



## BioLogic DuoFlow™ Valves

Expand Your Application Potential with  
Versatile and Reliable BioLogic DuoFlow Valves



# Making the Valve Connection

With Flexible Setup, Expanding Applications, and Software-Automated Functions



The numerous valve options available for the BioLogic DuoFlow Chromatography Systems easily accommodate simple to complex laboratory experiments. Valves can be easily added and programmed to expand the BioLogic DuoFlow System's capabilities, ensuring that you will always be able to run the experiments you need.



### SV5-4 Valve

A 5-port, 4-way low-pressure valve for:

- Selecting from up to 4 buffers or solutions
- Autosample loading from up to 4 samples directly onto a column or into an injection loop



### SVT3-2 Valve

A 3-port, 2-way low-pressure valve for:

- Selecting from 2 buffers or solutions
- Loading from 2 samples directly onto a column or into an injection loop
- Diverting buffer flow



### AVR7-3 Sample Inject Valve

A 7-port, 3-way high-pressure valve for:

- Loading samples into a static loop or a DynaLoop™ Dynamic Loop for injection onto a column
- Reverse-flow applications
- Switching between 2 columns



### AVR9-8 Select Valve

A 9-port, 8-way high-pressure valve for:

- Selecting from up to 8 buffers or solutions
- Autosample loading from up to 8 samples directly onto a column or into an injection loop
- Switching among 8 columns
- Collection of large-volume fractions

## Flexible Setup Leads to Optimized Results

The flexible setup design of the BioLogic DuoFlow System allows you to position valves for easy access and improved results. For example, place the sample valve outlet close to the column inlet for improved sample resolution.

## Expanding Applications Through Valving Options

BioLogic DuoFlow Systems are easily upgraded to meet your changing applications with valve additions. A total of 6 valves (3 low pressure and 3 high pressure) can be used at one time, allowing you to run more complex methods. Additionally, by adding a BioLogic Maximizer™ Valve System, the number of valves is doubled to 12 (6 low pressure and 6 high pressure).

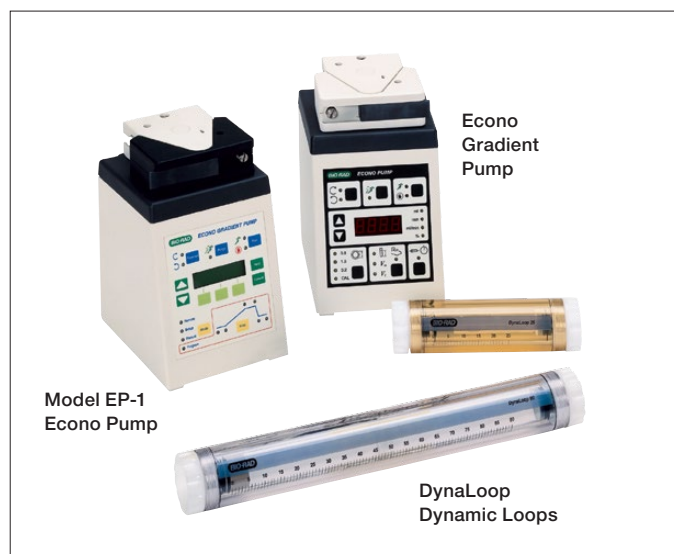
## Intuitive Software Automates Valve Functions

The easily programmed BioLogic DuoFlow Software incorporates valve functions into a method and allows each valve to be uniquely named to reflect its function.

# Valve Applications

## Flexible Sample Loading Options

- Basic AVR7-3 Sample Injection Valve operation using load, inject, and purge positions
- DynaLoop Dynamic Loops for large-volume sample loading or repetitive injections of smaller volumes
- Combination of a peristaltic pump, Model EP-1 Econo™ Pump, or Econo Gradient Pump with an inlet valve for:
  - Autosampling with automated loop fill and rinse
  - Sample loading directly onto a low-backpressure column
  - Repetitive loading of portions of large sample onto a column



## Multiple Buffer Selection

- Automated selection of multiple buffers, solutions, or samples

## Column Switching

- Automated switching between multiple columns
- Column scouting to compare separation characteristics within method

## Reverse Flow

- Elution of compounds bound to the top of a column by reverse flow

## Large-Volume Fraction Collection

- Collection of 2–8 large-volume sample fractions without a fraction collector

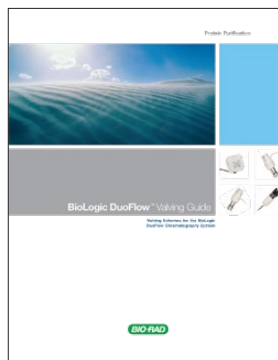
## Multidimensional Chromatography

- Automatic collection of selected fractions from one column and injection onto a second or third column

## The Make-Before-Break Advantage

The unique make-before-break design of the high-pressure valves rotates the inlet/outlet ports in a sequence that keeps the flow rate constant during port changes to:

- Prevent pressure spikes that cause baseline interference
- Prevent pump shutdowns from overpressure occurrences
- Protect fragile low-pressure columns and flow-sensitive detectors



Detailed schematics of valve-based applications are provided in the BioLogic DuoFlow Valving Guide (bulletin 2584 and on the Web at [bio-rad.com](http://bio-rad.com)) and the BioLogic DuoFlow Software Help screens.

