

## Monolith X Protocol MOX-P-104

# STAT3 – SD-36 (PROTAC) – Cereblon

A proteolysis targeting chimera (PROTAC) is a heterobifunctional small molecule composed of two active domains and a linker, capable of removing specific unwanted proteins. Rather than acting as a conventional enzyme inhibitor, a PROTAC works by inducing selective intracellular proteolysis. PROTACs consist of two covalently linked protein-binding molecules: one capable of engaging an E3 ubiquitin ligase, and another one that binds to a target protein meant for degradation. Cereblon forms an E3 ubiquitin ligase complex with damaged DNA binding protein 1 (DDB1), which then can ubiquitinate other proteins. Signal transducer and activator of transcription 3 (STAT3) is a transcription factor. SD-36 is a potent and efficacious STAT3 PROTAC degrader. It is composed of the STAT3 inhibitor SI-109, a linker, and an analog of Cereblon ligand Lenalidomide for E3 ubiquitin ligase.

protein – small molecule | PROTACs | ternary complex | Hook effect

### A1. Target/Fluorescent Molecule

Signal transducer and activator of transcription 3 (STAT3)

[uniprot.org/uniprot/P40763](https://uniprot.org/uniprot/P40763)

### A2. Molecule Class/Organism

Transcription factor

*Homo sapiens (Human)*

### A3. Sequence/Formula

GQANHPTAAV VTEKQQMLEQ HLQDVRKRVQ DLEQKMKVVE NLQDDDFDFNY KTLKSQGDMQ DLNGNNSQSVT RQKMQQLEQM  
 LTALDQMRRS IVSELAGLLS AMEYVQKTLT DEELADWKRR QQIACIGGPP NICLDRLLENW ITSLAESQLQ TRQIQKLEE  
 LQQKVSYSYKGD PIVQHRPMLE ERIVELFRNL MKSAFVVERQ PCMPMHPDRP LVIKTGVQFT TKVRLLVKFP ELNYQLKIKV  
 CIDKDSGDVA ALRGSRFNI LGTNTKVMNM EESNNGSLSA EFKHLTLREQ RCGNGGRANC DASLIVTEEL HLITFETEVY  
 HQGLKIDLET HSLPVVVISN ICQMPNAWAS ILWYNMLTNN PKNVNFFTKP PIGTWDQVAE VLSWQFSSTT KRGLSIEQLT  
 TLAEKLLGPG VNYSGCQITW AKFCKENMAG KGFSFWVWLD NIIDLKVKYI LALWNEGYIM GFISKERERA ILSTKPPGTF  
 LLRFSESSKE GGVTFTWVEK DISGKTQIQS VEPYTKQQLN NMSFAEIMG YKIMDATNIL VSPLVLYLPD IPKEEAFGKY  
 CRPESQEHPE ADPGSAAPYL KTKFICVTPG SGENLYFQGH HHHHH

### A4. Purification Strategy/Source

Recombinant

CreLux GmbH

Construct ID: DDO3, Lot-ID: PC10912

### A5. Stock Concentration/Stock Buffer

12.87 mg/mL | 195 µM

100 mM Tris-HCl pH 8.5, 100 mM NaCl, 2 mM DTT

## A6. Molecular Weight/Extinction Coefficient

66.1 kDa

84,340 M<sup>-1</sup>cm<sup>-1</sup> (ε<sub>280</sub>)

## A7. Dilution Buffer

20 mM HEPES pH 7.4, 150 mM NaCl, 0.5 mM TCEP, 0.01 % Pluronic® F-127

## A8. Labeling Strategy

Monolith Protein Labeling Kit RED – NHS 2nd Generation (MO-L014, NanoTemper Technologies GmbH)

