

# Plasma Ignition Troubleshooting For The 4100 MP-AES

Stephen Anderson, September 2012



# Protecting the torch during ignition

- Always watch plasma ignitions using the recommended position
- Any signs of Torch Overheating use the Emergency Plasma Off switch

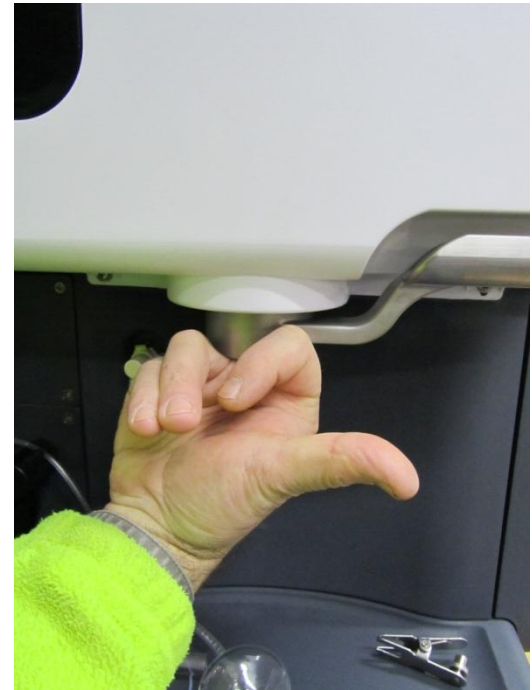


# Diagnosing Gas Flows

- Use your finger to confirm the presence of gas flows all the way to the torch connections
  - Plasma Outer Gas flow
  - Plasma Intermediate flow
  - Argon flow
  - Air Injection flow (if EGCM present)
- **DO NOT USE THIS TO CONFIRM IGNITER OPERATION**

# How to check gas flows to the torch loader

- Remove Plasma Torch.
- Using the Gas Control Diagnostics Tab, enable each gas flow to the Intermediate and Outer Flow, one at a time.
- Insert finger into the Torch Loader Assembly.
- Locate the flow channel (Outer/Upper or Intermediate/Lower) and place tool over the gas channel outlet. There should be distinct sense of PRESSURE, not just flow.



# Plasma Ignition Problems on the 4100 MP-AES

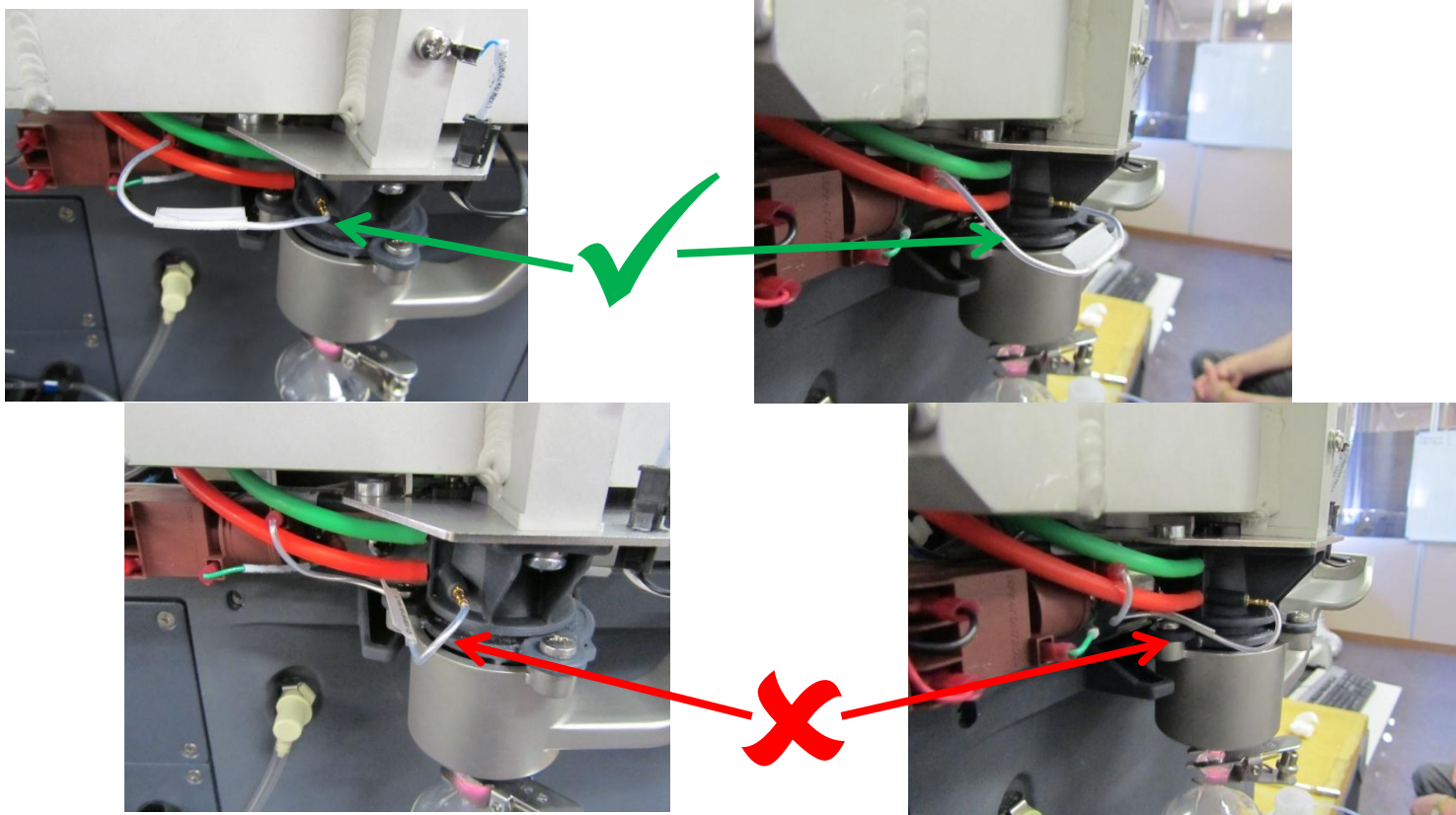
- Plasma Ignition on the 4100 MP-AES Requires
  - Spark
    - Generated by high voltage discharge
  - Argon Gas Flow
    - Required to transfer spark from igniter to the torch region inside the waveguide
    - Ar supplied by either small on board gas cylinder or connected to installation main Ar gas supply
    - Purity not critical (>99%, typically welding grade Ar acceptable)
  - Detection of plasma formation
    - Achieved using an LED located at the
  - Switch to N<sub>2</sub> gas

