



Guidelines for Trouble Shooting and Maintenance of AA Systems

Presented by Eric Vanclay,
Atomic Spectroscopy
Consumables Product
Manager

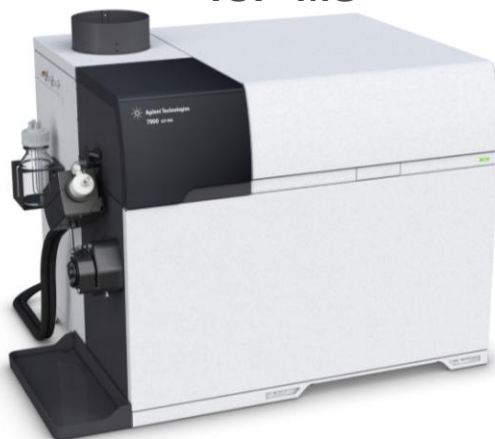
Today's Agilent: Atomic Spectroscopy

World's best, most complete atomic spectroscopy portfolio!

ICP-OES



ICP-MS



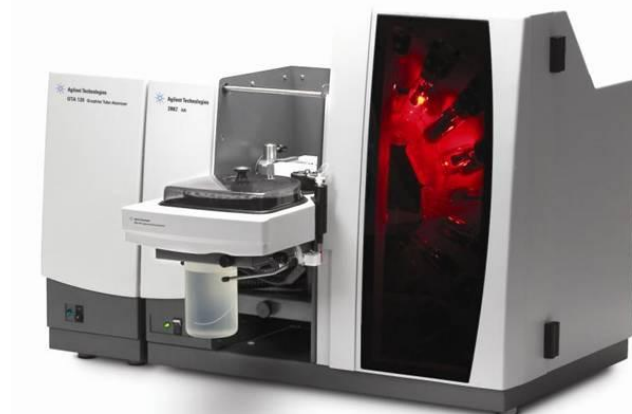
ICP-QQQ



Flame AAS



Graphite Furnace AAS



4200 MP-AES



New Agilent 7900 ICP-MS Performance Highlights

Rewriting the rules on ICP-MS

Better Analytical performance experience

- Ultra high matrix tolerance
- Superior sensitivity and lower background noise
- Wider dynamic range
- New Productivity Option (ISIS 3)
- Ultra fast scan speed for Single Nanoparticle analysis

Better Software experience

- ICP-MS MassHunter 4.1
- Method Wizard
- Mobile device support

Better Support experience

- Familiarization Tutorials/Videos
- Remote Advisor support



Next Generation 4200 MP-AES Performance Highlights

Change is in the Air!

Better Analytical performance experience

- 2nd generation of proven MP-AES technology
- Robust torch design for superior analytical performance with complex matrices
- Mass flow control of nebulizer gas and robust sample introduction
- Enhanced accuracy and long term stability with tough samples
- Expands the application range of MP-AES

Better Software experience

- MP Expert V1.2
- Intuitive software interface
- Advanced features in the 'PRO' pack

Better after sales experience

- No flammable gases or ongoing gas costs
- Runs from Nitrogen extracted from air using Agilent's nitrogen generator



Agilent's Atomic Spectroscopy Portfolio - Features

Flame AA



Lowest price

- Single element
- DLs typically ~100's ppb
- Fast (for 1 element)
- Good elemental coverage
- Low running cost

MP-AES



Lowest running cost

- Multi element
- DLs typically single to 10's ppb
- Faster
- Broader elemental coverage
- Lowest running cost

Graphite Furnace AA



Trace levels at lowest price

- Single element
- DLs typically 10's to 100's ppt
- Very slow
- Limited elemental coverage
- Moderate running cost

ICP-OES



Fastest measurement

- Multi element
- DLs typically single ppb
- Very fast
- Can measure most elements
- High running cost

ICP-MS



Broadest coverage

- Multi element
- DLs typically single or sub-ppt
- Fast
- Can measure almost all elements
- Highest running cost

Lowest

Selling Price

Highest



Agilent Technologies

Common AA Problems Reported by Customers

Sensitivity:

- Sensitivity is worse than it used to be
- I have a new application and I can't get the sensitivity I need
- How come I can't get the instrument to meet published detection limits?

Precision

- Sensitivity is acceptable but precision is terrible

High noise

- Can get the "right answers", but very noisy signal – this is also giving bad precision.

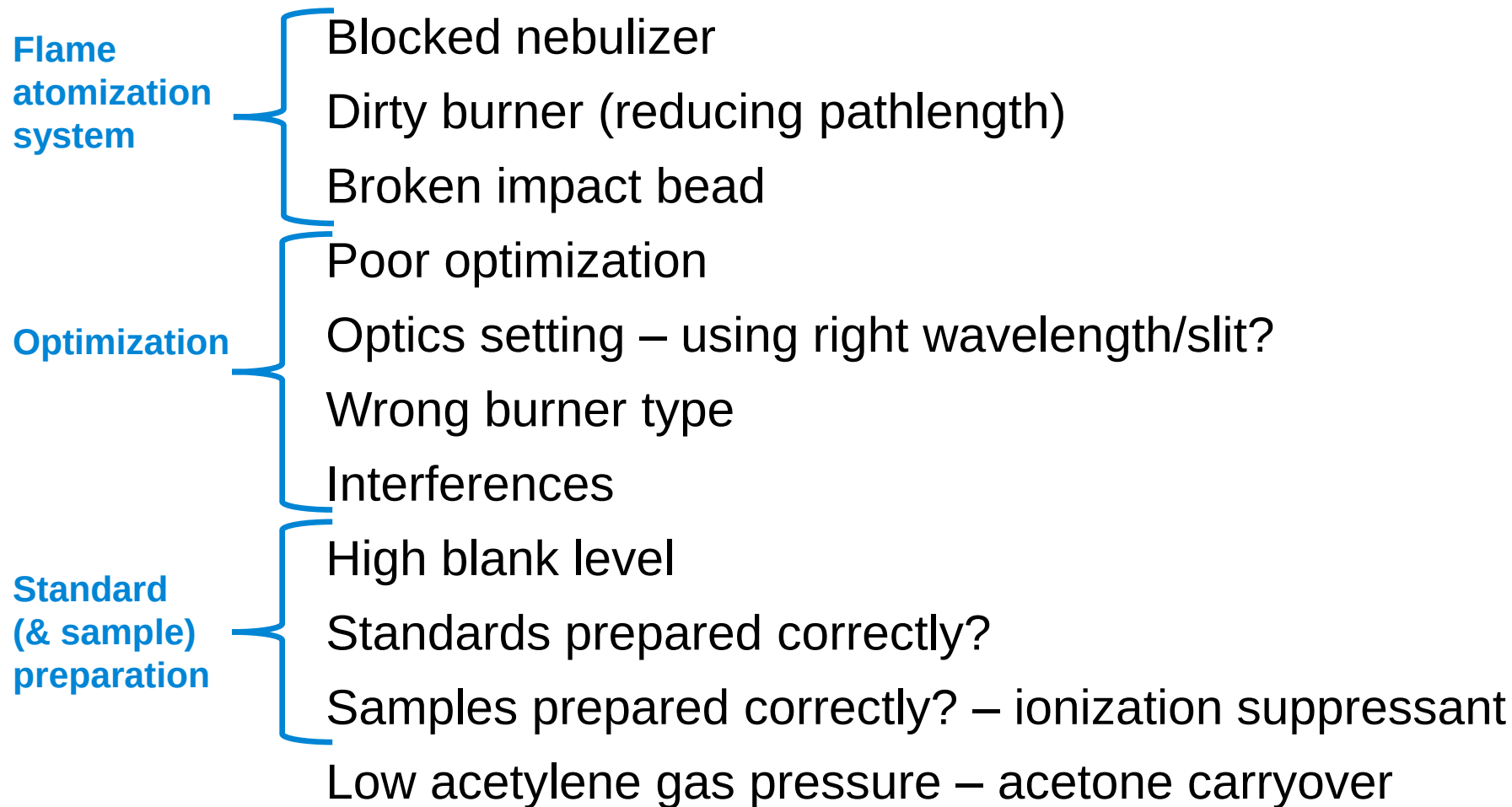
Accuracy

- Instrument does not give the "right" results.

