



## NGC Chromatography Systems

Comprehensive Solutions for Protein Purification



# DESIGNED BY YOU. BUILT BY BIO-RAD.

## NGC Medium-Pressure Chromatography Systems

The NGC instrument is an automated liquid chromatography system focused on biomolecule purification at the research, process development, and laboratory-scale levels. At the core of the NGC platform is a truly modular and scalable system combined with a single, intuitive software package for system control and evaluation. Together, the NGC Systems provide a total laboratory solution.



### ALIGNS

A single solution that aligns to your needs today and expands to support your future discoveries and throughput requirements.



### ADAPTS

A flexible system that adapts to your requirements and can be easily customized to suit your application needs.



### ENSURES

An intelligent design that ensures functional simplicity and guides you from experimental setup to analysis and support.





# ALIGNS

A single laboratory chromatography solution that aligns and scales to fit your throughput requirements

NGC Systems can be selected based on customer needs and can be further customized to fit changing customer requirements through the addition of more modules and capabilities.

### Capabilities Included in All NGC Systems

Choice of 10 ml/min or 100 ml/min system pumps, mixer module with multiple mixer barrel options (750  $\mu$ l, 2 ml, 5 ml, 12 ml), automated sample inject valve, ChromLab Software, and a touch screen.

### Enhancements Available for All Systems

Increase automation and functionality by adding modules for different phases of your purification scheme. All systems are compatible with the versatile, high-capacity NGC Fraction Collector and BioFrac Fraction Collector for automated fraction collection (analytical- to preparative-scale). See bulletin 6326 for more details.



### NGC Quest System

Designed for the easy, dependable, and all-purpose purification of biomolecules with accurate gradients and high-resolution separations.

#### Base system includes:

- Single-wavelength (UV) and conductivity detection
- ChromLab Software, for fast, easy automated and manual control — single platform compatible with all NGC Systems

### NGC Quest Plus System

Designed for the all-purpose purification of biomolecules and simultaneous detection of proteins, peptides, nucleic acids, and other chromogenic molecules.

#### Includes NGC Quest capability, plus:

- Multi-wavelength (UV/Vis) detection of up to 4 wavelengths simultaneously



### NGC Scout System

Designed for quick, reliable separation of proteins and peptides. Enables rapid scouting of protein purification conditions with automated gradients and buffer preparation.

#### Includes NGC Quest capability, plus:

- Buffer blending valve for automated inline buffer preparation
- pH valve to monitor buffer pH and separation by pH gradients

### NGC Scout Plus System

Designed for the simultaneous detection of proteins, peptides, nucleic acids, and other chromogenic molecules with expanded automation and scouting.

#### Includes NGC Scout capability, plus:

- Multi-wavelength (UV/Vis) detection of up to 4 wavelengths simultaneously



### NGC Discover System

Designed for higher throughput, rapid and secure methods, and process development. Provides expanded scouting options with the simultaneous detection of proteins, peptides, nucleic acids, and other chromogenic molecules.

#### Includes NGC Scout Plus capability, plus:

- Integrated sample pump, 100 ml/min
- Sample inlet valves
- Column switching valve, 10 ml or 100 ml options

### NGC Discover Pro System

Designed for higher throughput, rapid and secure methods, and process development.

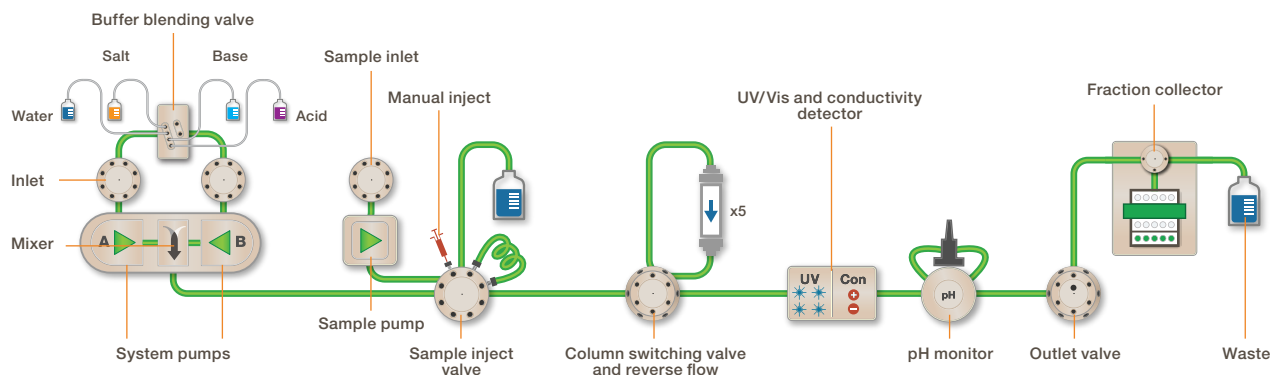
#### Includes NGC Discover capability, plus:

- Sample inlet valve
- Sample outlet valve

#### Options

- Tandem purification with additional column switching valve

# NGC SYSTEM CAPABILITIES



## System Pumps

Pump selection of up to 10 ml/min or 100 ml/min flow rates with the option to switch out pumps to meet your application requirements.

### F10 Pumps

- Flow rate of 0.001–10 ml/min at 3,650 psi (25.2 MPa)
- Ideal for small-scale preparative purifications
- Can also be used for analytical HPLC separations

### F100 Pumps

- Flow rate of 0.01–100 ml/min at 1,450 psi (10 MPa)
- Flexible flow rate range
- Ideal for scale-up applications

## Sample Pump

For automated sample application with the ability to load large sample volumes. Includes an integrated pressure sensor. Add a sample inlet valve or connect a third-party autosampler with a signal import module (SIM) for increased, automated sample loading capabilities.

## Mixer

Homogenizes buffers from two system pumps and can accommodate varying volumes (different sized barrels are available). Includes a mixer motor and integrated pressure sensor.

