

Monolith Protocol M0-P-011

Protein A – Herceptin

Protein A is a cell surface receptor found in the cell wall of the bacteria *Staphylococcus aureus*. It's frequent use in biochemical research derives from its ability to bind the Fc region of most immunoglobulins. Herceptin (Trastuzumab) is a monoclonal antibody used to treat breast cancer.

protein – protein interaction | antibody

A1. Target/Fluorescent Molecule

Protein A

uniprot.org/uniprot/P38507

A2. Molecule Class/Organism

Immunoglobulin-binding protein

Staphylococcus aureus

A3. Sequence/Formula

AQHDEAQQNA FYQVLNMPNL NADQRNGFIQ SLKDDPSQSA NVLGEAQKLN DSQAPKADAQ QNKFNKDQQS AFYEILNMPN
LNEEQRNGFI QSLKDDPSQS TNVLGEAKKL NESQAPKADN NFNKEQQNAF YEILNMPNLN EEQRNGFIQS LKDDPSQSAN
LLAEAKKLNE SQAPKADNKF NKEQQNAFYE ILHLPNLNEE QRNGFIQSLK DDPSQSANLL AEAKKLNDAAQ APKADNKFNK
EQQNAFYEIL HLPNLTEEQR NGFIQSLKDD PSVSKEILAE AKKLNDAAQAP KEEDNNKPGK EDGNKPGKED GNKPGKEDNK
KPGKEDGNKP GKEDNKKPGK EDGNKPGKED GNKPGKEDGN KPGKEDGNKP GKEDGNGVHV VKPGDTVNDI AKANGTTADK
IAADNKLADK NMIKPGQELV VDKKQPANHA DANKAQLPE T

A4. Purification Strategy/Source

Recombinant

anitkoerper-online

[ABIN1067599](https://uniprot.org/uniprot/P38507)

A5. Stock Concentration/Stock Buffer

5 mg/mL | 122 µM

H₂O

A6. Molecular Weight/Extinction Coefficient

42 kDa

6,930 M⁻¹cm⁻¹ (ε₂₈₀)

A7. Dilution Buffer

Phosphate-buffered saline (PBS, pH 7.4), 0.05% TWEEN® 20

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. Add 45 µL of Labeling Buffer NHS to 5 µL of 122 µM Protein A to obtain 50 µL of a 12 µM solution.
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 3 µL of the 600 µM dye solution with 47 µL of Labeling Buffer NHS to obtain 50 µL of a 36 µM dye solution (3x protein concentration).
4. Mix Protein A and dye in a 1:1 volume ratio (100 µL final volume, 3% 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 3 to the center of the column and let sample enter the bed completely.
9. Add 600 µL of dilution buffer after the sample has entered and discard the flow through.
10. Place column in a new collection tube, add 300 µL of dilution buffer and collect the eluate.
11. Keep the labeled protein A (~2 µM) on ice in the dark.

A10. Labeling Efficiency

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

Absorbance A ₂₀₅	3.70	Protein concentration	2.77 µM
Absorbance A ₆₅₀	0.499	Degree-of-labeling (DOL)	0.92

B1. Ligand/Non-Fluorescent Binding Partner

Herceptin (Trastuzumab)

B2. Molecule Class/Organism

Monoclonal antibody

B3. Sequence/Formula

Heavy chain

EVQLVESGGG LVQPGGSLRL SCAASGFNIK DTYIH^WVRQA PGKGLE^WVAR IYPTNGYTRY ADSVKGRFTI SADTSKNTAY
LQMNSLRAED TAVYYCSR^WG GDGFYAMDY^W GQGT^LLVTVSS ASTKGPSVFP LAPSSKSTSG GTAALGCLVK DYFPEPVTVS
^WNSGALTSGV HTFPAVLQSS GLYSLSSVVT VPSSSLGTQT YICNVNHKPS NTKVDKKVEP KSCDKTHTCP PCPAPELLGG
PSVFLFPPKP KDTLMISRTPEVTCVVDVS HEDPEVKFN^W YVDGVEVHNA KTKPREEQYN STYRVVSVLT VLHQD^WLNGK
EYKCKVSNKA LPAPIEKTIS KAKGQPREPQ VYTLPPSREE MTKNQVSLTC LVKGFYPSDI AVE^WESNGQP ENNYKTTPPV
LDSGDSFFLY SKLTVDKSR^W QQGNVFSCSV MHEALHNHYT QKSLSLSPG