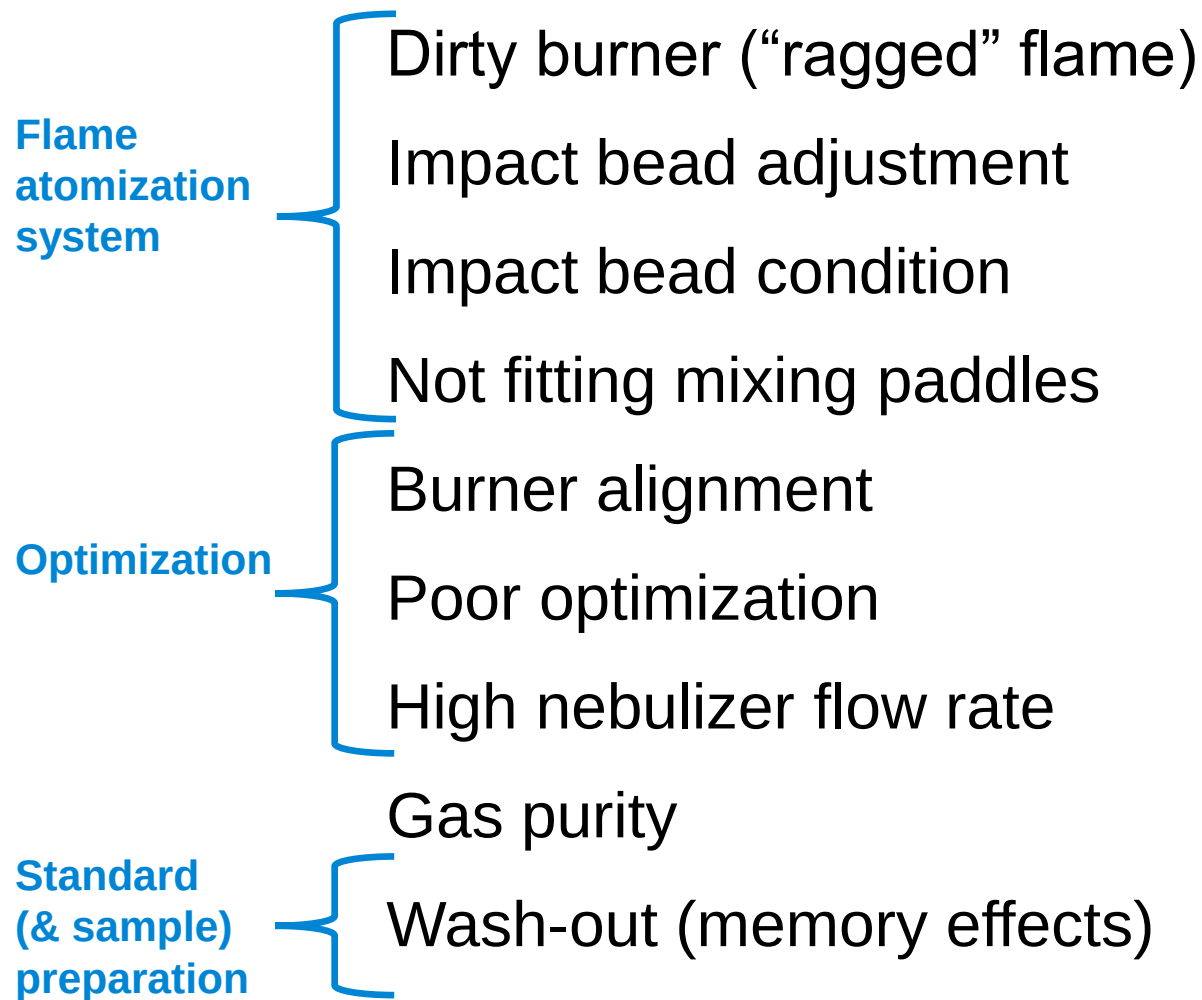
Poor Sample Throughput

- The instrument throughput needs to improve
- Burner blocks too quickly

Causes of Poor Flame AA Sensitivity



Causes of Poor Flame AA Precision



Causes of High Noise in Flame AA

Flame atomization system

Dirty burner (“ragged” flame)

Impact bead adjustment

Impact bead condition

Not fitting mixing paddles

Optimization

Burner alignment

Poor optimization – especially the HC lamp

Wrong lamp operating current

High nebulizer flow rate

Dirty optics

Gas purity

Standard (& sample) preparation

Incomplete digestion – particles in solution

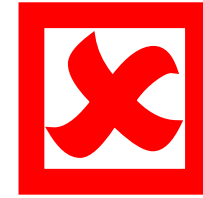


Flame Atomization System Tips



Do:

- Check optimization each analysis
- Check/monitor the nebulizer uptake
- Check/adjust the impact bead
- Check the blank reading
- Rinse between samples & at the end of the run
 - Rinse should match sample matrix
- Clean the burner/spray chamber regularly
 - Inspect condition of the impact bead
- Follow analytical recommendations in “cookbook”



Don't:

- Assume system is still optimized
- Assume nebulizer flow rate is the same
- Use a simple water blank
- Wait until you have blockage before cleaning



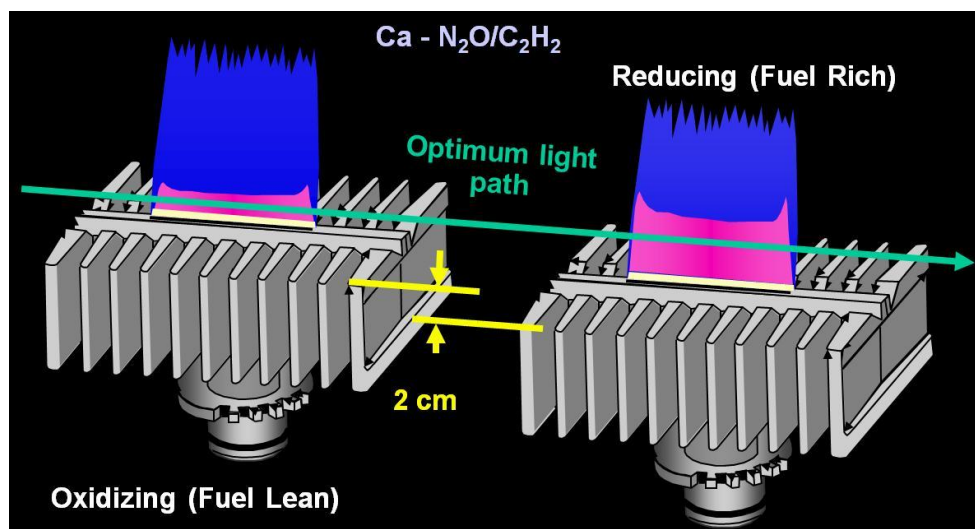
Burner Alignment Tips

Burner

- Burner position must be optimized (vertical, horizontal & rotational positioning)
- Use “target area” on burner alignment card to ensure light beam runs parallel to burner slot
- Adjust burner height while aspirating a standard and optimize for maximum signal
 - Optimum height varies depending on flame chemistry
- Burner type changes path length (for air/acetylene elements)