# Known Causes of Plasma Ignition Problems

- Faulty Plasma Igniter
- Incorrect looming of the HV lead of the Igniter
- Faulty Ar regulator supplied for the on-board Ar gas supply
- Ar leak at the barbed fitting on the Torch Loader assembly
- Faulty Torch Loader plastic moldings
- Contaminants on the soft rubber sections of the MP plasma Torch
- Blocked orifices in the MP torch
- Inadequate insulation of the HV Ignition Lead
- Leaking gas connections on the Main Gas Control System
- Compressed air supplied at greater than specified supply pressure to an EGCM connected to the Main Gas Control system

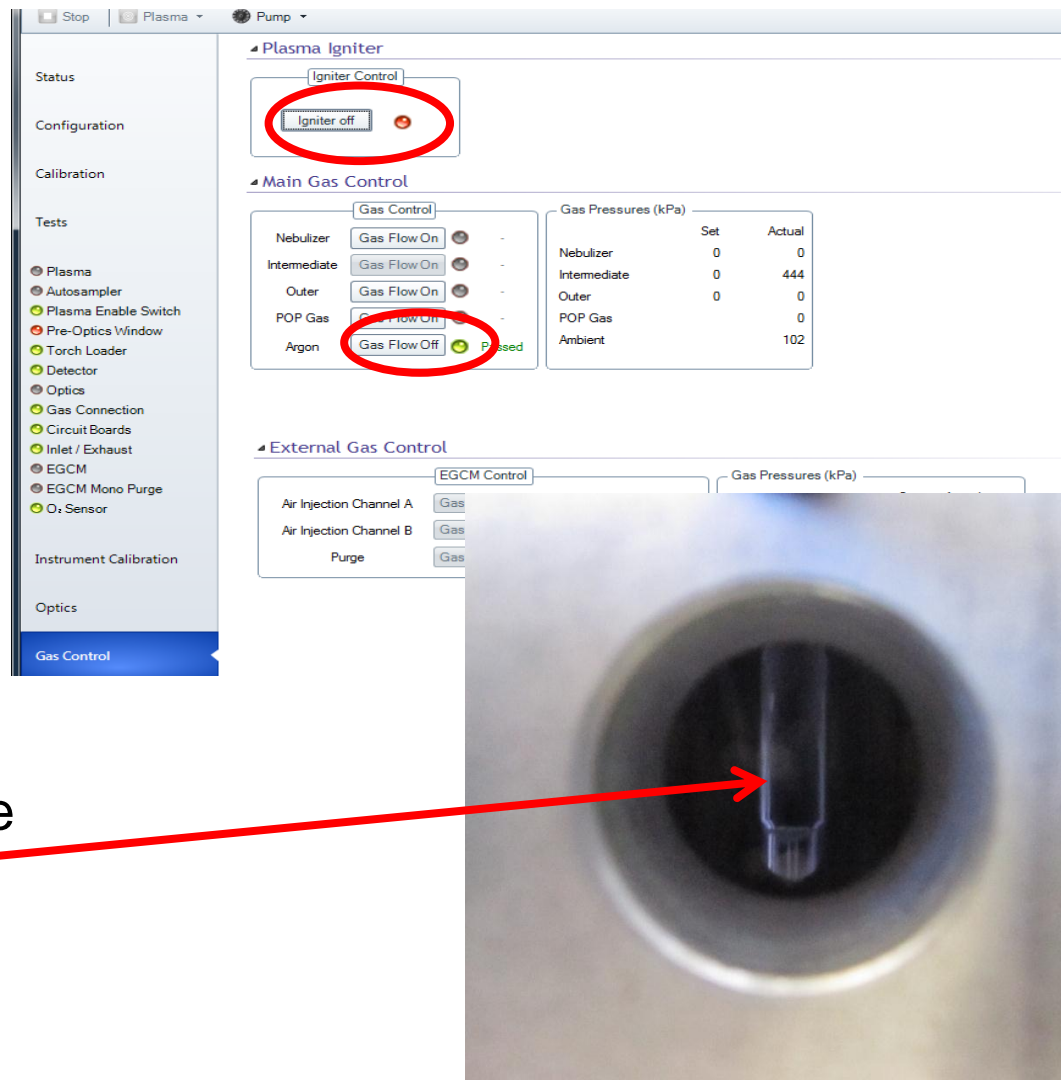
# Igniter HV Loom Location

- If the HV loom of the Igniter is located too close to an earth point the spark may not take the path into the torch
- Easily detected by removing the outer cover enclosing the Waveguide Assembly and inspecting the position of the loom
- Igniter loom can be inadvertently moved during the installation of the Waveguide Cover



# Confirm The Igniter Discharge Is Going To The Torch

- Remove the Waveguide Cover
- In service diagnostics:
  - Simultaneously Enable Ar flow and the Igniter
- Observe the Plasma Torch
  - An electric discharge should be visible inside the plasma torch



# Failure to Detect Igniter Discharge In the Plasma Torch

- Ar gas flow problem
  - Incorrect gas supplied to the Ar input
  - Confirm Ar flow using your finger to confirm flow is present to the Torch Loader assembly
- Faulty Igniter
