

## Monolith Protocol MO-P-048

# Mitochondrial Fission 1 Protein – 12 mer Peptide

Mitochondrial fission 1 protein (FIS1) is a component of the mitochondrial complex ARCosome, which promotes mitochondrial fission. FIS1 protein is involved in cell cycle and apoptosis and plays a role in several associated diseases, like diabetic cardiomyopathy, neurodegenerative diseases and cancer.

protein – peptide interaction | mitochondrial fission protein | FIS1

---

### A1. Target/Fluorescent Molecule

Mitochondrial fission 1 protein (FIS1)

[uniprot.org/uniprot/Q9Y306](https://uniprot.org/uniprot/Q9Y306)

### A2. Molecule Class/Organism

Cytosolic domain of mitochondrial outer membrane surface protein

*Homo sapiens (Human)*

### A3. Sequence/Formula

MEAVLNELVS VEDLLKFEKK FQSEKAAGSV SKSTQFEYAW CLVRSKYND D IRKGIVLLEE LLPKGSKEEQ RDYVFYLA VG  
 NYRLKEYEKA LKYVRGLLQT EPQNNQAKEL ERLIDKAMKK DGLVG

### A4. Purification Strategy/Source

Nickel chromatography followed by size-exclusion chromatography, > 95 %

Medical College of Wisconsin (John M. Egnér)

### A5. Stock Concentration/Stock Buffer

0.145 mg/mL | 10 µM

100 mM HEPES, pH 7.4, 200 mM NaCl, 1 mM DTT, 0.02 % sodium azide

### A6. Molecular Weight/Extinction Coefficient

14.5 kDa

15,930 M<sup>-1</sup>cm<sup>-1</sup> (ε<sub>280</sub>)

### A7. Dilution Buffer

100 mM HEPES, pH 7.4, 200 mM NaCl, 1 mM DTT, 0.02 % sodium azide, 0.05 % TWEEN® 20

### A8. Labeling Strategy

Monolith Protein Labeling Kit RED – NHS 2nd Generation (MO-L011, NanoTemper Technologies GmbH)

1\* Labeling Buffer NHS | 1\* Dye RED-NHS 2nd Generation (10 µg) | 1\* B-Column

## A9. Labeling Procedure

1. Prepare 100  $\mu\text{L}$  of a 10  $\mu\text{M}$  FIS1 solution.
2. Use the A-Column to perform a buffer exchange into Labeling Buffer NHS **supplemented** with 1 mM GSH.
  - a. Invert A-Column to suspend slurry and twist off bottom (twist slightly in both directions).
  - b. Loosen the cap of the column and place it in a 1.5 mL microcentrifuge collection tube.
  - c. Centrifuge at **1500  $\times$  g** for **1 min** to remove excess liquid.
  - d. Add 300  $\mu\text{L}$  of Labeling Buffer NHS with GSH and centrifuge at **1500  $\times$  g** for **1 min** (3x).
  - e. Place 100  $\mu\text{L}$  of the 10  $\mu\text{M}$  FIS1 solution in the center of the resin.
  - f. Place the sample in a **new** microcentrifuge collection tube and centrifuge at **1500  $\times$  g** for **2 min**.

The collected flow-through should yield around 100  $\mu\text{L}$  of  $\sim 5 \mu\text{M}$  FIS1 ( $\sim 50\%$  recovery).
3. Add 25  $\mu\text{L}$  of DMSO to Dye RED-NHS 2nd Generation (10  $\mu\text{g}$ ) to obtain a  $\sim 600 \mu\text{M}$  solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
4. Mix 10  $\mu\text{L}$  of the 600  $\mu\text{M}$  dye solution with 90  $\mu\text{L}$  of Labeling Buffer NHS with GSH to obtain 100  $\mu\text{L}$  of a 60  $\mu\text{M}$  dye solution (8x protein concentration).
5. Mix FIS1 and dye in a 1:1 volume ratio (200  $\mu\text{L}$  final volume).
6. Incubate for 30 minutes at room temperature in the dark.
7. In the meantime, remove the top cap of the B-Column and pour off the storage solution. Remove the bottom cap and place with adapter in a 15 mL tube.
8. Fill the column with dilution buffer and allow it to enter the packed resin bed completely by gravity flow. Discard the flow through collected. Repeat this step 3 more times.
9. Add 200  $\mu\text{L}$  of the labeling reaction from step 5 to the center of the column and let sample enter the bed completely.
10. Add 400  $\mu\text{L}$  of dilution buffer after the sample has entered and discard the flow through.
11. Place column in a new collection tube, add 500  $\mu\text{L}$  of dilution buffer and collect the eluate.
12. Keep the labeled FIS1 ( $\sim 1 \mu\text{M}$ ) on ice in the dark.

