

Monolith X Protocol MOX-P-117

HER2 – Herceptin ADCs

HER2 is a member of the human epidermal growth factor receptor (HER/EGFR/ERBB) family. Amplification or over-expression of this oncogene has been shown to play an important role in the development and progression of certain aggressive types of breast cancer. In recent years this protein has become an important biomarker and target of therapy for many breast cancer patients. Herceptin (Trastuzumab) is a monoclonal antibody used to treat breast cancer. Trastuzumab antibody-drug conjugates (ADCs) are a type of targeted cancer therapy that combines the specificity of monoclonal antibodies with the cytotoxic effects of chemotherapy drugs.

protein – protein | antibody | ADCs

A1. Target/Fluorescent Molecule

ErbB-2 (HER2)

uniprot.org/uniprot/P04626

A2. Molecule Class/Organism

Receptor tyrosine-protein kinase

Homo sapiens (Human)

A3. Sequence/Formula

TQVCTGDMK LRLPASPETH LDMLRHLYQG CQVVQGNLEL TYLPTNASLS FLQDIQEVQG YVLIAHNQVR QVPLQRLRIV
 RGTQLFEDNY ALAVLDNGDP LNNTTPVTGA SPGGLRELQL RSLTEILKGG VLIQRNPQLC YQDTILWKDI FHKNNQLALT
 LIDTNRSRAC HPCSPMCKGS RCWGESSEDC QSLTRTVGAG GCARCKGPLP TDCCHEQCAA GCTGPKHSDC LACLHFNHSG
 ICELHCPALV TYNTDTFESM PNPEGRYTFG ASCVTACPYN YLSTDVGSCT LVCPLHNQEV TAEDGTQRCE KCSKPCARVC
 YGLGMEHLRE VRAVTSANIQ EFAGCKKIFG SLAFLPESFD GDPASNTAPL QPEQLQVFET LEEITGYLYI SAWPDSLPLD
 SVFQNLQVIR GRILHNGAYS LTLQGLGISW LGLRSLRELG SGLALIHNT HLCFVHTVPW DQLFRNPHQA LLHTANRPED
 ECVGEGlach QLCARGHCWG PGPTQCVNCS QFLRGQECVE ECRVLQGLPR EYVNARHCLP CHPECQPQNG SVTCFGPEAD
 QCVACAHYKD PPFVARCPS GVKPDLSEMP IWKFPDEEGA CQPCPINCTH SCVDLDDKGC PAEQRASPLT

A4. Purification Strategy/Source

Recombinant, His-tagged (C-terminus)

Sino Biological

[10004-H08H](#)

A5. Stock Concentration/Stock Buffer

1 mg

Lyophilized (PBS, pH 7.4)

A6. Molecular Weight/Extinction Coefficient

71 kDa

65,000 M⁻¹cm⁻¹ (ε₂₈₀)

A7. Dilution Buffer

Phosphate-buffered saline (PBS), pH 7.4, 0.005% TWEEN® 20

A8. Labeling Strategy

Monolith Protein Labeling Kit RED – NHS 2nd Generation (MO-L011, NanoTemper Technologies GmbH)
 1* Dye RED-NHS 2nd Generation (10 µg) | 1* B-Column

A9. Labeling Procedure

1. Resuspend 1 mg HER2 in 500 µL ddH₂O to obtain a 2 mg/mL (28.2 µM) solution.
2. Prepare 10 µL aliquots and store at -80°C.
3. Mix 7.1 µL of 28.2 µM HER2 with 12.9 µL of dilution buffer to obtain 20 µL of 10 µM HER2.
4. Add 12.5 µL of DMSO to 10 µg RED-NHS 2nd Generation dye to obtain a ~1.2 mM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
5. Mix 0.5 µL of the 1.2 mM dye solution with 19.5 µL of dilution buffer to obtain 20 µL of a 60 µM dye solution (3x protein concentration).
6. Mix HER2 and dye in a 1:1 volume ratio (40 µL final volume, ~1.25% final DMSO concentration).
7. Incubate for 30 minutes at room temperature in the dark.
8. 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.
9. 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.
10. Add 40 µL of the labeling reaction from step 6 to the center of the column and let sample enter the resin bed completely.
11. Add 560 µL of dilution buffer after the sample has entered and discard the flow through.
12. Place column in a new collection tube, add 500 µL of dilution buffer and collect the eluate.
13. Prepare 20 µL aliquots of the labeled HER2 (~0.4 µM) and store at -80°C.

