

AUTOMATED ELECTROPHORESIS

Experion™ System Software

Quick Installation Guide

Windows XP Operating System

- 1 Insert the Experion software CD into the CD-ROM drive. Installation will begin automatically. (If not, follow the instructions printed on the CD booklet.)
- 2 Connect one end of the Experion system USB2 cable to the computer and the other end to the electrophoresis station. The USB2 cable is shipped with the electrophoresis station.
- 3 Connect the power cable to the electrophoresis station, and plug the other end into an appropriate outlet.
- 4 Power on the electrophoresis station by pressing the green button in front. Installation is a two-part process with installation of the I/O firmware driver first, followed by installation of the system device driver.
- 5 A dialog box appears when the installation begins (Figure 1). Select “No, not this time” and “Next” to continue the installation.

Part 1. Installation of the Experion System I/O Firmware Load Device Driver

- 6 A dialog box prompts you to begin installation of the device drivers for the Experion system (Figure 2). Select “Next”.
- 7 A warning dialog box appears (Figure 3). Disregard the warning and select “Continue Anyway”.
- 8 The next dialog box acknowledges completion of the first part of device driver installation (Figure 4). Select “Finish”.



Fig. 1.



Fig. 2.



Fig. 3.



Fig. 4.

Part 2. Installation of the Experion System Device Driver

- 9 A dialog box prompts you to begin installation of the Experion system device driver (Figure 5). Select “Next”.
- 10 A second warning dialog box appears (Figure 6). Again, disregard the warning and select “Continue Anyway”.
- 11 The next dialog box acknowledges completion of the second part of device driver installation (Figure 7). Select “Finish”.
- 12 Open the Experion software application by double-clicking on the Experion icon.
- 13 Verify that the electrophoresis station and software are communicating and fully functional by checking for a green dot and a 6-digit serial number in the bottom right-hand corner of the software application screen.

Installation is complete!



Fig. 5.



Fig. 6.



Fig. 7.

Ordering Information

Catalog #	Description
700-7050	Experion Software , system operation and standard data analysis tools, PC
700-7051	Experion Validation Kit , 3 test chips, qualification procedures, dongle, PC
700-7052	Experion Security Edition Software , system operation and standard and 21 CFR Part 11 data analysis tools, 3 test chips, qualification procedures, dongle, PC



LabChip and the LabChip logo are trademarks of Caliper Life Sciences, Inc. Bio-Rad Laboratories, Inc. is licensed by Caliper Life Sciences, Inc. to sell products using the LabChip technology for research use only. These products are licensed under US Patent Nos. 5,863,753, 5,658,751, 5,436,134, and 5,582,977, and pending patent applications, and related foreign patents, for internal research and development use only in detecting, quantitating, and sizing macromolecules, in combination with microfluidics, where internal research and development use expressly excludes the use of this product for providing medical, diagnostic, or any other testing, analysis, or screening services, or providing clinical information or clinical analysis, in any event in return for compensation by an unrelated party.

Windows and Windows XP are trademarks of Microsoft Corporation.