

Monolith X Protocol MOX-P-102

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

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

Von Hippel-Lindau E3 Ligase (VHL)

uniprot.org/uniprot/P40337

Elongin B (EloB)

uniprot.org/uniprot/Q15370

Elongin C (EloC)

uniprot.org/uniprot/Q15369

A2. Molecule Class/Organism

E3 protein ligase complex

Homo sapiens (Human)

A3. Sequence/Formula

```
MPRRAENWDE AEVGAEAGV EEYGPEDGG EESGAEESGP EESGPEELGA EEEMEAGRPR PVLRSVNSRE PSQVIFCNRS
PRVVLPPVWLN FDGEPQPYPT LPPGTGRRIH SYRGHLWLFK DAGTHDGLLV NQTELFVPSL NVDGQPIFAN ITLPVYTLKE
RCLQVVRSLV KPENYRRLDI VRSLYEDLED HPNVQKDLER LTQERIAHQK MGD
MDVFLMIRRH KTTIFTDAKE SSTVFELKRI VEGILKRPPD EQRLYKDDQL LDDGKTLGEC GFTSQTARPK APATVGLAFR
ADDTFEALCI EPFSSPELPEL DVMK
MYVKLISSDG HEFIVKREHA LTSGTIKAML SGPGQFAENE TNEVNFREIP SHVLSKVCMY FTYKVRYTNS STEIPEFPIA
PEIALELLMA ANFLDC
```

A4. Purification Strategy/Source

CreLux GmbH

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

A5. Stock Concentration/Stock Buffer

10.3 mg/mL | 206 µM

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

A6. Molecular Weight/Extinction Coefficient

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

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 – MALEIMIDE 2nd Generation¹ (MO-L014, NanoTemper Technologies GmbH)
 1* Labeling Buffer Maleimide | 1* Dye RED-MALEIMIDE 2nd Generation (10 µg) | 1* A-Column | 1* B-Column

A9. Labeling Procedure

1. Prepare 50 µL of 25 µM VCB (VHL/Elongin B/Elongin C) by mixing 6.1 µL of 206 µM VCB with 43.9 µL of Labeling Buffer Maleimide.
2. Use the A-Column to perform a buffer exchange into Labeling Buffer Maleimide.
 - 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 Labeling Buffer Maleimide and centrifuge at **1500 × g** for **1 min** (3x).
 - e. Place 50 µL of the 25 µM VCB 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 50 µL of ~20 µM VCB (~80% recovery).
3. Add 22 µL of DMSO to Dye RED-MALEIMIDE 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 45 µL of Labeling Buffer Maleimide to obtain 50 µL of a 60 µM dye solution (3x protein concentration).
5. Mix VCB and dye in a 1:1 volume ratio (100 µL final volume, 5% final DMSO concentration).
6. Incubate for 45 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 100 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. Centrifuge the eluate at 15,000 rpm and 4°C for 20 minutes to remove aggregates. Then, carefully transfer the supernatant into a fresh tube and mix well by pipetting.
13. Prepare 8 µL aliquots of the labeled VCB (~2.5 µM) and immediately store at -80°C.

¹ While Cys162 of VHL is buried at the interface between VHL and EloC, Cys77 is solvent exposed and adjacent to the native substrate binding site. Cys89 of EloB is also solvent exposed.

A10. Labeling Efficiency

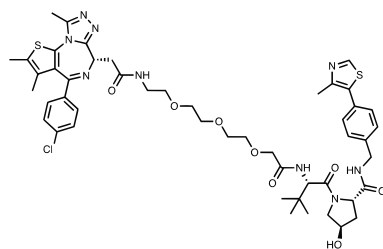
Measurement of protein concentration and degree of labeling (DOL) using a NanoDrop™:

nanotempertech.com/dol-calculator

Absorbance A_{280}	0.097	Protein concentration	2.51 μ M
Absorbance A_{650}	0.797	Degree-of-labeling (DOL)	1.63

