

AFFINITY PURIFICATION

Profinia™ Protein Purification System

Installation Instructions

The Profinia protein purification system is shipped in three boxes. The largest box contains the instrument enclosed in protective foam packaging. The smaller boxes contain the Profinia accessory and starter kits.

1 Profinia System Shipment



2 Remove Straps on Instrument Box

Cut and remove plastic straps. Remove staples (if present).



3 Remove Box Lid

Remove upper lid of box by using side handles.

4 Unpack Instrument

Remove protective foam packaging covering the instrument. The Profinia system instruction manual and quick guide are packed with the instrument.



5 Remove Instrument

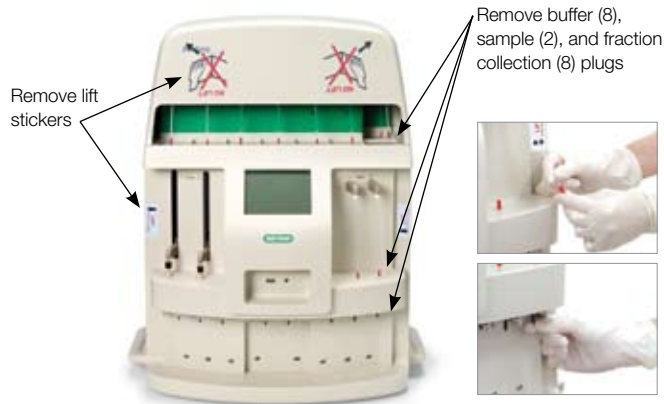
Hold lift handles and bottom of instrument to remove instrument from box.



BIO-RAD

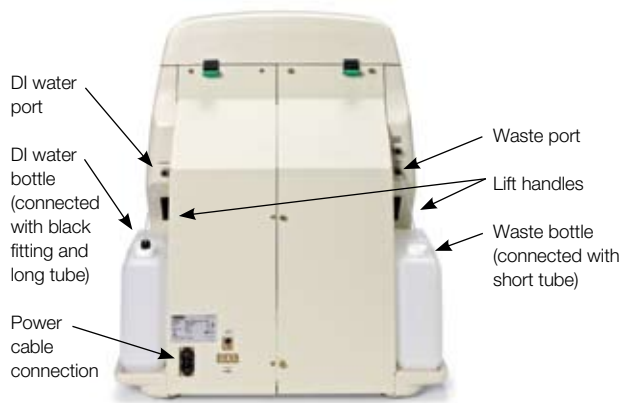
6 Remove Stickers and Plugs

Place instrument on lab bench. Remove plastic bag. Remove four lift instruction stickers from instrument top cover and side panels. Gently remove 18 plugs from tubing (as shown). Store plugs for later use.



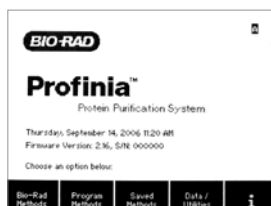
7 Install Water and Waste Bottles and Connect Power Cable

Remove plugs (2) from the deionized (DI) water and waste ports at the back of the instrument. Store plugs for later use. Fill and install the DI water bottle with black fitting and long tube and the waste bottle with white cap and short tube, and connect to the labeled ports above the bottles (as shown). Connect the power cable to the back of the instrument. Components are located in the accessory kit box.



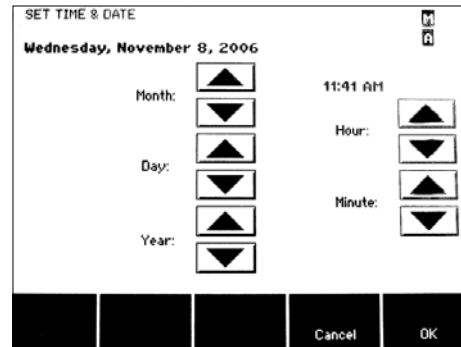
8 Turn Instrument On

Press the power button located on the lower right side of instrument to the on position. The Home screen will display. Locate stylus in accessory box. Touch the Data/Utilities button on the screen.



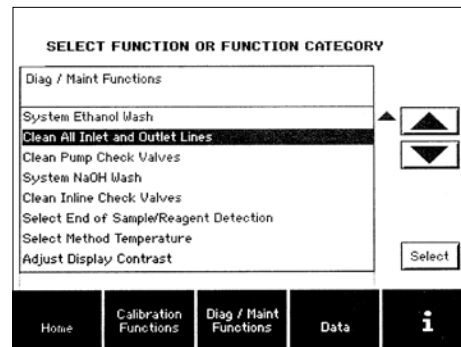
9 Set Time and Date

Touch the Diag/Maint Functions button. Use the arrow keys to highlight the Set Time and Date function. Touch the Select button. Set the time and date, then touch OK.



