	250.0	<input checked="" type="radio"/> 250.0																														
<b>Carrier Gas</b>	Helium	<input type="checkbox"/> Hydrogen																														
Enter one Setpoint		<input checked="" type="radio"/> Unlock																														
Head Pressure, psi	7.843	<input type="radio"/> 7.078																														
Flow Rate, mL/min	1	<input checked="" type="radio"/> 0.8347																														
Outlet Velocity, cm/sec	Very large	Very large																														
Average Velocity, cm/sec	36.14	<input type="radio"/> 60.23																														
Hold-up Time, min	1.38365	<input type="radio"/> 0.55346																														
Outlet Pressure (absolute), psi	0	<input checked="" type="checkbox"/> 0																														
Ambient Pressure (absolute), psi	14.696	<input type="checkbox"/> 14.696																														
<b>Oven Temperature</b> 1-ramp Program																																
	<table border="1"> <thead> <tr> <th>Ramp Rate</th> <th>Final Temp.</th> <th>Final Time</th> </tr> <tr> <th><math>^{\circ}\text{C}/\text{min}</math></th> <th><math>^{\circ}\text{C}</math></th> <th>min</th> </tr> </thead> <tbody> <tr> <td></td> <td>50</td> <td>2</td> </tr> <tr> <td>Ramp 1</td> <td>10</td> <td>300</td> </tr> <tr> <td></td> <td>300</td> <td>5</td> </tr> </tbody> </table>	Ramp Rate	Final Temp.	Final Time	$^{\circ}\text{C}/\text{min}$	$^{\circ}\text{C}$	min		50	2	Ramp 1	10	300		300	5	<table border="1"> <thead> <tr> <th>Ramp Rate</th> <th>Final Temp.</th> <th>Final Time</th> </tr> <tr> <th><math>^{\circ}\text{C}/\text{min}</math></th> <th><math>^{\circ}\text{C}</math></th> <th>min</th> </tr> </thead> <tbody> <tr> <td></td> <td>50</td> <td>0.800</td> </tr> <tr> <td></td> <td>25.000</td> <td>300</td> </tr> <tr> <td></td> <td>300</td> <td>2.000</td> </tr> </tbody> </table>	Ramp Rate	Final Temp.	Final Time	$^{\circ}\text{C}/\text{min}$	$^{\circ}\text{C}$	min		50	0.800		25.000	300		300	2.000
Ramp Rate	Final Temp.	Final Time																														
$^{\circ}\text{C}/\text{min}$	$^{\circ}\text{C}$	min																														
	50	2																														
Ramp 1	10	300																														
	300	5																														
Ramp Rate	Final Temp.	Final Time																														
$^{\circ}\text{C}/\text{min}$	$^{\circ}\text{C}$	min																														
	50	0.800																														
	25.000	300																														
	300	2.000																														
<b>Sample Information</b> None																																

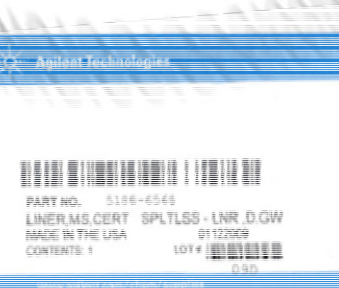
# Barcode Scanning Input



Columns



Liners



Barcode in the columns, liners, and syringes right into the Agilent data system method



