

High-throughput analysis of protein thermal stability

Protein Thermal Shift software and reagent kits

- Use Applied Biosystems™ real-time PCR instrumentation to analyze protein thermal stability
- Cost-effective method to screen for ligand–protein binding, optimal buffer conditions, or stability changes
- Simple and rapid method, with results in as little as 0.5–2 hours

The Applied Biosystems™ Protein Thermal Shift™ solution, which includes Protein Thermal Shift Software and reagents, offers an alternative application for real-time PCR systems—the analysis of protein thermal stability, including high-throughput screening of ligand–protein binding, optimization of buffer conditions that promote protein stability, and the effect of mutations or modifications on a protein’s thermal stability.

Proteins are the key molecules studied as targets for new drugs as part of the lead generation process of drug discovery. Drug discovery involves high-throughput screening of thousands of small molecules and ligands with a variety of different assays, requiring many months of time as part of a lead generation program. Protein targets are also a challenge to work with due to their susceptibility for degradation and aggregation—



requiring the addition of protein stability screening without ligand as well. Beyond drug discovery, protein thermal stability screening, which is performed with protein melting techniques, is also employed in many other research programs that utilize native proteins throughout academia and industry. For example, the identification and use of ligands and/or solution (buffer) conditions that maximally stabilize a protein are utilized as part of protein purification, crystallization, and functional characterization.

Protein Thermal Shift assay technology can help improve the success rates of protein purification and crystallization, and help reduce the cost of drug discovery—offering a rapid, cost-effective, high-throughput



screening method compared to other available technologies. The Protein Thermal Shift application, run on our real-time PCR systems, enables you to perform the following inexpensively and efficiently:

- Screen samples for relative protein thermal stability
- Screen samples for ligand binding to proteins in a high-throughput manner
- Screen for small-molecule and fragment-library drug candidates that bind to your protein target

- Screen for antibodies that bind to your protein target
- Screen samples for stability changes after protein point mutations
- Systematically identify suitable and optimal buffer conditions to measure protein–ligand interactions
- Identify optimal buffer conditions to improve protein purification and preparation
- Screen buffers to identify conditions for successful crystallization
- Screen buffers to identify optimal storage conditions for proteins

How does Protein Thermal Shift technology work?

Protein stability is dependent on buffer pH, salt content, and the presence of various cofactors in a protein’s storage or reaction buffer environment. A real-time melt experiment with a protein-binding dye such as the Applied Biosystems™ Protein Thermal Shift™ dye, run on any real-time PCR system (including the Applied Biosystems™ QuantStudio™ systems) yields a fluorescence profile specific to the protein of interest in a given test

buffer environment. Variations in the pH, salt content, or test buffer components are reflected in relative changes to this fluorescence profile and the T_m (melting temperature) calculated from it. The binding of a ligand to a protein also has a stabilizing effect on the protein’s thermal stability, thus leading to a measurable difference in the protein’s fluorescence profile.

Protein Thermal Shift assay

The Protein Thermal Shift reagents enable a protein melt assay, an efficient screening tool, to measure protein thermal stability, identify suitable buffer conditions, and measure protein–ligand interactions. This Protein Thermal Shift assay allows for systematic identification of optimal buffer conditions and ligands that stabilize proteins.

The Protein Thermal Shift assay is very easy to set up and run, and the entire workflow typically takes from 0.5–2 hours, depending on the system and run conditions used (Figure 1). The adjustable ramp rate and thermal accuracy of our real-time PCR systems allow complete flexibility to enable shorter run times

Steps in a Protein Thermal Shift assay

- Mix protein, buffer, ligand, if applicable, and Protein Thermal Shift dye
- Run a melt curve experiment on a real-time PCR instrument
- The protein unfolds as it is heated
- The Protein Thermal Shift dye binds to exposed hydrophobic regions and fluoresces
- Transfer *.eds file to the Protein Thermal Shift Software for analysis
- The melting temperature (T_m) is calculated from the melt curve
- Changes in T_m are correlated to changes in protein stability or ligand binding

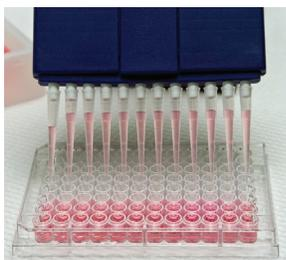
and accommodate the requirements of screening workflows.