Optimum viewing height for Ca

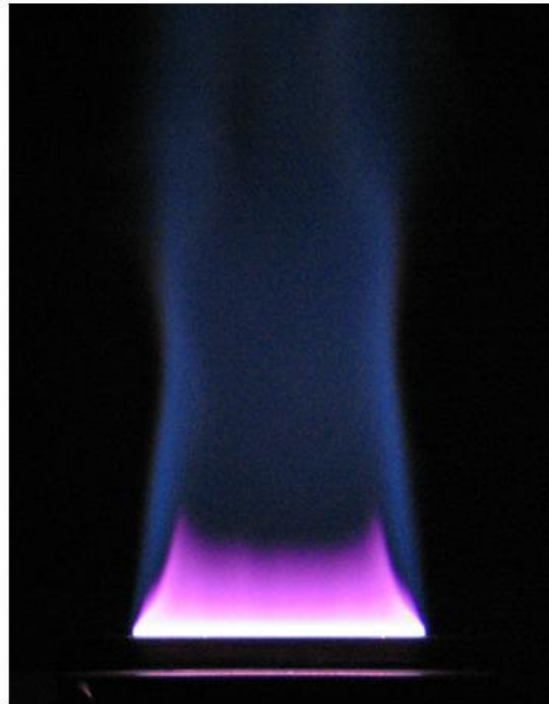


Optimization of the Nitrous Oxide/Acetylene Flame



Lean

Reduce acetylene flow



Stoichiometric



Rich

Add extra acetylene



Cleaning the Atomization System

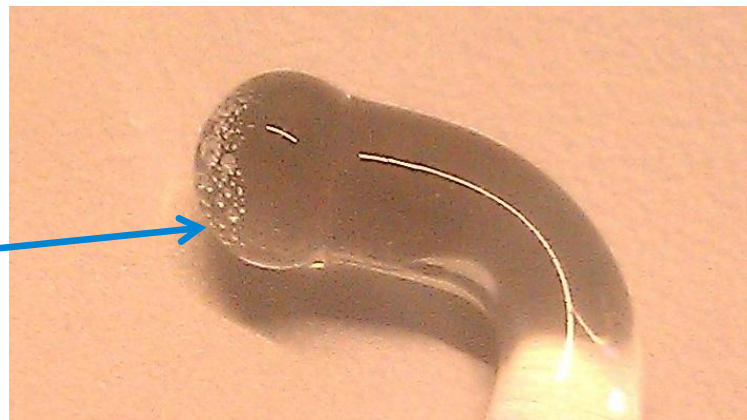
Cleaning the burner:

- Polish with metal polish like “Brasso”
 - Watch the video (#4)



Cleaning the spray chamber:

- Wash in detergent solution
- Check the impact bead condition & setting
 - Replace if badly pitted



Removing nebulizer blockage:

- Disassemble completely and wash in detergent
- Reassemble and test



Impact Bead Setting

Loosen bead securing screw

Turn external adjuster fully clockwise

Set the bead position

- 2 critical settings
 - Bead should be centred in front of venturi
 - Distance between bead & venturi should be $\sim 0.1\text{mm}$ (thickness of sheet of paper)

Tighten adjuster to lock bead in place



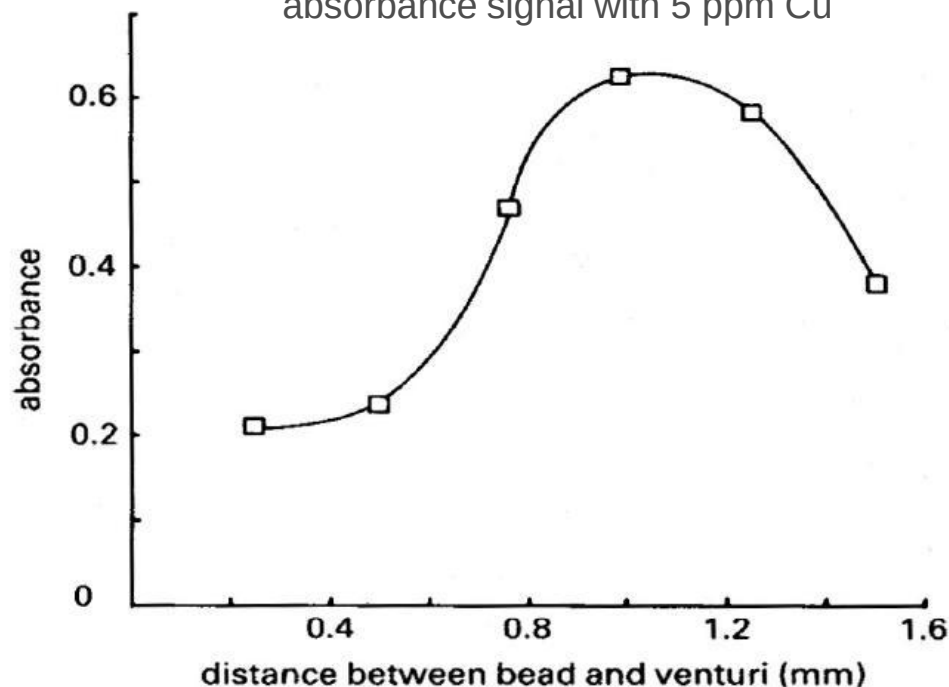
How Impact Bead Setting Impacts Sensitivity?

Use external bead adjuster

- Rotate this adjuster
 - Clockwise adjustment brings bead closer to nebulizer
- “Tune” performance to suit your application
 - For most applications, adjust for best mix of signal + precision
 - For best sensitivity, position bead away from nebulizer (anticlockwise)
 - For higher TDS samples, adjust for max. sensitivity – then rotate adjuster at least $\frac{1}{2}$ turn clockwise (closer to nebulizer)



Effect of impact bead position for absorbance signal with 5 ppm Cu



Flame Atomization – Rec. Settings

For most flame AA applications:

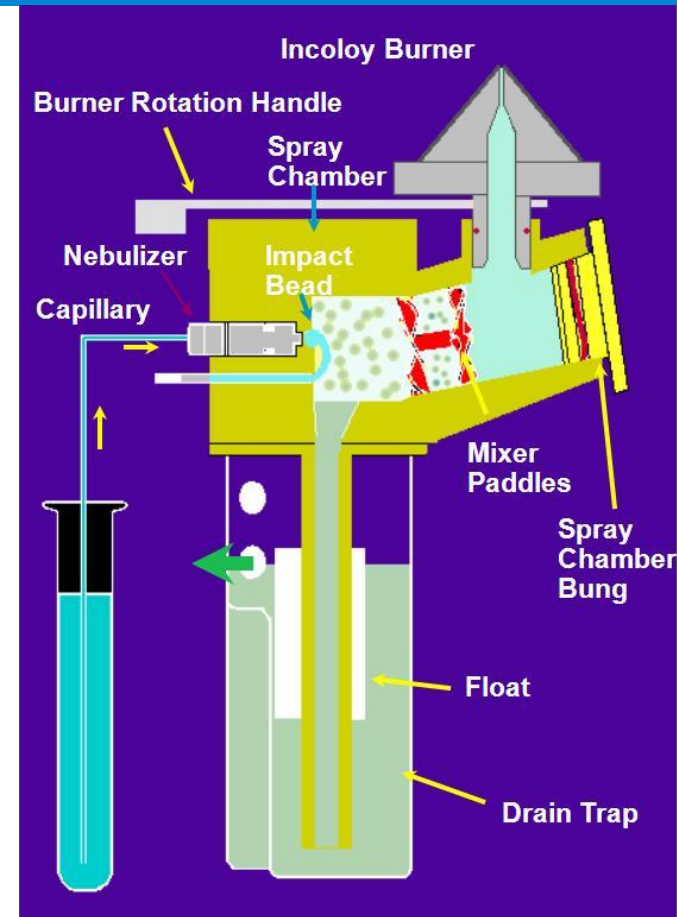
- Fit the mixing paddles
- Optimize the impact bead for best sensitivity
- Use narrow bore capillary tubing

For highest flame AA sensitivity:

- Remove mixing paddles
- Adjust impact bead further away from the nebulizer
- Use wide bore capillary tubing (highest uptake rate)

For higher TDS samples:

- Fit the mixing paddles
- Adjust impact bead $\frac{1}{2}$ to 1 turn clockwise (towards nebulizer) from optimum sensitivity position
- Use wide bore capillary tubing (to reduce chance of blockage)