  - Igniter Module
  - HV Igniter cable too close to a ground point
- Torch Problem
  - Check the plasma torch for blockages

# Faulty On-board Ar Gas Regulator

- A number of possible fault states
  - With the Ar regulator firmly attached to the on-board Ar cylinder and Swagelock connection in place, partially and slowly release the Swagelock connection at the Main Gas Control system and confirm that Ar is flowing.
  - In Service Diagnostics, remove the plasma torch and enable Ar flow and confirm pressure at the Intermediate output port of the torch loader assembly
  - In Service Diagnostics, install the torch and enable Ar flow and Igniter, no sparks should be visible in the vicinity of the torch loader assembly.

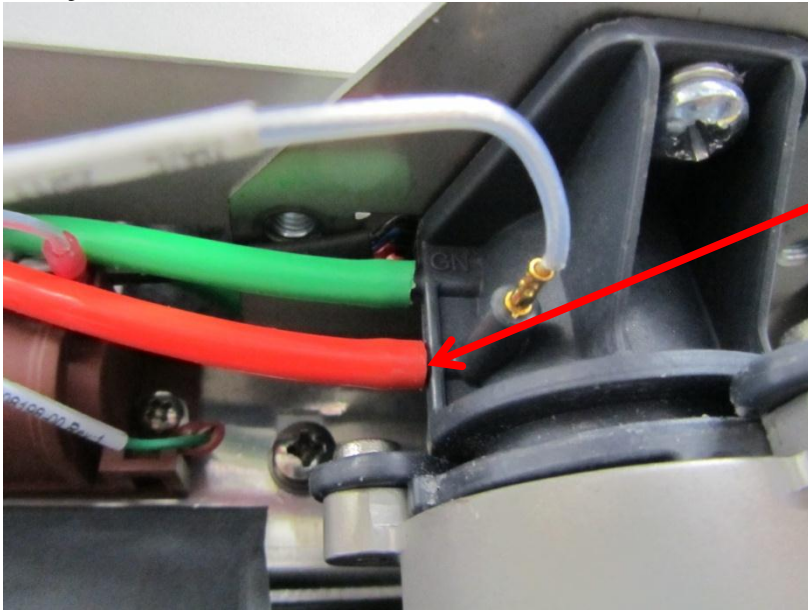


# Reduced Argon Purity Through Diffusion of Atmospheric Gases

- Typical Indication
  - Difficulty igniting after plasma has been off for an extended time ( > 1-2 hrs)
  - After successful plasma ignition plasma can be extinguished will reliably ignite if ignited immediately after being extinguished
  - Plasma reliably ignites if the Ar gas supply is purged immediately prior to plasma ignition (using diagnostics)
- Typical Problem
  - The use of Polymer hoses for Ar supply
  - Allows ingress of atmospheric gases via diffusion
  - If instrument is off for an extended time atmospheric gases can permeate through the Polymer Hoses and reduce Ar purity
  - Long Polymer supply hose are particularly prone to this problem