Light chain

DIQMTQSPSS LSASVGDRVT ITCRASQDVN TAVA^WYQQKP GKAPKLLIYS ASFLYSGVPS RFGSGRSGTD FTLTISSLQP
EDFATYYCQQ HYTTPPTFGQ GTKVEIKRTV AAPS^VFI^FPP SDEQLKSGTA SVVCLLN^NFY PREAKVQ^WKV DNALQSGNSQ
ESVTEQDSKD STYLSLSTLT LSKADYEKHK VYACEVTHQG LSSPVT^KSFN RGEC

B4. Purification Strategy/Source

N/A

B5. Stock Concentration/Stock Buffer

120 mg/mL | 825 μ M

B6. Molecular Weight/Extinction Coefficient

145.5 kDa

B7. Serial Dilution Preparation

1. Add 145 μ L of dilution buffer to 20 μ L of 825 μ M Herceptin to obtain 165 μ L of a 100 μ M solution.
2. Mix 2 μ L of 100 μ M Herceptin with 198 μ L of dilution buffer to obtain 200 μ L of a 1 μ M Herceptin solution.
3. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μ L of the 1 μ M Herceptin solution into tube **1**. Then, transfer 10 μ L of dilution buffer into tubes **2** to **16**.
4. Prepare a 1:1 serial dilution by transferring 10 μ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10 μ L from tube **16** to get an equal volume of 10 μ L for all samples.
5. Mix 2 μ L of labeled protein A with 38 μ L of dilution buffer to obtain 40 μ L of ~100 nM protein A.
6. Mix 4 μ L of ~100 nM labeled protein A with 196 μ L of dilution buffer to obtain 200 μ L of ~2 nM protein A.
7. Add 10 μ L of labeled protein A (~2 nM) to each tube from **16** to **1** and mix by pipetting.
8. Incubate for 5 minutes¹ at room temperature in the dark before loading capillaries.

¹ Longer incubation can lead to aggregation due to the multiple IgG-binding sites of Protein A.

BB. SD-Test

1. When checking for specificity of ligand-induced changes in initial fluorescence with the SD-Test, do **not** centrifuge² the remainder of tubes **1** to **3** and **14** to **16**.
2. Prepare the SD-mix: Dilute 400 μL of 10% SDS and 40 μL of 1 M DTT in 560 μL water to obtain a solution containing 4% SDS and 40 mM DTT.
3. Transfer 7 μL of the SD-mix to six PCR tubes.
4. Add 7 μL from the remainder of tubes **1** to **3** and **14** to **16** to the tubes containing 7 μL SD-mix. Mix well by pipetting.
5. Place samples on a heat block set to 95°C for 5 minutes to denature the protein, then allow to cool at 25°C for 10 minutes before loading into capillaries.

