

See Technical Manual #TM389 available at [www.promega.com/protocols/](http://www.promega.com/protocols/) for detailed information.



**CAUTION!** The packaged HSM 3.0 Instrument weighs approximately 38.8lb (~17.6kg) and should be handled by two people.



**CAUTION!** Strong magnet. Can be harmful to pacemaker wearers.



**WARNING!** Magnet may interfere with pacemakers and other devices. Pacemaker wearers stay back 30cm (12in).

1. Remove the HSM 3.0 Instrument, Power Supply and accessories from the shipping containers.

**Important:** Save the packaging material in case the equipment needs to be returned for service or repair at a later date.

2. Check that all parts have been included. Refer to the HSM 3.0 Instrument Technical Manual #TM389 for a list of parts.

3. The HSM 3.0 unit is shipped affixed to a wooden support, which must be removed before use. Carefully place the unit upside down and remove the 4 shipping screws with the supplied 6.0mm Hex Wrench (Figure 1). Then insert the Instrument Feet into the threaded holes on the bottom of the instrument.

4. Set the HSM 3.0 Instrument on a flat, level surface in a dust-free location with reasonable air circulation.



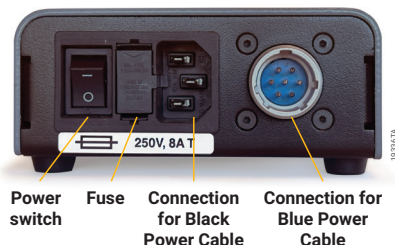
**Figure 1.** Place the instrument upside down on a stable surface. Remove the 4 screws holding the HSM 3.0 Instrument to the plywood.

5. Ensure that the power switch is in the off position. The power switch is located on the back of the Power Supply (Figure 2).
6. Connect the HSM 3.0 Instrument to the Power Supply using the Blue Power Cable. Plug the cable into the back of instrument (Figure 3) and the back of the Power Supply (Figure 2).

- The HSM 3.0 Instrument requires an external computer for operation, either a PC with HSM 3.0 Application Software installed (see Technical Manual #TM389 for instructions) or a PC controlling a liquid handler that has the appropriate software installed. Using the Serial Cable and USB to Serial Converter, attach the HSM 3.0 Instrument to the computer (see Figure 3 for the location of the RS-232 serial port).

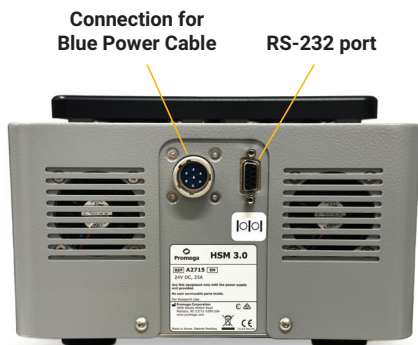
**Important:** Place the Power Supply in a dry location away from possible liquid spills.

- Plug the Black Power Cord into the back of the Power Supply (Figure 2). Plug the Power Cord into a grounded wall outlet.
- Turn on the power switch located on the back of the Power Supply (Figure 2).



**Figure 2.** Rear view of the Power Supply showing the main power switch, fuse and power connections.

- There are three LED lights on the front of the instrument. When the instrument is turned on, the green LED will light up, indicating that the power is on. The instrument will then perform a self-diagnostic test that will take about one minute. During that time you will see and hear the instrument following the diagnostic routine. If an error is detected, the red LED will light up. If there is no red light, the instrument has passed its self-diagnostic test.
- Once all diagnostic checks have passed, the HSM 3.0 Instrument is ready for operation.



**Figure 3.** Rear view of the HSM 3.0 Instrument showing the connections for the Blue Power Cable and the RS-232 Cable port.

### Ordering and Technical Information

www.promega.com • Phone 608-274-4330  
Toll free in USA 800-356-9526 • Fax 608-277-2601



AF9FB134 1224FB134  
© 2024 Promega Corporation.  
All Rights Reserved. Created in USA.  
Revised 12/24. Part #FB134