

Monolith Protocol MO-P-060

Lysozyme – NAG₃ (no column purification)

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. Tri-N-acetyl-D-glucosamine (NAG₃) is an inhibitor of lysozyme.

protein – small molecule | carbohydrate

A1. Target/Fluorescent Molecule

Lysozyme

uniprot.org/uniprot/B8YK79

A2. Molecule Class/Organism

Glycoside hydrolase

Gallus gallus (Chicken)

A3. Sequence/Formula

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

A4. Purification Strategy/Source

Sigma-Aldrich GmbH

[L6876](#)

A5. Stock Concentration/Stock Buffer

128 µg lyophilized powder

A6. Molecular Weight/Extinction Coefficient

14.3 kDa

37,970 M⁻¹cm⁻¹ (ε₂₈₀)

A7. Dilution Buffer

Phosphate-buffered saline (PBS), pH 7.4, 0.01% Pluronic® F-127

A8. Labeling Strategy

Monolith Protein Labeling Kit RED – NHS 2nd Generation (MO-L011, NanoTemper Technologies GmbH)
1* Dye RED-NHS 2nd Generation (10 µg)

A9. Labeling Procedure

1. Resuspend 128 µg of lysozyme in 43 µL of dilution buffer to obtain a ~208 µM lysozyme solution.
2. Add 25 µL of DMSO to Dye RED-NHS 2nd Generation (10 µg) to obtain a ~600 µM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
3. Transfer 1 µL of the ~600 µM dye solution into a new PCR tube.
4. Add 29 µL of the ~208 µM lysozyme solution to the tube and mix well (30 µL final volume, ~3% final DMSO concentration, ~200 µM final lysozyme concentration, ~20 µM final dye concentration¹).
5. Incubate for 20 minutes at room temperature in the dark.

A10. Labeling Efficiency

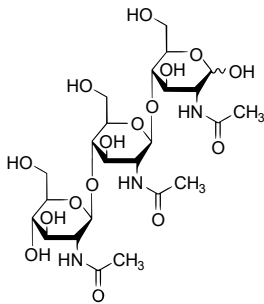
N/A²

¹ The excess of protein over dye assures that a maximum of dye can react with protein before hydrolysis.

² No purification from unreacted dye.

B1. Ligand/Non-Fluorescent Binding Partner

Tri-N-acetyl-D-glucosamine (NAG₃)



B2. Molecule Class/Organism

Carbohydrate

B3. Sequence/Formula

C₂₄H₄₁N₃O₁₆

B4. Purification Strategy/Source

Sigma-Aldrich GmbH

[T2144](#)