## A10. Labeling Efficiency

Measurement of protein concentration and degree of labeling (DOL) using a NanoDrop™:

[nanotempertech.com/dol-calculator](https://nanotempertech.com/dol-calculator)

Absorbance $A_{280}$	0.021	Protein concentration	1.1 $\mu\text{M}$
Absorbance $A_{650}$	0.096	Degree-of-labeling (DOL)	0.45

## B1. Ligand/Non-Fluorescent Binding Partner

12 mer peptide (pep2)

## B2. Molecule Class/Organism

Peptide

## B3. Sequence/Formula

N/A

#### B4. Purification Strategy/Source

Genscript peptide synthesis

#### B5. Stock Concentration/Stock Buffer

5.66 mM

100 mM HEPES, pH 7.4, 200 mM NaCl, 1 mM DTT, 0.02 % sodium azide

#### B6. Molecular Weight/Extinction Coefficient

N/A

#### B7. Serial Dilution Preparation

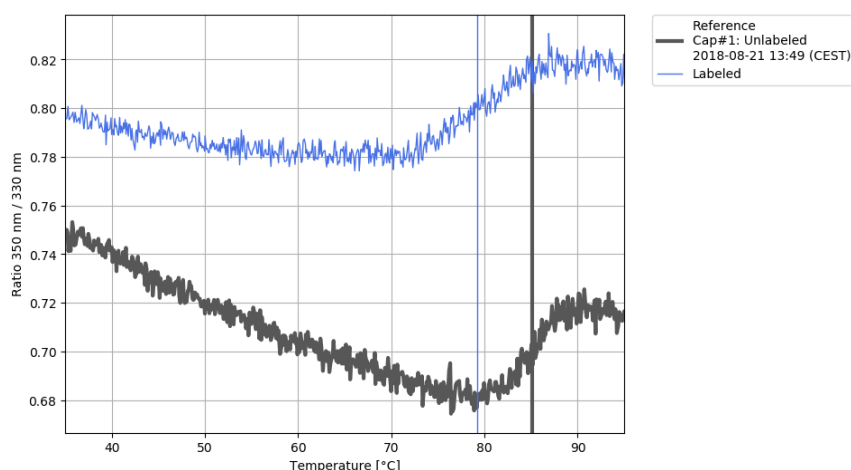
1. Prepare a PCR-rack with 16 PCR tubes. Transfer 7  $\mu\text{L}$  of the 5.66 mM pep2 stock solution into tube **1** and mix with 13  $\mu\text{L}$  of dilution buffer to obtain 20  $\mu\text{L}$  of a 2 mM solution. Then, transfer 10  $\mu\text{L}$  of dilution buffer into tubes **2** to **16**.
2. Prepare a 1:1 serial dilution by transferring 10  $\mu\text{L}$  from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10  $\mu\text{L}$  from tube **16** to get an equal volume of 10  $\mu\text{L}$  for all samples.
3. Mix 2  $\mu\text{L}$  of labeled FIS1 with 198  $\mu\text{L}$  of dilution buffer to obtain 200  $\mu\text{L}$  of  $\sim 10$  nM FIS1.
4. Add 10  $\mu\text{L}$  of labeled FIS1 ( $\sim 10$  nM) 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.

### C. Applied Quality Checks

Validation of structural integrity of labeled FIS1 using Tycho NT.6:

[nanotempertech.com/tycho](https://nanotempertech.com/tycho)

Unlabeled	10 $\mu\text{L}$ of FIS1 in dilution buffer ( $\sim 5$ $\mu\text{M}$ )	$T_i = 85.1^\circ\text{C}$
Labeled	10 $\mu\text{L}$ of B-Column eluate ( $\sim 1.1$ $\mu\text{M}$ )	$T_i = 79.2^\circ\text{C}$



## D1. MST System/Capillaries

Monolith NT.115<sup>Pico</sup> 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/>

