

# **Applied Biosystems PCR System 9700, Firmware v3.11 Release Notes**

## **Contents:**

1. Overview
2. Features
3. Downloading Firmware
4. Upgrading Firmware
5. Known Issues
6. Legal section

---

### **1. Overview**

#### **Firmware v3.11 for GeneAmp® PCR System 9700**

The GeneAmp® PCR System 9700 delivers the proven performance and reliability of earlier generation Applied Biosystems GeneAmp® PCR Systems.

New features in this firmware include:

- Support for replacement on-board battery
- Improved control for Aluminum Blocks

### **2. Features**

No new features were added in this firmware release.

### **3. Downloading Firmware:**

Please visit Applied Biosystems' support website to download the latest version of the firmware.

<http://www.appliedbiosystems.com/support/software/>

### **4. Upgrading Firmware:**

Using the N8051327 cable, there are settings that need to be done when upgrading using the HyperTerminal program in Windows XP.

#### **Part 1: Configurations of the GeneAmp 9700**

Step 1: On the control panel select "UTIL"

Step 2: select "DIAG"

Step 3: select "STOP" & "F3" buttons at the same time, which will lead user into the hidden menu.

Step 4: select "PORT MODE"  
Step 5: ensure that "TERM" is the selected mode  
Step 6: exit to main menu  
Step 7: select "UTIL"  
Step 8: select "CONFIG"  
Step 9: select "MORE"  
Step 10: select "BAUD RATE", note down the figure that is being displayed. This is the baud rate setting which we will be using for the HyperTerminal program later.

(The baud rate setting is determined by the block type and the baud rate setting on the HyperTerminal program must be of the same setting as that displayed on the 9700 unit)

Step 11: Select "More" twice, and ensure that "Get IP" is set to OFF  
Step 12: Select "Accept"  
Step 13: Select "Diag"  
Step 14: Select "Upgrad"  
Step 15: Select "Serial"  
Step 16: Select "Update"

## **Part 2: On your computer: Settings of HyperTerminal**

Note: HyperTerminal can be found in :

Start/Accessories/Communication/HyperTerminal in Windows XP.

Step 1: Name File, Select COM1 or COM2, depending on which port the cable is attached  
Set the "Bits per second" to the same setting as displayed in "BAUD RATE" found on the 9700 unit.  
Step 2: Ensure that the rest of port settings is configured (refer to Figure 1), Select Apply  
Step 3: Under File, Select Properties, Select Settings, Select "ASCII Setup"  
Step 4: Select both of the "ASCII Sending" settings (refer to Figure 2)  
Step 5: Under Transfer, Select Send File  
Step 6: In the Filename field, locate the upgrade file with the name under which you saved the file  
Step 7: In the protocol field, select Kermit, Select Send