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

## A9. Labeling Procedure

1. Mix 2 µL of 195 µM STAT3 with 38 µL of dilution buffer to obtain 40 µL of ~10 µM STAT3.
2. Use the A-Column to perform a buffer exchange into dilution buffer.
  - 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 2 mL microcentrifuge collection tube.
  - c. Centrifuge at **1500 × g** for **1 min** to remove excess liquid.
  - d. Add 300 µL of dilution buffer and centrifuge at **1500 × g** for **1 min** (3x).
  - e. Place 40 µL of the 10 µM STAT3 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 40 µL of ~5 µM STAT3 (~50% 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 2 µL of the 600 µM dye solution with 38 µL of dilution buffer to obtain 40 µL of a 30 µM dye solution (6x protein concentration).
5. Mix STAT3 and dye in a 1:1 volume ratio (80 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 80 µL of the labeling reaction from step 5 to the center of the column and let sample enter the resin bed completely.
10. Add 500 µL of dilution buffer after the sample has entered and discard the flow through.
11. Place column in a new collection tube, add 400 µL of dilution buffer and collect the eluate.
12. Prepare 8 µL aliquots of the labeled STAT3 (~0.5 µM) and immediately store at -80°C.

## A10. Labeling Efficiency

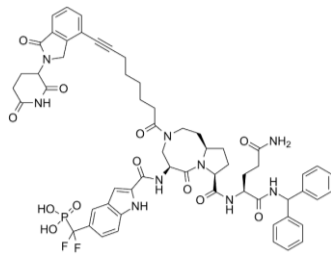
Measurement of protein concentration and degree of labeling (DOL) using a NanoDrop™:

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

Absorbance A <sub>280</sub>	0.059	Protein concentration	0.66 µM
Absorbance A <sub>650</sub>	0.086	Degree-of-labeling (DOL)	0.67

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

SD-36



Cereblon E3 ligase complex (CRBN/DDB1)

CRBN

[uniprot.org/uniprot/Q96SW2](https://uniprot.org/uniprot/Q96SW2)

DDB1

[uniprot.org/uniprot/Q16531](https://uniprot.org/uniprot/Q16531)

### B2. Molecule Class/Organism

PROTAC

E3 protein ligase complex

*Homo sapiens (Human)*

### B3. Sequence/Formula

C<sub>59</sub>H<sub>62</sub>F<sub>2</sub>N<sub>9</sub>O<sub>12</sub>P

```

MGSSHHHHHH SSGENLYFQG SMEAKKPNII NFDTSLPTSH TYLGADMEEF
HGRTLHDDDS CQVIPVLPQV MMILIPGQTL PLQLFHPQEV SMVRNLIQKD
RTFAVLAYSN VQEREAQFGT TAEIYAYREE QDFGIEIVKV KAIGRQRFKV
LELRTQSDGI QQAKVQILPE CVLPSTMSAV QLESLNKCQI FPSKPVSRED
QCSYKWWQKY QKRKFHCANL TSWPRWLYSL YDAETLMDRI KKQLREW DEN
LKDDSLPSNP IDFSYRVAAC LPIDDLVLIQ LLKIGSAIQR LRCELDIMNK
CTSLCCKQCQ ETEITTKNEI FSLSLCGPMA AYVNPHGYVH ETLTVYKACN
LNLIGRPSTE HSWFPGYAWT VAQCKICASH IGWKFATKK DMSPQKFWGL
TRSALLPTIP DTEDEISPK VILCL

MSYNYVVTAQ KPTAVNGCVT GHFTSAEDLN LLIKNTRLE IYVVTAEGLR
PVKEVGMYGK IAVMELFRPK GESKDLLFIL TAKYNACILE YKQSGESIDI
ITRAHGNVQD RIGRPSETGI IGIIDPECRM IGLRLYDGLF KVIPLDRDNK
ELKAFNIRLE ELHVIDVKFL YGCQAPTICF VYQDPQGRHV KTYEVSLEK
EFNKGPKWQKE NVEAEASMI AVPEPFGGAI IIGQESITYH NGDKYLAIA
PIIKQSTIVC HNRVDPNGSR YLLGDMEGRL FMLLLEKEEQ MDGVTTLKDL
RVELLGETSI AECLTYLDNG VVFGSRLGD SQLVKLNVD SNEQGSYVAM
ETFTNLGPIV DMCVVDLERQ GQGQLVTC SG AFKEGSLRII RINGIGIHEHA
SIDLPGIKGL WPLRSDPNRE TDDTLVLSFV GQTRVLMNG EEEVETELMG
FVDDQQTFFC GNAHQQLIQ ITSASVRLVS QEPKALVSEW KEPQAKNISV
ASCNSSQVV AVGRALYYLQ IHPQELRQIS HTEMEHEVAC LDITPLGDSN
GLSPLCAIGL WTDISARILK LPSFELLHKE MLGGEIIPRS ILMTTFESSH
YLLCALGDGA LFYFGLNIET GLLSDRKKVT LGTQPTVLR FRSLSTTNV
ACSDRPTVIY SSNHKLVFSN VNLKEVNYMC PLNSDGYPDS LALANNSTLT
IGTIDEIQKL HIRTVPLYES PRKICYQEV S QCFGVLS SRI EVQDTSGGTT
ALRPSASTQA LSSSVSSSKL FSSSTAPHET SFGEEVEVHN LLIIDQHTFE
VLHAHQFLQN EYALSLV SCK LGKDPNTYFI VGTAMVY PEE AEPKQGRIVV
FQYSDGKLQT VAEKEVKGAV YSMVEFNGKL LASINSTVRL YEWTTKEELR
TECNHYNIM ALYLKTKGDF ILVGDLMRSV LLLAYKPMEG NFEEIARDFN
PNWMSAVEIL DDDNFLGAEN AFNLFVCQKD SAATTDEERQ HLQEVGLFHL
GEFVNVFCHG SLVMQNLGET STPTQGSVLF GTVNGMIGLV TSLSES WYNL
LLDMQNRLNK VIKSVGKIEH SFWRSFHTER KTEPATGFID GD LIESFLDI
SRPKMQEVVA NLQYDDGSGM KREATADLI KVVEELTRIH