Favorite	Part Number	Description	Length, m	Diameter, µm	Flow Path, mm	Inner Dia, µm	Phase Ratio	Max Temp, °C	Min Temp, °C	New Temp, °C	Port Temp, °C	Port Factor	Comments	Time Stamp
<input type="checkbox"/>	19091A-012	ULTRA 1	25	320	6.17	463.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-013	ULTRA 1	25	320	6.17	463.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-013.7M	ULTRA 1	25	320	6.17	463.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-013	ULTRA 1	30	320	6.17	463.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-016	ULTRA 1	30	320	6.17	463.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-051	ULTRA 1	12	200	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-051R	ULTRA 1	12	200	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-051.7M	ULTRA 1	12	200	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-052	ULTRA 1	25	320	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-052R	ULTRA 1	25	320	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-052.7M	ULTRA 1	25	320	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-051	ULTRA 1	30	320	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-052	ULTRA 1	30	320	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-052.7M	ULTRA 1	30	320	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-058	ULTRA 1	17	200	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-058.7M	ULTRA 1	17	200	6.33	355.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-112	ULTRA 1	25	320	6.52	353.1	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-112R	ULTRA 1	25	320	6.52	353.1	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-112.7M	ULTRA 1	25	320	6.52	353.1	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-111	ULTRA 1	30	320	6.52	353.1	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-112R	ULTRA 1	30	320	6.52	353.1	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-052	ULTRA 2	25	300	6.11	453.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-052R	ULTRA 2	25	300	6.11	453.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-052.7M	ULTRA 2	25	300	6.11	453.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-051	ULTRA 2	30	300	6.11	453.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-051R	ULTRA 2	30	300	6.11	453.8	40	325	300	Perish				16/07/12 13:00:37
<input type="checkbox"/>	19091A-051.7M	ULTRA 2	30	300	6.11	453.8	40	325	300	Perish				16/07/12 13:00:37

\*Also syringes


Parts Finder

Agilent Technologies  
**PARTS FINDER**


FIND PARTS

ALS

7693A ALS




7650A Automatic Liquid Samp




GC


Agilent 7890B



Agilent 7890A




Agilent 6890N



MS

5975 MSD



PROMOTIONS

Select the instrument

FIND PARTS

PARTS LIST **3**

FAVORITES

UPDATES

HELP

PROMOTIONS

Ultra Inert Split Liners		
Add To Parts List   Add To Favorites		
<input type="checkbox"/>	Part Number	Part Description
<input type="checkbox"/>	<a href="#">5190-2294</a>	Liner, ultra inert, split, straight, glass wool
<input checked="" type="checkbox"/>	<a href="#">5190-3164</a>	Liner, Ultra Inert deactivation, split, straight, glass wool, 5/pk
<input checked="" type="checkbox"/>	<a href="#">5190-3168</a>	Liner, Ultra Inert deactivation, split, straight, glass wool, 25/pk
<input type="checkbox"/>	<a href="#">5190-2295</a>	Liner, ultra inert, universal, low pressure drop, GW
<input checked="" type="checkbox"/>	<a href="#">5190-3165</a>	Liner, Ultra Inert, universal, low pressure drop, glass wool, 5/pk
<input type="checkbox"/>	<a href="#">5190-3169</a>	Liner, Ultra Inert, universal, low pressure drop, glass wool, 25/pk
<input type="checkbox"/>	<a href="#">5190-4048</a>	Liner, GC, Ultra Inert, straight, 0.75 mm id, recommended for SPME injections

Click Add to Parts List

# 7890A Enhancements Kit

## 7890A Logic Board Enhancement Kit

- Makes a 7890A GC ready for most (but not all\*) 7890B GC features

- **“Integrated Intelligence” features:**

- Early Maintenance Feedback (EMF), Sleep/Wake modes
- Method Translation Calculator
- Enable 2-way GC/5977A MSD communication

- **Agilent Serial Port:**

- 2<sup>nd</sup> communication port for Remote Advisor
- Optional Bar Code Reader connects directly to GC  
[also a version for USB/PC]