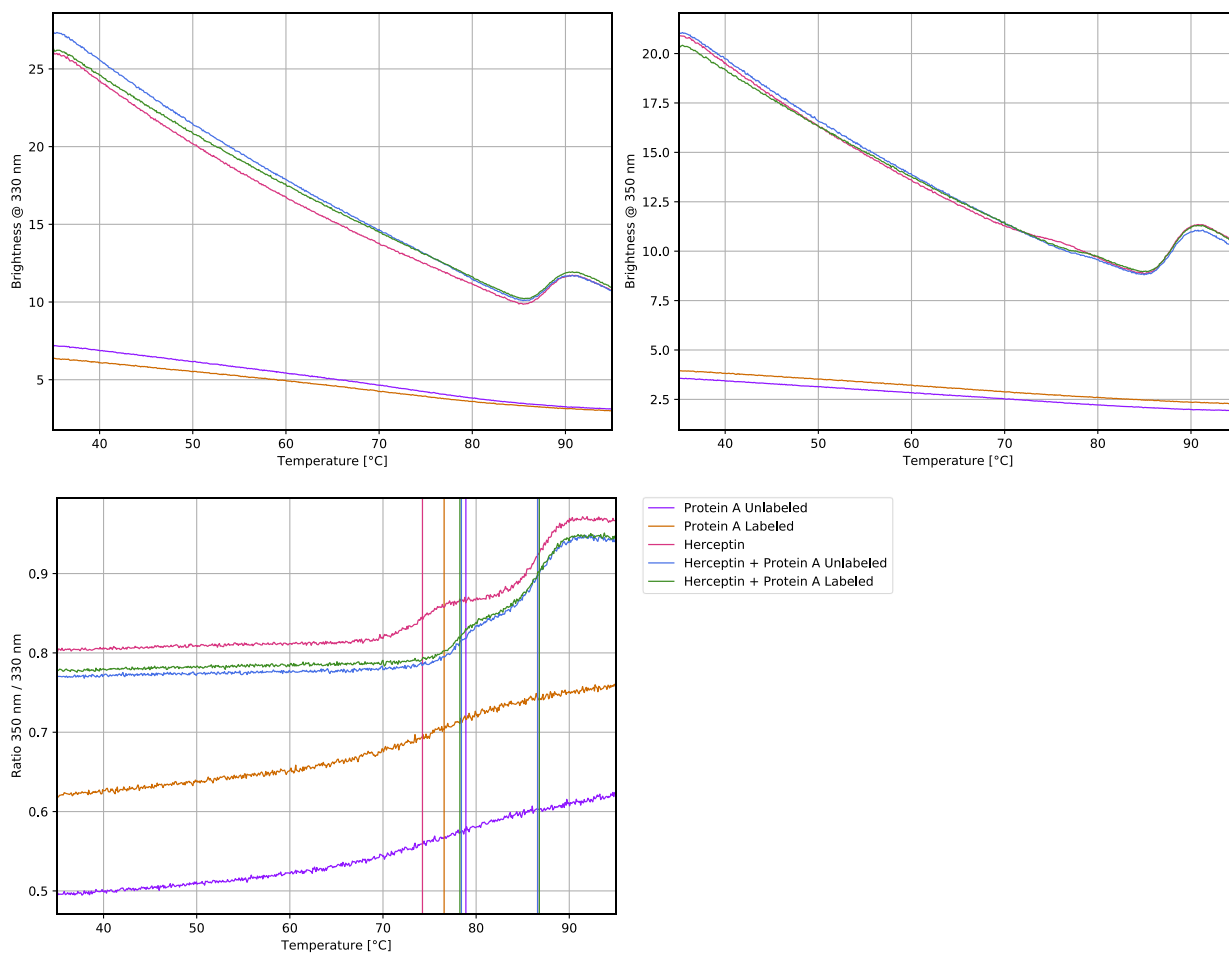
² Centrifugation will cause larger complexes of Herceptin and Protein A to precipitate out of the aqueous solution.

