

Monolith Protocol MO-P-036

Calmodulin – Ca²⁺

Calmodulin (CaM) is the primary receptor for calcium ions in human cells, acting as a mediator of calcium functions and regulating many cellular processes. Depending on its conformational state, CaM binds or regulates a multitude of protein targets. On binding to calcium, CaM undergoes a large conformational change.

protein – ion interaction | calcium

A1. Target/Fluorescent Molecule

Calmodulin bovine (CaM)

uniprot.org/uniprot/P62157

A2. Molecule Class/Organism

Calcium-binding messenger protein

Bos taurus (Bovine)

A3. Sequence/Formula

MGSSHHHHHH SSSLVPRGSH MADQLTEEQI AEFKEAFSLF DKDGDGTITT KELGTVMRSL GQNPTEAELQ DMINEVDADG
 NGTIDFPEFL TMMARKMKDT DSEEEIREAF RVFDKDGNGY ISAAELRHVM TNLGEKLTDE EVDEMIREAD IDGGQVNYE
 EFVQMMTAK

A4. Purification Strategy/Source

Recombinant, expressed in E. coli, His₆-tagged

Sigma Aldrich GmbH

[C4874](#)

A5. Stock Concentration/Stock Buffer

1 mg/mL | 52.6 μM

A6. Molecular Weight/Extinction Coefficient

19.0 kDa

2,980 M⁻¹cm⁻¹ (ε₂₈₀)

A7. Dilution Buffer

20 mM HEPES, pH 7.4, 150 mM NaCl, 0.005% TWEEN® 20

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. Prepare a 100 μM EDTA solution in dilution buffer.
2. Mix 12.1 μL of dilution buffer with 2.9 μL of 52.6 μM CaM to obtain 15 μL of a 10 μM CaM solution.
3. Mix 4 μL of dilution buffer with 4 μL of 10 μM CaM and 2 μL of 100 μM EDTA to obtain 10 μL of a 4 μM CaM solution in dilution buffer containing 20 μM EDTA (5-fold excess¹).
4. Incubate for 5 minutes to assure that all Ca^{2+} -ions are removed from CaM by EDTA.
5. Add 90 μL of dilution buffer to the solution of step 3 to obtain 100 μL of a 400 nM CaM solution in dilution buffer containing 2 μM EDTA.
6. Suspend 125 pmol RED-tris-NTA Dye 2nd Generation in 25 μL of dilution buffer to obtain a 5 μM dye solution.
7. Mix 123 μL of dilution buffer with 2 μL dye (5 μM) to obtain 125 μL of a 80 nM dye solution.
8. Add 100 μL of the dye solution (80 nM) to the solution of step 5 to obtain 200 μL of a 200 nM CaM, 40 nM dye solution in dilution buffer containing 1 μM EDTA².
9. Incubate for 20 minutes at room temperature in the dark.

A10. Labeling Efficiency

N/A

B1. Ligand/Non-Fluorescent Binding Partner

Calcium ion (Ca^{2+})

B2. Molecule Class/Organism

Divalent metal ion

B3. Sequence/Formula

Ca^{2+}

B4. Purification Strategy/Source

Calcium chloride dihydrate ($\text{CaCl}_2 \cdot 2\text{H}_2\text{O}$)

Carl Roth GmbH

[T885.2](#)

¹ A 5-fold excess of EDTA is used, as CaM can bind up to 4 Ca^{2+} -ions.

² High concentrations of EDTA are not recommended when using the His-Tag Labeling Kit, as EDTA can chelate Ni^{2+} -ions from tris-NTA. At a concentration of only 1 μM EDTA, however, labeling of proteins remains stable over several hours. Furthermore, the bound His-tag helps to 'protect' the Ni^{2+} -ions on tris-NTA from removal by EDTA.

