

Monolith X Protocol MOX-P-101

BRD2^{BD2} – MZ1 (PROTAC) – VCB (VHL) E3 Ligase

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. The von Hippel-Lindau E3 Ligase (VHL) and the elongin BC proteins form together the functional VCB complex which is able to use the PROTAC MZ1 to bind to the bromodomain of BRD2.

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

A1. Target/Fluorescent Molecule

BRD2^{BD2}

uniprot.org/uniprot/P25440

A2. Molecule Class/Organism

Bromo-domain containing protein

A3. Sequence/Formula

MHHHHHSGV DLGTENLYFQ SMGKLSEQLK HCNGILKELL SKKHAAYAWP FYKPVDSAL GLHDYHDIK HPMDLSTVKR
KMENRDYRDA QEFAADVRLM FSNCYKYNPP DHDVVAMARK LQDVFEFRYA KMPD

A4. Purification Strategy/Source

Crelux GmbH

Construct ID: CRN1, Lot-ID: PC16670

A5. Stock Concentration/Stock Buffer

5.72 mg/mL | 362 μ M

25 mM HEPES pH 7.5, 100 mM NaCl, 5% glycerol, 1 mM DTT

A6. Molecular Weight/Extinction Coefficient

15.8 kDa

17,420 $M^{-1}cm^{-1}$ (ϵ_{280})

A7. Dilution Buffer

50 mM Tris-HCl, pH 7.8, 150 mM NaCl, 10 mM MgCl₂, 2 mM GSH, 0.01 % Pluronic® F-127

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 50 µL of 25 µM BRD2^{BD2} by mixing 3.45 µL of 362 µM BRD2^{BD2} with 46.5 µL of Labeling Buffer NHS.
2. 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.
3. Mix 5 µL of the 600 µM dye solution with 45 µL of Labeling Buffer NHS to obtain 50 µL of a 60 µM dye solution (2.4x protein concentration).
4. Mix BRD2^{BD2} and dye in a 1:1 volume ratio (100 µL final volume, 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 100 µL of the labeling reaction from step 4 to the center of the column and let sample enter the resin bed completely.
9. Add 550 µL of dilution buffer after the sample has entered and discard the flow through.
10. Place column in a new collection tube, add 400 µL of dilution buffer and collect the eluate.
11. Centrifuge the eluate 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.
12. Prepare 8 µL aliquots of the labeled BRD2BD2 (~2.5 µM) and immediately store at -80°C.

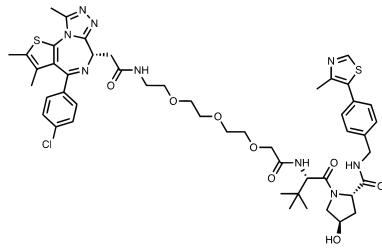
A10. Labeling Efficiency

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

Absorbance A ₂₈₀	0.047	Protein concentration	1.44 µM
Absorbance A ₆₅₀	0.461	Degree-of-labeling (DOL)	1.44

B1. Ligand/Non-Fluorescent Binding Partner

MZ1



Von Hippel-Lindau E3 Ligase (VHL)

uniprot.org/uniprot/P40337

Elongin B (EloB)

uniprot.org/uniprot/Q15370

Elongin C (EloC)

uniprot.org/uniprot/Q15369

B2. Molecule Class/Organism

PROTAC

E3 protein ligase complex

Homo sapiens (Human)

B3. Sequence/Formula

C₄₉H₆₀ClN₉O₈S₂

```
MPRRAENWDE AEVGAEAGV EEYGPEEDGG EESGAEESGP EESGPEELGA
EEEMEAGRPR PVLRSVNSRE PSQVIFCNRS PRVVLPVWLN FDGEPQPYPT
LPPGTGRRIH SYRGHLWLFV DAGTHDGLLV NQTELFVPSL NVDGQPIFAN
ITLPVYTLKE RCLQVVRSLV KPENYRRLDI VRSLYEDLED HPNVQKDLER
LTQERIAHQV MGD
MDVFLMIRRH KTTIFDAKE SSTVFELKRI VEGILKRPPD EQRLYKDDQL
LDDGKTLGEC GFTSQTARPQ APATVGLAFR ADDTFEALCI EPFSSPPELP
DVMK
MYVKLISSDG HEFIVKREHA LTSGTIKAML SGPGQFAENE TNEVNFREIP
SHVLSKVCMI FTYKVRVYNS STEIPEFPPIA PEIALELLMA ANFLDC
```

