

# Software development for improved sensitivity of mass spectrometry-based thermal shift assays (MS-TSA) for target engagement and drug discovery

Amanda M. Figueroa-Navedo<sup>a</sup>, Clifford Phaneuf<sup>ab</sup>, Konstantin Aizikov<sup>c</sup>, Alexander R. Ivanov<sup>a</sup>

<sup>a</sup>Department of Chemistry and Chemical Biology, Barnett Institute of Chemical and Biological Analysis, Northeastern University, Boston, MA, <sup>b</sup>Translational Sciences, Sanofi, Cambridge, MA, <sup>c</sup>Thermo Fisher Scientific, Bremen, Germany

## Overview

Thermal shift assays are designed to follow the rate of change in solubility of a protein with increasing temperatures

Thermal Proteome Profiling (TPP) is the first software implemented for mass spectrometry-based assay

-Identifies stabilized protein-drug interactions using melting temperature change ( $\Delta T_m$ )

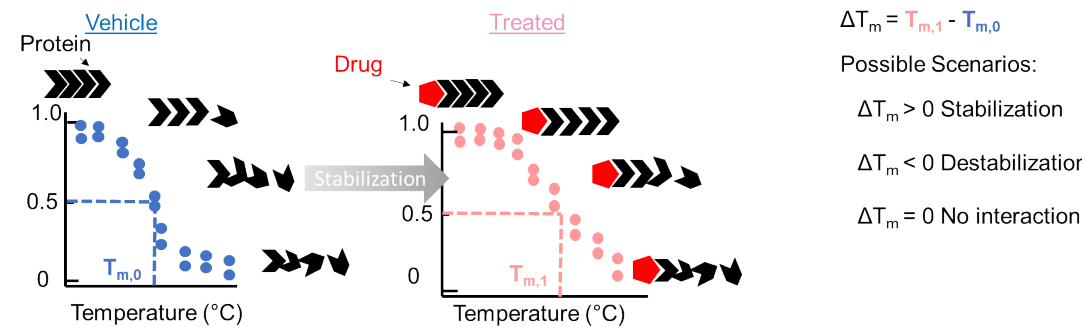
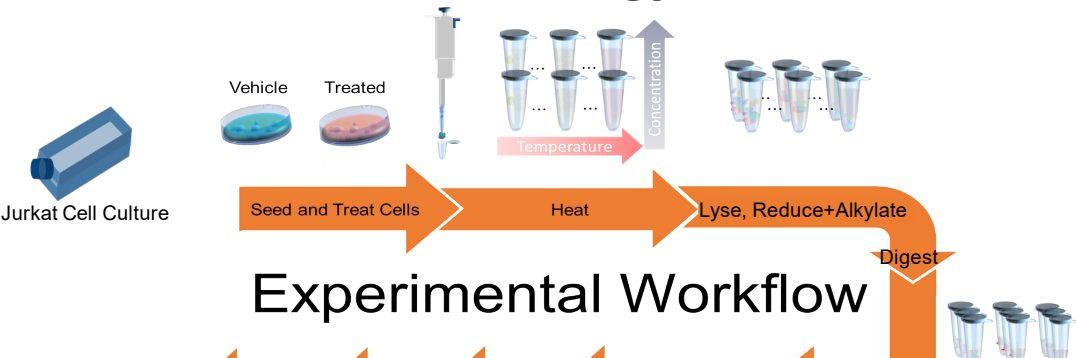
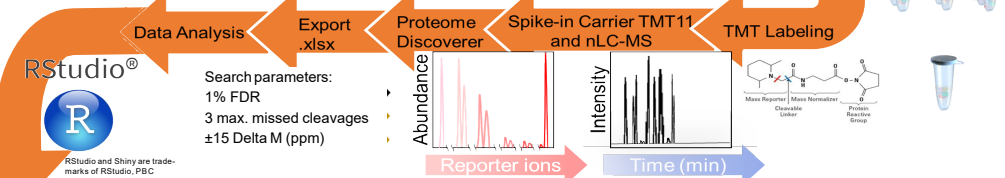


