

Monolith X Protocol MOX-P-115

DNA Hybridization - Thermodynamics

The stability of DNA duplexes depends on a fine balance of hydrogen bonds and base-stacking interactions. Factors affecting stability are ionic strength and temperature. Monitoring the shift in the dissociation constant (K_d) of the hybridization interaction over a range of temperatures can be used to calculate the free enthalpy (ΔH) and entropy (ΔS) for the hybridization reaction.

nucleic acid – nucleic acid | DNA | hybridization | thermodynamics | van't Hoff analysis | enthalpy | entropy

A1. Target/Fluorescent Molecule

Cy5-GB1

A2. Molecule Class/Organism

DNA

A3. Sequence/Formula

5' Cy5 GGA CTT CAG G 3'

A4. Purification Strategy/Source

metabion international AG

A5. Stock Concentration/Stock Buffer

0.36 mg/ml | 100 μ M
ddH₂O

A6. Molecular Weight/Extinction Coefficient

3602 Da
98,000 M⁻¹cm⁻¹ (ϵ_{260})

A7. Dilution Buffer

40 mM sodium phosphate buffer, pH 7.2, 200 mM NaCl, 0.1% TWEEN® 20¹

A8. Labeling Strategy

Cy5

¹ 2X assay buffer

A9. Labeling Procedure

1. Prepare the dilution buffer by dissolving 236 mg of Na_2HPO_4 (MW: 177.99 g/mol), 105 mg of NaH_2PO_4 (MW: 156.01 g/mol) and 584 mg of NaCl (MW: 58.44 g/mol) in 49.5 mL of ddH₂O. Then, add 500 μL of 10% TWEEN® 20.
2. Dilute 1.5 μL of 100 μM Cy5-GB1 in 15 mL of dilution buffer to obtain a 10 nM Cy5-GB1 solution.
3. Prepare 180 μL aliquots and store at -80°C .

A10. Labeling Efficiency

HPLC-purified, 100% labeled oligonucleotide

B1. Ligand/Non-Fluorescent Binding Partner

aGB1

B2. Molecule Class/Organism

DNA

B3. Sequence/Formula

5' CCT GAA GTC C 3'

B4. Purification Strategy/Source

metabion international AG

B5. Stock Concentration/Stock Buffer

0.60 mg/ml | 200 μ M
ddH₂O

B6. Molecular Weight/Extinction Coefficient

2988 Da
91,200 M⁻¹cm⁻¹ (ϵ_{260})

B7. Serial Dilution Preparation

1. Prepare 180 μ L of 10 nM Cy5-GB1.
2. Prepare a PCR-rack with 16 PCR tubes. Mix 2 μ L of 200 μ M aGB1 with 18 μ L of ddH₂O in tube **1**. Then, transfer 10 μ L of ddH₂O into tubes **2** to **16**.
3. 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.
4. Add 10 μ L of 10 nM Cy5-GB1 to each tube from **16** to **1** and mix by pipetting.
5. Load capillaries².
6. Set the Monolith capillary tray temperature to the desired value and wait at least 5 minutes before starting the measurement to ensure sufficient equilibration of the system.

² When performing multiple measurements at different temperatures, seal capillaries using the Capillary Sealing Paste Prometheus Series (PR-P001, NanoTemper Technologies GmbH).

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)

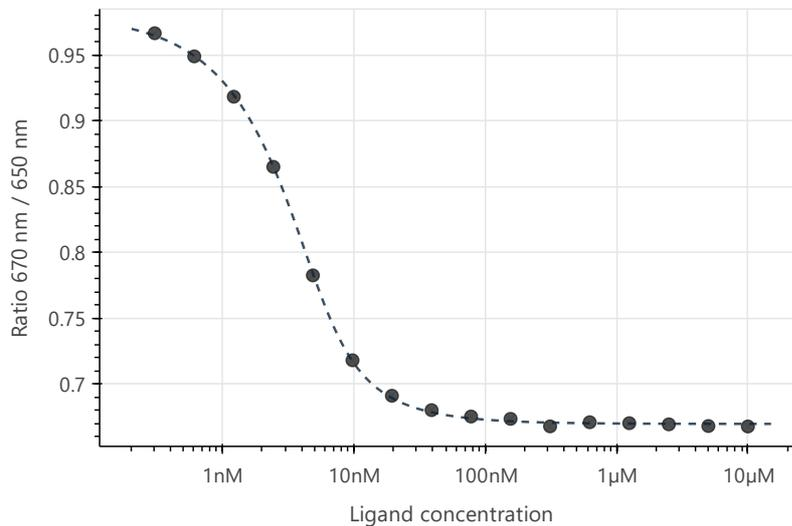
nanotempertech.com/monolith-mo-control-software

