

Nafion Dryer

Installation Manual

Version 1.0

March 2008

1.	Installation on an Air Server series 1.....	2
1.1.	Standard Air Server series 1	2
1.2.	Inverted Air Server series 1	4
2.	Installation on CIA 8	7
3.	Installation on an Air Server series 2.....	8

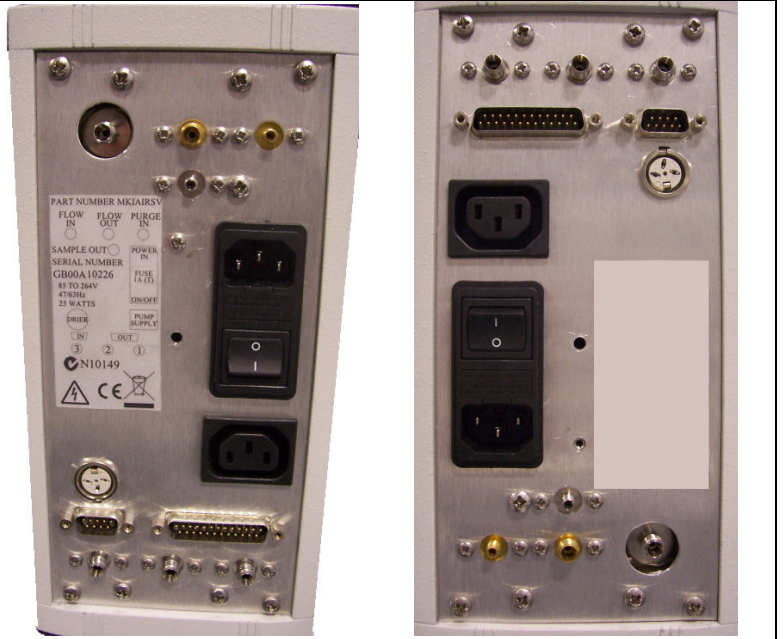
1. Installation on an Air Server series 1

Switch off power to the Air Server.

Disconnect the link tube if connected.

If the rear panel looks like the picture on the left follow the installation instructions in Section 1.1.

If the rear panel looks like the picture on the right then refer to Section 1.2

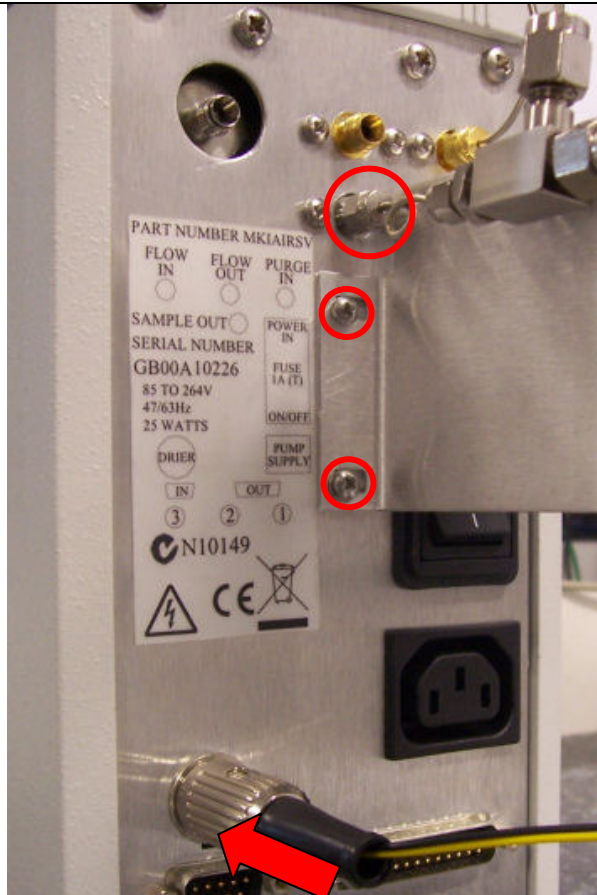


1.1. Standard Air Server series 1

Mount the nafion dryer assembly using the 2 screws highlighted.

