

Monolith Protocol MO-P-026

mRNA – sRNA₄₁

Methanosarcina mazei is a model organism for methanoarchaea growing under strictly anaerobic conditions. Next to regulatory proteins, also small noncoding RNAs (sRNAs) are involved in nitrogen (N)-regulation, such as sRNA₄₁, which is highly expressed under N-sufficiency. One of its targets is '5'UTR MM2089 mRNA, a polycistronic mRNA encoding for acetyl-CoA-decarbonylase/synthase complexes (ACDS).

RNA – RNA interaction | Nitrogen (N)-regulation

A1. Target/Fluorescent Molecule

'5'UTR MM2089 (mRNA)

A2. Molecule Class/Organism

Polycistronic mRNA Methanosarcina mazei Gö1

A3. Sequence/Formula

5' Cy3 AGU AGC UGA AGC UUU CAA GAA UCG UCG GUA AAG UGA CUU UAA GUA UAU UUA GGA GGU AAA GCU AAA AUG AGC AAA UUA AC 3'

A4. Purification Strategy/Source

Eurofins Genomics GmbH

A5. Stock Concentration/Stock Buffer

1 μM 2x Structure Buffer (Ambion, Thermo Fisher Scientific)

A6. Molecular Weight/Extinction Coefficient

26.1 kDa

A7. Labeling Strategy

5' Cy3 labeled

A8. Assay Buffer

1x Structure Buffer (Ambion, Thermo Fisher Scientific), 0.02% TWEEN® 20



A9. Labeling Procedure

N/A

A10. Labeling Efficiency

HPLC-purified, 100% labeled RNA

B1. Ligand/Non-Fluorescent Binding Partner

 $sRNA_{41}$

B2. Molecule Class/Organism

Noncoding RNA Methanosarcina mazei Gö1

B3. Sequence/Formula

5' GTC TAG CGA ACA GAC GTA AAA ATG GGA AAC GCC TCC TAA ATA TAT TGT ACA AAC TTG CAA TGC TGG AAT TCC CCA CCT GTG GTG CAA GGC GGT GTG GAG TTT CAG GTG CAA GGG AGT CAC 3'

B4. Purification Strategy/Source

In vitro transcription using TranscriptAid T7 High Yield Transcription kit (Thermo Fisher Scientific, Waltham, MA, USA)

B5. Stock Concentration/Stock Buffer

300 μM DEPC-H₂O

B6. Molecular Weight/Extinction Coefficient

39.3 kDa



B7. Serial Dilution Preparation

- 1. Prepare a PCR-rack with 16 PCR tubes. Transfer 10 μ L of 2x Structure Buffer and 5 μ L of DEPC-H₂O into tube **1**. Then add 5 μ L of 300 μ M sRNA₄₁ and mix by pipetting to obtain 20 μ L of a 75 μ M sRNA₄₁ solution in 1x Structure Buffer.
- 2. Transfer 10 μ L of 1x Structure 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 6 μ L of 1 μ M mRNA with 6 μ L of DEPC-H₂O to obtain 12 μ L of 500 nM mRNA in 1x Structure Buffer.
- 5. Add 188 μL 1x Structure Buffer to obtain 200 μL of 30 nM mRNA.
- 6. Incubate all tubes for 1 min at 95°C followed by 5 min at 4°C.
- 7. Add 10 μL of 30 nM mRNA to each tube from 16 to 1 and mix by pipetting.
- 8. Incubate for 15 minutes at 37°C in the dark before loading capillaries.

D1. MST System/Capillaries

Monolith NT.115 Green (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)

1x Structure buffer, 0.02% TWEEN® 20 15 nM mRNA | 37.5 μ M - 1.14 nM sRNA_{41} | 25°C | high MST power | 100% excitation power





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

D5. Reference Results/Supporting Results

 $K_{d} = 631 \pm 133 \text{ nM}$

Microscale Thermophoresis Buddeweg et al., Molecular Microbiology 107(5), 595–609 (2018)

E. Contributors

Anne Buddeweg¹, Teresia Hallström²

¹ Institute for General Microbiology, University of Kiel, Kiel, Germany

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