

SV5-4 SELECT VALVE

CATALOG NUMBER 750-0415

The SV5-4 Select Valve is a low pressure, 5-port, 4-position valve for stream selection with either the BioLogic HR or Duo-Flow System. Typically, it is used before the Gradient pump for buffer/solvent selection. Other uses could include stream selection for an Auxiliary pump. The valve is rated at 30 psi (max.). Refer to the BioLogic Instruction Manual for discussion of software control of the valve.

Connection to the BioLogic Rack and Workstation

The SV5-4 mounts to the System Rack using the clamp and Allen key provided with the valve. The valve is powered and controlled by the BioLogic Workstation. It plugs into any of the Workstation's rear panel connector ports identified as Solenoid Valves, ports 1, 2, or 3.

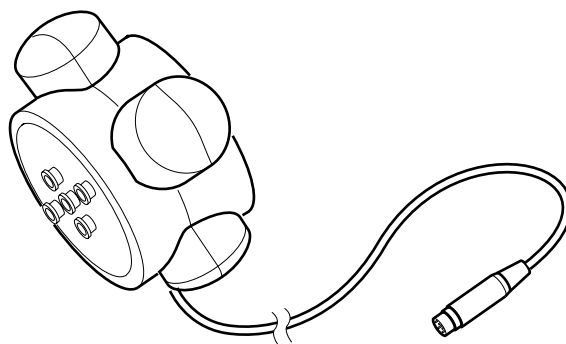


Figure 1. SV5-4 Select Valve

Example 1. Plumbing as a Gradient Pump Inlet Valve

To plumb the SV5-4 valve as a Gradient Pump Inlet valve, refer to the discussion below and the illustration on the following page:

1. Using 1/4-28 flat bottom fittings, connect a length of 1/8" OD (0.062" ID) PTFE tubing to any of ports 1 through 4 of the SV5-4.
Note: Only attach tubing to those ports that will deliver buffer. Block unused ports with a 1/4-28 plug.
2. Using 1/4-28 flat bottom fittings, connect a length of 1/8" OD (0.062" ID) PTFE tubing between port 5 (Common) of the SV5-4 and the inlet port of Pump A or Pump B of the Gradient Pump.
3. From the BioLogic Controller Manual screen, select each valve position by clicking on the appropriate position number and manually prime each line from the valve to the Gradient Pump.

Note: Always manually prime all pre-pump valves before initiating a run. Failure to displace air with buffer could stall the BioLogic Gradient Pump and/or damage the column during a chromatography run.

If the Workstation Purge feature is used, ensure that the sample Inject valve (AV7-3, V7-3, or AVR7-3) is also in the Purge position to bypass any column that would otherwise be in-line.

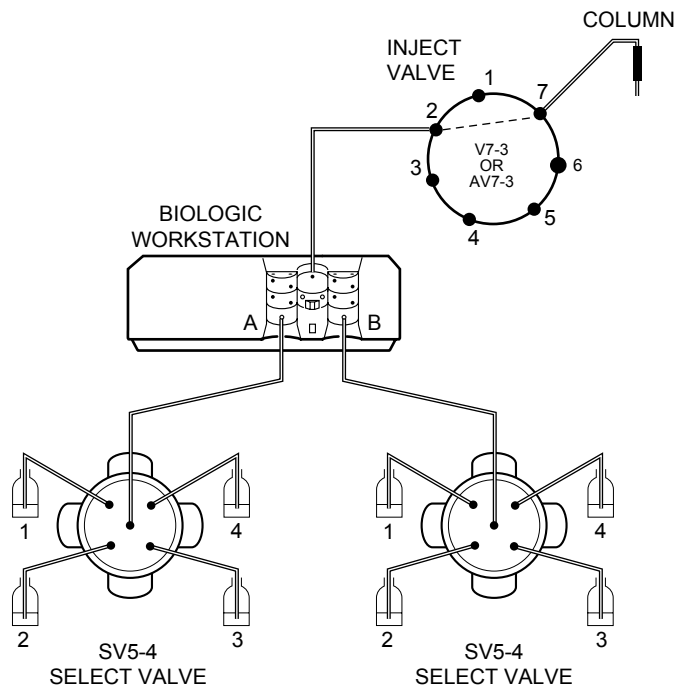


Figure 2. Configuration of the SV5-4 as a Gradient Pump Inlet Valve for buffer selection

Note that smaller diameter system tubing (1/16" OD, 0.020" ID) may also be used with this valve for other applications but the larger bore tubing **MUST** be used when the valve is used as a Gradient pump inlet valve.

Valve maintenance and repair

The Model SV5-4 valve should always be rinsed with water following use with aqueous buffers, or 20% ethanol following use with non-aqueous solutions. Never allow salt solutions to dry inside the valve.

Technical Assistance

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

Ordering Information

750-0415 Diverter Valve, SV5-4, BL
 750-0603 1/8" OD/0.062" ID Tefzel tubing, 15 feet