



Guidelines for Troubleshooting and Maintenance of AA Systems

Presented by Eric Vanclay,
Atomic Spectroscopy
Consumables Product
Manager



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

Flame atomization system

- Blocked nebulizer
- Dirty burner (reducing pathlength)
- Broken impact bead

Optimization

- Poor optimization
- Optics setting – using right wavelength/slit?
- Wrong burner type
- Interferences

Standard (& sample) preparation

- High blank level
- Standards prepared correctly?
- Samples prepared correctly? – ionization suppressant
- Low acetylene gas pressure – acetone carryover

Causes of Poor Flame AA Precision

Flame atomization system

- Dirty burner (“ragged” flame)
- Impact bead adjustment
- Impact bead condition
- Not fitting mixing paddles

Optimization

- Burner alignment
- Poor optimization
- High nebulizer flow rate

Standard (& sample) preparation

- Gas purity
- Wash-out (memory effects)

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

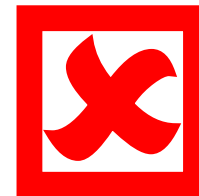
Rinse between samples & at the end of the run

- 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

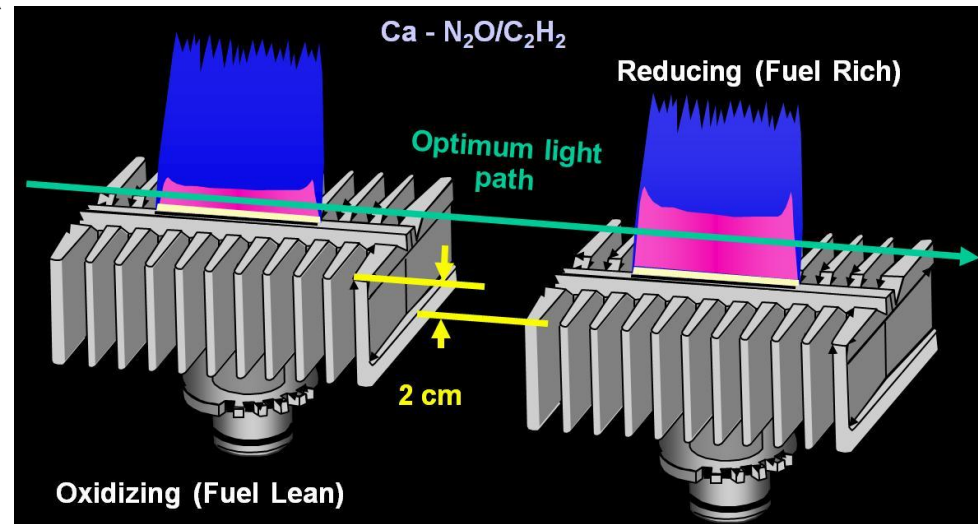
Burner Alignment Tips

Burner

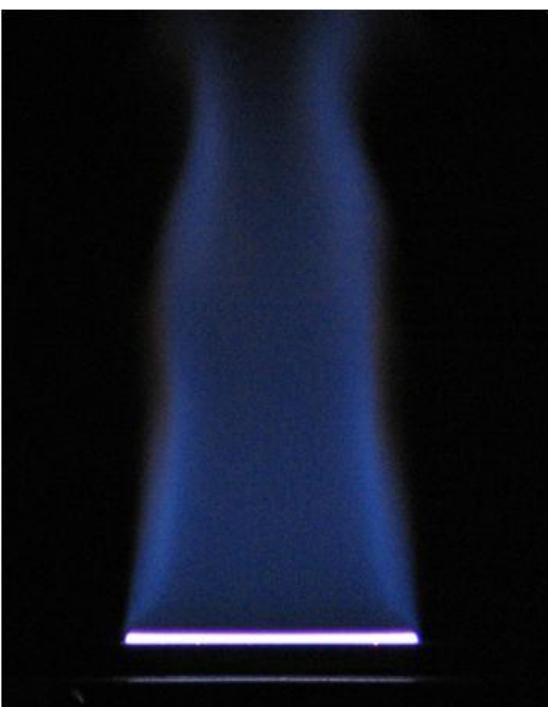
- Burner position must be optimized (vertical, horizontal & rotational positioning)
- Use “target area” on burner alignment card to ensure light beam runs parallel to burner slot
- Adjust burner height while aspirating a standard and optimize for maximum signal
 - Optimum height varies depending on flame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

