

Monolith Protocol MO-P-014

HSP90 – ADP

Heat shock protein 90 (Hsp90) is an ATP-dependent, ubiquitously expressed and highly abundant chaperone that aids in folding, stabilizing and degrading various other proteins, including kinases, steroid hormone receptors and transcription factors. By repeated ATP hydrolysis-coupled rounds of substrate binding and release, it helps both newly synthesized and misfolded proteins to find their native fold. Since a lot of Hsp90-dependent proteins are strongly upregulated in tumor cells, mammalian Hsp90 remains a promising target for anti-cancer drugs.

protein – small molecule interaction | chaperone

A1. Target/Fluorescent Molecule

Heat shock protein HSP90-alpha (Hsp90)

uniprot.org/uniprot/P07900

A2. Molecule Class/Organism

Heat shock protein 90 (Hsp90 family)

Homo sapiens (Human)

A3. Sequence/Formula

GSMPEETQTQ DQPMEEEEVE TFAFQAEIAQ LMSLIINTFY SNKEIFLREL ISNSSDALDK IRYESLTDPS KLD SGKELHI
NLIPNKQDRT LTIVDTGIGM TKADLINNLG TIAKSGTKAF MEALQAGADI SMIGQFGVGF YSAYLVAEKV TVITKHNDDE
QYAWESSAGG SFTVRTDTGE PMGRGTKVIL HLKEDQTEYL EERRIKEIVK KHSQFIGYPI TLFVEKERDK EVSDDEAE EK
EDKEEEKEKE EKESEDKPEI EDVGSDEEEE KKGDKKKKK KIKEYIDQE ELNKTPIWT RNPDDITNEE YGEFYKSLTN
DWEHLAVKH FSVEGQLEFR ALLFVPRRAP FDLFENRKKK NNIKLYVRRV FIMDNCEELI PEYLNFI RGV VDS EDLPLNI
SREMLQQSKI LKVIRKNLVK KCLELFTELA EDKENYKKFY EQFSKNIKLG IHEDSQNRKK LSELLRYYS ASGDEM VSLK
DYCTRMKENQ KHIYYITGET KDQVANS AFV ERLRKHGLEV IYMI EPIDEY CVQQLKEFEG KTLVSVTKEG LELPEDEEEK
KKQEEKTKTF ENLCKIMKDI LEKKVEKVVV SNRLVTSPCC IVTSTYGWTA NMERIMKAQA LRDNSTMGYM AAKKHLEINP
DHSIIETLRQ KAEADKNDKS VKDLVILLYE TALLSSGFSL EDPQTHANRI YRMIKLG LGI DEDDPTADDT SAAVTEEMP
LEGDDDTSRM EEVD

A4. Purification Strategy/Source

Expressed in *E. coli* RIL, His₆-tagged

Crelux GmbH

A5. Stock Concentration/Stock Buffer

4.82 mg/mL | 56.8 µM

20 mM Tris-HCl, pH 8.0, 250 mM NaCl, 1 mM TCEP

A6. Molecular Weight/Extinction Coefficient

84.8 kDa

59,250 M⁻¹cm⁻¹ (ε₂₈₀)

A7. Dilution Buffer

50 mM Tris-HCl, pH 7.8, 150 mM NaCl, 10 mM MgCl₂, 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* A-Column | 1* Dye RED-NHS 2nd Generation (10 µg) | 1* B-Column

A9. Labeling Procedure

1. Add 75 µL of Labeling Buffer NHS to 25 µL of 56.8 µM Hsp90 to obtain 100 µL of a 14.2 µM solution.
2. Use the A-Column to perform a buffer exchange into Labeling Buffer NHS.
 - 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 × g** for **1 min** to remove excess liquid.
 - d. Add 300 µL of Labeling Buffer NHS and centrifuge at **1500 × g** for **1 min** (3x).
 - e. Place 100 µL of the 14.2 µM Hsp90 solution in the center of the resin.
 - f. Place the sample in a **new** microcentrifuge collection tube and centrifuge at **1500 × g** for **2 min**.