## Detectors

Ensure accurate detection of biomolecules such as proteins, peptides, nucleic acids, and chromophores. Include an integrated conductivity monitor (0.01–999 mS/cm) and an optional pH monitor (pH 1–14).

### Single-Wavelength (UV) Detector

For the detection of standard proteins (280 nm) or nucleic acids (255 nm).

### Multi-Wavelength (UV/Vis) Detector

For greater sensitivity and flexible detection of any biomolecules and chromophores (190–800 nm). Simultaneous multi-wavelength (UV/Vis) detection of up to four wavelengths.

Connect external detectors to the NGC System via the signal import module.

## Air Sensor

Detects end of buffer and sample to protect against column damage. Air sensor extension enables use of up to four additional air sensors (eight total).

## Valves

### Sample Inject Valve

For accurate sample loading (µl to L volumes) with a low internal volume for minimal sample loss.

### Buffer Blending Valve

For fast pH scouting with automated inline buffer preparation and the ability to double the fluid output to 20 ml/min or 200 ml/min.

### pH Valve

For accurate inline pH monitoring (pH 1–14). Includes integrated bypass valve and calibration port for in situ calibration.

### Buffer Inlet Valve

Automated switching between buffers (up to eight inlets per valve) for accelerated method development, column cleaning, and regeneration. Option to include two inlet valves, one for each system pump.

### Column Switching Valve and Reverse Flow

Automated column/media scouting of up to five columns without replumbing. Includes reverse flow for rapid elution, sample concentration, and column cleaning. Internal bypass allows automated system priming and cleaning with integrated pressure sensors that measure pre- and delta-column pressures.

### Outlet Valve

For enhanced automated fraction collection of large volume fractions with up to 12 vessels.

## Accessories

### NGC Fraction Collector (catalog #17002070)

Provides automated collection options for discovery to small-scale batch production at flow rates up to 200 ml/min. It supports multiple rack and vessel collection combinations from microplates and tubes to bottles and carboys.

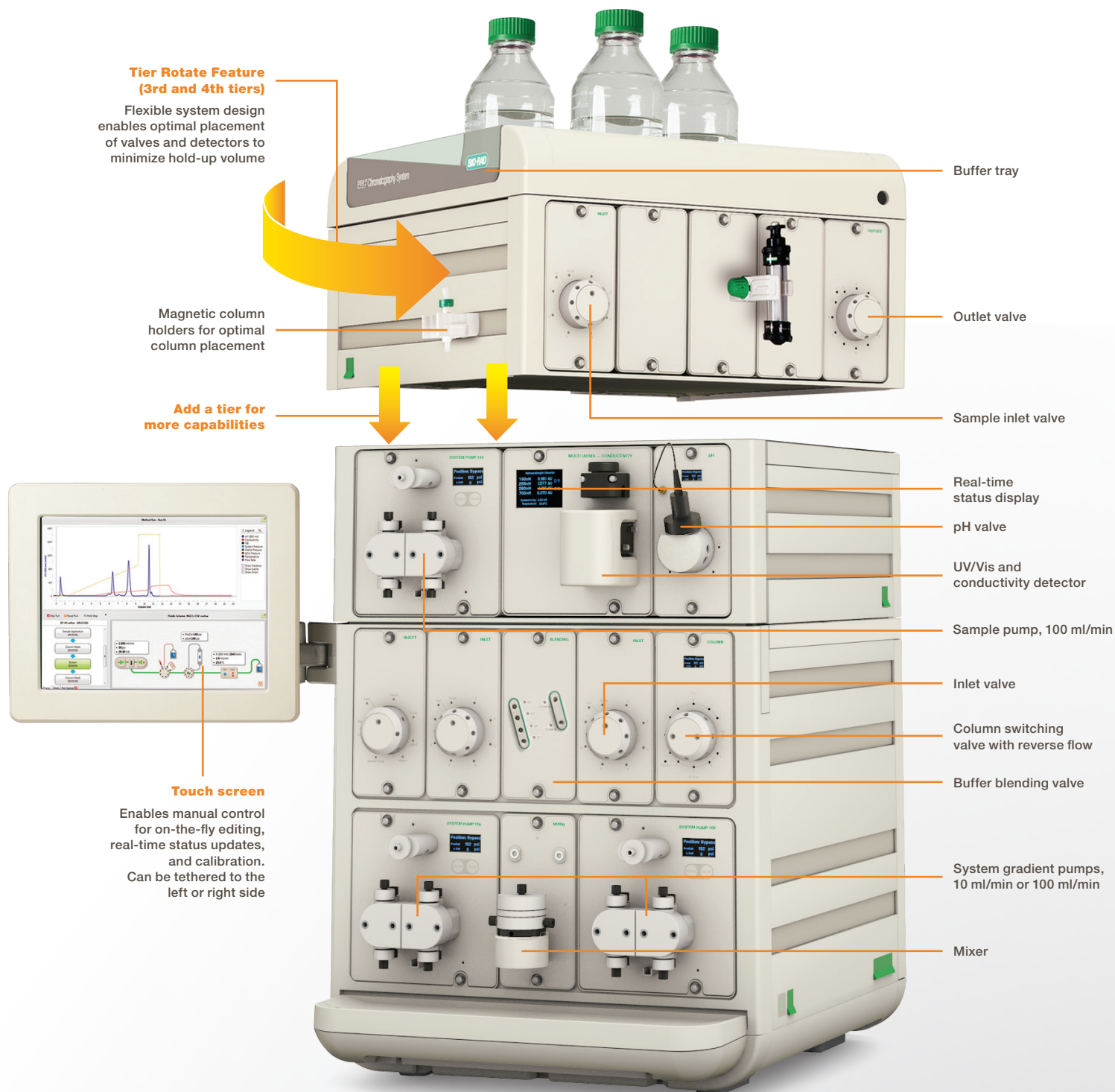


### BioFrac Fraction Collector (catalog #7410002)

Reliable fraction collection from analytical to preparative scale with versatile capability to collect from 96-well plates to 30 mm tubes.



