

## **Instructions for Mounting the Fan Control Kit for the BioFocus 3000**

1. The kit consists of:
  - 1 Fan Control Assembly, 800-5692
  - 1 Cable, 800-5693
  - 2 Mounting screws, 8-32 x 1/2 in. with captive lockwasher
  - 1 Cable tie
2. Turn off the BioFocus 3000 and remove the power connector.
3. Remove the top cover and the front lower door (to allow removal of the left side cover).
4. Remove the left side cover (the one opposite the reservoir).
5. Remove the back panel.
6. Remove the two screws below and to the sides of the exhaust fan which hold the TCU to the floor of the instrument.
7. Slip the Fan Control Assembly behind the exhaust fan.
8. Use the two mounting screws to install the Fan Control Assembly where the screws were removed in step 6.
9. On the left side of the instrument, unplug the waste heat fan cable to P129 in the middle of the TCU PCB.
10. Slide the 28 Volt power supply out a little. Carefully cut the cable tie which holds the waste heat fan cable to the other cables and remove the cable tie. Connect J129 to the Fan Control Assembly at the small connector.
11. Connect the new Cable, 800-5693 from the large connector on the Fan Control Assembly to P129 on the TCU PCB. Put a new cable tie at the place where the old one was removed.
12. Unscrew TB1-7. Slip the white wire from the cable into the terminal with the red wire, and retighten.
13. Reinstall all covers and test the operation. The fan control will react slowly, taking about 2 seconds to come to full speed when the sample area temperature control is turned on, and several minutes to react fully to a change in the temperature.