```

#### B4. Purification Strategy/Source

SD-36  
 MedChemExpress  
 HY-129602

Crelex GmbH  
 Construct ID: DLS7/CXU4, Lot-ID: PH15311

#### B5. Stock Concentration/Stock Buffer

11.6 mg/mL | 10 mM  
 DMSO

6.55 mg/mL | 37.2  $\mu$ M  
 10 mM HEPES pH 7.0, 240 mM NaCl, 1 mM TCEP

#### B6. Molecular Weight/Extinction Coefficient

1158.15 Da

48.8 kDa (CRBN), 127.0 kDa (DDB1)  
 166,510  $\text{M}^{-1}\text{cm}^{-1}$  ( $\epsilon_{280}$ )

## B7. Serial Dilution Preparation

1. Dissolve 5 mg of SD-36 in 432  $\mu\text{L}$  of DMSO to obtain a 10 mM stock solution.
2. Mix 2  $\mu\text{L}$  of 10 mM SD-36 with 98  $\mu\text{L}$  of DMSO to obtain 100  $\mu\text{L}$  of a 200  $\mu\text{M}$  SD-36 solution.
3. Mix 4  $\mu\text{L}$  of 200  $\mu\text{M}$  SD-36 with 196  $\mu\text{L}$  of dilution buffer to obtain 100  $\mu\text{L}$  of a 4  $\mu\text{M}$  SD-36 solution.
4. Mix 20  $\mu\text{L}$  of DMSO with 980  $\mu\text{L}$  of dilution buffer to obtain 1 mL of dilution buffer with 2% DMSO.
5. Mix 20  $\mu\text{L}$  of labeled STAT3 ( $\sim 0.5 \mu\text{M}$ ) with 480  $\mu\text{L}$  of dilution buffer to obtain 500  $\mu\text{L}$  of 20 nM STAT3.
6. Centrifuge at 15,000 rpm and 4°C for 30 minutes to remove aggregates. Then, carefully transfer the supernatant into a fresh tube and mix well by pipetting.
7. Mix 5.9  $\mu\text{L}$  of 37.2  $\mu\text{M}$  CRBN with 104.1  $\mu\text{L}$  of dilution buffer to obtain 110  $\mu\text{L}$  of 2  $\mu\text{M}$  CRBN.