Figure 1. Schematic showing the melting temperatures ( $T_m$ ) in a stabilizing event between vehicle and treated samples

## Methodology



## Experimental Workflow



## Data Analysis Workflow

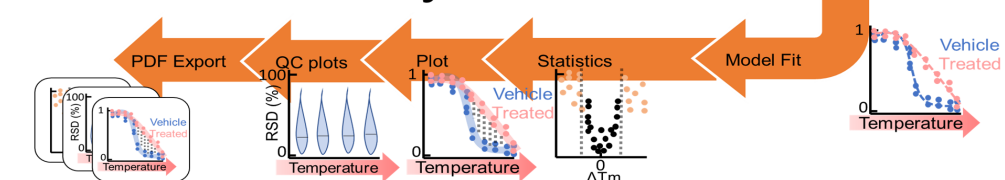


Figure 2. Experimental and data analysis workflows

## Thermal shifts detected by hypothesis testing using residual-sum-of-squares (RSS)

$H_0$ : No thermal shift identified, only one spline fitted curve for vehicle and treated

$H_1$ : Thermal shift identified, two fitted splines curves needed

No shift :  $RSS_0 \approx RSS_1$

Thermal shift:  $RSS_1 \ll RSS_0$

$$F = \frac{d_2}{d_1} \frac{RSS_0 - RSS_1}{RSS_1}$$

$d_1, d_2$  : effective degrees of freedom

## Results

Additional peptide filtering applied in the data analysis pipeline increases the number of overlaps in fitted curves for proteins with unique peptides

