

Monolith Protocol M0-P-071

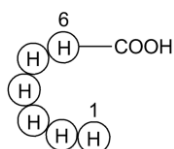
Zn(II)-Mediated Dimerization of His-Tags

The tris-NTA/His-tag system comprises one of the smallest high-affinity recognition elements known to date. This interaction is based on the capacity of the histidine's imidazole groups to form coordinative bonds with transition metal ions such as Ni(II). The presence of high concentrations of other transition metals such as Co(II), Cu(II) or Zn(II), however, can interfere with the tris-NTA/His-tag interaction. Zn(II), for example, has been described to mediate the dimerization of His-tagged proteins at low micromolar concentrations.

peptide – ion interaction | dimerization | His₆-tag | competitive assay

A1. Target/Fluorescent Molecule

His₆-peptide



A2. Molecule Class/Organism

Peptide

A3. Sequence/Formula

HHHHHH

A4. Purification Strategy/Source

APExBIO

A6006

A5. Stock Concentration/Stock Buffer

5 mg

Lyophilized powder

A6. Molecular Weight/Extinction Coefficient

840.8 Da

45,100 M⁻¹cm⁻¹ (ε₂₀₅)¹

¹ Calculated from amino acid sequence (Anthis et al., Protein Science 2013, 22, 851–858).

A7. Dilution Buffer

20 mM HEPES, pH 7.4, 150 mM NaCl, 0.01% Pluronic® F-127

A8. Labeling Strategy

Monolith His-Tag Labeling Kit RED-tris-NTA 2nd Generation (MO-L018, NanoTemper Technologies GmbH)
1* 125 pmol RED-tris-NTA Dye 2nd Generation

A9. Labeling Procedure

1. Suspend 125 pmol RED-tris-NTA Dye 2nd Generation in 25 μL of dilution buffer to obtain a 5 μM dye solution.
2. Prepare a 10 μM His₆-peptide stock solution by dissolving the lyophilized peptide into ddH₂O. Verify the concentration spectroscopically using an extinction coefficient of 45,100 $\text{M}^{-1}\text{cm}^{-1}$ (ϵ_{205}).
3. Mix 4 μL of 10 μM His₆-peptide and 2 μL of the RED-tris-NTA Dye 2nd Generation (5 μM) with 194 μL of dilution buffer to obtain 200 μL of a 200 nM His₆-peptide, 50 nM dye solution.
4. Incubate for 30 minutes at room temperature in the dark.

A10. Labeling Efficiency

N/A

B1. Ligand/Non-Fluorescent Binding Partner

Zn^{2+}

B2. Molecule Class/Organism

Divalent cation

B3. Sequence/Formula

ZnCl_2

B4. Purification Strategy/Source

Carl Roth GmbH

[T887.1](#)

B5. Stock Concentration/Stock Buffer

Powdered

B6. Molecular Weight/Extinction Coefficient

136.28 Da

B7. Serial Dilution Preparation

1. Dissolve 0.34 g of ZnCl_2 in 10 mL of ddH₂O to obtain a 250 mM ZnCl_2 stock solution. Add 4 drops of HCl (1.2 M) to acidify the solution and completely dissolve all ZnCl_2 .
2. Mix 4 μL of the 250 mM ZnCl_2 stock with 496 μL of ddH₂O to obtain a 500 μL of a 2 mM ZnCl_2 solution.
3. Prepare a PCR-rack with 16 PCR tubes. Mix 2 μL of the 2 mM ZnCl_2 solution with 18 μL of ddH₂O in tube **1**. Then, add 10 μL of ddH₂O to tubes **2** to **16**.
4. 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.
5. Add 10 μL of labeled His₆-peptide (200 nM) to each tube from **16** to **1** and mix by pipetting.
6. Incubate for 30 minutes at room temperature in the dark before loading capillaries.