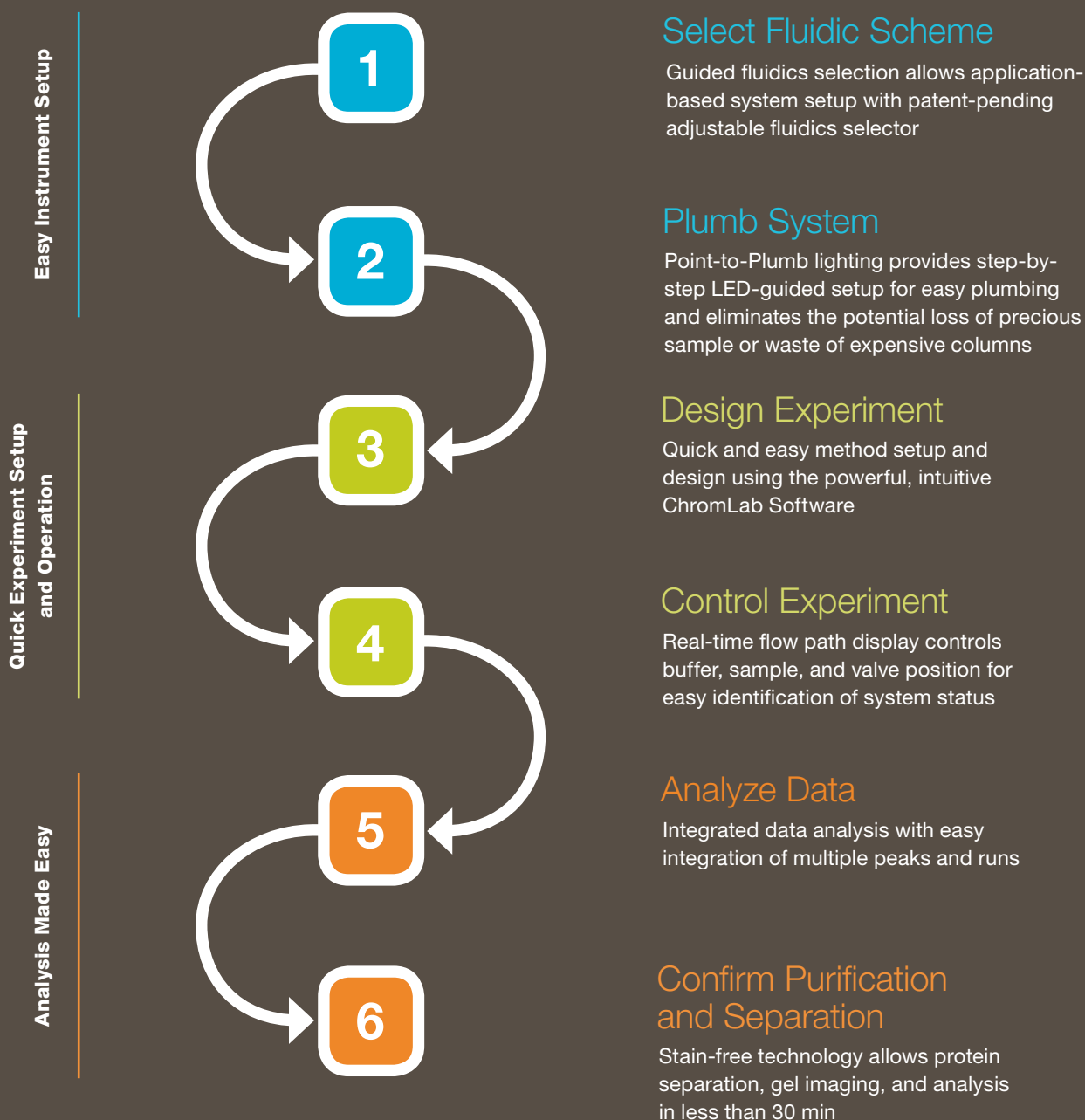
# PERSONALIZE AND EXPAND YOUR SYSTEM CAPABILITIES TO SUIT YOUR APPLICATION NEEDS AND WORKFLOW







Powerful ChromLab Software control, transferable across all NGC Systems, enables minimal training and fast setup to analysis.



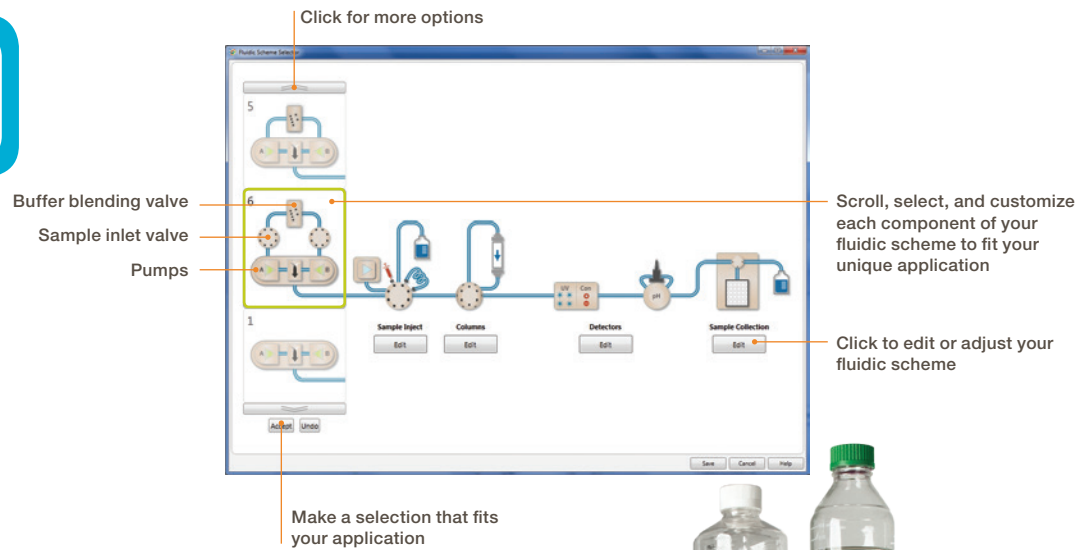
# EASY INSTRUMENT SETUP

1

## Select Fluidic Scheme

Select the fluidic scheme that best fits your application, set a default path, and optimize your module placement

