

Re-entering Lost Calibration Values on the GeneAmp® PCR System 2700

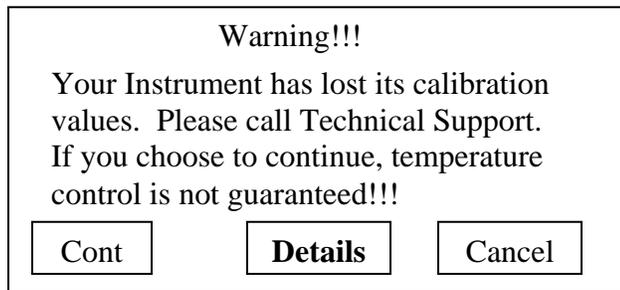
Calibration values can occasionally be lost from the GeneAmp® PCR System 2700 due to a variety of factors. If the calibration values have been lost, an error message will appear when attempting a run.

This document will explain the steps required to obtain and re-enter the lost calibration values for the GeneAmp® PCR System 2700.



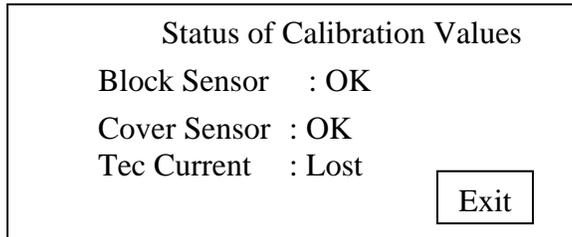
Directions for entering all the calibration values into a GeneAmp® PCR System 2700.

When calibration values have been lost an error message will be displayed as follows:



F1 F2 **F3** F4 F5

To obtain and re-enter calibration values, perform the following procedure. Go to the "Details" display by pressing the **F3** key to determine which calibration values have been lost.



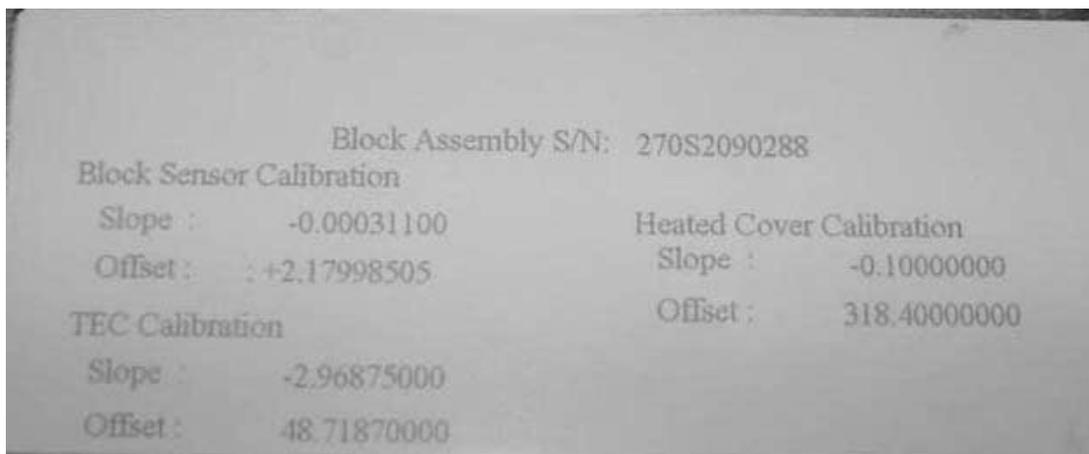
F1 F2 F3 F4 F5

The three types of calibration values that can be lost are:

- [Block Sensor values](#)
- [Cover Calibration values](#)
- [Tec Current values](#)

In the example above, the TEC current values have been lost. First, obtain your instrument serial number. The serial number of the instrument is located on the back of the instrument near the power on/off switch.

GeneAmp® PCR System 2700s with a serial number starting at 270S1102903 have the calibration values located on the underside of the machine. Write down these values, which contain a number for the **slope** and **offset** for each type of calibration value.



If the values on the calibration sticker are expressed in exponential form (Ex. 2.075410e-4), convert them to decimal numbers to enter into the instrument.

(Ex. 2.075410e-4 = .0002075410)

When entering calibration values, the instrument is limited to 7 characters. If your calibration value has more than 7 digits, enter only the first seven. In this example, you would enter 0002075

For instruments with a serial number below 270S1102903, the calibration values are **not** located on the instrument and you will need to contact your local Applied Biosystems support group to obtain these values.

For customers in the United States and Canada, please send and email to technical support at <http://fags.appliedbiosystems.com>. Please include in the email:

- the serial number of the machine
- your name
- phone number
- fax number
- email address
- your address

Once this information is received, your calibration values will be resent to you via e-mail or Fax.

For customers outside of the U.S. and Canada, please contact your local support group. You may find the contact information at: <http://www.appliedbiosystems.com/support/contact/internationalsupport.cfm>

Once the calibration values have been obtained, continue with the following pages for directions on entering the calibration values.