### STAT3 – SD-36 (binary complex)

8. Mix 100  $\mu\text{L}$  of 20 nM STAT3 with 100  $\mu\text{L}$  of dilution buffer to obtain 200  $\mu\text{L}$  of 10 nM STAT3.
9. Take a fresh 0.5 mL tube and mix 160  $\mu\text{L}$  of 10 nM STAT3 with 160  $\mu\text{L}$  of dilution buffer containing 2% DMSO to obtain 320  $\mu\text{L}$  of a 5 nM STAT3 solution.
10. Prepare a PCR-rack with 16 PCR tubes. Transfer 20  $\mu\text{L}$  of the 5 nM STAT3 solution into tubes **2** to **16**. Then, mix 20  $\mu\text{L}$  of the 10 nM STAT3 solution with 20  $\mu\text{L}$  of the 4  $\mu\text{M}$  SD-36 solution in tube **1**.
11. Prepare a 1:1 serial dilution by transferring 20  $\mu\text{L}$  from tube to tube. Mix carefully by pipetting up and down.
12. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

### STAT3 & high concentration of CRBN – SD-36 (ternary complex, saturation)

13. Mix 100  $\mu\text{L}$  of 20 nM STAT3 with 100  $\mu\text{L}$  of 2  $\mu\text{M}$  CRBN to obtain 200  $\mu\text{L}$  of 10 nM STAT3, 1  $\mu\text{M}$  CRBN.
14. Take a fresh 0.5 mL tube and mix 160  $\mu\text{L}$  of 10 nM STAT3, 1  $\mu\text{M}$  CRBN with 160  $\mu\text{L}$  of dilution buffer containing 2% DMSO to obtain 320  $\mu\text{L}$  of a 5 nM STAT3, 500 nM CRBN solution.
15. Prepare a PCR-rack with 20 PCR tubes. Transfer 20  $\mu\text{L}$  of the 5 nM STAT3, 500 nM CRBN solution into tubes **2** to **16**. Then, mix 20  $\mu\text{L}$  of 4  $\mu\text{M}$  SD-36 with 20  $\mu\text{L}$  of 10 nM STAT3, 1  $\mu\text{M}$  CRBN in tube **1**.
16. Prepare a 1:1 serial dilution by transferring 20  $\mu\text{L}$  from tube to tube. Mix carefully by pipetting up and down.
17. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