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

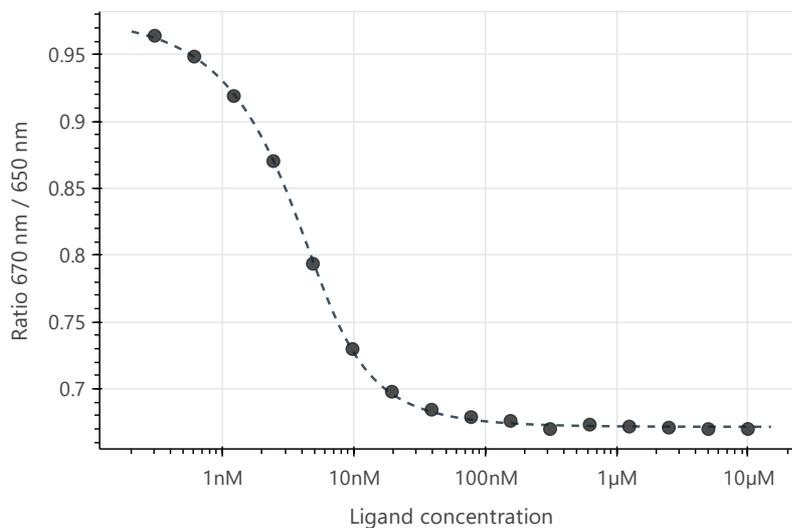
20 mM sodium phosphate buffer, pH 7.2, 100 mM NaCl, 0.05% TWEEN® 20
 5 nM Cy5-GB1 | 10 μ M – 305 pM aGB1 | 20°C – 37°C | 100% excitation power

D4. Monolith Results (Dose Response)

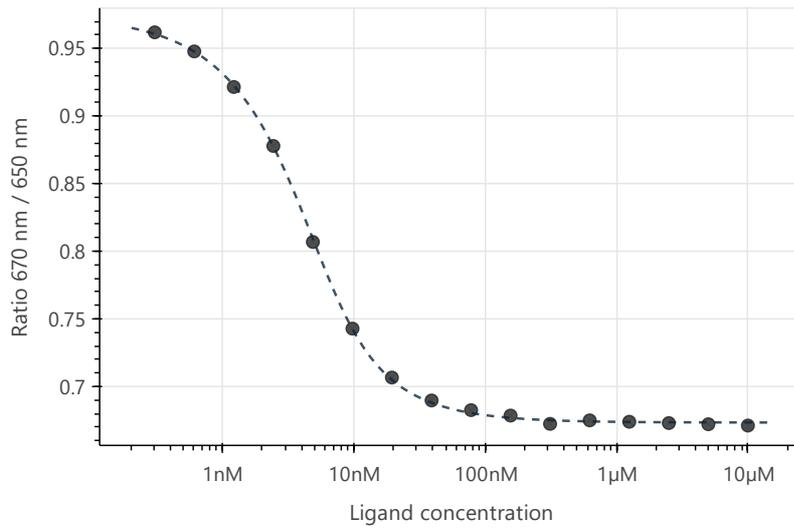
20°C | $K_d = 1.00 \pm 0.03$ nM (S/N = 203.9)



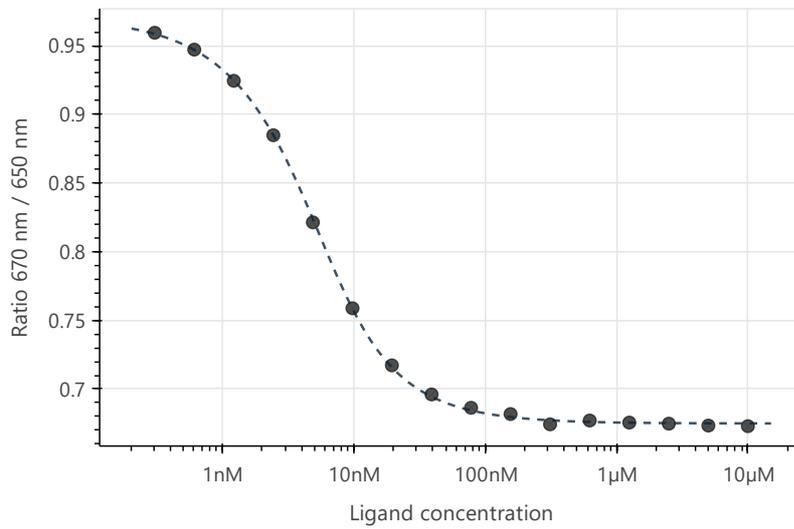
21°C | $K_d = 1.31$ nM \pm 0.04 (S/N = 187.3)



