

## Monolith Protocol MO-P-009

# Lysozyme – NAG<sub>3</sub>

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 (NAG<sub>3</sub>) is an inhibitor of lysozyme.

protein – small molecule interaction | carbohydrate

### A1. Target/Fluorescent Molecule

Lysozyme

[uniprot.org/uniprot/B8YK79](https://uniprot.org/uniprot/B8YK79)

### A2. Molecule Class/Organism

Glycoside hydrolase

*Gallus gallus* (Chicken)

### A3. Sequence/Formula

KVFGRCELAA AMKRHGLDNY RGYSLGNVVC AAKFESNFNT QATNRNTDGS TDYGILQINS RWWCNDGRTP GSRNLCNIPC  
SALLSSDITA SVNCAKKIVS DNGMNAWVA 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> (ε<sub>280</sub>)

### 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\* Labeling Buffer NHS | 1\* Dye RED-NHS 2nd Generation (10 µg) | 1\* B-Column

## A9. Labeling Procedure

1. Resuspend 32 µg lysozyme in 112 µL of Labeling Buffer NHS to obtain a 20 µM solution.
2. Transfer 80 µL of the 20 µM lysozyme solution into a new tube.
3. Add 25 µL of DMSO to 10 µg RED-NHS 2nd Generation dye to obtain a ~600 µM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
4. Mix 10 µL of the 600 µM dye solution with 70 µL of Labeling Buffer NHS to obtain 80 µL of a 75 µM dye solution (~4x protein concentration).
5. Mix lysozyme and dye in a 1:1 volume ratio (160 µL final volume, ~6% final DMSO concentration).
6. Incubate for 20 minutes at room temperature in the dark.
7. In the meantime, remove the top cap of the B-Column and pour off the storage solution. Remove the bottom cap and place with adapter in a 15 mL tube.
8. Fill the column with dilution buffer and allow it to enter the packed resin bed completely by gravity flow. Discard the flow through collected. Repeat this step 3 more times.
9. Add 160 µL of the labeling reaction from step 5 to the center of the column and let sample enter the resin bed completely.
10. Add 500 µL of dilution buffer after the sample has entered and discard the flow through.
11. Place column in a new collection tube, add 400 µL of dilution buffer and collect the eluate.
12. Keep the labeled lysozyme (~4 µM) on ice in the dark.

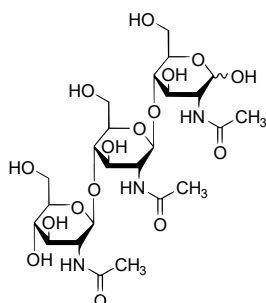
## A10. Labeling Efficiency

Measurement of protein concentration and degree of labeling (DOL) using a NanoDrop™:  
[nanotempertech.com/dol-calculator](https://nanotempertech.com/dol-calculator)

Absorbance A <sub>280</sub>	0.117	Protein concentration	~2.57 µM
Absorbance A <sub>650</sub>	0.490	Degree-of-labeling (DOL)	~0.98

## B1. Ligand/Non-Fluorescent Binding Partner

Tri-N-acetyl-D-glucosamine (NAG<sub>3</sub>)