Figure 1

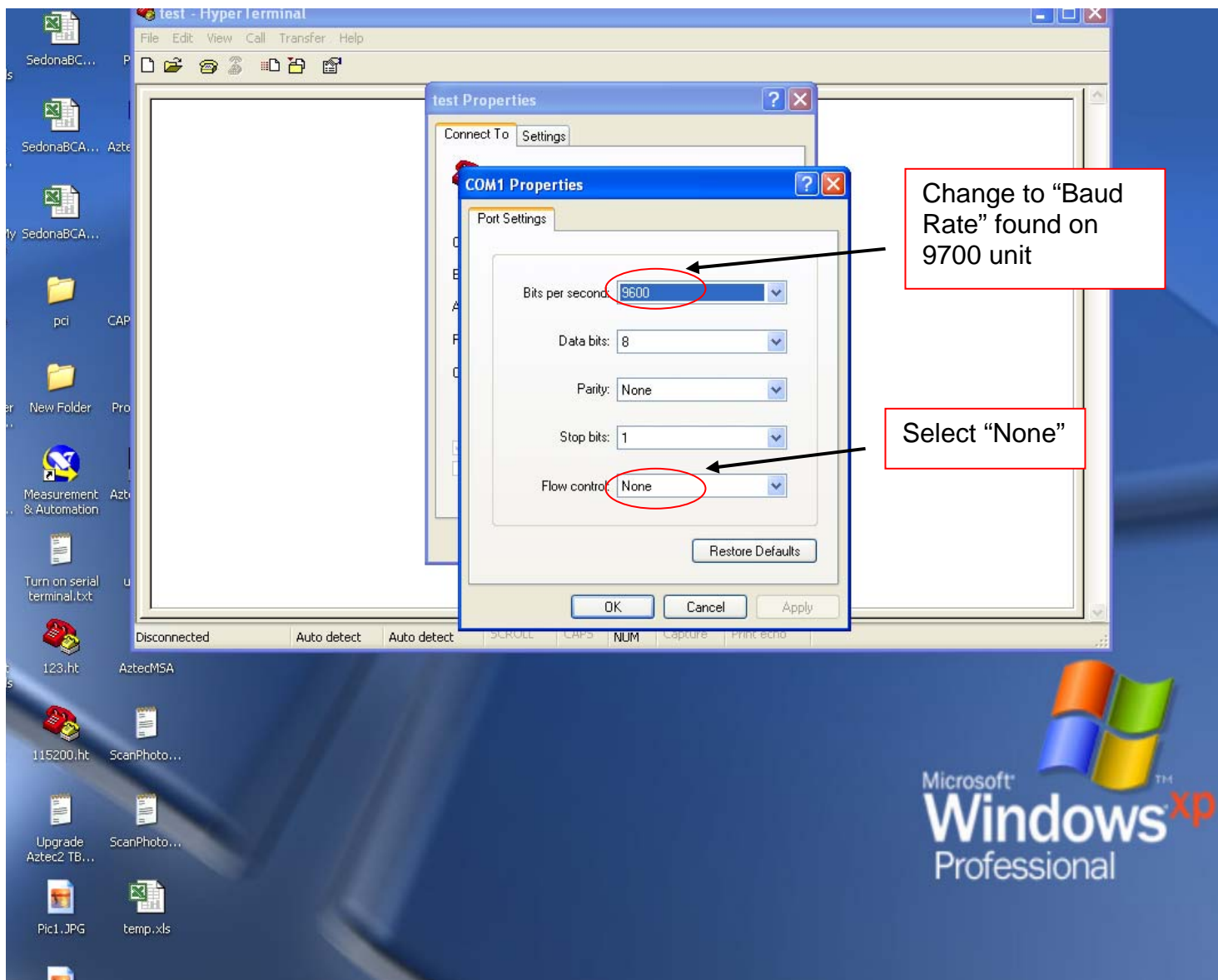
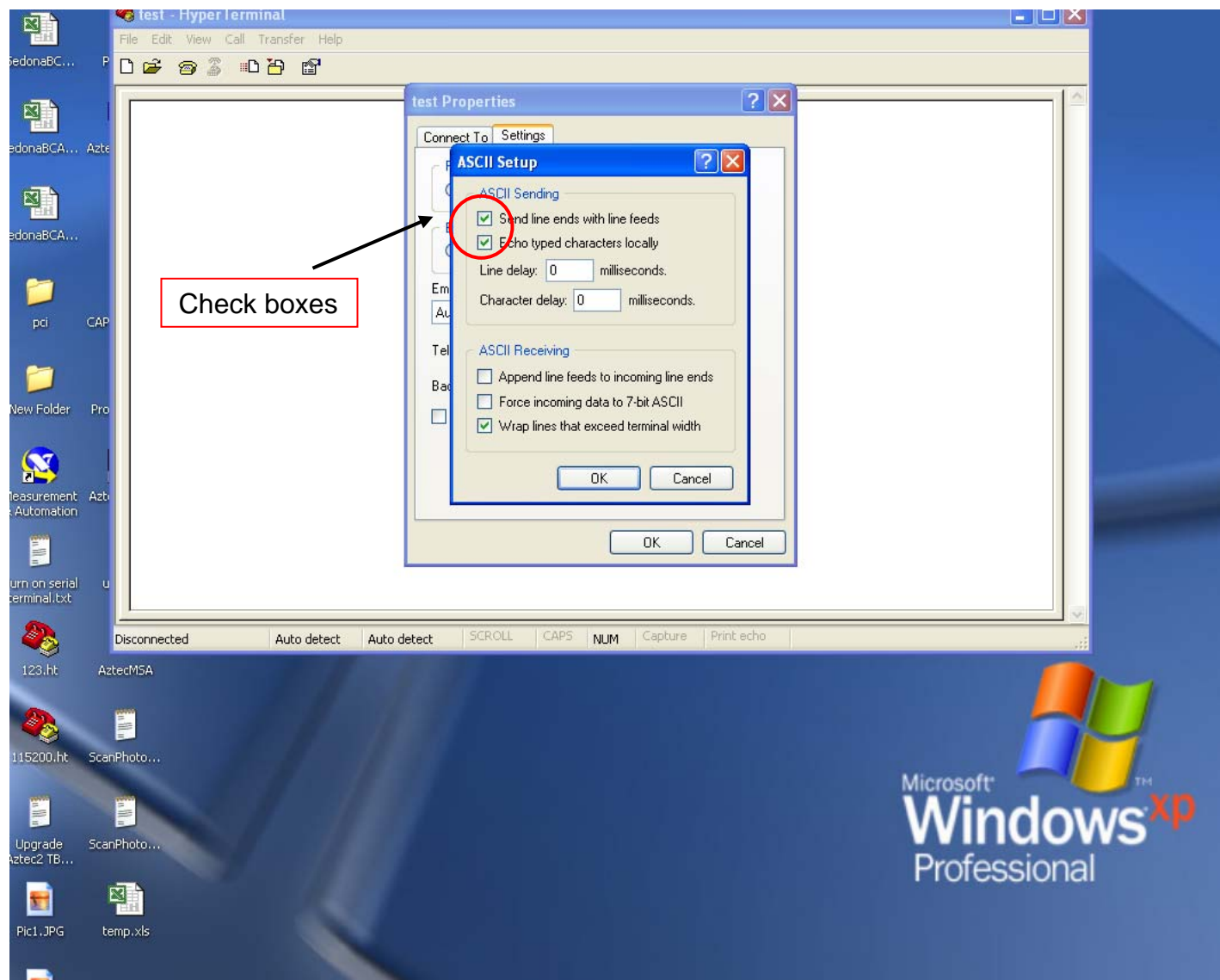


Figure 2



## 5. Known Issues

- Run does not return to exact point of the protocol at which the power loss occurred.

## 6. Legal Section

NOTICE TO PURCHASER: Label License

The GeneAmp PCR System 9700 is covered by patents owned by Applied Biosystems. No right is conveyed expressly, by implication, or by estoppel to apparatus, reagents, kits, or methods such as the patented 5' nuclease methods. For further information contact the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Applied Biosystems, AB (Design), and GeneAmp are registered trademarks of Applied Biosystems or its subsidiaries in the US and/or certain other countries.

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

**© Copyright 2009 Applied Biosystems. All Rights Reserved.**