To view the complete module library see bulletin 6326 or the NGC System Tour at [bio-rad.com/NGCSystems](http://bio-rad.com/NGCSystems)



**Choose** new modules to add to your system capabilities (for example, add a sample pump for loading large sample volumes)

**Change** module locations to adjust to your application and achieve optimal results

2

## Plumb System

Point-to-Plumb intuitive graphical indicators for simple, guided LED plumbing setup



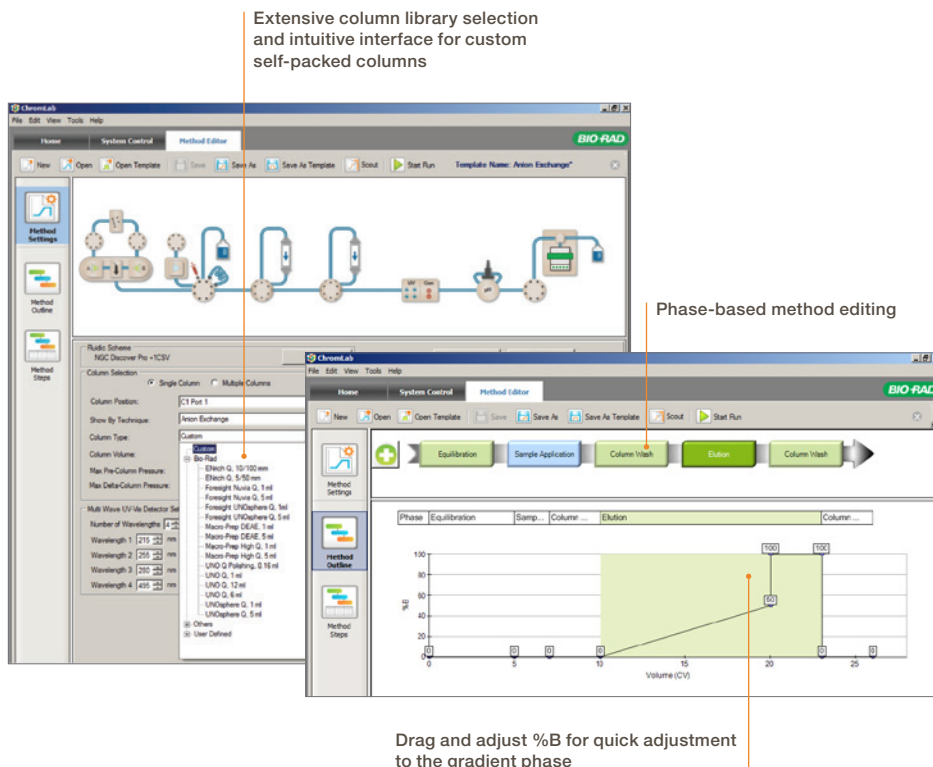
Click on each step in the flow path to guide system plumbing. Then, appropriate LEDs will light up to guide plumbing (as shown above).

# QUICK EXPERIMENT SETUP AND OPERATION

3

## Design Experiment

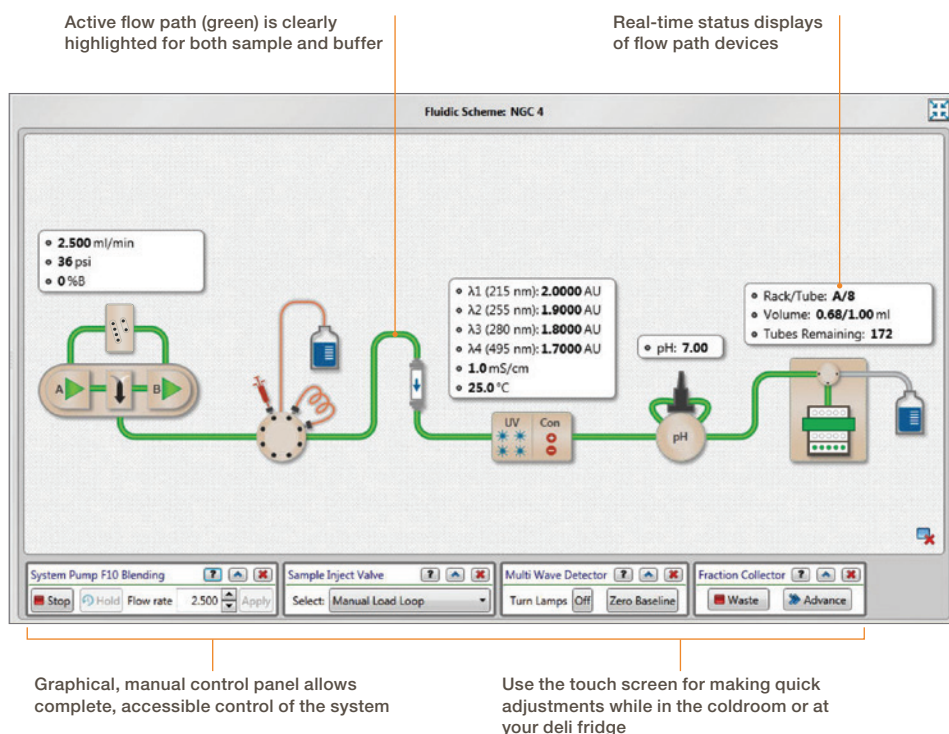
The ChromLab Method Editor enables confident, automated walk-away purification