Recommended Procedures at End of the day

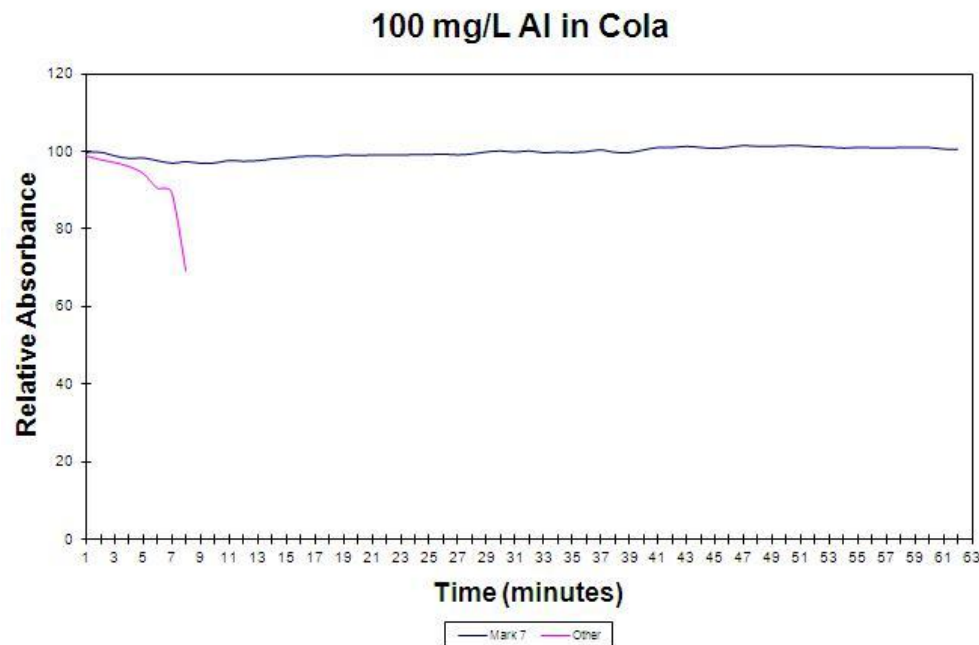
1. Aspirate distilled water for a few minutes before shutting off flame
2. Allow burner to cool
3. Remove burner and clean by running water through it
4. Dry burner by shaking
5. Pour 500 mL of water into spray chamber, through burner socket
6. Replace burner
7. Empty waste vessel

Agilent Flame AA Performance – Benefits

Flexibility, Ease of Use & Superior Flame Performance

Tunable performance means...

- Highest flame sensitivity: > 0.9 Abs. for 5 mg/L Cu
- Best precision: < 0.5 % RSD using 10 x 5 s readings
- Extended operation with difficult samples
- No loose gas hoses and no tools required for gas connection
- Fast change-over to furnace operation (manual - < 30 s)



Causes of Poor Furnace AA Sensitivity

Furnace workhead

- Missed injection
- Aged (or damaged) tube in use
- Wrong electrodes fitted

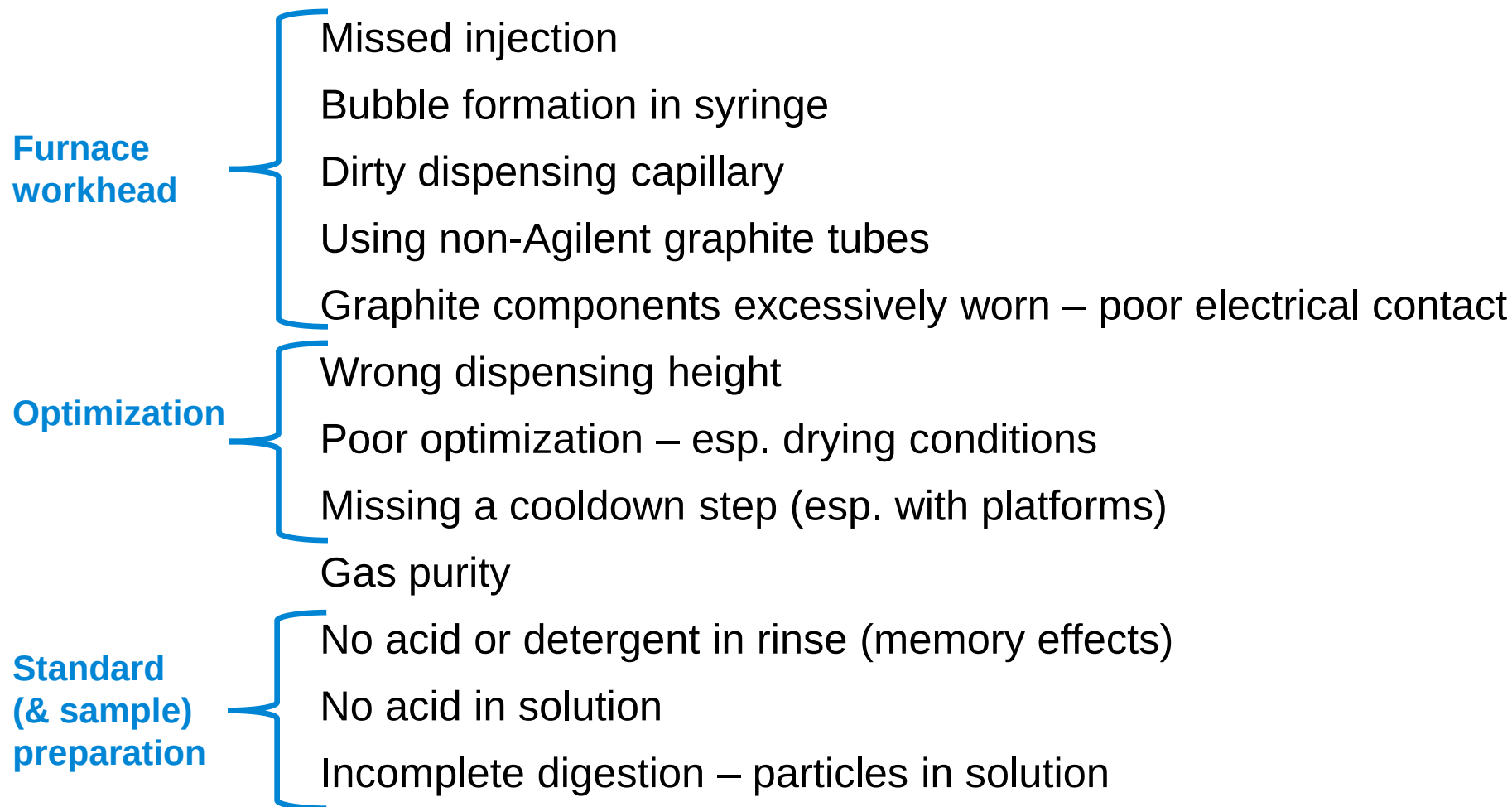
Optimization

- Poor optimization – esp. drying conditions
- Workhead incorrectly aligned
- Optics setting – using right wavelength/slit?
- No modifier (or incorrect) modifier used
- Use of nitrogen as inert gas