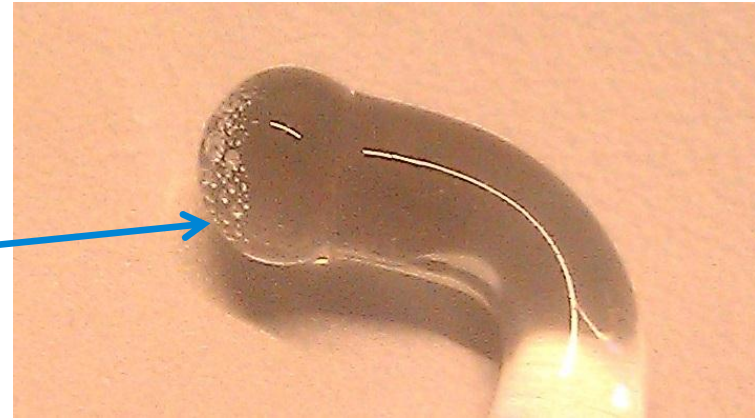
Cleaning the burner:

- 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 centered in front of venturi
 - Distance between bead & venturi should be ~ 0.1 mm (thickness of sheet of paper)

Tighten adjuster to lock bead in place



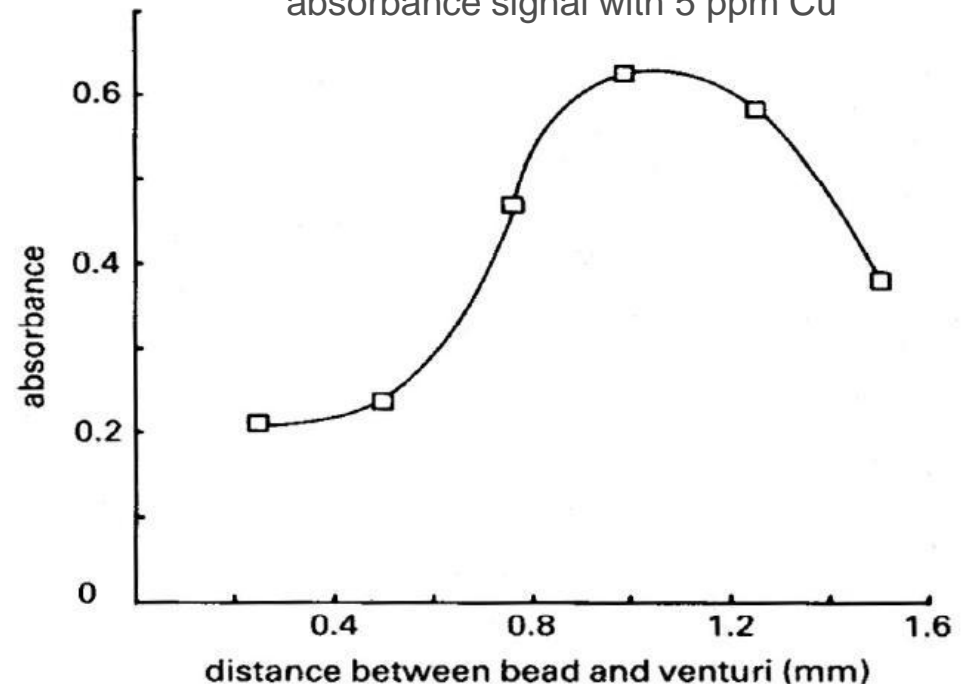
How Impact Bead Setting Impacts Sensitivity?