Connect the 1/16" stainless steel (SS) fitting highlighted and tighten using an 8 mm spanner.

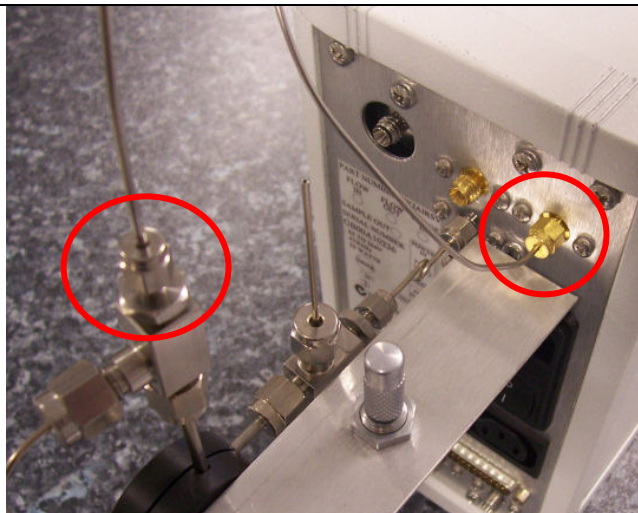
Plug the solenoid connector into the back panel.



The brass 1/16" connection from the link tube connects to the brass 1/16" union on the back panel.

The stainless steel 1/16" connection connects to the output on the nafion dryer as shown.

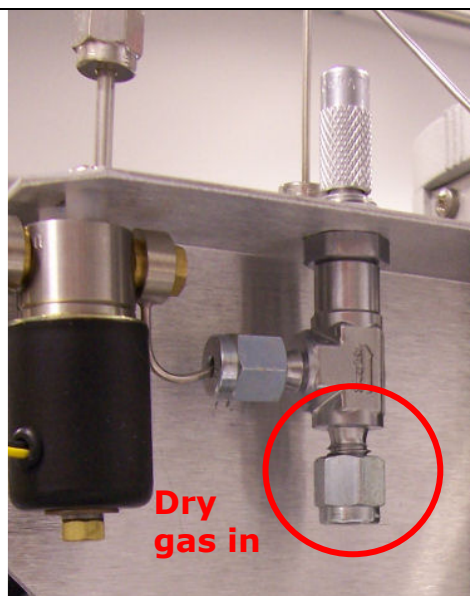
Caution: Do not rotate the fitting on the nafion dryer, use a 7/16" spanner to secure the fitting while tightening.



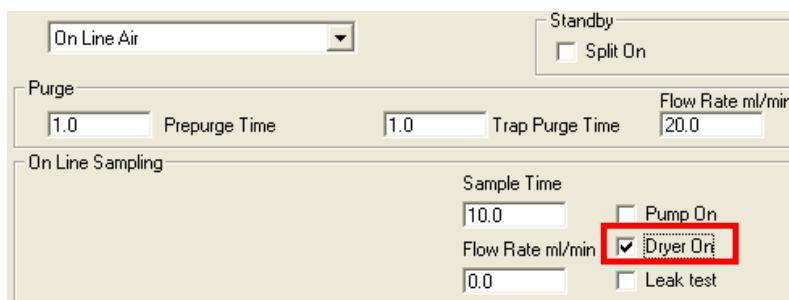
Dry gas (typically nitrogen) is connected to the 1/8" fitting on the nafion dryer as shown.

Note: The supply pressure should be below 50 psi.

The flow is controlled by the needle valve, once activated in the software.

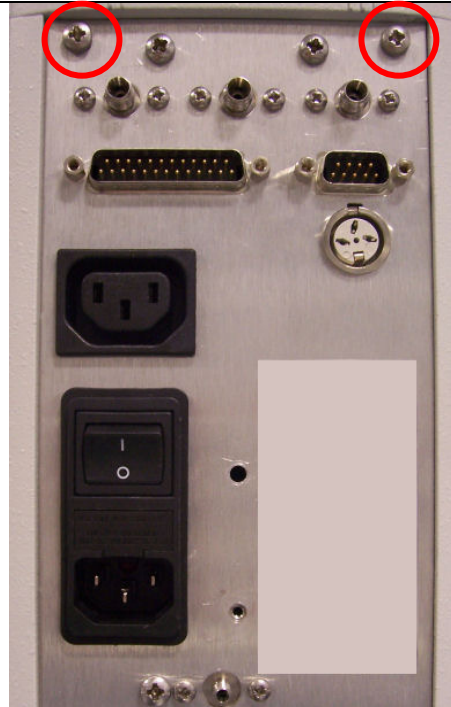


The dryer can now be selected in an 'On Line Air' method.

