



Dear Western Processor owner,

Please find enclosed a set of tubing that will prolong the life and improve the performance of your Western Processor. The tubing is installed on the inlet of each reagent valve. Remove the tubing that draws from the reagents into the valve. Place one piece of each new tubing on each valve connecting the end marked with the purple tape collar to the valve.

The increased diameter of the tubing reduces the vacuum drawn over the valve, which reduces stress on the valve, and improves ease of calibration by equalizing the performance of all six valves.

Bio-Rad is committed to the continuing improvement of our products and wanted to share this simple change in order to improve your long term satisfaction with this product.

***Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules, Ca 94553***