B1. Ligand/Non-Fluorescent Binding Partner

MZ1



BRD2^{BD2}

uniprot.org/uniprot/P25440

B2. Molecule Class/Organism

PROTAC

Bromo-domain containing protein

B3. Sequence/Formula

C₄₉H₆₀ClN₉O₈S₂

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

B4. Purification Strategy/Source

opnme.com
[MZ1](#)

CreLux GmbH
 Construct ID: CRN1, Lot-ID: PC16670

B5. Stock Concentration/Stock Buffer

10 mg/mL | 10 mM
 DMSO

5.72 mg/mL | 362 μM
 25 mM HEPES pH 7.5, 100 mM NaCl, 5% glycerol, 1 mM DTT

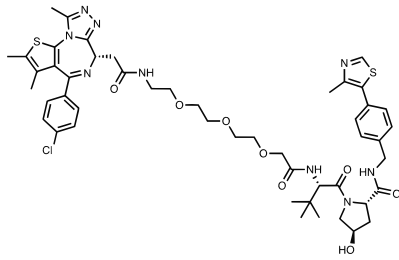
B6. Molecular Weight/Extinction Coefficient

1002.64 Da

15.8 kDa
 17,420 M⁻¹cm⁻¹ (ε₂₈₀)

B1. Ligand/Non-Fluorescent Binding Partner

MZ1



BRD2^{BD2}

uniprot.org/uniprot/P25440

B2. Molecule Class/Organism

PROTAC

Bromo-domain containing protein

B3. Sequence/Formula

C₄₉H₆₀ClN₉O₈S₂

MHHHHHSGV DLGTENLYFQ SMGKLSQLK HCNGILKELL SKKHAAYAWP FYKPVDASAL GLHDYHDIK HPMDLSTVKR
 KMENRDYRDA QEFAADVRLM FSNCYKYNPP DHDVVAMARK LQDVFEFRYA KMPD

B4. Purification Strategy/Source

opnme.com

MZI

Crelux GmbH

Construct ID: CRN1, Lot-ID: PC16670

B5. Stock Concentration/Stock Buffer

10 mg/mL | 10 mM

DMSO

5.72 mg/mL | 362 μM

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

B6. Molecular Weight/Extinction Coefficient

1002.64 Da

15.8 kDa

17,420 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 100 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 VCB ($\sim 2.5 \mu\text{M}$) with 496 μL of dilution buffer to obtain 500 μL of $\sim 20 \text{ nM}$ VCB.
6. Prepare a 4 μM BRD2^{BD2} solution by mixing 2 μL of 362 μM BRD2 with 179 μL of dilution buffer.

VCB – MZ1 (binary complex)

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

VCB – cis MZ1 (Control)