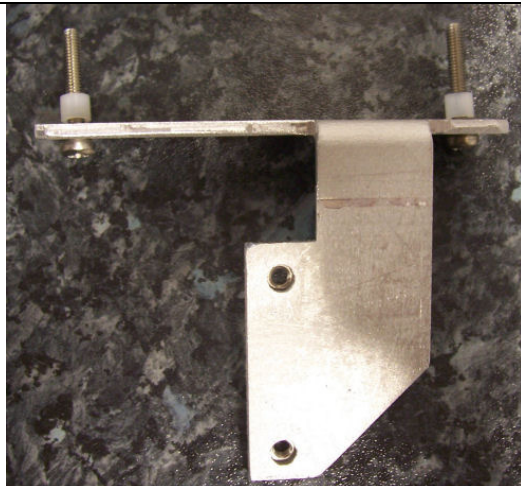


1.2. Inverted Air Server series 1

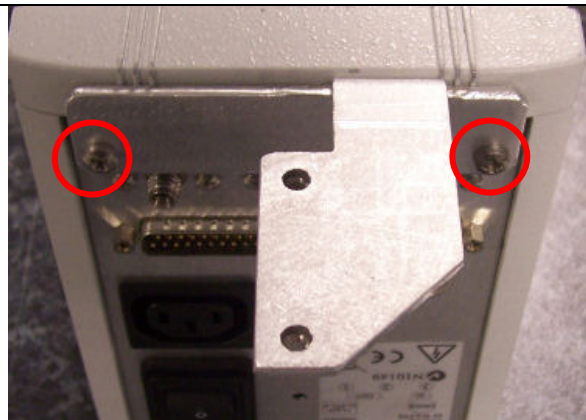
Remove the 2 screws from the top corners of the back panel.



Locate the M4 x 25 screws, spacers, washers and bracket and assemble as shown.



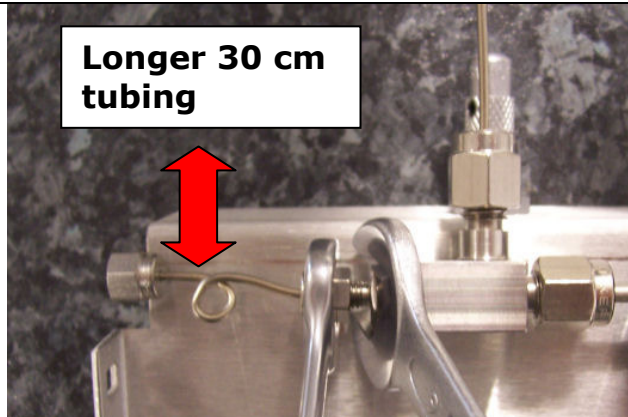
Mount the assembled bracket to the 2 holes at the top corners of the back panel.



Remove the short 1/16" SS tubing using 8 mm and 7/16" spanners to avoid rotating the fitting.