B5. Stock Concentration/Stock Buffer

25 µg powder

B6. Molecular Weight/Extinction Coefficient

627.59 Da

B7. Serial Dilution Preparation

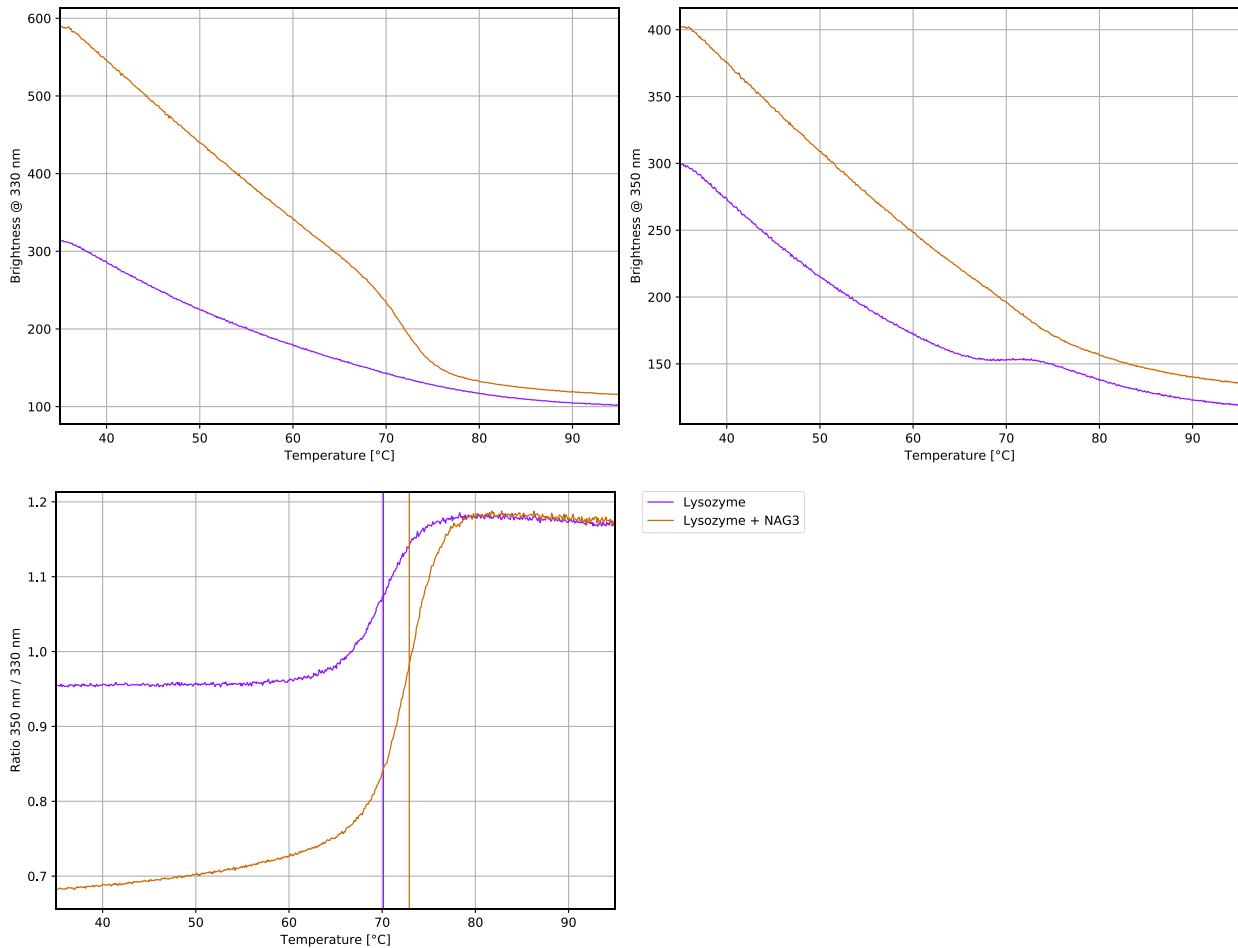
1. Resuspend NAG₃ in 25 µL of dilution buffer to obtain a 1.6 mM solution.
2. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 µL of the 1.6 mM NAG₃ solution into tube **1**. Then, transfer 10 µL of dilution buffer into tubes **2** to **16**.
3. Prepare a 1:1 serial dilution by transferring 10 µL from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10 µL from tube **16** to get an equal volume of 10 µL for all samples.
4. Mix 1 µL of labeled lysozyme (~200 µM) with 199 µL of dilution buffer to obtain 200 µL of ~1 µM labeled lysozyme solution.
5. Add 10 µL of labeled lysozyme (~1 µM) to each tube from **16** to **1** and mix by pipetting.
6. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

C. Applied Quality Checks

Validation of structural integrity and functionality of lysozyme using Tycho NT.6:

nanotempertech.com/tycho

Lysozyme	6 μ L of \sim 208 μ M lysozyme + 4 μ L of dilution buffer	$T_i = 70.1^\circ\text{C}$
Lysozyme + NAG3	6 μ L of \sim 208 μ M lysozyme + 4 μ L of 1.6 mM NAG ₃	$T_i = 72.9^\circ\text{C}$



D1. MST System/Capillaries

Monolith NT.115 Red (NanoTemper Technologies GmbH)

Capillaries Monolith NT.115 (MO-K022, NanoTemper Technologies GmbH)

D2. MST Software

MO.Control v1.6 (NanoTemper Technologies GmbH)

