

GeneAmp[®] PCR System 2400 Coolant Replacement Procedure

The GeneAmp[®] PCR System 2400 Thermal Cycler requires routine maintenance which can be performed by the end user. This routine maintenance will improve the performance of the instrument and may prevent more costly in-house repairs. The procedure described here is for the periodic checking of the coolant in the instrument and the procedure for re-filling the coolant, should the need arise.

Applied Biosystems service personnel can perform this procedure should you be reluctant to do so. This service requires the instrument to be shipped to our repair facility and the user would be charged all fees associated with this procedure. To request a service repair of your instrument please call our Service Call Center at 1-800-831-6844.

Applied Biosystems is not responsible for any damage to the machine, or its parts, during the performance of this procedure.

GeneAmp[®] PCR System 2400 Coolant Service Procedure



Instructions for checking the coolant level, cleaning the coolant reservoir filter, and re-filling the coolant reservoir for the GeneAmp[®] PCR System 2400 thermal cycler.

Step 1. Preparing the instrument:

Turn off and unplug the instrument. Unscrew and remove the four screws located on the sides of the instrument. Two screws are on the right side and two screws are on the left side.



Side Screws

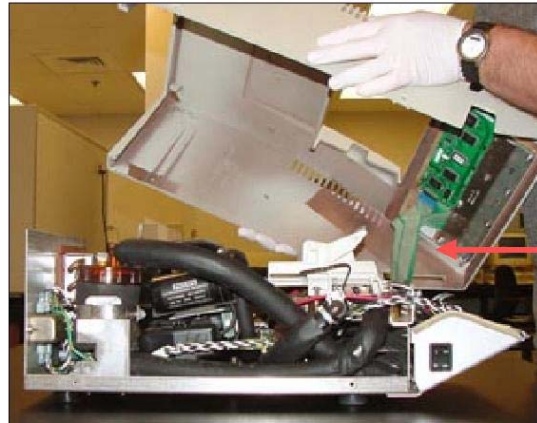
Step 2. Opening the instrument cover:

Carefully remove the cover of the instrument by lifting straight up.

NOTE: Use caution when lifting the cover. A ribbon cable is connected to the lid and the base of the instrument. Damage to the cable can result in the costly replacement of parts.

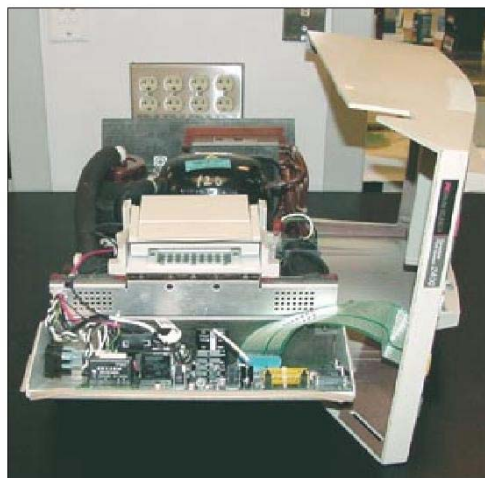


After the cover has cleared the instrument base, start to bring the cover to the side.



**Ribbon
Cable**

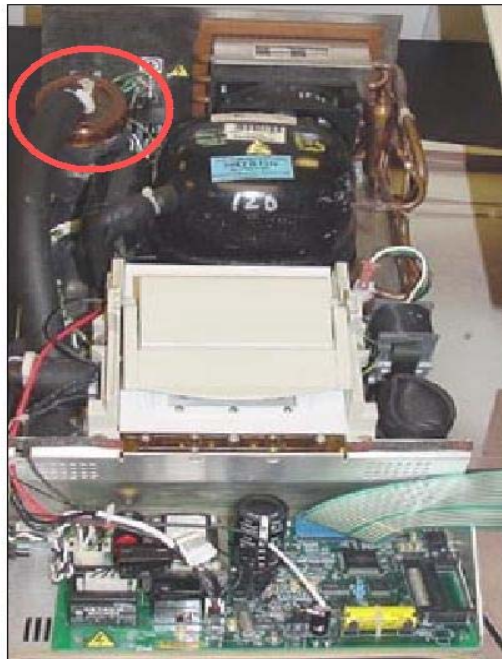
The ribbon cable is long enough that the cover can be placed on the benchtop next to the instrument.



Step 3. Locating the coolant reservoir:

After the cover is removed locate the coolant reservoir. The coolant reservoir is located in the back left-hand corner of the instrument.

**Coolant
Reservoir**

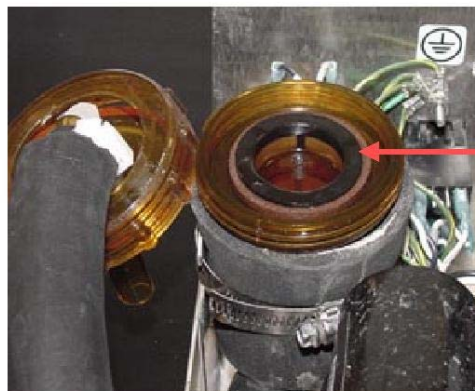


Step 4. Opening the coolant reservoir:

Unscrew the coolant reservoir lid.



Check the coolant level – the coolant reservoir should be full with coolant. If the coolant level is low please continue to step 5. If the coolant reservoir is full, please skip to step 6.



**Full
Line**

Step 5. Filling the coolant reservoir:

If the coolant level is low, prepare approximately 700 ml of 50% distilled water and 50% Prestone® 5/150 Extended Life Antifreeze. Pour the coolant (distilled water/antifreeze mixture) into the coolant reservoir until the reservoir is full. Properly dispose of the excess coolant.



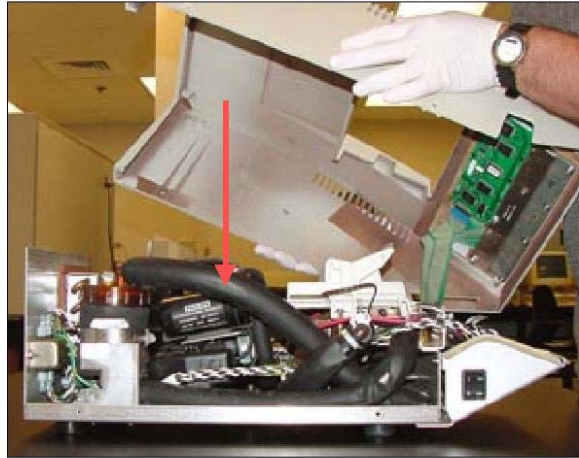
Step 6. Cleaning the coolant reservoir filter:

Pull out the basket filter located inside the coolant reservoir. Clean the filter with a scrub brush and distilled water. After cleaning, place the filter back in the coolant reservoir and screw on the coolant reservoir.



Step 7. Re-Assembling the Instrument:

When finished with previous steps, carefully place the cover back on- Please be careful not to crimp the ribbon cable. Replace the 4 screws removed in Step 1.



For Research Use Only. Not for use in diagnostic procedures.

Applied Biosystems is a registered trademark of Applied Biosystems Corporation or its subsidiaries in the U. S. and certain other countries.

GeneAmp is a registered trademark of Roche Molecular Systems, Inc.

All other trademarks are the sole property of their respective owners.