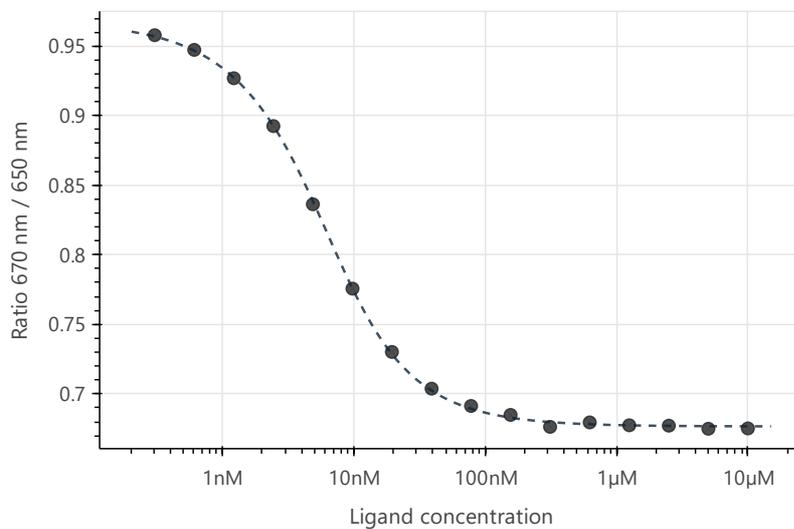
22°C | $K_d = 1.76 \pm 0.05$ nM (S/N = 192.4)



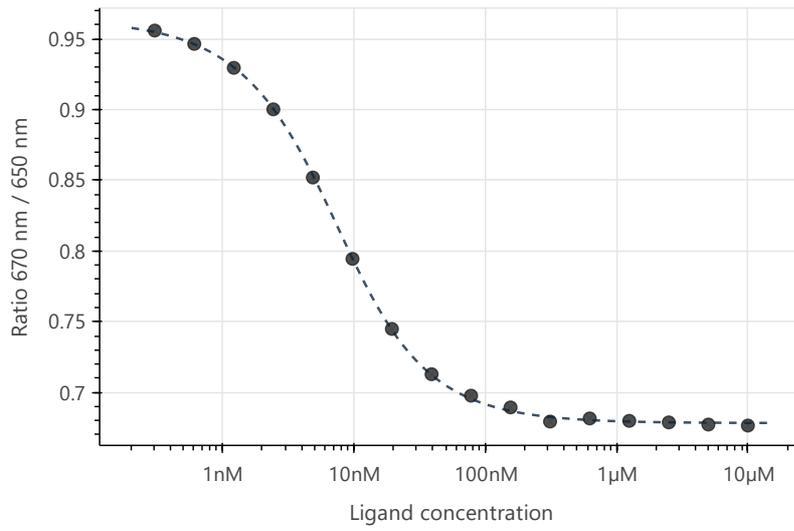
23°C | $K_d = 2.44 \pm 0.06$ nM (S/N = 187.8)



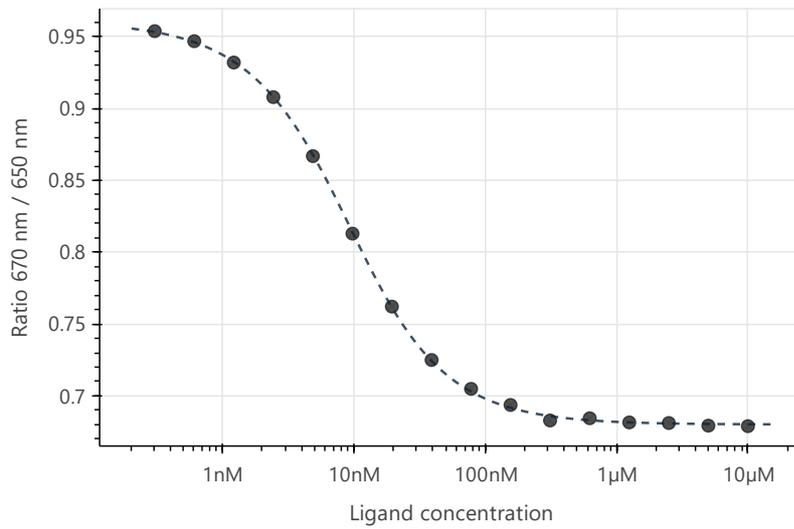
24°C | $K_d = 3.35 \pm 0.08$ nM (S/N = 181.6)