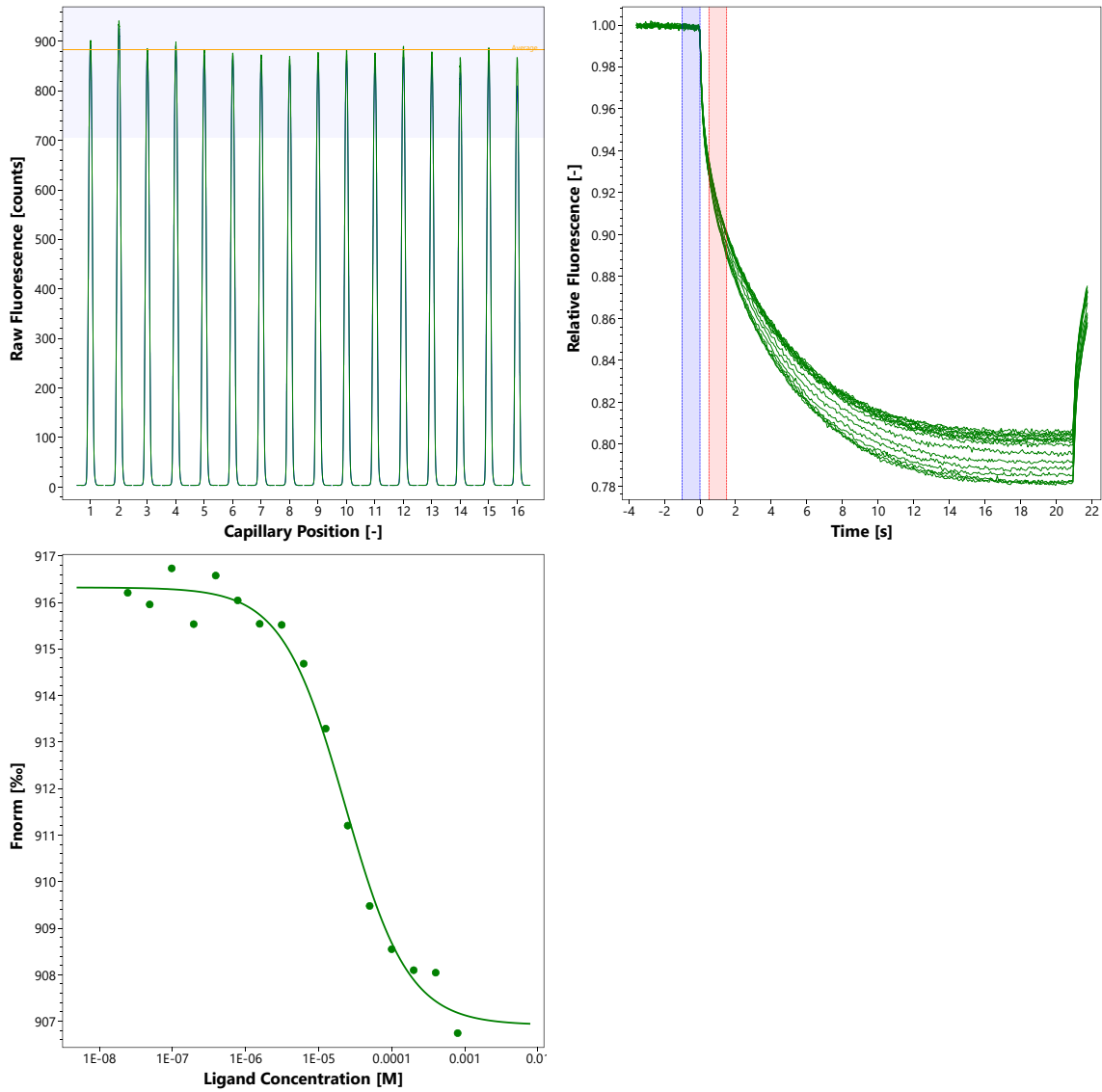
nanotempertech.com/monolith-mo-control-software

D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

Phosphate-buffered saline (PBS), pH 7.4, 0.01% Pluronic® F-127
 500 nM lysozyme (~50 nM dye) | 800 μ M – 24 nM NAG₃ | 22°C | medium MST Power | 20% excitation power

D4. MST Results (Capillary Scan/Time Traces/Dose Response)

$K_d = 23 \mu$ M



D5. Reference Results/Supporting Results

- $K_d = 39.8 \mu\text{M}$ NanoESI
[Jecklin et al., J Am Soc Mass Spectrom 19 \(2008\) 332–343](#)
- $K_d = 39 \mu\text{M}$ Microchip-ESI-MS
[Svobodova et al., Electrophoresis 31 \(2010\) 2680–2685](#)

E. Contributors

Andreas Langer³

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