

Monolith Protocol MO-P-037

DNA Aptamer – AMP (displacement assay)

The DNA aptamer for adenosine is a highly conserved sequence that is a widely used model aptamer for biosensor development. It also binds ADP and ATP, and with slightly weaker affinity AMP. The affinity between unlabeled aptamer and AMP can be determined by a displacement assay with a short, fluorescently labeled ssDNA (cDNA) complementary to parts of the aptamer sequence.

DNA – small molecule interaction | aptamer | displacement assay

A1. Target/Fluorescent Molecule

AMP aptamer

A2. Molecule Class/Organism

DNA aptamer

A3. Sequence/Formula

5' ACC TGG GGG AGT ATT GCG GAG GAA GGT 3'

A4. Purification Strategy/Source

metabion international AG

A5. Stock Concentration/Stock Buffer

17.1 nmol

A6. Molecular Weight/Extinction Coefficient

8485 Da

273,300 M⁻¹cm⁻¹ (ε₂₆₀)

A7. Dilution Buffer

20 mM Tris-HCl, pH 7.8, 300 mM NaCl, 5 mM MgCl₂, 0.05% TWEEN® 20

A8. Labeling Strategy

Competitive binding assay with Cy5 labeled complementary ssDNA (cDNA, metabion international AG)

5' Cy5 ACC TTC CTC C 3'

A9. Labeling Procedure

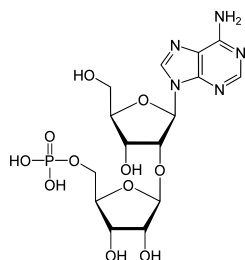
1. Dissolve 26.1 nmol cDNA in 261 μL ddH₂O to obtain a 100 μM cDNA solution.
2. Dissolve 17.1 nmol AMP aptamer in 171 μL ddH₂O to obtain a 100 μM AMP aptamer solution.
3. Add 98 μL of dilution buffer to 2 μL of 100 μM cDNA to obtain 100 μL of a 2 μM cDNA solution.
4. Add 96 μL of dilution buffer to 4 μL of 100 μM AMP aptamer to obtain 100 μL of a 4 μM AMP aptamer solution.
5. Mix 196 μL of dilution buffer with 2 μL of 2 μM cDNA and 2 μL of 4 μM AMP aptamer to obtain 200 μL of a 20 nM cDNA, 40 nM AMP aptamer solution.¹

A10. Labeling Efficiency

HPLC-purified, 100% labeled DNA

B1. Ligand/Non-Fluorescent Binding Partner

Adenosine monophosphate (AMP)



B2. Molecule Class/Organism

Nucleotide monophosphate

B3. Sequence/Formula

C₁₀H₁₄N₅O₇P

B4. Purification Strategy/Source

Sigma-Aldrich GmbH

01930

B5. Stock Concentration/Stock Buffer

17.4 mg/mL | 50 mM

20 mM Tris-HCl, pH 7.8, 300 mM NaCl, 5 mM MgCl₂, 0.05% TWEEN® 20

¹ As the K_d between cDNA and AMP aptamer is ~8 nM (c.f. section D4), a final concentration of 20 nM AMP aptamer results in a sufficient complex formation between cDNA and AMP aptamer.

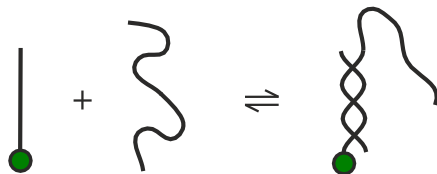
B6. Molecular Weight/Extinction Coefficient

347.22 Da

B7. Serial Dilution Preparation

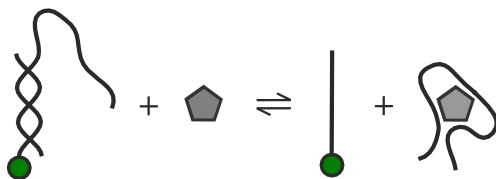
cDNA – AMP aptamer

1. Add 198 μL of dilution buffer to 2 μL of 2 μM cDNA to obtain 200 μL of a 20 nM solution.
2. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of 4 μM AMP aptamer 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. Add 10 μL of cDNA (20 nM) to each tube from **16** to **1** and mix by pipetting.
5. Incubate for 20 minutes at room temperature in the dark before loading capillaries.



