

Monolith X Protocol MOX-P-103

Cereblon E3 Ligase – SD-36 (PROTAC) – STAT3

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

Cereblon E3 ligase complex (CRBN/DDB1)

CRBN

DDB1

uniprot.org/uniprot/Q96SW2

uniprot.org/uniprot/Q16531

A2. Molecule Class/Organism

E3 protein ligase complex

Homo sapiens (Human)

A3. Sequence/Formula

```

MGSSHHHHHH SSGENLYFQG SMEAKKPNII NFDTSLPTSH TYLGADMEEF HGRTLHDDDS CQVIPVLPQV MMILIPGQTL
PLQLFHPQEV SMVRNLIQKD RTFAVLAYSN VQEREAQFGT TAEIYAYREE QDFGIEIVKV KAIGRQRFKV LELRTQSDGI
QQAKVQILPE CVLPSTMSAV QLESLNKCQI FPSKPVSRD QCSYKWWQKY QKRKFHCANL TSWPRWLYSL YDAETLMDRI
KKQLREWDEN LKDDSLPSNP IDFSYRVAAC LPIDDLVRIQ LLKIGSAIQR LRCELDIMNK CTSLCCKQCQ ETEITTKNEI
FSLSLCGPMA AYVNPBGYVH ETLTVYKACN LNLIGRPSTE HSWFPGYAWT VAQCKICASH IGWKFATKK DMSPQKFWGL
TRSALLPTIP DTEDEISPKD VILCL

MSYNYVVTAQ KPTAVNGCVT GHFTSAEDLN LLIAKNTRLE IYVVTAEGLR PVKEVGMYGK IAVMELFRPK GESKDLFLIL
TAKYNACILE YKQSGESIDI ITRAHGNVQD RIGRPSETGI IGIIDPECRM IGLRLYDGLF KVIPLDRDNK ELKAFNIRLE
ELHVIDVKFL YGCQAPTICF VYQDPQGRHV KTYEVSLREK EFNKGPWKQE NVEAEASMVI AVPEPFGGAI IIGQESITYH
NGDKYLAIAP PIIKQSTIVC HNRVDPNGSR YLLGDMEGRL FMLLLEKEEQ MDGTVTLKDL RVELLGETSI AECLTYLDNG
VVFVGSRLGD SQLVKLVNDS NEQGSYVVM ETFTNLGPIV DMCVVDLERQ GQGQLVTCSG AFKEGSLRII RNGIGIHEHA
SIDLPGIKGL WPLRSDPNRE TDDTLVLSFV GQTRVLMNG EEEVEELMG FVDDQQTFFC GNVAHQQLIQ ITSASVRLVS
QEPKALVSEW KEPQAKNISV ASCNSSQVV AVGRALYYLQ IHPQELRQIS HTEMEHEVAC LDITPLGDSN GLSPLCAIGL
WTDISARILK LPSFELLHKE MLGGEIIPRS ILMTTFESSH YLLCALGDGA LFYFGLNIET GLLSDRKKVT LGTQPTVLR
FRSLSTTVNF ACSDRPTVIY SSNHKLVSFN VNLKEVNYMC PLNSDGYPDS LALANNSTLT IGTIDEIQKL HIRTVPLYES
PRKICYQEVS QCFGVLSRI EVQDTSGGTT ALRPSASTQA LSSSVSSSKL FSSSTAPHET SFGEEVEVHN LLIIDQHTFE
VLHAHQFLQN EYALSLVSK LGKDPNTYFI VGTAMVYPEE AEPKQGRIVV FQYSDGKLQT VAEKEVKGAV YSMVEFNGKL
LASINSTVRL YEWTTEKELR TECNHYNNIM ALYLKTKGDF ILVGDLMRSV LLLAYKPMEG NFEEIARDFN PNWMSAVEIL
DDDNFLGAEN AFNLFVCQKD SAATTDEERQ HLQEVGLFHL GEFVNVFCHG SLVMQNLGET STPTQGSVLF GTVNGMIGLV
TSLSESWYNL LLDQMRLNKK VIKSVGKIEH SFWRSFHTER KTEPATGFID GDLESFLDI SRPKMQEVVA NLQYDDGSGM
KREATADDLI KVVEELTRIH