### STAT3 & low concentration of CRBN – SD-36 (ternary complex, Hook effect)

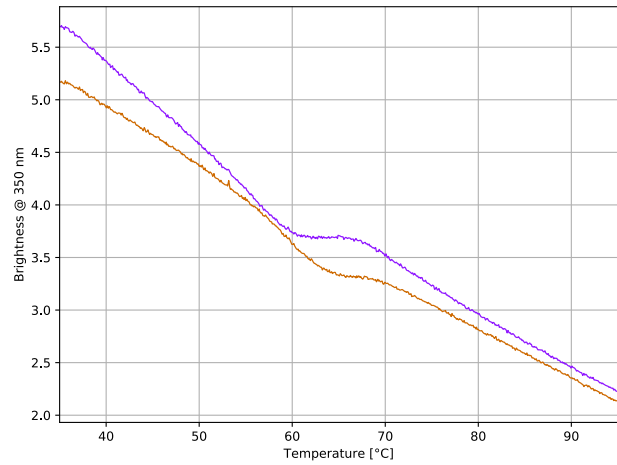
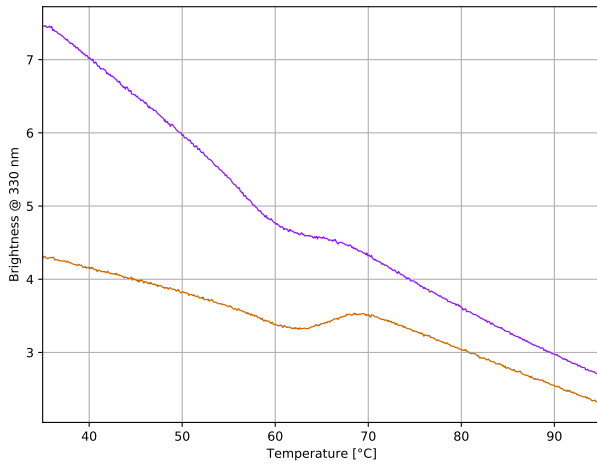
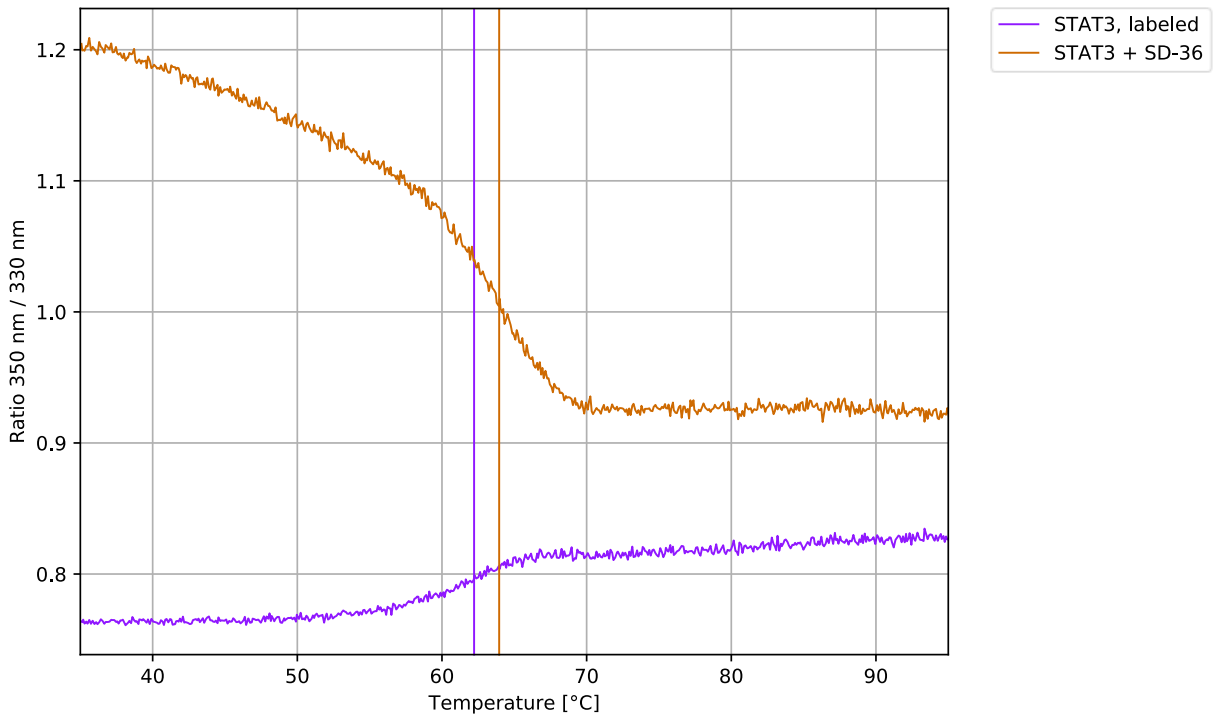
18. Mix 100  $\mu\text{L}$  of 20 nM STAT3 with 8  $\mu\text{L}$  of 2  $\mu\text{M}$  CRBN and 92  $\mu\text{L}$  of dilution buffer to obtain 200  $\mu\text{L}$  of 10 nM STAT3, 80 nM CRBN.
19. Take a fresh 0.5 mL tube and mix 160  $\mu\text{L}$  of 10 nM STAT3, 80 nM CRBN with 160  $\mu\text{L}$  of dilution buffer containing 2% DMSO to obtain 320  $\mu\text{L}$  of a 5 nM STAT3, 40 nM CRBN solution.
20. Prepare a PCR-rack with 20 PCR tubes. Transfer 20  $\mu\text{L}$  of the 5 nM STAT3, 40 nM CRBN solution into tubes **2** to **16**. Then, mix 20  $\mu\text{L}$  of 4  $\mu\text{M}$  SD-36 with 20  $\mu\text{L}$  of 10 nM STAT3, 80 nM CRBN in tube **1**.
21. Prepare a 1:1 serial dilution by transferring 20  $\mu\text{L}$  from tube to tube. Mix carefully by pipetting up and down.
22. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

