

Monolith X Protocol MOX-P-118

# KRAS G12D (biotinylated) – BI-2852 & MRTX1133

BI-2852 is a potent inhibitor for in vitro use that directly targets GTP-bound KRAS, which is the major form present in cancer cells carrying KRAS mutations. MRTX1133 is a noncovalent, potent, and selective KRAS G12D inhibitor which optimally fills the switch II pocket and extends three substituents to favorably interact with the protein. It selectively inhibits KRAS G12D mutant, but not KRAS wild-type, tumor cells. MRTX1133 has single digit nanomolar activity in cellular assays and marked in vivo efficacy in tumor models harboring KRAS G12D mutations.

protein – small molecule | biotinylated target | competition assay | thermodynamics | van't Hoff analysis | enthalpy | entropy

---

## A1. Target/Fluorescent Molecule

KRAS<sup>G12D</sup>

[uniprot.org/uniprot/P01116](https://uniprot.org/uniprot/P01116)

## A2. Molecule Class/Organism

GTPase

Human (*Homo sapiens*)

## A3. Sequence/Formula

MTE<sup>Y</sup>KLVVVG ADG<sup>V</sup>GKSALT IQLIQNHFVD EYDPTIEDS<sup>Y</sup> RKQVVIDGET CLLDILD<sup>T</sup>AG QEE<sup>Y</sup>SAMRDQ YMRTGEGFLC  
VFAINNTKSF EDIH<sup>Y</sup>H<sup>Y</sup>RE<sup>Y</sup>QI KRVK<sup>D</sup>SE<sup>D</sup>V<sup>P</sup> MVLVGNKCDL PSRTVDTKQA QDLARS<sup>Y</sup>GIP FIETSAKTRQ GVDDAF<sup>Y</sup>TLV  
REIRKHKEK

## A4. Purification Strategy/Source

N-terminal His-tag and Avi-tag, biotinylated

Sino Biological

[12259-H56E-B](#)

## A5. Stock Concentration/Stock Buffer

0.25 mg/mL | 10.5 μM

50 mM Tris-HCl, pH 8.0

## A6. Molecular Weight/Extinction Coefficient

23.8 kDa

13,410 M<sup>-1</sup>cm<sup>-1</sup> (ε<sub>280</sub>)

## A7. Dilution Buffer

20 mM HEPES, pH 7.4, 150 mM NaCl, 2 mM MgCl<sub>2</sub>, 0.5 mM TCEP, 0.01 % TWEEN® 20<sup>1</sup>

## A8. Labeling Strategy

Biotinylated Target Labeling Kit (Cat# NT-L020)  
1\* 20 pmol labeling dye (50 µL, 400 nM)

## A9. Labeling Procedure

1. Mix 1.9 µL of 10.5 µM KRAS with 18.1 µL of dilution buffer to obtain 20 µL of a 1 µM KRAS solution.
2. Mix 1 µL of labeling dye with 99 µL of dilution buffer to obtain 100 µL of a 4 nM dye solution.

### Affinity of dye to KRAS

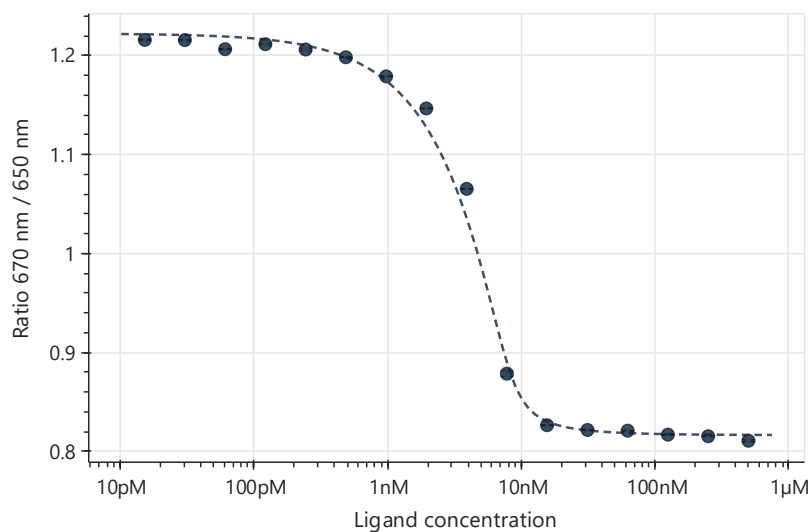
3. Take a PCR tube and mix 80 µL of 4 nM dye with 80 µL of dilution buffer to obtain 160 µL of a 2 nM dye solution.
4. Prepare a PCR-rack with 16 PCR tubes. Transfer 10 µL of the 2 nM dye solution into tubes **2** to **16**. Then, mix 10 µL of 1 µM KRAS with 10 µL of 4 nM dye in tube **1**.
5. Prepare a 1:1 serial dilution by transferring 10 µL from tube to tube. Mix carefully by pipetting up and down.
6. Incubate for 30 minutes at room temperature in the dark before loading capillaries.

