TmTool[™] Application User Guide



Table of Contents

P	REFAC	CE	2
1	OVI	ERVIEW OF THE ANALYSIS TOOL	2
2	НО	W TO OBTAIN THE TOOL	2
3	INS	TALL THE MATLAB RUNTIME ENVIRONMENT	2
4	WO	RKING WITH TMTOOL TM	4
	4.1	STARTING THE TOOL	4
	4.2	IMPORTING A SOURCE FILE FOR ANALYSIS	6
	4.3	SELECTING THE INSTRUMENT TYPE	7
	4.4	IDENTIFYING REFERENCE WELLS	8
	4.5	IDENTIFYING WELLS TO OMIT FROM ANALYSIS	. 10
	4.6	REGION OF ANALYSIS: DEFAULT SELECTION	. 11
	4.7	CHOOSING A LOCATION TO STORE THE RESULTS	.12
	4.8	SELECTING PLOT OPTIONS	.12
	4.9	CALCULATE TM	. 13
	4.10	REVIEWING THE RESULTS	. 13
	4.11	DEFINING CUSTOM REGION OF ANALYSIS FOR WELLS WHERE NECESSARY	. 18
	4.12	Designating Reference Wells, generating $\Delta T m$ and Plots with References .	.20
5	TRO	DUBLESHOOTING	. 22
6	DEF	FINITIONS	. 26
A	PPEND	DIX	.27

Preface

The TmToolTM application analyzes fluorescent readings from melt experiments performed using Applied Biosystems Real-time PCR Instruments¹ and generates a table of the melting temperatures based on the Boltzmann equation. The tool is designed for melt profiles where the fluorescence readings rise with temperature.

The tool is a suite of scripts written in MATLAB and compiled into an executable. MATLAB is not required to run the tool, but MATLAB runtime environment is required.

The TmToolTM application was developed and tested on a Windows® XP 32 bit system with Service Pack 3. A minimum of 10G available disk space, 1G RAM available memory and a 2.4GHz CPU is recommended. TmToolTM application works best at a screen resolution of 1280 by 800 pixels.

Text conventions:

- • Note Indicates information that may be of interest or help but is not critical to the use of the TmTool.
- Courier font Indicates code or command line code.
- Italicized text -- Used for emphasis.

1 Overview of the analysis tool

The TmToolTM application directly imports instrument run files from the StepOnePlusTM and the 7500 Fast Applied Biosystems Real-time PCR platforms. It can use text exports of multi-component data from the 7900 HT systems.

The TmToolTM application enables you to:

- a. calculate a melting temperature (Tm) from their melt curve data using the Boltzmann equation,
- b. designate reference samples,
- c. compare samples to the reference samples by computing a Δ Tm, defined as a shift in Tm value of their sample from the average Tm of reference samples,
- d. visualize the melt curve of a sample with or without a reference curve in a picture format.

2 How to obtain the tool

Contact your local Life Technologies support representative for more information.

3 Install the MATLAB runtime environment

The following files are supplied as part of the installation:

MCRInstaller.exe TmToolv1_0.ctf TmToolv1_0.exe UserGuidev1_0.pdf

¹ For further information on experiment run protocols, refer to TmToolTM Quick Set up Guide.

QuickSetupGuide.pdf

?Note the installation does *not* create a shortcut to the $TmTool^{TM}$ application on the desktop.

- Double click and install the MCRInstaller.exe to set up the MATLAB run time environment. The installer prompts you to choose the installation location. The default location is: C:\Program Files\MATLAB\MATLAB Component Runtime\
- 2. Add the following to the environment PATH variable:
 - a. From the desktop, right-click "My Computer", then select "Properties".
 - b. In the "System Properties" window, select the "Advanced" tab then click "Environment Variables".

