

# p38- $\alpha$ – SB 203580 (purification-free)

Mitogen-activated protein kinase 14 (also called p38- $\alpha$ ) is an enzyme that has been implicated in the regulation of many proinflammatory pathways. It is a drug target for many diseases such as rheumatoid arthritis, endotoxic shock and osteoporosis. SB 203580 is a specific and potent inhibitor of kinases that binds p38- $\alpha$  with nM affinity. Using His-tag labeling, the interaction can be measured purification-free in cell lysate.

protein – small molecule interaction | kinase | inhibitor | His-tag | cell lysate | purification-free

## A1. Target/Fluorescent Molecule

Mitogen-activated protein kinase 14 (p38- $\alpha$ )

[uniprot.org/uniprot/Q16539](https://uniprot.org/uniprot/Q16539)

## A2. Molecule Class/Organism

p38 mitogen-activated protein kinase (MAP kinase)

*Homo sapiens* (Human)

## A3. Sequence/Formula

MSQERPTFYR QELNKTIWEV PERYQNLSPV GSGAYGSVCA AFDTKTGLRV AVKKLSRPFQ SIIHAKRTYR ELRLKHKMH  
ENVIGLLDVF TPARSLEEFN DVYLVTHLMG ADLNNIVKCQ KLTDDHVQFL IQILRGLKY IHSADIIHRD LKPSNLAVNE  
DCELKILDFG LARHTDDEMT GYVATRWYRA PEIMLNWMHY NQTVDIWSVG CIMAELLTGR TLFPGTDHID QLKLIILRLVG  
TPGAELLKKI SSESARNYIQ SLTQMPKMNF ANVFIGANPL AVDLLEKMLV LDSDKRITAA QALAHAYFAQ YHDPDDEPVA  
DPYDQSFESR DLLIDEWKS L TYDEVISFVP PPLDQEEMES

Constructs: p38- $\alpha$ -mNeonGreen-His<sub>6</sub> | mNeonGreen-His<sub>6</sub> (negative control)

## A4. Purification Strategy/Source

### Cloning

Using In-Fusion Cloning technology (Takara Bio USA, Inc.), p38- $\alpha$  coding sequence was cloned in a pcDNA3.1 mammalian expression vector behind a CMV promoter and separated from the mNeongreen-His<sub>6</sub>-tag by the linker sequence ESGSGS. A pcDNA3.1 vector coding for only mNeongreen-His<sub>6</sub> was used as a control.

### Expression

The two plasmids expressing mNeongreen-His<sub>6</sub> with and without the p38- $\alpha$  sequence were transfected into  $3 \times 10^6$  HeLa cells using separate T-75 flasks. Cells were grown for 24 h reaching approximately  $10^7$  cells.

### Lysate Preparation

Cells were pelleted by centrifugation and resuspended in 1 mL PBS-T buffer<sup>1</sup> (137 mM NaCl, 2.5 mM KCl, 10 mM Na<sub>2</sub>HPO<sub>4</sub>, 2 mM KH<sub>2</sub>PO<sub>4</sub>, pH 7.4, 0.05% TWEEN® 20), supplemented with protease inhibitors. At this step, the cells were disrupted using a Dounce homogenizer and centrifuged again at  $14\,000 \times g$  for 30 min at 4°C to remove cell debris. Obtained supernatant was diluted 1:10 in PBS-T and supplemented with protease inhibitors (PMSF)<sup>2</sup>.

<sup>1</sup> RIPA buffer was tested but did not work for this assay.

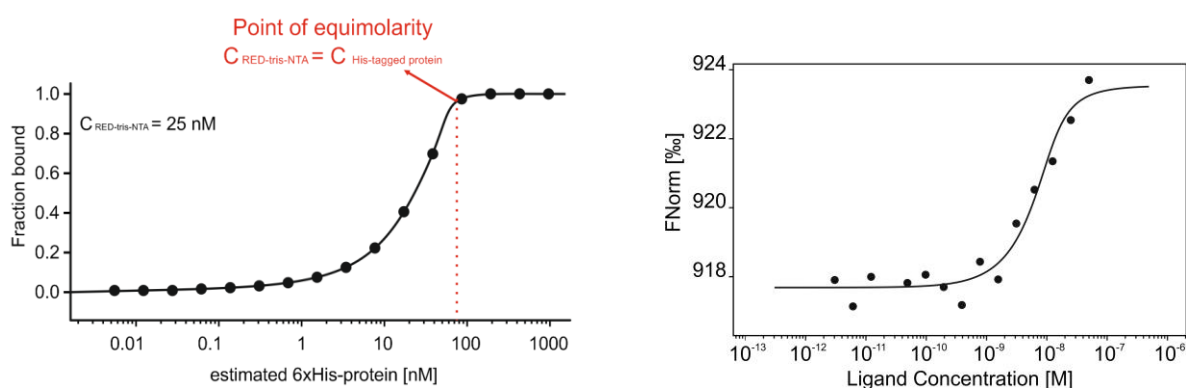
<sup>2</sup> It is recommended to work with freshly prepared cell lysates when performing MST experiments.