C. Applied Quality Checks

Validation of structural integrity and functionality of labeled protein A and Herceptin using Tycho NT.6:

nanotempertech.com/tycho

Protein A (UL)	1 μ L of 122 μ M protein A + 40 μ L of dilution buffer	$T_i = 78.9^\circ\text{C}$
Protein A (L)	10 μ L of B-Column eluate ($\sim 2.9 \mu\text{M}$)	$T_i = 76.7^\circ\text{C}$
Herceptin	5 μ L of 1 μ M Herceptin + 5 μ L of dilution buffer	$T_{i,1} = 74.2^\circ\text{C} \mid T_{i,2} = 86.6^\circ\text{C}$
Herceptin + protein A (UL)	5 μ L of 1 μ M Herceptin + 5 μ L of 2.9 μ M protein A	$T_{i,1} = 78.4^\circ\text{C} \mid T_{i,2} = 86.6^\circ\text{C}$
Herceptin + protein A (L)	5 μ L of 1 μ M Herceptin + 5 μ L of B-Column eluate ($\sim 2.9 \mu\text{M}$)	$T_{i,1} = 78.3^\circ\text{C} \mid T_{i,2} = 86.8^\circ\text{C}$



D1. MST System/Capillaries

Monolith NT.115^{Pico} Red (NanoTemper Technologies GmbH)

Premium Capillaries Monolith NT.115 (MO-K025, NanoTemper Technologies GmbH)

D2. MST Software

MO.Control v1.6 (NanoTemper Technologies GmbH)

nanotempertech.com/monolith-mo-control-software

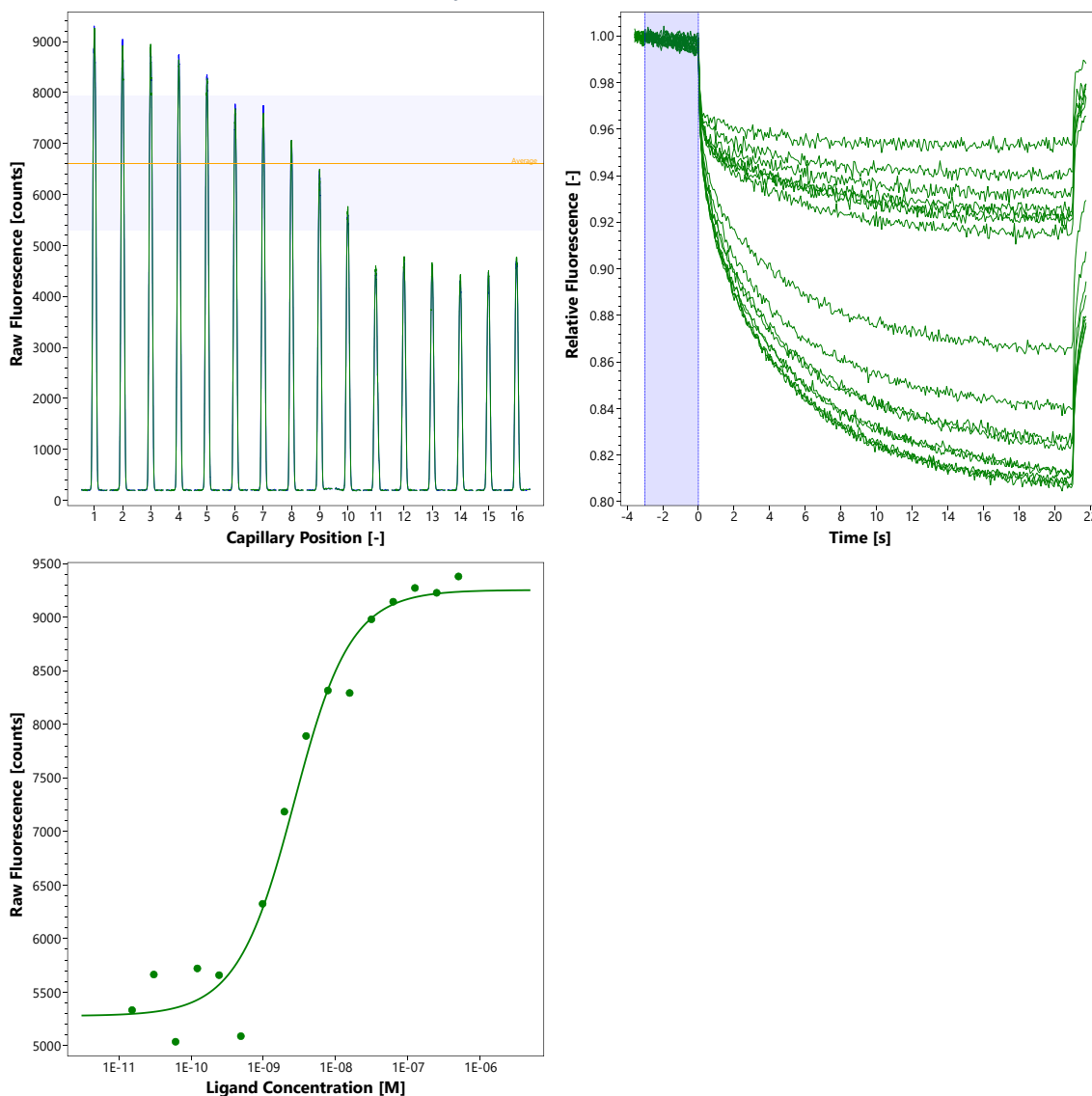
D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