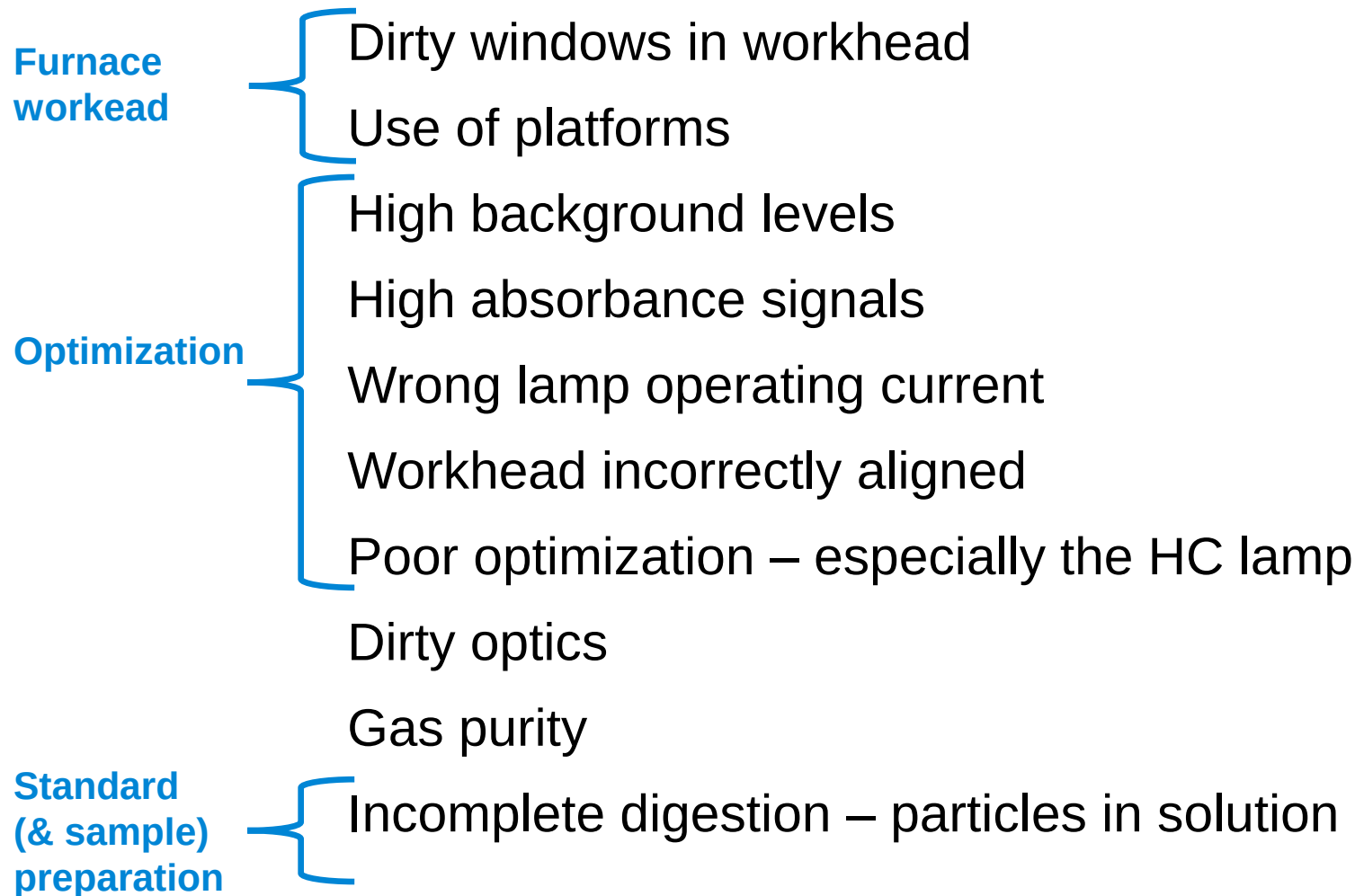
Standard (& sample) preparation

- No acid in solution
- High blank level
- Standards prepared & stored correctly?
- Samples prepared correctly – digestion?

Causes of Poor Furnace AA Precision



Causes of High Noise in Furnace AA



Furnace AA System Tips



Do:

Check optimization each analysis

Check/monitor the dispensing height

Ensure the rinse solution has 10 drops
conc. HNO_3 + 5 drops Triton X-100

Remove residue from the dispensing
capillary

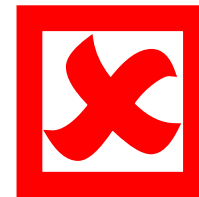
Check/monitor the graphite tube

Check the blank reading

Clean the workhead regularly

- Inspect condition of the graphite components

Follow analytical recommendations in
“cookbook”



Don't:

Assume system is still optimized

Assume dispensing height is the same

Use a simple water rinse

Start analysis with a dirty capillary tip

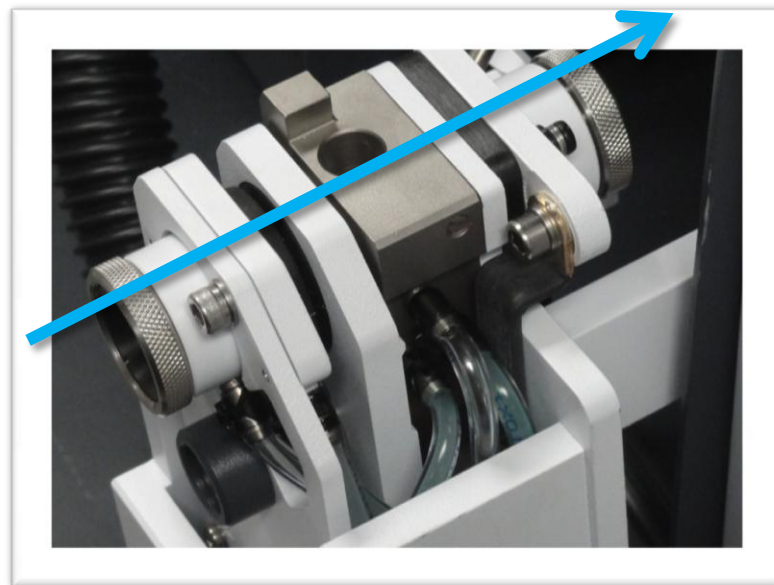
Start analysis with a tube near the end of
it's life



Furnace AA System Tips - Workhead Alignment

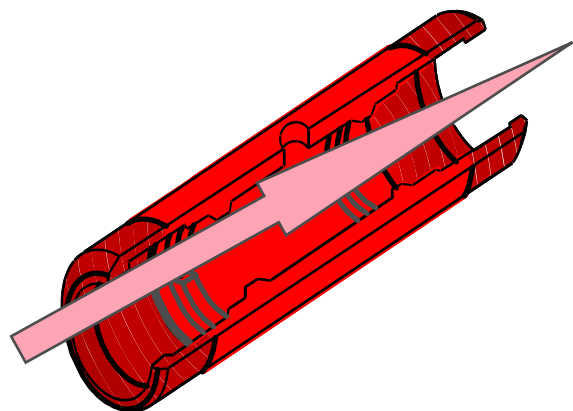
Furnace workhead

- Workhead position must be optimized (want light beam to pass through centre of graphite tube)
 - Align lamp first (no workhead), then place workhead in position and align



Sample Dispenser settings

- Carefully adjust injection depth – easy with the furnace camera



**Light Beam Aligned
Through Center of
Graphite Tube**



Furnace AA System Tips - Setting Injection Depth

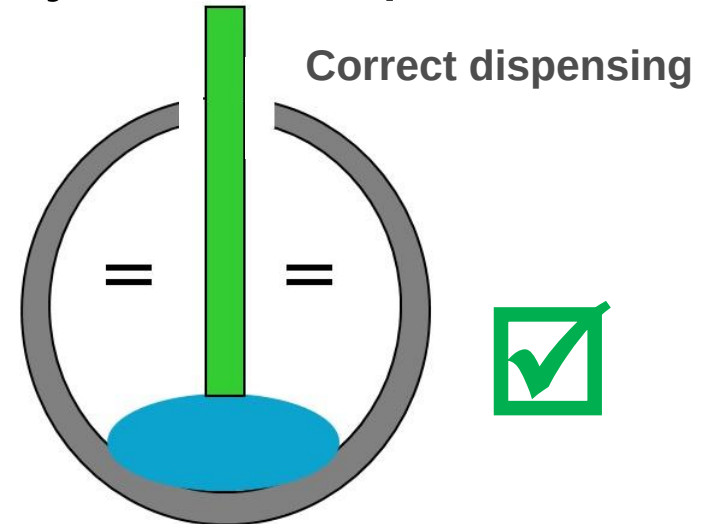
Capillary tip must remain in contact with solution during dispensing

Reduce dispensing height if sample spreads due to low surface tension

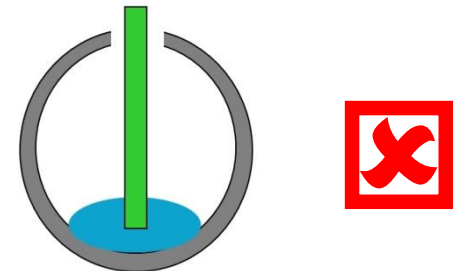
Ensure there is no liquid on the outside of the capillary after dispensing

Ensure there is no liquid inside the capillary tip after dispensing

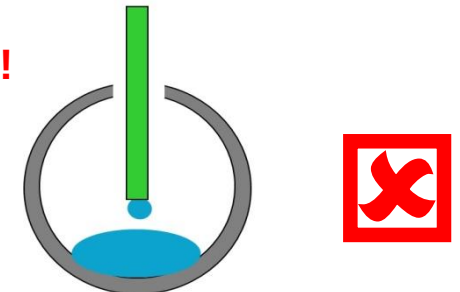
Sample should remain as a drop in the centre of the tube



Pipet tip too low!



Pipet tip too high!



Furnace AA System Tips – Tube Conditioning

Why condition the tubes before use?

- Helps remove residual contamination
- Gently “beds” a new tube in
 - Important when determining concentrations near detection limit
 - Also important with some complex matrices
- Critical when using modifiers
 - Helps to build up coating inside the tube
 - Improves efficiency of the modifier
- Improves reproducibility



Recommended process

- Use “Tube Condition” facility (or otherwise, fire tube 5-10 times using either reagent blank or representative sample)



Furnace AA System Tips – Method Parameters

What to Check?