## B2. Molecule Class/Organism

Carbohydrate

## B3. Sequence/Formula

C<sub>24</sub>H<sub>41</sub>N<sub>3</sub>O<sub>16</sub>

## 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

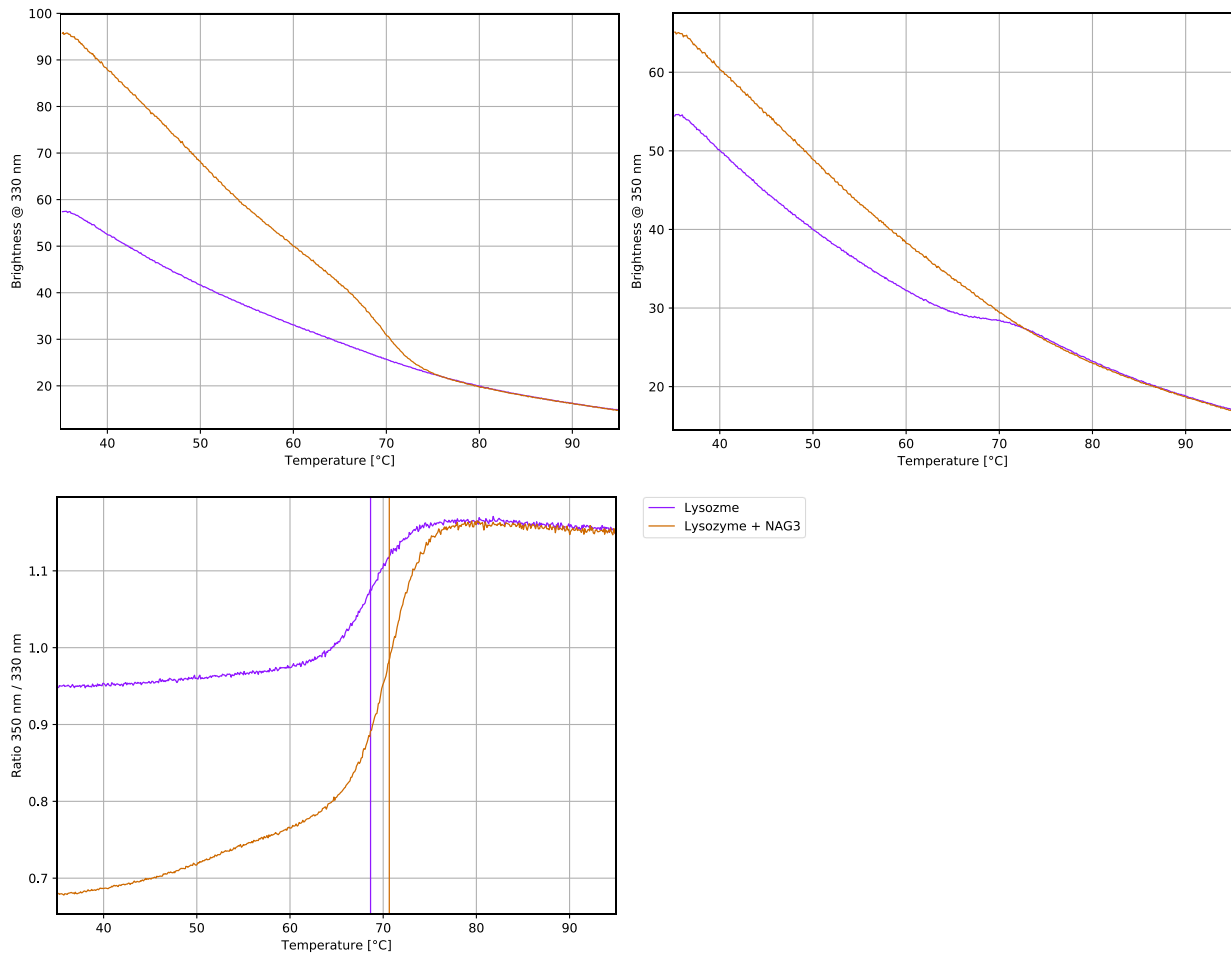
1. Resuspend NAG<sub>3</sub> in 25 µL of dilution buffer to obtain a 1.6 mM solution.
2. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 µL of the 1.6 mM NAG<sub>3</sub> solution into tube **1**. Then, transfer 10 µL of dilution buffer into tubes **2** to **16**.
3. Prepare a 1:1 serial dilution by 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 4 µL of labeled lysozyme (~4 µM) with 196 µL of dilution buffer to obtain 200 µL of a ~100 nM lysozyme solution.
5. Add 10 µL of labeled lysozyme (~100 nM) 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](https://nanotempertech.com/tycho)

Lysozyme	8 $\mu$ L of 20 $\mu$ M lysozyme + 2 $\mu$ L of dilution buffer	$T_i = 68.6^\circ\text{C}$
Lysozyme + NAG <sub>3</sub>	8 $\mu$ L of 20 $\mu$ M lysozyme + 2 $\mu$ L of 1.6 mM NAG <sub>3</sub>	$T_i = 70.7^\circ\text{C}$