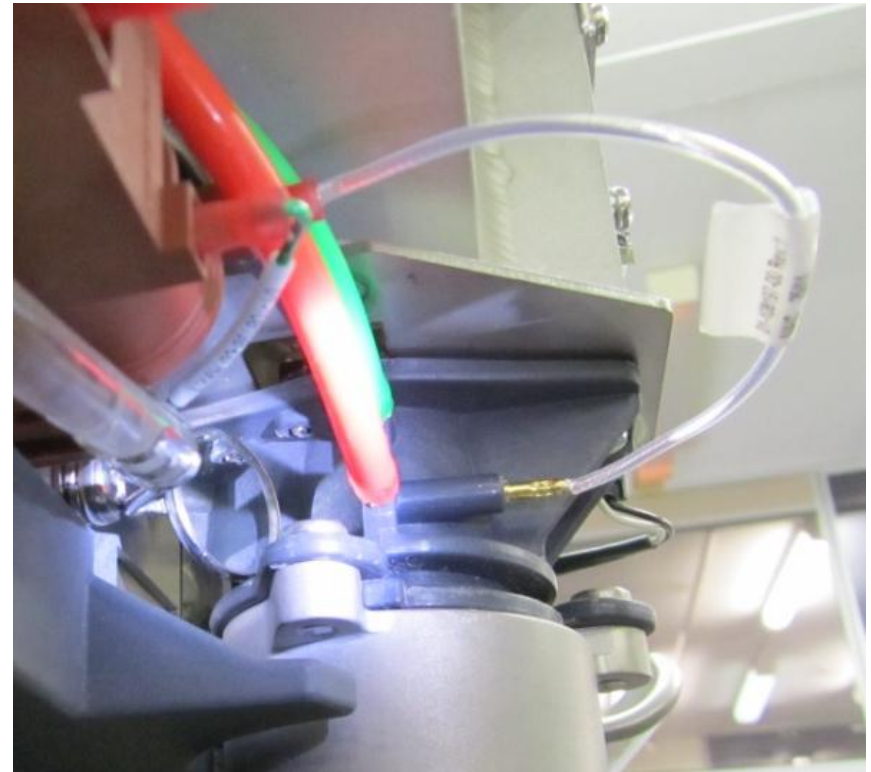
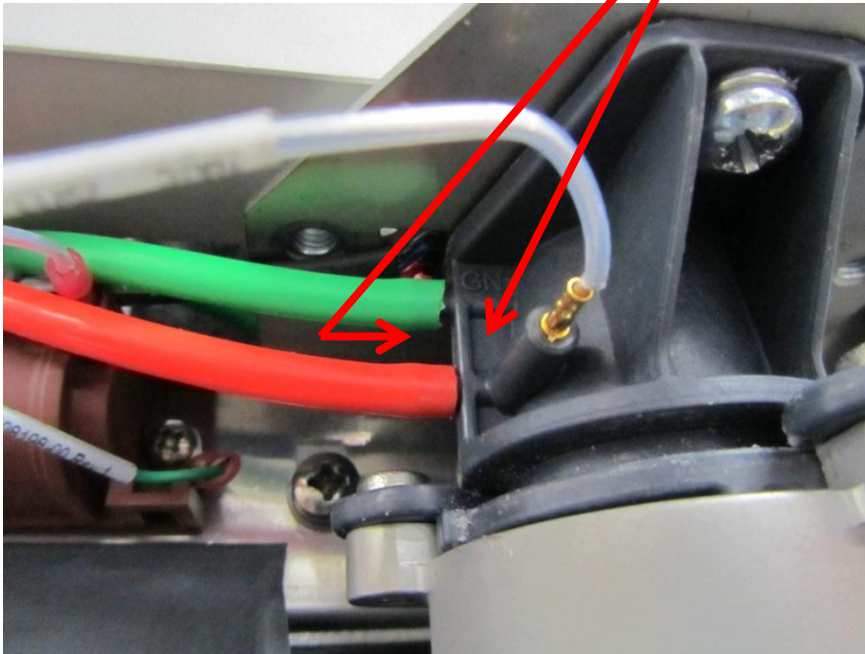
# Ar leak at the barbed fitting on the Torch Loader assembly

- Typically results in HV arcing to earth points around the Torch Loader
- If a leak is suspected, carefully remove the red hose (intermediate flow) from the barbed fitting, cut off the end of the hose and reattach
- It is recommended to immerse the newly trimmed end of the hose in very hot water to soften before reattaching.



