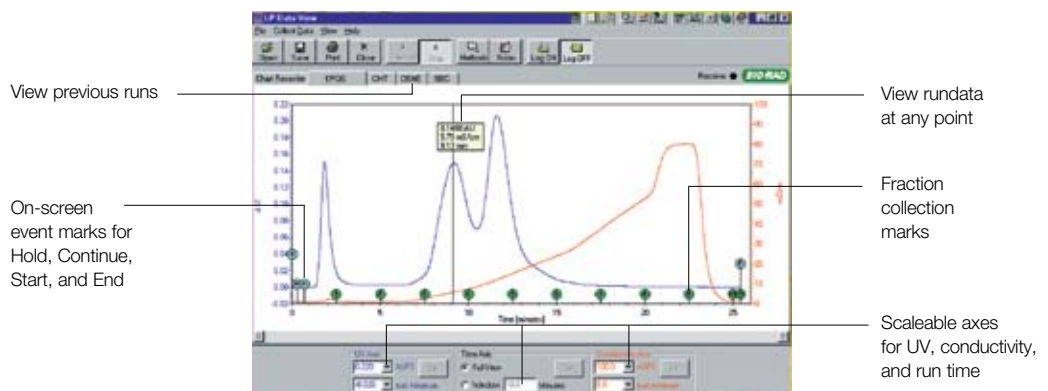


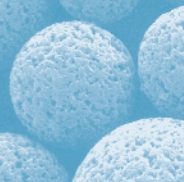
Off the Chart
Flexibility

BioLogic™ LP Data View™ Software

Low Pressure Chromatography for Protein Purification Just Got Better

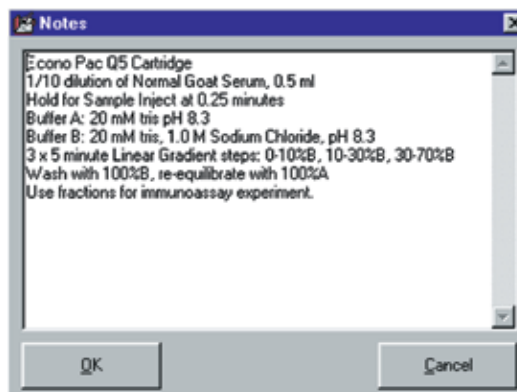
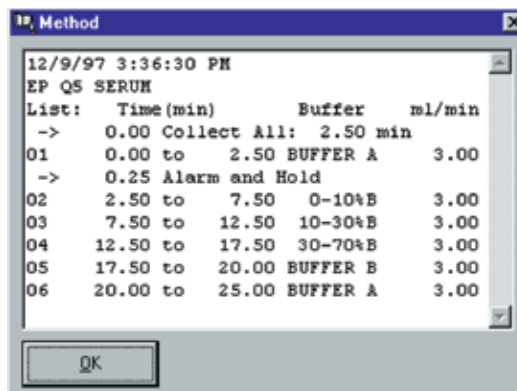
LP Data View software is a simple, easy-to-use software that transforms any computer running Windows 95, 2000, XP, or NT Version 4.0 into a powerful electronic chart recorder for the BioLogic LP system. Far more flexible than a traditional chart recorder, LP Data View software allows total freedom to rescale chromatogram axes both during and after a run, to multi-task 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 95, 2000, XP, or NT Version 4.0
- All three axes, UV data, conductivity data, and time, can be re-scaled 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
731-8365	LP Data View Software for the BioLogic LP System , includes software CD, 25' serial cable, port adaptor

Windows 95, 2000, XP, Windows NT, and Excel are trademarks of Microsoft Corporation.

BIO-RAD**Bio-Rad
Laboratories, Inc.**Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
 Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050
 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723
 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000