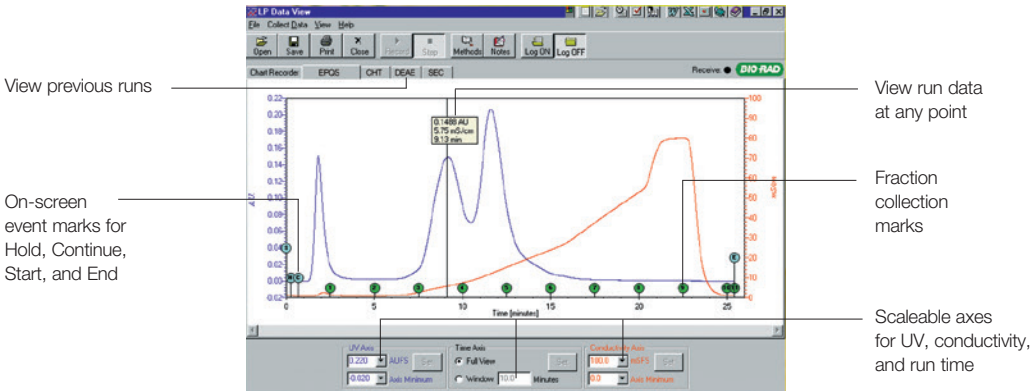


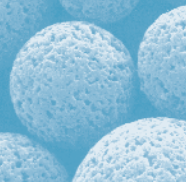
Off the Chart Flexibility

BioLogic LP Data View Software

Low-Pressure Chromatography for Protein Purification Just Got Better

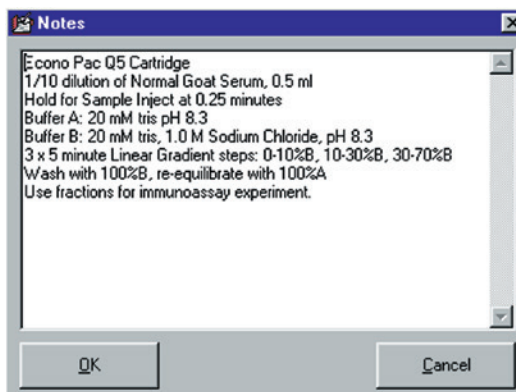
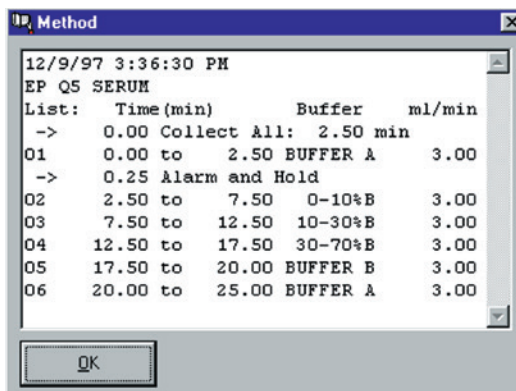
LP Data View Software is a simple, easy-to-use software that transforms any computer running Windows 10 into a powerful electronic chart recorder for the BioLogic LP System. Far more flexible than a traditional chart recorder, LP Data View Software allows total freedom to rescale chromatogram axes both during and after a run, to multitask during chromatogram data capture, and to print using any dedicated or networked printer. For more information, contact your local Bio-Rad representative.





LP Data View Software

- Automatically collects UV and conductivity data from the BioLogic LP System using any PC running Windows 10
- All three axes — UV data, conductivity data, and time — can be rescaled independently, both during and after data are collected, using the keyboard or the mouse
- Automatically records Method information for each run and allows Notes to be recorded with the data
- Automatically records run events such as **Start, End, Fraction Advance, Pause, Hold, Continue,** and **User Event Mark**
- Includes data logging feature, which can be used to record all data rather than only chromatogram run data and prevents loss of information in the event of a power failure
- Saves data on disk for future reference
- Selected parts of either run data or logged data can be saved in independent data files
- Prints customized reports, which can include the chromatogram, Method information, and/or Notes
- Exports data to other applications such as Microsoft Excel
- No special hardware required. Uses only a standard serial cable, which is provided (the PC must have one available serial port)
- Saves money on expensive chart recorder paper and ink pens



LP Data View Software creates a powerful research tool for the BioLogic LP Liquid Chromatography System. LP Data View Software automatically prints your Method and allows you to input Notes so you never lose track of important details about your separation runs.

Ordering Information

Catalog #	Description
7318365	LP Data View Software for the BioLogic LP System , includes link to download software from Bio-Rad website, installation and user guide, 25' serial cable, and serial to USB adaptor.

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** 47 0 233 841 30 **Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **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