### KRAS labeling

7. Mix 4 µL of 400 nM dye with 0.8 µL of KRAS and 195.2 µL of dilution buffer to obtain 200 µL of a 40 nM KRAS, 8 nM dye solution. Keep labeled KRAS on ice in the dark.

## A10. Labeling Efficiency

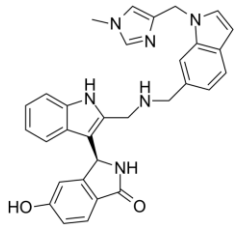
Affinity between labeling dye and biotinylated KRAS –  $K_d = 277 \pm 107$  pM (S/N = 38.2)



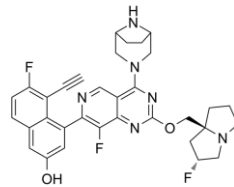
<sup>1</sup> To prepare the dilution buffer, mix 100 µL of HEPES (1 M, pH 7.4) with 250 µL of NaCl (3 M), 5 µL of TCEP (500 mM), 10 µL of MgCl<sub>2</sub> (1 M), 5 µL of TWEEN® 20 (10%) and 4.6 mL of ddH<sub>2</sub>O.

## B1. Ligand/Non-Fluorescent Binding Partner

BI-2852



MRTX1133



## B2. Molecule Class/Organism

KRAS inhibitors

## B3. Sequence/Formula

$C_{31}H_{28}N_6O_2$

$C_{33}H_{31}F_3N_6O_2$

## B4. Purification Strategy/Source

opnme.com  
BI-2852

MedChemExpress  
HY-134813

## B5. Stock Concentration/Stock Buffer

1 mg/mL | 2 mM  
DMSO

1.2 mg/mL | 2 mM  
DMSO

## B6. Molecular Weight/Extinction Coefficient

516.59 Da

600.63 Da

## B7. Serial Dilution Preparation

### BI-2852

1. Prepare a 2 mM stock solution of BI-2852 in DMSO.
2. Mix 2  $\mu$ L of 2 mM BI-2852 with 48  $\mu$ L of dilution buffer to obtain 48  $\mu$ L of an 80  $\mu$ M BI-2852 solution.
3. Mix 20  $\mu$ L of DMSO with 480  $\mu$ L of dilution buffer to obtain 500  $\mu$ L of dilution buffer with 4% DMSO.
4. Take a PCR tube and mix 40  $\mu$ L of dilution buffer with 40  $\mu$ L of 40 nM KRAS, 8 nM dye to obtain 80  $\mu$ L of 20 nM KRAS, 4 nM dye.
5. Take a fresh PCR tube and mix 60  $\mu$ L of dilution buffer containing 4% DMSO with 60  $\mu$ L of 20 nM KRAS, 4 nM dye to obtain 120  $\mu$ L of a 10 nM KRAS, 2 nM dye solution.
6. Prepare a PCR-rack with 12 PCR tubes. Transfer 10  $\mu$ L of the 10 nM KRAS, 2 nM dye solution into tubes 2 to 12. Then, mix 10  $\mu$ L of 80  $\mu$ M BI-2852 with 10  $\mu$ L of 20 nM KRAS, 4 nM dye in tube 1.
7. Prepare a 1:1 serial dilution by transferring 10  $\mu$ L from tube to tube. Mix carefully by pipetting up and down.
8. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