The collected flow-through should yield around 100 µL of ~10 µM Hsp90 (70 – 80% recovery).
3. Add 25 µL of DMSO to Dye RED-NHS 2nd Generation (10 µg) to obtain a ~600 µM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
4. Mix 5 µL of the 600 µM dye solution with 95 µL of Labeling Buffer NHS to obtain 100 µL of a 30 µM dye solution (3x protein concentration).
5. Mix Hsp90 and dye in a 1:1 volume ratio (200 µL final volume, 2.5% final DMSO concentration).
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 µL of the labeling reaction from step 5 to the center of the column and let sample enter the bed completely.
10. Add 400 µL of dilution buffer after the sample has entered and discard the flow through.
11. Place column in a new collection tube, add 500 µL of dilution buffer and collect the eluate.
12. Keep the labeled Hsp90 (~2 µM) on ice in the dark.

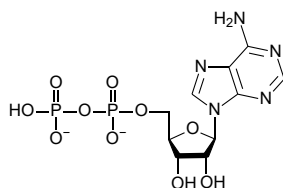
A10. Labeling Efficiency

Measurement of protein concentration and degree of labeling (DOL) using a NanoDrop™:
nanotempertech.com/dol-calculator

Absorbance A ₂₈₀	0.140	Protein concentration	2.03 µM
Absorbance A ₆₅₀	0.490	Degree-of-labeling (DOL)	1.24

B1. Ligand/Non-Fluorescent Binding Partner

Adenosine diphosphate (ADP)



B2. Molecular Class/Organism

Nucleotide diphosphate

B3. Sequence/Formula

$C_{10}H_{15}N_5O_{10}P_2$

B4. Purification Strategy/Source

Sigma-Aldrich GmbH

[A2754](#)