4

## Control Experiment

Manual controls, conveniently located for quick and easy access, provide total graphical user control of the NGC System with a coldroom-compatible touch screen or a computer



For further details see the NGC System Tour at [bio-rad.com/NGCSystems](http://bio-rad.com/NGCSystems)



# ANALYSIS MADE EASY

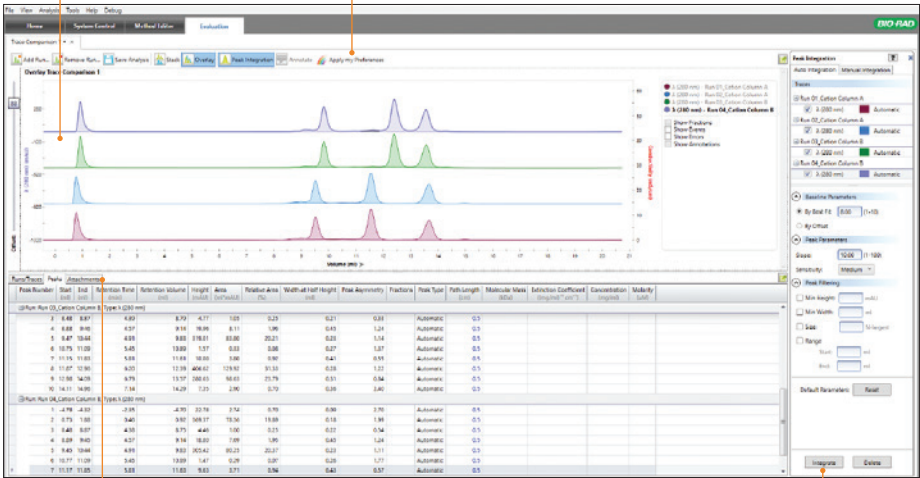
5

## Analyze Data

Comprehensive data analysis that enables fast, accurate data comparison

Ability to overlay and zoom in on multiple chromatograms

Enhanced user-specific preferences for personalized analysis



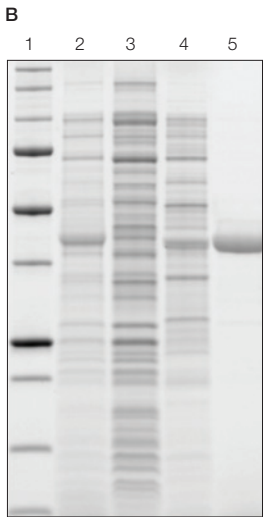
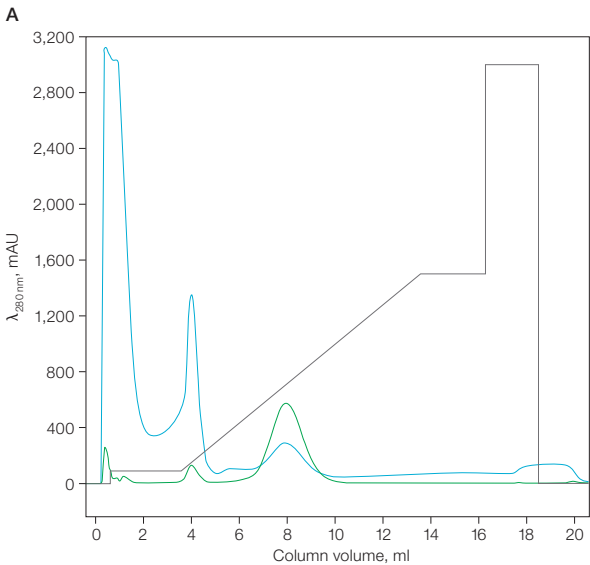
Keep all relevant purification data together like fraction gel images

Simple peak integration across multiple runs with single button autointegration

6

## Confirm Purification and Separation

Stain-free technology allows protein separation, gel imaging, and analysis in less than 30 min

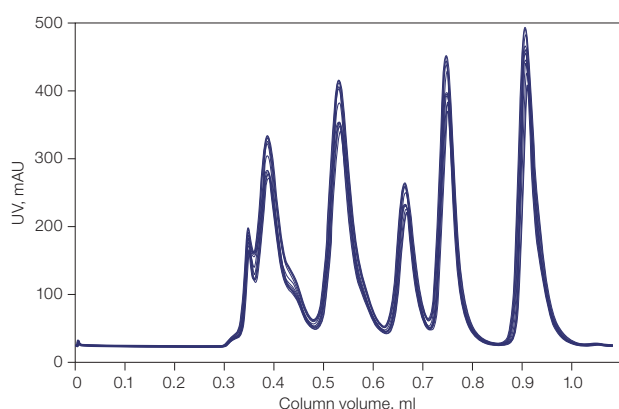


**Visual confirmation of chromatography results using stain-free gels and imaging. A**, isolation of a histidine-tagged green fluorescent protein (GFP) from a crude *Escherichia coli* lysate by affinity chromatography using an IMAC column; **B**, purification was confirmed by SDS-PAGE using a Criterion TGX Stain-Free Gel run for 20 min and directly visualized on the Gel Doc EZ Imaging System without the need for Coomassie staining. Samples in lanes 2 (crude *E. coli* lysates), 3 (flowthrough from the IMAC column), 4 (10% imidazole column wash), and 5 (purified histidine-tagged GFP) were compared against Precision Plus Protein Unstained Standards (lane 1).

## Intelligent design that guides your setup and operation

### Preplumbed System

Quality control (QC)–validated performance optimized for low hold-up volume translates to more reproducible results and sharper peaks.



**High-quality results with reproducible separations.** Eleven overlaid separations of a Bio-Rad size exclusion standard — composed of thyroglobulin,  $\gamma$ -globulin, ovalbumin, myoglobin, and vitamin B<sub>12</sub> — performed on the NGC Quest System with a 10 x 300 mm size exclusion column.