- **FPD Plus**

- \* **7890B Only**

- 2<sup>nd</sup> MMI on GC, 3<sup>rd</sup> Detector as uECD



# Agilent's Response - 7890B GC & 5977A MSD - OpenLAB & MassHunter

## Economic - Routine -

- Productivity
- Investment Protection
- Cost of Ownership
- Ease-of-Use

**GC/MSD Communications**  
**GC/MSD Communications**  
**Gas/Power Resource Management**  
**OpenLAB, MassHunter**

## Economic - Method - Development

- Time = Money
- Skilled Resources  
... Harder to Find

**Integrated Intelligence**

## Sensitivity

- Trend – Lower LoD's
- "Imperfect" Samples

**5977A MSD**  
**GC FPD Plus**

**Inert Flow Path**

# **Agilent 5977 MSD** New GC/MSD Introduction **Overview**

January 2013



# New Agilent 5977 Series GC/MSD

The 5977 is the new Agilent GC/MSD platform and replaces the 5975C and 5975E systems

- Improves performance with high sensitivity **Extractor Ion Source**
- **MassHunter GC/MS Acquisition**
- **MassHunter** and **MSD ChemStation** data analysis
- Integration of **new Agilent Vacuum Pumps** brings intelligent operation
- **Direct Communication** between 7890B and 5977A for more productivity



# New Agilent 5977 Series GC/MSD

What is the same as 5975?

- ✓ Quality
- ✓ Reliability
- ✓ Proven core technology since 5973/5975
  - Inert Ion Source
  - Heatable Gold Quartz Quadrupole
  - Triple Axis Detector
  - High Energy Diode
  - Triple Channel EM Detector
  - Productivity Software  
(MSD ChemStation Data Analysis)



# New Agilent 5977 Series GC/MSD

What is new with 5977?

- ✓ Better Performance
  - High Sensitive Extractor Ion Source
  - Improved Agilent Vacuum Pump
- ✓ More Productivity
  - Direct Communication 7890B ↔ 5977A
    - Protect system
    - Vent the system faster
    - Power save (Sleep/Wake)
  - MassHunter Software
    - Data Acquisition (Method/Sequence)
    - Data Analysis



## New Extractor Ion Source for 5977

Why the Extractor EI Source?



