

Prometheus Protocol PR-P-018

Incremental Cycling of Trastuzumab

Thermal unfolding of a lot of proteins can be partly reversible as long as the temperature is not increased beyond a critical value that induces non-reversible changes in the protein structure. Finding this temperature requires a different heating program than the widely used linear temperature ramp that is commonly applied to measure T_m values. In this protocol, the temperature of non-reversible unfolding of Trastuzumab (Herceptin) is determined using the *Incremental Cycle* mode of the *PR.TimeControl* software. Herceptin is a monoclonal antibody used to treat breast cancer.

thermal cycling | refolding | non-reversibility curve

A1. Target/Fluorescent Molecule

Herceptin (Trastuzumab)

A2. Molecule Class/Organism

Monoclonal antibody

A3. Sequence/Formula

Heavy chain

EVQLVESGGG LVQPGGSLRL SCAASGFNIK DTYIH^WVRQA PGKGLE^WVAR IYPTNGYTRY ADSVKGRFTI SADTSKNTAY
 LQMNSLRAED TAVYYCSR^WG GDGFYAMDY^W GQGTLVTVSS ASTKGPSVFP LAPSSKSTSG GTAALGCLVK DYFPEPVTVS
^WNSGALTS^WGV HTFPAVLQSS GLYSLSSVVT VPSSSLGTQT YICNVNHKPS NTKVDKKVEP KSCDKTHTCP PCPAPELLGG
 PSVFLFPPKP KDTLMISRTP EVTCVVDVVS HEDPEVKFN^W YVDGVEVHNA KTKPREEQYN STYRVVSVLT VLHQD^WLNGK
 EYKCKVSNKA LPAPIEKTIS KAKGQPREPQ VYTLPPSREE MTKNQVSLTC LVKGFYPSDI AVE^WESNGQP ENNYKTTTPPV
 LDSDGSFFLY SKLTVDKSR^W QQGNVFSCSV MHEALHNHYT QKSLSLSPG

Light chain

DIQMTQSPSS LSASVGDRVT ITCRASQDVN TAVAW^WYQQKP GKAPKLLIYS ASFLYSGVPS RFSGSRSGTD FTLTISLQP
 EDFATYYCQQ HYTTPPTFGQ GTKVEIKRTV AAPS^WVIFPP SDEQLKSGTA SVVCLLN^WFY PREAKVQ^WKV DNALQSGNSQ
 ESVTEQDSKD STYLSLSTLT LSKADYEKHK VYACEVTHQG LSSPVTKSFN RGEC

A4. Purification Strategy/Source

N/A

A5. Stock Concentration/Stock Buffer

120 mg/mL | 825 μ M

A6. Molecular Weight/Extinction Coefficient

145.5 kDa

A7. Dilution Buffer

Phosphate buffered saline (PBS, pH 7.4), 0.05% TWEEN® 20

D1. nanoDSF System/Capillaries

Prometheus NT.48 (NanoTemper Technologies GmbH)

High Sensitivity Capillaries Prometheus NT.48 nanoDSF Grade (PR-C006, NanoTemper Technologies GmbH)

D2. nanoDSF Software

PR.TimeControl v1.0.2 (NanoTemper Technologies GmbH)

