

Replacing the Dual Plasma Controller Printed Circuit Board

Accessory G6600-67000

Purpose

To replace the printed circuit board (PCB) in a Dual Plasma Controller.

Scope

This procedure describes how to replace the PCB in a Dual Plasma Controller for use with either the 355 SCD or the 255 NCD.

Safety

Make sure all devices are powered off while performing the installation.

Follow ESD precautions while performing the procedure.

Materials

- A Dual Plasma Controller printed circuit board kit
- The Dual Plasma Controller
- A Philips screwdriver

Procedure

1. Make sure all devices are powered off (Controller, pump, and GC).
2. Remove all connections from the back of the Controller.
3. Take ESD precautions; put on an ESD wrist strap before proceeding.
4. Remove the cover from the Controller.
5. Remove all electrical connections (8 total) on the PCB (see **Figure 1**). Make note of where each quick connect attaches to the PCB, as these connections must be properly re-established in the new PCB later.



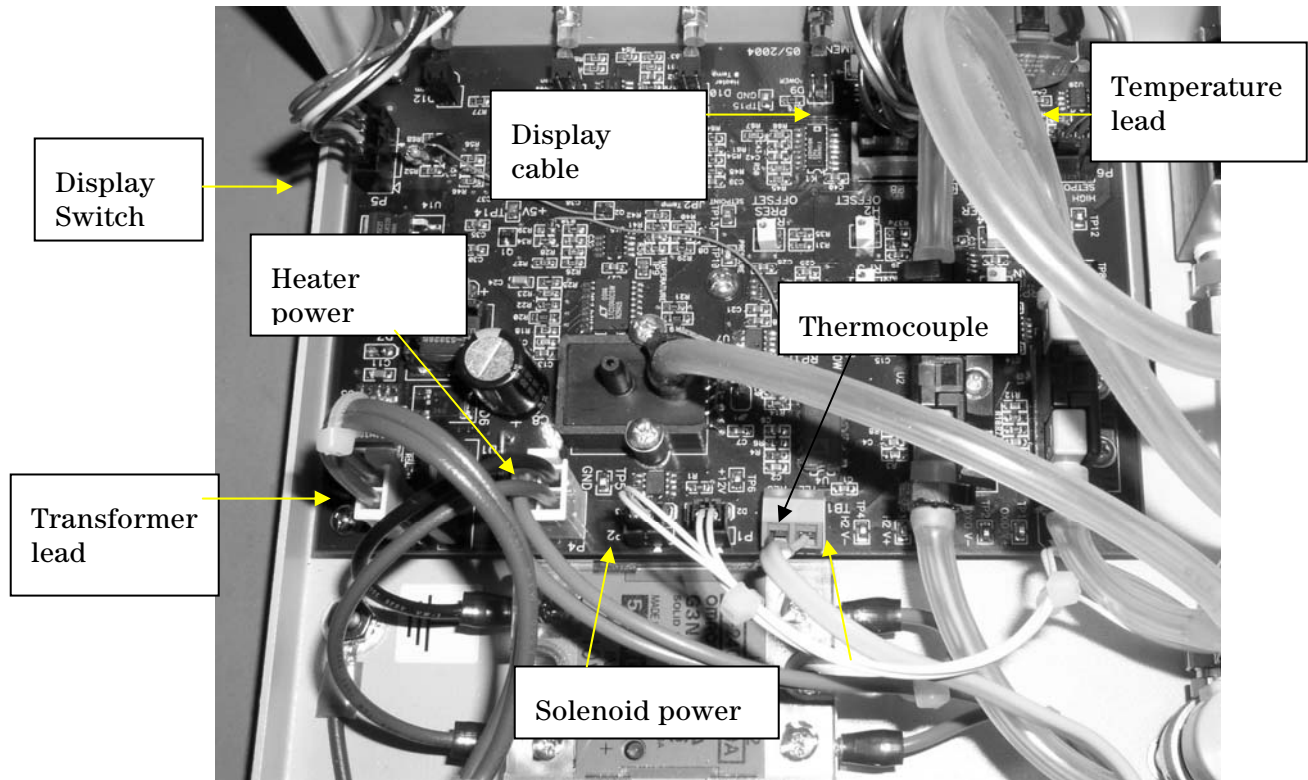


Figure 1 Electrical connections on the PCB

6. Remove all fluidics connections (5 total) from the PCB by loosening the hose clamp and gently pulling the hose free of the barb. Make note of where each line attaches; you may want to attach a small piece of tape to the tubing to identify the connections.

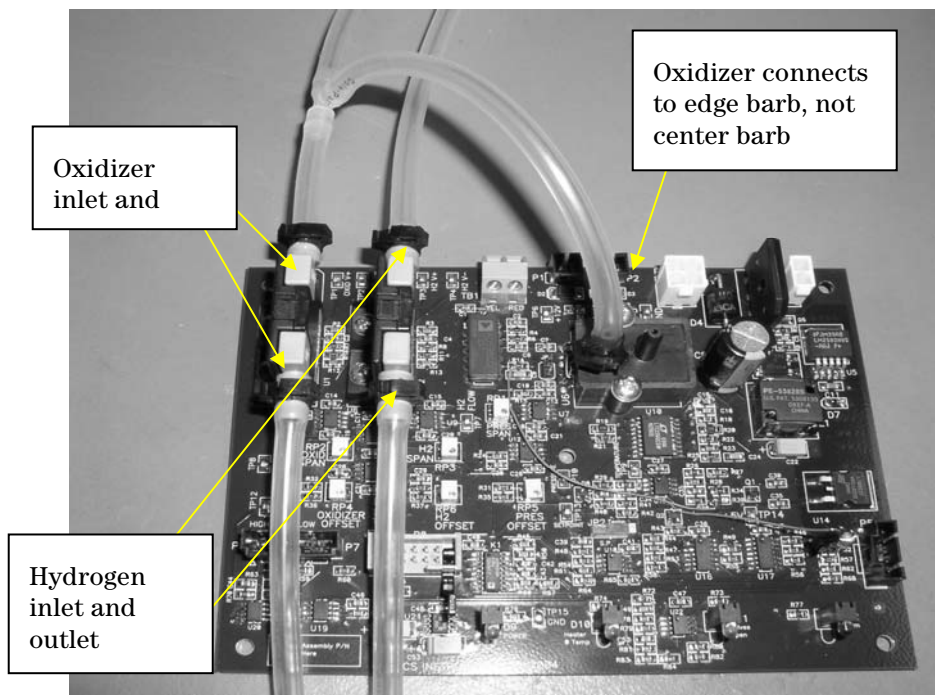


Figure 2 Fluidics connections on the PCB

7. Remove the PCB from the Controller by loosening the five screws (one at each corner and one in center) and pulling the board free.
8. Remove the new PCB from the packaging.
9. Attach the new PCB to the bottom of the case by securing the five screws.
10. Reconnect each of the five fluidics connections to the appropriate barb and reattach the hose clamp.
11. Reconnect the electrical connections to the PCB.
12. Replace the Controller cover and secure properly.
13. Restore all connections to the back of the Controller.

Copyright © 2007
Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610
Printed in the USA Feb 2007



G6600-90033