

Prometheus Protocol PR-P-019

# Chemical Unfolding of a Monoclonal Antibody

Herceptin (Trastuzumab) is a monoclonal antibody used to treat breast cancer. In this protocol, the chemical stability is determined using a chemical unfolding measurement with PR.ChemControl.

chemical unfolding | guanidine hydrochloride |  $\Delta G$  |  $C_{50}$

---

## A1. Target/Fluorescent Molecule

Herceptin (Trastuzumab)

## A2. Molecule Class/Organism

Monoclonal antibody

## A3. Sequence/Formula

Heavy chain

EVQLVESGGG LVQPGGSLRL SCAASGFNIK DTYIH<sup>W</sup>VRQA PGKGLE<sup>W</sup>VAR IYPTNGYTRY ADSVKGRFTI SADTSKNTAY  
 LQMNSLRAED TAVYYCSR<sup>W</sup>G GDGFYAMDY<sup>W</sup> GQGT<sup>L</sup>LVTVSS ASTKGPSVFP LAPSSKSTSG GTAALGCLVK DYFPEPVTVS  
<sup>W</sup>NSGALTS<sup>G</sup>V HTFPAVLQSS GLYSLSSVVT VPSSSLGTQT YICNVNHKPS NTKVDKKVEP KSCDKTHTCP PCPAPELLGG  
 PSVFLFPPKP KDTLMISRTP EVTCVVDVDS HEDPEVKFN<sup>W</sup> YVDGVEVHNA KTKPREEQYN STYRVVSVLT VLHQD<sup>W</sup>LNGK  
 EYKCKVSNKA LPAPIEKTIS KAKGQPREPQ VYTLPPSREE MTKNQVSLTC LVKGFYPSDI AVE<sup>W</sup>ESNGQP ENNYKTTTPV  
 LDSDGSFFLY SKLTVDKSR<sup>W</sup> QQGNVFSCSV MHEALHNHYT QKSLSLSPG

Light chain

DIQMTQSPSS LSASVGDRTV ITCRASQDVN TAVAW<sup>Y</sup>QQKP GKAPKLLIYS ASFLYSGVPS RFSGSRSGTD FTLTISLQP  
 EDFATYQCQQ HYTTPPTFGQ GTKVEIKRTV AAPS<sup>V</sup>IFPP SDEQLKSGTA SVVCLLN<sup>N</sup>FY PREAKVQ<sup>W</sup>KV DNALQSGNSQ  
 ESVTEQDSKD STYLSLSTLT LSKADYEKHK VYACEVTHQG LSSPVTKSFN RGEK

## A4. Purification Strategy/Source/Batch-No.

N/A

## A5. Stock Concentration/Stock Buffer

120 mg/mL | 825  $\mu$ M

## A6. Molecular Weight/Extinction Coefficient

145.5 kDa  
 225,000  $M^{-1}cm^{-1}$  ( $\epsilon_{280}$ )

## A7. Dilution Buffer

Phosphate-buffered saline (PBS), pH 7.4

## D1. nanoDSF System/Capillaries

Prometheus NT.48 (NanoTemper Technologies GmbH)

Standard Capillaries Prometheus NT.48 nanoDSF Grade (PR-C002, NanoTemper Technologies GmbH)

## D2. nanoDSF Software

PR.ChemControl v1.4 (NanoTemper Technologies GmbH)

[nanotempertech.com/prometheus-software](http://nanotempertech.com/prometheus-software)

## D3. nanoDSF Experiment

1. Add 145  $\mu\text{L}$  of 1x PBS to 20  $\mu\text{L}$  of 825  $\mu\text{M}$  Herceptin to obtain 165  $\mu\text{L}$  of a 100  $\mu\text{M}$  Herceptin solution.
2. Prepare 1 mL of a 8 M solution of Guanidine Hydrochloride (GuHCl) in ddH<sub>2</sub>O.
3. Prepare two 1.5 mL tubes and label them with **A** and **B**.
4. Solution **A** – 1280  $\mu\text{L}$  of a 5  $\mu\text{M}$  Herceptin solution in **5 M GuHCl**, 1x PBS:  
*Mix 800  $\mu\text{L}$  of 8 M GuHCl with 200  $\mu\text{L}$  of a 5x PBS stock solution. Then, add 216  $\mu\text{L}$  of 1x PBS and 64  $\mu\text{L}$  of the 100  $\mu\text{M}$  Herceptin solution. Mix carefully using a pipette.*
5. Solution **B** – 1280  $\mu\text{L}$  of a 5  $\mu\text{M}$  Herceptin solution in **0.2 M GuHCl**, 1x PBS:  
*Mix 32  $\mu\text{L}$  of 8 M GuHCl with 768  $\mu\text{L}$  of ddH<sub>2</sub>O and 200  $\mu\text{L}$  of a 5x PBS stock solution. Then, add 216  $\mu\text{L}$  of 1x PBS and 64  $\mu\text{L}$  of the 100  $\mu\text{M}$  Herceptin solution. Mix carefully using a pipette.*
6. Prepare 48 small PCR tubes according to the following table:

Tube #	1	2	3	4	5	6	7	8	9	10	11	12
Solution A ( $\mu\text{L}$ ):	48	47	46	45	44	43	42	41	40	39	38	37
Solution B ( $\mu\text{L}$ ):	0	1	2	3	4	5	6	7	8	9	10	11
Final GuHCl (M):	5	4.9	4.8	4.7	4.6	4.5	4.4	4.3	4.2	4.1	4.0	3.9

Tube #	13	14	15	16	17	18	19	20	21	22	23	24
Solution A ( $\mu\text{L}$ ):	36	35	34	33	32	31	30	29	28	27	26	25
Solution B ( $\mu\text{L}$ ):	12	13	14	15	16	17	18	19	20	21	22	23
Final GuHCl (M):	3.8	3.7	3.6	3.5	3.4	3.3	3.2	3.1	3.0	2.9	2.8	2.7

Tube #	25	26	27	28	29	30	31	32	33	34	35	36
Solution A ( $\mu\text{L}$ ):	24	23	22	21	20	19	18	17	16	15	14	13
Solution B ( $\mu\text{L}$ ):	24	25	26	27	28	29	30	31	32	33	34	35
Final GuHCl (M):	2.6	2.5	2.4	2.3	2.2	2.1	2.0	1.9	1.8	1.7	1.6	1.5

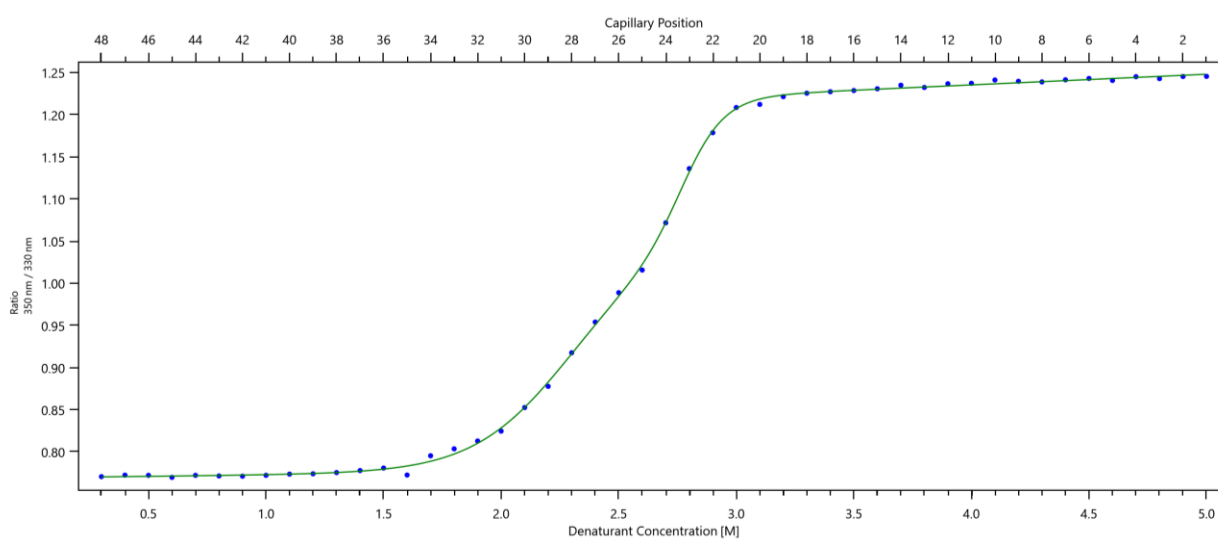
Tube #	37	38	39	40	41	42	43	44	45	46	47	48
Solution A ( $\mu\text{L}$ ):	12	11	10	9	8	7	6	5	4	3	2	1
Solution B ( $\mu\text{L}$ ):	36	37	38	39	40	41	42	43	44	45	46	47
Final GuHCl (M):	1.4	1.3	1.2	1.1	1.0	0.9	0.8	0.7	0.6	0.5	0.4	0.3

7. Mix well and incubate for at least 48 hours<sup>1</sup> at room temperature to ensure equilibrium.
8. After incubation, load 48 capillaries from the tubes and place them on positions 1 to 48 of the capillary tray according to the vial number. Place the magnetic lid to fix the capillaries.
9. Start a new session of the *PR.ChemControl* software.
10. Go to 'Plan' and prepare a run with the following settings:
  - a. Start concentration: 5 M
  - b. End concentration: 0.3 M
11. Start the measurement.
12. Use the 'Three-State' fit to analyze the data.

#### D4. nanoDSF Results

$\Delta G_1 = 28.0 \text{ kJ/mol}$  |  $c_{50_1} = 2.3 \text{ M}$

$\Delta G_2 = 74.1 \text{ kJ/mol}$  |  $c_{50_2} = 2.8 \text{ M}$



#### D5. Reference Results/Supporting Results

N/A

#### E. Contributors

Andreas Langer<sup>2</sup>

<sup>1</sup> Large biomolecules such as monoclonal antibodies often require a very long incubation time to equilibrate sufficiently and at least overnight incubation (better 24 – 72 hours) is recommended. Performing the measurement too early will lead to incorrect  $\Delta G$  and  $c_{50}$  values.

<sup>2</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](http://nanotempertech.com)