Protein Thermal Shift Software

Protein Thermal Shift Software has been developed to analyze protein melt fluorescence readings directly from real-time PCR instrument files.

The complete Protein Thermal Shift solution workflow typically takes 0.5–2 hours

1 Prepare assays



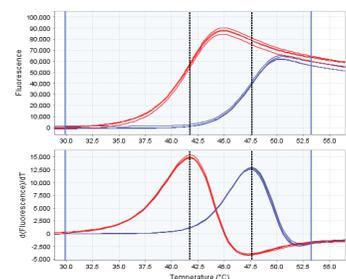
Mix protein, buffer, ligand (if applicable), and Protein Thermal Shift dye.

2 Run melt assay on real-time PCR system (QuantStudio 3, 5, 6, 7, and 12K Flex systems)



The protein unfolds with increasing temperature, and the Protein Thermal Shift dye binds to exposed hydrophobic regions and fluoresces.

3 Analyze data with Protein Thermal Shift Software



The melting temperature (T_m) is calculated from the melt curve; changes in T_m are correlated to changes in protein stability or ligand binding.

Figure 1. Protein Thermal Shift workflow.

Different proteins will have different thermal shift profiles, each with a unique melt curve shape, slope, signal-to-noise ratio, and temperature melt range. The Protein Thermal Shift Software generates one or multiple melting temperatures (T_m values) from these curves by the following two methods: Boltzmann-derived T_m and derivative curve-determined T_m . The Boltzmann T_m values are taken from the inflection point of the fluorescence melt curve plot (Figure 2, top panel) and the derivative T_m values are taken from the top of the peak in the derivative plot (Figure 2, bottom panel).

The Boltzmann method fits the data within an automatically (or manually) identified melt region to the two-state Boltzmann model to generate the T_m . Typically, for proteins with a single melt domain, the Boltzmann method is used. However, for proteins with multiple melt domains, the derivative method can be utilized to determine up to six T_m values per sample. The derivative method uses a numerically computed second derivative of the raw data to estimate the temperatures where up to six

peaks (local maxima) may occur in the derivative profile. An empirically derived threshold on the signal-to-noise ratio is used to determine which local maxima will be detected.

Case study: protein–ligand binding

Protein researchers use a variety of methods to study protein stability and screen for ligands, including biosensors and other high-throughput screening (HTS) methods, as well as lower-throughput methods such as calorimetry and circular dichroism systems. These methods can yield a great level of detail for the protein under study, but range from being very slow and requiring large quantities of sample protein (calorimetry and circular dichroism) to fast but expensive (biosensors). Protein Thermal Shift technology fills the need for fast and inexpensive screening of samples in a high-throughput fashion to quickly narrow down the number of candidates that merit more detailed studies with other technologies. The flexibility, speed, and ease-of-use of the Protein Thermal Shift solution makes it an

excellent option for protein researchers who need to understand how the stability of their proteins is affected at all stages of their research.

Figure 2 demonstrates the identification of a protein-binding ligand, as shown by the increase in T_m of the protein when bound to a ligand. The red curves represent replicate measurements of the protein in solution, while the blue curves represent replicate measurements of the same protein after incubation with a ligand that binds to the protein. For this sample the T_m shifted from $\sim 41.5^\circ\text{C}$ to 47.5°C , indicating that the protein stability increased upon ligand binding.

Summary

Protein researchers use a number of methods to study protein stability and screen for ligands. These methods are generally very slow and tedious to perform and require large quantities of sample protein. Some fast methods exist, but they are very expensive to run compared to the Protein Thermal Shift assay.