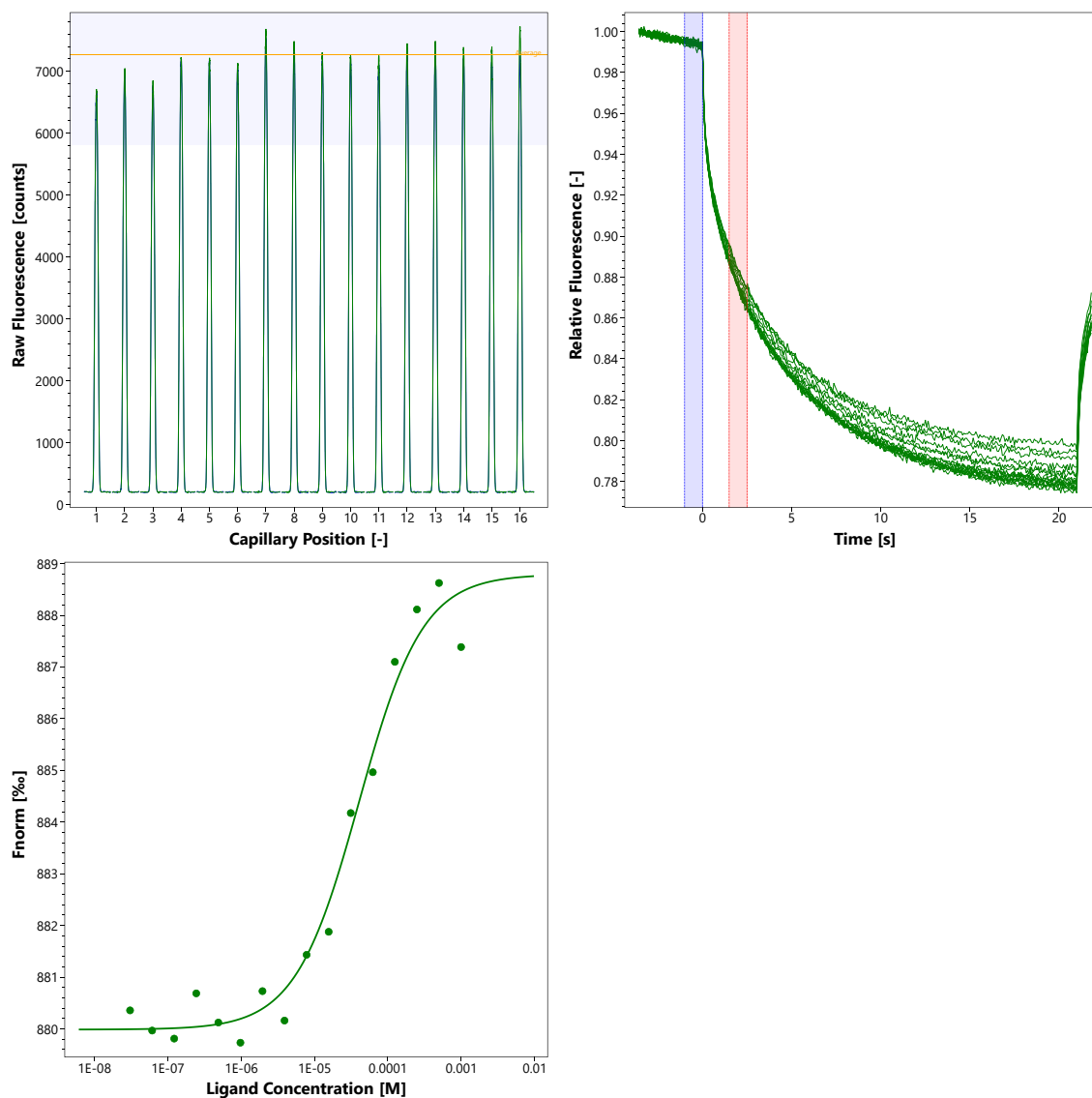
## D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

100 mM HEPES, pH 7.4, 200 mM NaCl, 1 mM DTT, 0.02 % sodium azide, 0.05 % TWEEN® 20

5 nM FIS1 | 1 mM – 30.5 nM pep2 | 25°C | medium MST power | 20% excitation power

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

$K_d = 40.6 \mu\text{M}$



## D5. Reference Results/Supporting Results

$K_d = 200 \mu\text{M}$       2D-protein detected NMR titration experiments  
[Inhouse-measurements at Medical College of Wisconsin](#)

---

## E. Contributors

John M. Egner<sup>1</sup>, Blake Hill<sup>1</sup>, Tanja Bartoschik<sup>2</sup>

---

<sup>1</sup> Medical College of Wisconsin, Milwaukee, United States

<sup>2</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](https://nanotempertech.com)