

# 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



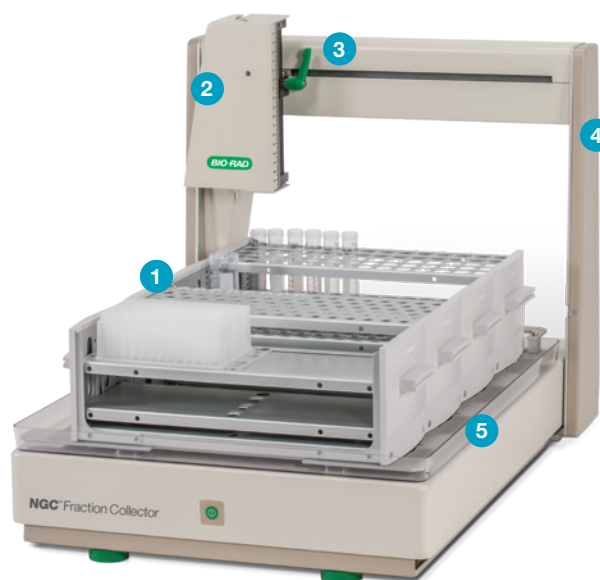
The NGC Fraction Collector paired with the NGC Chromatography System has the capability to produce purified biomolecules for discovery and small-scale batch production on one hardware and software platform.

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 4 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



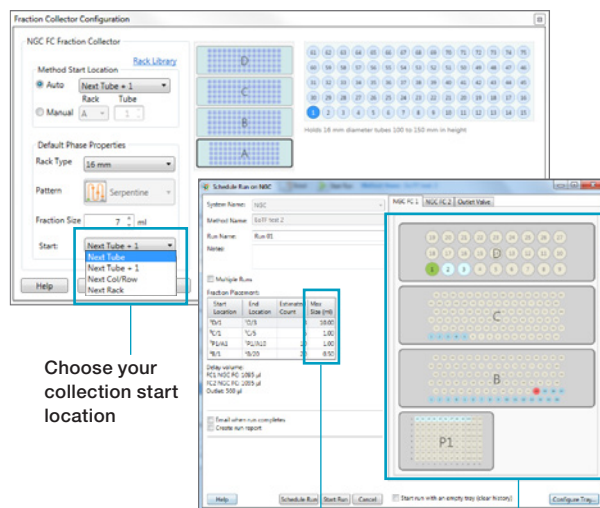
### Prep Rack Adaptor

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.

## 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	<b>NGC Fraction Collector</b> , 100/240 V, includes 2 racks for 13 mm tubes
7880020IQOQ	<b>NGC System with NGC Fraction Collector IQ/OQ</b>
7880030IQOQ	<b>NGC System with NGC Fraction Collector IQ/OQ</b>
12003752	<b>Rack for 13 mm Tubes</b> , for 96 x 13 mm tubes
12003755	<b>Rack for 16 mm Tubes</b> , for 75 x 16 mm tubes
12003756	<b>Rack for 18 mm/15 ml Tubes</b> , for 70 x 18 mm tubes
12003757	<b>Rack for 30 mm/50 ml Tubes</b> , for 27 x 29 mm tubes
12003754	<b>Rack for Deep Well Microtiter Plates</b> , for 2 x 24-, 48-, or 96-deep well plates
12003753	<b>Rack for 1.5–2 ml Capless Microtubes</b> , for 96 tubes
12003760	<b>Plate Rack Riser</b>
12009848	<b>NGC Fraction Collector Enclosure</b> , includes cover and frame, 56 x 79 x 66 cm (22 x 31 x 26 in.) (W x D x H)

### Kits

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

Visit [bio-rad.com/FractionCollector](http://bio-rad.com/FractionCollector) 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**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Website** [bio-rad.com](http://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

