



User Manual PR.ThermControl Software

Content

1. Technical Information	3
1.1. System Requirements	3
1.2. Backward Compatibility	3
2. Home Screen and General Usage.....	4
3. Discovery Scan	4
4. Melting Scan	6
5. Annotations and Results.....	8
6. Data Export	10
6.1. Discovery Scan	10
6.2. Melting Scan	10
6.3. Annotations and Results	11

NanoTemper® Technologies' Prometheus™ instruments detect changes in the fluorescence of the amino acid tryptophan (and fluorophores with equivalent spectroscopic properties) over a wide range of temperatures using nanoDSF™, an advanced Differential Scanning Fluorimetry technology. The instruments can be used to induce thermal unfolding of proteins and to determine thermal unfolding transition temperatures. Furthermore, the instruments are equipped to investigate chemical unfolding and the free energy of unfolding ΔG in an extraordinarily straightforward and fast manner.

The PR.ThermControl software is dedicated to running and analyzing thermal unfolding (and optionally refolding) experiments on Prometheus instruments.

1. Technical Information

1.1. System Requirements

If the necessary licenses have been purchased, PR.ThermControl software can be installed on additional computers for convenient data analysis. The computers have to meet the following requirements:

Operating system:	Windows 7 64 Bit or higher
CPU:	Intel Core i5 or better
RAM:	8 GB or more
Hard disk:	20 GB or more free disk space available
Display resolution:	1600 x 900 or better
Software:	Microsoft .NET 4.6.2 framework (included in installer of PR.ThermControl software)
Operating system language:	English or German

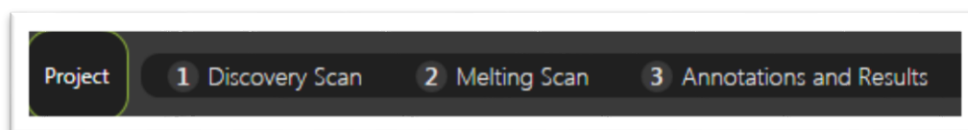
An external computer mouse is necessary to access all software features.

1.2. Backward Compatibility

Measurements collected with all previous versions of PR.ThermControl (previously called PR.Control) software are compatible with this version.

NanoTemper Technologies provides support for all current and previous versions of PR.ThermControl.

2. Home Screen and General Usage



To perform a new measurement, start the PR.ThermControl software, which will show the *Project* home screen. Click *Create New Project* to enter a file name and location. The file will be saved in .prc format. Optionally, enter relevant information into the *Project Name* and *Comments* fields.

Alternatively, previous project files can be loaded by clicking *Browse* to analyze previous experiments or to add additional measurements. Recently loaded files are listed chronologically on the right.

Use the *Save Changes* button at any time to save modifications of the file. An asterisk in the software title bar indicates unsaved changes in the open file. Closing the software will trigger a dialogue box asking whether you want to save the changes.

Three tabs guide the user through running and analyzing nanoDSF measurements:

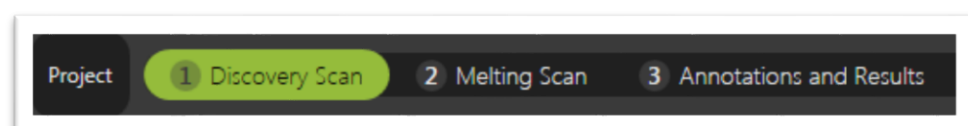
1. Discovery Scan
2. Melting Scan
3. Annotations and Results

Proceed from 1 to 3 to set up, run and analyze a nanoDSF measurement. As soon as the measurement is started, navigate between all three tabs freely to re-analyze, modify or annotate. More details on each tab follow below.

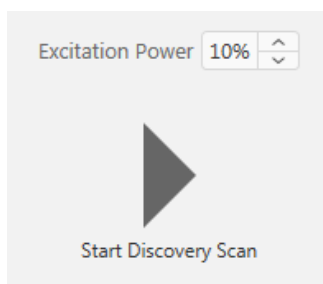
The keyboard shortcut ctrl + z will undo any action, while ctrl + y will redo.

