

BioLogic DuoFlow™ Chromatography System

Congratulations!

Enclosed is the BioLogic DuoFlow system software – version 4.0.

Enclosed you should find:

1. Software version 4.0 CD with installation instructions inside the CD case.
2. The New Features summary includes solutions to known discrepancies. This information is also found in your Read Me file.
3. BioLogic DuoFlow Instruction Manual
4. Windows 2000 software
5. USB Bitbus Communicator with USB and bus communication cables

If you are missing any items, please call the Technical Support number listed below.

Installation of 4.0 Software

- Installation instructions are found on the inside cover of the CD.
- Remember to save all of your current data either to disk or to your hard drive before installing your new software. For details of the copy-out/copy-in functions see pages 6-3 and 6-4 of the new instruction manual.

If you have additional questions, please call Bio-Rad Technical Support at 1 800 4-BIORAD, 1 800 424-6723.