

ECONO GRADIENT PUMP

QUICK START

1) Install pump tubing.

- Tubing selection and installation is described in Section 4.
- For typical applications, ensure the platen adjusting screw is tightened 3 to 5 turns clockwise.

2) Plumb the tubing as needed for the application.

- The right side of the pump is the inlet when the pump is run in the default clockwise direction.

3) Connect peripherals.

- Connect valves and mixer to connectors on the back panel. Refer to Sections 3 and 6.

4) Turn on the Econo Gradient Pump.

- Ensure the correct input voltage is selected (Section 2.2).
- The power switch is underneath the pump.
- When the Econo Gradient Pump is first turned on, the control panel opens up in the Manual mode (Section 5.3).

5) Calibrate the pump flow rate.

- Pump flow rate calibration is done from Screen 2 in Setup mode.
- Set up other parameters in Setup mode.

6) Program a run.

- Press the Mode key until the Program LED is lit. Write a six-step program; one program is saved in non-volatile memory.

7) Start the pump by pressing the Run key.

- Press the Run key to start and stop the pump.
- When the pump is stopped, pressing the Purge key will run the pump at full speed (25 RPM) to rapidly purge air from the tubing.

More sophisticated controls and features are accessed through the Program, Setup and Remote modes.

ECONO GRADIENT PUMP SETUP MODE