## 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](https://nanotempertech.com/monolith-mo-control-software)

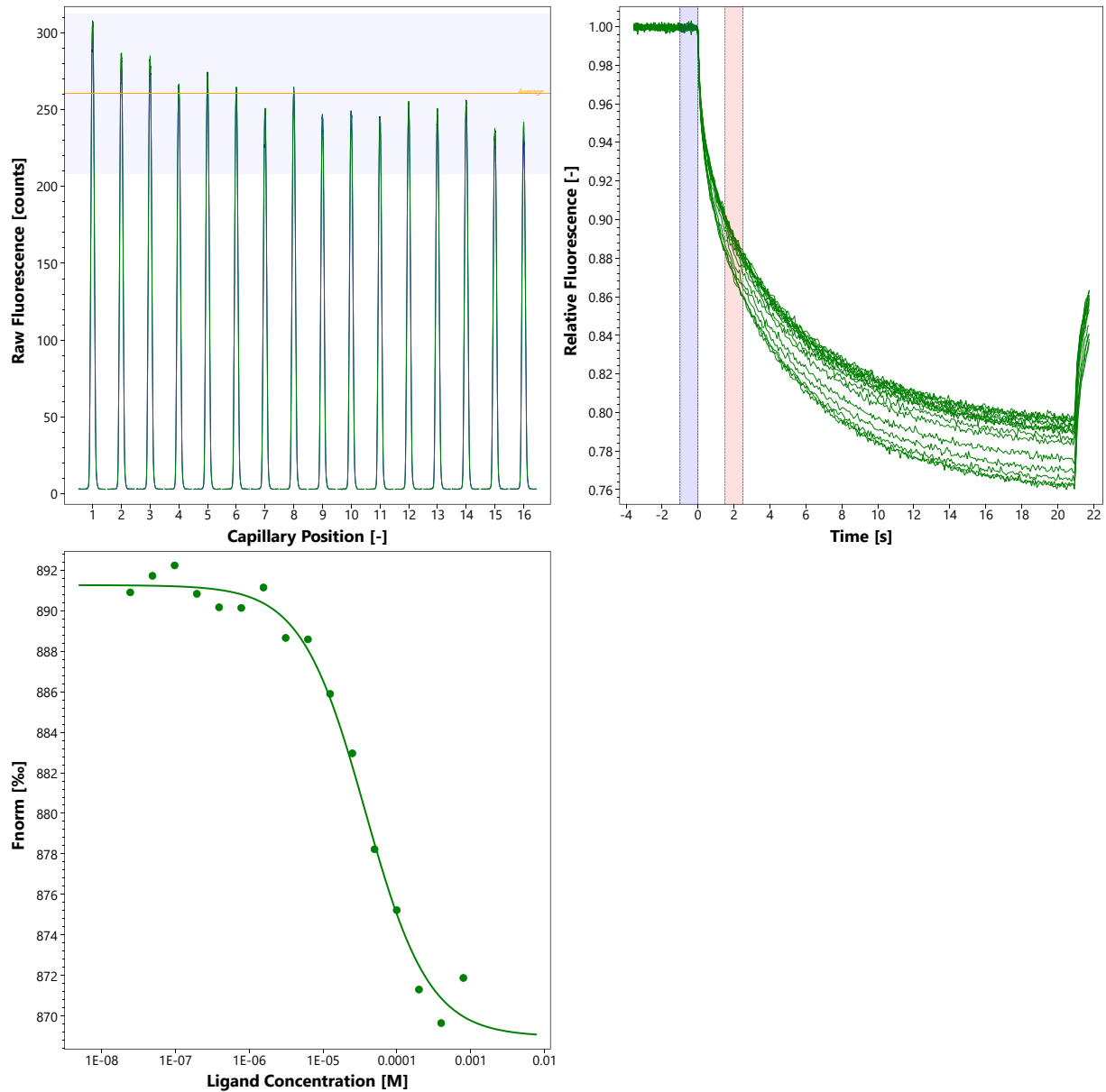
### D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

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

50 nM lysozyme | 800  $\mu$ M – 24 nM NAG<sub>3</sub> | 22°C | medium MST Power | 20% excitation power

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

$K_d = 37.3 \mu$ M



## D5. Reference Results/Supporting Results

$K_d = 39.8 \mu\text{M}$	NanoESI <a href="#">Jecklin et al., J Am Soc Mass Spectrom 19 (2008) 332–343</a>
$K_d = 39 \mu\text{M}$	Microchip-ESI-MS <a href="#">Svobadova et al., Electrophoresis 31 (2010) 2680–2685</a>

## E. Contributors

Andreas Langer<sup>1</sup>

---

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