Increase in ion count

- ✓ True increase in sensitivity

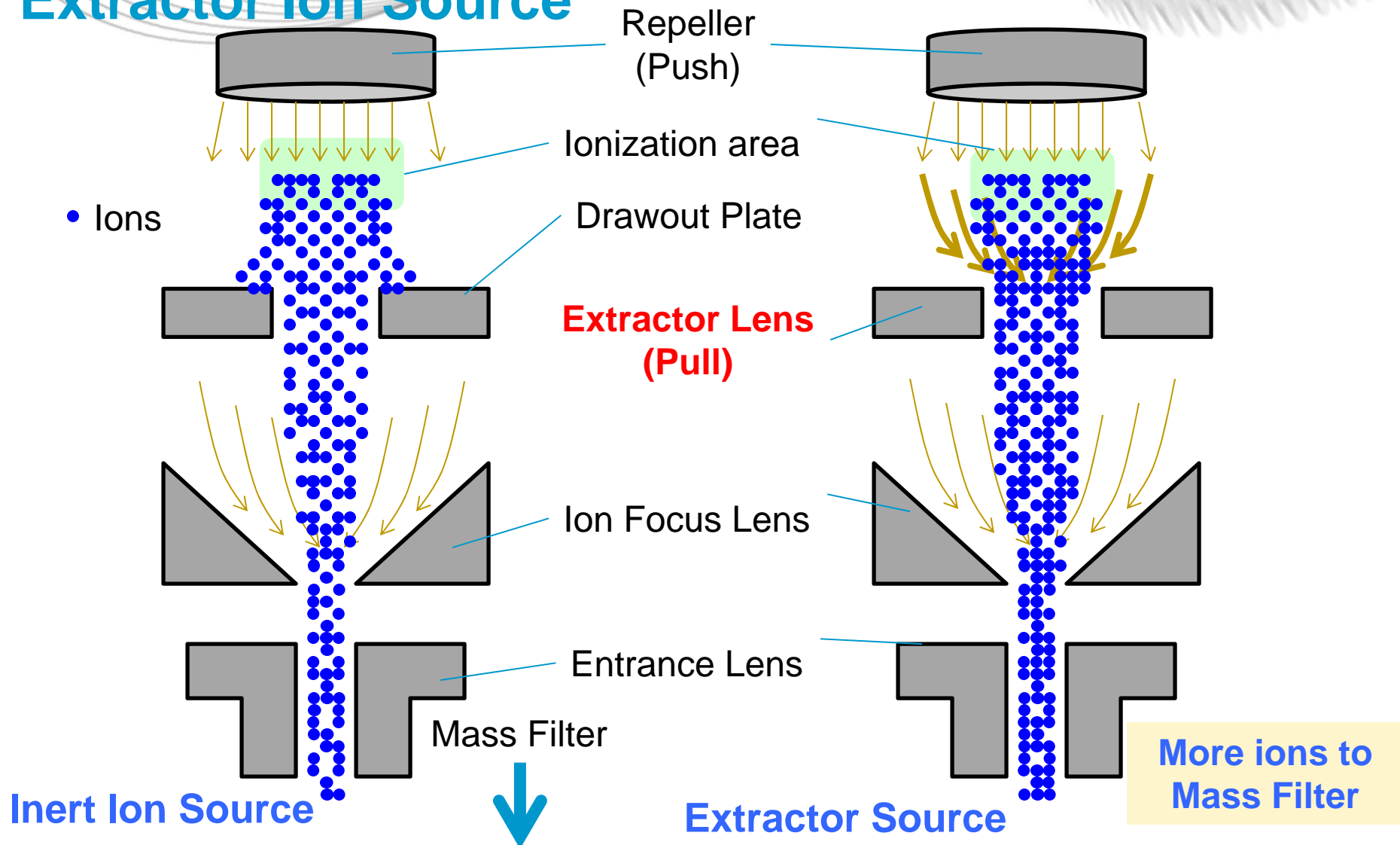


Better trace level precision

- ✓ Lower IDL and MDL

**Proven sensitivity increase demonstrated with  
over 1500 7000 Series GC/TQ!**