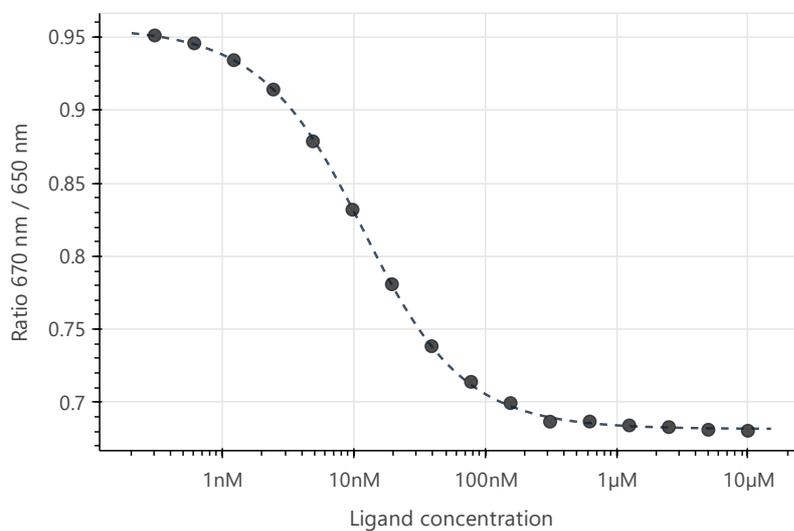
25°C | $K_d = 4.69 \text{ nM} \pm 0.1$ (S/N = 192.0)



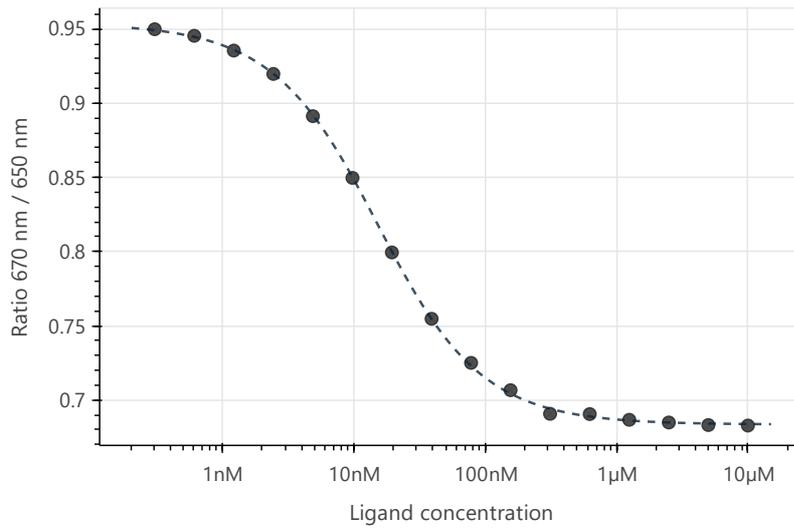
26°C | $K_d = 6.53 \text{ nM} \pm 0.12$ (S/N = 213.2)



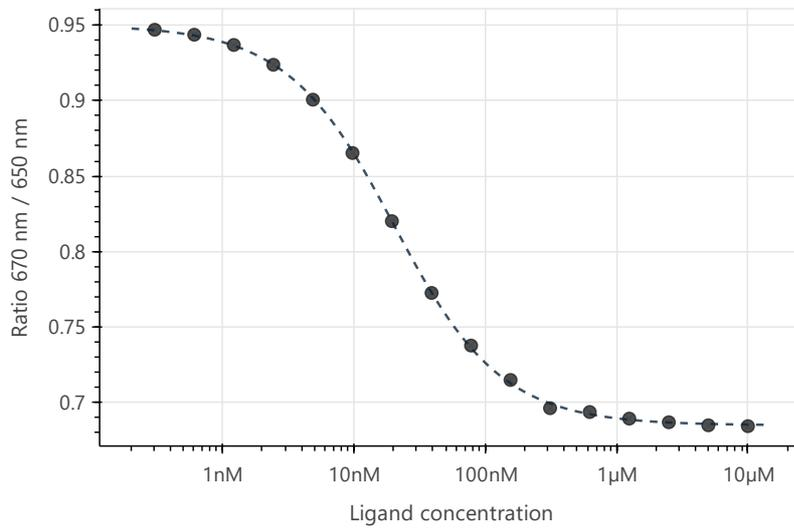
27°C | $K_d = 9.11 \text{ nM} \pm 0.16$ (S/N = 208.8)