Any graph displayed in the software can be exported by clicking the *Export* button. Options are to copy the graph to the clipboard, to save it as an image (.png or .svg file formats), or to export the raw data needed to recreate the graph in third-party software (.xlsx file format).

3. Discovery Scan



After creating a new project or loading a previous project file, perform a *Discovery Scan* (1) to determine optimal settings for the unfolding experiment. You can vary the excitation power between 1 % and 100 %. Re-scan with different excitation power settings if necessary.



The discovery scan is used to detect the fluorescence intensity and position of each capillary along the entire length of the capillary tray. Each capillary will be visible as a combination of two peaks, which represent the fluorescence intensity at 330 nm and 350 nm, respectively (**Figure 1**). If the Prometheus instrument is equipped with Aggregation Optics, the scattering signal is also displayed for each capillary.

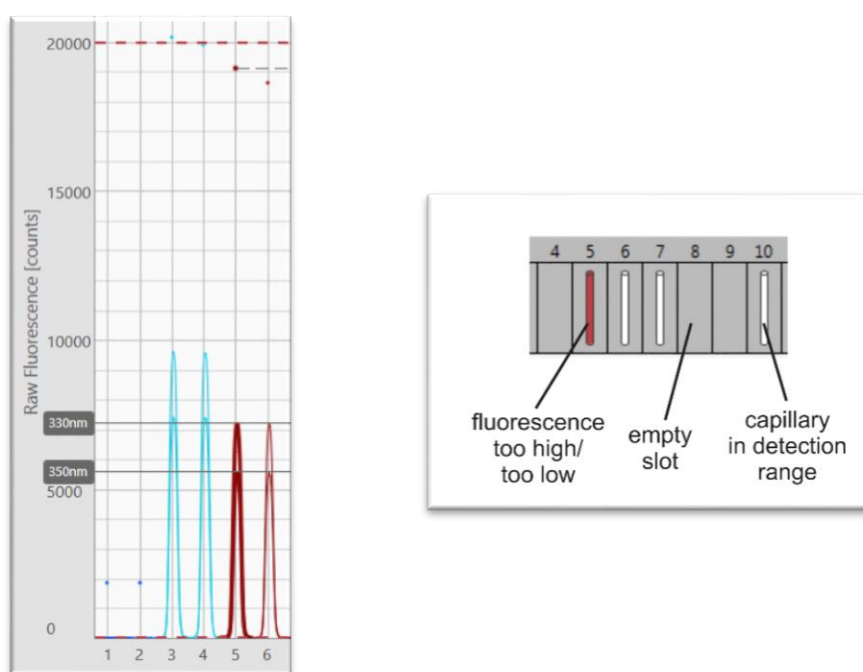


Figure 1: Discovery Scan. The discovery scan provides information about the fluorescence intensities of the samples, and is used to determine optimal measurement settings and the positions of occupied capillary slots. (Left) The fluorescence intensities of the 330 nm channel and 350 nm channel are displayed as two separate, nested peaks for each capillary. Clicking on a specific peak or capillary will highlight this capillary's fluorescence at 330 nm and 350 nm, and also display its values for Peak Fluorescence and Integrated Fluorescence on the right of the screen. Both wavelength peaks should be between the upper and lower detection limits (dotted red lines). If the Prometheus instrument is equipped with Aggregation Optics, the scattering signal is plotted on the right y-axis and represented in the graph by a dot symbol for each capillary. Capillaries with non-aggregated samples typically exhibit a scattering signal of 90-110 mAU, while maximum aggregation reaches values of up to 600 mAU. (Right) Capillaries with fluorescence intensities within the dynamic detection range are colored white; capillaries with too high fluorescence intensities are colored red.

The upper and lower limits of detection are highlighted by dotted red lines in the discovery scan profile. The upper limit is 20,000 fluorescence counts (Peak Fluorescence, meaning the height of the capillary peak). Please note that the lower detection limit is dynamically adjusted to the excitation power settings and thus may vary between experiments with different settings.