B4. Purification Strategy/Source

opnme.com

[MZ1](#)

CreLux GmbH

Construct ID: ESA2, Lot-ID: PH14288-1

B5. Stock Concentration/Stock Buffer

10 mg/mL | 10 mM

DMSO

10.3 mg/mL | 206 μM

20 mM Bis-Tris/HCl, 150 mM NaCl, 1 mM DTT, pH 7.0

B6. Molecular Weight/Extinction Coefficient

1002.64 Da

24 kDa (VHL), 13.1 kDa (EloB), 12.8 kDa (EloC)
25,900 M⁻¹cm⁻¹ (ε₂₈₀)

B7. Serial Dilution Preparation

1. Dissolve 4.9 mg of MZ1 in 489 μL of DMSO to obtain a 10 mM stock solution.
2. Mix 2 μL of 10 mM MZ1 with 98 μL of DMSO to obtain 200 μL of 200 μM MZ1.
3. Mix 4 μL of 200 μM MZ1 with 196 μL of dilution buffer to obtain 200 μL of 4 μM MZ1.
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 BRD2^{BD2} (~2.5 μM) with 496 μL of dilution buffer to obtain 500 μL of ~20 nM BRD2^{BD2}.
6. Prepare a 4 μM VCB solution by mixing 2 μL of 206 μM VCB with 101 μL of dilution buffer.

BRD2^{BD2} – MZ1 (binary complex)

7. Mix 100 μL of 20 nM BRD2^{BD2} with 100 μL of dilution buffer to obtain 200 μL of 10 nM BRD2^{BD2}.
8. Take a fresh 0.5 mL tube and mix 160 μL of 10 nM BRD2^{BD2} with 160 μL of dilution buffer containing 2% DMSO to obtain 320 μL of a 5 nM BRD2^{BD2} solution.
9. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of the 5 nM BRD2^{BD2} solution into tubes **2** to **16**. Then, mix 20 μL of 4 μM MZ1 with 20 μL of 10 nM BRD2^{BD2} 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.

BRD2^{BD2} & high concentration of VCB – MZ1 (ternary complex, saturation)

12. Mix 100 μL of 20 nM BRD2^{BD2} with 50 μL of 4 μM VCB and 50 μL of dilution buffer to obtain 200 μL of 10 nM BRD2^{BD2}, 1 μM VCB.
13. Take a fresh 0.5 mL tube and mix 160 μL of 10 nM BRD2^{BD2}, 1 μM VCB with 160 μL of dilution buffer containing 2% DMSO to obtain 320 μL of a 5 nM BRD2^{BD2}, 500 nM VCB solution.
14. Prepare a PCR-rack with 20 PCR tubes. Transfer 20 μL of the 5 nM BRD2^{BD2}, 500 nM VCB solution into tubes **2** to **16**. Then, mix 20 μL of 4 μM MZ1 with 20 μL of 10 nM BRD2^{BD2}, 1 μM VCB 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.

BRD2^{BD2} & low concentration of VCB – MZ1 (ternary complex, Hook effect)

17. Mix 100 μL of 20 nM BRD2^{BD2} with 2.5 μL of 4 μM VCB and 97.5 μL of dilution buffer to obtain 200 μL of 10 nM BRD2^{BD2}, 50 nM VCB.
18. Take a fresh 0.5 mL tube and mix 160 μL of 10 nM BRD2^{BD2}, 50 nM VCB with 160 μL of dilution buffer containing 2% DMSO to obtain 320 μL of a 5 nM BRD2^{BD2}, 25 nM VCB solution.
19. Prepare a PCR-rack with 20 PCR tubes. Transfer 20 μL of the 5 nM BRD2^{BD2}, 25 nM VCB solution into tubes **2** to **16**. Then, mix 20 μL of 4 μM MZ1 with 20 μL of 10 nM BRD2^{BD2}, 50 nM VCB in tube **1**.
20. Prepare a 1:1 serial dilution by transferring 20 μL from tube to tube. Mix carefully by pipetting up and down.
21. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

BRD2^{BD2} & low concentration of VCB – cis MZ1 (Control)

