

BioLogic LP Diverter/Column Bypass Valve

Catalog Number 731-8322

The Model SV-3 valve may either be used as a Diverter Valve or a Column Bypass Valve. Fluid enters the valve through the "common" port and is diverted to either the "Bypass/Waste" or "Column/Collect" ports as commanded by the BioLogic LP and EGP Controller. Although the valve is the same, its function depends on which receptacle (at the rear of the BioLogic LP and EGP Controller) is selected. Used as a Diverter Valve, the SV-3 serves to either divert the eluant stream to waste or to collect it into fractions. Used as a Column Bypass Valve, the SV-3 serves to remove or maintain the chromatography column in line. This is useful for purging the system lines.

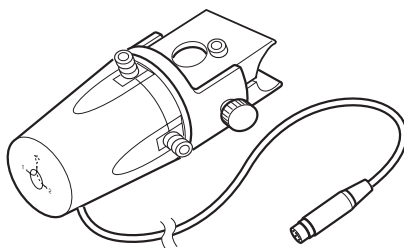


Figure 1. SV-3 Valve

Using the SV-3 as a Collect/Waste Diverter Valve

1. To install the valve onto the BioLogic LP and EGP Rack:
 - a. Ensure that the clamping screw is loose - loosen with a 1/4" Allen wrench, if necessary.
 - b. Clip the attaching bracket to a convenient place on the rack.
 - c. Position the valve, as desired, then tighten the clamping screw.
2. Plug the connector into the receptacle marked "Diverter Valve" on the rear panel of the BioLogic LP and EGP case.
3. Using the provided barb-to-male luers, connect an appropriate length of tubing between the conductivity flowcell and the "Common" port of the diverter valve (see figure 2). Match the appropriate luer connector with tubing ID, and ensure a firm connection but do not overtighten. Similarly, connect the "Collect" port to a fraction collector drophead and the "Waste" port to a waste container.

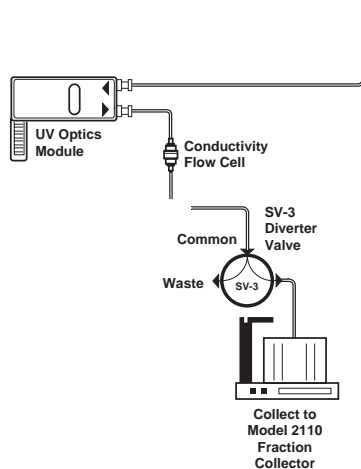


Figure 2. Using the SV-3 as a Diverter Valve

- To **manually operate** the LP Divert Valve, press the "Manual" key on the front panel of the BioLogic LP and EGP Controller then press the "Valve" key located near the bottom edge of the front panel. You may now select "Divert" or "Collect" using the soft keys located under the BioLogic LP and EGP Controller display.
- During **automated operation**, if a Method includes fraction collection, then the SV-3 diverter valve will automatically send fluid to the fraction collector during "collect" periods and divert fluid to waste when not collecting.

Using the SV-3 as a Column Bypass Valve

- To install the valve onto the BioLogic LP and EGP Rack:
 - Ensure that the clamping screw is loose - loosen with a 1/4" Allen wrench, if necessary.
 - Clip the attaching bracket to a convenient place on the rack.
 - Position the valve, as desired, then tighten the clamping screw.
- Plug the connector into the receptacle marked "Bypass Valve" on the rear panel of the BioLogic LP and EGP case.
- Using the provided barb-to-male luers, connect an appropriate length of tubing between the BioLogic LP and EGP Injection Valve and the "Common" port of the bypass valve (see figure 3). Match the appropriate luer connector with tubing ID, and ensure a firm connection but do not overtighten. Similarly, connect the "Column" port to the column inlet and the "Bypass" port to a waste container.
- To operate the Bypass Valve from the BioLogic LP and EGP Controller, press the "Manual" button on the Controller, then select the "Valve" key located near the bottom edge of the front panel. You may now select "Column" or "Bypass" with the soft keys below the BioLogic LP and EGP Controller display. Note that bypass valve operation is manual only - it cannot be programmed into a method.

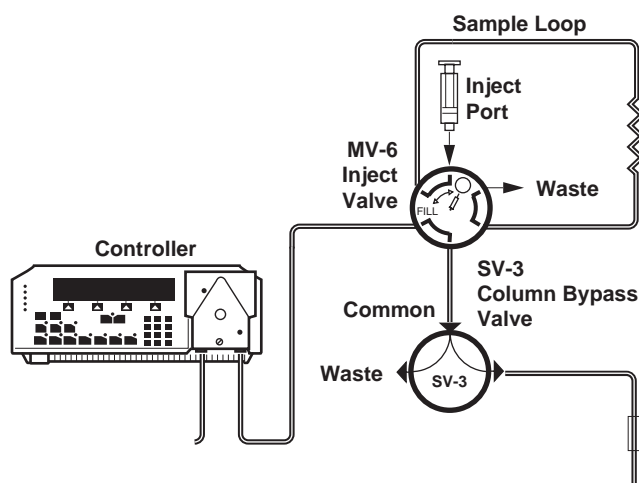


Figure 3. Using the SV-3 as a Column Bypass Valve

Valve Maintenance

After use, flush the valve with water, being sure to open both ports during the flush procedure. Valves allowed to stand for long period of time with high salt fluids may be damaged when the salts crystallize inside the valve. For long-term storage, store the valve in a 20% EtOH solution and do **not** allow the valve to dry out.

Technical Assistance

For additional help, contact your local Bio-Rad representative. In the United States, call Technical Service at 1-800-4BIORAD.

BIO-RAD

**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 2527 3454
Canada (905) 712-2771 China (86-21) 63052255 Czech Republic + 420 2 41 43 05 32 Denmark 44 52 10 00 Finland 09 804 22 00
France 01 47 95 69 65 Germany 089 318 84-0 Hong Kong 852-2789-3300 Hungary 36 1 455 8800 India (91-124)-6398112/113/114, 6450092/93
Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 55-52-00-05-20
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 7 095 721 1404 Singapore 65-6415 3188 South Africa 00 27 11 4428508 Spain 34 91 590 5200 Sweden 08 555 12 700
Switzerland 061 717-9555 Taiwan (8862) 2578-7189/2578-7241 United Kingdom 020 8328 2000