## Concentration Determination

Concentration of His<sub>6</sub>-tagged p38- $\alpha$  in cell lysate was determined as described below. A 16-point serial dilution of cell lysate expressing p38- $\alpha$  was prepared using non-transfected HeLa cell lysates in 10  $\mu$ L of final volume. 10  $\mu$ L of 50 nM RED-tris-NTA dye were added to each reaction tube, resulting in a final dye concentration of 25 nM. The reaction mixture was incubated for 30 min at room temperature and loaded into Premium Capillaries Monolith NT.115 (MO-K025, NanoTemper Technologies GmbH).

The MST experiment was carried out using 50% Excitation and medium MST power. MST data were evaluated after 10 s MST-on time. Since the affinity of the RED-tris-NTA dye towards the His<sub>6</sub>-tag is in the low nM range (2.1 nM as measured for the purified protein) and therefore well below the concentration of RED-tris-NTA, the binding curve displays a 'kink' at the point where equimolar concentrations of dye and His-tagged protein are present (simulated dose response, left panel).

As there is a single binding site for the RED-tris-NTA dye, one can assume that dye and protein concentrations are equivalent at the 'kink', so that the concentration of the protein at the start of the dilution can be calculated. In this experiment, saturation of the binding curve was not reached (right panel) suggesting that the concentration of the protein in the lysate was less than 2-fold above the concentration of the dye (less or equal to 50 nM).



## A5. Stock Concentration/Stock Buffer

50 nM (p38- $\alpha$ -mNeonGreen-His<sub>6</sub>)

Phosphate-buffered saline (PBS, pH 7.4), 2 mM PMSF, 0.05% TWEEN® 20

## A6. Molecular Weight/Extinction Coefficient

44.7 kDa

50,100 M<sup>-1</sup>cm<sup>-1</sup> ( $\epsilon_{280}$ )

## A7. Dilution Buffer

Phosphate-buffered saline (PBS, pH 7.4), 0.05% TWEEN® 20

### **A8. Labeling Strategy**

Monolith His-Tag Labeling Kit RED-tris-NTA (MO-L008, NanoTemper Technologies GmbH)  
1\* PBS-T | 1\* 250 pmol RED-tris-NTA Dye

### **A9. Labeling Procedure**

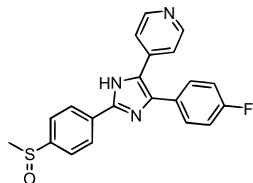
1. Prepare a 100 nM dye solution by mixing 2  $\mu\text{L}$  of dye (5  $\mu\text{M}$ ) and 98  $\mu\text{L}$  PBS-T.
2. Add 100  $\mu\text{L}$  of lysate containing p38 $\alpha$ -mNeonGreen-His<sub>6</sub>.
3. Incubate for 30 minutes at room temperature in the dark.
4. Centrifuge the reaction mixture for 10 min at 15,000  $\times$  g.

