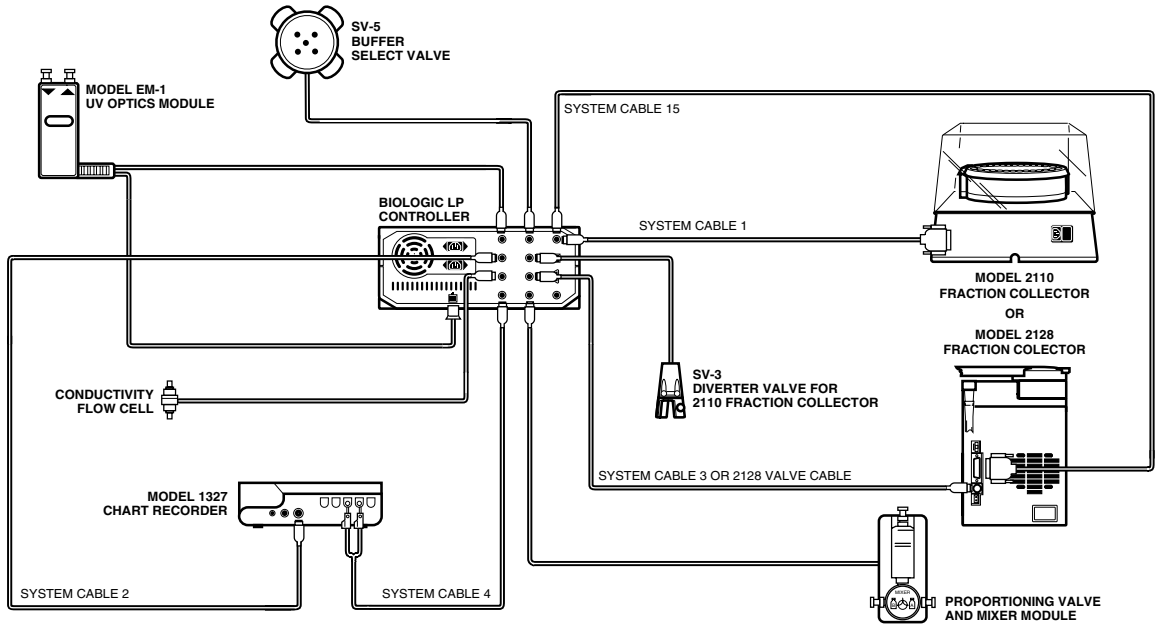


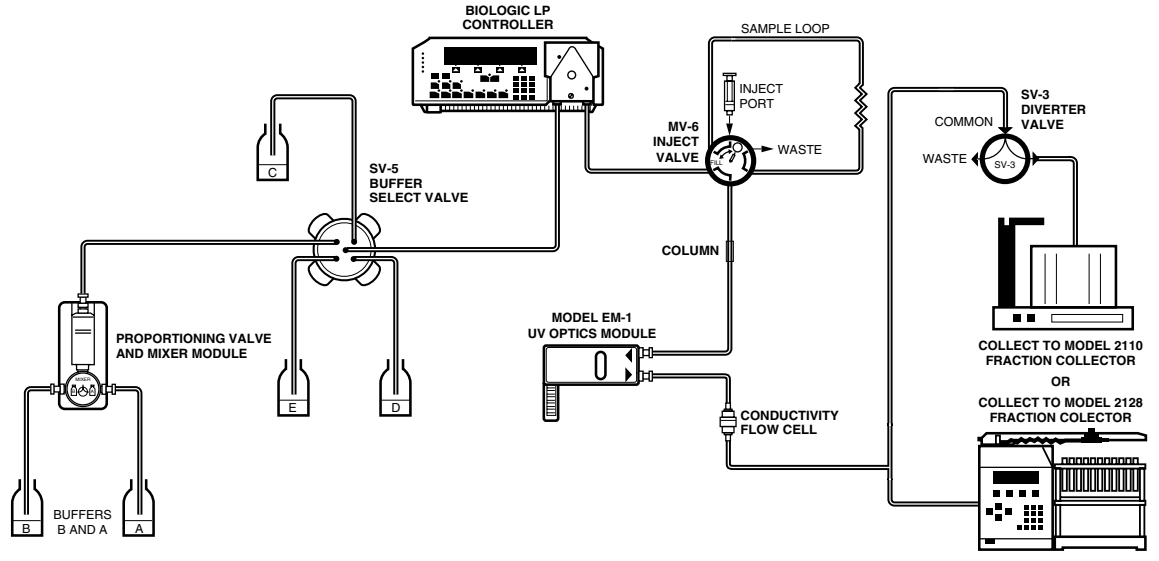
BioLogic LP Quick Start Tutorial

The BioLogic LP Starter Kit (Catalog # 731-8350) allows separation of a standard anion exchange mixture using the Econo Pac High Q cartridge. Cabling and plumbing, platen adjustment, UV setup, and conductivity min and max settings are to be completed prior to programming; instructions can be found in the BioLogic LP Starter Kit Instruction Manual, page 3-9. The illustrations below show the suggested cabling and tubing connections for the BioLogic LP system.

