

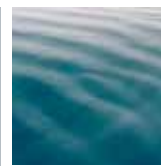


BioLogic DuoFlow™ Systems

A family of high-resolution chromatography systems.



BioLogic DuoFlow Systems



The BioLogic DuoFlow family of chromatography systems offers flexibility with multiple system configurations, many optional upgrades, and a common software platform that is intuitive and easy to follow. Key features of all systems include:

- Friendly, intuitive software
- Onscreen help menu that is easily accessible and provides extensive explanations, diagrams, and a system and application troubleshooting guide
- Modular hardware components that allow for optimal placement of valves, detectors, and columns to minimize dead volumes; modularity makes the systems easily upgradable to fit all applications and budgets
- Flow rate flexibility — the workstation is configured with an F10 or F40 pump to accommodate different flow rates; optional pump kits provide convenient reconfiguration at any time
- Make-Before-Break (MBB) technology in all automated valves to provide uninterrupted flow (when valve is switching positions), which eliminates baseline interference
- Endless valving options to perform diverse applications: reverse-flow chromatography, column switching, tandem chromatography, multidimensional chromatography, buffer scouting, and large sample loading
- A starter kit and UNO® Q1 column — easy-to-follow instructions enable the user to learn how to run a sample under controlled conditions
- Scouting wizard that provides quick and easy setup to optimize a purification method of a target molecule by changing a variable (e.g., buffer pH, flow rate, or length of step) over multiple runs automatically
- Method templates that allow easy method creation with predefined chromatography methods for all common chromatography applications, including affinity, ion exchange, chromatofocusing, hydrophobic interaction, hydroxyapatite, size exclusion, and multidimensional chromatography
- Trace Compare function that permits overlay of any number of different chromatograms for comparison of runs performed under different conditions (e.g., pH, flow rate, or buffer gradients)
- Ability to connect auxiliary pump for large sample loading (e.g., Econo™ gradient pump or Model EP-1 Econo pump)
- Ability to integrate up to 2 external detectors (e.g., refractive index or fluorescence detectors)
- Optional EZLogic™ integration software that offers additional data analysis functions to the BioLogic DuoFlow software
- Optional BioFrac™ ice bath/microplate rack to keep samples cool
- Coldroom compatibility



BioLogic DuoFlow Systems

BioLogic DuoFlow 10 System

10 ml/min flow rate at 3,500 psi
(233 bar, 23.3 MPa)

BioLogic DuoFlow 40 System

40 ml/min flow rate at 1,000 psi
(66 bar, 6.6 MPa)

Both systems include:

- Workstation with dual-piston pumps and valving connections — able to connect up to 3 low-pressure and 3 high-pressure valves
- Gradient mixer to ensure accurate gradient formation
- AVR7-3 automated injection valve with MBB technology
- Standard UV detector with fixed 254 and 280 nm filters; additional drop-in expansion filters (214, 365, 405, and 436 nm) are available
- Conductivity monitor that monitors salt concentration to ensure reliable gradient formation and pump functions
- BioFrac fraction collector that offers flexibility with 15 different rack options to collect microliter to liter sample volumes



BioLogic DuoFlow QuadTec™ Systems

BioLogic DuoFlow QuadTec 10 System

10 ml/min flow rate at 3,500 psi
(233 bar, 23.3 MPa)

BioLogic DuoFlow QuadTec 40 System

40 ml/min flow rate at 1,000 psi
(66 bar, 6.6 MPa)

Both systems include:

- Workstation with dual-piston pumps and valving connections — able to connect up to 3 low-pressure and 3 high-pressure valves
- Gradient mixer to ensure accurate gradient formation
- AVR7-3 automated injection valve with MBB technology
- BioLogic QuadTec™ UV/Vis detector that enables simultaneous monitoring of 4 wavelengths for greater sensitivity without background interference

- Conductivity monitor that ensures reliable gradient formation and pump functions
- BioFrac fraction collector that offers flexibility with 15 different rack options to collect microliter to liter sample volumes



BioLogic DuoFlow Maximizer™ Systems

BioLogic DuoFlow Maximizer 20 System

20 ml/min flow rate at 3,500 psi
(233 bar, 23.3 MPa)

BioLogic DuoFlow Maximizer 80 System

80 ml/min flow rate at 1,000 psi
(66 bar, 6.6 MPa)

Both systems include:

- Maximizer valve system that enables automated buffer blending, which eliminates manual buffer titrations; doubles flow rates for F10 and F40 pumpheads; and doubles valving capacity (6 low-pressure and 3 high-pressure valves)
- AVR7-3 automated injection valve with MBB technology
- Gradient mixer with barrel extenders to ensure accurate gradient formation at higher flow rates