# Problems with the Torch Loader Molding Assembly

- By removing the Torch Loader Assembly from the Waveguide Assembly or by using a mirror and torch, inspect the Torch Loader molding for holes.
- Inspect both front and back sides



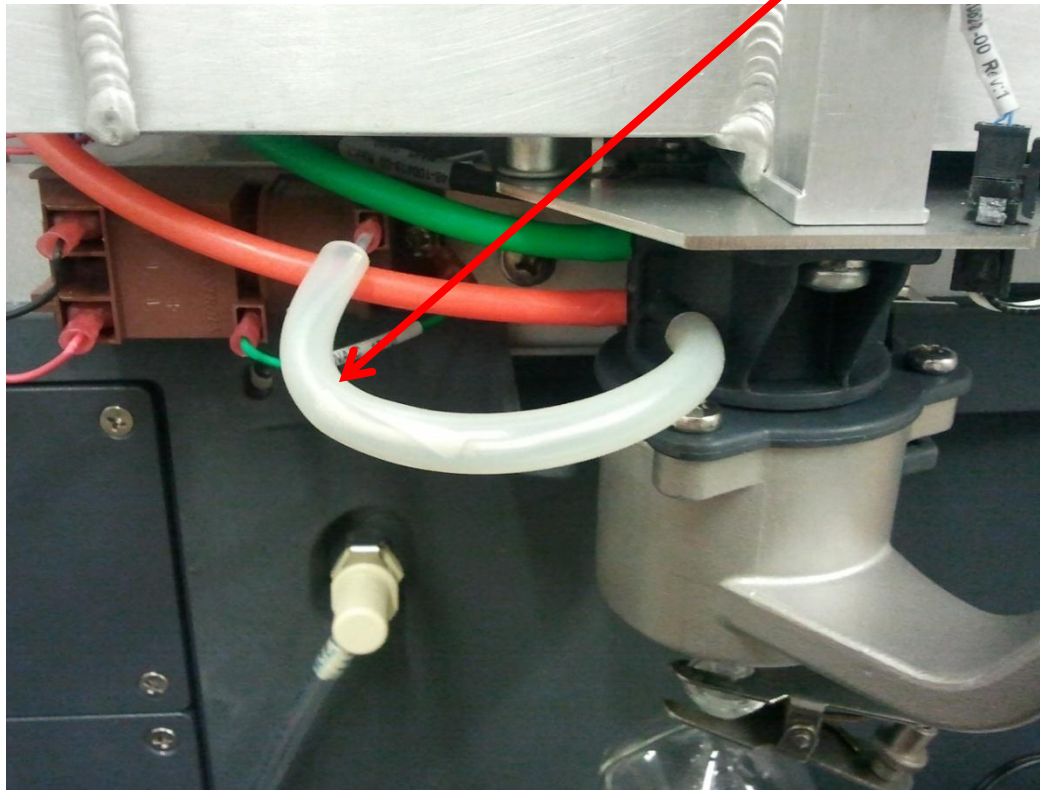
# Particulates on the Gas Sealing Surfaces or Blocked Orifices of the MP Plasma Torch

- Checked for blocked orifices
  - Visual Inspection
  - Compressed Gas Flow
- Check for Particulates on the Gas sealing surfaces
- Consider replacing with a New torch



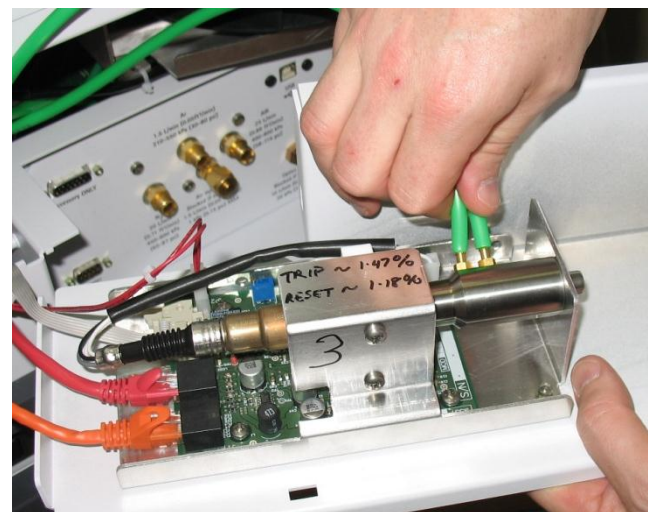
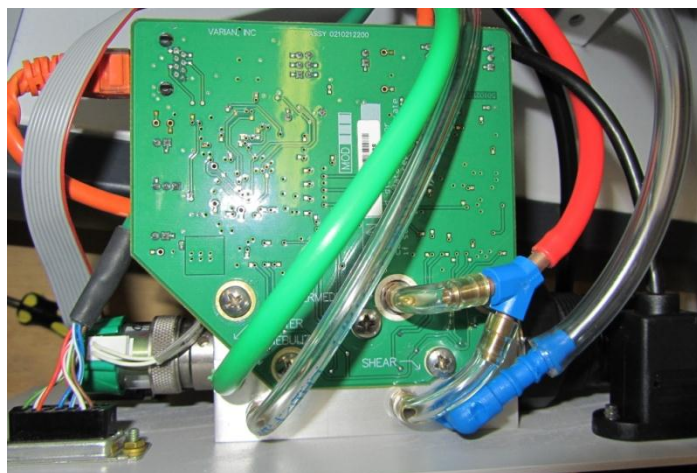
# Improved Insulation on the HV Igniter Lead