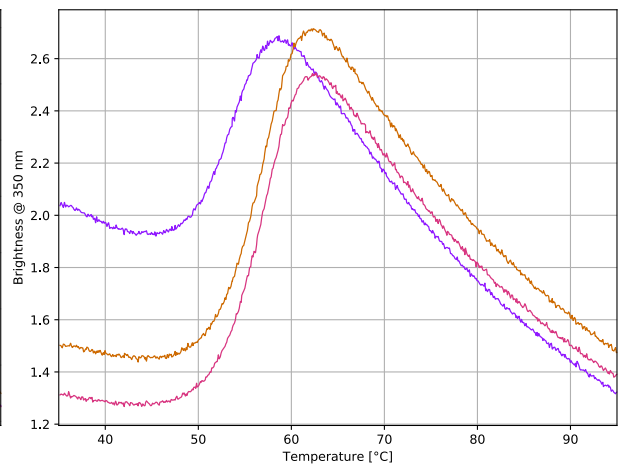
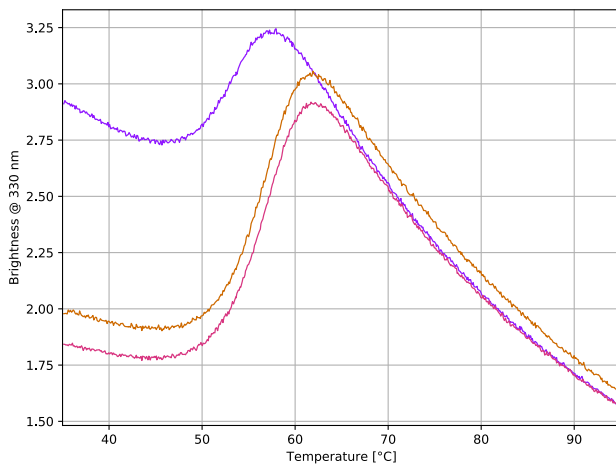
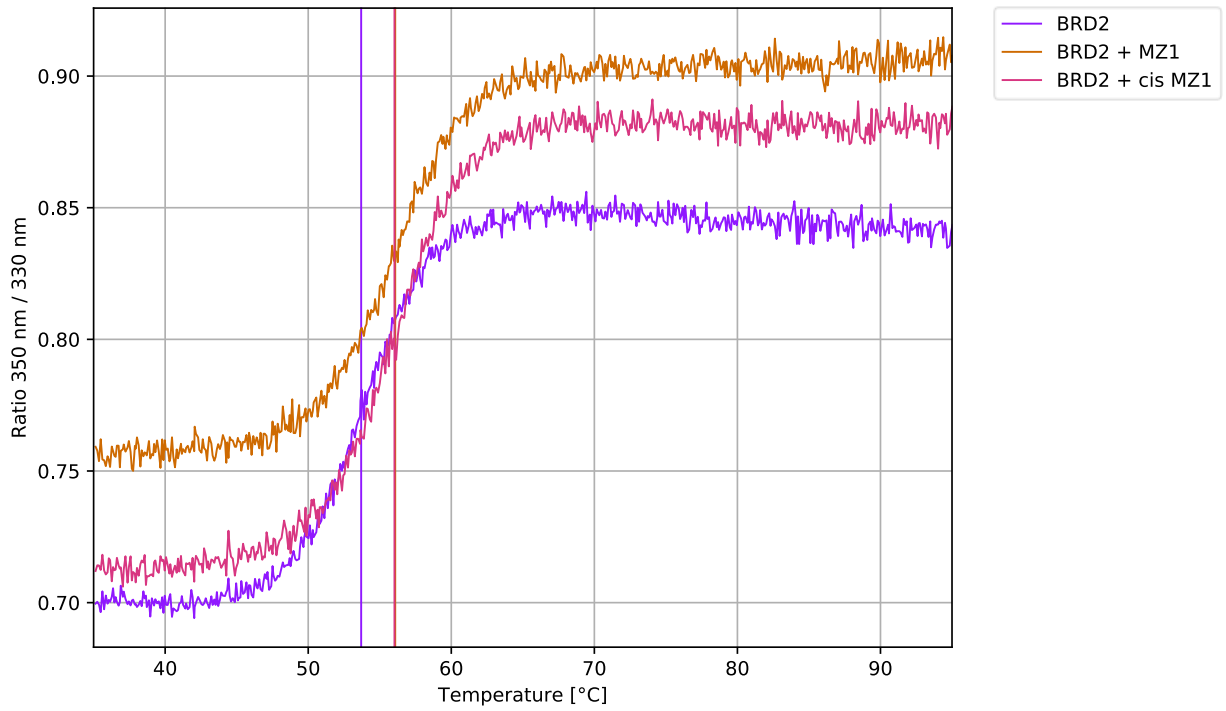
22. Mix 100 μL of 20 nM BRD2^{BD2} with 2.5 μL of 4 μM VCB and 97.5 μL of dilution buffer to obtain 200 μL of 10 nM BRD2^{BD2}, 50 nM VCB.
23. Take a fresh 0.5 mL tube and mix 160 μL of 10 nM BRD2^{BD2}, 50 nM VCB with 160 μL of dilution buffer containing 2% DMSO to obtain 320 μL of a 5 nM BRD2^{BD2}, 25 nM VCB solution.
24. Prepare a PCR-rack with 20 PCR tubes. Transfer 20 μL of the 5 nM BRD2^{BD2}, 25 nM VCB solution into tubes **2** to **16**. Then, mix 20 μL of 4 μM cis MZ1 with 20 μL of 10 nM BRD2^{BD2}, 50 nM VCB in tube **1**.
25. Prepare a 1:1 serial dilution by transferring 20 μL from tube to tube. Mix carefully by pipetting up and down.
26. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