System Properties	Environment Variables	X
System Restore Automatic Updates Remote General Computer Name Hardware Advanced	User variables for majumdn1	
	Variable Value	
You must be logged on as an Administrator to make most of these changes.	TEMP %USERPROFILE%\Local Settings\Temp	
Performance	TMP %USERPROFILE%\Local Settings\Temp	
Visual effects, processor scheduling, memory usage, and virtual memory		
Settings	New Edit Delete]
User Profiles		
Desktop settings related to your logon	System variables	
	Variable Value	
Settings	FP_NO_HOST_C NO	
	NUMBER_OF_P 4	
Startup and Recovery	Path C:\Applied Biosystems\StepOne Softwa	
System startup, system failure, and debugging information	PATHEXT .COM;.EXE;.BAT;.CMD;.VBS;.VBE;.JS;	
Settings	New Edit Delete]
Environment Variables Error Reporting	OK Cancel	
OK Cancel Apply	Edit System Variable	
	Variable name: Path	
	Variable value: : : Program Files MATLAB \R 2007b \bin \win64	
	OK Cancel	

c. In the "System variables" section of the "Environment Variables" dialog box, select the "Path" variable, then click "Edit". Add or modify the path lines with the paths you wish the computer to access. Each different directory is separated with a semicolon. To access the tool, add:

 $\label{eq:linstallDirectoryv77binwin32} where installDirectory is the installation location you chose in step 1.$

4 Working with TmTool[™] Application

Run the experiment on the 7500 Fast, StepOnePlusTM, or the 7900 HT systems following the guidelines in the QuickSetupGuide.pdf. Analyze the raw data using the instrument system software, then save the experiment file (*.eds or *.sds).

For the 7900 HT system, export the multi-component data from the ROXTM dye filter to a *.txt file to analyze the data using the TmToolTM application.

Note that no passive reference dye should be assigned to wells at the time of analysis. At this point you can exit from the instrument software.

4.1 Starting the tool

To start the TmTool application, double click the TmToolv1_0.exe. This file is present in the installationDirectory. The first time this tool is run, it will bring up the end user license agreement. Click on accept. This will bring up the main interface as follows:

ioni v 1.0		
Source File	Please locate data file for analysis using the Search button	Browse
Instrument Type	Select	
Reference		
Omit		Clear
		Clear
Region Selection	O Use defaults	Browse
	C Use Custom	
Results Location	Please select a directory location to store results using the Search button	Browse.
Plot Options	🕑 Generate Plot	
	Display References	
	Show Boltzmann fit	

Figure 1: Main Interface of the TmToolTM Application

4.2 Importing a source file for analysis

The TmToolTM application can import data from *.eds files generated by the StepOnePlusTM and 7500 Fast Real-Time PCR Systems. For the 7900 HT system, the multi-component data from the ROX filter must be first exported from the SDS run file for use with the TmToolTM application.

1. In the TmToolTM application, click Browse to navigate to the folder that contains the data files.

📣 Tm Tool v1.0		
Source File Instrument Type	Please locate data file for analysis using the Search button.	Browse
Reference		Clear
Omit		Clear
Region Selection	 O Use defaults ○ Use custom 	Browse
Results Location	Please select a directory location to store results using the Search button.	Browse
Plot Options	 Generate Plot Display References Show Boltzmann fit Export Plot Data 	
		Calculate Tm

2. Select the data file, then click **Open:**

By default, the *.eds files are listed.

• To analyze files from the StepOnePlusTM or 7500 Fast systems, select an *.eds file, then click **Open:**

Select input da	ta file				? 🛛
Look in:	Completed Ru	ins	• (=	1 💣 🎫	
à	example 1.eds				
My Recent Documents					
Desktop					
My Documents					
My Computer					
(
My Network	File name:			•	Open
Places	Files of type:	Step One Plus or 7500 Fast of	lata file(*.eds)	•	Cancel

• To analyze a text export file from the 7900HT system, select from the "Files of Type" dropdown list, select a *.txt file, then click **Open**.

Select input da	ta file	? 🛛
Look in:	7900 💌 🗢	⊨ 🗈 📸 •
à	🗐 example1.txt	
My Recent Documents		
Desktop		
1		
My Documents		
My Computer		
		>
My Network Places	File name: example1.txt	▼ Open
	Files of type: Multicomponent text export for 7900(*.bx	t) Cancel

4.3 Selecting the instrument type

From the Instrument $\ensuremath{\mathbb{T}ype}$ dropdown list, select the instrument type used to generate the data.

You need to select a source file and the instrument type to enable the Reference, Omit, and Calculate Tm buttons. The instrument type is validated with the source file at the time when the Calculate Tm button is pressed.