- Ignition problems have also be resolved by improving the insulation on the HV igniter lead
  - Installation of Silicone hosing over the HV lead



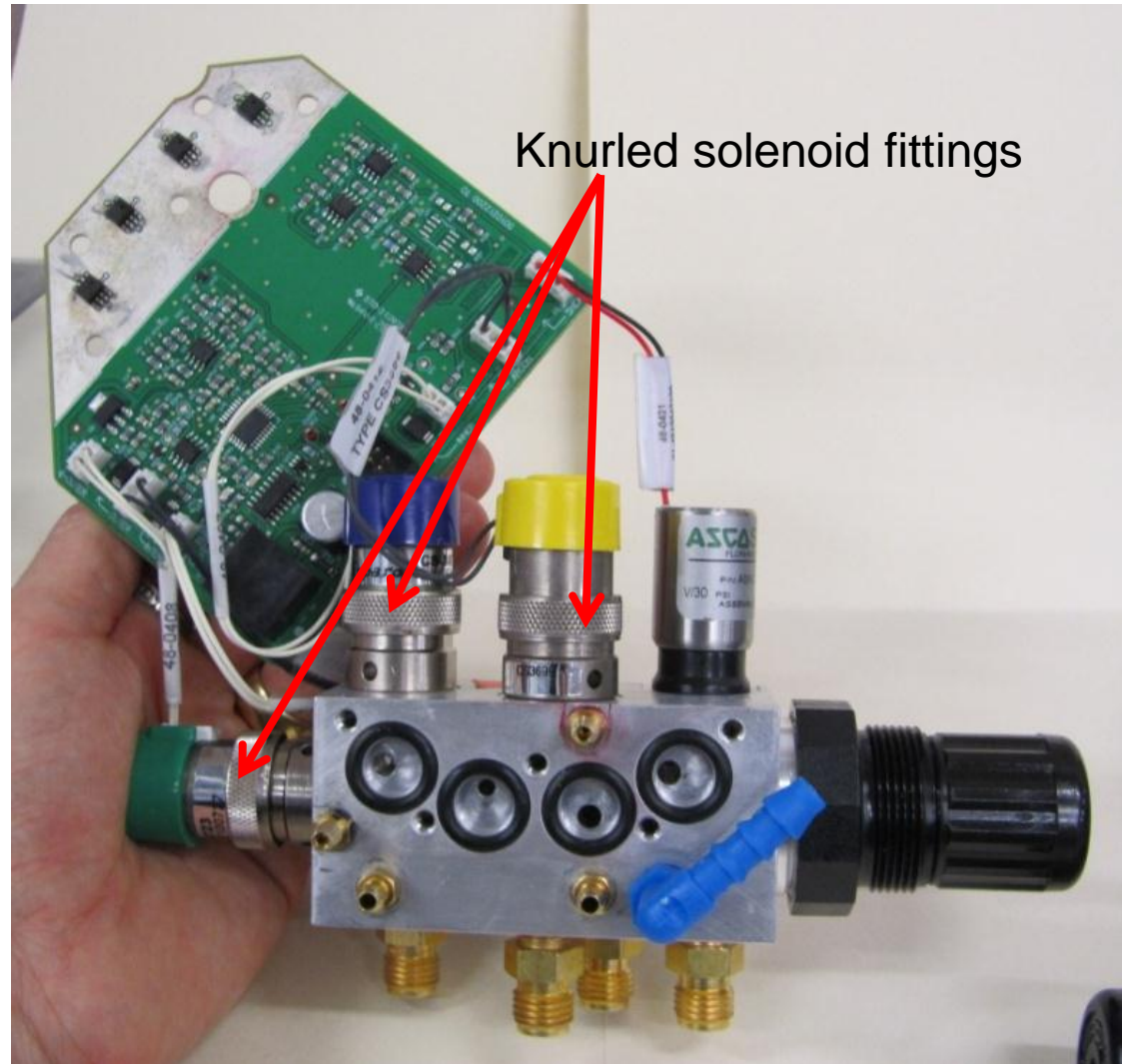
# Leaking Gas Connections on the Main Gas Control System and Oxygen Sensor Connections

- Confirm gas connections to the Main Gas Control System are firm
- Confirm gas connections to the Oxygen Sensor are firm
- Remove any suspect connection, trim hose ends and refit
  - Consider immersing the cut, trimmed hoses in hot water before reconnecting