C. Tycho

Validation of structural integrity and functionality of BRD2^{BD2} using Tycho NT.6:

nanotempertech.com/tycho

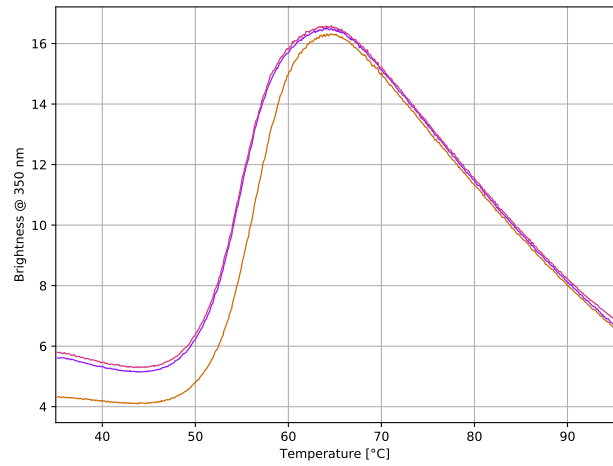
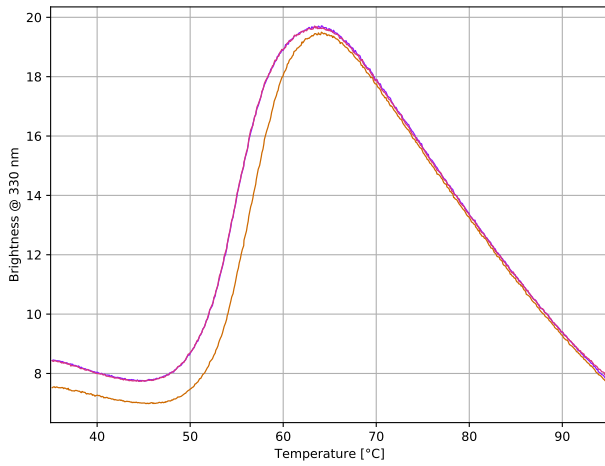
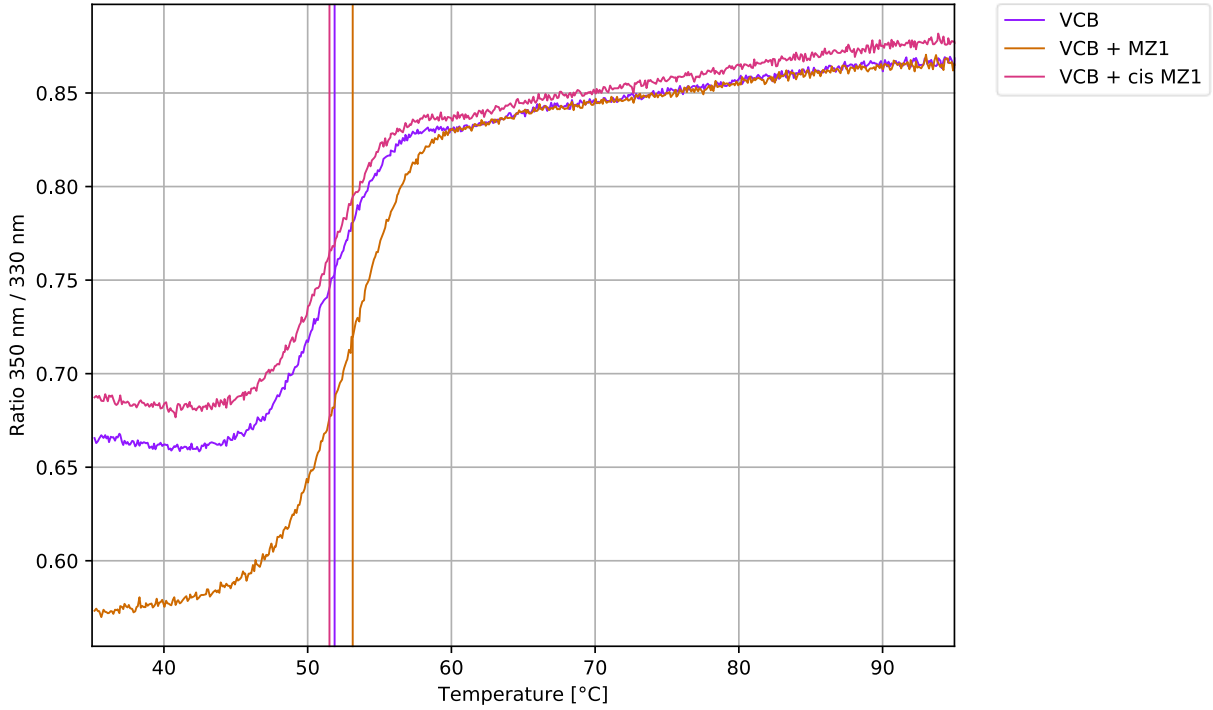
BRD2	5 μ L of 4 μ M BRD2 ^{BD2} + 5 μ L of dilution buffer containing 2% DMSO	T _i = 53.7°C
BRD2 + MZ1	5 μ L of 4 μ M BRD2 ^{BD2} + 5 μ L of 4 μ M MZ1	T _i = 56.1°C
BRD2 + cis MZ1	5 μ L of 4 μ M BRD2 ^{BD2} + 5 μ L of 4 μ M cis MZ1	T _i = 56.0°C