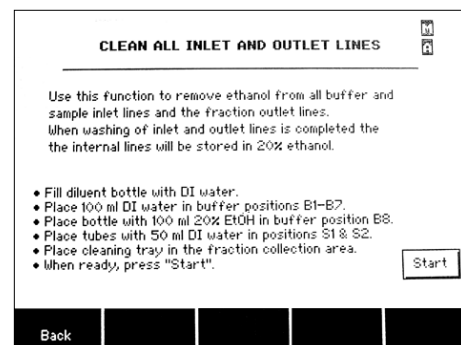
10 Clean All Inlet and Outlet Lines

Use the arrow keys to highlight the Clean All Inlet and Outlet Lines function. Touch the Select button.



11 Perform Cleaning Function Setup

Read screen instructions to guide you through the setup. See steps 12–15 for detailed instructions on completing the procedures for the Clean All Inlet and Outlet lines function.



12 Fill and Place Buffer Bottles

Fill seven buffer bottles (in accessory kit box) with 100 ml of DI water and place in positions 1–7. Fill one buffer bottle with 100 ml of 20% ethanol and place in position 8 (see step 14). Place buffer lids on bottles to prevent dust and other contaminants from entering the system.



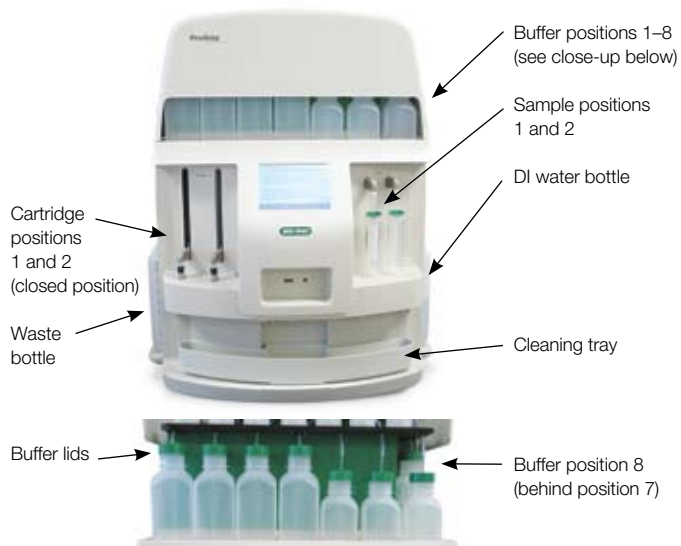
13 Fill and Place Sample Tubes

Fill two conical sample tubes with 50 ml of deionized water. Place lids on tubes, then place in sample positions 1 and 2 (see step 14).



14 Place Cleaning Tray Into Position

Keep cartridge holders in the closed position. Instrument is now ready for cleaning.



15 Touch the Start Button

16 Repeat Clean All Inlet and Outlet Lines

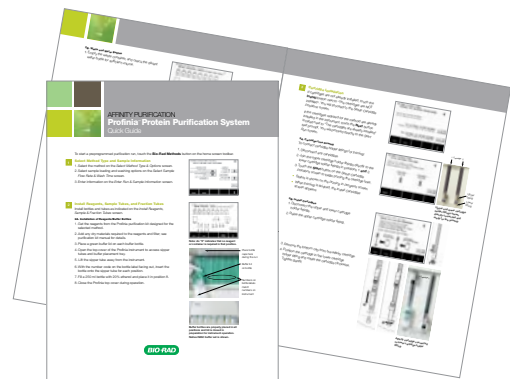
It is recommended that this function be performed a second time prior to first use to remove any air that might have entered the system.

17 Remove Components

Remove cleaning tray, sample tubes, buffer bottles (positions 1-7) before performing a purification run.

18 Perform Purification Run

Use the Profina starter kit to validate instrument. See the Profina system quick guide or refer to section 6 in the instruction manual for instructions on running a Bio-Rad method. (For software installation instructions, see next page.)



1 Open Profinia Software Box



2 Connect USB Cable

For real-time data collection, connect USB cable first to instrument, then to computer, before installing software (USB port is located on the back of the Profinia instrument).



3 Insert Software CD Into Computer



4 Install Software

Follow screen instructions to install Profinia software.



5 Begin Operation

Profinia software is now installed and ready to collect real-time data or view data files from the Profinia system.

Minimum Computer Requirements

- Windows 2000 or Windows XP operating system
- 2 GHz Pentium 4 processor
- 256 MB RAM
- 1,024 x 768 screen resolution
- 40 GB hard drive
- CD-ROM drive
- USB 2.0 port

Pentium is a trademark of Intel Corporation. Windows and Windows XP are trademarks of Microsoft Corporation.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
 Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050
 Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723
 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000