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

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

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

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

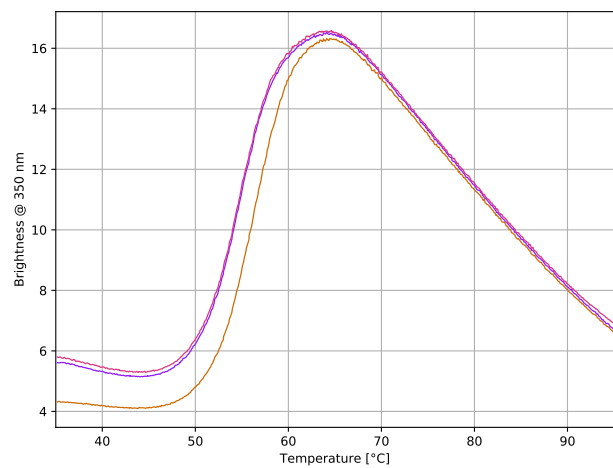
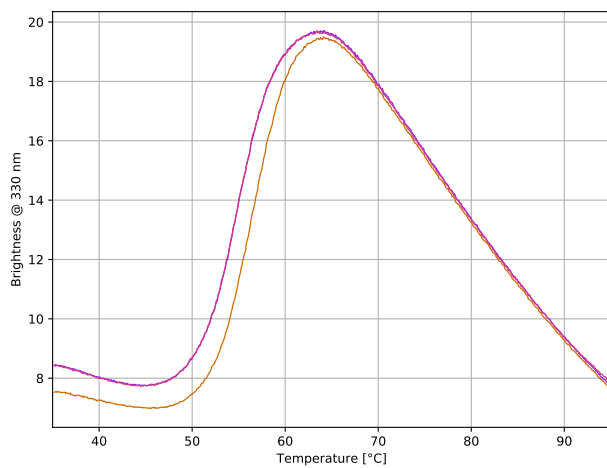
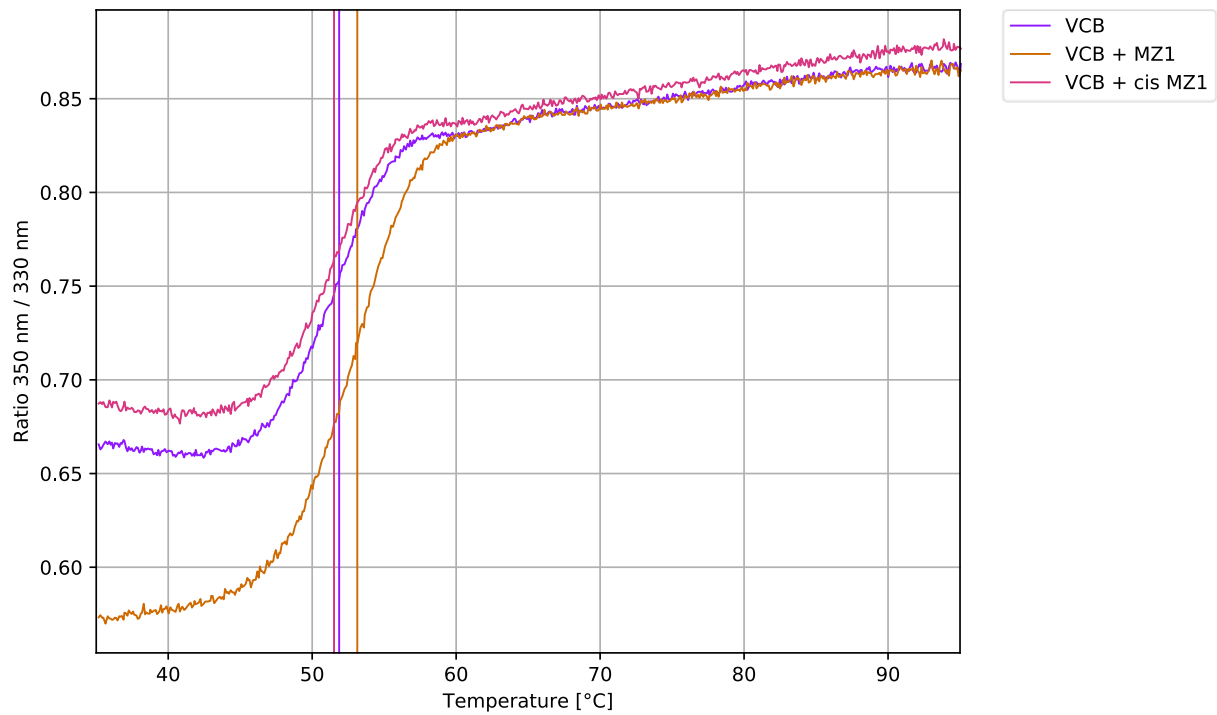
22. Mix 100 μL of 20 nM VCB with 2.5 μL of 4 μM BRD2^{BD2} and 97.5 μL of dilution buffer to obtain 200 μL of 10 nM VCB, 50 nM BRD2^{BD2}.
23. Take a fresh 0.5 mL tube and mix 160 μL of 10 nM VCB, 50 nM BRD2^{BD2} with 160 μL of dilution buffer containing 2% DMSO to obtain 320 μL of a 5 nM VCB, 25 nM BRD2^{BD2} solution.
24. Prepare a PCR-rack with 20 PCR tubes. Transfer 20 μL of the 5 nM VCB, 25 nM BRD2^{BD2} solution into tubes **2** to **16**. Then, mix 20 μL of 4 μM MZ1 with 20 μL of 10 nM VCB, 50 nM BRD2^{BD2} 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 VCB using Tycho NT.6:

nanotempertech.com/tycho

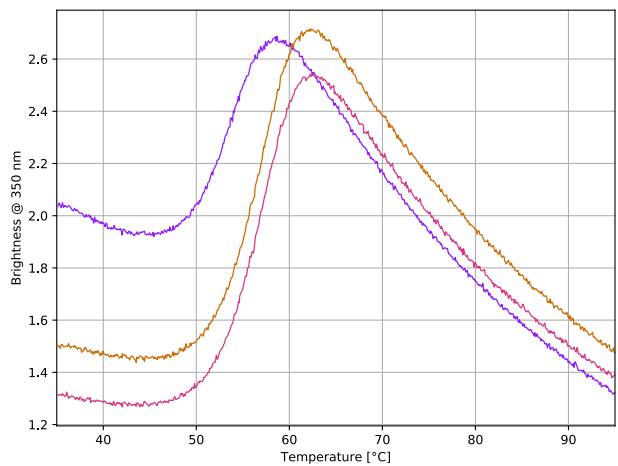
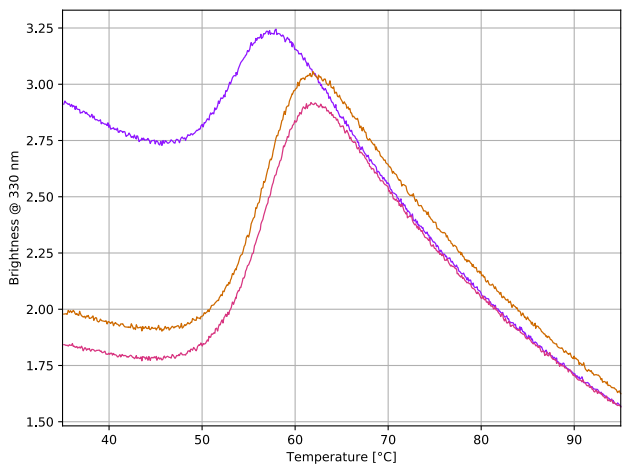
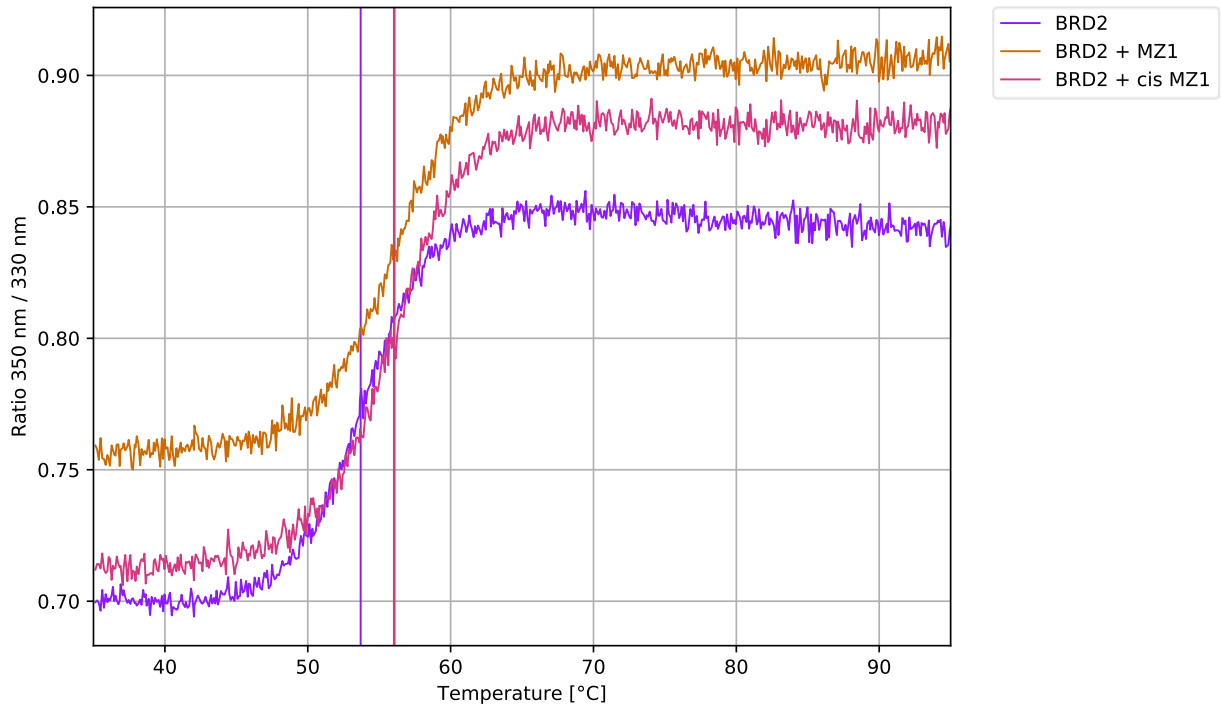
VCB	5 μ L of 4 μ M VCB + 5 μ L of dilution buffer containing 2% DMSO	$T_i = 51.9^\circ\text{C}$
VCB + MZ1	5 μ L of 4 μ M VCB + 5 μ L of 4 μ M MZ1	$T_i = 53.1^\circ\text{C}$
VCB + cis MZ1	5 μ L of 4 μ M VCB + 5 μ L of 4 μ M cis MZ1	$T_i = 51.5^\circ\text{C}$