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

nonotempertech.com/tycho

VCB	5 μ L of 4 μ M VCB + 5 μ L of dilution buffer containing 2% DMSO	T _i = 51.9°C
VCB + MZ1	5 μ L of 4 μ M VCB + 5 μ L of 4 μ M MZ1	T _i = 53.1°C
VCB + cis MZ1	5 μ L of 4 μ M VCB + 5 μ L of 4 μ M cis MZ1	T _i = 51.5°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)

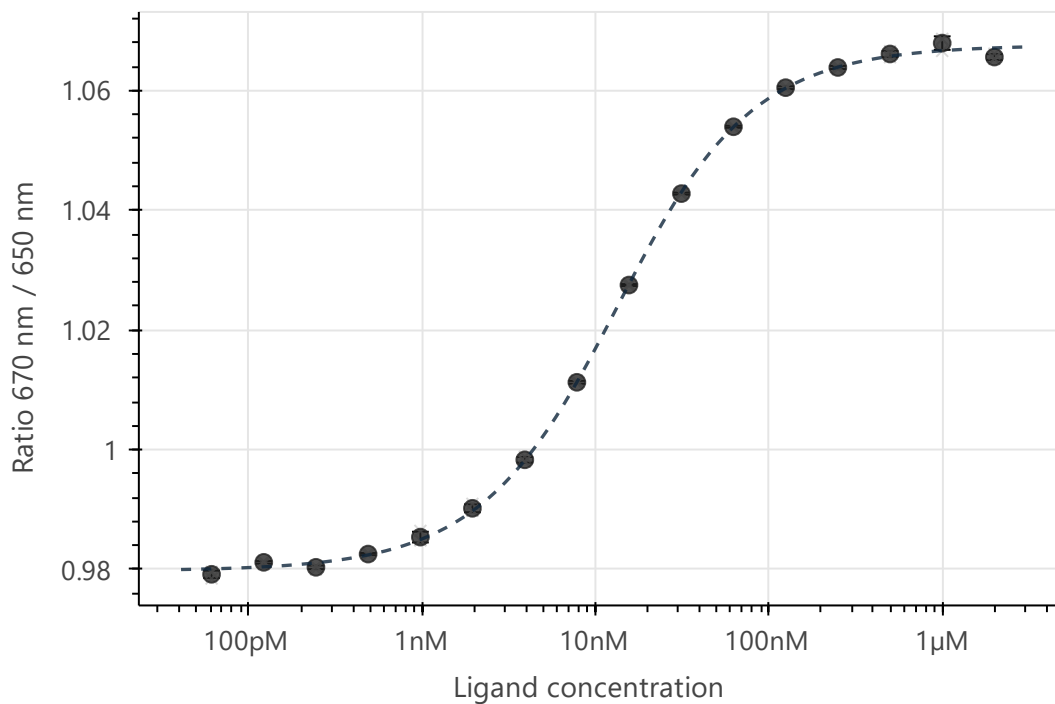
nanotempertech.com/monolith-mo-control-software