### MRTX1133 (competitive assay)

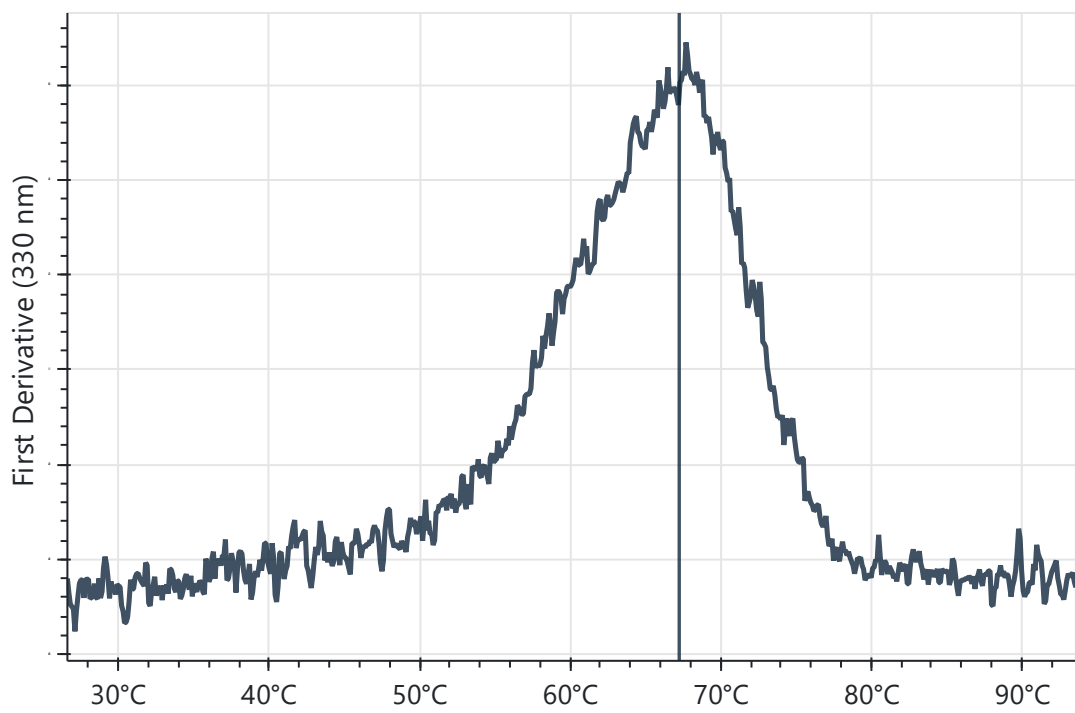
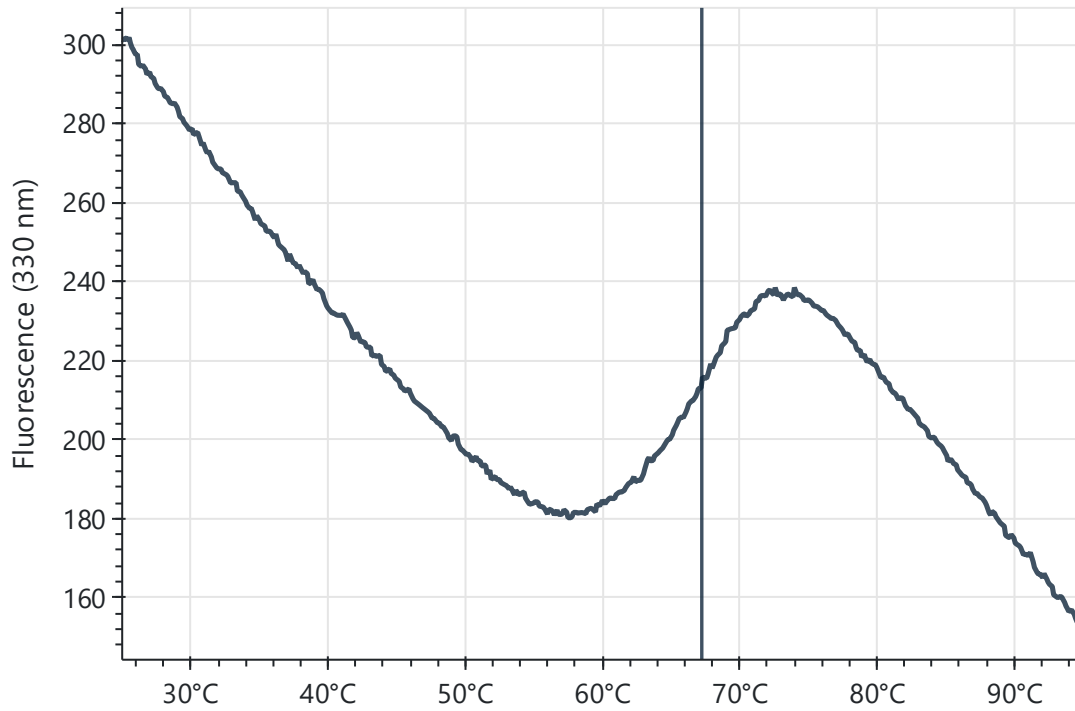
1. Prepare a 2 mM stock solution of MRTX1133 in DMSO.
2. Mix 1  $\mu\text{L}$  of 2 mM MRTX1133 with 999  $\mu\text{L}$  of dilution buffer to obtain 1 mL of 2  $\mu\text{M}$  MRTX1133.
3. Mix 4  $\mu\text{L}$  of 2 mM BI-2852 with 46  $\mu\text{L}$  of dilution buffer. Then, add 50  $\mu\text{L}$  of 40 nM KRAS, 8 nM dye to obtain 100  $\mu\text{L}$  of a 20 nM KRAS, 4 nM dye, 80  $\mu\text{M}$  BI-2852 solution.
4. Take a fresh PCR tube and mix 80  $\mu\text{L}$  of dilution buffer with 80  $\mu\text{L}$  of 20 nM KRAS, 4 nM dye, 80  $\mu\text{M}$  BI-2852 to obtain 160  $\mu\text{L}$  of a 10 nM KRAS, 2 nM dye, 40  $\mu\text{M}$  BI-2852 solution.
5. Prepare a PCR-rack with 16 PCR tubes. Transfer 10  $\mu\text{L}$  of the 10 nM KRAS, 2 nM dye, 40  $\mu\text{M}$  BI-2852 solution into tubes **2** to **16**. Then, mix 10  $\mu\text{L}$  of 2  $\mu\text{M}$  MRTX1133 with 10  $\mu\text{L}$  of 20 nM KRAS, 4 nM dye, 80  $\mu\text{M}$  BI-2852 in tube **1**.
6. Prepare a 1:1 serial dilution by transferring 10  $\mu\text{L}$  from tube to tube. Mix carefully by pipetting up and down.
7. Incubate for 15 minutes at room temperature in the dark before loading capillaries.

