

BioLogic DuoFlow™ Chromatography System Software Version 5.0

Enclosed is the BioLogic DuoFlow software 5.0. There are three software package configurations to support the various computer operating system options.

If you ordered the BioLogic DuoFlow software 5.0 only (Catalog # 760-2051) you will find:

1. BioLogic DuoFlow software 5.0 CD with installation instructions on the inside of the CD case.
2. A "New Features" summary. This information is also found in the Read Me file.
3. BioLogic DuoFlow Software 5.0 instruction manual.

If you ordered the BioLogic DuoFlow Software 5.0 and Windows 2000 Upgrade Kit (Catalog # 760-2052) you will find:

1. BioLogic software version 5.0 CD with installation instructions on the inside of the CD case.
2. A "New Features" summary. This information is also found in the Read Me file.
3. BioLogic DuoFlow Software 5.0 instruction manual.
4. Windows 2000 software
5. USB Bitbus Communicator with USB and bus communication cables

If you ordered the BioLogic DuoFlow Version 5.0 Software Upgrade Kit for New Windows 2000 Computers (Catalog # 760-2053) you will find:

1. BioLogic DuoFlow Software 5.0 CD with installation instructions on the inside of the CD case.
2. A "New Features" summary. This information is also found in the Read Me file.
3. BioLogic DuoFlow Software 5.0 instruction manual.
4. USB Bitbus Communicator with USB and bus communication cables.

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

BioLogic DuoFlow Software 5.0 Installation

Installation instructions are found on the inside cover of the installation CD. Remember to save all of your current data either to your hard drive or a CD before installing the new software. For details of the copy-out/copy-in functions see Section 6.1 of the BioLogic DuoFlow 5.0 instruction manual.

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