A10. Labeling Efficiency

N/A

B1. Ligand/Non-Fluorescent Binding Partner

Herceptin (Trastuzumab)

Trastuzumab emtansine (Kadcyla)

Trastuzumab deruxtecan (Enhertu)

Disitamab vedotin (RC48)

B2. Molecule Class/Organism

Monoclonal antibody, IgG1

B3. Sequence/Formula

Heavy chain

EVQLVESGGG LVQPGGSLRL SCAASGFNIK DTYIHVVRQA PGKGLEWVAR IYPTNGYTRY ADSVKGRFTI SADTSKNTAY
 LQMNSLRAED TAVYYCSRWG GDGFYAMDYW GQGTLVTVSS ASTKGPSVFP LAPSSKSTSG GTAALGCLVK DYFPEPVTVS
 WNSGALTSKV HTFPAVLQSS GLYSLSSVVT VPSSSLGTQT YICNVNHKPS NTKVDKKVEP KSCDKTHTCP PCPAPELLGG
 PSVFLFPPPK KDTLMISRTP EVTCVVDVS HEDPEVKFNW YVDGVEVHNA KTKPREEQYN STYRVVSVLT VLHQDWLNGK
 EYKCKVSNKA LPAPIEKTIS KAKGQPREPQ VYTLPPSREE MTKNQVSLTC LVKGFYPSDI AVEWESNGQP ENNYKTTTPV
 LSDSGSFFLY SKLTVDKSRW QQGNVFSCSV MHEALHNHYT QKSLSLSPG

Light chain

DIQMTQSPSS LSASVGRVT ITCRASQDVN TAVAWYQQKPK GKAPKLLIYS ASFLYSGVPS RFGSRSRGTD FTLTISSLQP
 EDFATYYCQQ HYTTPPTFGQ GTKVEIKRTV AAPSVEIFPP SDEQLKSGTA SVVCLLNNFY PREAKVQWKV DNALQSGNSQ
 ESVTEQDSKD STYLSLSTLT LSKADYEKHK VYACEVTHQG LSSPVTKSFN RGEK

B4. Purification Strategy/Source

MedChemExpress

[HY-P9921](#)

[HY-138298A](#)

[HY-P9985](#)

B5. Stock Concentration/Stock Buffer

5 mg/mL | 34.3 μ M

5 mg/mL | 34.3 μ M

5 mg/mL | 34.3 μ M

B6. Molecular Weight/Extinction Coefficient

145.5 kDa

225,000 $M^{-1}cm^{-1}$ (ϵ_{280})

