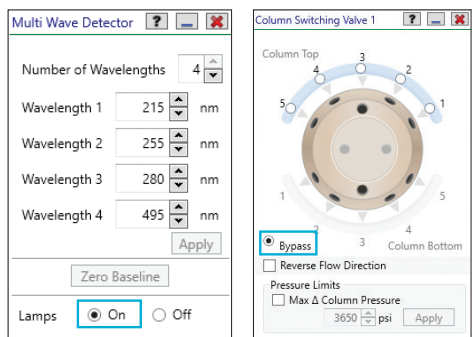


NGC Liquid Chromatography System Quick Guide

Starting Your NGC Chromatography System

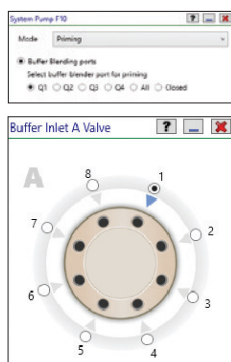
- 1 Turn on the system and detector to allow lamps to warm up.**
Note: If the system is in a cold room or box, the system should remain powered on to prevent condensation from forming inside.
- 2 Ensure all column switching valves are set to Bypass.**



Priming the Pumps

All lines should be primed with filtered and degassed water using a syringe at the pump head priming port.

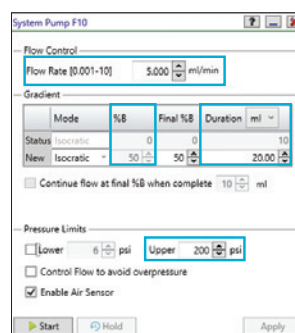
- 1 Insert a syringe onto the priming port, rotate the priming port one half turn, and use the syringe to pull 15 ml.**



- If using a buffer blending valve, change the mode to **Priming** and select the buffer blending port you want to prime
- If using an inlet valve, select the port you want to prime

Purging the System

- 1 Set the pumps to flush 20 ml of filtered and degassed water through the system.**



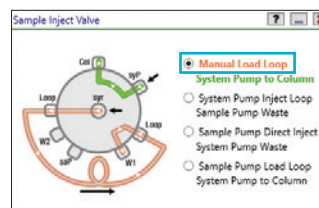
- Set Flow Rate
- Set %B
- Set Duration
- Set Pressure Limits

- 2 If a column is stored in 20% ethanol, be sure to flush the column with water.**
 - If the system contains a column switching valve, select the correct column position on the column switching valve
 - In the pump control window, enter the recommended flow rate, adequate duration to flush the column, and the upper pressure limit of the column in line
 - LED lights will indicate flow to the column position

- 3 Repeat priming steps using the appropriate filtered and degassed buffer for the run.**

Sample Application Using the Sample Loop

- 1 Load sample in Manual Load Loop mode.**



- Flush the loop with water and then equilibration buffer before introducing the sample
- Load a minimum of 200 µl greater than the loop volume to ensure air has been purged from the loop

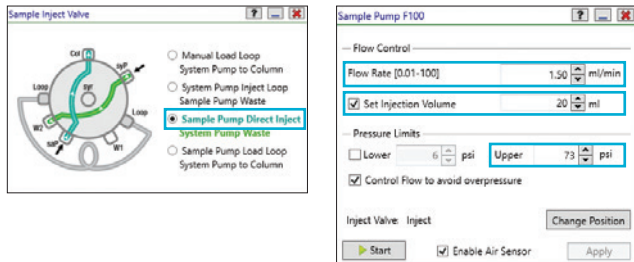
2 Select System Pump Inject Loop on the Sample Inject Valve to start sample injection with the system pumps to the column.

- Use 2.5x the loop volume to ensure all sample has been loaded to the column

Sample Application Using the Sample Pump

1 Prime the sample pump with water and then equilibration buffer, followed by sample.

2 Select Sample Pump Direct Inject on the Sample Inject Valve.



- Enter the flow rate
- Set the Injection Volume. If equipped with sample air sensing, enable air sensor and set the injection volume to greater than the sample volume
- Set the Column Pressure Limits
- Start the pump

Cleaning the System

1 After priming the pumps and ensuring the column switching valves are in bypass, flush the system with 30 ml water followed by 30 ml 0.5 M NaOH, incubate for 30 minutes to an hour, and then flush the system with 30 ml of water. Collect all 90 ml using the fraction collector to ensure the dispenser is cleaned.

- To clean the sample loop, set the injection valve to **System Pump Inject Loop**

2 If the sample pump was used, repeat the procedure in step 1 using the Sample Pump while the Sample Inject Valve is in the Sample Pump Waste position.

System Shutdown

- Turn off lamps and reset column switching valve to Bypass.
- Prime the pumps and purge the system with water to remove high salt buffers.
- For storage, prime the pumps and purge the system with 20% ethanol.

Maintenance

- Change pump wash solution (20% ethanol) weekly.
- Power cycle every 2 weeks.
 - Go to **Shutdown** on the touch screen, shut the system down, wait 10 seconds, and turn it back on
 - If your system is in a cold room or box, do not leave the system off for more than 10 minutes
- Clean system after each use to prevent damage to flow cell. For more information about system maintenance, see Appendix A of the NGC Chromatography Systems and ChromLab Software Instrument Guide (1000049091).

Need More Help?

Tutorial videos are in ChromLab Software.

- For priming pumps, go to **Help > Tutorials > Priming the Pumps**
- For priming the system and sample application, go to **Help > Tutorials > Performing a Manual Run**

Visit bio-rad.com/NGC for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