Source File	Please locate data file for analysis using the Search button.	Browse
Instrument Type	Select	
Reference	StepOnePlus TM 7500 Fast 7900 HT 384 wells 7900 HT 96 wells	Clear
Omit_		Clear
Region Selection	I Use defaults O Use custom	Browse
Results Location	Please select a directory location to store results using the Search button	Browse
Plot Options	Generate Plot Display References Show Boltzmann fit Export Plot Data	

4.4 Identifying reference wells

If you designate one or more reference wells, the difference between the Tm of the sample in a given well and the average Tm value across the reference wells is calculated and reported as the Δ Tm. You can choose to have the data from the reference wells superimposed on the data from the sample well in the well data plots that the tool generates.

To identify reference wells:

- 1. Click "Reference" to view a list of available wells. The list of available wells is dependent on the type of instrument you selected in the previous step. You need to select the instrument type to enable the "Reference" button.
- 2. Select the wells that contain reference samples.
 - Press the ctrl key while you click to select non-contiguous wells.
 - Press the shift key while you click to select contiguous wells.
 - Click Select all to select all wells in the list.
- 3. Press "OK" when you are finished identifying the reference wells.

🛃 Tra Teal ví), 7		(= (= 💌	4 🔳 🛛 🛛
			Select wells (multiple by using
Source File	exampie Leds	Browse	A1 🔨
instrument Type	Step One Flus 📉		A2 A3
			A4
Deference	▶		A6
Reference		Class	A7 A8
			A9 A10
Omit			A11 A12
		Clear	B1 B2
Region Selection	© Lise detaults		B3
	O Lise clistom	Srowse	B4 B5
			B6 B7
Results Location	Peace select a directory location to store results using the Search button	Browse]	B8
Plot Options	☑ C-recal- Phil		B10
	 Display References 		B11
	□ Srtw intermentin		Select all
	The Export Mot Data		
	Caler	ulate Tm	OK Cancel

4.5 Identifying wells to omit from analysis

You may want to omit certain wells from analysis. Figure 2 shows selection of wells A2 and A9 to A12 for omission.

To identify wells to omit from analysis:

- 1. Click Omit to view a list of available wells. The list of available wells is dependent on the type of instrument you selected in the previous step. You need to select the instrument type to enable the Omit button.
- 2. Select the wells to omit from analysis.
 - Press the ctrl key while you click to select non-contiguous wells.
 - Press the shift key while you click to select contiguous wells.
 - Click Select all to select all wells in the list.
- 3. Press OK when you are finished identifying the wells to omit.

🖬 Tra Tasi yil 2		[1] [1] [2]	
Searce File Instrument Type	inampintade Step Cea Flat	frase.	Select wells (multiple by using A1 A2 A3 A4
Reference		Clear	A5 A6 A7 A8 A9 A10
Omit		Gear	A11 A12 B1 B2
Region Selecter	di Crix deta da O une natione	Drawn.	B3 B4 B5 B6
Results Longitor	Now sold a contry locks is downards very lefter that to	Boste.	87 88 89
Piet Option	2 Development (Mat Dependent trans Development (Mathematical) Development (Mathematical)		B10 B11 Select all
		Calasiane Tan	OK Cancel

• Note that if certain wells are selected for omission, they will not be available in the "Reference..." well selection window.

⁽²⁾Also, note that the axis limits of the data plots per well are set using the range of fluorescence values from the set of non-omitted wells.

4.6 Region of analysis: default selection

The region of analysis is automatically chosen such that there is a continuous rise in fluorescence values over that temperature range. If there are multiple regions in the melt profile that meet this definition, the steepest of these regions is selected for analysis. Given below is a typical melt profile. The regions between the black dashed lines indicate the region of analysis.



It is recommended that the default settings for region of analysis be used for the first analysis of every file.

Note that the Tm detection process is impacted by the appropriate selection of the region of analysis. For fluorescence profiles with multiple phases of melting, the ability to choose the region of analysis allows the flexibility of analyzing multiple regions of interest in the melt profile in separate runs of the TmToolTM application.

4.7 Choosing a location to store the results

From the TmToolTM application, you can select a directory location for storing analysis results. The TmToolTM creates a folder at that location with the same name as the name of the file given for analysis.

