

Prometheus Protocol PR-P-003

Folding Kinetics of Lysozyme

Lysozyme is an enzyme that prevents bacterial infections by attacking peptidoglycan, a component of certain bacterial cell walls. Peptidoglycan is composed of the repeating amino sugars N-acetylglucosamine (NAG) and N-acetylmuramic acid (NAM), which are crosslinked by peptide bridges. Lysozyme hydrolyzes the bond between NAG and NAM, increasing the bacteria's permeability and causing the bacteria to burst. It is widely distributed in plants and animals. The majority of the lysozyme used in research is purified from hen egg whites.

[chemical unfolding](#) | [guanidine hydrochloride](#) | [unfolding kinetics](#) | [refolding kinetics](#) | ΔG | c_{50}

A1. Target/Fluorescent Molecule

Lysozyme

uniprot.org/uniprot/B8YK79

A2. Molecule Class/Organism

Glycoside hydrolases

Gallus gallus (Chicken)

A3. Sequence/Formula

KVFGRCELAA AMKRHGLDNY RGYSLGNWVC AAKFESNFNT QATNRNTDGS TDYGILQINS RWWCNDGRTP GSRNLCNIPC
SALLSSDITA SVNCAKKIIVS DGNMNAWVA WRNRCKGTDV QAWIRGCRL

A4. Purification Strategy/Source

Sigma-Aldrich GmbH

[L6876](#)

A5. Stock Concentration/Stock Buffer

32 μg lyophilized powder

A6. Molecular Weight/Extinction Coefficient

14.3 kDa

37,970 $\text{M}^{-1}\text{cm}^{-1}$ (ϵ_{280})

A7. Dilution Buffer

50 mM Tris-HCl, pH 7.8, 150 mM NaCl, 10 mM MgCl_2 , 0.05% TWEEN® 20

D1. nanoDSF System/Capillaries

Prometheus NT.48 (NanoTemper Technologies GmbH)

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

or

Prometheus NT.Plex (NanoTemper Technologies GmbH)

Standard 24-Capillary Chips Prometheus NT.Plex nanoDSF Grade (PR-AC002, NanoTemper Technologies GmbH)

D2. nanoDSF Software

PR.TimeControl v1.0.2 (NanoTemper Technologies GmbH)

nanotempertech.com/prometheus-software

D3. nanoDSF Experiment

1. Dissolve 764 mg of guanidine hydrochloride (Gua-HCl, MW 95.53) in 1 mL of dilution buffer to obtain an 8 M solution. Heat the solution until all Gua-HCl is dissolved.

Unfolding

2. Prepare 12 small PCR tubes according to the following table:

Tube	1	2	3	4	5	6	7	8	9	10	11	12
8 M Gua-HCl (μL):	140	130	120	110	100	90	80	70	65	60	55	50
Dilution buffer (μL):	10	20	30	40	50	60	70	80	85	90	95	100

3. Resuspend 32 μg lysozyme in 32 μL dilution buffer to obtain a 1 mg/mL solution.
4. Prepare 7 new tubes with 2 μL of lysozyme solution in each.
5. Start a new session of the *PR.TimeControl* software.
6. Go to 'Measurement Scan' and prepare a run with the following settings:
 - a. Only capillary 1 selected
 - b. Isothermal
 - c. 25°C
 - d. 10 min
 - e. 100% excitation power

The following steps should happen **as fast as possible**:

7. Add 30 μL of the solution from tube **1** of step 2 to one of the lysozyme vials from step 4 to obtain 32 μL of a **7 M** GuaHCl solution (this will start the **unfolding** process). Quickly mix 4 – 5 times by pipetting.¹
8. Load one capillary from this solution and place it on position 1 of the Prometheus capillary tray.
9. Place the magnetic lid to fix the capillary.
10. Start the measurement.
11. Repeat these steps for tubes **2 – 7** to obtain unfolding curves for **6.5 M, 6 M, 5.5 M, 5 M, 4.5 M, 4 M** Gua-HCl.

¹ When using the Prometheus NT.Plex, GuaHCl solution can be added to multiple tubes at once using a multichannel pipette. The solutions can then be loaded into a capillary chip and all unfolding/refolding kinetics can be recorded simultaneously.

Refolding

12. Resuspend new 32 μg lysozyme in 32 μL of the 8 M Guanidine hydrochloride solution from step 1 to obtain a 1 mg/mL solution and **incubate** for **10 minutes**.
13. Prepare 5 new tubes with 2 μL of lysozyme solution in each.
14. Go to 'Measurement Scan' and prepare a run with the following settings:
 - a. Only capillary 1 selected
 - b. Isothermal
 - c. 25°C
 - d. 10 min
 - e. 100% excitation power