Zoom into the graph using the scrolling function of the mouse wheel. Holding down shift or ctrl while scrolling will zoom horizontally or vertically, respectively. Holding the mouse wheel also allows to move the graph. To reset the view, click *Zoom Extent*.

Note: The optimal detection range is between 2,000 and 18,000 counts. Some proteins might require limiting the maximal fluorescence counts to 15,000, since the unfolding might result in an atypical increase in fluorescence intensity.

Note: Photobleaching effects are negligible even at high excitation intensities due to the rapid on-the-fly measurement mode.

Note: If the capillary fluorescence exceeds the upper limit, the capillary position will be indicated in red and excluded from the measurement. Capillaries with fluorescence intensities below the measurement limit may not be recognized by the software but can be manually chosen in the Melting Scan submenu (see below).

4. Melting Scan



Once the optimal excitation power settings are determined, continue with the *Melting Scan* (2). The PR.ThermControl software automatically identifies the capillaries for unfolding analysis from the discovery scan. By default, only capillaries that lie in the dynamic detection range are pre-selected for the unfolding experiment. However, you can also add capillaries that fall below the detection limits.

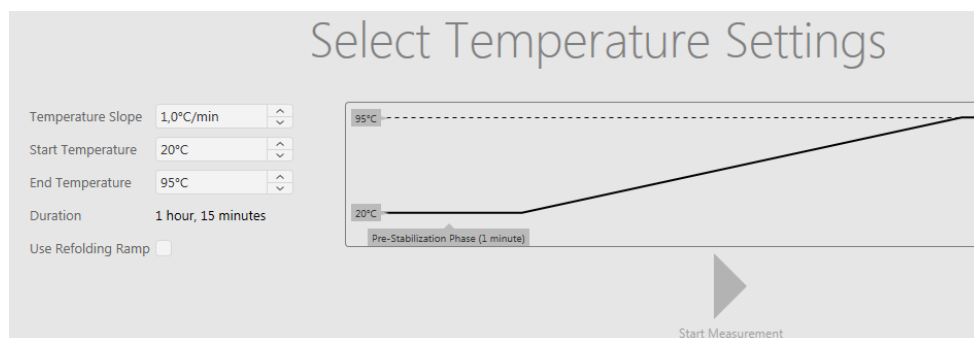
Note: You can multi-select capillaries using shift or ctrl + left click.




Next, define the “Start” and “End” temperature for the thermal unfolding experiment, and set the “Temperature Ramp” to the desired heating rate.

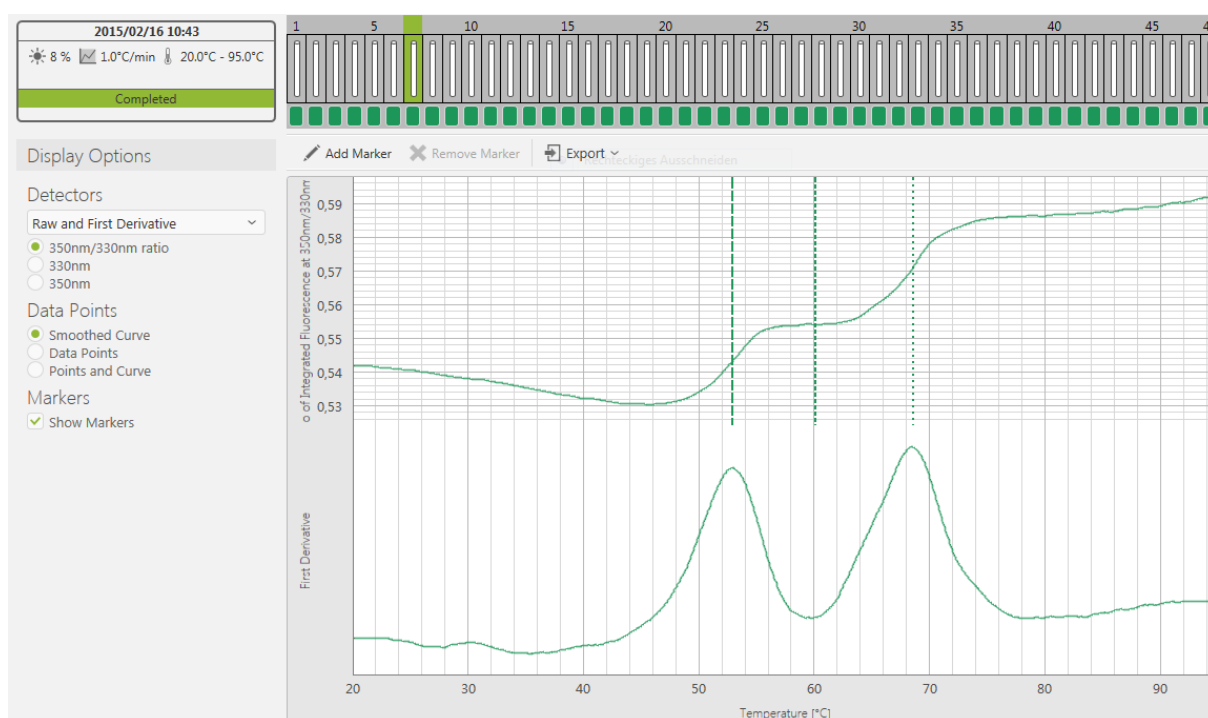
Activate the “Use Refolding Ramp” checkbox to include a refolding experiment with temperature slopes identical to the unfolding experiment.