AMP aptamer – AMP

1. Prepare a new PCR-rack with 16 PCR tubes. Transfer 20 μL of the 50 mM AMP solution into tube **1**. Then, transfer 10 μL of dilution buffer into 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. Add 10 μL of the 20 nM cDNA, 40 nM AMP aptamer solution to each tube from **16** to **1** and mix by pipetting.
4. Incubate for 20 minutes at room temperature in the dark before loading capillaries.



D1. MST System/Capillaries

Monolith NT.115^{PICO} 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)

cDNA – AMP aptamer

20 mM Tris, pH 7.8, 300 mM NaCl, 5 mM MgCl₂, 0.05% TWEEN® 20
 10 nM cDNA | 2 μM AMP – 61 pM aptamer | 25°C | medium MST power | 3% excitation power

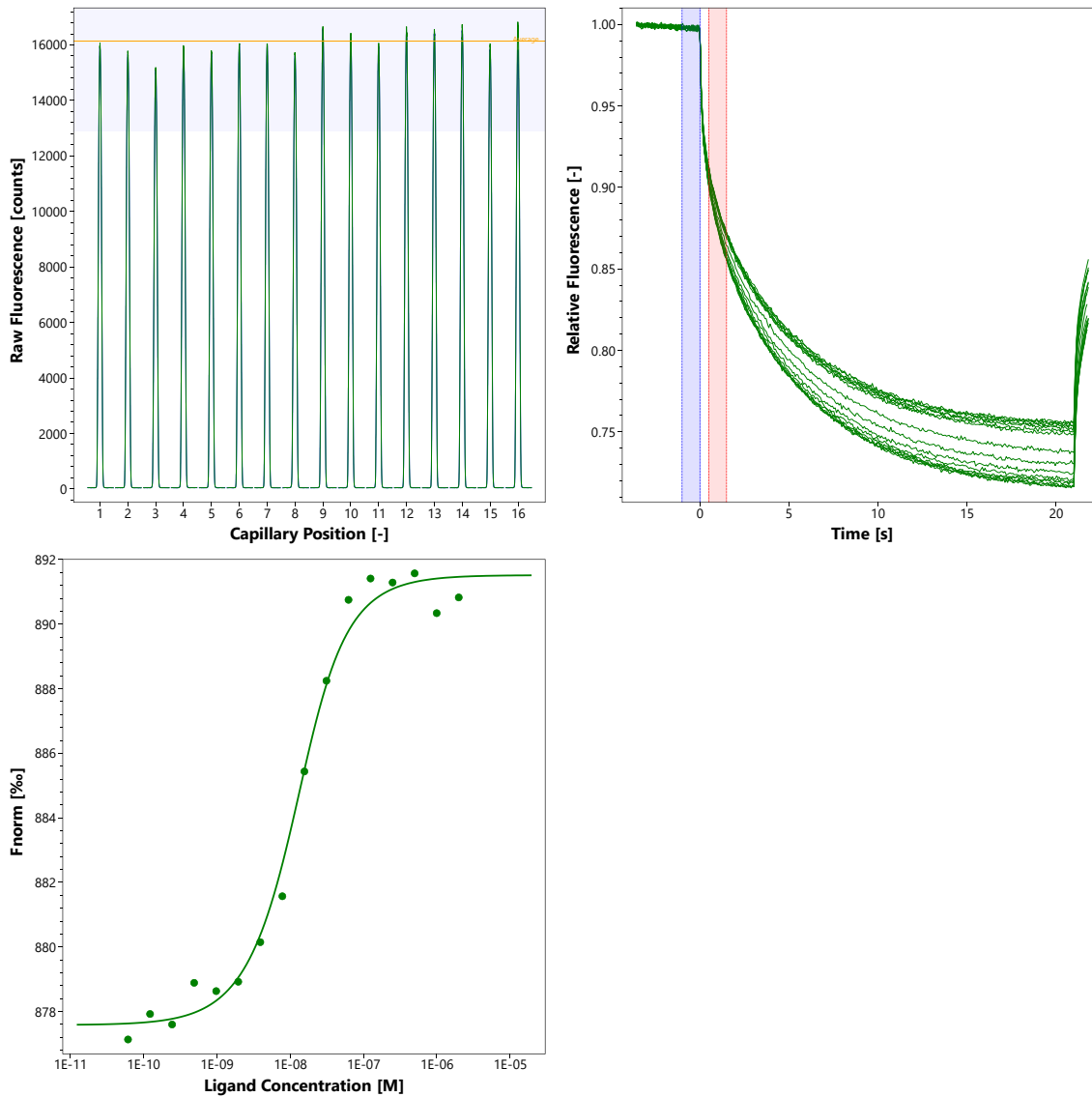
AMP aptamer – AMP

20 mM Tris, pH 7.8, 300 mM NaCl, 5 mM MgCl₂, 0.05% TWEEN® 20
 10 nM cDNA, 20 nM AMP aptamer | 25 mM – 763 nM AMP | 25°C | medium MST power | 3% excitation power