B7. Serial Dilution Preparation

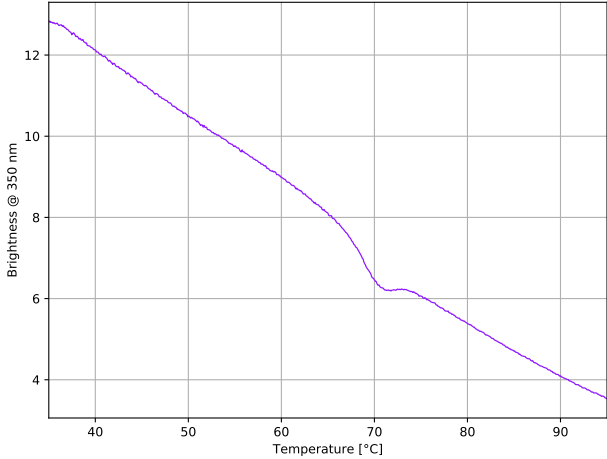
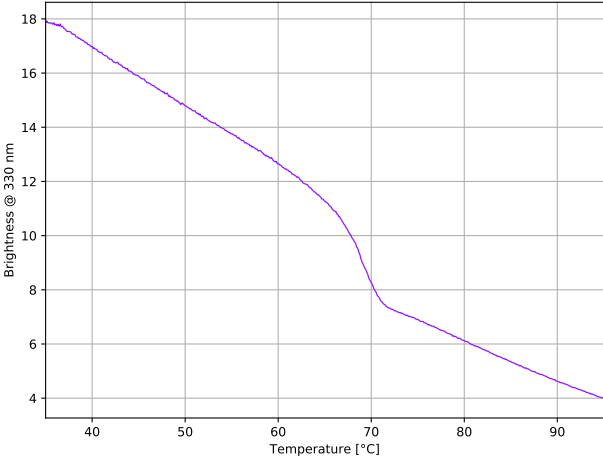
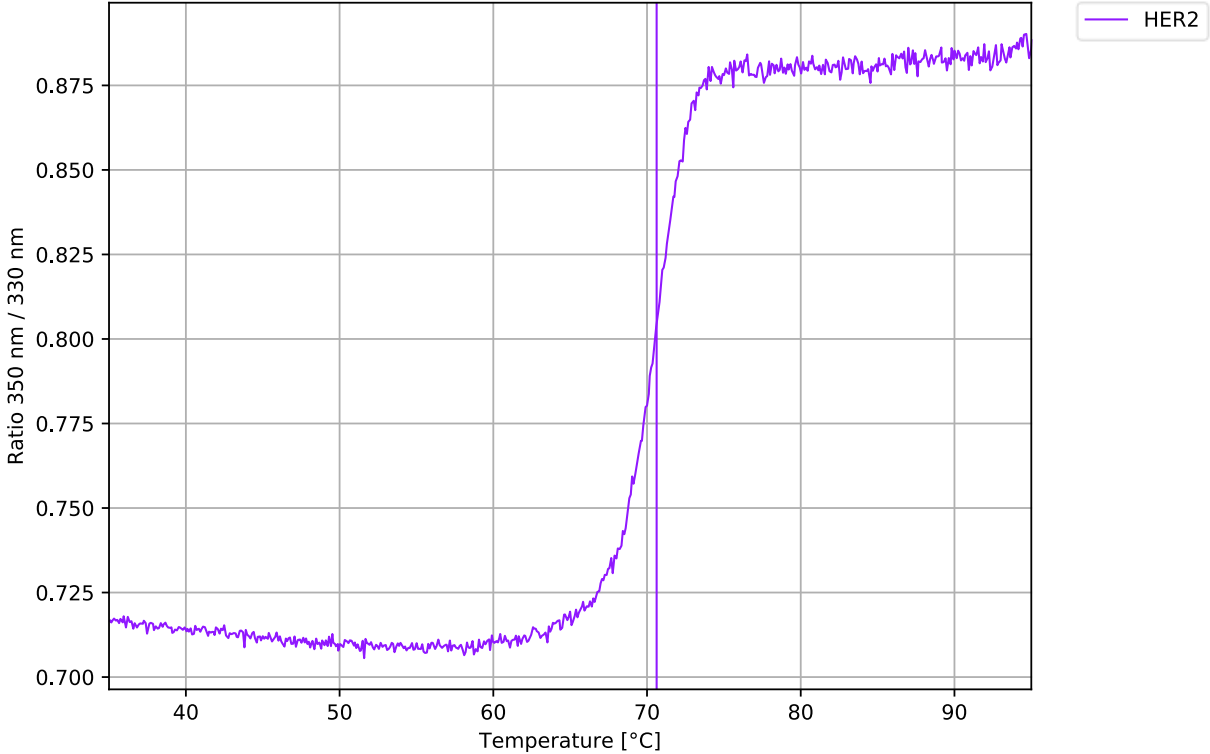
1. Mix 1 μ L of 34.3 μ M ADC with 170.5 μ L of dilution buffer to obtain a 200 nM ADC solution.
2. Mix 8 μ L of 400 nM labeled HER2 with 392 μ L of dilution buffer to obtain 400 μ L of an 8 nM HER2 solution.
3. Take a fresh PCR tube and mix 80 μ L of dilution buffer with 80 μ L of 8 nM HER2 to obtain 160 μ L of 4 nM HER2.
4. Prepare a PCR-rack with 16 PCR tubes. Transfer 10 μ L of 4 nM HER2 into tubes **2** to **16**. Then, mix 10 μ L of 8 nM HER2 with 10 μ L of 200 nM ADC in tube **1**.
5. Prepare a 1:1 serial dilution by transferring 10 μ L from tube to tube. Mix carefully by pipetting up and down.
6. Incubate for at least 30 minutes at room temperature in the dark before loading capillaries.