B5. Stock Concentration/Stock Buffer

42.7 mg/mL | 100 mM
ddH₂O

B6. Molecular Weight/Extinction Coefficient

427.20 Da

B7. Serial Dilution Preparation

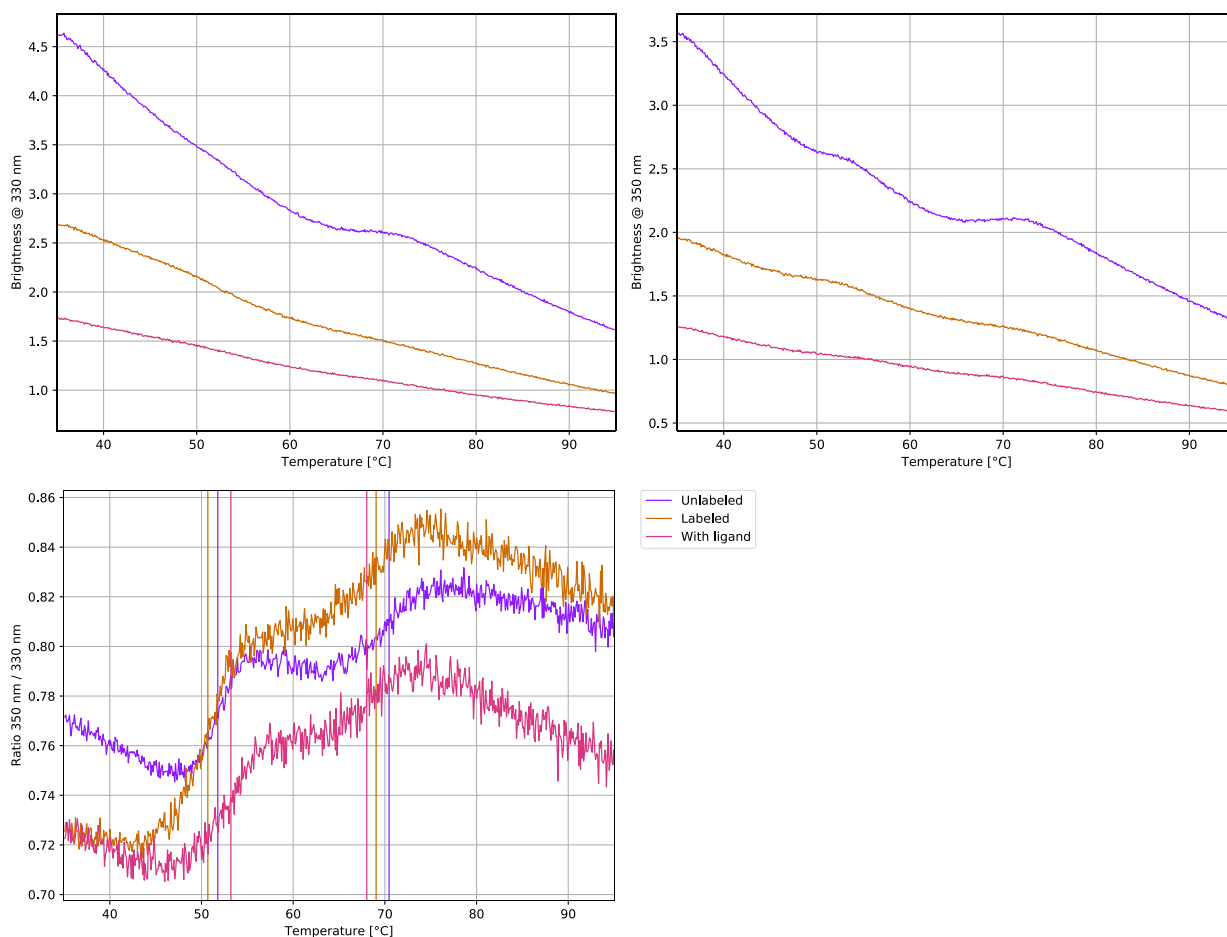
1. Add 27 μ L of dilution buffer to 3 μ L of the 100 mM ADP stock to obtain 30 μ L of a 10 mM solution.
2. Add 180 μ L of dilution buffer to 20 μ L of ddH₂O to obtain 200 μ L of diluted dilution buffer.
3. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μ L of the 10 mM ADP solution into tube **1**. Then, transfer 10 μ L of diluted dilution buffer into tubes **2** to **16**.
4. Prepare a 1:1 serial dilution by transferring 10 μ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10 μ L from tube **16** to get an equal volume of 10 μ L for all samples.
5. Mix 4 μ L of labeled Hsp90 (~2 μ M) with 196 μ L of dilution buffer to obtain 200 μ L of ~40 nM labeled Hsp90.
6. Add 10 μ L of labeled Hsp90 (~40 nM) to each tube from **16** to **1** and mix by pipetting.
7. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

C. Applied Quality Checks

Validation of structural integrity and functionality of labeled Hsp90 using Tycho NT.6:

nanotempertech.com/tycho

Unlabeled	10 μ L of 2 μ M Hsp90	$T_{i,1} = 51.8^{\circ}\text{C}$ $T_{i,2} = 70.5^{\circ}\text{C}$
Labeled	5 μ L of B-Column eluate ($\sim 2 \mu\text{M}$) + 5 μ L of dilution buffer	$T_{i,1} = 50.7^{\circ}\text{C}$ $T_{i,2} = 69.0^{\circ}\text{C}$
With ligand	5 μ L of B-Column eluate ($\sim 2 \mu\text{M}$) + 5 μ L of 4 mM ADP	$T_{i,1} = 53.2^{\circ}\text{C}$ $T_{i,2} = 68.0^{\circ}\text{C}$



D1. MST System/Capillaries

Monolith NT.115 Red (NanoTemper Technologies GmbH)