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

nonotempertech.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



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)

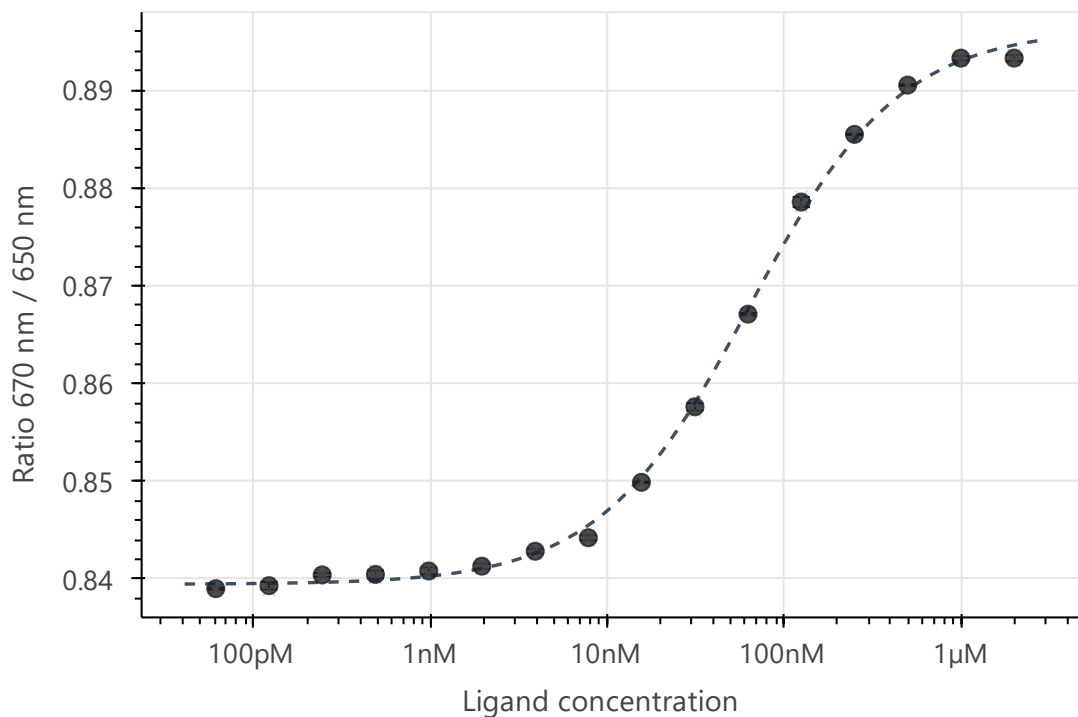
nanotempertech.com/monolith-mo-control-software