B5. Stock Concentration/Stock Buffer

Powdered

B6. Molecular Weight/Extinction Coefficient

40 Da

B7. Serial Dilution Preparation

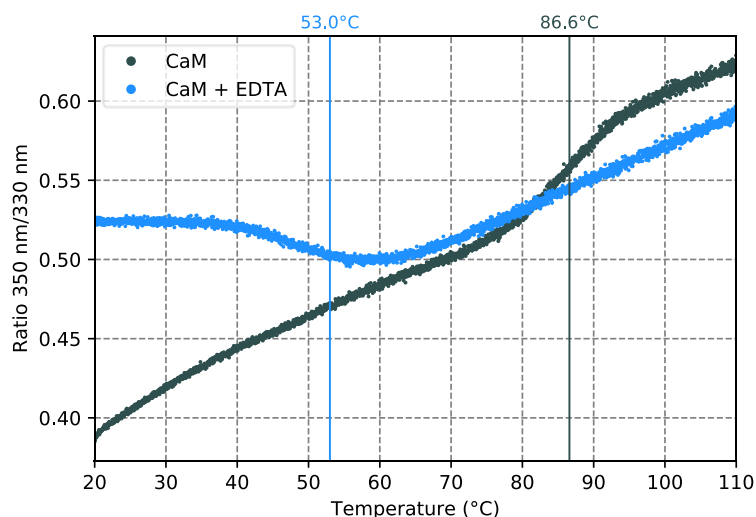
1. Dissolve 73.5 mg of calcium chloride dihydrate (MW: 147,02 g/mol) in 10 mL of dilution buffer to obtain a 50 mM CaCl₂ stock solution.
2. Mix 2 μL of the 50 mM CaCl₂ stock with 198 μL of dilution buffer to obtain 200 μL of a 500 μM CaCl₂ solution.
3. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of the 1 mM CaCl₂ solution into tube **1**. Then, transfer 10 μL of dilution buffer into 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 CaM (~200 nM) to each tube from **16** to **1** and mix by pipetting.
6. Incubate for 5 minutes in the dark before loading capillaries.

C. Applied Quality Checks

Validation of structural integrity and functionality of CaM via a thermal shift assay (1°C/min) with EDTA:

nanotempertech.com/prometheus

CaM	5 μL of 10 μM CaM + 5 μL of dilution buffer	T _m = 86.6°C
CaM + EDTA	5 μL of 10 μM CaM + 5 μL of 100 μM EDTA	T _m = 53.0°C ³



³ As CaM is already complexed with Ca²⁺ in storage buffer, it needs to be removed prior to the MST experiment (c.f. A9, step 4). Since CaM does **not** contain any Tryptophans, unfolding of the Ca²⁺-free form shows a 'blue-shift', while the conformational change upon unfolding of the Ca²⁺-bound form results in a 'red-shift'.

D1. MST System/Capillaries

Monolith NT.115^{PICO} Red (NanoTemper Technologies GmbH)

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

D2. MST Software

MO.Control v1.6 (NanoTemper Technologies GmbH)