Entering Block Sensor Calibration Values on the GeneAmp® PCR System 2700

Press F4(Util)

```

08:00 AM      09/25/00  25.0°C
GeneAmp® PCR System 2700
Version
User: adrian
Run  Create  Edit  Util  User
F1    F2    F3    F4    F5
    
```

Press F1(Diag)

```

Utilities
Diag   - Instrument diagnostics
TmCalc - Calculate melting temp
Config - Instrument configuration
Diag  TmCalc Config More Exit
F1    F2    F3    F4    F5
    
```

Hold down the Stop key and press F3 (TempVer)

```

Diagnostics
Hard   - Hardware Diagnostics
System - System Performance Tests
TmpVer - Temperature Verification
Upgrad - Firmware Upgrade
Hard  System TempVer Util User
F1    F2    F3    F4    F5
    
```

Press the down arrow key twice to highlight Calibrations, then press F1(Run)

```

Service Functions
Service Info  Block Params  Final Test
Base Serial#  Manual Ctrl   Manual TEC
Calibrations  Port Mode     Other Diags
Init BRAM     Stability     Log
Run           User
F1    F2    F3    F4    F5
    
```

Press F2 to get to the block sensor calibration page and enter the slope and offset

```

Calibrations
Elect - Electronic Calibration
Block - Block sensor calibration
Tec I - Amp. Cur. Calibration
Print - Print Tec I Slope/Offset
Elect Block Tec I Print Exit
F1    F2    F3    F4    F5
    
```

```

Block Sensor Calibration
Resist  Temp °C  Slope  Offset
Xxxx   xx.xx  -x.xxxxxx +x.xxxxxx
Accept Calib +/- Default Exit
F1    F2    F3    F4    F5
    
```

Enter the new slope and offset. When complete, press F1 (Accept). Press F5 (Exit) **four** times to return to the main menu

Entering Cover Calibration Values on the GeneAmp® PCR System 2700

Press F4(Util)

```

08:00 AM      09/25/00  25.0°C
GeneAmp® PCR System 2700
Version
User: adrian
Run  Create  Edit  Util  User
F1   F2   F3   F4   F5
    
```

Press F1 (Diag)

```

Utilities
Diag   - Instrument diagnostics
TmCalc - Calculate melting temp
Config - Instrument configuration
Diag  TmCalc Config More Exit
F1    F2    F3    F4    F5
    
```

Hold down the Stop key and press F3 (TempVer)

```

Diagnostics
Hard   - Hardware Diagnostics
System - System Performance Tests
TmpVer - Temperature Verification
Upgrad - Firmware Upgrade
Hard  System TempVer Util User
F1    F2    F3    F4    F5
    
```

Press the right arrow key once to highlight Block Params, then press F1(Run)

```

Service Functions
Service Info  Block Params  Final Test
Base Serial#  Manual Ctrl   Manual TEC
Calibrations  Port Mode     Other Diags
Init BRAM     Stability      Log
Run           User
F1    F2    F3    F4    F5
    
```

Press F4 (Next) twice to Cover Calibrations

```

Block Parameter
Block Type : 0
Block F/W version : 53
Block assembly S/N : 096
Block is configured : yes
Accept Prev Next Cancel
F1    F2    F3    F4    F5
    
```

```

Cover Calibration
Slope : -x.xxxxxx
Offset : +xxx.xxxx
Accept Prev Print Next Cancel
F1    F2    F3    F4    F5
    
```

Enter the new calibration values. When complete, press F1 (Accept). Press F5 (Exit) **three** times to return to the main menu

Entering Tec Current Calibration Values on the GeneAmp® PCR System 2700

Press F4(Util)

```

08:00 AM      09/25/00  25.0°C
GeneAmp® PCR System 2700
Version
User: adrian
Run  Create  Edit  Util  User
F1   F2   F3   F4   F5
    
```

Press F1 (Diag)

```

Utilities
Diag  - Instrument diagnostics
TmCalc - Calculate melting temp
Config - Instrument configuration
Diag  TmCalc  Config  More  Exit
F1    F2    F3    F4    F5
    
```

Hold down the Stop key and press F3 (TempVer)

```

Diagnostics
Hard  - Hardware Diagnostics
System - System Performance Tests
TmpVer - Temperature Verification
Upgrad - Firmware Upgrade
Hard  System  TempVer  Util  User
F1    F2    F3    F4    F5
    
```

Press the right arrow key once to highlight Block Params, then press F1(Run)

```

Service Functions
Service Info  Block Params  Final Test
Base Serial#  Manual Ctrl   Manual TEC
Calibrations  Port Mode     Other Diags
Init BRAM     Stability      Log
Run           User
F1    F2    F3    F4    F5
    
```

Press F4 (Next) **four** times to Tec Current Calibrations

```

Block Parameter
Block Type : 0
Block F/W version : 53
Block assembly S/N : 096
Block is configured : yes
Accept  Prev           Next  Cancel
F1    F2    F3    F4    F5
    
```

```

Tec Current Calibration
Slope : -x.xxxxxx
Offset : +xx.xxxx
Accept  Prev  Print  Next  Cancel
F1    F2    F3    F4    F5
    
```

Enter the new calibration values. When complete, press F1 (Accept). Press F5 (Exit) **three** times to return to the main menu

104GU11-01

4370997 Revision B