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

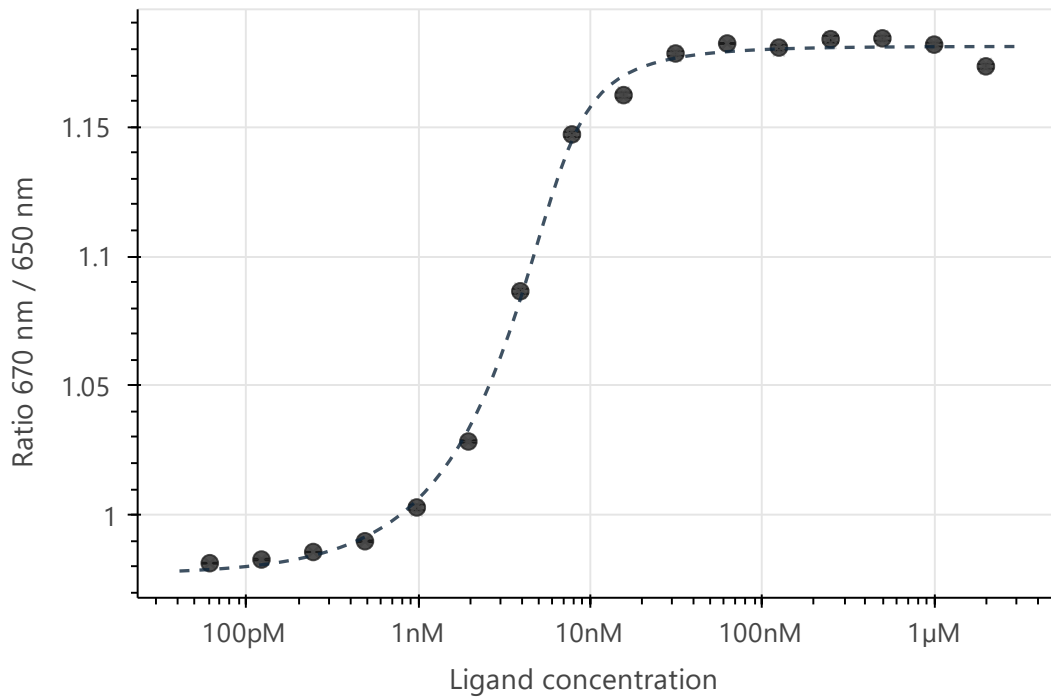
50 mM Tris-HCl, pH 7.8, 150 mM NaCl, 10 mM MgCl₂, 2 mM GSH, 0.01 % Pluronic® F-127, 1% DMSO
 5 nM BRD2^{BD2} | 2 μM – 61 pM MZ1 | 20°C | 100% Excitation Power

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

BRD2^{BD2} – MZ1 (binary complex) | $K_d = 10.7 \pm 0.2$ nM (S/N = 132.3)