### **A10. Labeling Efficiency**

N/A

## B1. Ligand/Non-Fluorescent Binding Partner

SB 203580



## B2. Molecule Class/Organism

p38 MAPK inhibitor

## B3. Sequence/Formula

C<sub>21</sub>H<sub>16</sub>FN<sub>3</sub>OS

## B4. Purification Strategy/Source

Sigma-Aldrich GmbH

[S8307](#)

## B5. Stock Concentration/Stock Buffer

1 mg/mL | 2.65 mM

DMSO

## B6. Molecular Weight/Extinction Coefficient

377.43 Da

## B7. Serial Dilution Preparation

1. Mix 2 µL of the 2.65 mM SB 203580 solution with 98 µL of dilution buffer to obtain 100 µL of a 53 µM SB 203580 solution in dilution buffer containing 2% DMSO.
2. Mix 4 µL of DMSO with 196 µL of dilution buffer to obtain 200 µL of a 2% DMSO solution.
3. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 µL of the 53 µM SB 203580 solution into tube **1**. Then, transfer 10 µL of the 2% DMSO solution into tubes **2** to **16**.
4. 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.
5. Add 10 µL of labeled p38α-mNeonGreen-His<sub>6</sub> to each tube from **16** to **1** and mix by pipetting.
6. Incubate for 20 minutes in the dark and then centrifuge for 10 min at 14000 × g before loading capillaries.

## 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](https://nanotempertech.com/monolith-mo-control-software)

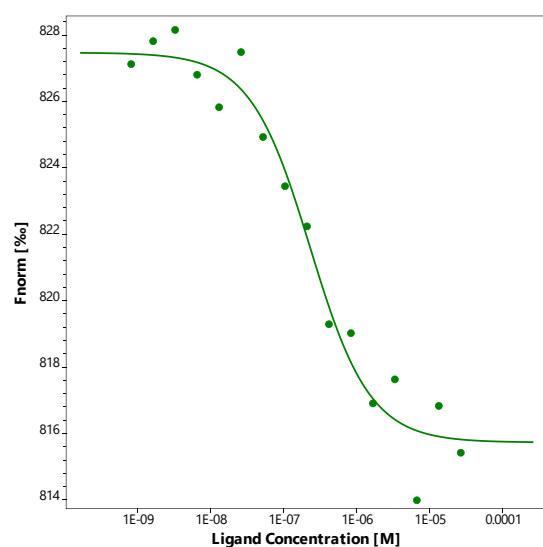
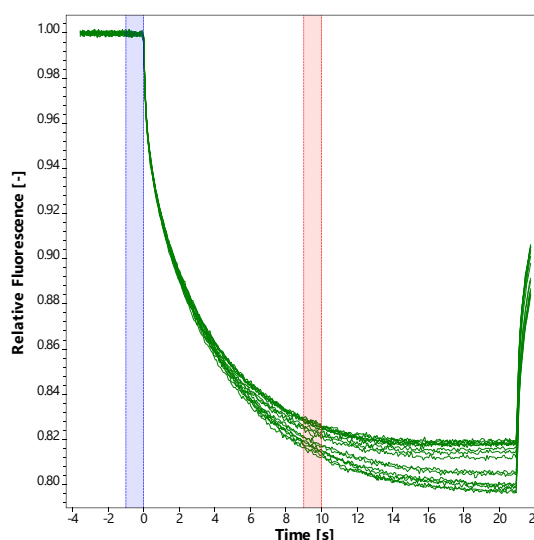
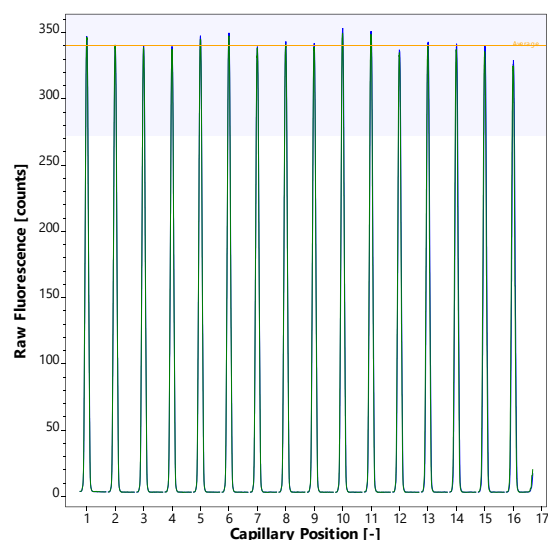
## D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