SYSTEM CABLING



SYSTEM TUBING



1. Buffer and Sample Preparation.

- Buffer "A": Empty the contents of the bottle labeled Buffer "A" into a 500 ml graduated cylinder. Add degassed, deionized water to 500 ml volume. Stir briefly. Label container as Buffer "A".
- Buffer "B": Prepare Buffer "B" and label container as described above.
- Sample: Dissolve the anion exchange standard in Buffer "A" using the syringe to measure 6.5 ml of buffer into the bottle of protein standards. Replace the stopper and shake gently to dissolve.

2. Press the Program mode key; select New Method.

3. Select Time programming mode.

4. Program the Pump.

- Press ADD.
- Select Buffer "A" with the Previous/Next keys; press OK.
- Enter the step length of 1 minutes; press OK.
- Enter flow rate of 3.0 ml/min; press OK.
- You have now entered the first step of the method (Buffer "A", step length of 1 minute, flow rate of 3.0 ml/min.) Enter the remaining steps.
Step 2. Gradient 0% to 50% "B", step length 5 minutes, flow rate 3.0 ml/minute.
Step 3. Buffer "B", step length 3 minutes, flow rate 3.0 ml/minute.
Step 4. Buffer "A", step length 3 minutes, flow rate 3.0 ml/minute.
- After entering Step 4; press OK.

5. Set the Alarm.

- Press the Alarm soft key.
- Press ADD.
- Enter a time of 1 minute for Alarm 1 and check that Hold Methods is set to NO.
- Press OK twice.

The alarm sounds after 1 minute to remind you to turn the valve back to the load position. If left in the Inject position, the gradient will flow through the loop first instead of going directly to the column.

6. Enter the fraction collector program.

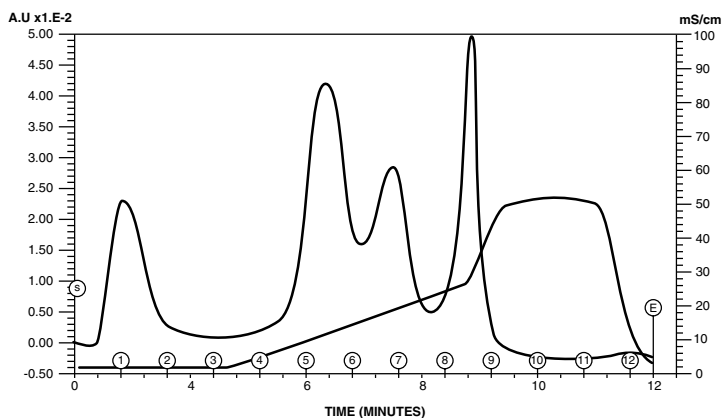
- Press the Frac Coll soft key.
- Select ALL.
- Enter a fraction size of 0.5 minute.
- Press OK twice, then select DONE.
- Press SAVE, name the method "DEMO 1", and press DONE.

7. Load Sample loop with protein standard.

- Turn the MV-6 Injector valve knob counter-clockwise as far as it will go.
- Draw 2 ml of the protein standard mixture into the syringe.
- Insert the syringe in the top port of the MV-6 injector valve, and fill the sample loop. Leave the syringe in the port when you have injected the sample — this prevents the sample from siphoning out of the loop.

8. Start Method.

- Press the Run mode key. The system will count down 10 seconds, then start the method.
- When the method starts, turn the injector valve knob to the right as far as it will go. When the alarm sounds, turn the valve knob to the left.
- If you are using a Bio-Rad 2128 fraction collector, press "Park", then "Yes" when method is complete.



RUN CONDITIONS

Column: Econo Pac High Q Cartridge (Catalog #732-0028)
Buffer A: 25 mM Tris HCl, pH 8.1
Buffer B: 25 mM Tris HCl, pH 8.1 + 0.5 M NaCl
Standard: Anion Exchange Std. (Catalog #125-0561)

Equine myoglobin is not retained on the column — it elutes in the void volume (first peak). The remaining three proteins bind to the column and elute separately as the salt concentration increases.



**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 21 507 6191
Canada (905) 712-2771 China 86-10-8201-1366/68 Denmark 45 44 52-1000 Finland 358 (0)9 804 2200 France 01 47 95 69 65 Germany 089 318 84-177
Hong Kong 852-2789-3300 India (91-124) 6398112/113/114 Israel 03 951 4124 Italy 34 91 590 5200 Japan 03-5811-6270 Korea 82-2-3473-4460
Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 47-23-38-41-30
Russia 7 095 979 98 00 Singapore 65-2729877 Spain 34-91-590-5200 Sweden 46 (0)8-55 51 27 00 Switzerland 061-717-9555 United Kingdom 0800-181134