

Installation and Operation Guide

LP 'Data View Software

for the

BioLogic LP System

Catalog Number 731-8365

installation and Operation of **LP** Data View software for the **BioLogic LP** System

Note: Due to the ease of use of the **LP** Data View software, a conventional hard copy instruction manual is not included. This Installation and Operation guide will enable you to install **LP** Data View on your computer and to become familiar with its operation. See the on-line **HELP** function of **LP** Data View for additional information about features of this program.

Computer Requirements.

PC with one available serial port and with Windows **95** or Windows NT version **4.0** operating systems.

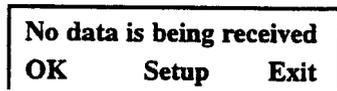
Section I installation

1. Ensure that both the **BioLogic LP** system and the computer are turned off.
2. Connect the serial cable (supplied with the software) from the **COMM** port on the rear of the **BioLogic LP** System to either **COM1**, **COM2** or **COM3** on the computer. You may need to use the supplied port adapter.

Note that **25** feet is the maximum **serial cable** length supported by the **BioLogic LP** system and **LP** Data View at this time.

3. Turn on the computer and install the **LP** Data View software as follows.
 - i. For both Windows **95** and NT users, insert diskette number **1**.
 - ii. Go to **Start, Settings, Control Panel, Add/Remove Programs** and follow the screen prompts to install **LP Data View**.
4. Turn on the **BioLogic LP** System and then launch the **LP** Data View application from the computer.
5. Data View uses **COM 2** as the default port.

If another **COM port** is used, you will see a dialog box reading:



6. Click on **Setup**. Another dialog box will ask you to select the correct **COM** port. Make your selection and click **OK**.
7. You should now see the main **Chromatogram** Run screen with the **Receive** indicator light green.
8. Congratulations! You are now ready to record your data.

Sections **2**, and **3** will provide a brief explanation of how to use Data View. A, more comprehensive explanation can be found by selecting the on-screen **HELP** function.

Section **2**. Using Data View as an electronic chart recorder

To Monitor Pre-Run Column Equilibration Steps

1. Click on **Record**. **LP** Data View will act just like a chart recorder and record **UV** and conductivity data from the **BioLogic LP** System. You may scale the axes at any time. To stop recording, click on **Stop**.
2. To clear the screen, click on **Clear**.

Recording Method Steps

1. If you wish to view and print the Method before it is run, use the following procedure.
 - i. Close any Run you are viewing and use the **Clear** button or **Chart Recorder** tab to obtain a blank chart recorder screen.
 - ii. Click on the **Record** button.

- iii. Go to the **BioLogic LP System** and press the **Program** hardkey.
 - iv. Select **List of Methods**. Select your Method of interest and open it using the **Open** softkey.
 - v. Press the **Program** hardkey again (the key light will remain on).
 - vi. Press the **Export** softkey to send the Method to Data View.
 - vii. The Method may now be viewed in Data View by clicking the **Method** button. Click the **Print button** to access print dialog box where you can choose to print just the Method.
2. After a Method has been run, the **BioLogic LP System** automatically sends the list of Method steps to the **LP Data View** application where it may be viewed by clicking on the **Method** button.

To Monitor a Run

1. If necessary, click the **Clear** button or the **Chart Recorder** tab to obtain a blank chart recorder screen.
2. When a Method is started from the **BioLogic LP System**, **LP Data View** will start automatically and will record all W and conductivity data as well as event marks such as fraction collection events. **Start, Hold, Continue,** and **Pause events** are automatically recorded.

To Save a Run

1. At the end of a Run, click on the **Save** button. You will be asked to name the Run. Runs are saved as **Data Files**.
2. **LP Data View** will automatically ask if you wish to save a Run if you try to clear the chart recorder screen or exit the application without a prior Save.

To Add Notes to a Run

1. Click on the **Notes** button.

To View The Method After The Run is Finished

1. Click on the **Method** button.

To Print The Run Chromatogram, Method and Notes

1. Open the Run of interest and **re-scale** axes as desired. Click on the **Print** button. A dialog box will ask you to select what Run parameters are printed.

Exporting Data to a Spreadsheet Program

1. Open the run of interest and **re-scale** the time axis as required.
2. From the **File** menu select **Export Data**. A dialog box will ask you to select what run parameters should be exported as a text file. This text file may then be opened in a spreadsheet program such as Microsoft® Excel.