Phosphate-buffered saline (PBS, pH 7.4), 0.05% TWEEN® 20, 1% DMSO

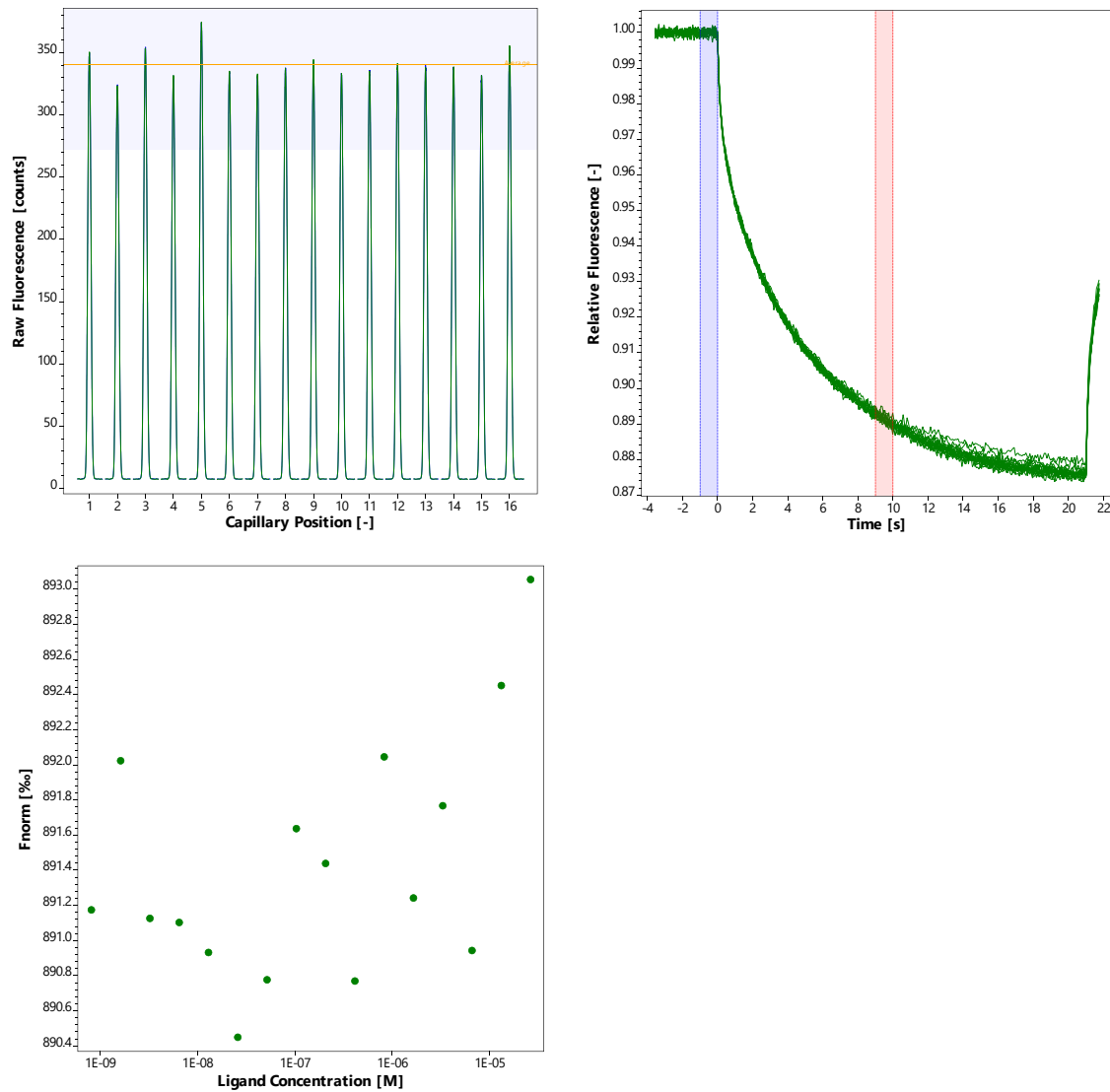
25 nM p38- $\alpha$  | 26.5  $\mu$ M – 809 pM SB 203580 | 25°C | high MST power | 20% excitation power

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

$K_d = 116$  nM (p38- $\alpha$ -mNeonGreen-His<sub>6</sub> – SB 203580)



$K_d = \text{N/A}$  (mNeonGreen-His<sub>6</sub> – SB 203580, negative control)



## D5. Reference Results/Supporting Results

$K_d = 116 \text{ nM}$       Microscale Thermophoresis  
[Bartoschik et al., Scientific Reports 8 \(2018\) 4977](#)

## E. Contributors

Katarzyna Walkiewicz<sup>3</sup>

<sup>3</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](http://nanotempertech.com)