

Installing the NGC Fraction Collector

Introduction

The NGC Fraction Collector provides versatile options for preparative and analytical chromatography applications. The high-capacity and flexible rack options enable fractions ranging in size from milliliters to liters to be collected and accessed freely during a purification run.

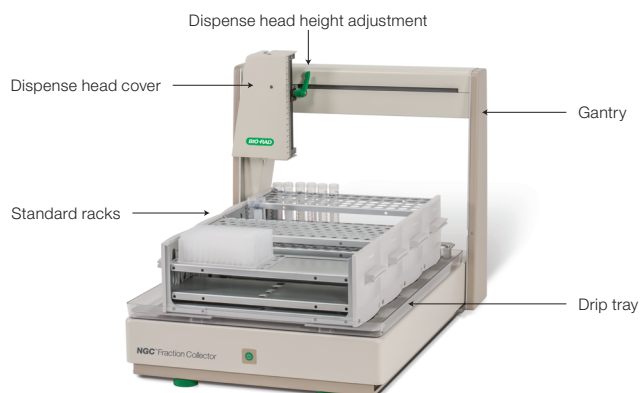
Setting Up the NGC Fraction Collector

This quick guide outlines the process of installing and plumbing the NGC Fraction Collector to your NGC System. The fraction collector (FC) can be installed on a laboratory bench or in a coldroom or chromatography refrigerator.

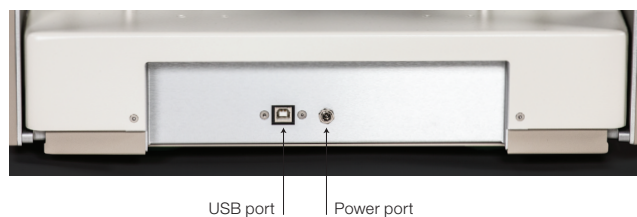
- 1 Open the shipping box and remove the foam packaging. Carefully lift the FC from the box. **Do not** lift the FC by the gantry; reach down and lift it from the base. Grip the base of the FC by the sides to lift from the shipping box.

Check the box for the tubing kit, external power supply, and USB cable. The tubing kit contains:

- 92 cm green polyetheretherketone (PEEK) tubing from NGC union to dispense head for F100 systems (fittings on both ends)
- 92 cm orange PEEK tubing for F10 systems (fittings on both ends)
- 92 cm clear fluorinated ethylene propylene (FEP) waste line from dispense head to union (fittings on both ends)
- 150 cm clear FEP waste line from union to waste bottle (fitting on one end)
- Black magnetic union to be attached to side of NGC System to act as anchor point for FC tubing; position as necessary to avoid kinking any of the tubes

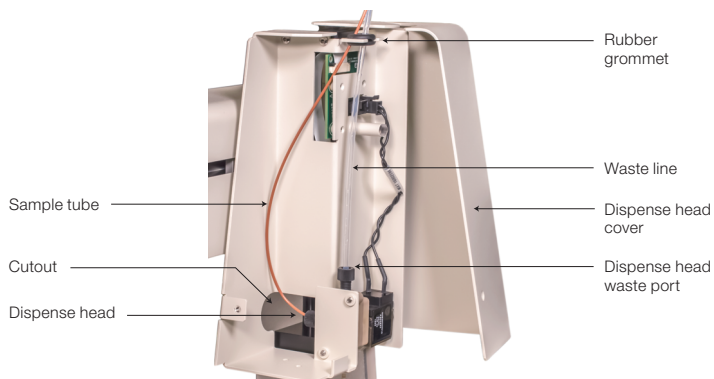


- 2 Connect the power cable and USB to the back of the FC. Plug in the power supply and connect the USB cable into a USB port on the NGC Chromatography System.



Plumbing the NGC Fraction Collector

Use the supplied tubing and union to plumb from the NGC System to the dispense head on the fraction collector. Plumb the waste line from the dispense head to the waste bottle via the union.



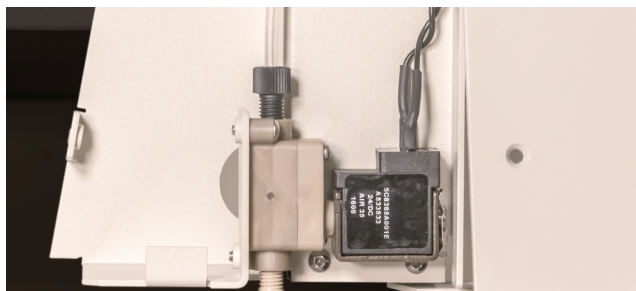
- 1 Open the dispense head cover using a small Phillips screwdriver.