```

A4. Purification Strategy/Source

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

A5. Stock Concentration/Stock Buffer

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

A6. Molecular Weight/Extinction Coefficient

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

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-L011, NanoTemper Technologies GmbH)
1* Dye RED-NHS 2nd Generation (10 μ g) | 1* B-Column

A9. Labeling Procedure

1. Mix 5 μ L of 37.2 μ M CRBN/DDB1 with 15 μ L of dilution buffer to obtain 20 μ L of ~10 μ M CRBN/DDB1.
2. Add 25 μ L of DMSO to Dye RED-NHS 2nd Generation (10 μ g) to obtain a ~600 μ M solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
3. Mix 1 μ L of the 600 μ M dye solution with 19 μ L of dilution buffer to obtain 20 μ L of a 30 μ M dye solution (3x protein concentration).
4. Mix CRBN/DDB1 and dye in a 1:1 volume ratio (40 μ L final volume, 2.5% final DMSO concentration).
5. Incubate for 30 minutes at room temperature in the dark.
6. 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.
7. 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.
8. Add 40 μ L of the labeling reaction from step 4 to the center of the column and let sample enter the resin bed completely.
9. Add 550 μ L of dilution buffer after the sample has entered and discard the flow through.
10. Place column in a new collection tube, add 400 μ L of dilution buffer and collect the eluate.
11. Prepare 8 μ L aliquots of the labeled CRBN/DDB1 (~0.5 μ M) and immediately store at -80°C.

A10. Labeling Efficiency

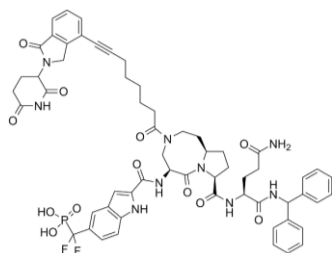
Measurement of protein concentration and degree of labeling (DOL) using a NanoDrop™:

nanotempertech.com/dol-calculator

Absorbance A_{280}	0.100	Protein concentration	0.57 μ M
Absorbance A_{650}	0.118	Degree-of-labeling (DOL)	1.06

B1. Ligand/Non-Fluorescent Binding Partner

SD-36



Signal transducer and activator of transcription 3 (STAT3)

uniprot.org/uniprot/P40763

B2. Molecule Class/Organism

PROTAC

Transcription factor

Homo sapiens (Human)

B3. Sequence/Formula

C₅₉H₆₂F₂N₉O₁₂P

```

GQANHPTAAV VTEKQQMLEQ HLQDVRKRVQ DLEQMKVVE NLQDDDFDNY
KTLKSQGMQ DLNGNQSVT RQKMQQLEQM LTALDQMRRS IVSELAGLLS
AMEYVQKTLT DEELADWKRR QQIACIGGPP NICLDRLLENW ITSLAESQLQ
TRQQIKKLEE LQQKVSYKGD PIVQHRPMLE ERIVELFRNL MKSAFVVERQ
PCMPMHPDRP LVIKTGVQFT TKVRLLVKFP ELNYQLKIKV CIDKDSGDVA
ALRGSRKFN I LGTNTKVMNM EESNNGLSA EFKHLTLREQ RCGNGGRANC
DASLIVTEEL HLITFETEVY HQGLKIDLET HSLPVVVISN ICQMPNAWAS
ILWYNMLTNN PKNVNFFTP PIGTWDQVAE VLSWQFSSTT KRGLSIEQLT
TLAEKLLGPG VNYSQCITW AKFCKENMAG KGFSFWVWLD NIIDLVKKYI
LALWNEGYIM GFISKERERA ILSTKPPGTF LLRFSESSKE GGVFTFWVEK
DISGKTQIQS VEPYTKQQLN NMSFAEIIMG YKIMDATNIL VSPLVLYLPD
IPKEEAFGKY CRPESQEHP E ADPGSAAPYL KTKFICVTPG SGENLYFQGH HHHHH
    
```

B4. Purification Strategy/Source

SD-36
MedChemExpress
[HY-129602](#)

Recombinant
CreLux GmbH

B5. Stock Concentration/Stock Buffer

11.6 mg/mL | 10 mM
DMSO

12.87 mg/mL | 195 µM
100 mM Tris-HCl pH 8.5, 100 mM NaCl, 2 mM DTT

B6. Molecular Weight/Extinction Coefficient

1158.15 Da

66.1 kDa

84,340 M⁻¹cm⁻¹ (ε₂₈₀)