C. Tycho

Validation of structural integrity of HER2 using Tycho NT.6:

nanotempertech.com/tycho

HER2	10 µL of 1 µM HER2	T _i = 70.6°C
------	--------------------	-------------------------



D1. Monolith System/Capillaries

Monolith X (NanoTemper Technologies GmbH)

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

D2. Monolith Software

MO.Control v2.6.3 (NanoTemper Technologies GmbH)

nanotempertech.com/monolith-mo-control-software

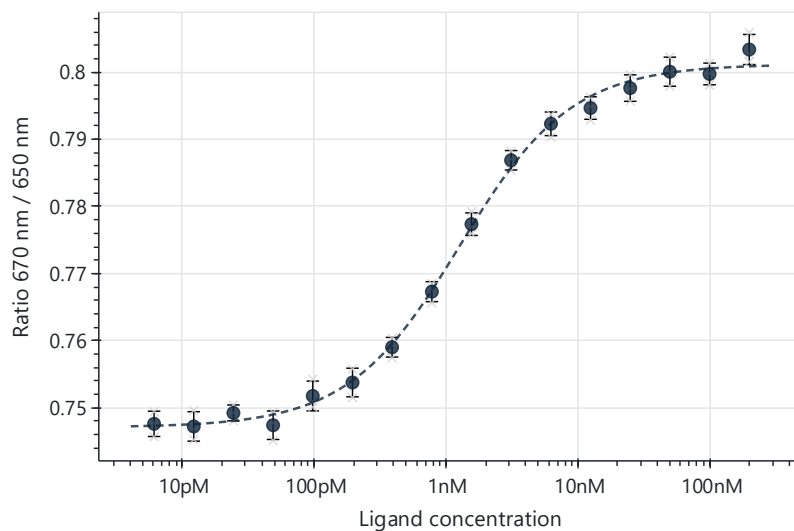
D3. Monolith Experiment (Assay Buffer/Concentrations/Temperature/Excitation Power)

Phosphate-buffered saline (PBS), pH 7.4, 0.005% TWEEN® 20

4 nM HER2 | 100 nM¹ – 3.0 pM Trastuzumab ADC | 25°C | 100% excitation power

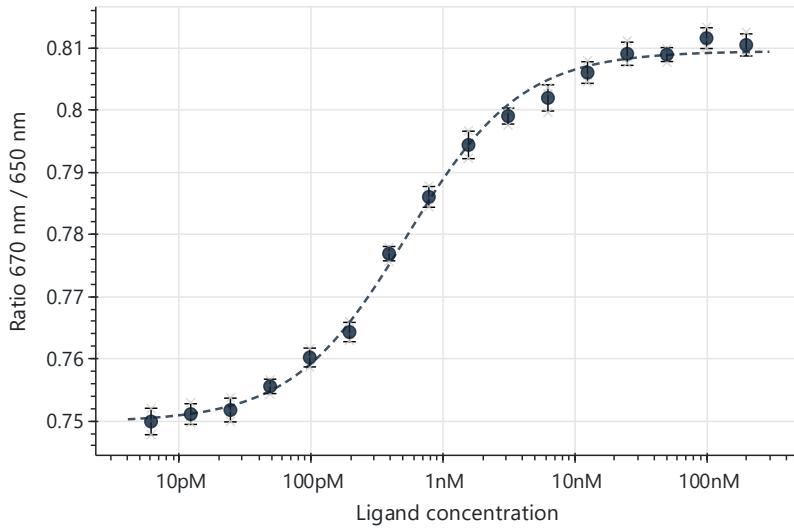
D4. Monolith Results (Dose Response)