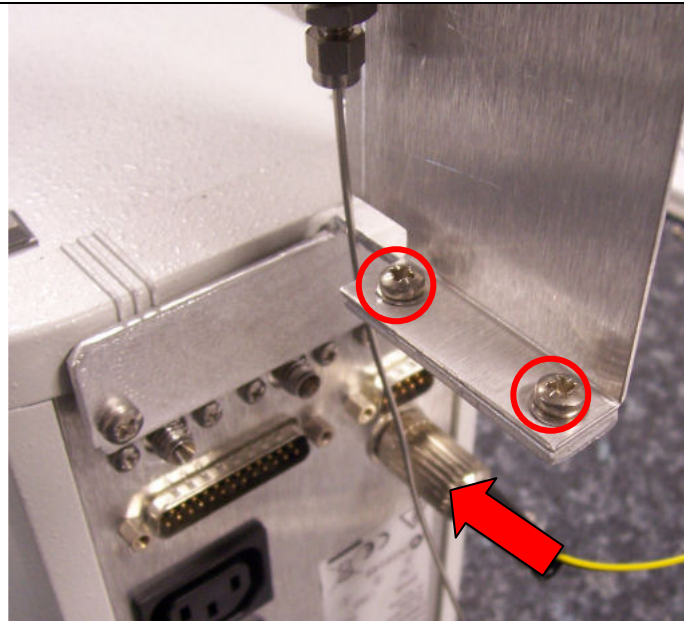
Attach the nuts and ferrules to the 30 cm length of SS tubing supplied. Fit the 1/16" nut first then the back ferrule and front ferrule.

Attach the SS tubing to the nafion dryer using 2 spanners to avoid rotating the fitting.



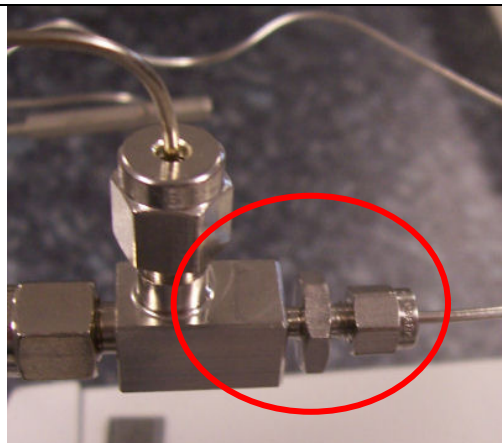
Attach the nafion dryer assembly to the mounting bracket using M4 x 12 screws and washers.

Plug the solenoid connection into the back panel.



Attach the 1/16" SS connection on the link tube to the output of the nafion dryer as shown.

Caution: Do not rotate the fitting on the nafion dryer, use a 7/16" spanner to secure the fitting while tightening.



Connect the 30 cm SS line to the SS 1/16" connector on the back panel.

Connect the brass connector on the link tube to the 1/16" brass connector on the back panel.



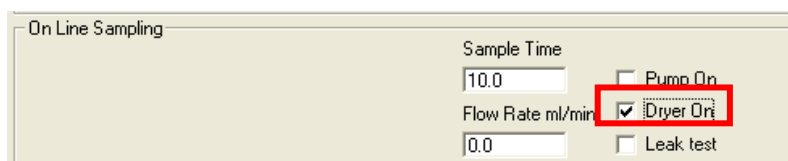
Dry gas (typically nitrogen) is connected to the 1/8" fitting on the nafion dryer as shown.

Note: The supply pressure should be kept below 50 psi.

The flow is controlled by the needle valve, once activated in the software.



The dryer can now be selected in an 'On Line Air' method.

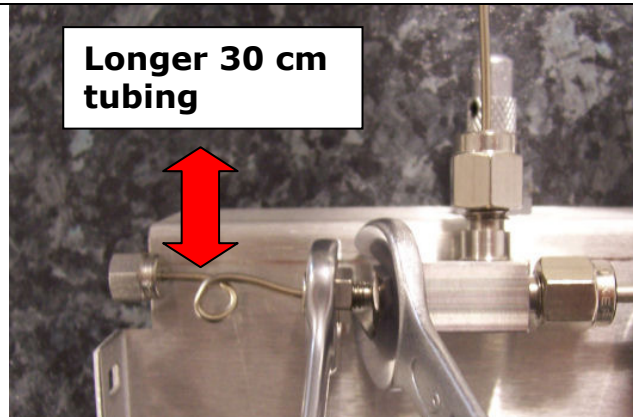


2. Installation on CIA 8

Remove the short 1/16" SS tubing from the nafion dryer assembly using 8 mm and 7/16" spanners to avoid rotating the fitting.

