

Clearing User-Protected Stored Files from the Applied Biosystems GeneAmp® PCR Thermal Cyclers

The GeneAmp® PCR System 2400, GeneAmp® PCR System 9600, GeneAmp® PCR System 9700, and GeneAmp® PCR System 2700 have the option to protect methods using a customer selected PIN number. With time, users may leave the laboratory leaving many of these programs protected by unknown PIN numbers. This document will provide instructions for clearing user-protected files.

Warning: Clearing user-protected files will permanently remove the file.

Clearing User Protected files from the GeneAmp® PCR System 2400, GeneAmp® PCR System 9700, and GeneAmp® PCR System 2700 systems

In order to delete user protected files, the customer selected PIN number must be known. However, if this original PIN number is not known, the GeneAmp® PCR System 2400, GeneAmp® PCR System 9700, and GeneAmp® PCR System 2700 have a back-up PIN number that can be used. This back-up PIN number is the date in which you are using this PIN written in the format of **mm/dd**. For example, if today is July 4th, the PIN number would be 0704. When prompted for the PIN number, enter this back-up PIN.

Clearing User Protected files from the GeneAmp® PCR System 9600

The following instructions will allow you to delete protected files from the GeneAmp® PCR System 9600.

Using the touch pad on the GeneAmp® PCR System 9600 return to the **Main Menu** by pressing **STOP** either one or two times

Select Option 9600

RN-CREATE-EDIT-UTIL

Move the Cursor to **UTIL** by pressing the **OPTION** key

Select Function

DIR-CONFIG-DIAG-DEL

The cursor will be under the **DIR** Function
Press **Enter**

Directory

PROGRAM-TYPE-USER-PRINT

Press the following keys in the following order (do not be concerned with the display):

"MORE 1,2,3"

DO NOT PRESS ENTER

A complete program directory, including user numbers, is now accessible. Use the **STEP** key to view each program beginning with #1. For those programs that have a user's number, the number will be listed on the display

Write down each user number as you progress through the list of programs

Return to the **Main Menu** by pressing **STOP** either one or two times

Select Option 9600

RN-CREATE-EDIT-**UTIL**

Move the cursor to **UTIL** by pressing the **OPTION** key

Select Function

DIR-CONFIG-DIAG-DEL

When the Utilities Menu appears, move the cursor to **DEL** by pressing the **OPTION** key

The Delete Function Menu offers three choices

Delete

PROGRAM-**USER**-ALL

Move the cursor to **USER** by pressing the **OPTION** key. Press **ENTER**. Enter the user numbers of the user-protected programs that are to be deleted. Press **ENTER** after each entry

Programs that are not user-protected can simply be deleted as a group by then moving the cursor to **ALL** and pressing **ENTER**

Programs that are not user-protected can be deleted individually by moving the cursor to **PROGRAM** and pressing **ENTER**. As prompted, enter the number of the individual programs to be deleted.

To Verify that the programs have been deleted, return to the **Main Menu**. Move the cursor to **UTIL**. Press **ENTER**. The cursor should be under **DIR**. Press **ENTER**. Move the cursor to **USER** and press **ENTER**.

A message will indicate that there are no files in the directory.

104GU18-01

4371086 RevB