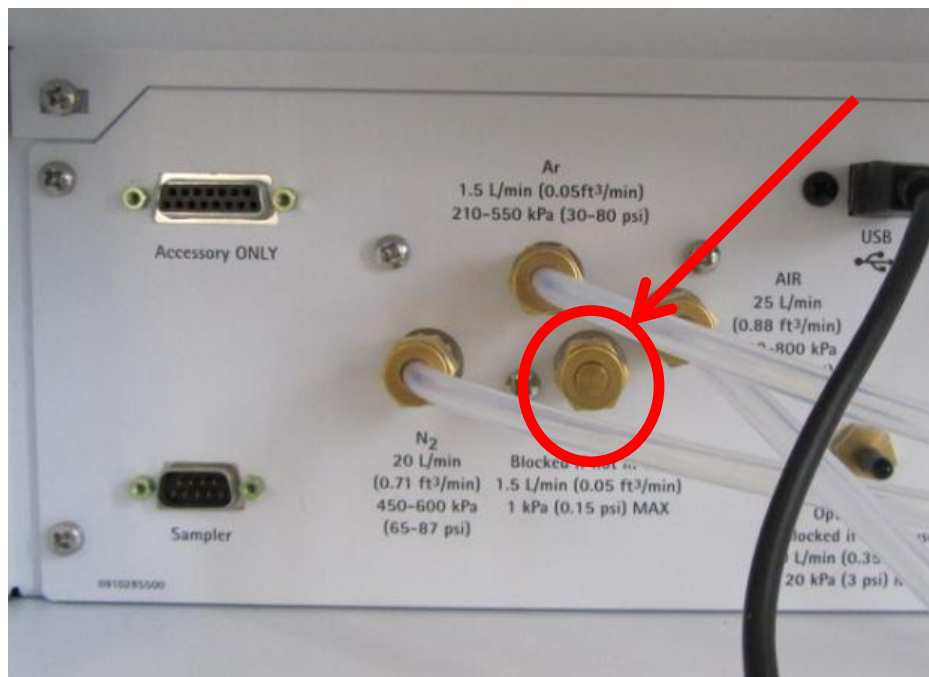
# Gas Leaks Due to Loose Solenoid Valve Assembly

- Confirm the knurled fittings holding the solenoid valve assembly together is NOT loose
- Confirm the Valve is NOT loosely attached to the Manifold



# Air Injection Port

- If not in use ensure the Air Injection port is capped
  - If left uncapped, known to cause
  - Performance problems
  - Plasma Ignition problems
  - Torch melt problems
- The correct Swagelok fitting should be used to ensure proper sealing



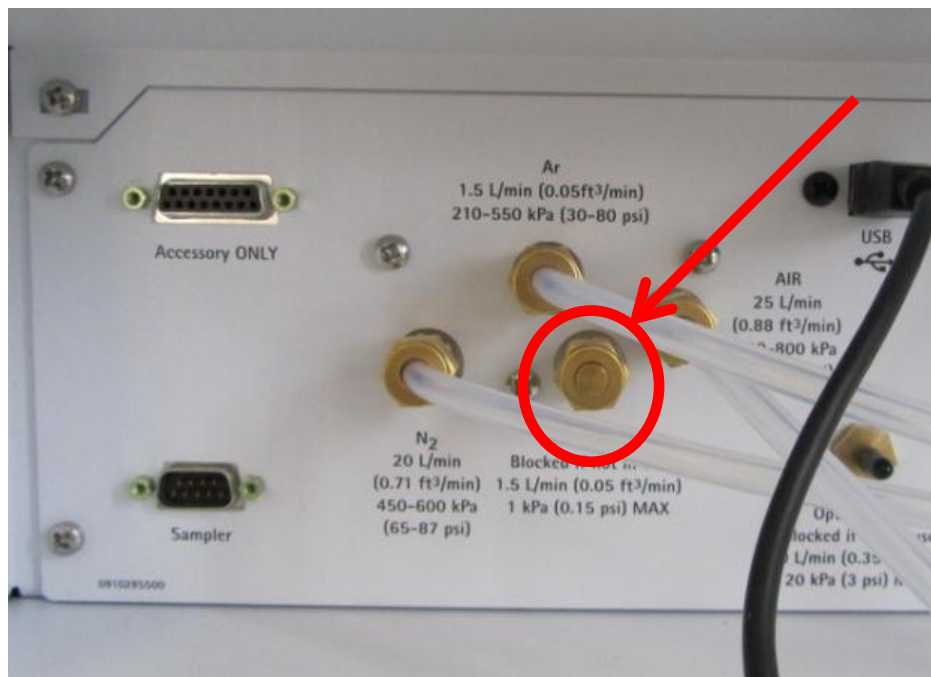
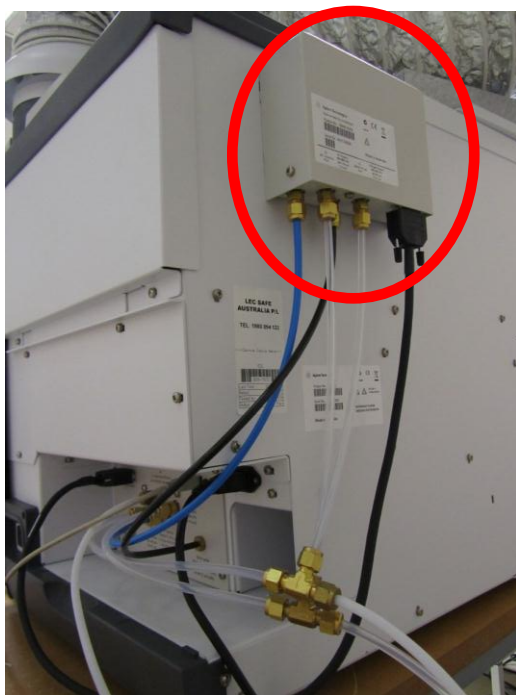
# Air Injection via the EGCM