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

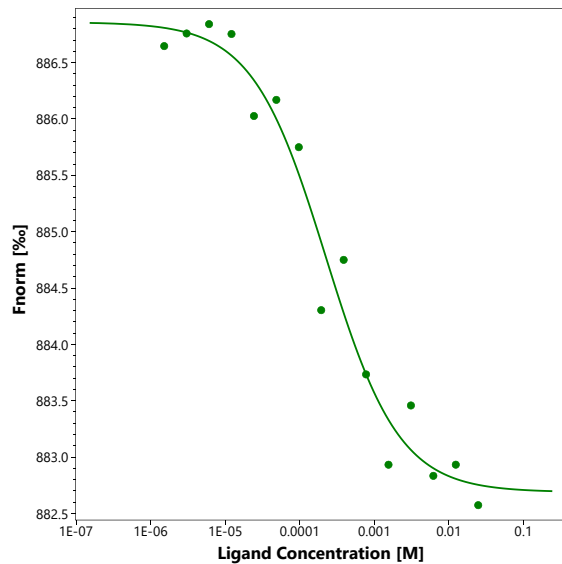
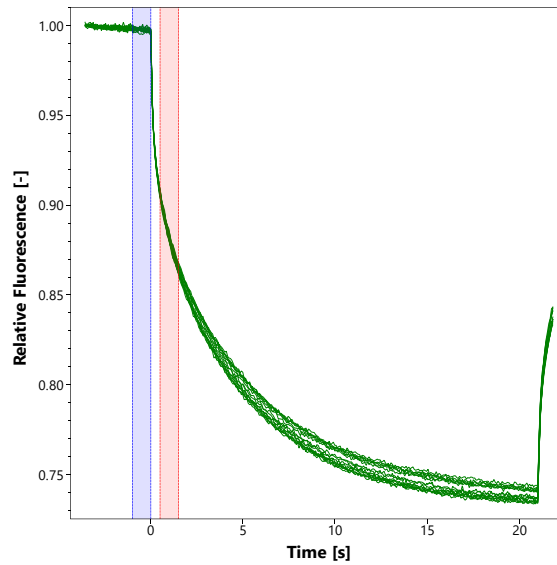
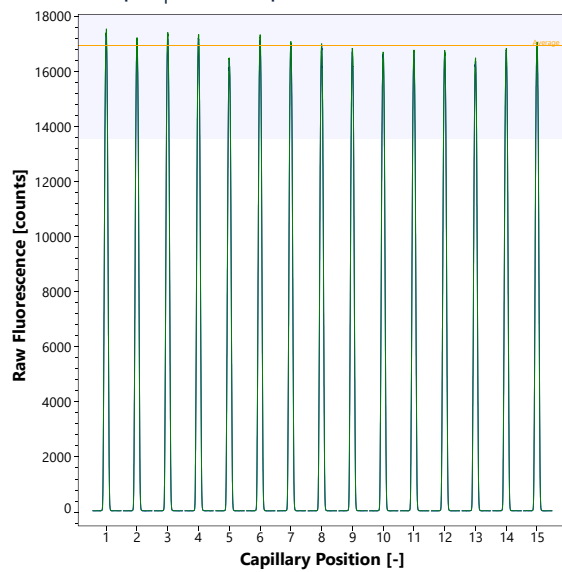
cDNA – AMP aptamer

$K_d = 7.61$ nM



AMP optamer – AMP

$EC_{50} = 226 \mu\text{M}$ | $K_i = 60.6 \mu\text{M}^2$



D5. Reference Results/Supporting Results

$K_d = 58 \pm 2 \mu\text{M}$ Frontal chromatography analysis
[Deng et al., Anal Chem 73 \(2001\) 5415-5421](#)

E. Contributors

Andreas Langer³

² For calculation of K_i , see the NanoTemper 'FAQ Competitive Binding Assay'.

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