• Note that if you do not select a location for storing the results, the default location for storing analysis results is the current working directory where the TmToolTM has been installed.

M Tm Tool v1.0			
			ĸ
Source File	example1.eds		Browse
Instrument Type	StepOnePlus TM		
Reference			Clear
	A1 A8 A10 A11 A12	Browse For Folder	
Omit			
		Select Directory to Open	Clear
Region Selection	 O Use defaults O Use custom 		Browse
Results Location	Please select a directory location to store results using the Se		Browse
Plot Options	Generate Plot		
	Display References	Folder: Desktop	
	Show Boltzmann fit	Make New Folder OK Cancel	
	Export Plot Data		
			Calculate Tm

4.8 Selecting plot options

You can select plot options in the main window of the TmTool[™] application.

- Select the Generate Plot checkbox to create a data plot for each well that is not omitted.
- Select the Display References checkbox if you designated reference wells (as per Section 4.4) and to superimpose reference well data on top of the plot of the data from the well (see Figure 2)
- Select the Show Boltzmann Fit checkbox to superimpose the fit data on the sample data from the well. The Boltzmann curve display can be used to visually check the quality of the calculated fit (see Figure 2).
- Select the Export Plot Data checkbox to create an output file of the raw data in a tab delimited format within the results location (see Figure 2).

🛃 Tm Tool v0.7		
Source File	example1.eds	Browse
Instrument Type	StepOnePlus TM	
Reference		Clear
Omit	A1,A9,A10,A11,A12,	Class
Region Selection	 O Use defaults ○ Use custom 	Browse
Results Location	Please select a directory location to store results using the Search button.	Browse
Plot Options	Generate Plot Display References	
	 Show Boltzmann fit Export Plot Data 	
		Calculate Tm

Figure 2: Plot Options

4.9 Calculate Tm

Press the Calculate Tm button (see Figure 2) to start calculating the Tm values. A progress bar appears to indicate that the processing is in progress.



When analysis is complete, the plots per well and the table of Tm values are generated. Results are stored in the folder indicated in the results location in section 4.7.

4.10 Reviewing the Results

4.10.1 Data Plots

After one round of execution, the main interface of the TmToolTM application as seen in Figure 1 is restored.

Plots of data per well are generated in the results location. An example data plot generated for well A2 is shown in Figure 3.

- The green line shows the Boltzmann fit produced by the tool.
- The dotted vertical lines delimit the region of analysis.
- The blue vertical line indicates the Tm value produced by the fitting.
- The blue line showing the melt curve is the fluorescence data from the sample in that well.
- The position where warnings or error flags appear, if any are generated, is also indicated in the figure. This example has no flags associated with the result.



Figure 3: A sample plot of data from a well

4.10.2 Table of Tm Values

The $TmTool^{TM}$ application generates a table of melting point temperatures for each well. The results are stored in a tab-delimited file with the naming convention "<filename>_Tm" where <filename> is the name of the input data file. This table has the following columns:

- Wells: well name
- Tm Boltzmann (deg C): Tm from fitting data in the region of analysis to the Boltzmann Equation.
- Delta Tm: If reference wells were indicated, the difference between Tm from the given well and the average Tm of the reference wells is reported as the delta Tm.
- Warning: Warning/error flags that may apply to the result. For a complete list of error/warning messages see <u>Table 1</u>,
- Start T for Boltzmann Fit: Start temperature of the region of analysis
- End T for Boltzmann Fit: End temperature of the region of analysis

The output file "<filename>_Tm.txt" can be opened using Microsoft Excel software. An example of the output file is shown in Figure 4.