Section 3 Using Data View with the Data Logging function

1. **LP Data View** includes a Data Logging feature which can be turned on and off using the **Log On/Log Off** buttons. The advantage of using this feature is that, in the event of a power failure during a Run, the temporary hidden file is not deleted and the Run data may be recovered.
2. With **Log On** selected, all data sent from the **BioLogic LP System** is automatically stored to the computer hard drive in a temporary hidden file. Data is stored even if the **LP Data View** application is not active as a chart recorder.
3. Log Data may be saved as a discrete Log File using the **Save Log** command under the **File** menu. A **Log File may be** opened using **Open Log** under the **File** menu. User-selected parts of a file may then be viewed, **re-named** and stored using the **Save selected data** as command under the **File** menu. The selected data is then stored as a Data File and may be viewed by clicking the **Open** button.
4. If the Log File is not saved by the user, then the temporary hidden file is deleted whenever the **LP Data View** application is exited. This prevents the hard drive of your computer from being filled with unnecessary data.

See the **HELP** feature for a more detailed description of **Data Logging**.

LP Data View Version 1.0 for the BioLogic LP system

Every effort has been made to ensure that the LP Data View software is as trouble-free as possible. However, as in any software application program, certain “bugs” are found after the software is released. This document lists the known bugs in version 1.0 and ways to work around them. These bugs will be fixed in a future release of LP Data View. please monitor our web site at www.bio-rad.com for information concerning software upgrades.

	Bug	Work-around
1. International versions of Windows' 95	The LP Data View program does not work correctly with international versions of Windows where the comma (,) and the period (.) characters used in formatting numbers are reversed from the US version of Window?.	Change the number formatting characters to match the US convention. This is done by opening the “Regional Settings” program within the Windows 95 “Control Panel”. Select the “Number” Tab and change the “Decimal symbol” to a period (.) and change the “Digit grouping symbol” to a comma (,). Restart your computer after making these changes. If this does not work for any reason, you must install the US version of Windows@95.
2. Manual mode fraction collection	The LP Data View program does not automatically record data when fraction collection is started in manual mode of the BioLogic LP system and does not reset the fraction number counter correctly in manual mode.	Click the “record” button on the toolbar of LP Data View to record data in manual mode before starting fraction collection and click on the “stop” button when fraction collection has been stopped. To reset the fraction number counter without running a method on the BioLogic LP you must exit Data View and then start the program again.
3. Time Axis Window Text Box	If an incorrect entry, for example a blank entry or a decimal point, is made in the Time Axis Window Text Box and the user clicks on “Set” or presses the “Enter” key, then the program will crash display the message “Run time error ‘13’. Type mismatch” and crash. Any unsaved data will be lost, but if Data Logging was ON the Log file can be used to recover the data when the program is started again. To prevent this problem, always ensure that the correct numerical entry is made beforehand.	If this situation arises, then PAUSE the run from the BioLogic LP Controller. Re-boot the LP Data View program, save any log file which will contain the pre-crash run data and click on “record”. Then resume the run from the BioLogic LP Controller.

4. Saving a Log file as a Data file	If a Log file is saved as a Data file without entering a different name for the file, the original. Log file remains but the Data file is not saved. Data View also incorrectly presents a warning about “replacing an existing file”. When this problem occurs, no data is lost.	To prevent the problem, enter a different name when saving a Log file as a Data file.
5. Setting axes	If multiple entries are made in the Axes Settings Text Boxes for different axes before clicking on the “Set!” button or pressing the “Enter” key, you must click on each “Set” button for all entries to take effect. The “Enter” key will only activate the “Set” button for the axis Text Box which has the focus.	Click on the “Set” button as appropriate for each axis.
6. Viewing multiple open applications	When data recording is started automatically as the BioLogic LP System starts a run, Data View takes the computer focus away from other applications and does not return it even if LP Data View is minimized.	The user should click on the program they were using previously or use “ Alt+Tab ” to send the focus back to the desired program.
7. Scroll bars active/inactive	While recording -data and using Window mode for the Time Axis, the scroll bar is not active.	The solution is to use the Full View setting for the Time Axis to see all data and then use the mouse again to “zoom in” on a new region.
8. Viewing runs of less than one minute duration	LP Data View does not use the scroll bar correctly when using the Window setting for the Time Axis when less than 1 minute of data is present.	The Full View setting will show all data and the mouse can be used to rubberband the data if necessary.
9. Printing in Landscape format	The program does not print in Landscape mode in Windows NT.	No work-around for this release.