Attach the nuts and ferrules to the 30 cm length of SS tubing supplied. Fit the 1/16" nut first then the back ferrule and front ferrule.

Attach the SS tubing to the nafion dryer using 2 spanners to avoid rotating the fitting.



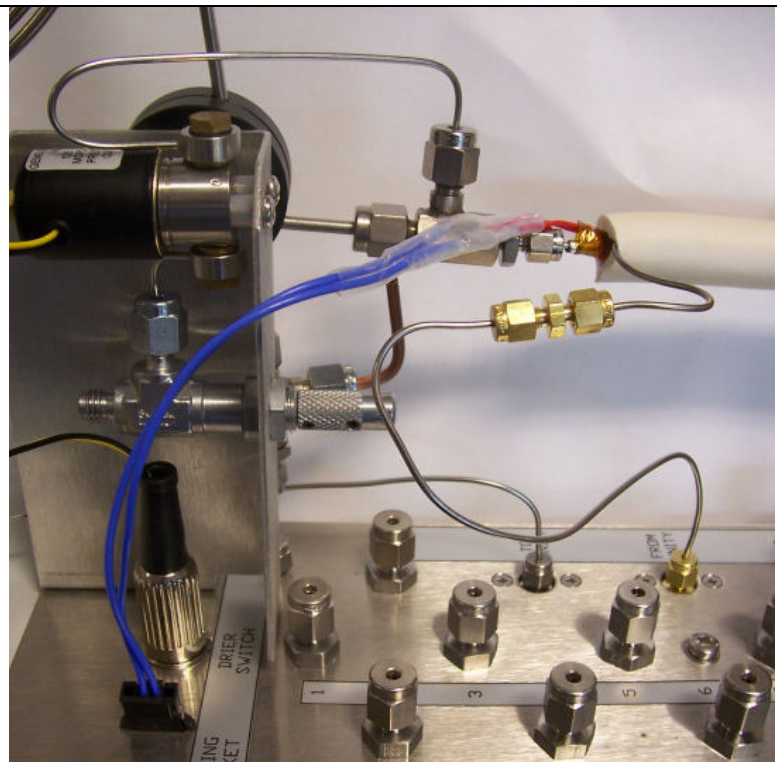
The dryer mounts with a single screw to the 'DRIER MOUNT' point.

The 30 cm SS tubing connects to the 'TO UNITY' union.

The SS fitting on the heated transfer link connects to the nafion output.

The brass fitting on the heated transfer link connects to the 'FROM UNITY' union via a 1/16" union and 20 cm length of SS tubing.

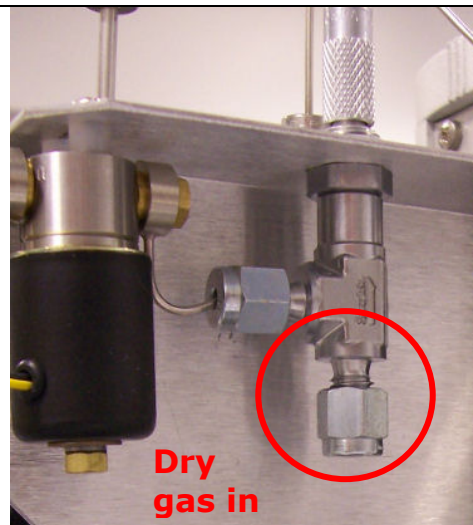
Plug the solenoid connection into the 'DRIER SWITCH' position.



Dry gas (typically nitrogen) is connected to the 1/8" fitting on the nafion dryer as shown.

Note: The supply pressure should be kept below 50 psi.

The flow is controlled by the needle valve, once activated in the software.



The dryer can now be selected in an 'On Line Air' method.

The screenshot shows a software interface for 'On Line Sampling'. It contains the following elements:

- Sample Time:** A text input field containing the value '10.0'.
- Flow Rate ml/min:** A text input field containing the value '0.0'.
- Control Options:** Three checkboxes are listed:
 - Pump On
 - Dryer On (This checkbox is highlighted with a red rectangular box)
 - Leak test

3. Installation on an Air Server series 2

Please refer to the Air Server series 2 installation manual