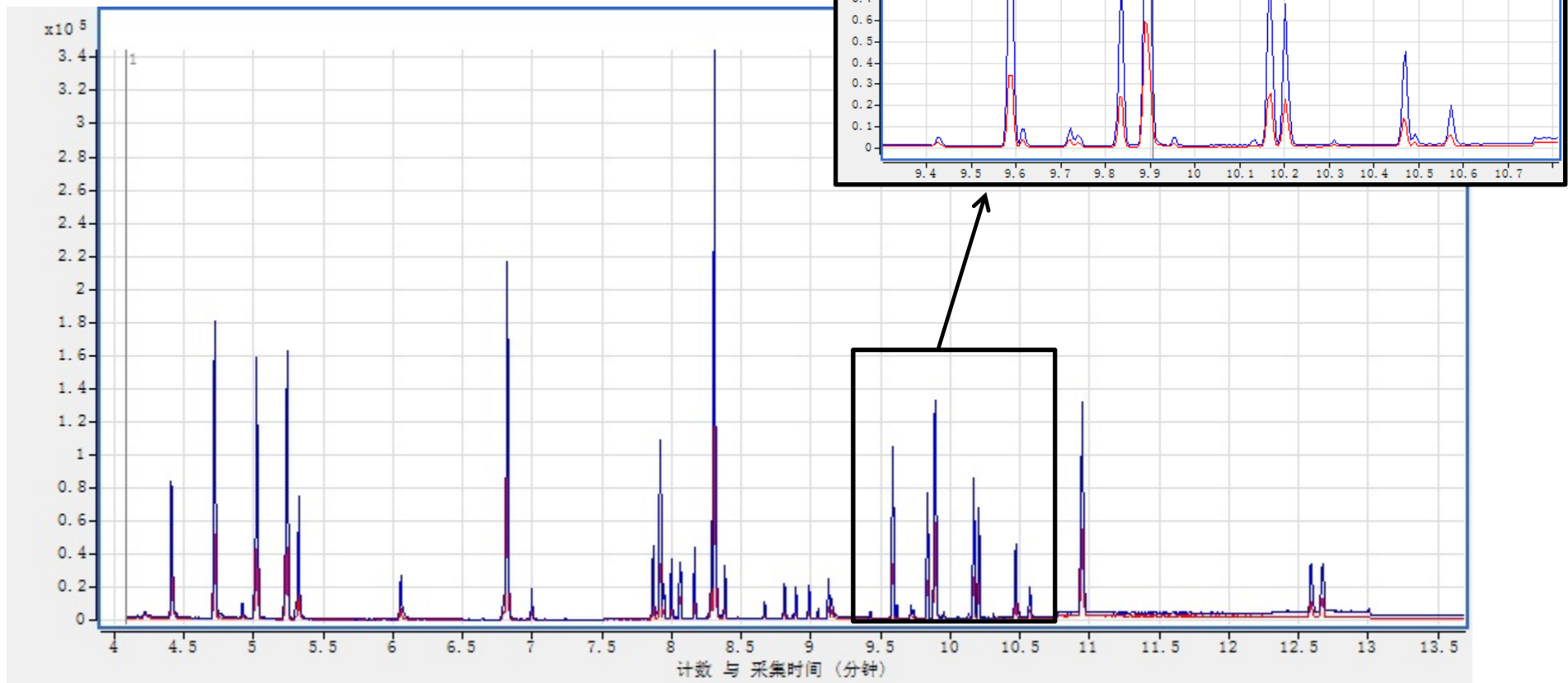
# Comparison between Inert Ion Source and Extractor Ion Source



# How Extractor Ion source with “Etune” be seen in the Chromatogram?

Blue – Extractor 40ppb solution (SVOC)

