

BioLogic™ LP System Unique Specifications

The BioLogic LP chromatography system includes the LP controller housing the peristaltic pump, the BioLogic rack, and the BioLogic LP accessory kit (MV-6 inject valve, proportioning valve/mixer, UV optics, conductivity cell, tubing and fittings kit, column and conductivity cell holder, user manual, starter kit). The BioLogic LP low-pressure chromatography system may also include a fraction collector (either Model 2110 or BioFrac fraction collector), BioLogic LP Data View software a dual-pen chart recorder (Model 1327), and extra valves (buffer select valve, column divert/bypass valve).

Unique Features of the BioLogic LP

Stores Up to 50 Methods. Each method can include up to 50 pump steps (flow rate, buffer selection, and step duration), 50 collection windows (with or without threshold values), and three alarms (with or without programmed hold).

Alphanumeric Naming Protocol. Identify the saved method with a name up to 12 characters (letters, numbers, spaces) in length.

Buffer Selection. Program up to five buffers with the addition of the buffer select valve.

Variable Programming Modes. Program the method in progress and edit future steps.

On Line Help. Pressing a dedicated HELP key on front panel of BioLogic LP immediately brings up the help screens relevant to the task in process.

Single-Point Control. All system components and functions are controlled from the BioLogic LP controller, including the fraction collector and chart recorder.

Flow Rates to 40 ml/min.

Conductivity Monitoring. Conductivity flow cell monitors gradient progress and is included with every LP system.

LP Dataview™ Software. UV and conductivity data can be transferred to any PC. Saves, stores, and prints methods and run data. Replaces expensive chart recorders and provides real-time continuous peak scaling.

Automatic Valve Sensing. Valves are automatically sensed when plugged into the BioLogic LP; therefore, no manual entry of system configuration is required.

Range Setting. UV and conductivity monitors have unique user-set range settings in addition to discrete range settings. The user-set range settings allow automatic 0–100% scaled output to the chart recorder.

Programmable Methods Cycling. Program a method to repeat up to 999 times.

Small Volume Mixing. 750 µl dynamic mixer for accurate, responsive, reproducible gradients.

Sample Injection Flexibility. Automated and manual sample injection options.

Portable Optics Module. Minimizes dead volume between column and UV detector for more precise protein separations.

Precalibrated Peristaltic Pump. Peristaltic pump flow rates are precalibrated for three common tubing sizes (0.8, 1.6, 3.2 mm ID tubing).

Flexible, Space-Saving, Modular, and Stackable Design. The compact design maximizes available workspace in the coldroom and on the lab bench.

Safety Certification. The BioLogic LP is certified to meet the I.E.C. 1010 safety standard for safety of laboratory equipment. (I.E.C. 1010 is an internationally accepted electrical safety standard for laboratory equipment.)