- Furnace parameters
 - Set appropriate drying temperature and time (2-3 sec/uL of solution injected)
 - Optimize ashing temperature using ashing study – use SRM optimization
 - Ensure inner gas flow “off” just prior to atomization



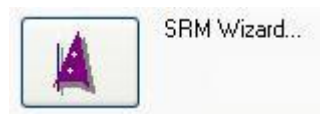
Does the sample sizzle or splatter during the dry stage?

- Listen for the sound
- Use the mirror or furnace video to monitor the sample drying



Furnace AA System Tips – Method Optimization

- SRM “Wizard” automates furnace optimization
- Optimizes absorbance as a function of ashing and atomization temperature
- Automatically creates a method using recommended conditions
- Reduces training requirements for new users



SRM Wizard...

SRM (Surface Response Methodology) Wizard

Set up the Ash and Atomiser start and stop steps, temperature and change temperature to create the values for the furnace methods. When finished, press 'Next'.

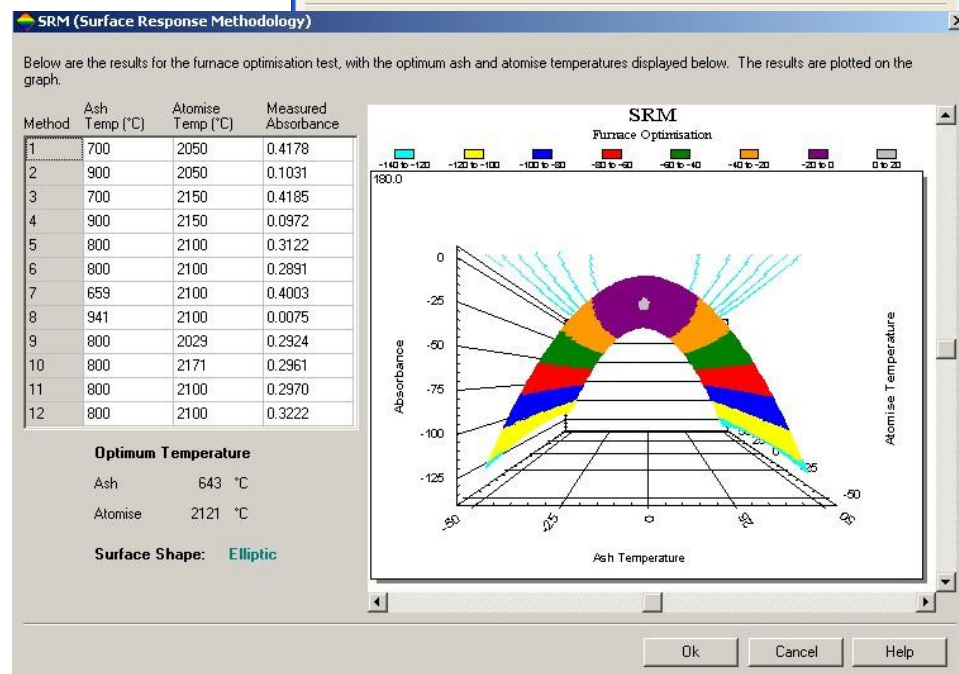
Ash

Start Step: 4
Stop Step: 6
Temperature °C: 400
Change °C: 100

Atomise

Start Step: 9
Stop Step: 11
Temperature °C: 2100
Change °C: 50

Step	Temp (°C)	Method	Ash Temp	Atomise Temp
1	85	1	300	2050
2	95	2	500	2050
3	120	3	300	2150
4	400	4	500	2150
5	400	5	400	2100
6	400	6	400	2100
7	200	7	259	2100
8	200	8	541	2100
9	2100	9	400	2029
10	2100	10	400	2171
11	2100	11	400	2100
12	2100	12	400	2100



Optimization results for Pb determination using phosphate modifier



Furnace AA System Tips – Reducing Sensitivity

May be required due to sensitivity of this technique:

- Switch to an alternate wavelength
 - Select a less sensitive wavelength (if available)
- Reduce sample volume
- Use slower ramp rate to atomization
 - Aim to broaden the peak during atomization
- Use low gas flow during the atomization step

Furnace AA System Tips – Factors Influencing Tube Lifetimes

Graphite components excessively worn – poor electrical contact

No tube preconditioning (or always using Tube Clean)

Sample matrix

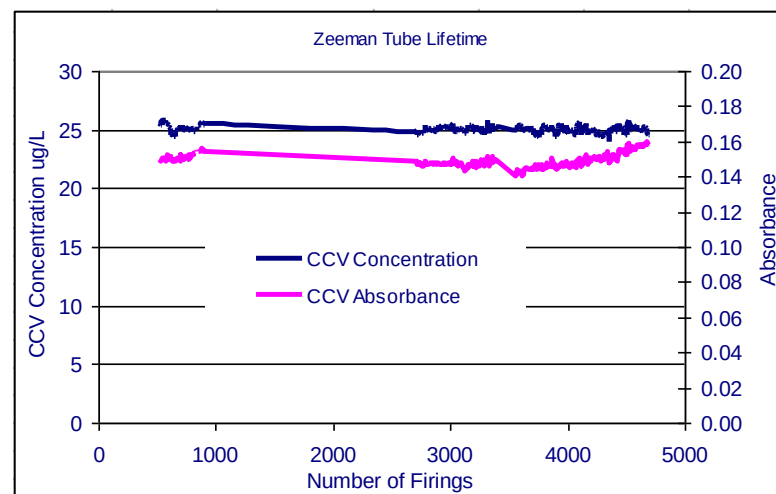
Inert gas used

- Argon gives longest life
- Nitrogen degrades tube faster due to oxygen presence

Temperature program used

Type of chemical modifiers used

- Powerful oxidizing agents degrade tube faster
 - Perchloric acid
 - Perchlorates
 - Sodium nitrate