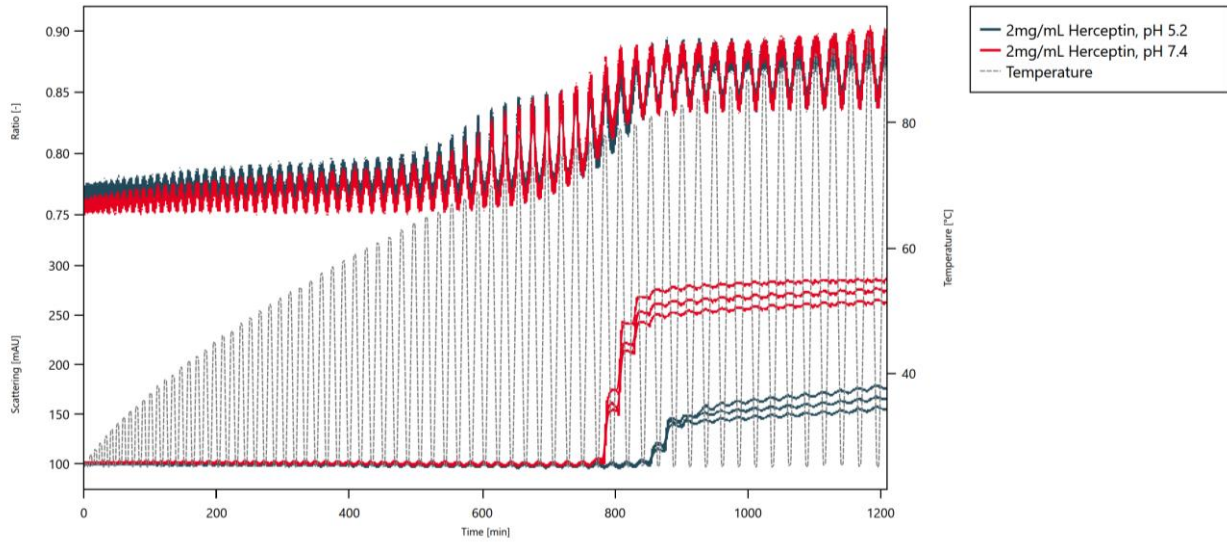
nanotempertech.com/prometheus-software

D3. nanoDSF Experiment

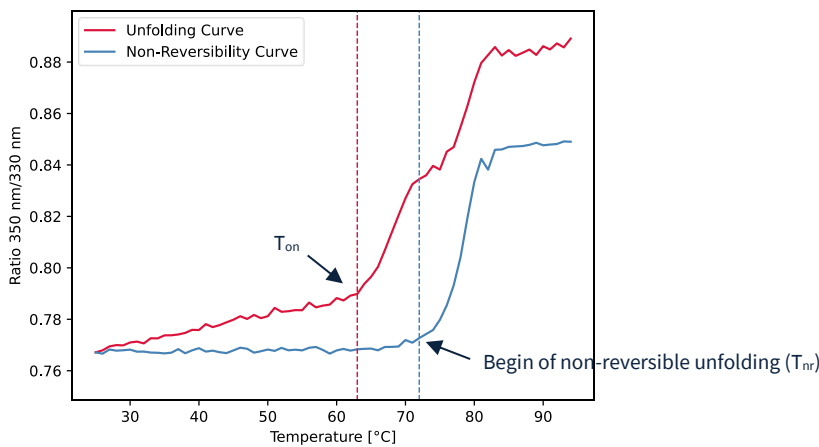
1. Prepare a 20 mg/mL solution of Trastuzumab in PBS buffer.
2. Mix 4 µL of the 20 mg/mL Trastuzumab solution with 36 µL of 20 mM Na-Acetate buffer, pH 5.2, to obtain 40 µL of a 2 mg/mL Trastuzumab pH 5.2 solution.
3. Then, mix 4 µL of the 20 mg/mL Trastuzumab solution with 36 µL of PBS buffer, pH 7.4, to obtain 40 µL of a 2 mg/mL Trastuzumab pH 7.4 solution.
4. Start a new session of the *PR.TimeControl* software.
5. Change the Thermostat Set Point to 25°C by using the touch display on the instrument and wait for the system to reach this temperature.
6. Go to 'Measurement Scan' and prepare a run with the following settings:
 - a. Capillaries 1 – 6 selected
 - b. Incremental Cycle
 - c. First Temperature 25°C
 - d. First Duration 2 min
 - e. Temperature Increase 1°C
 - f. Second Duration 2 min
 - g. Measurement Cycles 70
 - h. 5% excitation power
7. Load 3 capillaries from the 2 mg/mL Trastuzumab **pH 5.2** solution and place them on positions **1 – 3** of the Prometheus capillary tray.
8. Load 3 capillaries from the 2 mg/mL Trastuzumab **pH 7.4** solution and place them on positions **4 – 6** of the Prometheus capillary tray.
9. Place the magnetic lid to fix the capillary.
10. Start the measurement (duration: **~24 hours**).

D4. nanoDSF Results

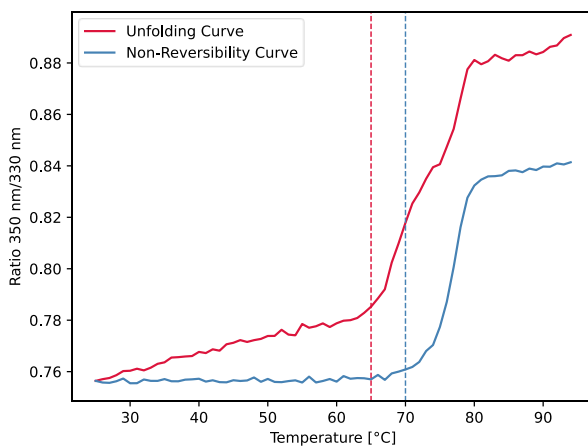
Data from *PR.TimeControl*:



Separating the data into ratio values at ‘hot’ (unfolding curve) and ‘cold’ (non-reversibility curve)¹:



pH 5.2



pH 7.4

¹ See also [Svilenov et al., Mol. Pharmaceutics 2020, 17, 7, 2638–2647](#) for more information on data analysis.

Trastuzumab	pH 5.2	pH 7.4
T_{on} (Onset temperature)	~63°C	~65°C
T_{nr} (Non-reversibility onset temperature)	~72°C	~70°C
T_{agg} (Aggregation onset temperature)	~81°C	~78°C

While Trastuzumab begins to unfold at a slightly lower temperature at pH 5.2 than at pH 7.4 (~63°C vs. ~65°C), the temperature of the first non-reversible unfolding is in fact higher (~72°C vs. ~70°C), which also correlates with the non-reversible aggregation information from the scattering signal (~81°C vs. ~78°C).

D5. Reference Results/Supporting Results

Svilenov et al., *Mol. Pharmaceutics* 2020, 17, 7, 2638–2647

E. Contributors

Andreas Langer²

² NanoTemper Technologies GmbH, München, Germany | nanotempertech.com