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 ½ 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:

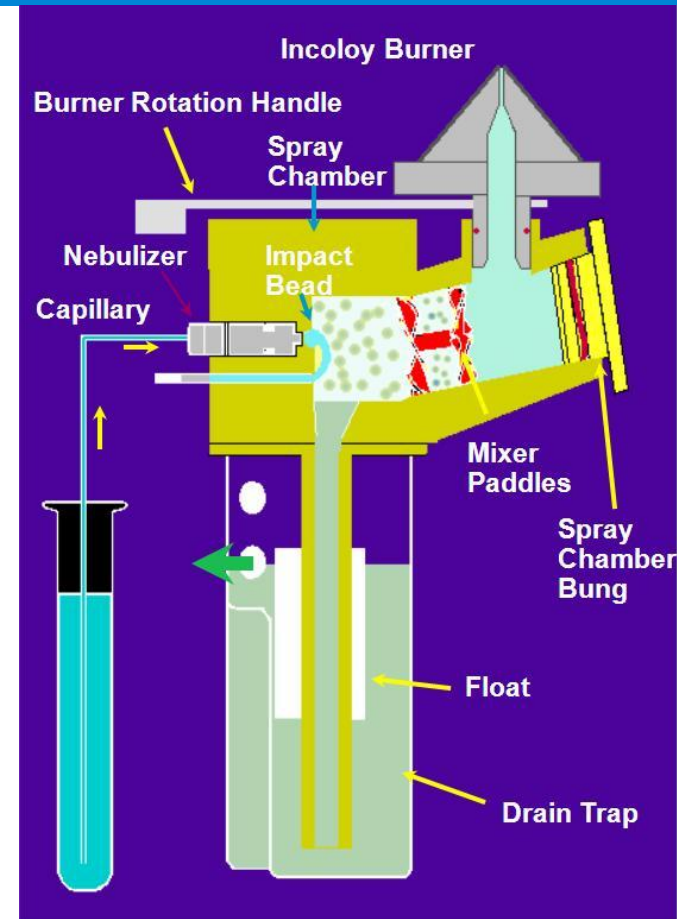
- Fit the mixing paddles
- Optimize the impact bead for best sensitivity
- Use narrow bore capillary tubing

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

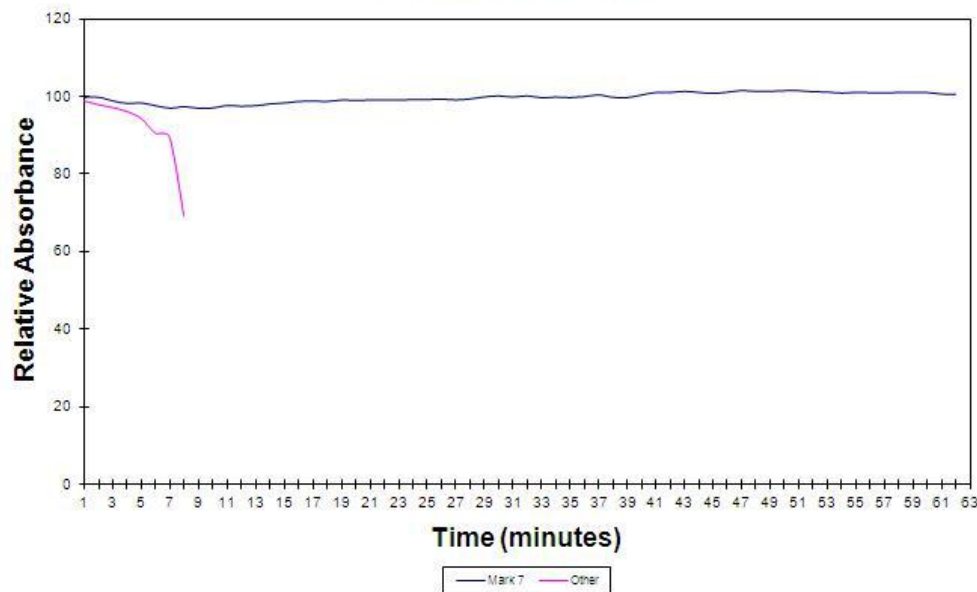
Flexibility, Ease of Use & Superior Flame Performance