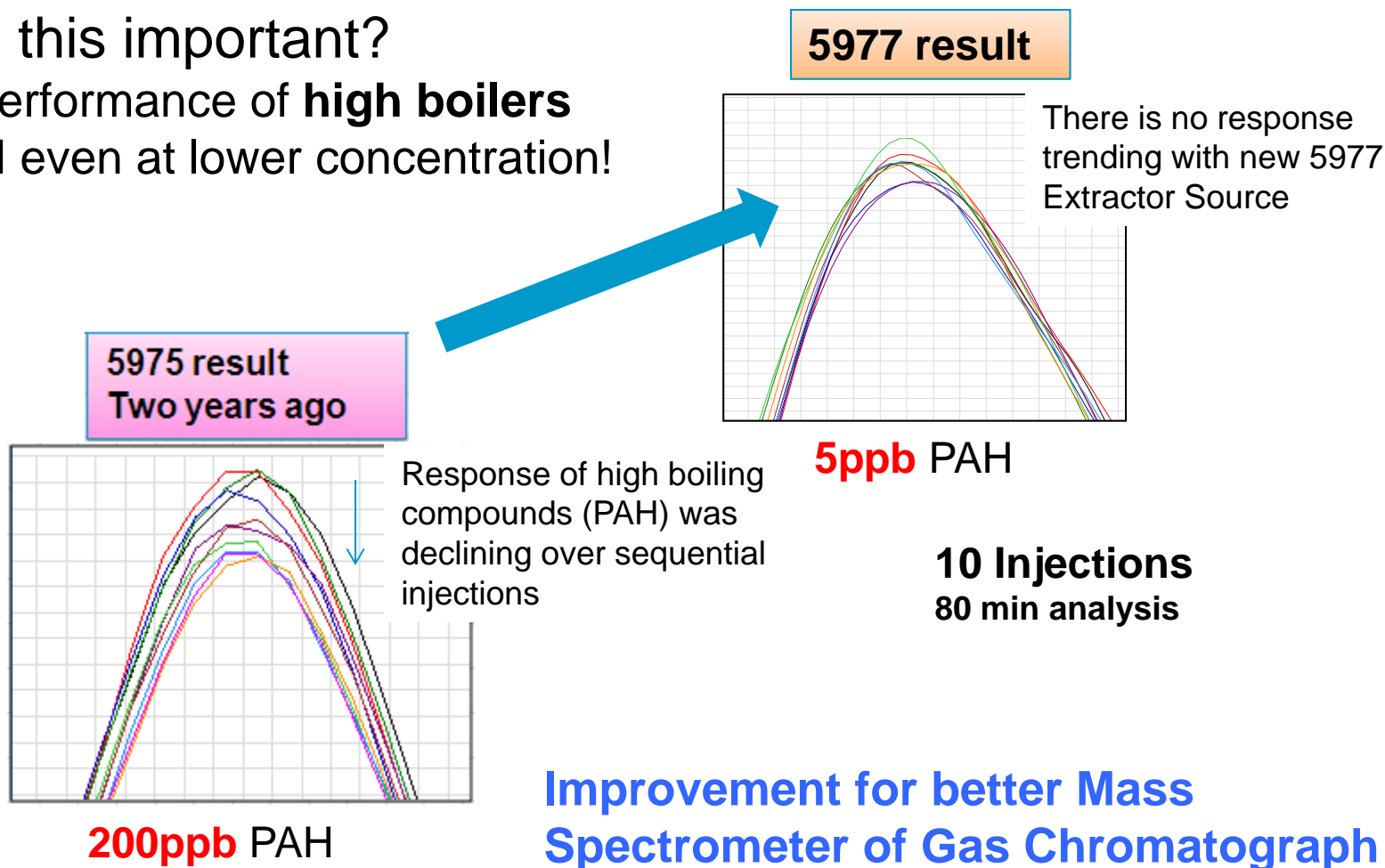
Red – Standard 40ppb solution (SVOC)



# More improvements on 5977's Extractor Ion Source: Improved Thermal Profile

Why is this important?

Better performance of **high boilers**  
**i.e. PAH** even at lower concentration!



## Sensitivity Specification of 5977

Signal-to-noise (S/N) specification has been used as the sensitivity specification of GC/MS for long years.

Agilent 5977 has highest S/N (1500:1) in Extractor Ion Source among Single Quadrupole GC/MS system.

Agilent 5977A introduces *new* sensitivity specification of Instrument Detection Limit (IDL), which shows the detectability of trace compounds of the GC/MS system.



## Why IDL specification is useful for 5977?

- S/N **can be** higher when the Noise is too low, such as SIM acquisition.
- S/N can only show the sensitivity performance of Single-shot injection. It doesn't indicate reproducibility of system performance.
- IDL uses statistic calculation to estimate the detection limit of the instruments (the same calculation of Method Detection Limit; MDL) both sensitivity and reproducibility.

You will get more accurate sensitivity performance of GC/MS system from IDL than single-shot S/N.



## 5977A GC/MSD IDL Specificaiton

IDL is represented in the concentration of the compound with the confidence level.

For example, Agilent 7890B/5977A with Auto Liquid Sampler (ALS) has following specification at 99% confidence level with 8 repeated injections.

Ion Source type	High Vacuum pump type	IDL of OFN by 8 replicates of ALS
Extractor Source	Turbo Pump	10fg
Inert Source	Turbo Pump	24fg
	Diffusion Pump	30fg

## Meaning of IDL

10 fg IDL specification of 5977A Extractor Ion Source with Turbo Pump means that the GC/MSD with ALS will detect 10fg OFN or lower at 99% confidence from 8 repeated injections.

For accurate estimation of IDL, accurate sample injection to GC/MSD system is critical.