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

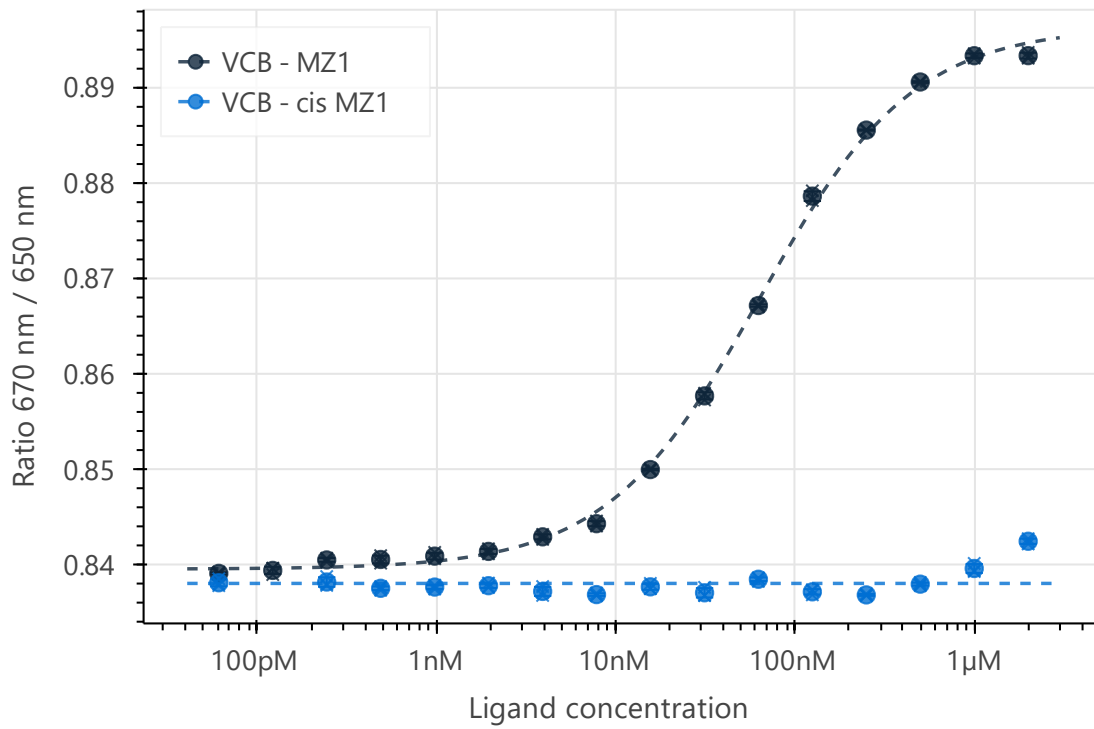
50 mM Tris-HCl, pH 7.8, 150 mM NaCl, 10 mM MgCl₂, 2 mM GSH, 0.01 % Pluronic® F-127, 1% DMSO
5 nM VCB | 2 μM – 61 pM MZ1 | 20°C | 100% Excitation Power

