



NGC Fraction Collector

From Discovery to Scale-Up, the NGC Fraction Collector Has the Flexibility and Collection Capacity Needed for Any Application



Flexible and Customizable

Adapts for Small- to Process-Scale Collection with Multivessel Collection Modes



The NGC Fraction Collector paired with the NGC Chromatography System has the capability to produce purified biomolecules for discovery and small-scale batch production on one hardware and software platform.

Access your fractions when, how, and in the collection vessel you need. With automated collection options for analytical to preparative chromatography applications, the NGC Fraction Collector:

- Ensures maximum flexibility within a method by collecting into any combination of four racks for 96-well plates, tubes, and bottles
- Collects based on customized trigger options for each phase: slope, %B, pH, and detector signals
- Enables faster downstream analysis with front-to-back dispensing that allows easy access to fractions before method completion

Setting a New Standard for Industry

There may be bottlenecks in other aspects of your research, but the NGC Chromatography System and NGC Fraction Collector will increase your purification efficiency. Along with the customizability and flexibility of the NGC System, the high-capacity and versatile NGC Fraction Collector streamlines the purification process. Optimize collection with the ability to start fractionation into any vessel you designate and quickly change the starting location if needed. Additionally, the combination of high capacity with user-specified preferences for personalized analysis works perfectly in multi-user settings. Queue multiple methods requiring various collection parameters, press start, and walk away.



- 1 Collection racks
- 2 Dispense head cover
- 3 Dispense head height adjustment
- 4 Gantry
- 5 Drip tray

Versatile Rack Options

Ideal for Meeting Your Exact Fractionation Needs

Stain-resistant metal racks offer extreme durability and factory calibration provides accurate positioning. Mix and match racks as needed to accommodate all your collection needs.

Increase Capacity When Needed



Standard Racks

The NGC Fraction Collector improves high-throughput purification workflows with its high-capacity collection capabilities. Need even more capacity? Couple with outlet valves or a second NGC Fraction Collector to reach the capacity your laboratory requires.

Experience Limitless Flexibility



Prep Rack Adaptor

Scale up your purification from milligrams to grams on a single platform. Maximize versatility in batch production with a flow rate capacity up to 200 ml/min and a prep rack adaptor that will carry projects from discovery to small-batch production. The prep rack adaptor provides up to 80 unlimited volume fractions.

Maintain Sample Integrity



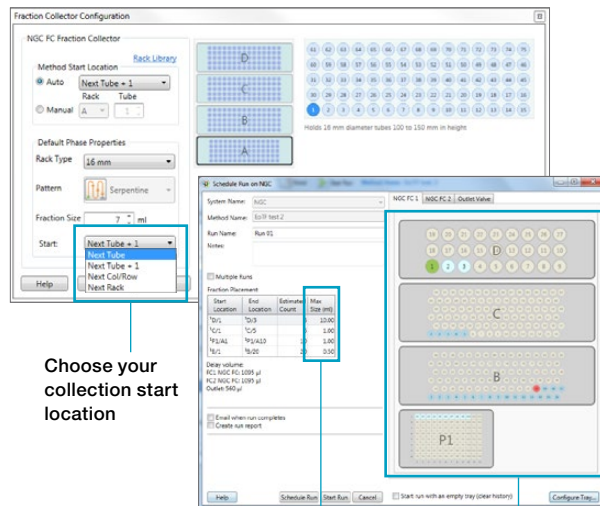
Peltier Cooling Modules

After all the hard work that goes into preparing samples for purification, you'll want to ensure temperature-sensitive biomolecules aren't degrading. Maximize valuable lab space and maintain sample integrity by integrating highly efficient benchtop Peltier cooling with optional enclosure to protect fractions from particulates.

NGC Fraction Collector Rack Capacity

Vessel Type/Size	Vessels/Rack	Fraction Collector Capacity
13 mm tube	96	384
16 mm tube	75	300
18 mm/15 ml tube	70	280
30 mm/50 ml tube	27	108
24-, 48-, or 96-deep well plate	2	8
1.5-2 ml microtube	96	384
250 ml bottle	4	16
Prep-20 rack	20	40

Get Flexible Collection Options with ChromLab Software



Choose your collection start location

Safeguard against going over your collection limit

Collect into different vessels and visualize where you fractionate through color coding by method phase

Specifications

Fractionation	Volume: 0.030–100,000 ml
Collection mode	Collect All, Threshold, Time/Volume windows
Peak detection	Intensity (UV, Conductivity, pH, %B), Slope (UV), Intensity or Slope, Intensity + Slope
Wetted parts	Polyetheretherketone (PEEK) Fluorinated ethylene propylene (FEP) Ethylene propylene diene monomer (EPDM) rubber
Weight	28.7 lb (13 kg)
Dimensions (W x D x H)	42 x 60 x 54.5 cm (16.5 x 23.6 x 21.5 in.)
Test tube change time	~200 ms with 13 mm tubes

Ordering Information

Catalog # Description

NGC Fraction Collector

17002070	NGC Fraction Collector , 100/240 V, includes 2 racks for 13 mm tubes
7880020IQOQ	NGC System with NGC Fraction Collector IQ/OQ
7880030IQOQ	NGC System with NGC Fraction Collector IQ/OQ
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 30 mm/50 ml Tubes , for 27 x 29 mm tubes
12003754	Rack for Deep Well Microtiter Plates , for 2 x 24-, 48-, or 96-deep well plates
12003753	Rack for 1.5–2 ml Capless Microtubes , for 96 tubes
12003760	Plate Rack Riser
12009848	NGC Fraction Collector Enclosure , includes cover and frame, 56 x 79 x 66 cm (22 x 31 x 26 in.) (W x D x H)

Peltier Cooling

12003747	Peltier Rack Module Microtiter Plate Adaptor , for 2 deep well microtiter plates
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
12003751	Insulated Rack for 15 ml Tubes , for 30 x 15 ml tubes
12003759	Insulated Rack for 50 ml Tubes , for 14 x 50 ml tubes
12005799	Peltier Cooling Kit , includes Peltier Rack Power Supply, 2 Peltier Rack Modules, and 2 insulated racks

Kits

12005822	Dispense Tip Replacement Kit , includes 2 tips
17002942	Prep-20 Preparative Rack Kit , includes funnel rack for 20 adaptors, tubing, and clamps

Visit bio-rad.com/web/FractionCollector for more information.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.

All trademarks used herein are the property of their respective owner.

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 01 877 89019 Belgium 32 03 710 53 00 Brazil 55 11 3065 7550
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 36 01 459 6192 Denmark 45 04 452 10 00 Finland 35 08 980 422 00 France 33 01 479 593 00
Germany 49 089 3188 4393 Hong Kong 852 2789 3300 Hungary 36 01 459 6190 India 91 124 4029300 Israel 972 03 963 6050 Italy 39 02 49486600
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 0 233 841 30
Poland 36 01 459 6191 Portugal 351 21 4727717 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 36 01 459 6193 Spain 34 091 49 06 580
Sweden 46 08 555 127 00 Switzerland 41 0617 17 9555 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300
United Kingdom 44 01923 47 1301