### C. Tycho

Validation of structural integrity and functionality of labeled STAT3 using Tycho NT.6:

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

STAT3, labeled	8 $\mu$ L of B-Column eluate ( $\sim 0.5 \mu$ M) + 2 $\mu$ L of dilution buffer with 2% DMSO	$T_i = 62.2^\circ\text{C}$
STAT3 + SD-36	8 $\mu$ L of B-Column eluate ( $\sim 0.5 \mu$ M) + 2 $\mu$ L of 4 $\mu$ M SD-36	$T_i = 64.0^\circ\text{C}$



## D1. Monolith System/Capillaries

Monolith X (NanoTemper Technologies GmbH)

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

## D2. Monolith Software

MO.Control v2.4.2 (NanoTemper Technologies GmbH)

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

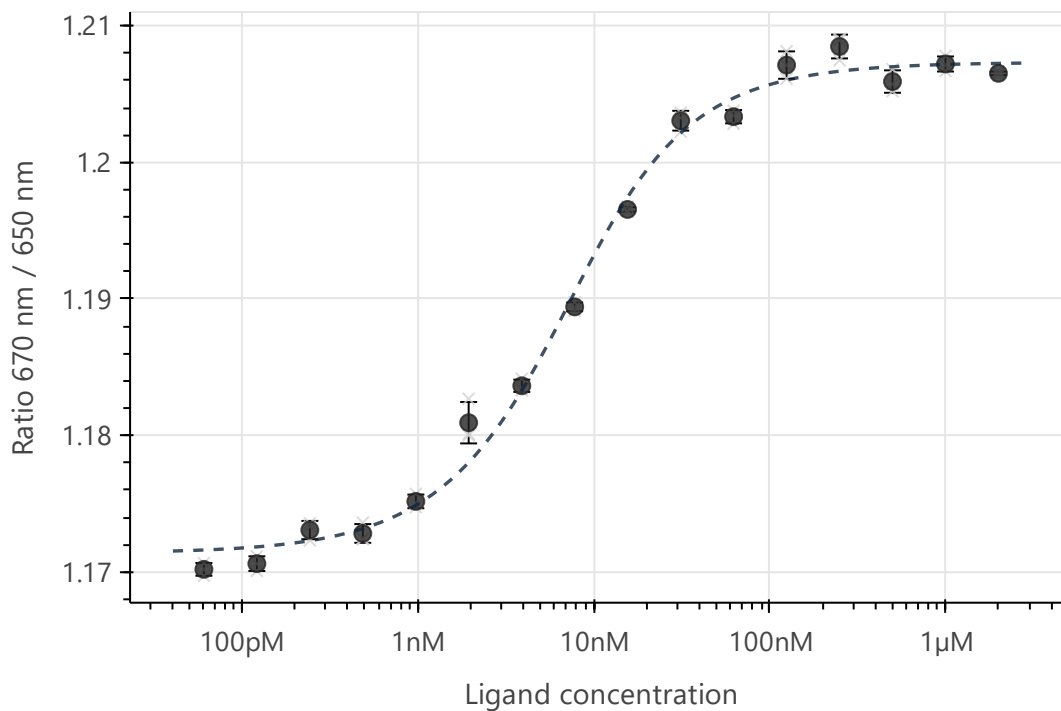
## D3. Monolith Experiment (Assay Buffer/Concentrations/Temperature/Excitation Power)