Start the thermal unfolding experiment by pressing “*Start Measurement*”. After a pre-melting phase in which the start-temperature is reached and the system is equilibrated, continuous scanning of the selected capillaries proceeds while a thermal ramp is applied. The thermal unfolding of proteins in each capillary can be followed in real time. For visualization purposes, the fluorescence emission values at 330 nm and 350 nm and the F350/F330 ratio can be displayed either together or separately.

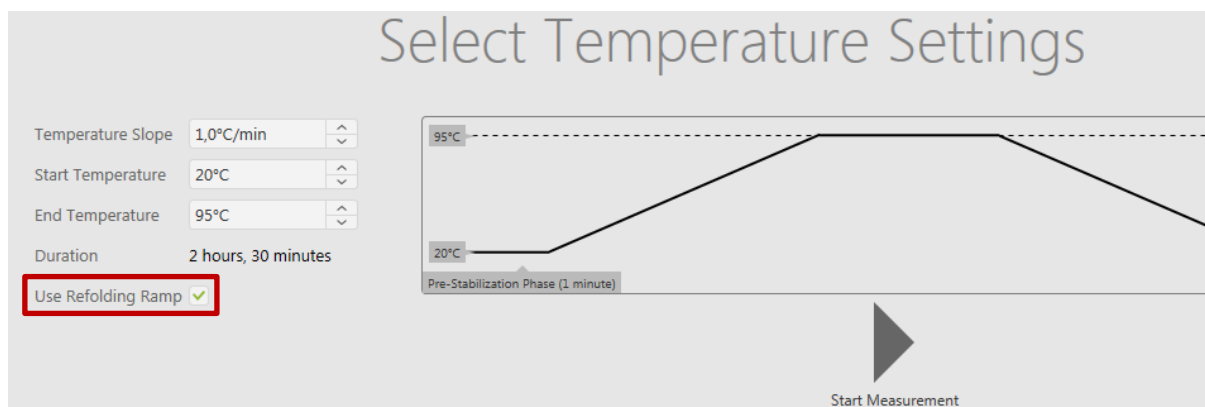
Note that the values plotted in the *Melting Scan* for 330 nm and 350 nm are the Integrated Fluorescence values (area of the capillary peak), not Peak Fluorescence values (height of the capillary peak).