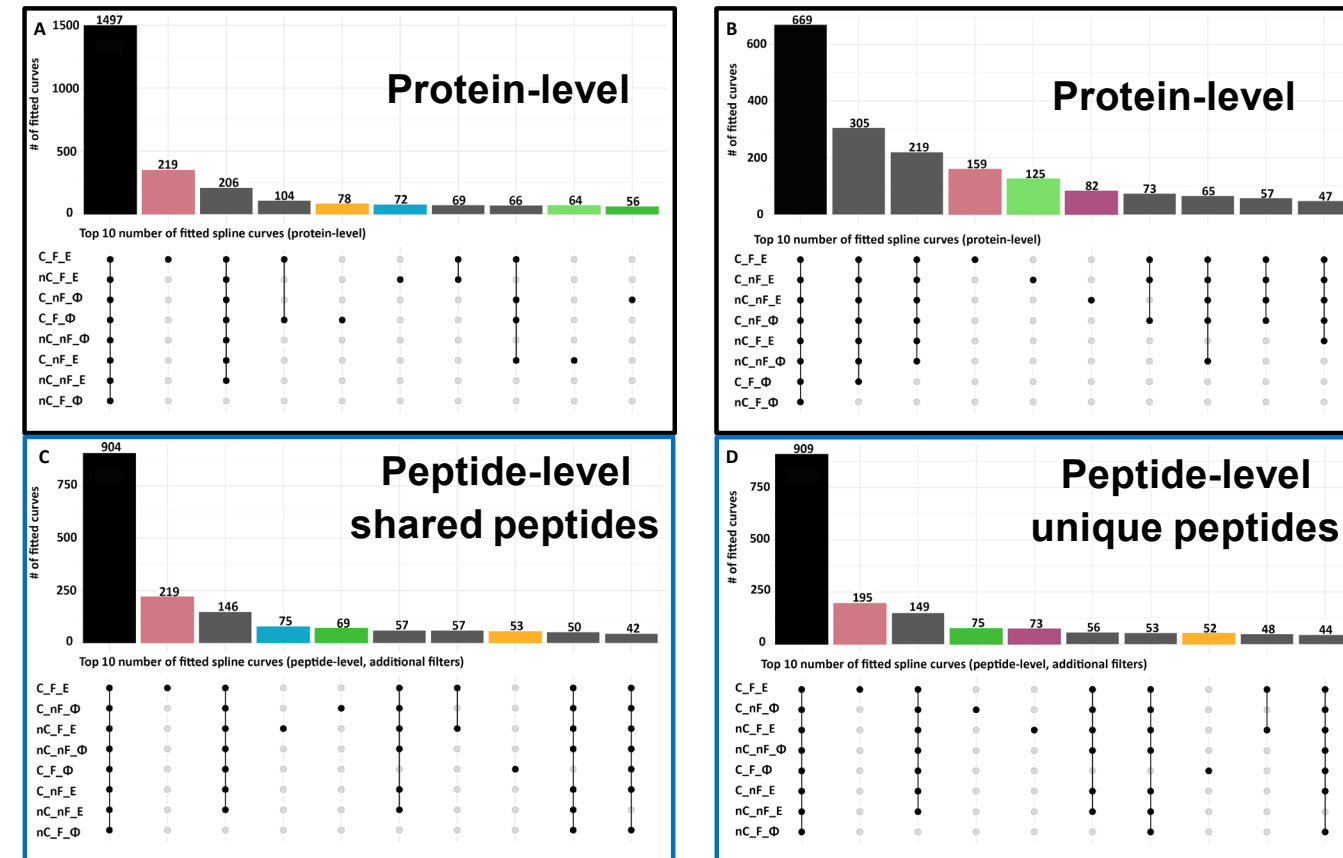


Figure 3. A) Shared peptide quantification results increase the number of overlaps in fitted melt curves B) Protein melt curves using unique peptides show a 65% decrease in overlaps C) Additional filtering on shared peptides decreases the overlaps in number of fitted curves by 40% D) Unique peptide filtering increases the melt curves by 27% in overlaps from 8 different approaches

## Quality Control

Missing values, signal-to-noise ratios and variability in reporter ion abundance channels are key components that affect melt curve quality

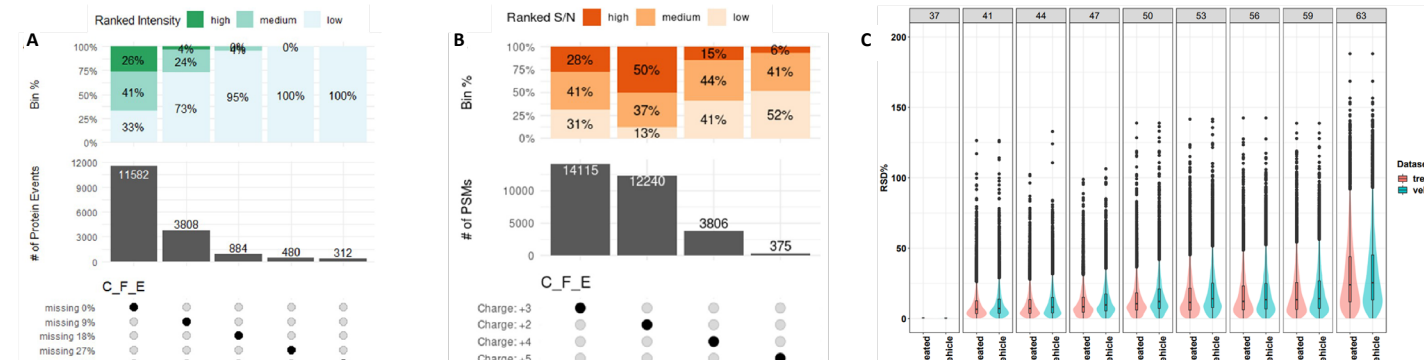


Figure 4. A) Upset plot shows that missing values in melt curves are more likely if the raw reporter ion abundance intensities are lower at the baseline temperature (37° C) B) Upset plot at the PSM-level show higher charge states have lower S/N values at the baseline temperature C) Relative standard deviation of the reporter ion abundances increase with higher temperatures due to a decrease in reporter ion abundance