B8. EDTA Control

1. Prepare a PCR-rack with 16 fresh PCR tubes. Mix 2 μL of the 2 mM ZnCl_2 solution with 18 μL of ddH₂O in tube **1**. Then, add 10 μL of ddH₂O to tubes **2** to **16**.
2. 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.
3. Mix 10 μL of a 1 mM buffered EDTA solution (pH ~7 – 8) with 184 μL of dilution buffer. Then, add 4 μL of 10 μM His₆-peptide and 2 μL of the RED-tris-NTA Dye 2nd Generation (5 μM) to obtain 200 μL of a 200 nM His₆-peptide, 50 nM dye, 50 μM EDTA solution.
4. Add 10 μL of this solution to each tube from **16** to **1** and mix by pipetting.
5. Incubate for 30 minutes at room temperature in the dark before loading capillaries.

D1. MST System/Capillaries

Monolith NT.115 Red (NanoTemper Technologies GmbH)

Premium Capillaries Monolith NT.115 (MO-K025, NanoTemper Technologies GmbH)

D2. MST Software

MO.Control v1.6 (NanoTemper Technologies GmbH)

nanotempertech.com/monolith-mo-control-software

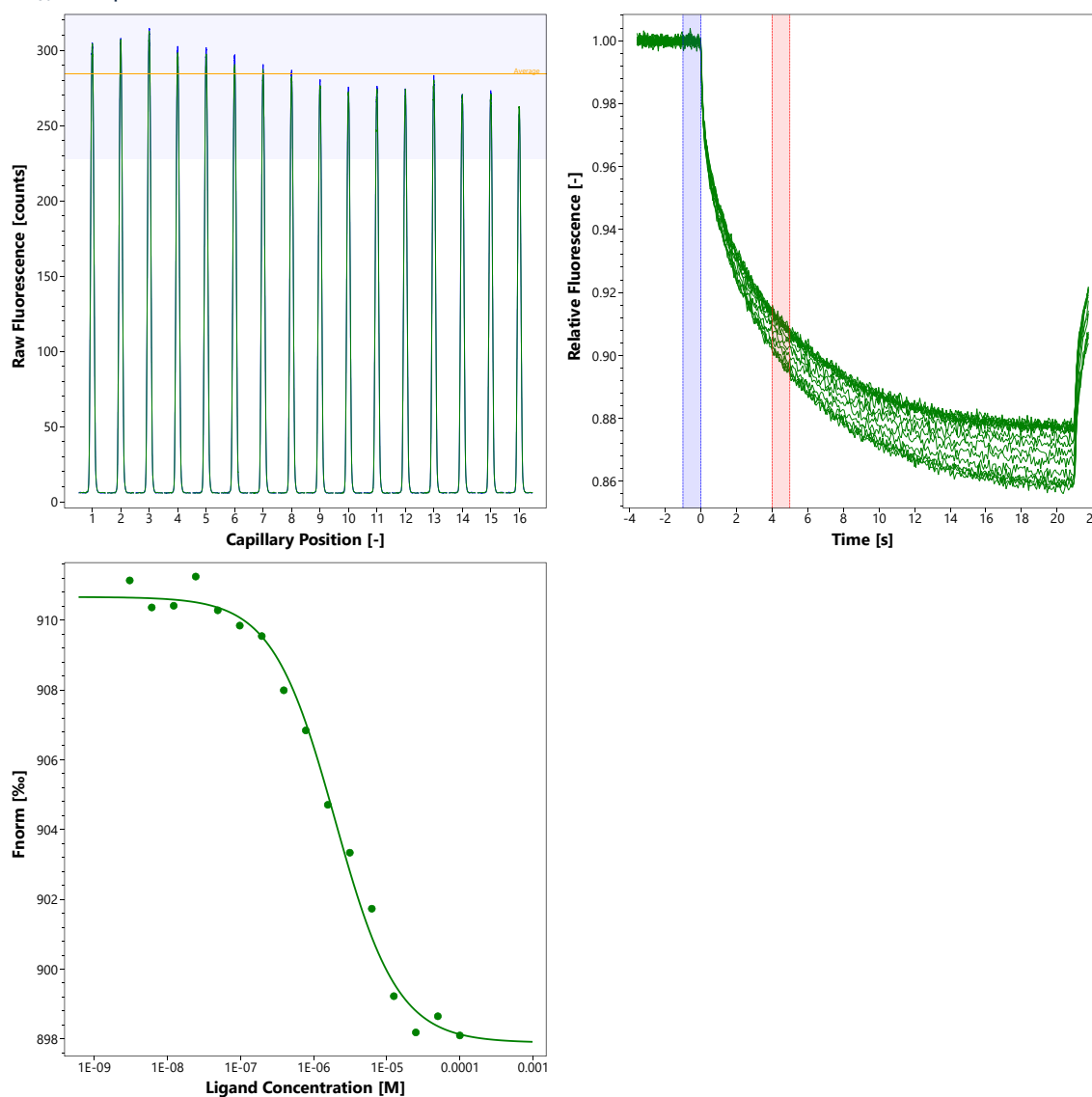
D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