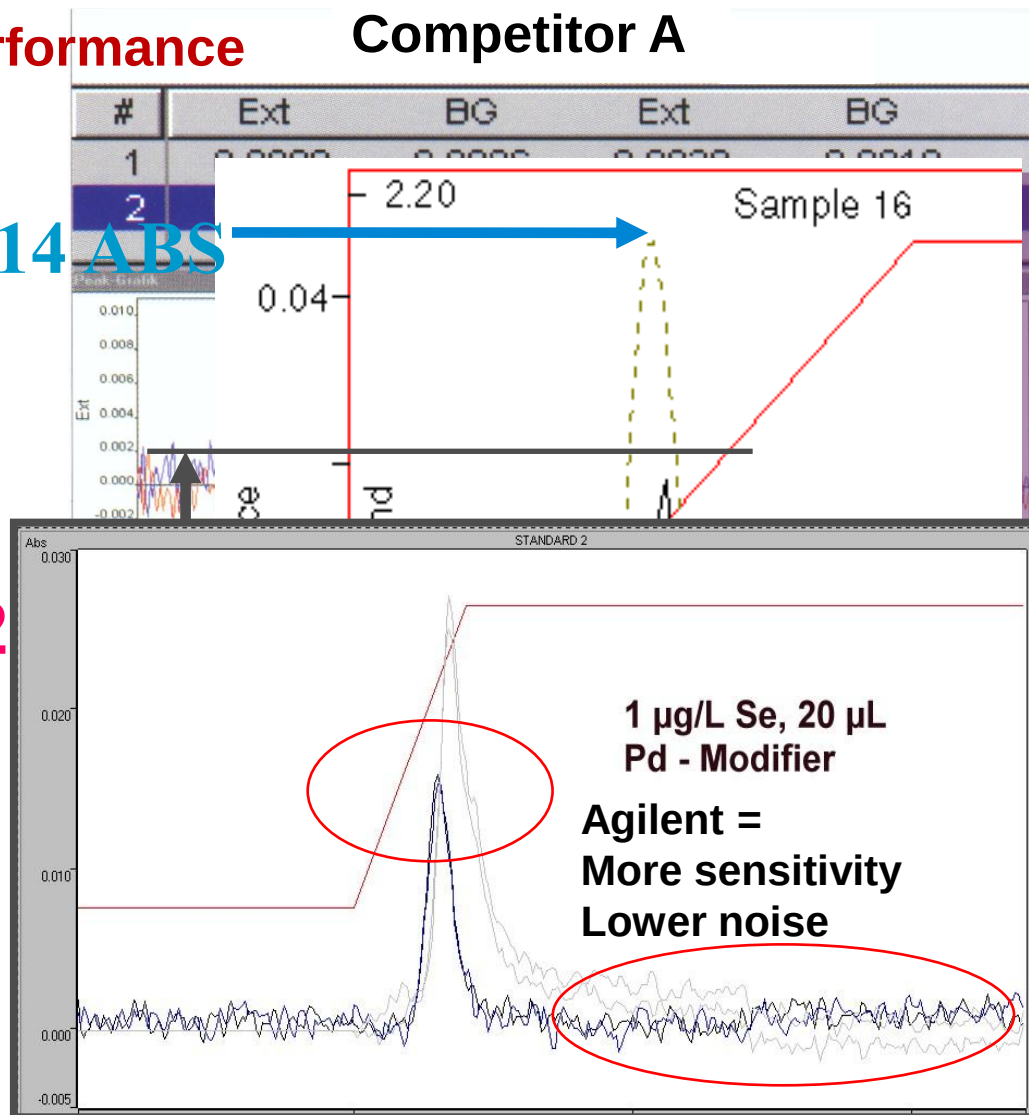
Agilent Furnace AA Systems – Benefits

Flexibility & Superior Furnace Performance

Competitor A

- Highest furnace sensitivity
- Best Zeeman correction capability:
< 2 % error at > 2 Abs. b'ground
 - Polynomial interpolation of the background
 - 100/120 readings every second
- Best capability to handle difficult samples

2.14 ABS



Tips to Improve Standard Preparation

- How are they prepared?
 - Ensure purchased standards are still within “Use By” date
 - Use calibrated pipettes and class ‘A’ volumetric flasks for dilutions
 - Periodically, check accuracy & reproducibility of your pipettes
 - Use de-ionized water (Type I - conductivity $\geq 18 \text{ M}\Omega/\text{cm}^3$)
 - Lower grades may have contamination
 - Use serial dilutions for preparing low concentrations from 1,000 ppm stock
 - Please don’t do large dilutions ($> 1:10,000$) in 1 step
- What concentration are they?
 - Low concentration standards have a finite life
 - Prepare ppb (ug/L) concentration standards daily from high conc. stock
 - Prepare low ppm (mg/L) concentration standards weekly
- How are they stored?
 - Plastic vessels ensure better stability
 - Stabilize with acid – low pH ensures better stability



Tips to Reduce Contamination

Contamination can come from anything that comes into contact with your sample during storage, digestion (dilution) and analysis



- Check reagent purity
 - Always buy the best reagents
 - Always check the certificate of analysis for elevated levels
 - Caution if buying in large quantities
 - Worst case – can use contaminated acid for cleaning
 - Ensure still within “use by” date
 - Reseal immediately after use

CERTIFICATE OF ANALYSIS																											
BASELINE [®] Nitric Acid																											
PRODUCT NUMBER: S020101							LOT NUMBER: 1211120							ASSAY (HNO ₃ , w/w): 68%													
1A		2A		3A		4A		5A		6A		7A															
3	Li	4	Be	5	B	6	C	7	N	8	O	9	F	10	Ne	11	Na	12	Mg	13	Al	14	Si	15	P	16	S
< 0.5		< 1		< 10												< 10											
19	K	20	Ca	21	Sc	22	Ti	23	V	24	Cr	25	Mn	26	Fe	27	Co	28	Ni	29	Cu	30	Zn	31	Ga	32	Ge
< 1		< 5		< 1		< 5		< 0.5		< 10		< 1		< 10		< 1		< 10		< 1		< 1		< 0.1		< 0.1	
37	Rb	38	Sr	39	Y	40	Zr	41	Nb	42	Mo	43	Tc	44	Ru	45	Rh	46	Pd	47	Ag	48	Cd	49	In	50	Sn
< 1		< 1		< 0.1		< 0.5		< 1		< 1			< 1		< 1		< 5		< 0.5		< 0.5		< 0.1		< 0.1		< 5
55	Cs	56	Ba	57	La	58	Ce	59	Pr	60	Nd	61	Pm	62	Sm	63	Eu	64	Gd	65	Tb	66	Dy	67	Ho	68	Er
< 0.05		< 1		< 0.05		< 0.05		< 0.05		< 1		< 1					< 1		< 1		< 1		< 10		< 0.05		< 0.05
ALL VALUES ARE REPORTED IN PARTS PER TRILLION (PPT)																											

- Other common contamination sources
 - Reagent water
 - Clean glassware?
 - Airborne dust in the lab.
 - Pipette tips
 - Don't insert pipette tips into your acids
 - Use natural tips – colored tips may increase contamination (Cu, Fe, Zn, Cd)
 - Powdered gloves (esp. for Zn)



Tips to Improve Accuracy of Results

- Sample preparation
 - Is the most appropriate digestion being used?
 - Are all of the analytes being quantitatively (and reproducibly) extracted and dissolved?
 - Many digestions are only partial extracts – efficiency will vary with the sample matrix
 - Some volatile analytes may be “lost” during digestion
 - Confirm by taking a solid certified reference material through your preparation and analysis procedure
- Is the digest stable – or are you seeing any precipitates or a suspension?
- Do you see any potential contamination from either reagents or the digestion equipment? e.g. especially with Si, B or Ca
 - Include a “Reagent Blank” with every sample batch to monitor



Tips for Cleaning Dirty Optics

Monitor the windows regularly

- Check lamp for fingerprints
- Check sample compartment windows for build-up of film/chemical residue



Smudges or chemical residue reduces light and increases noise

Cleaning the windows?

- Wipe clean with an optical tissue (as you would use to clean a camera lens)
- If necessary, use optical tissue moistened with ethanol



Cleaning end windows from furnace workhead

Flame AA – Recommended Maintenance Schedule

Daily:

- Check the gas delivery pressures & cylinder contents (esp. acetylene)
- Check exhaust system
- Check the nebulizer uptake rate and burner condition
- Clean burner compartment & instrument
- Empty waste container

Weekly:

- Clean burner (or earlier if required)
- Disassemble flame atomization system and clean
 - Check condition of O rings and impact bead (no pitting)

Monthly:

- Check windows in sample compartment (clean if necessary)

Furnace AA – Rec. Maintenance Schedule

Daily:

- Check the gas delivery pressures & cylinder contents
- Check exhaust system
- Check condition of the graphite tube – replace as necessary
 - When replacing the tube, check the condition of the electrodes
- Check dispensing capillary “free” and syringe
- Top up rinse reservoir as required
- Clean the workhead around the sample injection hole
- Empty waste container

Weekly:

- Check furnace workhead windows (clean if necessary)

Overview - Key Consumables for AA

All instruments:

- HC lamps
- AA standard solutions

Flame AA systems:

- Glass impact beads, burner cleaning strips, nebulizer components, capillary tubing, burners etc
- Ionization suppressant / buffer solutions
- With the SIPS dilution system – SIPS pump tubing and transfer tubing
- With an autosampler - sample tubes, racks, probes and transfer tubing

Graphite furnace AA systems:

- Graphite tubes
- Sample vials, dispensing capillary and syringe for autosampler
- Matrix modifiers

Vapor generation AA systems:

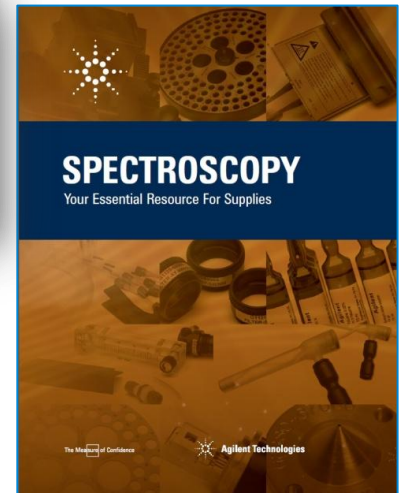
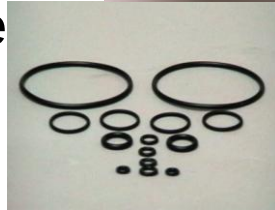
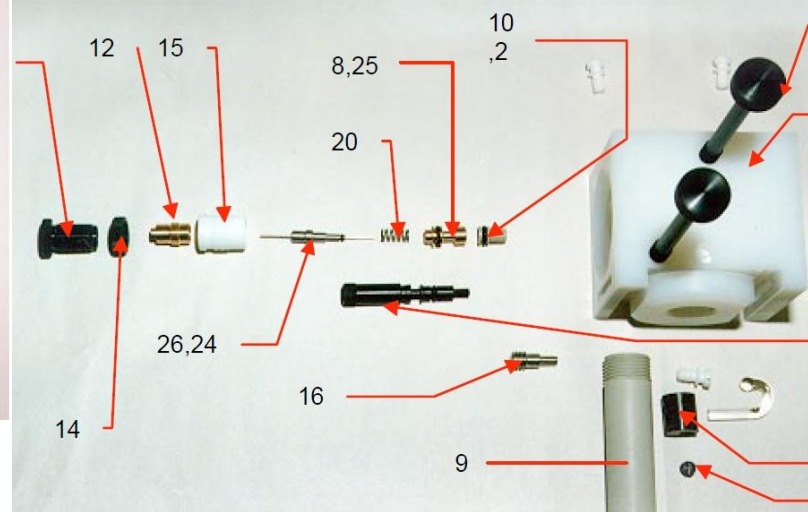
- Quartz atomization cells
- Peristaltic pump tubing
- Connecting tubing



Flame Atomizer - Mark 7 Design

Key consumable items requiring frequent replacement:

- O rings
- Glass impact beads
- Burner cleaning strips
- Nebulizer capillary kit
- Nebulizer venturi kit
- Capillary tubing
- Nebulizer cleaning wire
- Liquid trap assembly
- Mixing paddles
- Burners



Where to find ordering details?

- Agilent website – dedicated webpage: [Mark 7 Spares](#)
- Agilent Quick Reference Guide for AA
- Agilent Spectroscopy Supplies Catalogue

Agilent AA Consumable Kits

Part Number	Description	Content
190034100	Flame AA operating supplies kit (for Mark 7 atomization system)	Nebulizer venturi, capillary kit, nebulizer block (excl. integral nebulizer), Glass impact beads Capillary tubing O ring kit Mixing paddles Burner cleaning strips
190065400	SPS 3 Flame Autosampler operating supplies kit	0.8 mm ID inert probe 2 packs grey/grey 3 bridged pump tubing (12/pk) Connecting tubing, drain tubing and capillary tubing Rinse reservoir (10 L) 1 pack 16 mm OD polypropylene tubes (1000/pk) 3 sample racks for 30mm OD tubes (21 positions) 1 pack 30mm OD polypropylene sample tubes (500/pk)
190067900 (for GTA 120); or 190068000 (for GTA 120 Zeeman)	Graphite Furnace AA operating supplies kits	2 sets graphite electrodes Graphite shroud 5 packs Omega tubes (each 10/pk) 100 uL syringe for PSD 1 pack of capillary assemblies for PSD (5/pk) 1 pack of plastic beakers (5/pk) 2 packs 2 mL furnace vials (1000/pk)
190025200	VGA 77 Vapor Generation AA operating supplies kit	2 sets tubing and connector kits 2 packs sample pump tubes (12/pk) 2 packs reagent pump tubes (12/pk) 1 set replacement pump beds 1 replacement AA gas-liquid separator 1 Hg Flow Through Cell (1/pk) 2 packs hydride absorption cell (2/pk) 1 spare AA hydride module
190025400	SIPS Flame Dilution System operating supplies kit	2 ea 500 mL constant pressure vessel 1 x 1L diluent bottle 1 x 3 way tee piece assembly 1 Pack SIPS pump tubing (6/pk) 1 Pack Pump Bands (10/pk) 1 SIPS tubing kit



Where to Find the Right Consumable?

Analytical Consumables: Consumables & Supplies

1-800-227-9770 (Option 1,1)
www.agilent.com/chem/contactus

On-Line resources:

[Atomic Absorption Supplies](#)

[Mark 7 Sample Introduction Spares](#)

[AA FAQs](#)

[ICP-OES Parts & Supplies Portfolio](#)

[ICP-MS Supplies](#)

[Instrument Parts & Supplies](#)

[Atomic Spectroscopy Application Notes](#)

[Recorded Agilent e-Seminars](#)

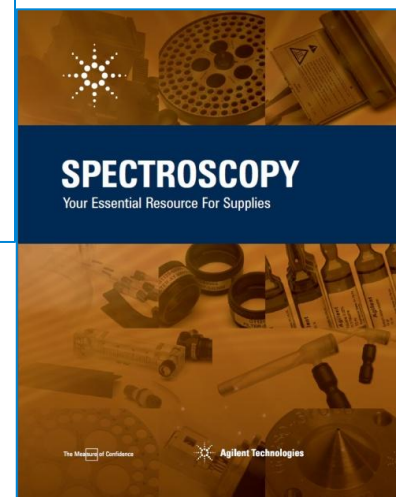
Agilent Quick Reference Guide for AA (pub. # [5990-9476EN](#))

Agilent Spectroscopy Supplies Catalogue (pub # [5991-1060EN](#))

Instrument User Manuals

Agilent Assist: Instrument Sales & Services

1-800-227-9770 (Option 1,3)
www.agilent.com/chem/contactus



Summary – To Achieve Quality Data

- Most “instrument” failures occur in the sample introduction area
 - Includes
 - Burner
 - Spray chamber
 - Nebulizer
 - All tubing
 - Drain Assembly
- Improper maintenance of this area can result in poor data quality
- Frequently less experienced analysts can fail to recognize problems resulting in productivity losses
- Establishing maintenance procedures can prevent problems



Questions?



Agilent MP-AES



Agilent AAS



Agilent ICP-OES



Agilent ICP-MS

The Market Leaders in Atomic Spectroscopy

Other Support Resources

Are you a subscriber to the Access Agilent newsletter?

Access Agilent is:

- A monthly e-newsletter newsletter tailored to your preferences
- Includes in-depth articles, new products, literature, offers and events

To register, use this link to the registration form on the Agilent website (or ask your Agilent representative):

<https://www.chem.agilent.com/en-US/layouts/Agilent/Security/ProfileSelectionPage.aspx>

[Receive this email as text only.](#)

 **Agilent Technologies**

Access Agilent

January 2014 [Update My Profile](#) | [Contact Us](#)

Dear Eric Vanclay,

Welcome to your new monthly newsletter, personalized to your selected profile. If you would like more information on how to select your information preferences and update your profile, [click here](#).

In-Depth Articles



Water, water everywhere... but is it fit to drink? [Read article](#)



Better USP <467> analysis of residual solvents [Read article](#)



Gain better food science insight from better data using the new Agilent ProPulse NMR System [Read article](#)



Reduce detection limits for pesticide residues in foods with the new Agilent 7000C Triple Quadrupole GC/MS System [Read article](#)



OpenLAB ELN makes it easy to capture, manage, and share scientific results [Read article](#)



Large Valve Oven GC system meets the petrochemical industry's growing demands [Read article](#)



Choosing the right filter for food and drink analysis made easy by Agilent [Read article](#)

Special Offers

Get up to 30% off for Bond Elut SPE or QuEChERS products Use promo code 9174 | Expires March 31, 2014

New Literature

Application Notes

- Determination of Ochratoxin A in Roasted Coffee According to DIN EN 14132
- Optimizing Sample Preparation for LC/MS/MS of Pesticide Residues in Herbal Teas
- Identification and Quantitation of PCB Aroclor Mixtures in a Single Run Using the Agilent 7000B Triple Quadrupole GC/MS
- Simultaneous determination of multiclass antibiotic residues in Nile Tilapia (*Oreochromis niloticus*) by LC-MS/MS
- Volatile Profiling of U.S. Cabernet Sauvignon Wines Using HS-SPME and the Agilent 5975 Series GC/MSD System: Relating the Chemical Profile to Sensory Properties
- Automated Sample Preparation in Quality Control of Eye-Drop Formulation