Stabilization ● Destabilized Shift ● Destabilized  $T_m$  ● Not Shifted ● Stabilized  $T_m$  ● Stabilized Shift

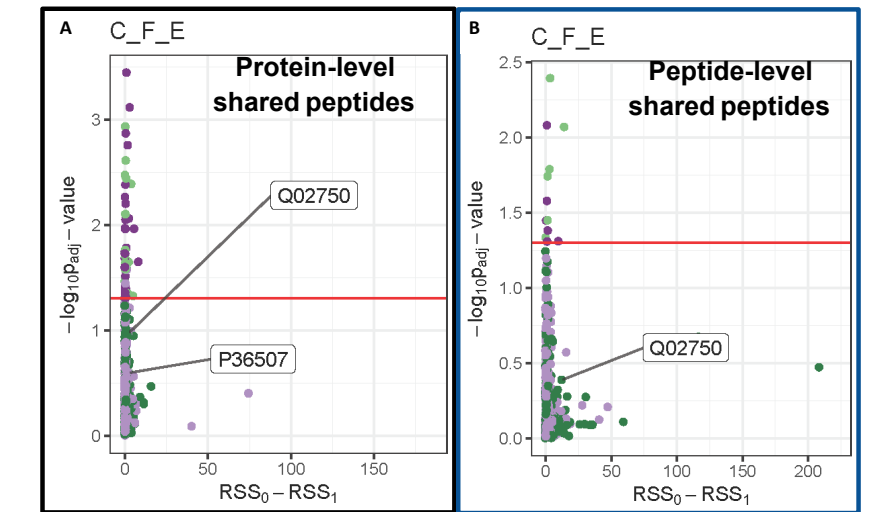


Figure 5. A) Protein-level data prior to employment of additional filters shows both stabilized (dark green) and destabilized (dark purple) shifts as proteins that have shown thermal shifting ability using the F-statistic and  $\Delta T_m$  values B) Peptide-level data using shared peptides after filtering shows that MEK2 was filtered out with the filters implemented as well as other possible targets MEK 1 and MEK 2 targets in this study are labeled and horizontal red lines depict  $-\log(0.05)$  value threshold

## Conclusions

The software pipeline developed can analyze from 1 to 8+ experiments to record overlaps as well as outliers for thermal shift assays

Novel peptide filtering approaches help increase the number of fitted curves for proteins with unique peptides by 27%

Quality control experiments aid in identification of missing values, signal-to-noise and variability to diagnose their effect on the number of fitted melt curves

## Future work

Analyze external data to evaluate potential target identification improvements

Develop receiver operating characteristic (ROC) curves to evaluate sensitivity using STRINGdb

Develop a software package in R to share with the scientific community

## References

Franken, H., Mathieson, T., Childs, D., Sweetman, G. M., Werner, T., Tögel, I.,... & Savitski, M. M. (2015). Thermal proteome profiling for unbiased identification of direct and indirect drug targets using multiplexed quantitative mass spectrometry. *Nature protocols*, 10(10), 1567-1593.

Childs, D.; Bach, K.; Franken, H.; Anders, S.; Kurzawa, N.; Bantscheff, M.; Savitski, M.; Huber, W., Non-parametric analysis of thermal proteome profiles reveals novel drug-binding proteins. *Molecular & Cellular Proteomics* 2019, mcp.TIR119.001481.

## Acknowledgements

NIH NIGMS  
R35GM1364211, R01GM120272 and R43GM117932;

NIH NCI R01CA218500; ASMS 2015 Research Award;

DFCI/NU Program in Cancer Drug Development Northeastern University, TIER1 Program;

Thermo Fisher Scientific

Alla Kloss

Harvey Lieberman