	A	В	С	D	E	F
1	Wells	Tm Boltzmann (deg C)	Delta Tm	Warning	Start T for Boltzmann fit (deg C)	End T for Boltzmann fit (deg C)
2	A1					
3	A2	40.8			28.25	52.66
4	A3	40.8			29.09	52.66
5	A4	40.3			28.25	51.82
6	A5	41.2			29.09	53.5
7	A6	40.9			29.09	53.5
8	A7	40.6			29.09	52.66
9	A8	40.4			28.25	52.66
10	A9					
11	A10					
12	A11					
13	A12					
14	B1	41.1			28.25	53.5
15	B2	40.9			28.25	52.66
16	B3	40.6			28.25	53.5
17	B4	40.3			28.25	52.66
18	B5	41.2			28.25	54.34
19	B6	41.1			28.25	54.34
20	B7	40.7			28.25	54.34
21	B8	40.5			28.25	54.34
22	B9	92.1		Low Signal	82.11	96.42
23	B10	94.3		Low Signa	85.48	98.95
24	B11	94.3		Low Signal	89.69	96.42
25	B12			Undetermin	ned	
26	C1	46.7			24.05	53.5
27	C2	46.7			30.78	54.34

Figure 4: Table <filename>_Tm.txt

²Entries corresponding to omitted wells are left blank in the <filename>_Tm.txt. In Figure 4, wells A1 and A9 to A12 were omitted from analysis.

Formatte Roman, 12

Formatte Roman, 12 grammar

Deleted:

Formatte Roman, 12

Formatte Roman, 12 grammar

Deleted:

• For quality purposes, it is recommended that you review the data plots for all Tm values generated by the tool.

4.10.3 Table of Region of Analysis

The TmToolTM application generates a tab-delimited text file called <filename>_ AnalysisRegion.txt. This table has three columns:

- Wells: well name
- Start T for Boltzmann Fit: Start temperature of the region of analysis
- End T for Boltzmann Fit: End temperature of the region of analysis

The <filename>_AnalysisRegion.txt can be opened using Microsoft Excel® software. An example is shown in Figure 5. The use of this file is explained in section 4.11.

	A	В	С
1	Wells	Start T for Boltzmann fit (deg C)	End T for Boltzmann fit (deg C)
2	A1		
3	A2	28.25	52.66
4	A3	29.09	52.66
5	A4	28.25	51.82
6	A5	29.09	53.5
7	A6	29.09	53.5
8	A7	29.09	52.66
9	A8	28.25	52.66
10	49		

Figure 5: Table <filename>_AnalysisRegion.txt

4.10.4 Table of Plotted Melt Curve Data (optional)

The TmToolTM application generates a tab-delimited text file called as <filename>_ SmoothedMeltCurves.txt. This table has the plotted data for each well along the rows.

4.11 Defining Custom Region of Analysis for wells where necessary

In some instances the default region of selection for analysis may not be appropriate to the data, as deemed by the closeness of fit of the Boltzmann fitting line that can be superimposed over the sample data. To define a custom region of analysis:

1. Locate the "<filename>_ AnalysisRegion.txt" in the results location and open it with Microsoft Excel software. Edit the start and end temperature for the region of analysis of each well as required (see Figure 6). For example, for well C1, the tool had selected 24.05 to 53.5, but the region was redefined from 35 to 53.5° C.

Formatte

Roman, 12 Formatte Roman, 12

	Δ	B	C	D	F		grammar
1	Wells	Start T for Boltzmann fit (deg C)	End T for	Boltzma	nn fit (dea C)		Deleted:
2	A1		2.1.0 1 101		(009 0)		
3	A2	28.25	52.66				
4	A3	29.09	52.66				
5	A4	28.25	51.82				
6	A5	29.09	53.5				
7	A6	29.09	53.5				
8	A7	29.09	52.66				
9	A8	28.25	52.66				
10	A9						
11	A10						
12	A11						
13	A12						
14	B1	28.25	53.5				
15	B2	28.25	52.66				
16	B3	28.25	53.5				
17	B4	28.25	52.66				
18	B5	28.25	54.34				
19	B6	28.25	54.34		Edit start te	mperature for region of analysis	
20	B7	28.25	54.34				
21	B8	28.25	54.34				
22	B9	82.11	96,42		East end ter	nperature for region of analysis	
23	B10	85.48	98.95				
24	B11	89.69	/ 96.42	/			
25	B12			<u> </u>			
26	<u>C1</u>	35	53.5				
27	C2	30.78	54.34				

Figure 6: Define Custom Range of Temperature for Analysis

2. Save this edited version of the analysis region file anywhere. Name the file as <any text> AnalysisRegion.txt.

Note that the file containing the custom region of analyses must be named as "*_AnalysisRegion.txt" for the TmToolTM application to locate the file in the next step. "*" in the file name can stand for any text.