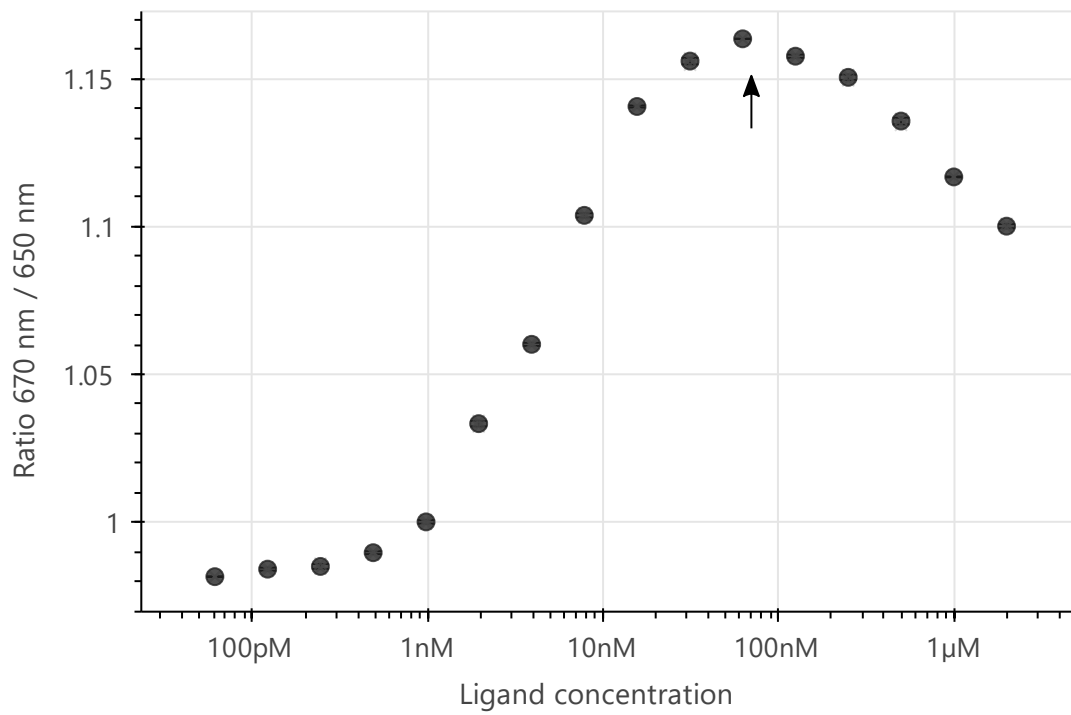


BRD2^{BD2} & VCB (500 nM) – MZ1 (ternary complex, saturation) | $K_d = 566 \pm 94$ pM (S/N = 54.4)