Agilent only have IDL specification for ALS system. (7693 Auto injector or CTC sampler)

The GC/MSD system without ALS has S/N specification.

# Better HW Features: Agilent Turbo Pump (V304)



Innovation for:

- **Higher pumping performance**
  - especially for He and H<sub>2</sub>
- **High reliability**
  - lower cost of operation

**Agilent Quality Standard**

**TwisTorr 304 FS**

Press release : <http://www.agilent.com/about/newsroom/presrel/2012/19dec-ca12077.html>

# Better HW Features: Agilent Rough Pumps



**New** DS42i

High performance  
**Energy saving** design  
“Inverter” mode not active with diff pump

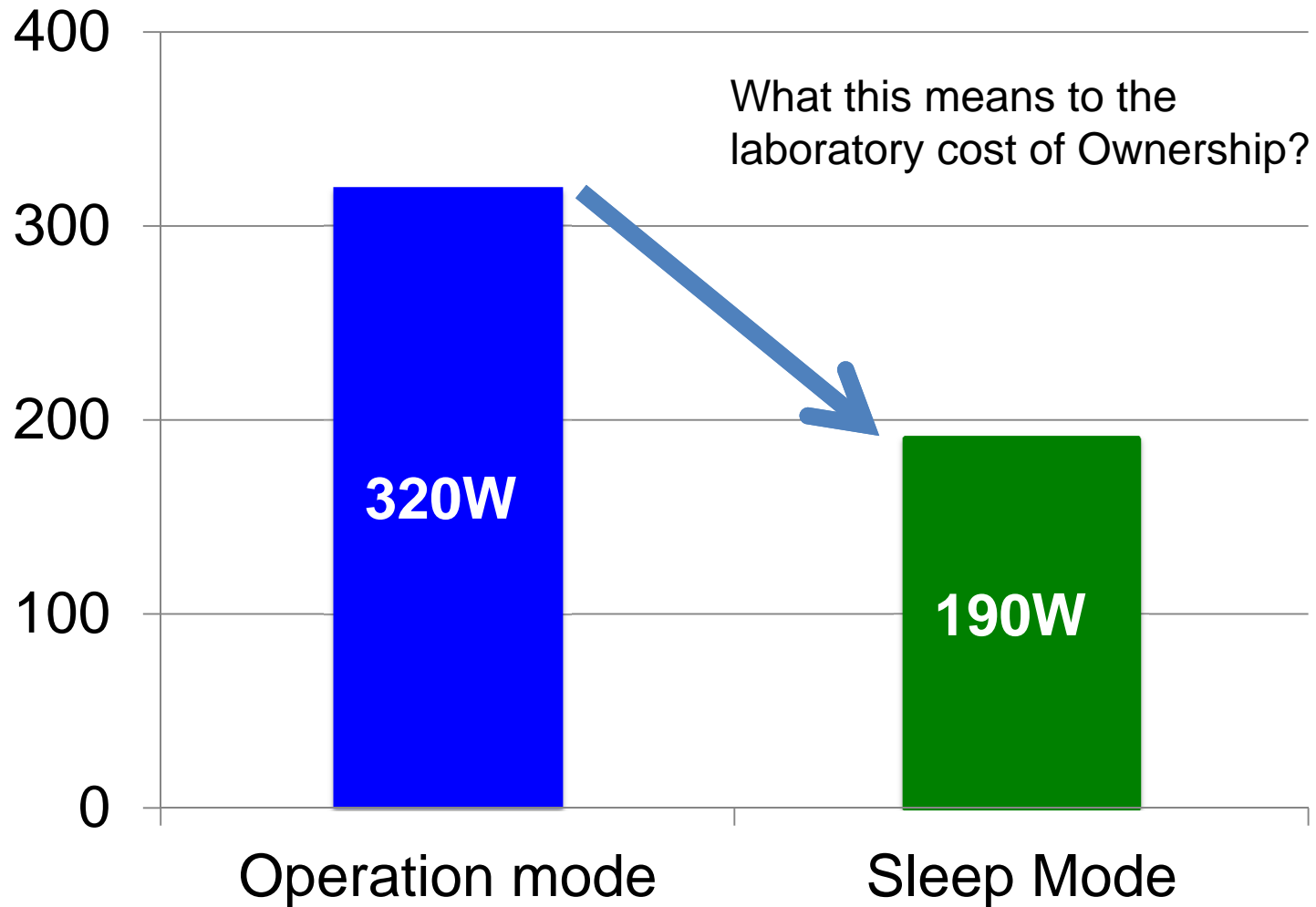


**New** IDP-3c

High performance  
Oil-free scroll pump  
NH<sub>3</sub> Cl compatible

**with Agilent Quality Standards**

# Better HW Features: Power consumption saving of MSD by Inverter Pump (DS42i)



# Example of Cost simulation of saving power with Agilent 7890B/5977A at Environmental Lab

## Assumptions

- GC/MSD cycle time: 40 minutes (30min for data acq./10 min for cool down & prep)
- Sample: 30 samples/day
- Working day: 5 days/week, no operating on Sat/Sun

GC/MSD **wake time** = 40 min/sample \* 30 samples \* 5 days/week  
= 6,000 min/week = **100 hours/week**  
= **32 kWh**

GC/MSD **sleep time** = 168 hours/week – 100 hours/week = **68 hours/week**  
= **13 kWh**

**Total Power usage (1 week) = 45 kWh (with Sleep mode)**

**Total Power Usage (1 week) = 54 kWh (without Sleep mode)**

**Weekly MS Power saving per instrument = 9kWh (17%)**

# Key features of “Direct Communication” between 7890B and 5977A



## Sleep/Wake mode

- In sleep mode, MSD will go to the Power save mode (320W → 190W)