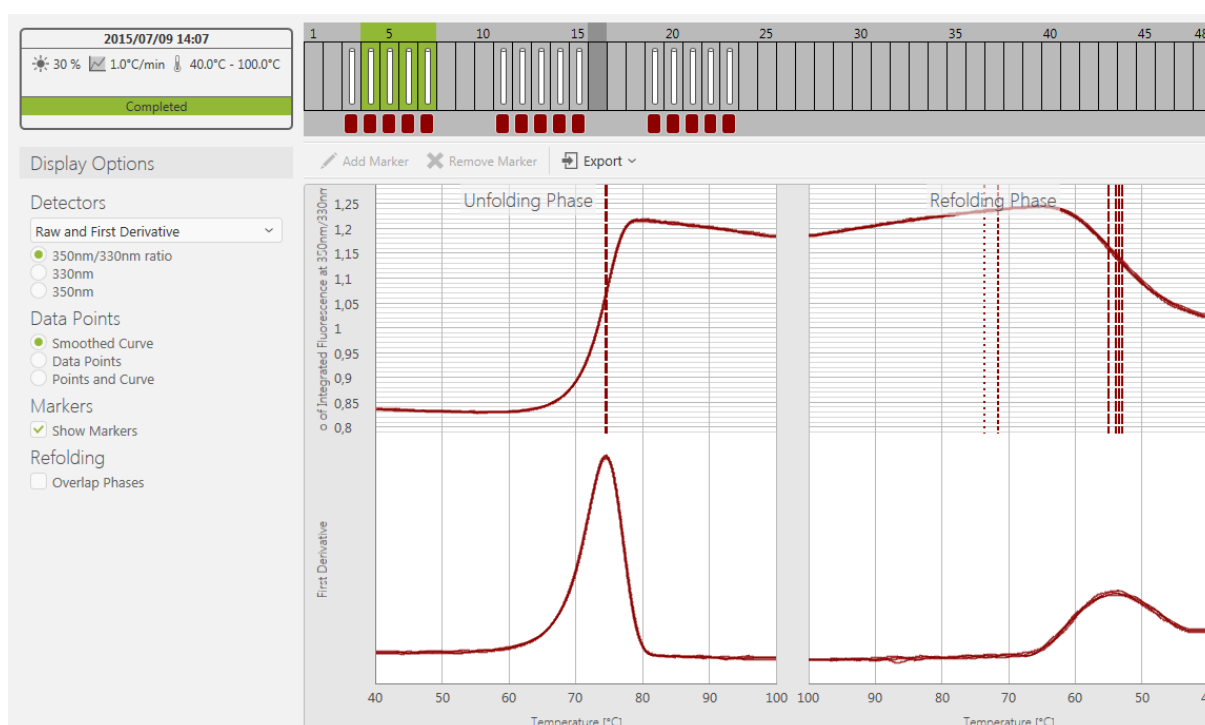
Once the final temperature is reached, unfolding transition midpoints  and unfolding onset temperatures  will be automatically calculated by the software and can be displayed in the melting scan window. The window will also display the unfolding curves and, if selected, the respective first derivative of the curves. In the first derivative view, each local maximum or minimum corresponds to a transition temperature. Transition temperatures and unfolding onsets can be manually added  or removed by clicking the *Add / Remove Marker* buttons. Colored buttons below the capillaries can be used to manually change the color of each curve.



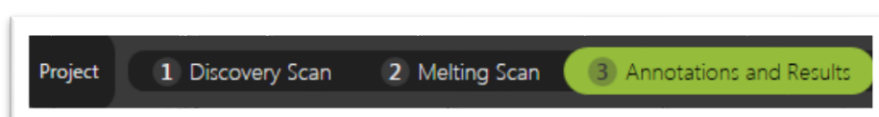
The calculated unfolding transition temperatures as well as manually annotated markers will be added to the annotations table. Raw fluorescence and processed data (raw fluorescence and first derivative values) as well as the corresponding calculated unfolding transition midpoints and annotations can be exported in Excel™ format using the *Export* button. Images can be exported as .svg or .png files.



Folding and refolding can either be shown by overlapping phases or side by side.



5. Annotations and Results



It is not required to enter sample annotations prior to the experiment. Annotations for each capillary can be entered at any time while the melting scan is running or after it is finished. Annotations can either be entered by simple copy-and-paste from a spreadsheet software like Excel™ or manually. Columns can be added or removed by using the *Create New Column* and *Remove Column* buttons on the right side of the screen:

Create New Column

Name

Create New Column

Remove Column

Buffer

pH

NaCl(mM)

Delete Selected Column

Annotations can be entered into multiple fields simultaneously after multi-selection, and subsequently sorted in ascending or descending manner by left-clicking on the column header.