Protein Thermal Shift reagent kits utilized on real-time PCR systems enable fast and inexpensive screening of samples in a high-throughput fashion to quickly narrow down the number of candidate ligands or buffer conditions for a wide range of applications that are impacted by protein stability.

Protein Thermal Shift Software allows researchers to quickly compare the shift in T_m (ΔT_m or ΔT_m) between different assay conditions or different ligands added to a sample, relative to a reference sample, thus providing a tool to screen and identify conditions that stabilize (or destabilize) a protein or to screen ligands that bind to the protein of interest.

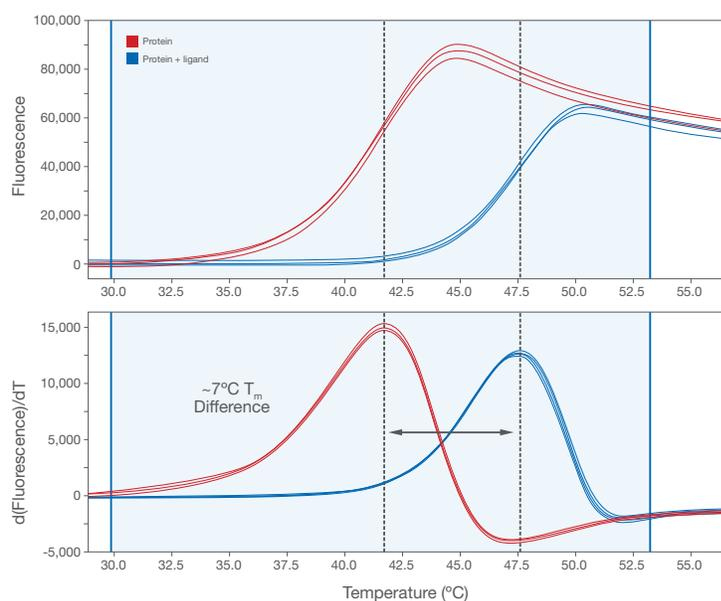


Figure 2. Protein stabilization upon ligand binding. The protein T_m increases after a ligand is added. **(Top)** Fluorescence melt curve plot. **(Bottom)** Derivative curve plot.

Compatible instruments

System	System features	Cat. No.
QuantStudio 3 Real-Time PCR System	<ul style="list-style-type: none"> Capacity: 96-well plates, 0.1 mL tubes or 0.2 mL tubes Fast block: yes Colors: 4 	A28136
QuantStudio 5 Real-Time PCR System	<ul style="list-style-type: none"> Capacity: 96-well plates, 0.1 mL tubes or 0.2 mL tubes, 384-well plates Fast block: yes Colors: 6 	A28138 A28140
QuantStudio 6 Flex Real-Time PCR System	<ul style="list-style-type: none"> Capacity: 96-well, 384-well Fast block: yes Colors: 5 	4485697
QuantStudio 7 Flex Real-Time PCR System	<ul style="list-style-type: none"> Capacity: 96-well, 384-well, TaqMan Array micro fluidic cards Fast block: yes Colors: 6 (21 filter combinations) 	4485698
QuantStudio 12K Flex Real-Time PCR System	<ul style="list-style-type: none"> Capacity: 96-well, 384-well, TaqMan Array micro fluidic cards, OpenArray plates Fast block: yes Colors: 6 (21 filter combinations) 	4471050

Ordering information

Product	Quantity	Cat. No.
Protein Thermal Shift Dye Kit This kit includes the following components: <ul style="list-style-type: none"> Protein Thermal Shift Dye (2,000 reactions) Protein Thermal Shift Buffer (2,000 reactions) 	1 kit	4461146
Protein Thermal Shift Starter Kit This kit includes the following components: <ul style="list-style-type: none"> Protein Thermal Shift Dye (2,000 reactions) Protein Thermal Shift Buffer (2,000 reactions) Protein Thermal Shift Control Ligand (100 reactions) Protein Thermal Shift Control Protein (100 reactions) 	1 kit	4462263
Related products	Quantity	Cat. No.
Protein Thermal Shift Software	1 kit—10 licenses 1 license	4466037 4466038