<https://nanotempertech.com/monolith-mo-control-software/>

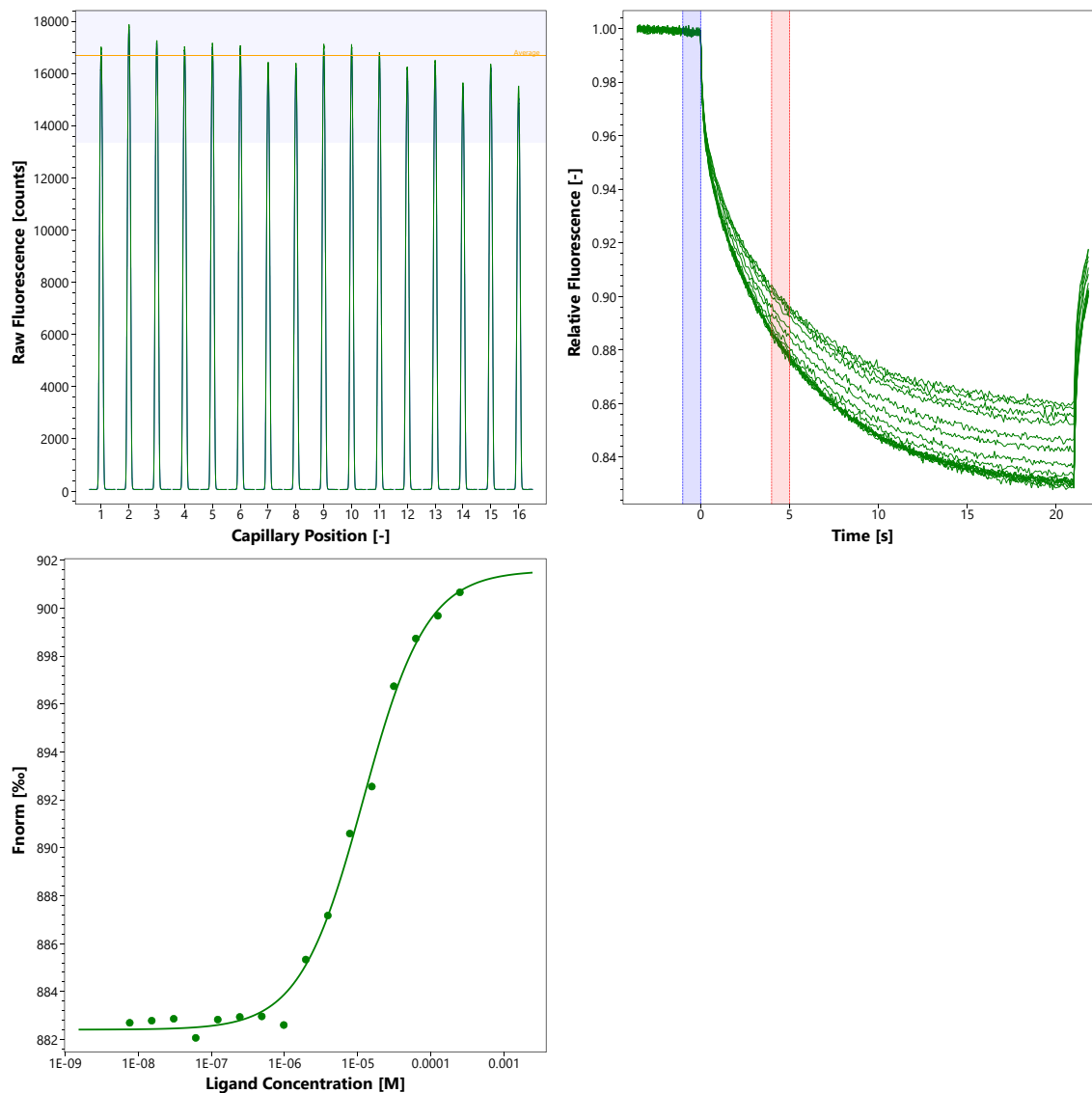
D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