20 mM HEPES pH 7.4, 150 mM NaCl, 0.5 mM TCEP, 0.01 % Pluronic® F-127, 1% DMSO

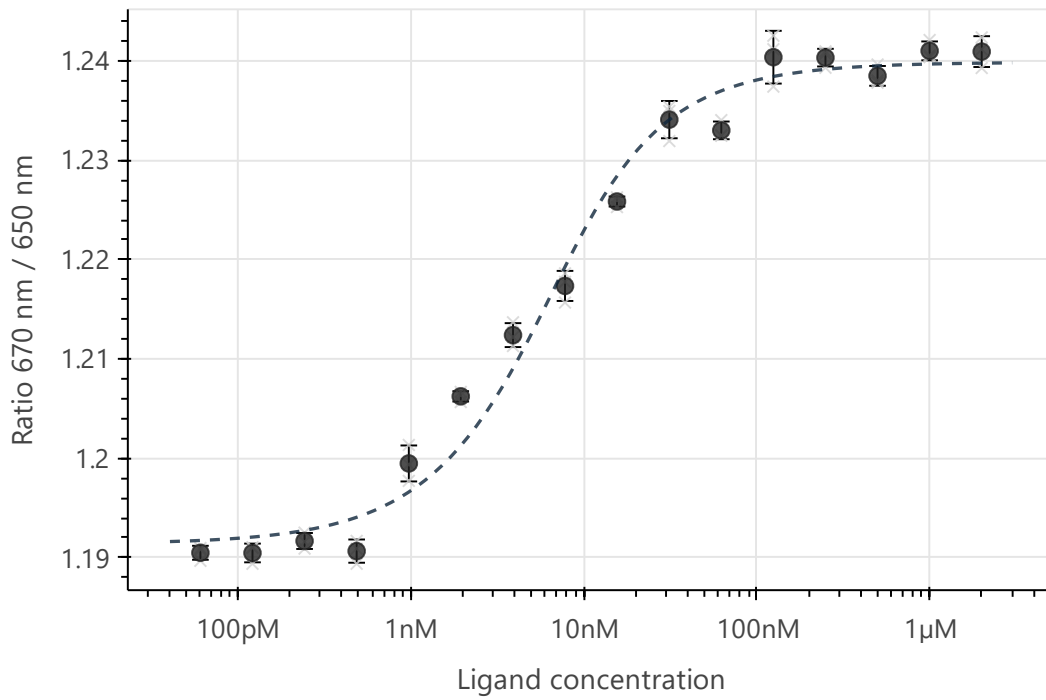
5 nM STAT3 | 2  $\mu$ M – 61 pM SD-36 | 20°C | 100% Excitation Power

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

STAT3 – SD-36 (binary complex) |  $K_d = 4.54 \pm 0.65$  nM (S/N = 28.5)