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

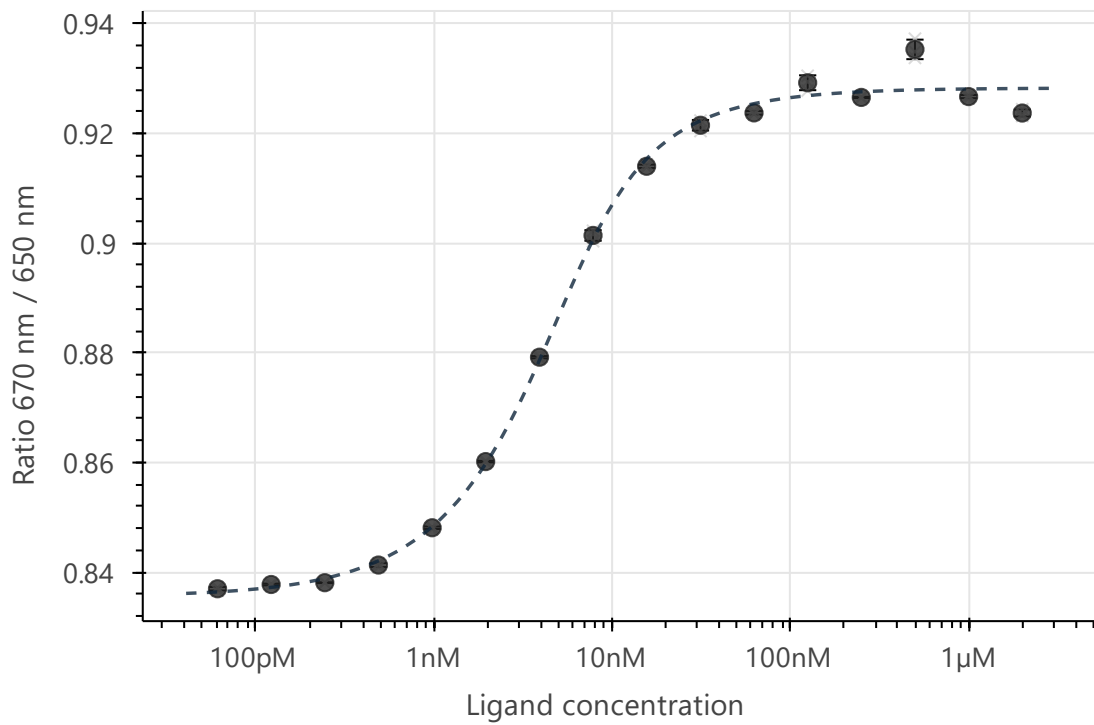
VCB – MZ1 (binary complex) | $K_d = 61.2 \pm 2.8$ nM (S/N = 76.1)