B7. Serial Dilution Preparation

1. Dissolve 5 mg of SD-36 in 432 μL of DMSO to obtain a 10 mM stock solution.
2. Mix 2 μL of 10 mM SD-36 with 98 μL of DMSO to obtain 100 μL of a 200 μM SD-36 solution.
3. Mix 4 μL of 200 μM SD-36 with 196 μL of dilution buffer to obtain 100 μL of a 4 μM SD-36 solution.
4. Mix 20 μL of DMSO with 980 μL of dilution buffer to obtain 1 mL of dilution buffer with 2% DMSO.
5. Mix 4 μL of labeled CRBN/DDB1 (~0.5 μM) with 96 μL of dilution buffer to obtain 200 μL of 20 nM CRBN/DDB1.
6. Mix 1 μL of 195 μM STAT3 with 121 μL of dilution buffer to obtain 122 μL of 1.6 μM STAT3.

CRBN – SD-36 (binary complex)

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

CRBN & STAT3 – SD-36 (ternary complex)

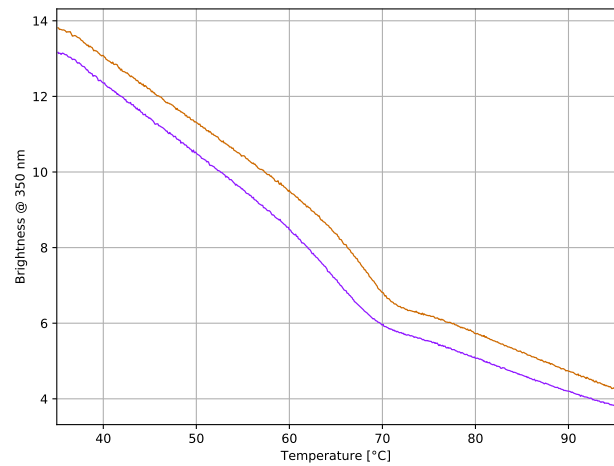