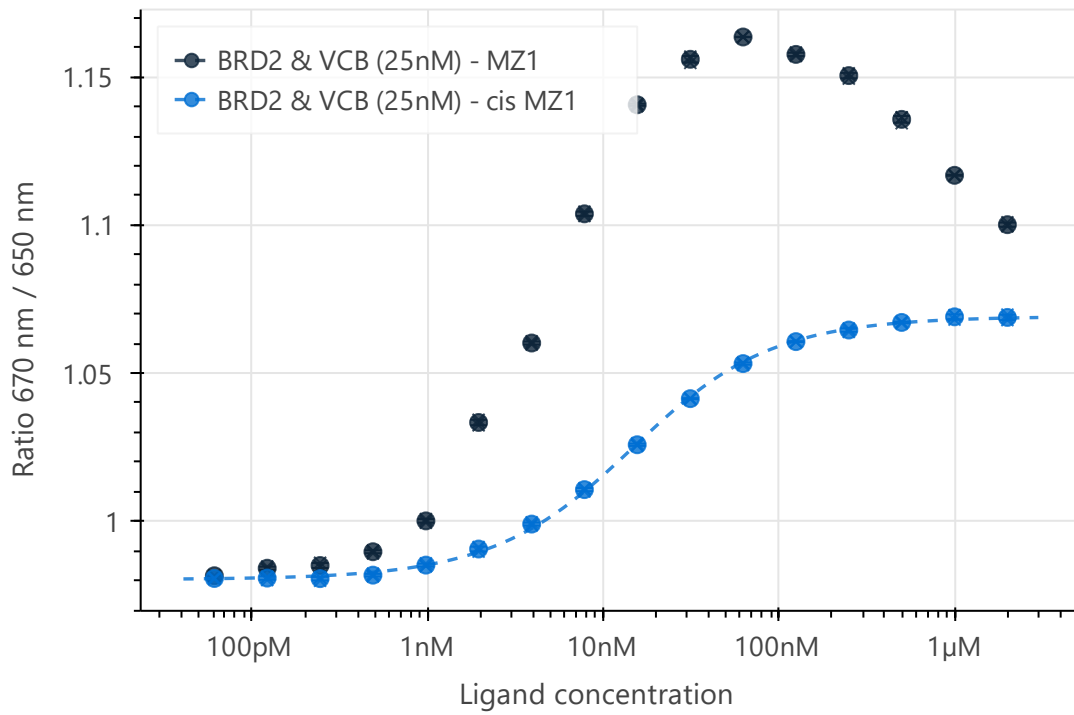


Cooperativity factor: $\alpha = \frac{10.7 \text{ nM}}{0.566 \text{ nM}} \approx 18.9$

BRD2^{BD2} & VCB (25 nM) – MZ1 (ternary complex, Hook effect)

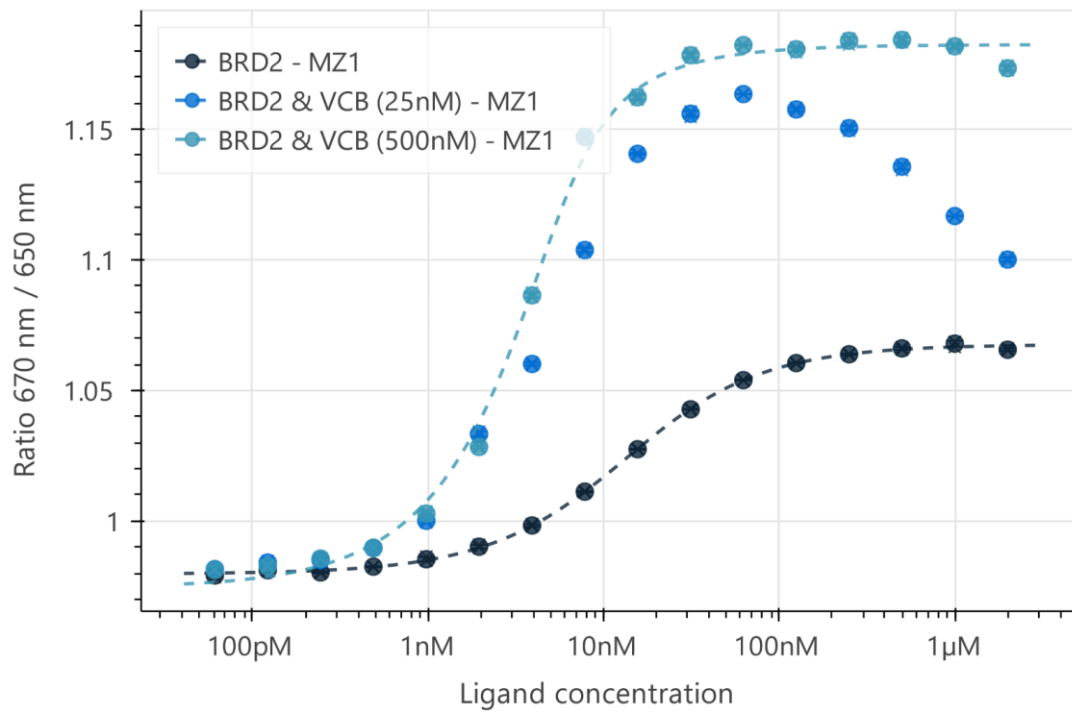


BRD2^{BD2} & VCB (25 nM) – cis MZ1 (control) | $K_d = 12.2 \pm 0.4$ nM (S/N = 124)



No Hook effect and ternary complex formation

Overlay plot:



D5. Reference Results/Supporting Results

PROTAC		+ target	Monolith X		ITC	
			K_d (nM)	α^1	K_d (nM)	α
MZ1	binary	-	10.7	-	60	-
MZ1	ternary	VCB	0.566	18.9	-	-

Isothermal Titration Calorimetry (ITC)

[Gadd et al., Nat Chem Biol 13, 514–521 \(2017\)](#)

E. Contributors

Andreas Langer²

¹ Cooperativity factor

² NanoTemper Technologies GmbH, München, Germany | nanotempertech.com