

- 7 An information window is displayed, indicating that the device driver is being installed (Figure 3).
- 8 An information window is displayed, indicating that the device driver installation is complete (Figure 4).
- 9 An information window is displayed, indicating that you can now use your Experion instrument (Figure 5).

Installation is complete!



Fig. 3.



Fig. 4.



Fig. 5.

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 Vista are trademarks of Microsoft Corporation.