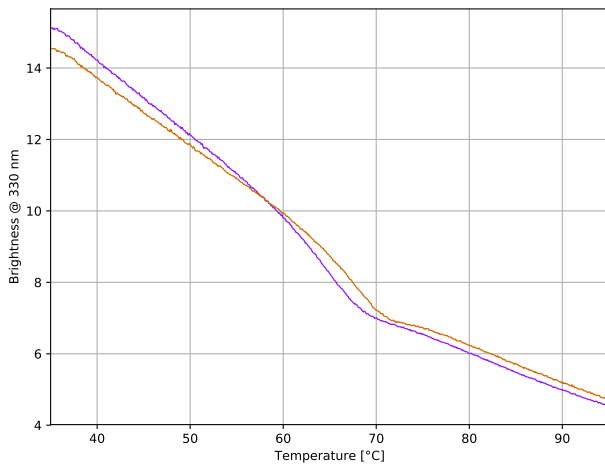
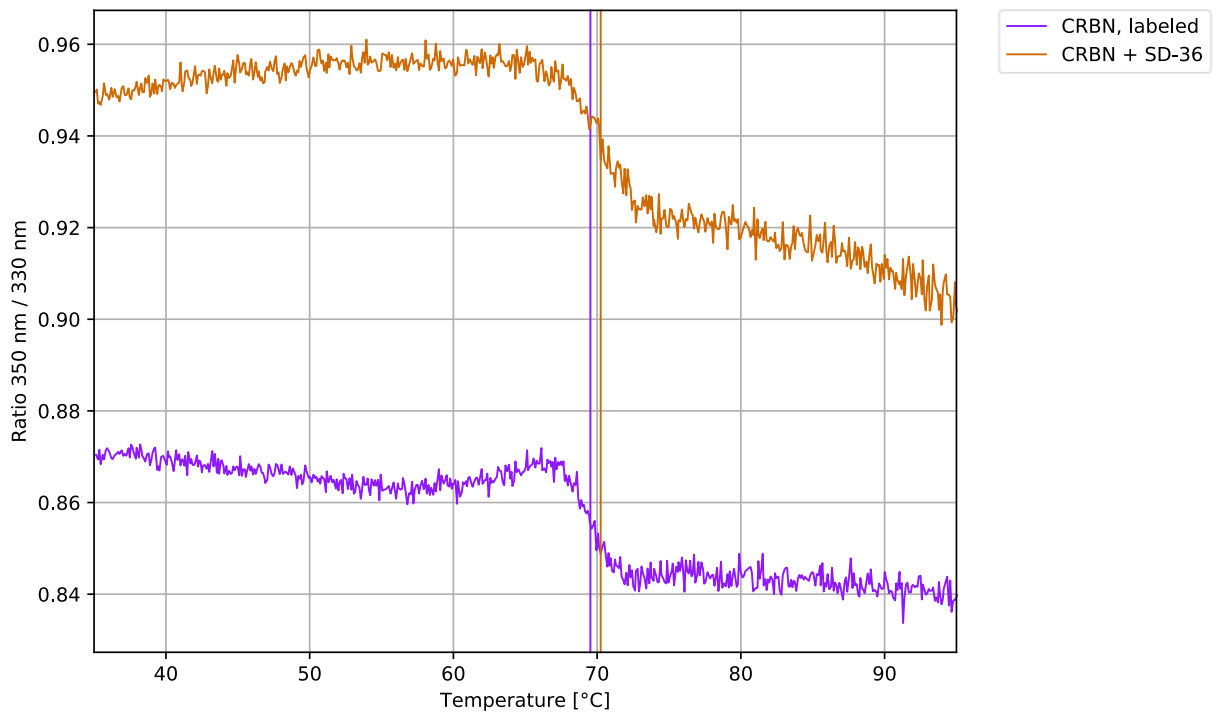
12. Mix 100 μL of 20 nM CRBN with 100 μL of 1.6 μM STAT3 to obtain 200 μL of 10 nM CRBN, 800 nM STAT3.
13. Take a fresh 0.5 mL tube and mix 160 μL of 10 nM CRBN, 800 nM STAT3 with 160 μL of dilution buffer containing 2% DMSO to obtain 320 μL of a 5 nM CRBN, 400 nM STAT3 solution.
14. Prepare a PCR-rack with 20 PCR tubes. Transfer 20 μL of the 5 nM CRBN, 400 nM STAT3 solution into tubes **2** to **16**. Then, mix 20 μL of 4 μM SD-36 with 20 μL of 10 nM CRBN, 800 nM STAT3 in tube **1**.
15. Prepare a 1:1 serial dilution by transferring 20 μL from tube to tube. Mix carefully by pipetting up and down.
16. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