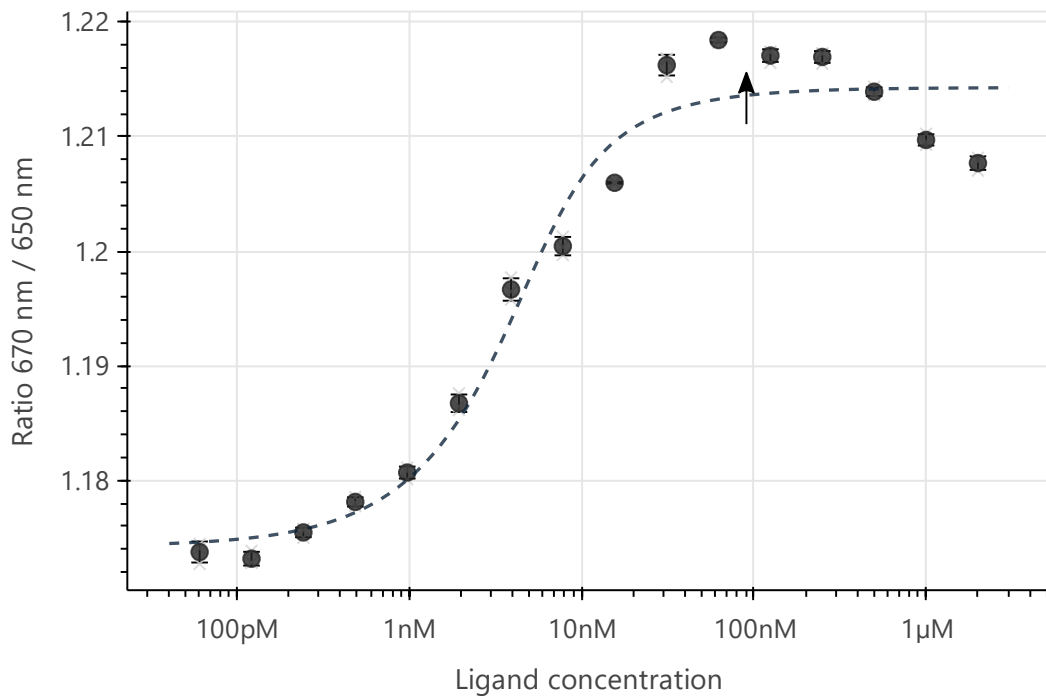


STAT3 & CRBN (500 nM) – SD-36 (ternary complex, saturation) |  $K_d = 3.64 \pm 0.82$  nM (S/N = 19.1)



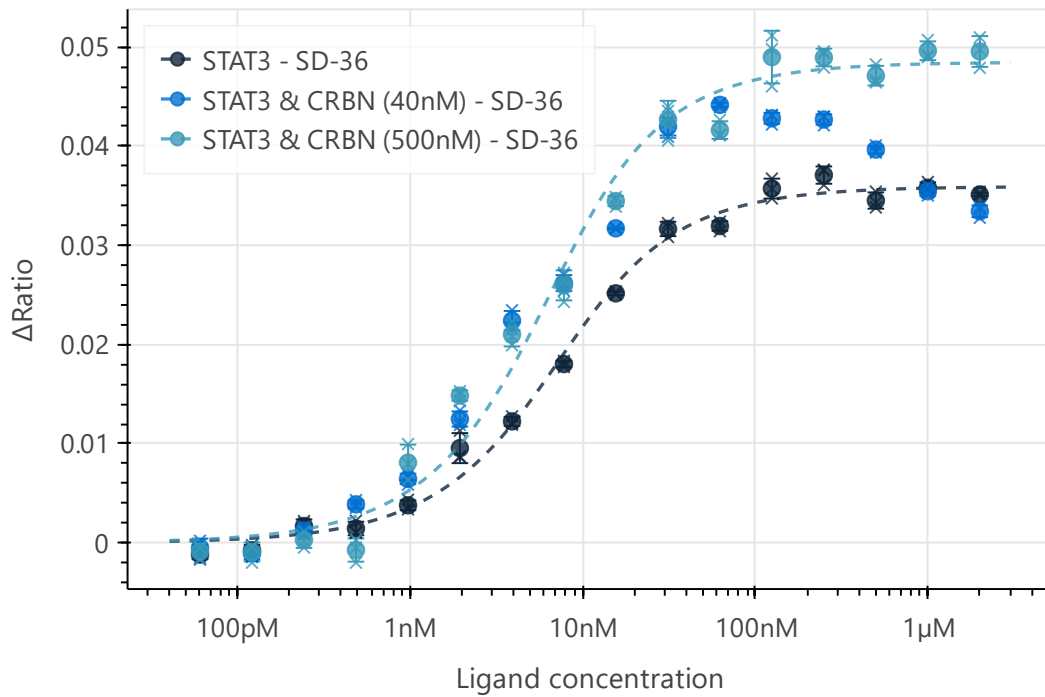
Cooperativity factor:  $\alpha = \frac{4.54 \text{ nM}}{3.64 \text{ nM}} \approx 1.2$

STAT3 & CRBN (40 nM) – SD-36 (ternary complex, Hook effect)





Overlay



### D5. Reference Results/Supporting Results

STAT3 - SD-36  $K_d = 44.4$  nM      Biolayer Interferometry (BLI)  
 $K_i = 9$  nM                      FP assay  
[Bai et al, Cancer Cell 36, 498-511 \(2019\)](#)

### E. Contributors

Andreas Langer<sup>1</sup>

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