

Prometheus Protocol PR-P-001

# Thermal Unfolding 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.

thermal unfolding | conformational stability | colloidal stability | aggregation |  $T_m$  |  $T_{on}$  |  $T_{agg}$

---

## A1. Target/Fluorescent Molecule

Lysozyme

[uniprot.org/uniprot/B8YK79](https://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 M<sup>-1</sup>cm<sup>-1</sup> ( $\epsilon_{280}$ )

## A7. Dilution Buffer

50 mM Tris-HCl, pH 7.8, 150 mM NaCl, 10 mM MgCl<sub>2</sub>, 0.05% TWEEN® 20

## D1. nanoDSF System/Capillaries

Prometheus NT.48 (NanoTemper Technologies GmbH)

Prometheus Aggregation Detection Optics (PR-AGO, NanoTemper Technologies GmbH)

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

## D2. nanoDSF Software

PR.ThermControl v2.1 (NanoTemper Technologies GmbH)

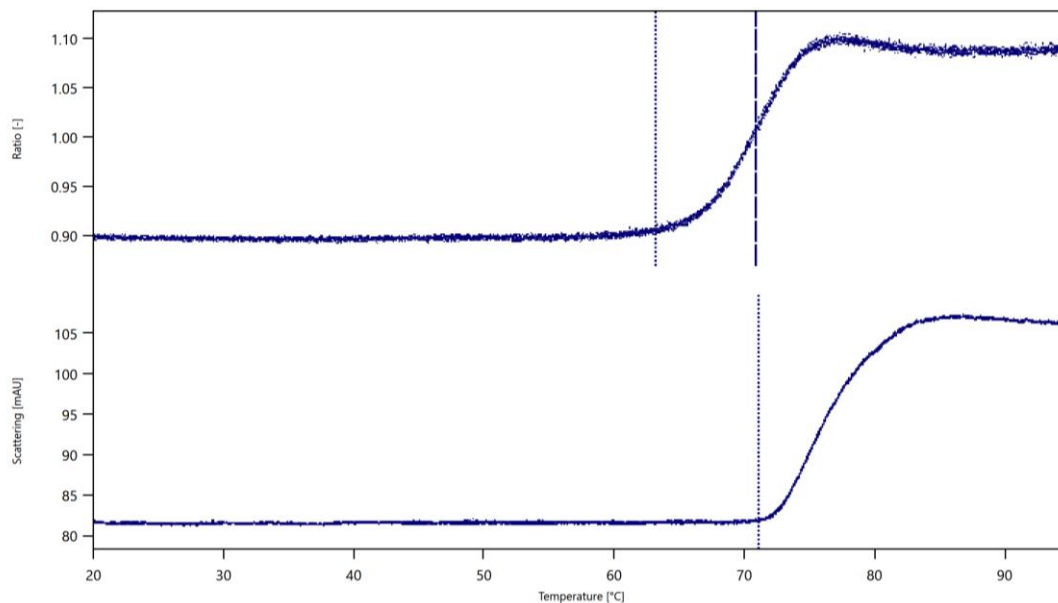
[nanotempertech.com/prometheus-software](http://nanotempertech.com/prometheus-software)

## D3. nanoDSF Experiment

1. Resuspend lysozyme in 32  $\mu\text{L}$  of dilution buffer to obtain a 1 mg/mL solution.
2. Completely fill one capillary from this solution, place it on position 1 of the capillary tray and place the magnetic lid to fix the capillary.
3. Start a new session of the *PR.ThermControl* software.
4. Go to 'Melting Scan' and prepare a run with the following settings:
  - a. Only capillary 1 selected
  - b. 1.0°C/min
  - c. 20°C – 95°C
  - d. 20% excitation power
5. Start the measurement.

## D4. nanoDSF Results

$T_m = 70.9^\circ\text{C}$  |  $T_{on} = 63.2^\circ\text{C}$  |  $T_{agg} = 71.1^\circ\text{C}$



## D5. Reference Results/Supporting Results

$T_m = 70.3^\circ\text{C}$	FTIR spectroscopy <a href="#">Venkataramani et al., J Pharm Bioallied Sci. 2 (2013) 148–153</a>
$T_m = 71.7^\circ\text{C}$	Differential Scanning Calorimetry (DSC) <a href="#">James et al., J Phys Chem B 116 (2012) 10182–10188</a>

---

## E. Contributors

Andreas Langer<sup>1</sup>

---

<sup>1</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](http://nanotempertech.com)