C. Tycho

Validation of structural integrity and functionality of labeled CRBN-DDB1 using Tycho NT.6:

nanotempertech.com/tycho

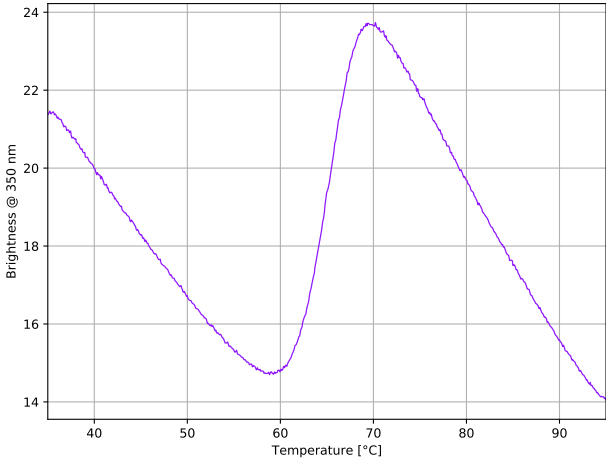
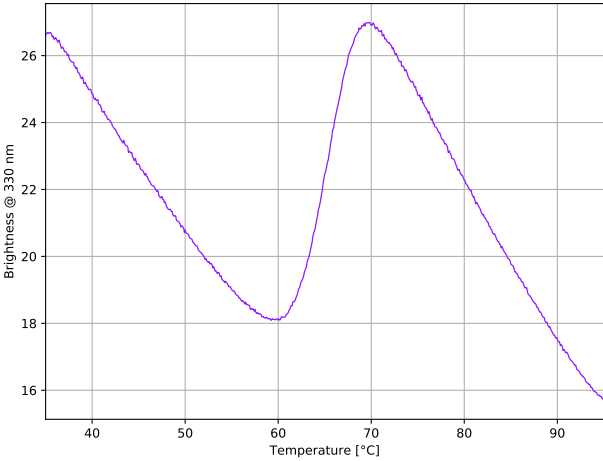
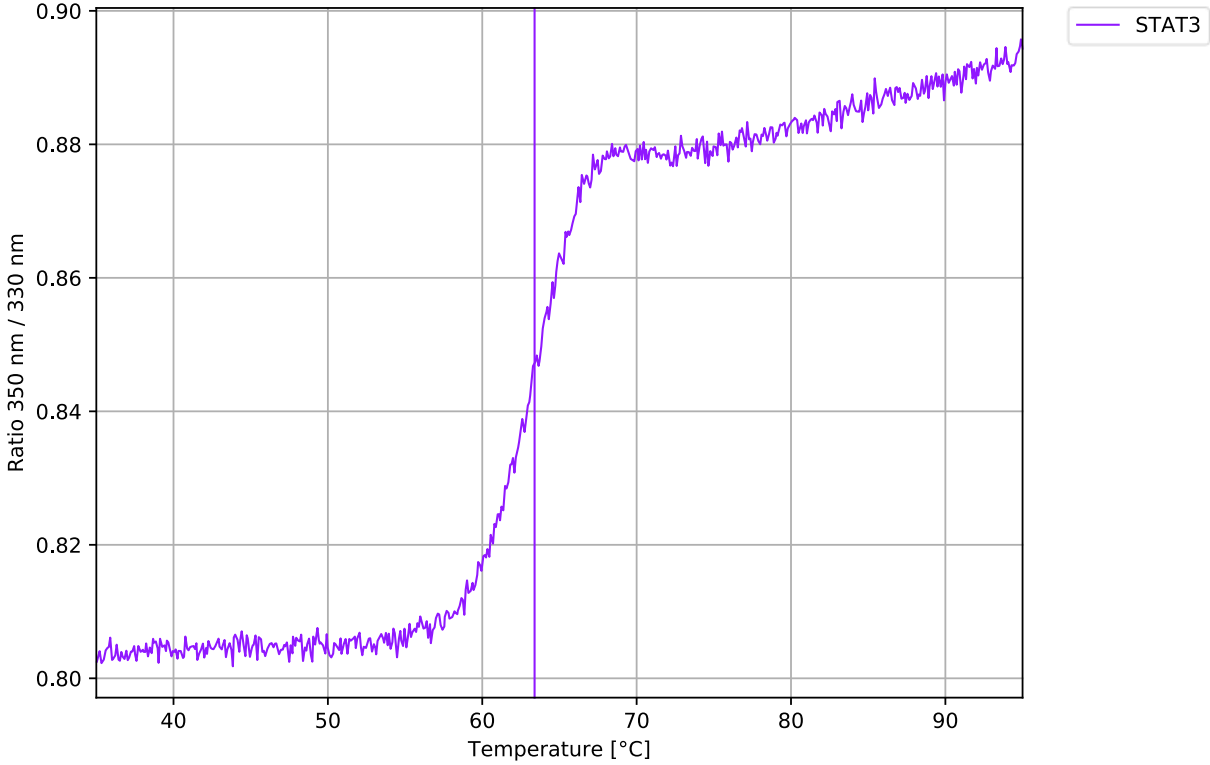
CRBN, labeled	8 μ L of B-Column eluate ($\sim 0.5 \mu$ M) + 2 μ L of dilution buffer with 2% DMSO	$T_i = 69.5^\circ\text{C}$
CRBN + SD-36	8 μ L of B-Column eluate ($\sim 0.5 \mu$ M) + 2 μ L of 4 μ M SD-36	$T_i = 70.2^\circ\text{C}$



Validation of structural integrity of STAT3 using Tycho NT.6:

nonotempertech.com/tycho

STAT3	10 μ L of 1.6 μ M STAT3	$T_i = 63.4^\circ\text{C}$
-------	---------------------------------	----------------------------



