



CHEF Variable Speed Pump

Catalog Number 170-3644

Temperature control is a key factor in achieving high resolution and reproducibility in pulsed field separations. Temperature regulation is closely linked to the buffer flow rate in the electrophoresis chamber. The CHEF Variable Speed Pump gives you a controllable and reproducible buffer flow rate. As a result, CHEF separations are more reproducible from run to run.

The variable speed pump is powered from the CHEF-DR[®] II Drive Module, CHEF-DR II Power Module, or CHEF Mapper[®] Power Module. Plug the pump into the socket marked "Pump" at the front or back of the module. Turn the pump on using the switch marked "Pump Switch". Adjust the flow rate by turning the dial on the variable speed pump. The setting of 100 on the dial is 100% of the pump's maximum flow rate, though the flow is modified by varying line voltage. A setting of 60 on the dial is recommended for most pulsed field electrophoresis experiments with the CHEF system. This corresponds to a flow rate of about 1 liter/minute.

If the pump does not run, check the fuse on the back side, located under the round gray lid. The pump uses a 0.5 Amp SB fuse.

Note: The variable speed pump is ground isolated for safety. It should not be plugged into a wall socket.