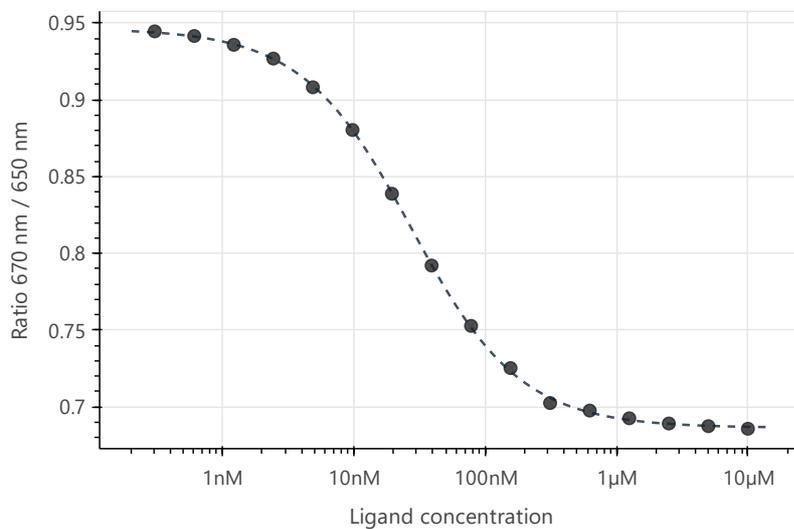
28°C | $K_d = 12.6 \pm 0.2$ nM (S/N = 193.2)



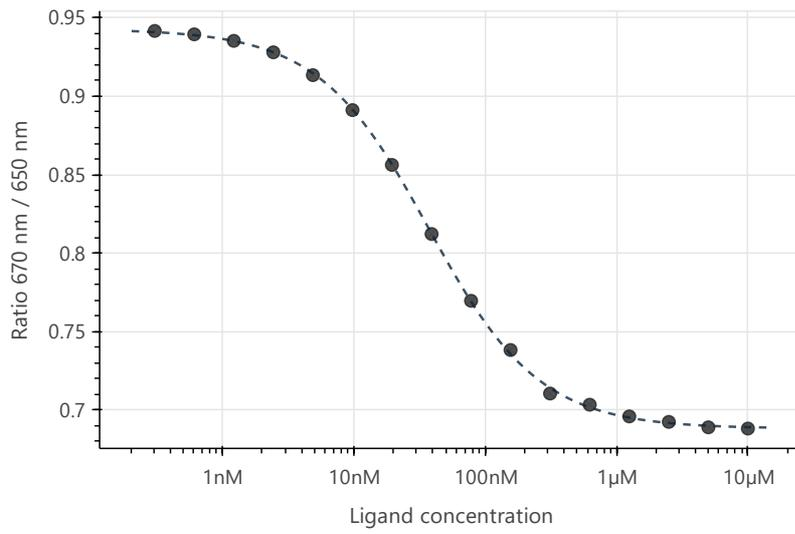
29°C | $K_d = 17.7 \pm 0.3$ nM (S/N = 207.1)



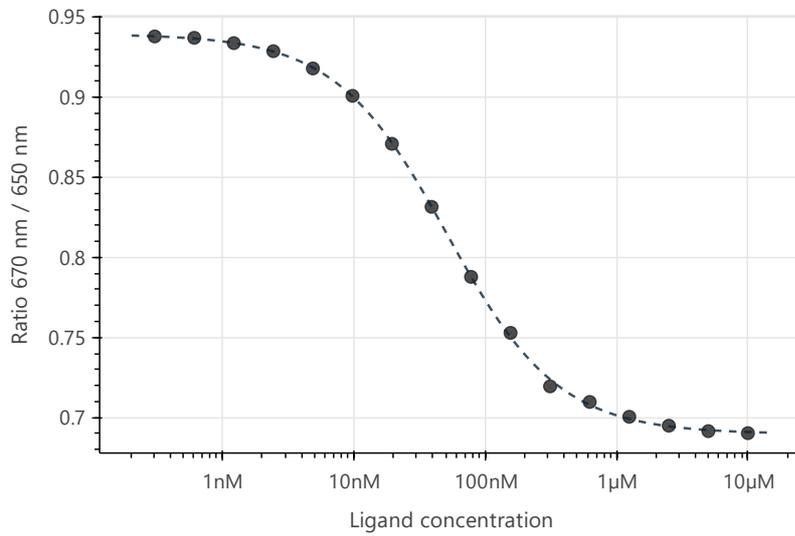
30°C | $K_d = 24.9 \pm 0.4$ nM (S/N = 204.1)