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

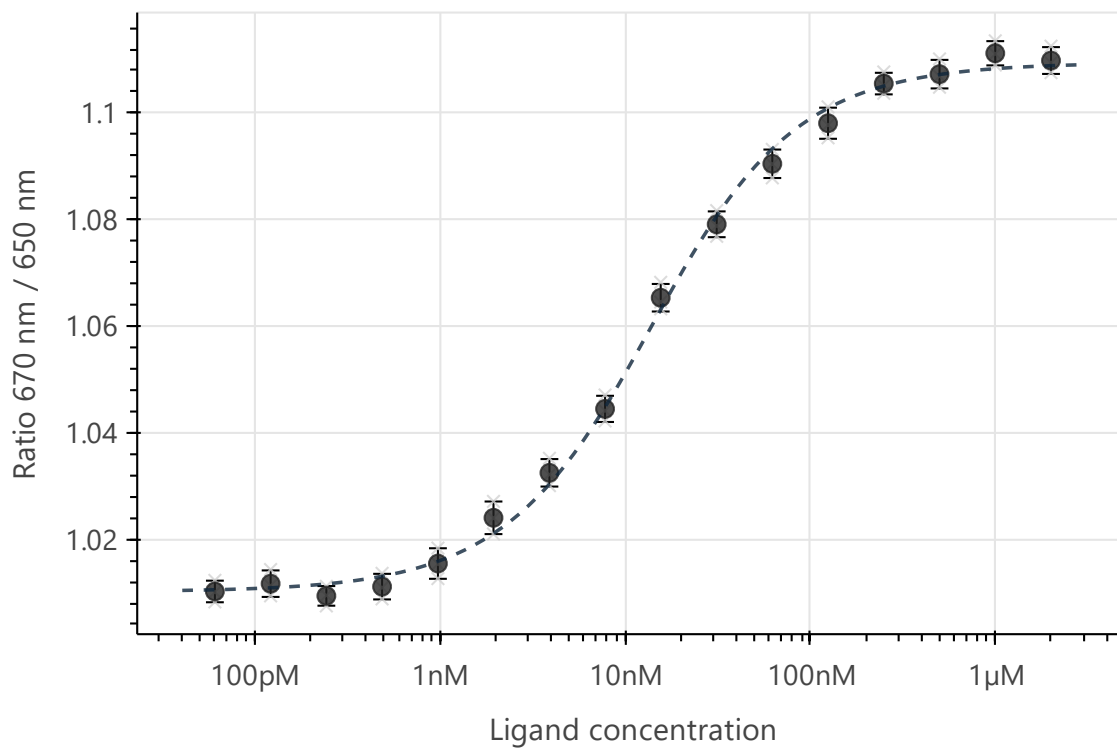
D3. Monolith Experiment (Assay Buffer/Concentrations/Temperature/Excitation Power)

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

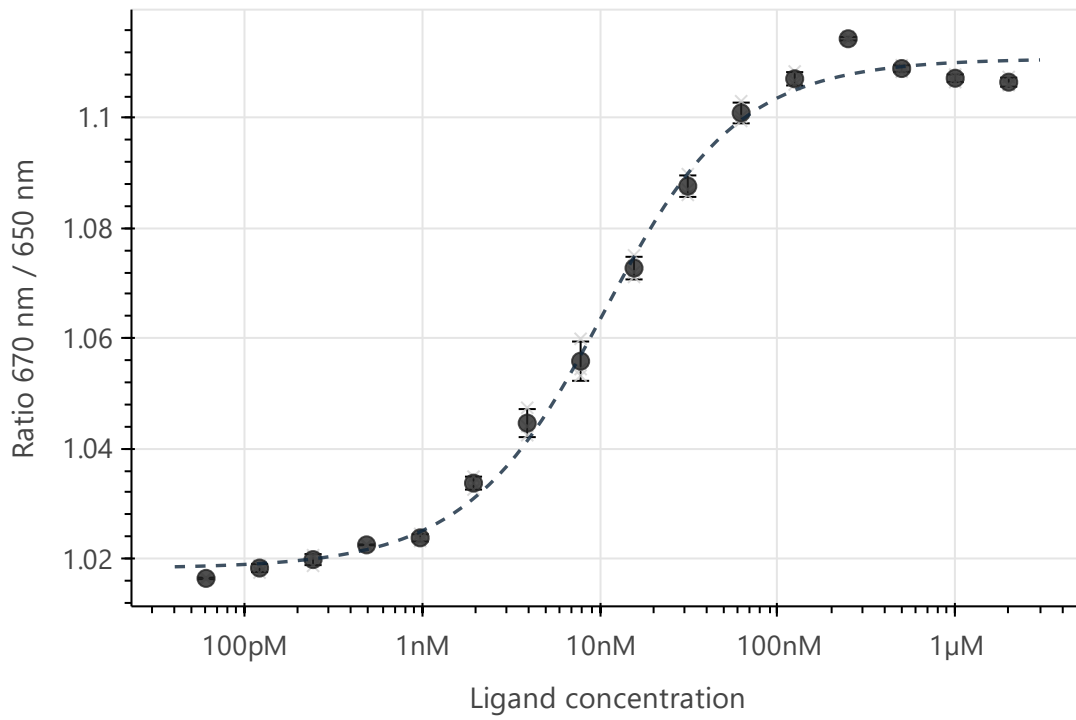
5 nM CRBN-DDB1 | 2 μ M – 61 pM SD-36 | 20°C | 100% Excitation Power

