

# Profinia™ pH Monitor Kit

**Catalog #**

620-0411

**Quantity**

Includes 1 x calomel pH probe, 1 x pH flow cell, 1 x tubing assembly, and 1 x mounting bracket and screw kit

**Description**

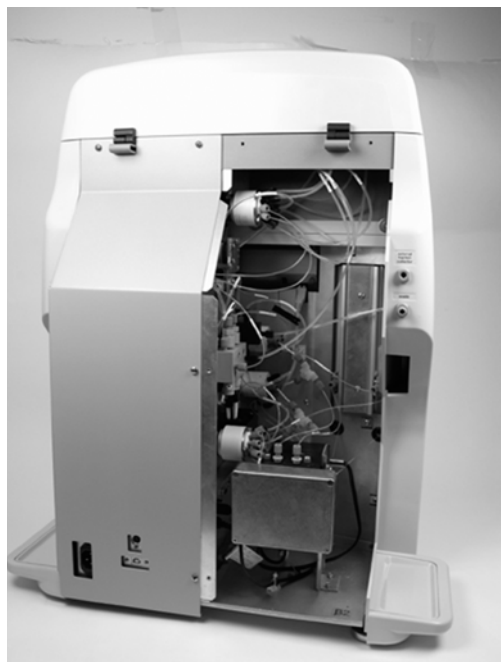
The optional pH monitor kit allows the Profinia instrument to monitor the pH of the fluid stream. The pH information is displayed on the instrument's screen and stored in the run data file that can be viewed on the Profinia software (catalog #620-0010).

**Instructions for Use**

Additional materials required are standard pH calibration buffers, pH 4.0 and 10.0, available from laboratory supply companies.

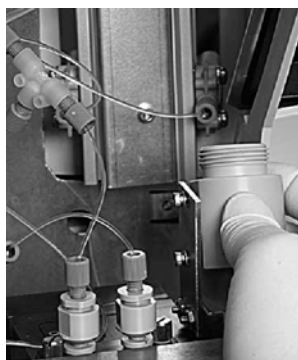
To install the pH flow cell and probe:

1. Use the power switch to turn the instrument off and unplug the power cord.
2. Use the handles on the back of the unit to turn it so the back of the unit is facing you.
3. Disconnect and remove the waste bottle (right side of the instrument).
4. Remove the thumbscrews connecting the rear access panel to the instrument; there are two thumbscrews at the top of the panel, and two near the center.
5. Lift off the access panel by pulling it straight toward you.

**Access panel removed****Bracket mounted in instrument**

pH flow cell  
mounting bracket  
holes and screws

6. Locate the two holes provided for mounting the pH flow cell.
7. Install the pH flow cell bracket using the two screws provided.
8. Find the tubing connecting the conductivity flow cell to the center of the outlet select valve. Unscrew the fitting at the outlet select valve.
9. Connect the fitting you disconnected in step 8 to the inlet of the pH flow cell.
10. Find the length of tubing supplied with the pH monitor kit. Connect one end to the outlet fitting of the pH flow cell.
11. Connect the other end of the tubing to the center port of the outlet select valve.
12. Add a small amount of water (2 ml) to the pH flow cell.
13. Connect the pH probe's electrical connector to the receptacle on the instrument.
14. Remove the plastic cap from the pH probe and install it on the pH flow cell; tighten the black collar until it is snug. If a small amount of liquid drips from the flow cell during this operation, wipe it off with a paper towel.
15. Replace the access panel, four thumbscrews, and the waste bottle, and turn the unit around so it is facing you.
16. Reconnect the waste bottle tube and power and USB cables to prepare the instrument for operation.
17. Turn the instrument power on.
18. On the home screen, touch the Data/Utilities button, then touch the Calibration Functions button.
19. Touch the arrow buttons to scroll through the list of calibration functions and highlight "Calibrate pH", then touch the select button.
20. Follow the instructions displayed on the screen to calibrate the pH probe.



**Installing the flow cell**



**Connecting the tubing**



**Connecting electrical connector**



**Installing flow cell**