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)



100 mg/L Al in Cola



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

C E R T I F I C A T E O F A N A L Y S I S

BASELINE® Nitric Acid

PRODUCT NUMBER: S020101 LOT NUMBER: 1211120 ASSAY (HNO₃, w/w): 68%

1A		2A		3A		4A		5A		6A		7A	
3	LI	4	Bb	5	B								
< 0.5	< 1	< 10		< 10									
11	Na	12	Mg	13	Al								
< 5	< 5	< 10		< 10									
19		20		21		22		23		24		25	
K	Ca	Sc	Ti	V	Cr	Mn	Fe	Co	Ni	Cu	Zn	Ga	Ge
< 1	< 5	< 1	< 5	< 0.5	< 10	< 1	< 10	< 1	< 10	< 1	< 1	< 0.1	< 10
37		38		39		40		41		42		43	
Rb	Sr	Y	Zr	Nb	Mo	Ru	Rh	Pd	Ag	Cd	In	Sn	Sb
< 1	< 1	< 0.1	< 0.5	< 1	< 1	< 1	< 1	< 5	< 0.5	< 0.1	< 0.1	< 5	< 5
55		56		57		58		59		60		61	
Cs	Ba	La	Hf	Ta	W	Re	Pt	Au	Hg	Tl	Pb	Bi	Po
< 0.05	< 1	< 0.05	< 0.05	< 10	< 1	< 0.1	< 1	< 1	< 10	< 0.05	< 0.5	< 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)

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:

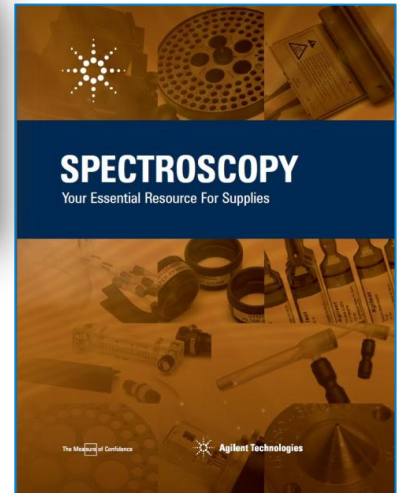
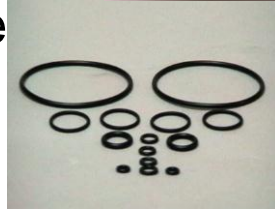
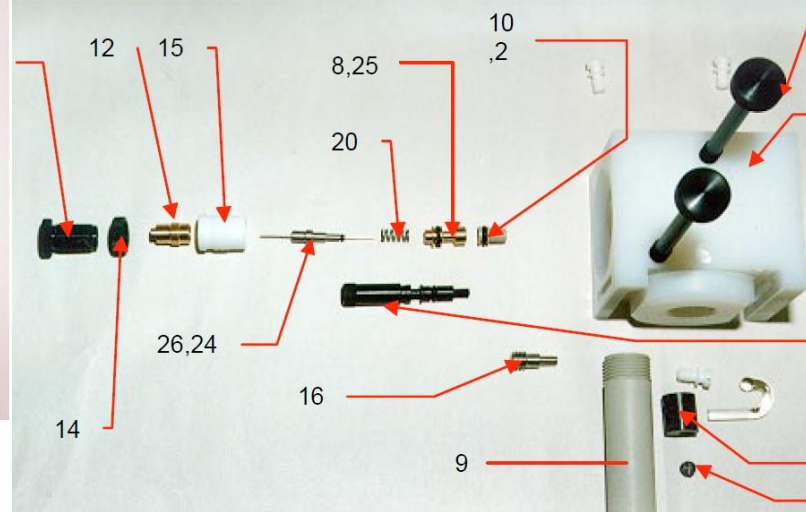
- Quartz atomization cells
- Peristaltic pump tubing
- Connecting tubing



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 30 mm OD tubes (21 positions) 1 pack 30 mm 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 µL 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 1 L 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

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