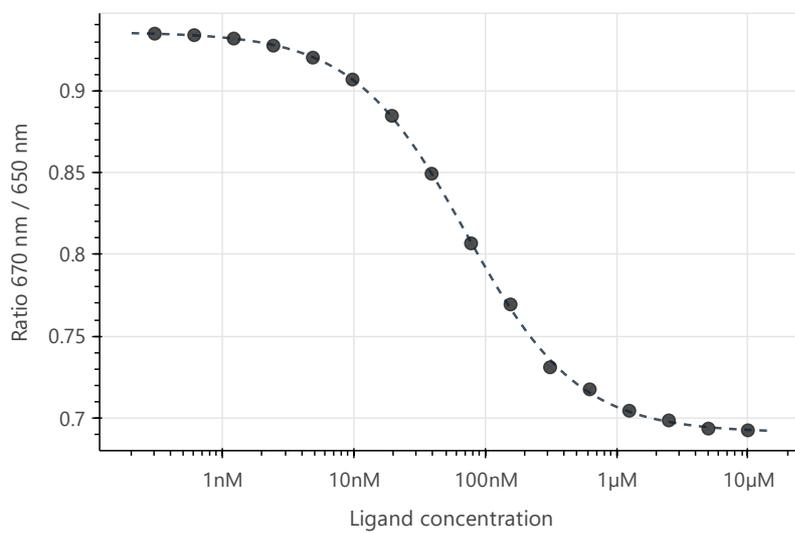
31°C | $K_d = 34.8 \pm 0.6$ nM (S/N = 185.5)



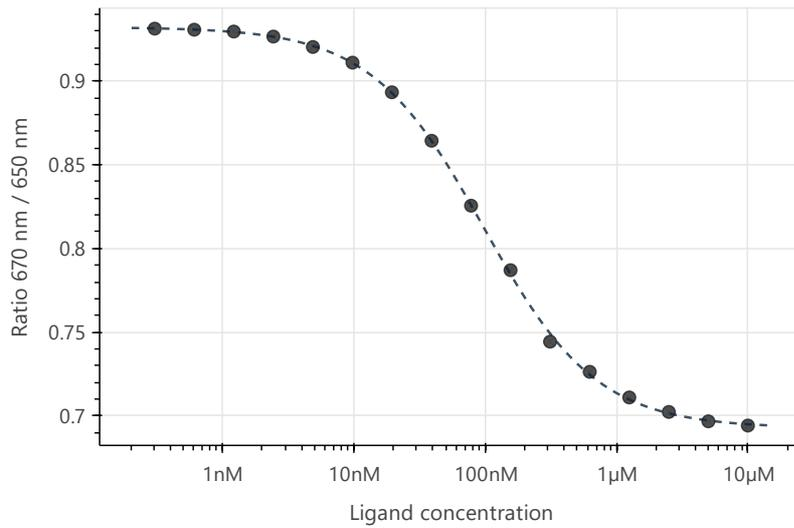
32°C | $K_d = 48.8 \pm 1$ nM (S/N = 166.5)



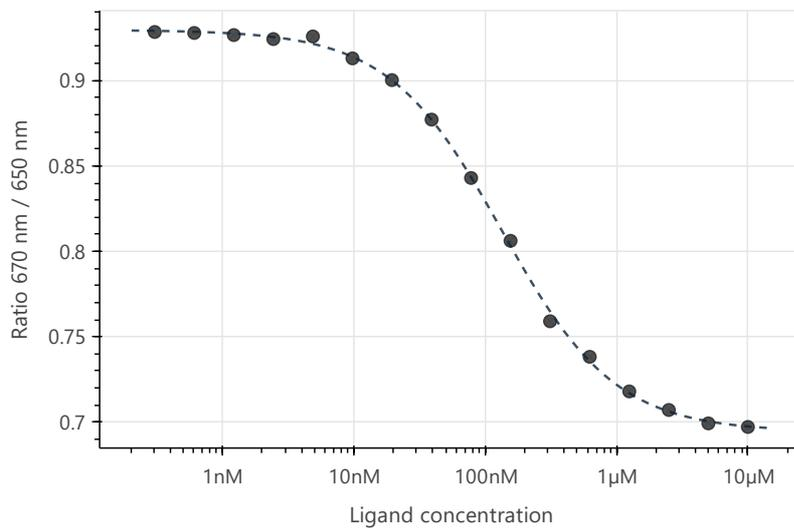
33°C | $K_d = 68.9 \pm 1.4$ nM (S/N = 158.1)