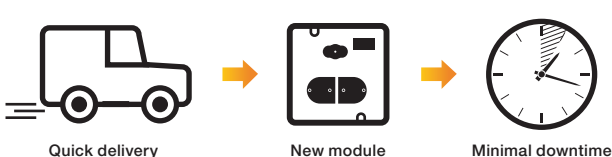
### Real-Time Status Displays

Provide immediate status of important parameters for clear diagnostics of key NGC instrument modules.



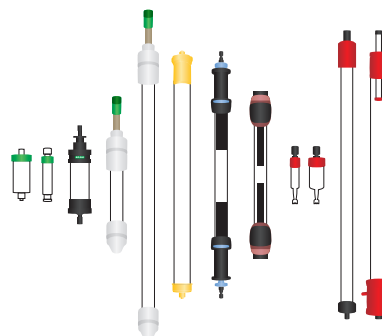
### Module Replacement Service

Directly replace plug and play modules — eliminate lengthy downtime and costly service visits.

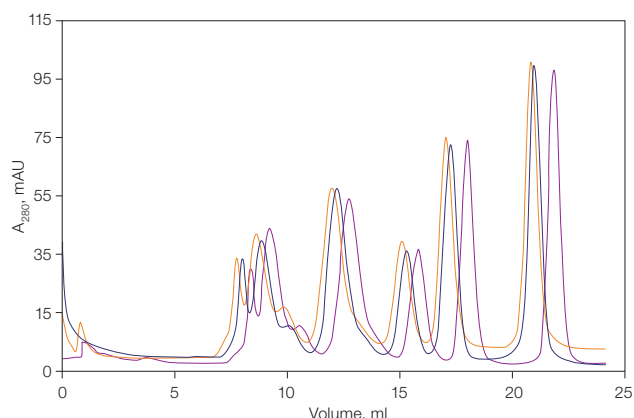


### Open Platform

Compatible with all medium-pressure columns and ChromLab Software includes method templates with column libraries.



Validated column applications on the NGC System.



**Completely transferable applications.** Identical comparisons of a Bio-Rad size exclusion standard (catalog #1511901) performed on a Superdex 200 10/300 GL Size Exclusion Column with separations performed on the NGC Quest (—), ÄKTApurifier (—), and ÄKTA avant (—) Systems.

# SELECTION GUIDE

## NGC Chromatography Systems

Catalog #	Product Description	NGC Quest 10 7880001	NGC Quest 10 Plus 7880003	NGC Quest 100 7880002	NGC Quest 100 Plus 7880004	NGC Scout 10 7880005	NGC Scout 10 Plus 7880007	NGC Scout 100 7880006	NGC Scout 100 Plus 7880008	NGC Discover 10 7880009	NGC Discover 100 7880010	NGC Discover 10 Pro 7880011	NGC Discover 100 Pro 7880012
7884002	NGC F10 Pump Module	●	●			●	●			●		●	
7884003	NGC F100 Pump Module			●	●			●	●		●		●
7884018	NGC Mixer Module	●	●	●	●	●	●	●	●	●	●	●	●
7884007	NGC Sample Inject Valve Module	●	●	●	●	●	●	●	●	●	●	●	●
7884008	NGC Single-Wavelength Detector Module, includes conductivity monitor	●		●		●		●					
7884009	NGC Multi-Wavelength Detector Module, includes conductivity monitor	○	●	○	●	○	●	○	●	●	●	●	●
7884010	NGC Buffer Blending Valve Module	○	○	○	○	●	●	●	●		●	●	●
7884011	NGC pH Valve Module, includes pH probe	○	○	○	○	●	●	●	●	●	●	●	●
7884004	NGC Sample Pump Module, integrated	○	○	○	○	○	○	○	○	●	●	●	●
7884006	NGC Inlet Valve Module	○	○	○	○	○	○	○	○	●	●	●	●
7884012	NGC Column Switching Valve Module, 10 ml	○	○	○	○	○	○	○	○	●	○	●	○
7884026	NGC Column Switching Valve Module, 100 ml	○	○	○	○	○	○	○	○	○	●	○	●
7884013	NGC Outlet Valve Module	○	○	○	○	○	○	○	○	○	○	●	●
12009390	ChromLab Software	●	●	●	●	●	●	●	●	●	●	●	●

- Standard
- Optional

**Note:** All NGC Systems include a touch screen and NGC Fittings Kit (catalog #7884017) and are compatible with the NGC Fraction Collector and BioFrac Fraction Collector.

Visit [bio-rad.com/web/NGCChromSystems](http://bio-rad.com/web/NGCChromSystems) for more information.



## Specifications

### System Specifications

<b>Control system</b>	ChromLab Software (compatible across all NGC Systems)
<b>Dimensions (W x D x H)</b>	49 x 61 x 56 cm (NGC Quest and NGC Scout Systems) 49 x 61 x 74 cm (NGC Discover System)
<b>Weight (excluding computer)</b>	41–46 kg (NGC Quest and NGC Scout Systems) 64 kg (NGC Discover System)
<b>Power supply</b>	100–240 V, 50–60 Hz
<b>Power consumption</b>	750 W maximum

### System Pump

<b>Pump type</b>	Reciprocating piston
<b>Flow rate setting</b>	<b>10 ml/min pumps:</b> 0.001 to 10 ml/min (normal range) <b>100 ml/min pumps:</b> 0.01 to 100 ml/min (normal range)
<b>Flow rate accuracy</b>	±2% (conditions: F10 pump — 0.1 to 10 ml/min, F100 pump — 1.0 to 100 ml/min; pressure: <600 psi [4.1 MPa, 41 bar]; viscosity: 0.5–3.7 cP)
<b>Pressure range</b>	<b>10 ml/min pumps:</b> 0 to 25.2 MPa (3,650 psi) <b>100 ml/min pumps:</b> 0 to 10 MPa (1,450 psi)
<b>Viscosity range</b>	0.5–10.8 cP (for 10 ml/min and 100 ml/min pumps)