  - Laboratory Operating Cost Reduction (see the previous slide for detail)

## Safety Control

- If 5977A MSD Pumps fail, 7890B will shut off the gas supply for protecting GC/MSD system.
  - Helium Carrier gas : This prevents wasting gas.
  - Hydrogen Carrier gas: This reduces the risk of explosion in MSD.

## Cool Down MSD Faster (FastVent)

- When MSD Vent Cycle started, GC will cool down MSD analyzer at the optimum gas flow to MSD.
- → Less time to wait for cooling down MSD analyzer (40 min → 25 min when cooling down from 280C to 100C)



# Expanded Capabilities: 7890B/5977A

## Improving the Industry Standards for GC/SQ

### Higher Ion Count

- Extractor EI source
  - Increased ion count
  - S/N > 1500:1
  - IDL < 10 fg (*industry first*)
  - Improved thermal profile

### Better Vacuum by Agilent Pump

- New Agilent V304 Turbo
  - TwisTorr
    - Molecular Drag technology
  - Improved pumping light gases
    - H<sub>2</sub> (220 L/s); He (255 L/s)
  - Agilent Floating Suspension
    - Longer life
    - Quiet operation



### Safety/Protection

- Vacuum failure → Shut off gas (He or H<sub>2</sub>)

### Save Time & Energy

- Sequence Scheduler
  - “Standby” post run
  - “Ready to run” from standby
  - “Bake Mode” as needed
- Factory backflush option
- Fast Vent capability
- DS42i rough pump “standby” power reduction

### Faster Scanning

- 20,000 u/s
  - But never equal to the 7200 Q-TOF speed or ‘scan’ sensitivity

### 7890B GC Features

- EMF, calculator, etc.

All of 5975’s features also built in 5977, such as “Gold Quad”, SIM/Scan, Gain tuning, Triple-Axis Detector, Trace Ion Detection, Auto CI, etc..



**RESOLVE**  
YOUR SEARCH  
FOR VALUE