Entering annotations

Capillary	Target	Comment
1		
2		
3		
4		
5		
6		
7		
8		
9		
10		
11		
12		
13		
14		

multiselect ctrl+left click



Capillary	Target	Comment
1	Antibody	
2	Antibody	
3		
4	Antibody	
5	Antibody	
6		
7	Antibody	
8	Antibody	
9		
10		
11		
12		
13		
14		

Sorting function

Capillary	Target	Comment	Number
1			1
2			
3			5
4			
5			2
6			
7			4
8			
9			3
10			
11			



left click on header

Capillary	Target	Comment	Number
3			5
7			4
9			3
5			2
1			1
29			
30			
31			
32			
33			
34			

Note: Multi-selections of capillaries applied in the Melting Scan or Annotations and Results tabs persist after switching tabs.

Capillaries can be selected and a single color can be assigned by clicking on the rectangle next to “Single Color”. Furthermore, capillaries can be colored by categories or gradients after selecting the cells that are to be used for categorizing. The colors will directly translate into the melting curves in the *Melting Scan* window. Color assignment will only apply to selected fields. Multi-selection by using shift / ctrl + left click is possible.

Assign Colors

Single Color:

Categories Colors:

Gradient Colors:

The annotations table includes separate columns for melting onsets (ON), unfolding transition midpoints (inflection points, IP) and manually annotated points (M). Arrows in the column headers indicate whether the values were determined during a heating ramp or a cooling ramp .

The thermal stability of a given protein is typically described by the thermal unfolding transition midpoint, also called inflection point (IP) or T_m [°C], at which half of the protein population is unfolded. It is determined automatically by the PR.ThermControl software via the derivative of the curve. This method circumvents the subjective determination of baseline levels and also allows for the determination of multiple unfolding transition midpoints, e.g. for antibody unfolding or more complex multi-domain proteins. The onset of unfolding (ON) is calculated using the transition midpoint and the slope of the unfolding signal.

Annotation tables can be exported in Excel™ format (.xlsx).

6. Data Export

Each measurement and analysis performed can be exported to be used in third-party software. Throughout this software manual, different export options are mentioned. This section aims to give an overview regarding the file format and the content of each export.

6.1. Discovery Scan

The Discovery Scan tab contains three export options.

Clicking on *Export Raw Data* will create an Excel™ file (.xlsx) containing two sheets. The first sheet gives the major information for each peak, whereas the second one contains the information to reproduce the graph.

The *Copy Chart to Clipboard* button will copy the graph displayed so you can paste it into another software.

The *Export Graph* button will create an image file (.png or .svg) of the zoom extended view displayed by the software.

6.2. Melting Scan

The export button in this tab will display four options.

The *Copy Chart to Clipboard* button will copy the graph exactly how it is displayed in the software (for example only showing selected capillaries, or showing both raw data and first derivative) so you can paste it into another software.

The *Export Graph* button will create an image file (.png or .svg) of the graph as described above.

Clicking on *Export Raw Data* will create an Excel™ file containing four sheets. The first sheet gives an overview of all capillaries contained in the run, including unfolding transition midpoints, unfolding onsets and manually set markers. The other sheets contain the fluorescence at 330 nm, 350 nm and the ratio as a function of temperature for all capillaries.

Using *Export Processed Data* will create the exact same Excel™ file but including also the data of the first derivative for each fluorescence channel and their ratio as separate sheets.

6.3. Annotations and Results

The export option is located at the bottom right of the screen and it will create an Excel™ file containing the table displayed in this tab.

Contact

NanoTemper Technologies GmbH

Floessergasse 4
81369 Munich
Germany

Phone: +49 (0)89 4522895 0
Fax: +49 (0)89 4522895 60

info@nanotemper-technologies.com
<http://www.nanotemper-technologies.com>

Prometheus™ and nanoDSF™ are trademarks.
NanoTemper® is a registered trademark.
NanoTemper® is registered in the U.S. Patent and
Trademark Office.

V06_2018-11-13