- Standard UV detector with fixed 254 and 280 nm filters; optional filters (214, 365, 405, and 436 nm) are available
- Conductivity monitor that ensures reliable gradient formation and pump functions
- pH monitor — high-precision calomel electrode, fully compatible with buffers such as Tris
- BioFrac fraction collector that offers flexibility with 15 different rack options to collect microliter to liter sample volumes



BioLogic DuoFlow Pathfinder™ Systems

BioLogic DuoFlow Pathfinder 20 System

20 ml/min flow rate at 3,500 psi
(233 bar, 23.3 MPa)

BioLogic DuoFlow Pathfinder 80 System

80 ml/min flow rate at 1,000 psi
(66 bar, 6.6 MPa)

Both systems include:

- Maximizer valve system that enables automated buffer blending, which eliminates manual buffer titrations; doubles flow rates for F10 and F40 pumpheads; and doubles valving capacity (6 low-pressure and 3 high-pressure valves)
- AVR7-3 automated injection valve with MBB technology
- Gradient mixer with barrel extenders to ensure accurate gradient formation at higher flow rates

- BioLogic QuadTec UV/Vis detector that enables simultaneous monitoring of 4 wavelengths for greater sensitivity without background interference
- Conductivity monitor that ensures reliable gradient formation and pump functions
- pH monitor — high-precision calomel electrode, fully compatible with buffers such as Tris
- BioFrac fraction collector that offers flexibility with 15 different rack options to collect microliter to liter sample volumes