Phosphate-buffered saline (PBS, pH 7.4), 0.05% TWEEN® 20

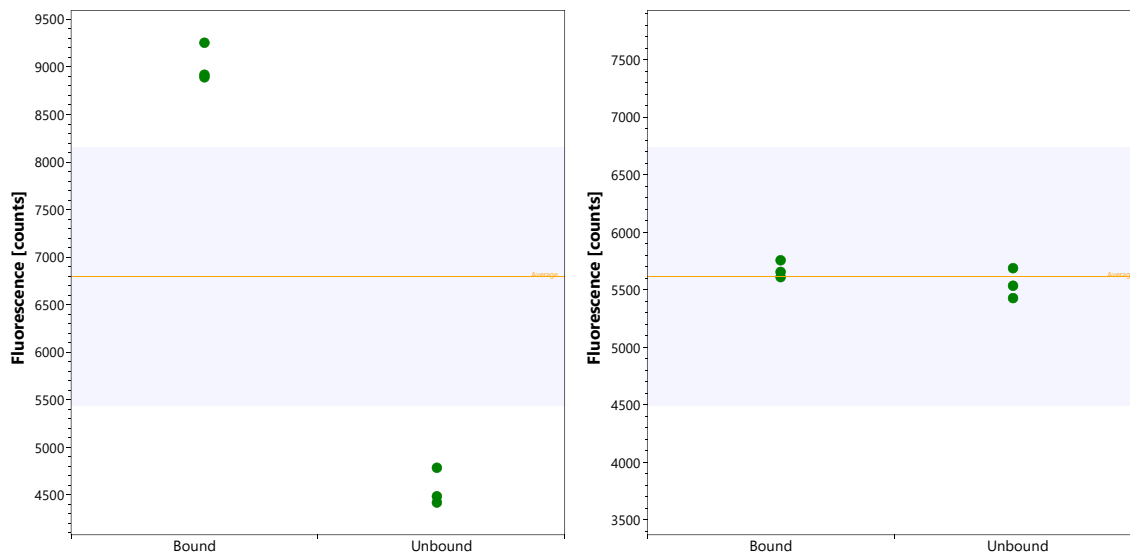
1 nM protein A | 500 nM – 15 pM Herceptin | 25°C | medium MST power | 20% excitation power

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

$K_d = 2.15$ nM Initial Fluorescence Analysis



SD-Test³:



D5. Reference Results/Supporting Results

$K_d = 7.1 \text{ nM}$ Surface Plasmon Resonance (SPR)
[Svensson et al., Eur J Biochem 258 \(2001\) 890-896](#)

E. Contributors

Andreas Langer⁴

³ Due to the ligand-dependent fluorescence changes, an SD-test was performed to confirm binding-dependent change. As the changes in initial fluorescence were concluded to be binding-specific, the initial fluorescence data was used for binding curve fit and K_d determination.

⁴ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com