
HydroTech™ Vacuum Pump

QUICK REFERENCE GUIDE

SETUP AND OPERATION



Warning

Turn the vacuum line's two-way valve to the closed position when the dryer is not in use; the vacuum pump then can continue operating without foaming.

1. **Set the Power On/Off Timer switch for the desired length of time.**
To select "continuous on," turn the knob counter-clockwise.
2. **Set the vacuum line two-way valve to the Closed position.**
This prevents foaming in the reservoir when the vacuum is running but the gel dryer is not sealed for vacuum.
3. **Position the gel in the gel dryer and close the lid.**
4. **Turn the vacuum line's two-way valve to the Open position.**
5. **When the gel is dry,**
 - a. **Lift the transparent Sealing Gasket to break the vacuum seal.**
 - b. **Turn the vacuum line's two-way valve to the Closed position.**
 - c. **Turn off the HydroTech pump.**

MAINTENANCE

The HydroTech pump's reservoir water should be changed at least once a month or more frequently if the water level in the reservoir has reached MAX. Regular reservoir maintenance helps maintain pump performance and reduces foaming. To drain the reservoir:

1. Turn off the Vacuum pump.
2. Connect the drain line to the Reservoir Drain Port.
3. Place the drain line into a large bucket or sink.
4. Turn on the pump. Draining the reservoir requires approximately 30 seconds.
5. Turn off the Pump and disconnect the drain line by pushing the button on top of the reservoir's drain port. The drain port will automatically close.
6. Refill the reservoir with water.