- If an EGCM is connected to the Air Injection port of the Main Gas Control system
  - Confirm the supplied air pressure to the EGCM is within recommended limits
  - Excessive air pressure to the EGCM can cause the solenoid on the EGCM to leak, injecting air into the plasma continuously



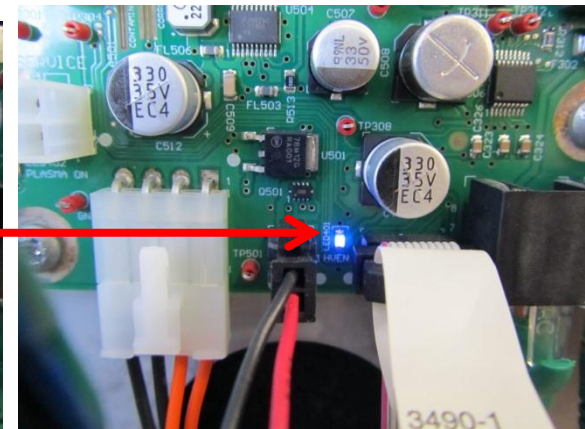
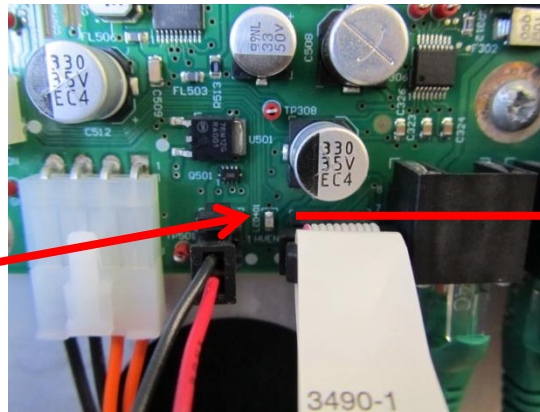
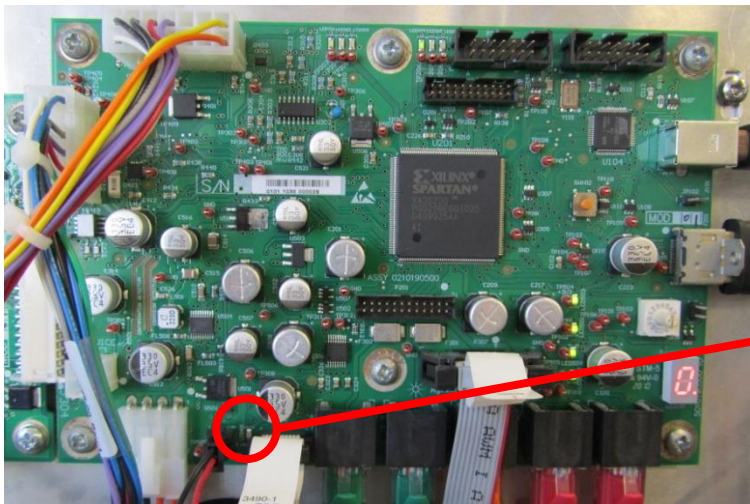
# Accessories

- If an EGCM is attached and the Air injection connected to the Instrument Services Panel disconnect/disable
  - Disconnection and removal ensures no possible interference of the accessory with the operation of the instrument
  - Ensure the air injection port is capped if not connected



# Main Control Board “Blue” LED401, HVEN (High Voltage Enabled)

- If all interlock conditions are met, (HVEN) should turn on during the gas purge process and stay on until the plasma is extinguished
- Failure of LED401 to turn on and stay on, can indicate a Main Control Board Problem
  - If LED401 does not remain ON and there is no indication of error state that will prevent the plasma ignition replace the main control board



# Check Electrical Connection to the HVPS

- Turn the off the 4100 MP-AES
- Disconnect and reconnect all electrical looms to the HVPS
  - High Voltage Loom to Magnetron
  - Communication and Interlocks
  - Mains Voltage and Fan Mains Control