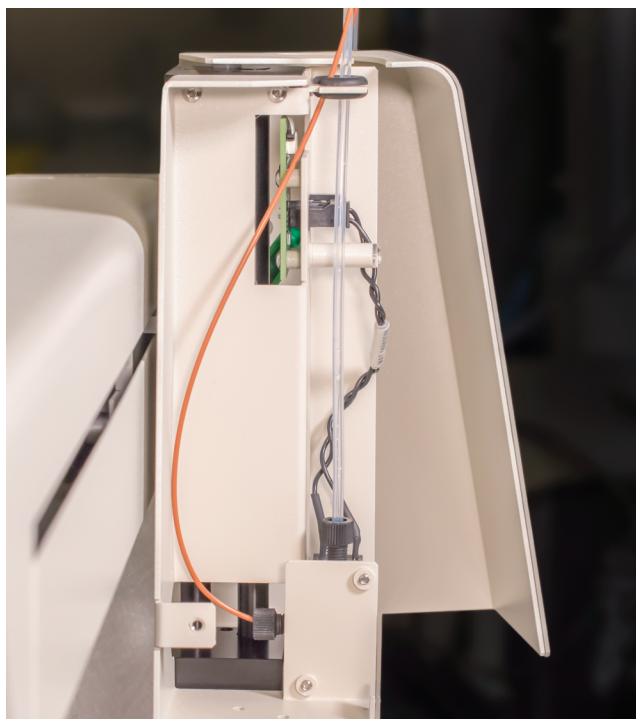
- 2 Connect the sample tube to the dispense head. Use the cutout in the back of the cover to insert a fittings tightener and tighten.



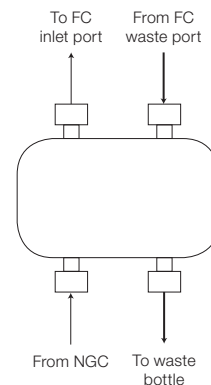
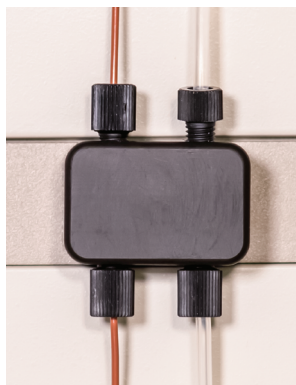
- 3 Attach the waste line from the dispense head waste port to the union on the side of the NGC System.



- 4 Route both tubes through the top of the dispense head cover by sliding them into the rubber grommet. Finally, close the dispense head cover and fasten with the screw.



- 5 Attach the tubes to the union so the direction of flow is as shown below.



Completing NGC Fraction Collector Installation

After plumbing the NGC Fraction Collector, complete the installation and adjustments.

- 1 Place the plastic drip tray on the FC with the notched corner facing the left rear of the FC platform.



- 2 Place the metal rack tray into position, ensuring it is aligned correctly on the locating pins.



- 3 Turn on the fraction collector using the power button on the front.

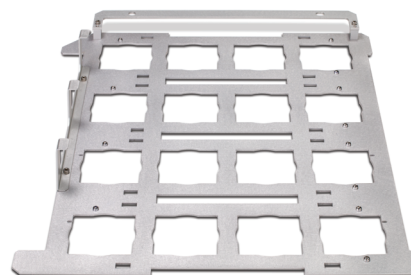


- 4 The gantry and dispense head will move to the home position when the NGC is powered up and the fraction collector is detected by the system.

Note: The gantry will move quickly from the home position to the first collection tube at the start of a run. Ensure there are no objects in the way before starting collection.



- 5 Place the racks on the FC, ensuring they are correctly aligned on the raised locating pins. Identifier labels on the racks should face right.



- 6 Before using the fraction collector, adjust the z-height of the dispense head to sit within 10 mm of the collection vessels.



Ordering Information

Catalog #	Description
12003748	Peltier Rack Power Supply , 4-channel power supply for up to 4 Peltier Rack Modules
12003749	Peltier Rack Module , 1 insulated rack per module
12003759	Insulated Rack for 50 ml Tubes , for 14 x 50 ml tubes
12003751	Insulated Rack for 15 ml Tubes , for 30 x 15 ml tubes
12003747	Peltier Rack Module Microtiter Plate Adaptor , for 2 deep well microtiter plates
12003752	Rack for 13 mm Tubes , for 96 x 13 mm tubes
12003755	Rack for 16 mm Tubes , for 75 x 16 mm tubes
12003756	Rack for 18 mm/15 ml Tubes , for 70 x 18 mm tubes
12003757	Rack for 50 ml Tubes , for 27 x 29 mm tubes
12003754	Rack for Deep Well Microtiter Plates , for 2 x 24-, 48-, or 96-well plates
12003753	Rack for 1.5–2 ml Capless Microtubes , for 96 tubes
12003760	Plate Rack Riser

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

PEEK is a trademark of Victrex plc. Phillips is a trademark of Phillips Screw Co.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00 **France** 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026 674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

