
LP Data View Software, Version 2.0, for the BioLogic LP System

Installation and Operation Guide

Catalog number 7318365

BIO-RAD

Table of Contents

Section 1	Introduction	1
Section 2	Installation	1
Section 3	Using LP Data View Software as an Electronic Chart Recorder	2
	Monitoring Prerun Column Equilibration Steps	2
	Recording Method Steps	2
	Monitoring a Run	3
	Saving a Run	3
	Adding Notes to a Run	3
	Viewing the Method after the Run Is Finished.	3
	Printing the Run Chromatogram, Method, and Notes.	3
	Exporting Data to a Spreadsheet Program.	3
Section 4	Using LP Data View Software with the Data Logging Function	4
Section 5	Troubleshooting Guide	5

Section 1 Introduction

This installation and operation guide will enable you to install LP Data View Software on your computer and to become familiar with its operation. See the online HELP function of LP Data View for additional information about features of this program.

A PC with an available USB port and/or serial port running Windows 10 is required.

Note: Due to the ease of use of the LP Data View Software, a conventional hard copy instruction manual is not included.

Section 2 Installation

1. Ensure both the BioLogic LP System and computer are turned off.
2. Connect the supplied serial cable to the COM port on the rear of the BioLogic LP System. Note that 25 feet is the maximum serial cable length supported.
3. Connect the other end of the serial cable to your computer. If your computer has one or more serial ports, connect it to one of the serial ports. If not, then use the serial to USB adaptor to connect to an available USB port on the computer. Once connected restart the computer.
4. Install the LP Data View Software as follows:
 - a. If there is a previous version of LP Data View Software, then uninstall it via the control panel.
 - b. Go to bio-rad.com and download the LP Data View Software.
 - c. Unzip the LP Data View Software installer zip file.
 - d. Double click the LP Data View installer and follow the instructions.
5. Turn on the BioLogic LP System and launch the LP Data View application from the computer. First you will need to select a COM port.
 - a. When you open the software, an Enter BioLogic LP System Name dialog box will appear.
 - Enter the Identifier you will use for the BioLogic LP System that will be connected at the COM port being configured. If you are connecting multiple BioLogic LP Systems, make sure each Identifier is unique so you can distinguish between data from each system port
 - Press **OK** once you have entered the Identifier to save it
 - The Identifier may be edited by pressing **CollectData\LP System Name** in the menu

- b. Press **Setup** in the dialog box.
 - Select the first available (active) COM port
 - Press **OK**
 - To view the COM port setting, press **CollectData\COM Port** in the menu
 - c. To use LP Data View Software to collect data from a second BioLogic LP System, right click the LP Data View application icon and start a second instance of the application. Follow steps 5a and 5b to set the LP System Name and COM port for the second BioLogic LP System. This procedure may be used to connect up to ten BioLogic LP Systems to LP Data View Software.
6. You should now see the main Chromatogram Run screen with the Receive indicator light green.
 7. LP Data View Software is now ready to record data.

Section 3

Using LP Data View Software as an Electronic Chart Recorder

Monitoring Prerun Column Equilibration Steps

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

Recording Method Steps

1. To view and print the Method before it is run, use the following procedure.
 - a. Close any Run you are viewing and click **Clear** or **Chart Recorder** to obtain a blank chart recorder screen.
 - b. Click **Record**.
 - c. Go to the BioLogic LP System and press the **Program** hardkey.
 - d. Select **List of Methods**. Select your Method of interest and open it using the **Open** softkey.
 - e. Press the **Program** hardkey again (the key light will remain on).
 - f. Press the **Export** softkey to send the Method to Data View.
 - g. The Method may now be viewed in Data View by clicking **Method**. Click **Print** to access the 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 **Method**.

Monitoring a Run

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

Saving a Run

1. At the end of a run, click **Save**. You will be asked to name the Run. Runs are saved as Data Files.
2. LP Data View Software 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.

Adding Notes to a Run

Click **Notes**.

Viewing the Method after the Run Is Finished

Click **Method**.

Printing the Run Chromatogram, Method, and Notes

Open the Run of interest and rescale the axes as needed. Click **Print**. A dialog box will ask you to select which Run parameters to print.

Exporting Data to a Spreadsheet Program

1. Open the Run of interest and rescale the time axis as needed.
2. In the File menu, select **Export Data**. A dialog box will ask you to select which 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 4

Using LP Data View Software with the Data Logging Function

1. LP Data View Software 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 are automatically stored to the computer hard drive in a temporary hidden file. Data are 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 in the File menu. A Log File may be opened using **Open Log** in the File menu. User-selected parts of a file may then be viewed and stored using the **Save selected data** command in the File menu. The selected data are then stored as a Data File and may be viewed by clicking **Open**.
4. If the Log File is not saved by the user, then the temporary hidden file is deleted when the LP Data View application is exited. This prevents the hard drive of your computer from being filled with unnecessary data.

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

Section 5

Troubleshooting Guide

Every effort has been made to ensure that 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. Known bugs in version 2.0 and ways to work around them are listed in the following table. These bugs will be fixed in a future release of LP Data View Software. Go to bio-rad.com for information about software upgrades.

Feature	Bug	Work-Around
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 Record in the toolbar of LP Data View to record data in manual mode before starting fraction collection and click Stop when fraction collection has been stopped. To reset the fraction number counter without running a method on the BioLogic LP System, you must exit LP Data View and then start the program again.
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 Set or presses Enter , then the program will 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. Reboot the LP Data View program, save any log file that contains the precrash run data, and click Record . Then resume the run from the BioLogic LP Controller.

Feature	Bug	Work-Around
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 are lost.	To prevent the problem, enter a different name when saving a Log file as a Data file.
Setting axes	If multiple entries for different axes are made in the Axes Settings Text Boxes before clicking Set or pressing Enter , click each Set button for all entries to take effect. The Enter key will activate the Set button only for the axis Text Box that has the focus.	Click Set as appropriate for each axis.
Viewing multiple open applications	When data recording is started automatically as the BioLogic LP System starts a run, LP Data View takes the computer focus away from other applications and does not return it, even if LP Data View is minimized.	Click the program previously used or use Alt+Tab to send the focus back to the desired program.
Scroll bars active/inactive	While recording data and using Window mode for the Time Axis, the scroll bar is not active.	Use the Full View setting for the Time Axis to see all data. Then use the mouse again to zoom in on a new region.
Viewing runs of less than 1 min duration	LP Data View does not use the scroll bar correctly when using the Window setting for the Time Axis when less than 1 min of data are present.	The Full View setting will show all data and the mouse can be used to capture the data if necessary.
Printing in Landscape format	The program does not print in Landscape mode.	There is no work-around for this release.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.

All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00
Brazil 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00
Finland 35 08 980 422 00 **France** 33 01 479 593 00 **Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190
India 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670
The Netherlands 310 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 470 233 841 30 **Poland** 36 01 459 6191 **Portugal** 351 21 472717
Russia 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580 **Sweden** 46 08 555 127 00
Switzerland 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