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

CRBN – SD-36 (binary complex) | $K_d = 11.3 \pm 0.8$ nM (S/N = 51.2)

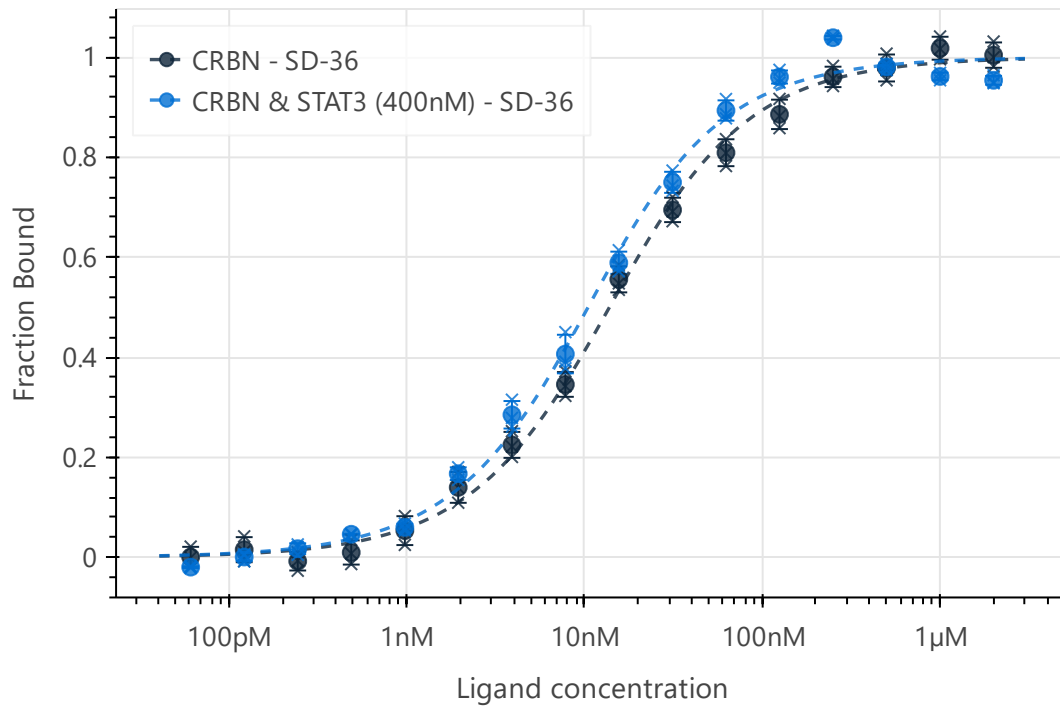


CRBN & STAT3 – SD-36 (ternary complex) | $K_d = 7.98 \pm 0.87$ nM (S/N = 34.3)



Cooperativity factor: $\alpha = \frac{11.3 \text{ nM}}{7.98 \text{ nM}} \approx 1.4$

Overlay



D5. Reference Results/Supporting Results

N/A

E. Contributors

Andreas Langer¹

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