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

D2. MST Software

MO.Control v1.6 (NanoTemper Technologies GmbH)

nanotempertech.com/monolith-mo-control-software

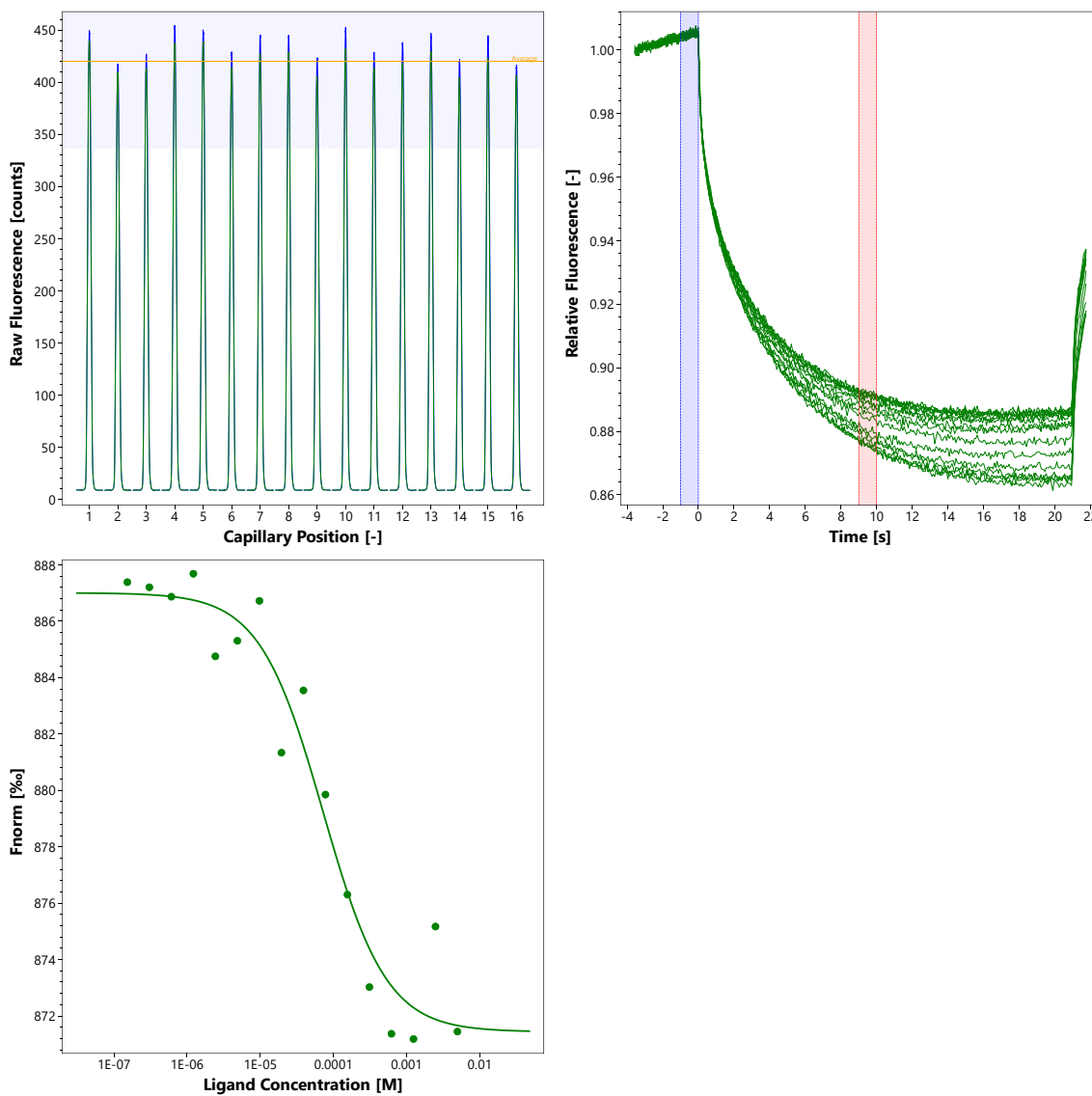
D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

50 mM Tris-HCl, pH 7.8, 150 mM NaCl, 10 mM MgCl₂, 0.05% TWEEN® 20

20 nM Hsp90 | 5 mM ADP – 153 nM | 25°C | medium MST power | 80% excitation power

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

$K_d = 72.9 \mu\text{M}$



D5. Reference Results/Supporting Results

$K_d = 29 \mu\text{M}$

Isothermal Titration Calorimetry (ITC)

[Prodromou et al., Cell 90\(1\), 65-75 \(1997\)](#)

E. Contributors

Timm Hassemer¹

¹ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com