VCB – cis MZ1 (control) | $K_d > 10 \mu\text{M}$

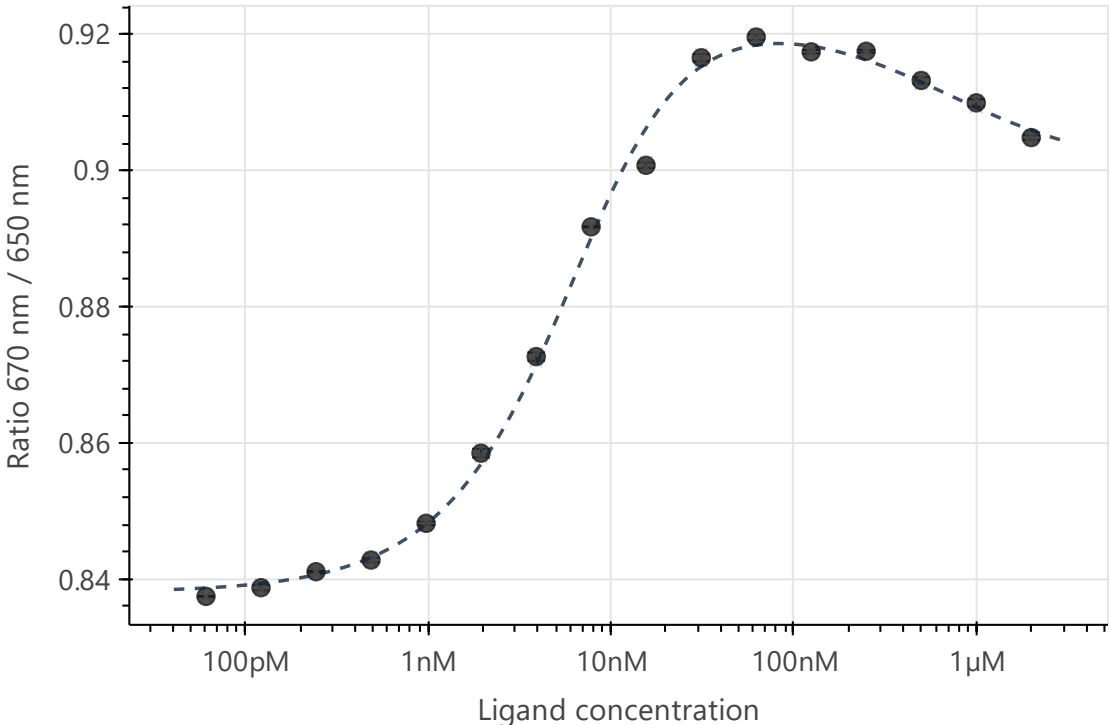


VCB & BRD2^{BD2} (500 nM) – MZ1 (ternary complex, saturation) | $K_d = 1.83 \pm 0.25 \text{ nM}$ (S/N = 37.5)

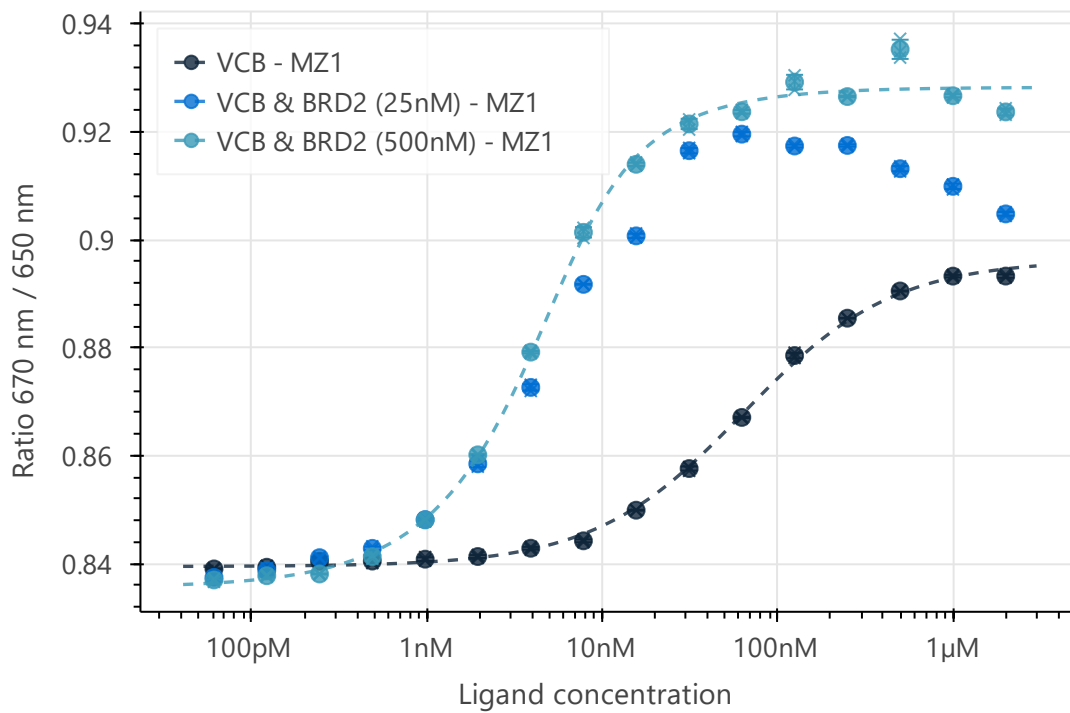


Cooperativity factor: $\alpha = \frac{61.2 \text{ nM}}{1.83 \text{ nM}} \approx 33.4$

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



Overlay



D5. Reference Results/Supporting Results

PROTAC		+ target	Monolith X		SPR		ITC	
			K _d (nM)	α ²	K _d (nM)	α	K _d (nM)	α
MZ1	binary	-	61.2	-	29	-	66	-
MZ1	ternary	BRD2 ^{BD2}	1.83	33.4	0.9	32	28	2.3

Isothermal Titration Calorimetry (ITC)

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

Surface Plasmon Resonance (SPR)

[Roy et al., ACS Chem. Biol. 14, 361–368 \(2019\)](#)

E. Contributors

Andreas Langer³

² Cooperativity factor

³ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com