

Profinia™ Software Version 1.0

Catalog #

620-0010

Quantity

1 software CD and 1 USB cable

Description

Profinia software comes with a USB cable to connect your Profinia protein purification system directly to a computer for real-time data monitoring.

Storage

Store at ambient temperature. Keep away from direct heat sources.

Instructions for Use

For real-time data monitoring, connect the USB cable to the USB port located on the back of the Profinia instrument. Connect the opposing end of USB cable to the USB slot on the computer.

Software installation instructions:

1. Place the CD into the CD drive; installation will start automatically. Follow the screen instructions for Profinia software installation.
2. If installation does not run automatically, follow these instructions:
 - Double-click on the My Computer desktop icon
 - Select the CD drive by double-clicking
 - Double-click on the file setup.exe
 - Follow the screen instructions for Profinia software installation



Profinia software