3. Select the "Use Custom" option from the TmToolTM application as indicated in <u>Figure</u> <u>7</u>, The "Browse" button is activated. Click Browse to navigate to the location where you saved the analysis region file (step 2).

Tm Tool v1.0							
Source File	example1.eds						Browse
Instrument Type	StepOnePlus TM						
							_
Reference		Select tab delin	nited text file	to specify analysis region	for each well	? 🗙	Char
Omit	A1,A8,A10,A11,A12,	Look in:	istrib	ed_AnalysisRegion.txt	▼ ← È d* III	•	Clear
Region Selection	O Use defaults ⊙ Use custom ➤	Documents Documents Desktop					Clear Browse
		My Documents					
Results Location	C:\Documents\	My Computer					Browse
	 Display Referer Show Boltzmar 	My_Network	File name:	example 1_modified_Analysis Re	gion.txt 💌	Open	
	✓ Export Plot Data	Places	Files of type:	text	•	Cancel	
							Calculate Tm

4. Run the analysis again using the "Calculate Tm" button.

Figure 7: Select file with custom analysis regions defined

If the results location was not changed, a new directory called <filename> 2 is created and the new results files are saved there. Figure 8 shows the data plot from well C1 with the region of analysis as defined in the previous steps.

grammar

Deleted:



Figure 8: Well C1 with custom region of analysis

4.12 Designating Reference Wells, generating ∆Tm and Plots with References

It is also possible to estimate the degrees of shift in Tm values between all the wells and wells C1 and C2.

- 1. Import a .eds file (see section 4.2).
- 2. Select the instrument type (see section 4.3).
- 3. Designate reference wells (see section 4.4).
- 4. Run the tool with the plot option Display References checkbox selected.
- 5. Use custom region of analysis defined in the <filename>_AnalysisRegion.txt. The analysis settings for this scenario are shown in Figure 9.

Martin Tool v1.0	
Source File Instrument Type	example1.eds Browse StepOnePlus TM
Reference Omit	C1,C2, C1ear
Region Selection	O Use custom
Results Location Plot Options	Z \Software Browse Ø Generate Plot Display References Ø Show Boltzmann fit Export Plot Data
	Calculate Tm

Figure 9: Using reference wells to generate ΔTm

An example of the plot generated with "Display References" selected is given in Figure 10. A Δ Tm of 5.9 degrees C is observed between the average Tm of the reference wells and the Tm generated from data in well A3.



Figure 10: A sample plot of data from a well with reference wells

The result file $< \texttt{filename} _ \texttt{Tm.txt}$ now has the column $\Delta \texttt{Tm}$ populated. If you open this file using Microsoft Excel software, you can sort the wells by the $\Delta \texttt{Tm}$ value to identify the well with the largest shift in Tm. The buffer condition for that well can then be identified as the buffer in which the sample proved to be maximally stable.

5 Troubleshooting

- Observation: The Microsoft Excel software has difficulty with opening the tab-delimited result files
 Possible cause: The path and file name combinations of the tab delimited result files may exceed 270 characters.
 Solution: Rename the tab-delimited results files, rename the file using a shorter name and then use MS Excel software to open it.
- Observation: The TmToolTM application is not able to locate the file where you saved the modified regions of analysis.

Possible cause: The path and file name combinations of the tab delimited result files may exceed 270 characters. Or, you did not save the file using (*_AnalysisRegion.txt) as the naming convention.

Solution: Rename the file with the analysis regions as *AnalysisRegion.txt file where the path and filename combination is under 270 characters.

Observation: If there are multiple melt phases (Figure 11), the auto selected region of analysis selects the entire region over which a rise in fluorescence value was observed. Possible cause: The tool is programmed to select the *largest* contiguous region of rise in fluorescence until the point where the slope becomes negative. Solution: If this is not the desired behavior, it is possible to define a different region of analysis following the instructions in section 4.11.



Figure 11: Multiple phases of melt, or, initial melt phase slowly rising

• Observation: If there are two melt phases, with one rising to a higher magnitude than the other (Figure 12), the auto selected region of analysis will still select the region of the steeper rise, even if it has a smaller magnitude.