20 mM HEPES, pH 7.4, 150 mM NaCl, 0.5 μ M EDTA, 0.005% TWEEN[®] 20

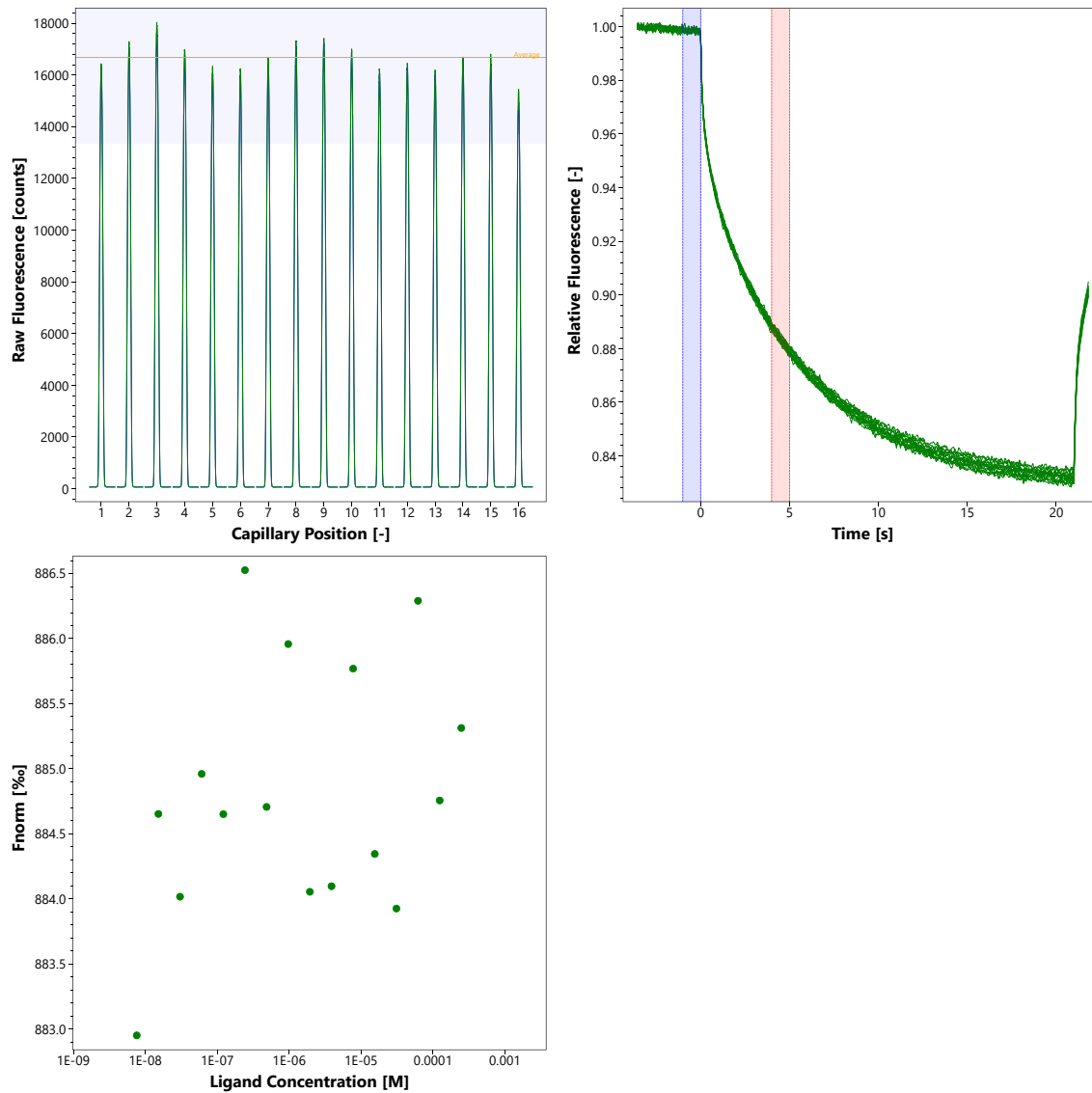
100 nM CaM | 250 μ M – 7.6 nM CaCl₂ | 25°C | medium MST power | 5% excitation power

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

$K_d = 11.8 \mu$ M



Negative control: MgCl₂ instead of CaCl₂



D5. Reference Results/Supporting Results

$K_d = 2 - 50 \mu\text{M}$ Tyrosine fluorescence changes
 Tsuruta and Sano, *Biophysical Chemistry*, 35, 75-84 (1990)

$T_{1/2} = 55^\circ\text{C}$ (w/ EDTA), Circular Dichroism
 $T_{1/2} > 90^\circ\text{C}$ (w/ Ca²⁺) Brzeska et al., *FEBS Letters* 153 (1), 169-173 (1983)

E. Contributors

Andreas Langer⁴

⁴ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com