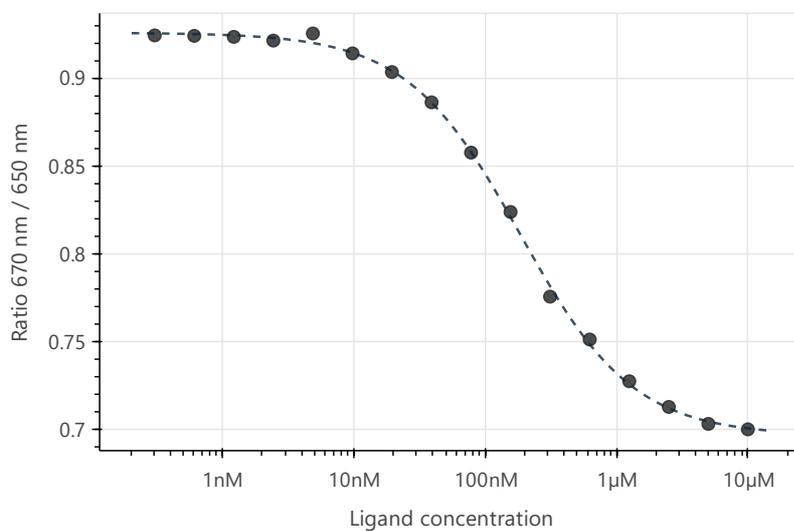
34°C | $K_d = 96.6 \pm 2 \text{ nM}$ (S/N = 151.1)



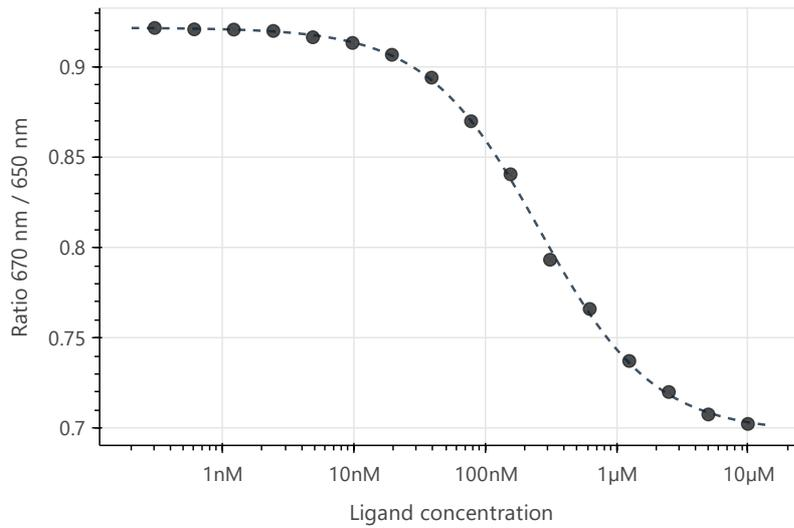
35°C | $K_d = 134 \pm 4 \text{ nM}$ (S/N = 103.4)



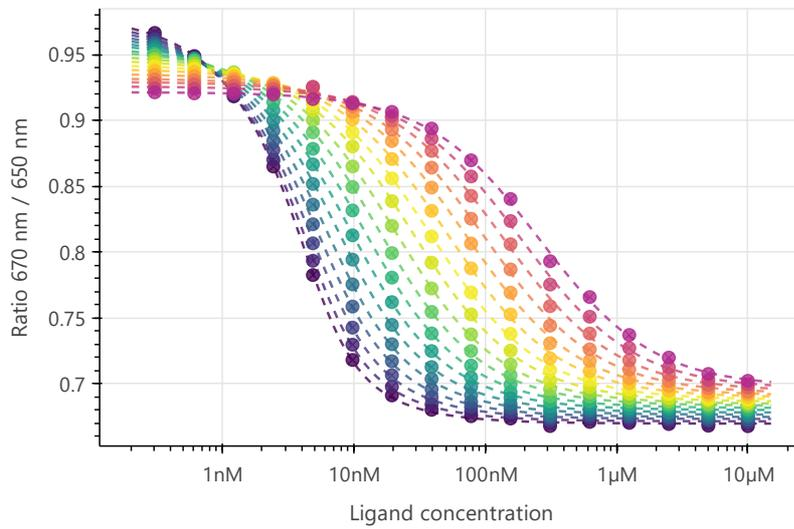
36°C | $K_d = 182 \pm 7 \text{ nM}$ (S/N = 90.7)