Ordering Information

Catalog #	Description	Catalog #	Description
BioLogic DuoFlow Systems*			
760-0037	BioLogic DuoFlow 10 System, 100/120 V, includes Dell controller and monitor, USB Bitbus communicator, workstation, MX-1 mixer, 3-tray rack, AVR7-3 sample inject valve, fittings kit, UV detector with 5 mm flow cell and 254/280 nm filters, conductivity monitor, starter kit, UNO Q1 column, instructions	760-2246	BioLogic DuoFlow Maximizer 80 System, 100/120 V, for Japan and Korea only
760-0036	BioLogic DuoFlow 10 System, 100/120 V, for Japan and Korea only	760-2248	BioLogic DuoFlow Maximizer 80 System, 220/240 V
760-0038	BioLogic DuoFlow 10 System, 220/240 V	BioLogic DuoFlow Pathfinder Systems*	
760-4037	BioLogic DuoFlow 40 System, 100/120 V	760-2257	BioLogic DuoFlow Pathfinder 20 System, 100/120 V, includes Maximizer valve unit, Maximizer mixer, pH monitor, Dell controller and monitor, USB Bitbus communicator, workstation, MX-1 mixer, 3-tray rack, AVR7-3 sample inject valve, fittings kit, QuadTec UV/Vis detector with 3 mm PEEK flow cell, system cable 25 (RS-232), conductivity monitor, starter kit, UNO Q1 column, BioFrac fraction collector, diverter valve, two F1 racks, instructions
760-4036	BioLogic DuoFlow 40 System, 100/120 V, for Japan and Korea only	760-2256	BioLogic DuoFlow Pathfinder 20 System, 100/120 V, for Japan and Korea only
760-4038	BioLogic DuoFlow 40 System, 220/240 V	760-2258	BioLogic DuoFlow Pathfinder 20 System, 220/240 V
760-0047	BioLogic DuoFlow 10 System With BioFrac Fraction Collector, 100/120 V, includes same as 760-0037 with diverter valve and two F1 racks	760-2267	BioLogic DuoFlow Pathfinder 80 System, 100/120 V
760-0046	BioLogic DuoFlow 10 System With BioFrac Fraction Collector, 100/120 V, for Japan and Korea only	760-2266	BioLogic DuoFlow Pathfinder 80 System, 100/120 V, for Japan and Korea only
760-0048	BioLogic DuoFlow 10 System With BioFrac Fraction Collector, 220/240 V	760-2268	BioLogic DuoFlow Pathfinder 80 System, 220/240 V
760-4047	BioLogic DuoFlow 40 System With BioFrac Fraction Collector, 100/120 V	Pump Kits	
760-4046	BioLogic DuoFlow 40 System With BioFrac Fraction Collector, 100/120 V, for Japan and Korea only	760-0110	F10 Pump Kit, converts F40 workstation to F10 pumps to enable flow rates as low as 0.01 ml/min; includes 2 fully assembled pumpheads, 4 piston assemblies, F10 tubing kit, tools, instructions
760-4048	BioLogic DuoFlow 40 System With BioFrac Fraction Collector, 220/240 V	760-0180	F40 Pump Kit, expands pumping capabilities to 40 ml/min; includes 2 fully assembled pumpheads with seals and check valves installed, 4 piston assemblies, mixer barrel extender, 2 mm UV flow cell, F40 tubing kit, tools, instructions
BioLogic DuoFlow QuadTec Systems*			
760-1137	BioLogic DuoFlow QuadTec 10 System, 100/120 V, includes Dell controller and monitor, USB Bitbus communicator, workstation, MX-1 mixer, 3-tray rack, AVR7-3 sample inject valve, fittings kit, QuadTec UV/Vis detector with 3 mm PEEK flow cell, instrument control module (ICM), system cables 25 and 26 (RS-232 and ICM power), conductivity monitor, starter kit, UNO Q1 column, instructions	BioLogic Maximizer™ Valve System	
760-1136	BioLogic DuoFlow QuadTec 10 System, 100/120 V, for Japan and Korea only	760-2200	BioLogic Maximizer Kit, 100/120 V, for US, Japan, Korea; includes Maximizer base unit, pH monitor, Maximizer mixer, starter kit, tubing kit, system cable 30, US power cord
760-1138	BioLogic DuoFlow QuadTec 10 System, 220/240 V	760-2204	BioLogic Maximizer Kit, 220/240 V, same as 760-2200 without US power cord
760-4137	BioLogic DuoFlow QuadTec 40 System, 100/120 V	Detectors and Accessories	
760-4136	BioLogic DuoFlow QuadTec 40 System, 100/120 V, for Japan and Korea only	750-0200	BioLogic DuoFlow Detector Kit, includes UV optics module and conductivity monitor, 5 mm analytical flow cell
760-4138	BioLogic DuoFlow QuadTec 40 System, 220/240 V	760-1300	BioLogic QuadTec Detector Kit, includes QuadTec detector with 3 mm PEEK flow cell, instrument control module (ICM), system cables 25, 26, and 17 (QuadTec RS-232, ICM power, and bus communication), US power cord, instructions, 40 psi backpressure regulator
760-1147	BioLogic DuoFlow QuadTec 10 System With BioFrac Fraction Collector, 100/120 V, includes same as 760-1137 with diverter valve and two F1 racks	760-2040	BioLogic DuoFlow pH Monitor, includes SIM module, pH electrode, flow cell, and tubing
760-1146	BioLogic DuoFlow QuadTec 10 System With BioFrac Fraction Collector, 100/120 V, for Japan and Korea only	750-0502	Signal Import Module, includes 4' communication cable (system cable 17), required for external detectors
760-1148	BioLogic DuoFlow QuadTec 10 System With BioFrac Fraction Collector, 220/240 V	760-2034	Universal AC/DC Inline Adaptor for USB Bitbus Device, required when connecting pH monitor or other external detector through SIM module on BioLogic DuoFlow systems
760-4147	BioLogic DuoFlow QuadTec 40 System With BioFrac Fraction Collector, 100/120 V	BioLogic DuoFlow Valves	
760-4146	BioLogic DuoFlow QuadTec 40 System With BioFrac Fraction Collector, 100/120 V, for Japan and Korea only	760-0406	AVR7-3 Automated Sample Injection Valve, 7-port, 3-position high-pressure valve, 3,500 psi (233 bar) limit
760-4148	BioLogic DuoFlow QuadTec 40 System With BioFrac Fraction Collector, 220/240 V	760-0408	AVR9-8 Stream-Select Valve, 9-port, 8-position high-pressure valve, 3,500 psi (233 bar) limit
BioLogic DuoFlow Maximizer Systems*			
760-2237	BioLogic DuoFlow Maximizer 20 System, 100/120 V, includes Maximizer valve unit, Maximizer mixer, pH monitor, Dell controller and monitor, USB Bitbus communicator, F10 workstation, MX-1 mixer, 3-tray rack, AVR7-3 sample inject valve, fittings kit, UV detector with 5 mm flow cell and 254/280 nm filters, conductivity monitor, starter kit, UNO Q1 column, BioFrac fraction collector, diverter valve, two F1 racks, instructions	760-0410	SVT3-2 Diverter Valve, 3-port, 2-position solenoid valve, 30 psi (2 bar) limit
760-2236	BioLogic DuoFlow Maximizer 20 System, 100/120 V, for Japan and Korea only	750-0415	SV5-4 Select Valve, 5-port, 4-position solenoid valve, 30 psi (2 bar) limit
760-2238	BioLogic DuoFlow Maximizer 20 System, 220/240 V	* The 10 and 20 systems include F10 workstation; the 40 and 80 systems include F40 workstation. 100/120 V systems for Japan and Korea and 220/240 V systems do not include Dell monitor.	
760-2247	BioLogic DuoFlow Maximizer 80 System, 100/120 V	Dell is a trademark of Dell Laboratories. MBB is a trademark of Rheodyne, LLC. MBB technology is covered by US patent 4,506,558, issued to Rheodyne, LLC. PEEK is a trademark of Victrex PLC.	



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300/5013478 Israel 03 963 6050 Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 55 5200 05 20 The Netherlands 0318 540666 New Zealand 64 9415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 095 721 14 04 Singapore 65 6415 3188 South Africa 27 0861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189/2578 7241 United Kingdom 020 8328 2000