Trastuzumab | $K_d = 1.42 \pm 0.1$ nM (S/N = 49.1)

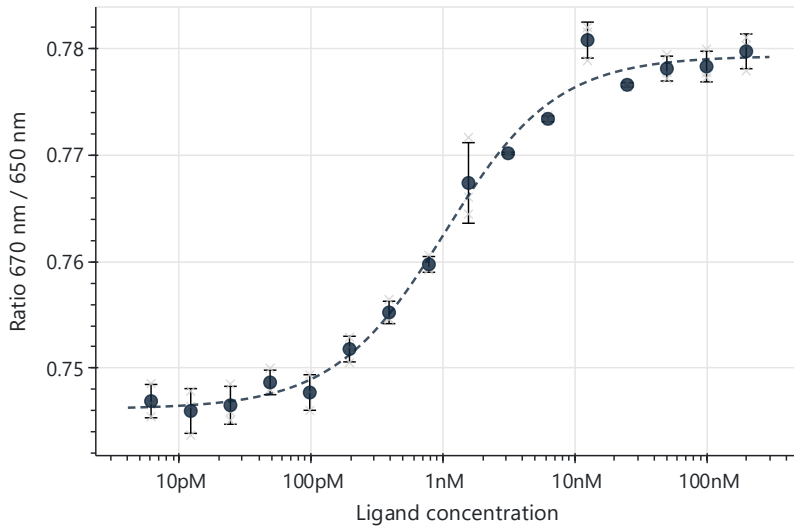


¹ 100 nM Trastuzumab corresponds to a ligand concentration of 200 nM (two antigen binding sites per antibody)

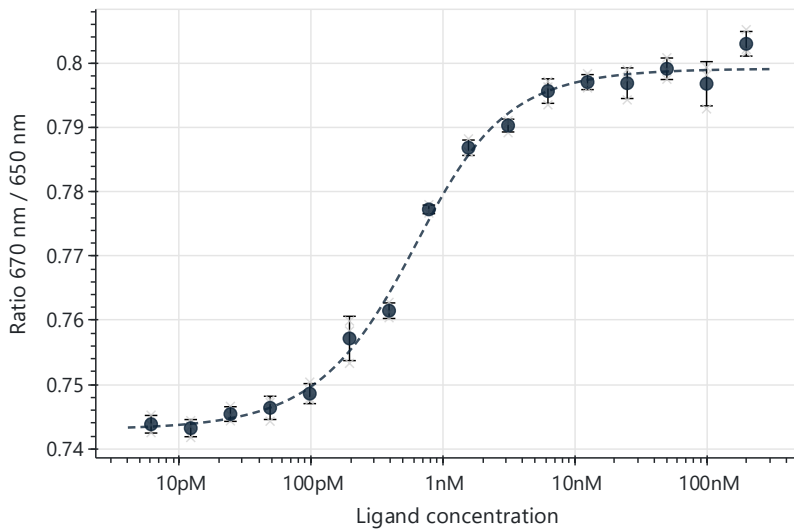
Trastuzumab emtansine | $K_d = 646 \pm 49 \text{ pM}$ (S/N = 42.6)



Disitamab vedotin | $K_d = 1.12 \pm 0.15 \text{ nM}$ (S/N = 24.6)



Trastuzumab deruxtecan | $K_d = 441 \pm 55 \text{ pM}$ (S/N = 33.0)



D5. Reference Results/Supporting Results

$K_d = 1.3 \text{ nM}$
($k_a = 1.2 \times 10^5 \text{ M}^{-1}\text{s}^{-1}$, $k_d = 1.6 \times 10^{-4} \text{ s}^{-1}$)

Surface Plasmon Resonance (SPR)
[Epa et al., PLOS One, 8\(3\), 2013](#)

$K_d = 0.4 \text{ nM}$
($k_a = 2.8 \times 10^5 \text{ M}^{-1}\text{s}^{-1}$, $k_d = 1.2 \times 10^{-4} \text{ s}^{-1}$)

Surface Plasmon Resonance (SPR)
[Pedersen et al., Mol Cancer Ther, 14\(3\) 2015](#)

E. Contributors

Andreas Langer²

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