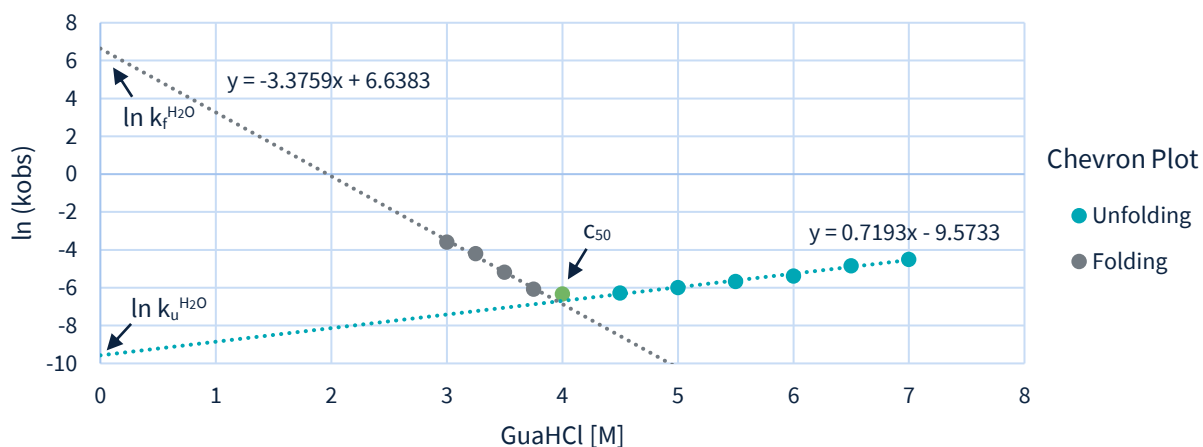
The following steps should happen **as fast as possible**:

15. Add 30 μL of the solution from tube **8** of step 2 to one of the lysozyme vials from step 13 obtain 32 μL of a **4 M** GuaHCl solution (this will start the **refolding** process). Quickly mix 4 – 5 times by pipetting.
16. Load one capillary from this solution and place it on position 1 of the Prometheus capillary tray.
17. Place the magnetic lid to fix the capillary.
18. Start the measurement.

Repeat these steps for tube **9 – 11** to obtain refolding curves for **3.75 M, 3.5 M, 3.25 M, 3 M** Gua-HCl.

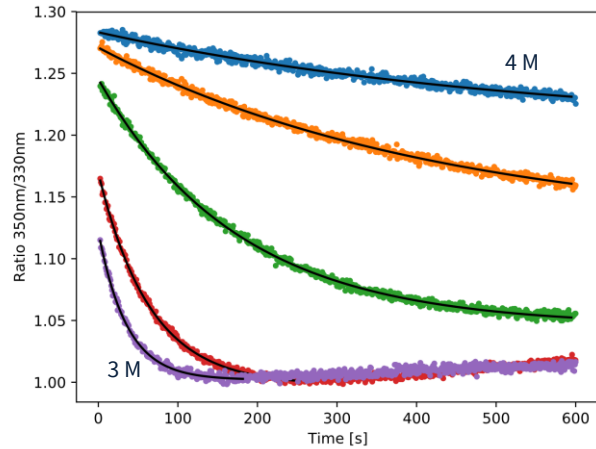
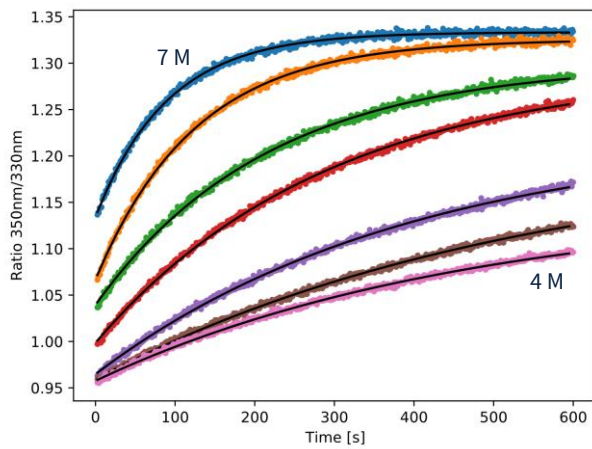
D4. nanoDSF Results

$$k_f^{\text{H}_2\text{O}} = 762 \text{ s}^{-1} \mid k_u^{\text{H}_2\text{O}} = 5.3 \cdot 10^{-5} \text{ s}^{-1} \mid \Delta G = RT \cdot [\ln(k_f^{\text{H}_2\text{O}}) - \ln(k_u^{\text{H}_2\text{O}})] = 40.8 \text{ kJ/mol} \mid c_{50} \sim 4 \text{ M}$$



The denaturation midpoint c_{50} is defined as the denaturant concentration at which both folded and unfolded states are equally populated at equilibrium.

Gua-HCl (M)	7	6.5	6	5.5	5	4.5	4	3.75	3.5	3.25	3
$k_{\text{obs}} (10^{-3} \text{ s}^{-1})$	11.1	8.0	4.6	3.4	2.5	1.9	1.8	2.3	5.7	15.1 ²	27.7 ¹



D5. Reference Results/Supporting Results

$\Delta G = 37.2 \text{ kJ/mol}$ | $c_{50} = 4.2 \text{ M}$

Absorbance measurements

[Ahmand et al., Biochem J 287 \(1992\) 481-485](#)

$k_f^{\text{H}_2\text{O}} = 447 \text{ s}^{-1}$ | $k_u^{\text{H}_2\text{O}} = 6.2 \cdot 10^{-7} \text{ s}^{-1}$

Tryptophan fluorescence

[Kiefhaber et al., PNAS 92 \(1995\) 9029-9033](#)

E. Contributors

Franziska Toppel³, Andreas Langer³

² Under strongly native conditions, lysozyme refolds on parallel pathways, resulting in double-exponential refolding kinetics ([Kiefhaber et al., PNAS 92 \(1995\) 9029-9033](#)).

³ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com