Possible cause: The tool is programmed to select the *steepest* contiguous region of rise in fluorescence.

Solution: If this is not the desired behavior, it is possible to define a different region of analysis following the instructions in section 4.11.



Figure 12: Multiple melt phases

Analysis tips

- **To generate plots of data from multiple wells:** The TmtoolTM application provides plot of data from a single well. To generate plots of data from multiple wells together, use the data output in example_SmoothedmeltCurves.txt produced when the "Export Plot Data" option is selected prior to the analysis.
- **Review data plots**: It is recommended that, for quality purposes, all Tm values generated by the tool be reviewed by observing the data plots.
 - For example, for well B9, there is a low signal flag raised. Well B9 data is presented in Figure 13: Well with warning flag "Low Signal". The signal was very low, compared to the highest magnitude of the signal seen in that group of wells analyzed together. This result should be rejected.





• Another example is presented in Figure 14. The automated region of selection in this case was too broad which resulted in a poor fitting to the data. Fitting can be improved by a good choice of the region of analysis. Look at section 4.11 to learn how to define custom region of analysis for more accurate results.



Figure 14: An example of where a custom region of analysis selection is recommended

• Flags: The software applies flags to a broad range of results. For flagged results, review the Tm value and the Boltzmann fit on the plot and if the results are reasonable, you may choose to ignore the flag.

6 Definitions

Warning	Definitions
Message	
Empty Well	The well contains no data.
Custom Region	User-designated region of analysis instead of using default region of analysis auto selected by the tool.
Low Signal	The ratio of average signal value in well to the absolute maximum seen in the set of data from wells being analyzed (excluding omitted wells) is low
Few Points In	The Boltzmann fitting is performed with less than 5 points because of a very
Region of	short region of positive slope observed in the data.
Analysis	
Fitting with	If a good convergence was not reached in the fitting process, this warning is
warning	supplied.
Custom	If a custom region of analysis was selected, but Boltzmann fitting did not
Region,	execute properly, this error flag is generated.
Boltzmann	
failed	
Undetermined	Additional errors such as a fitting error or a region of analysis selection error
	happened inside the algorithm for which a Tm generation failed for the well.

Table 1: Warning/Error Messages and their Definitions

Table 2: Acronyms/Terms and Definitions

Acronyms	Definitions			
NPC	No Protein Control			
LC	Ligand Control			
Tm	The inflection point in a temperature series data that is ideally a sigmoid.			
Reference	One or more data wells may be designated as a reference well. Each well will have its			
Well, ∆Tm	own Tm value based upon data obtained from that well. If one or more reference wells			
	are designated, an average of the Tm values corresponding to the reference wells is			
	calculated. Δ Tm, defined as difference between average Tm value of reference wells			
	and the Tm value of the given well will be calculated for each well.			
Undetermined If the data from a well is such that a reliable Tm cannot be generated, a T				
	Undetermined is returned. User can force the software to return a Tm value			
	nevertheless by selecting custom region of analysis.			
Region of	A region of the melt curve data is automatically selected by the software as the region			
analysis	of analysis. The data from the region of analysis is fit to the Boltzmann equation to			
	return a Tm value. The algorithm looks for a region of the curve with the steepest slope			
	increase and uses that as the region of analysis.			
	User may choose to override this auto selected region by selecting custom region of			

	analysis.
Instrument	The Applied Biosystem Real-time instrument on which the experiment was run needs
Туре	to be selected to ensure that software applies the algorithm parameters that are suitable for the chosen instrument
	for the chosen instrument.

Appendix

Boltzmann Fitting Method

The Boltzmann equation: $F(T) = F(pre) + \frac{[F(post) - F(pre)]}{1 + e^{\frac{(T_m - T)}{C}}}$

where:

- F(pre) is the fluorescence reading before the transition or melting at the start of the analysis region
- F(post) is the fluorescence reading after the transition or melting, at the end of the analysis region
- T_m is the melting temperature
- C is the enthalpy of the reaction

Data are fit to this equation to generate the T_m.

© 2010 Life Technologies Corporation. All rights reserved.

The trademarks mentioned herein are the property of Life Technologies Corporation unless otherwise noted. Windows and Excel are registered trademarks of Microsoft Corporation in the United States and other countries.