10 mM HEPES, pH 7.4, 75 mM NaCl, 0.005% Pluronic® F-127

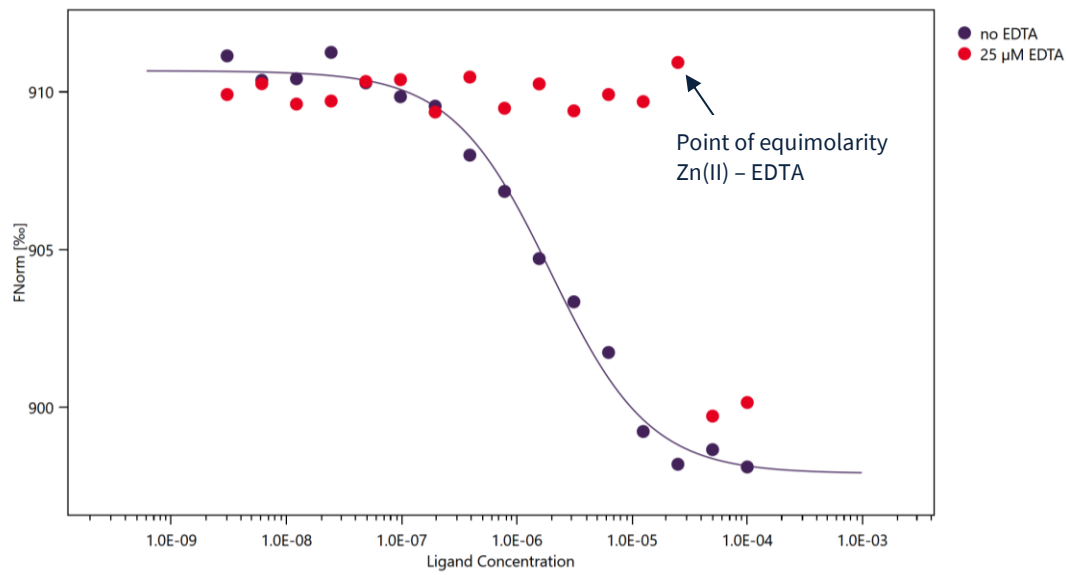
100 nM His₆-peptide | 100 μ M – 3.05 nM ZnCl₂ | 22°C | medium MST power | 40% excitation power

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

EC₅₀ = 1.9 μ M



EDTA control (25 μ M final concentration in the assay):



D5. Reference Results/Supporting Results

Zn(II)-induced dimerization of His-tags

$K_A = 7.7 \cdot 10^{10} \text{ M}^{-2}$ Fluorescence Spectroscopy

Evers et al., *Protein Engineering, Design & Selection* 21 (8), 529–536, 2008

E. Contributors

Andreas Langer²

² NanoTemper Technologies GmbH, München, Germany | nanotempertech.com