

# Installation Instructions for Redesigned CHEF Electrophoresis Cell

The CHEF Electrophoresis Cell has been redesigned with improved features. The installation instructions for the new cell may not have been incorporated into the system instruction manuals for the CHEF Mapper<sup>®</sup>, CHEF-DR<sup>®</sup> III, CHEF-DR II, or GenePath<sup>®</sup> system at the time this cell was introduced. Please follow these instructions for proper connection of your new cell to the other components of CHEF systems.

To connect the cell to the Cooling Module (may be optional), attach approximately 1–2 feet of 1/4 inch internal diameter Tygon tubing to both the Flow In and Flow Out ports on the Cooling Module, and secure the tubing with plastic clamps (provided). Connect a quick release connector (provided) to 2 feet of 3/8 inch internal diameter Tygon tubing. Attach the quick release connector to the left front port of the cell. Attach the other end of the 3/8 inch tubing to the 1/4 inch tubing from the Flow In of the Cooling Module using 3/8 to 1/4 inch reducer (provided). The pump should be positioned between the outlet of the Cooling Module

and the inlet (rear) of the electrophoresis cell. Connect the 1/4 inch tubing from the Flow Out of the Cooling Module to the inlet of the pump using a 3/8 to 1/4 inch reducer. Connect approximately two feet of 3/8 inch Tygon tubing to the outlet of the pump using a 3/8 to 3/8 straight connector (provided). Connect a quick release connector to the other end of the 3/8 inch tubing. Connect the quick release connector to the inlet of the cell.

Connect a quick release connector to a six inch of 3/8 inch Tygon tubing, and connect it to the right front port of the cell. This tube will be used to drain the buffer in the electrophoresis cell.

Connect the 9 pin gray temperature probe cable (included) from the back of the electrophoresis cell to the Remote Sensor port on the back of the Cooling Module.

Refer to the system instruction manuals for connection of the electrophoresis cell to various CHEF power modules.