37°C | $K_d = 257 \pm 8 \text{ nM}$ (S/N = 106.7)



Overlay

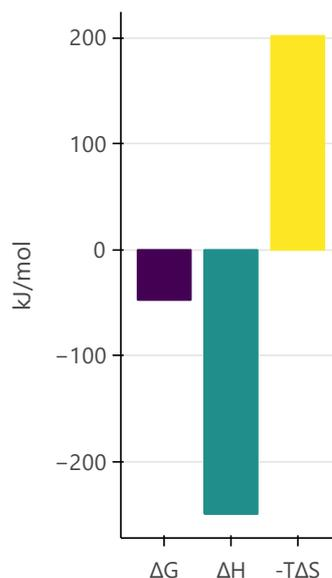
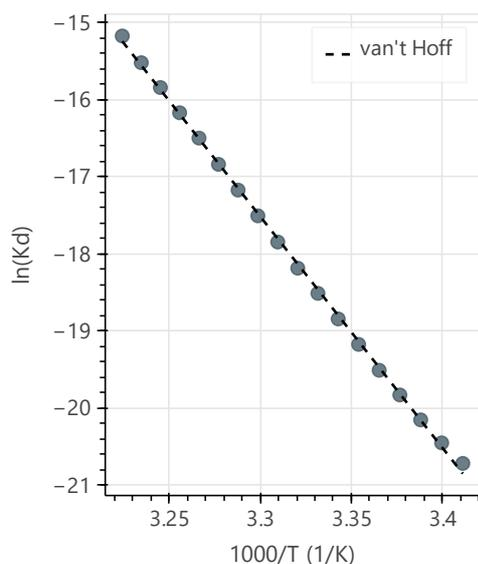


Overview of determined K_d values at different temperatures:

T (°C)	20	21	22	23	24	25	26	27	28
K_d (nM)	1.00	1.31	1.76	2.44	3.35	4.69	6.52	9.11	12.6

T (°C)	29	30	31	32	33	34	35	36	37
K_d (nM)	17.7	24.9	34.8	48.8	68.9	96.6	134	186	257

Van't Hoff analysis³:



$$\Delta H = \text{slope} \cdot R = -30096 \text{ K} \cdot 8.314 \frac{\text{J}}{\text{mol} \cdot \text{K}} = -250.2 \frac{\text{kJ}}{\text{mol}} = -59.8 \frac{\text{kcal}}{\text{mol}}$$

$$\Delta S = -\text{intercept} \cdot R = -81.804 \cdot 8.314 \frac{\text{J}}{\text{mol} \cdot \text{K}} = -680.1 \frac{\text{J}}{\text{mol} \cdot \text{K}} = -162.4 \frac{\text{cal}}{\text{mol} \cdot \text{K}}$$

D5. Reference Results/Supporting Results

$K_d = 11.6 \text{ nM}$ (23°C)
sequence CTC ACA ACA G

$K_d = 6.77 \text{ nM}$ (23°C)
sequence CTC ACA ACA G

$K_d = 3.9 \text{ nM}$ (25°C) | $\Delta H = -56.1 \text{ kcal/mol}$ | $\Delta S = -149.5 \text{ cal/(mol K)}$
sequence GAC GTG CGA AG

Surface Plasmon Resonance (SPR)
[Palau and Di Primo, Biochimie 94 \(2012\) 1891-1899](#)

switchSENSE
[Dynamic Biosensors, Application Note](#)

Isothermal Titration Calorimetry (ITC)
[Halvorsen et al., Anal Biochem 465 \(2014\) 127-133](#)

E. Contributors

Andreas Langer⁴

³ Calculations can be performed with Monolith X's Thermodynamics Measurement mode in MO.Control 2.7.0 and later versions. Plots were created outside of MO.Control 2 with temperature set at 25°C (298.15K) for ΔG calculation.

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