

Monolith Protocol MO-P-021

# DNA Aptamer – AMP

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.

DNA – small molecule interaction | aptamer

---

## A1. Target/Fluorescent Molecule

AMP aptamer

## A2. Molecule Class/Organism

DNA aptamer

## A3. Sequence/Formula

5' Cy5 ACC TGG GGG AGT ATT GCG GAG GAA GGT 3'

## A4. Purification Strategy/Source

metabion international AG

## A5. Stock Concentration/Stock Buffer

0.36 µg/ml | 40 nM

20 mM Tris-HCl, pH 7.8, 300 mM NaCl, 5 mM MgCl<sub>2</sub>, 0.05% TWEEN® 20

## A6. Molecular Weight/Extinction Coefficient

9019 Da

273,300 M<sup>-1</sup>cm<sup>-1</sup> (ε<sub>260</sub>)

## A7. Dilution Buffer

20 mM Tris-HCl, pH 7.8, 300 mM NaCl, 5 mM MgCl<sub>2</sub>, 0.05% TWEEN® 20

## A8. Labeling Strategy

5' Cy5 labeled

## A9. Labeling Procedure

N/A

---

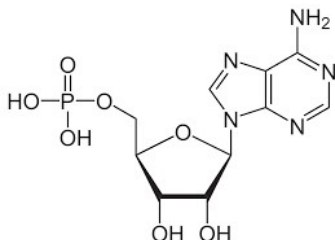
## A10. Labeling Efficiency

HPLC-purified, 100% labeled DNA

---

### B1. Ligand/Non-Fluorescent Binding Partner

Adenosine monophosphate (AMP)



### B2. Molecule Class/Organism

Nucleotide monophosphate

### B3. Sequence/Formula

$C_{10}H_{14}N_5O_7P$

### 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 MgCl<sub>2</sub>, 0.05% TWEEN® 20

### B6. Molecular Weight/Extinction Coefficient

347.22 Da

### B7. Serial Dilution Preparation

1. Prepare a PCR-rack with 16 PCR tubes. Transfer 20  $\mu$ L of the 50 mM AMP solution into tube **1**. Then, transfer 10  $\mu$ L of dilution buffer into tubes **2** to **16**.
2. Prepare a 1:1 serial dilution by transferring 10  $\mu$ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10  $\mu$ L from tube **16** to get an equal volume of 10  $\mu$ L for all samples.
3. Add 10  $\mu$ L of 40 nM AMP aptamer to each tube from **16** to **1** and mix by pipetting.
4. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

### D1. MST System/Capillaries

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

### D2. MST Software

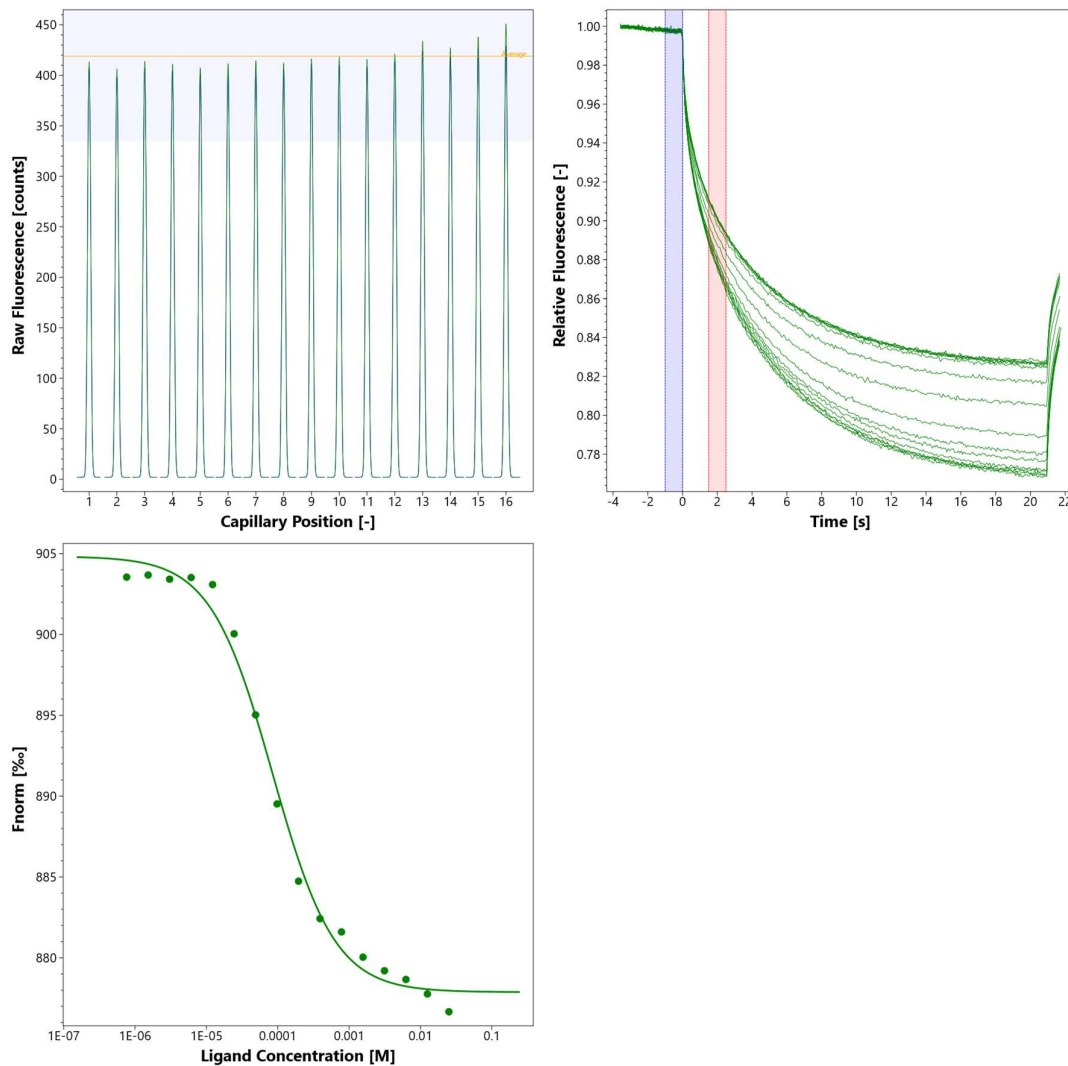
MO.Control v1.6 (NanoTemper Technologies GmbH)  
[nanotempertech.com/monolith-mo-control-software](http://nanotempertech.com/monolith-mo-control-software)

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

20 mM Tris, pH 7.8, 300 mM NaCl, 5 mM MgCl<sub>2</sub>, 0.05% TWEEN® 20  
 20 nM DNA aptamer | 25 mM AMP – 763 nM | 25°C | low MST power | 20% excitation power

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

$K_d = 85 \mu\text{M}$



## 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 Langer<sup>1</sup>

---

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