## Valve Specifications

Valve	Type	Ports	Pressure Rating	Fitting	Tubing size
<b>SV5-4</b>	Low-pressure solenoid, 4-way stream selection valve	5 ports 4 positions	30 psi (2 bar)	1/4–28	1/8" (3.2 mm) OD
<b>SVT3-2</b>	Low-pressure solenoid, 2-way flow diversion or stream selection valve	3 ports 2 positions	30 psi (2 bar)	1/4–28	1/8" (3.2 mm) OD
<b>AVR7-3</b>	High-pressure, 3-way (load, inject, and purge) valve	7 ports 3 positions	3,500 psi (233 bar)	1/4–28	1/16" (1.6 mm) OD
<b>AVR9-8</b>	High-pressure, 8-way stream selection valve	9 ports 8 positions	3,500 psi (233 bar)	1/4–28	1/16" (1.6 mm) OD

## Guaranteed Reliability

Thousands of BioLogic DuoFlow Valves are in daily use worldwide. Composed of biocompatible PEEK polymer with no metallic parts, the valves are compatible with a wide range of buffers, salt solutions, and solvents (1.0 M salt to 100% acetonitrile) to accommodate any application requirement.

## Ordering Information

Catalog #	Description	Catalog #	Description
<b>BioLogic DuoFlow Valves</b>		<b>BioLogic Dynamic (Sliding Piston) Loops</b>	
760-0406	<b>AVR7-3 Automated Sample Injection Valve</b> , 7-port, 3-position high-pressure valve, 3,500 psi (233 bar) limit	750-0450	<b>DynaLoop Parts Kit</b> , includes 4 end cap O-rings, 1 sliding seal O-ring, 1 filter, 4 nut fittings, four 1/8" ferrules, five 1/4–28 nuts and ferrules, 10' of 1/8" PTFE tubing
760-0401	<b>AVR7-3 Valve Rebuild Kit</b>	750-0451	<b>DynaLoop 25 Kit</b> , includes 25 ml DynaLoop Sliding Piston Loop, DynaLoop Parts Kit, instructions
760-0408	<b>AVR9-8 Stream-Select Valve</b> , 9-port, 8-position high-pressure valve, 3,500 psi (233 bar) limit	750-0452	<b>DynaLoop 90 Kit</b> , includes 90 ml DynaLoop Sliding Piston Loop, DynaLoop Parts Kit, instructions
760-0403	<b>AVR9-8 Valve Rebuild Kit</b>	750-0455	<b>DynaLoop Sliding Seal Replacement</b>
760-0410	<b>SVT3-2 Diverter Valve</b> , 3-port, 2-position solenoid valve, 30 psi (2 bar) limit	750-0475	<b>DynaLoop 25 Sample Loop</b> , 25 ml replacement
760-0411	<b>SVT3-2 Valve Rebuild Kit</b>	750-0476	<b>DynaLoop 90 Sample Loop</b> , 90 ml replacement
750-0415	<b>SV5-4 Select Valve</b> , 5-port, 4-position solenoid valve, 30 psi (2 bar) limit	<b>Econo Gradient Pump</b>	
<b>BioLogic Single Injection Sample Loops</b>		731-9001	<b>Econo Gradient Pump</b> , 100/120 V, includes tubing and fittings kit
750-0471	<b>Sample Injection Port</b>	731-9002	<b>Econo Gradient Pump</b> , 220/240 V, includes tubing and fittings kit
125-0224	<b>Injection Needle</b> , 22 gauge, blunt	Contact your local Bio-Rad representative or visit us on the Web at <a href="http://bio-rad.com">bio-rad.com</a> for more information.	
750-0490	<b>Small-Volume Sample Loop Kit</b> , includes 100, 250, and 500 µl PEEK Loops	PEEK is a trademark of Victrex PLC. Tefzel is a trademark of E.I. du Pont de Nemours and Co.	
750-0491	<b>Large-Volume Sample Loop Kit</b> , includes 1, 2, and 5 ml PEEK Loops		
750-0482	<b>25 µl Tefzel Sample Loop</b>		
750-0483	<b>50 µl Tefzel Sample Loop</b>		
750-0492	<b>100 µl PEEK Sample Loop</b>		
750-0493	<b>250 µl PEEK Sample Loop</b>		
750-0494	<b>500 µl PEEK Sample Loop</b>		
750-0495	<b>1 ml PEEK Sample Loop</b>		
750-0496	<b>2 ml PEEK Sample Loop</b>		
750-0497	<b>5 ml PEEK Sample Loop</b>		



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 Belgium 03 710 53 00 Brazil 55 11 3065 7550  
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00  
France 01 47 95 69 65 Germany 49 89 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666  
New Zealand 64 9 415 2280 Norway 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 08 555 12700 Switzerland 026674 55 05  
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000