### Sample Pump

<b>Pump type</b>	Piston pump, metering type
<b>Flow rate setting</b>	0.01 to 100 ml/min
<b>Flow rate accuracy</b>	±2%
<b>Pressure range</b>	0 to 10 MPa (1,450 psi)
<b>Viscosity range</b>	0.5–10.8 cP

### Mixer

<b>Mixing principle</b>	Chamber with magnetic stirrer
<b>Mixer volume</b>	263 µl (included), 750 µl (included), 2 ml, 5 ml (F10) 750 µl (included), 2 ml (included), 5 ml, 12 ml (F100)
<b>Gradient composition accuracy</b>	±0.5% (conditions: 3 to 97%B, 0.25 to 10 ml/min F10 pumps) ±0.8% (conditions: 5 to 95%B, 1 to 100 ml/min F100 pumps)

### Valves

<b>Type</b>	Rotary valves and rocker solenoid
<b>Number of valves</b>	1 inject valve, up to 2 x 8-port sample inlet valves, and 2 x 8-port buffer inlet, 2 x 12-port outlet, and 3 x 5-port column switching valves
<b>Functions</b>	Loop selection (PEEK Loop and DynaLoop offerings)

### Pressure Sensors

<b>Placement of sensors</b>	Standard: after system pump Options: precolumn, postcolumn
<b>Range</b>	0–3,650 psi
<b>Accuracy</b>	±2 psi or 2%, whichever is greater

### Inlet Valves

<b>Inlet A</b>	8 inlets
<b>Inlet B</b>	8 inlets
<b>Sample inlet</b>	8 inlets

**Note:** All NGC Systems include a touch screen and are compatible with the NGC Fraction Collector and BioFrac Fraction Collector.

### UV and UV/Vis Detectors

(each includes an integrated conductivity monitor)

<b>Wavelength</b>	<b>Single-wavelength:</b> 255 nm (nucleic acids) or 280 nm (proteins) <b>Multi-wavelength (up to 4):</b> 190–800 nm
<b>Absorbance range</b>	0 to >2.8 AU*
<b>Linearity</b>	0 to 2 AU within ±5%
<b>Operating pressure</b>	750 psi (5.2 MPa) for 2, 5, 10 mm flow cells
<b>Flow cells</b>	Preparative: 2 mm (cell volume: 20 µl) Analytical: 5 mm (cell volume: 16 µl) Analytical: 10 mm (cell volume: 18 µl)

\* For 5 mm and 10 mm flow cells.

### Conductivity Monitor

<b>Conductivity reading range</b>	0.01–999 mS/cm
<b>Accuracy</b>	±2%
<b>Operating pressure</b>	0–5.5 MPa (800 psi)
<b>Flow cell volume</b>	6 µl
<b>Temperature monitor range</b>	4–50°C
<b>Temperature monitor accuracy</b>	±2%

### pH Monitor

<b>pH reading range</b>	0 to 14
<b>Accuracy</b>	±0.1 pH unit within pH 2–12
<b>Operating pressure</b>	0 to 70 psi with pH probe inline and 0–500 psi in bypass mode
<b>Flow cell volume</b>	100 µl (210 µl including internal flow paths)

### NGC Fraction Collector

#### Collection modes

Collect All, Threshold, and Time/Volume windows

Flow rate	0.01–200 ml/min
-----------	-----------------

#### Collection rack options

(each NGC Fraction Collector can accommodate 4 racks)

	96 x 13 mm tubes, 75 x 16 mm tubes, 75 x 18 mm tubes, 27 x 50 ml tubes, 2 x deep well microplates (24-/48-/96-well), 96 x 1.5–2 ml capless tubes, 16 x 250 ml bottles, and 40 x unlimited volume prep-rack adaptors
<b>Peltier cooling option</b>	Yes
<b>Operating temperature</b>	4–40°C
<b>Dimensions (W x D x H)</b>	42 x 60 x 54.5 cm

### BioFrac Fraction Collector

#### Collection modes

Time	0.02–99,999 min
Volume	0.02–99,999 ml
Flow rate	0.01–100 ml/min

#### Collection rack options

	180 x 12–13 mm tubes, 120 x 15–16 mm tubes, 80 x 18–20 mm tubes, 168 x 1.5 ml microtubes, 24 x 30 mm tubes, 4 x 96-, 48-, 24-, or 12-position microplates, 4 x 250 ml bottles, and 20 x unlimited volume prep-rack adaptors
<b>Operating temperature</b>	4–40°C
<b>Dimensions (W x D x H)</b>	44.5 x 35.6 x 38.7 cm

### Column Switching

<b>Five-column valve</b>	Can connect up to 5 columns with forward and reverse flow and bypass capability
--------------------------	--

## Specifications (cont.)

### Buffer Blending

Standard in the NGC Scout and NGC Discover Systems

### Air Sensor Module

<b>Number of sensors</b>	Up to 8 total air sensors (1 for end of sample detection, remaining are buffer)
<b>Placement of built-in sensors</b>	End of buffer, end of sample
<b>Sensing principle</b>	Acoustic

## Ordering Information

### NGC Medium-Pressure Chromatography Systems

#### NGC Quest Chromatography Systems

For the all-purpose purification of biomolecules:

7880001	<b>NGC Quest 10 System</b>
7880003	<b>NGC Quest 10 Plus System</b>
7880002	<b>NGC Quest 100 System</b>
7880004	<b>NGC Quest 100 Plus System</b>

#### NGC Scout Chromatography Systems

For rapid scouting of proteins, peptides, and nucleic acids:

7880005	<b>NGC Scout 10 System</b>
7880007	<b>NGC Scout 10 Plus System</b>
7880006	<b>NGC Scout 100 System</b>
7880008	<b>NGC Scout 100 Plus System</b>

#### NGC Discover Chromatography Systems

For method development:

7880009	<b>NGC Discover 10 System</b>
7880011	<b>NGC Discover 10 Pro System</b>
7880010	<b>NGC Discover 100 System</b>
7880012	<b>NGC Discover 100 Pro System</b>

### NGC System Modules and Accessories

#### System Pumps

7884002	<b>NGC F100 Pump Module</b> , pkg of 1, includes 10 ml/min system pump kit with necessary tubing and fittings, for creating buffer gradients; for use with the buffer blending valve to generate flow rates of up to 20 ml/min
7884003	<b>NGC F100 Pump Module</b> , pkg of 1, includes 100 ml/min system pump kit with necessary tubing and fittings, for creating buffer gradients; for use with the buffer blending valve to generate flow rates of up to 200 ml/min

#### Sample Pump

7884004	<b>NGC Sample Pump Module</b> , pkg of 1, includes 100 ml/min sample pump kit with necessary tubing and fittings, for automated large-volume sample application via sample inject valve
---------	---

#### Detectors

7884008	<b>NGC Single-Wavelength Detector Module</b> , pkg of 1, includes UV/conductivity detector kit with necessary tubing and fittings, for nucleotide and protein detection, salt gradient generation
7884009	<b>NGC Multi-Wavelength Detector Module</b> , pkg of 1, includes UV/Vis and conductivity detector kit with necessary tubing and fittings, for simultaneous 4-wavelength monitoring of elution fractions between 190 and 800 nm and salt gradient generation

#### Mixers

7884018	<b>NGC Mixer Module</b> , pkg of 1, can be extended with 2, 5, and 12 ml barrels for efficient gradient mixing at higher flow rates, for use with all NGC Systems; does not include mixer base or barrels
7884019	<b>NGC F100 Mixer</b> , pkg of 1, 750 µl base and top assembly, included with all 100 ml/min NGC Systems
7884020	<b>NGC F10 Mixer</b> , pkg of 1, 263 µl base and top assembly, included with all 10 ml/min NGC Systems
7884021	<b>NGC F10 Mixer Barrel Kit</b> , pkg of 1, 750 µl extension barrel for F10 263 µl mixer, part of NGC Scout 10, NGC Discover 10 Systems
7884022	<b>NGC F10 Mixer Barrel Kit</b> , pkg of 1, 2 ml extension barrel for F10 263 µl mixer, optional part

7884028

**NGC F100 Mixer Barrel Kit**, pkg of 1, 2 ml extension barrel for F100 750 µl mixer, part of NGC Scout 100, NGC Discover 100 Systems

7884023

**NGC F100 Mixer Barrel Kit**, pkg of 1, 5 ml extension barrel for F100 750 µl mixer, optional part

7884024

**NGC F100 Mixer Barrel Kit**, pkg of 1, 12 ml extension barrel for 750 µl mixer, optional part

#### Valves

7884010

**NGC Buffer Blending Valve Module**, pkg of 1, kit includes necessary tubing and fittings for inline buffer preparation and generating pH gradients for quick pH scouting

7884006

**NGC Inlet Valve Module**, pkg of 1, kit includes necessary tubing and fittings for automated switching between multiple buffers and samples during method development

7884011

**NGC pH Valve Module**, pkg of 1, kit includes the pH valve kit, pH probe, tubing, and fittings for accurate inline pH measurement

7884012

**NGC Column Switching Valve Module (10 ml)**, kit includes the necessary tubing and fittings to accommodate the most common column types, holds 5 columns or sample loops; for use with F10 systems for quick column scouting, automated multicolumn, and reverse flow applications

7884026

**NGC Column Switching Valve Module (100 ml)**, kit includes the necessary tubing and fittings to accommodate the most common column types, holds 5 columns or sample loops; for use with F100 systems for quick column scouting, automated multicolumn, and reverse flow applications

7884013

**NGC Outlet Valve Module**, pkg of 1, kit includes necessary tubing and fittings for automated fraction collection of large-volume fractions with up to 12 vessels

7884016

**NGC Signal Import Module**, pkg of 1, enables analog to digital signal conversion and connection to third-party autosamplers and detectors

#### Air Sensors

7885017

**NGC Air Sensor Module**, pkg of 1, kit includes 2 large-bore air sensors for detection of end of buffer and sample to protect against air entering pumps and columns; supports up to 4 large- and small-bore air sensors

7885018

**NGC Air Sensor Extension Module**, pkg of 1, connects to the base air sensor module to support 4 additional air sensors; does not include any air sensors, optional part

7885020

**NGC Small Air Sensor**, pkg of 1, includes air sensor to exclude air from system and columns; detects air in small-diameter PEEK Tubing

7885021

**NGC Large Air Sensor**, pkg of 1, includes air sensor to exclude air from system and columns; detects air in large-diameter PTFE tubing

7885019

**NGC Air Sensor Extension Cable**, pkg of 1, for placement of air sensors outside air sensor module

#### Fraction Collectors

17002070

**NGC Fraction Collector**, 100/240 V, fraction collector compatible with all NGC Systems, includes power cord, rack set (two 13 mm tube racks), tubing, union

7410002

**BioFrac Fraction Collector**, 100/240 V, fraction collector compatible with all NGC Systems, includes power cord, rack set F1 (2 x flatpack, 13 mm), BioFrac Diverter Valve, fittings kit

7884025

**NGC Communication Adaptor**, pkg of 1, enables communication with Bio-Rad devices, such as the BioFrac Fraction Collector (#7410002), with the NGC System

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions. Precision Plus Protein Standards are sold under license from Life Technologies Corporation, Carlsbad, CA for use only by the buyer of the product. The buyer is not authorized to sell or resell this product or its components. All trademarks used herein are the property of their respective owner.



# DESIGNED BY YOU BUILT BY BIO-RAD

**BIO-RAD**



Available on the  
**App Store**



Download the NGC System Tour on the App Store  
and use this augmented reality (AR) target to visualize the  
NGC System in your lab.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

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