**C. Prometheus Panta**

Validation of structural integrity of KRAS using Prometheus Panta (heating ramp of 7°C/min):

[nanotempertech.com/prometheus](http://nanotempertech.com/prometheus)

KRAS	10 µL of 1 µM KRAS	T <sub>i</sub> = 67.25°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 (NanoTemper Technologies GmbH)

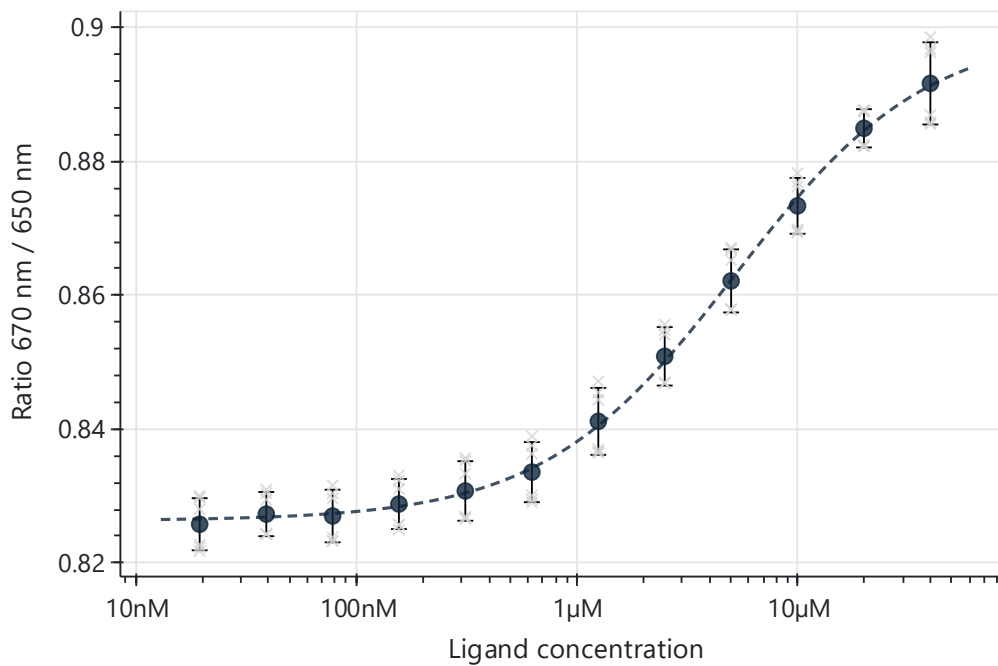
[nanotempertech.com/monolith-mo-control-software](https://nanotempertech.com/monolith-mo-control-software)

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

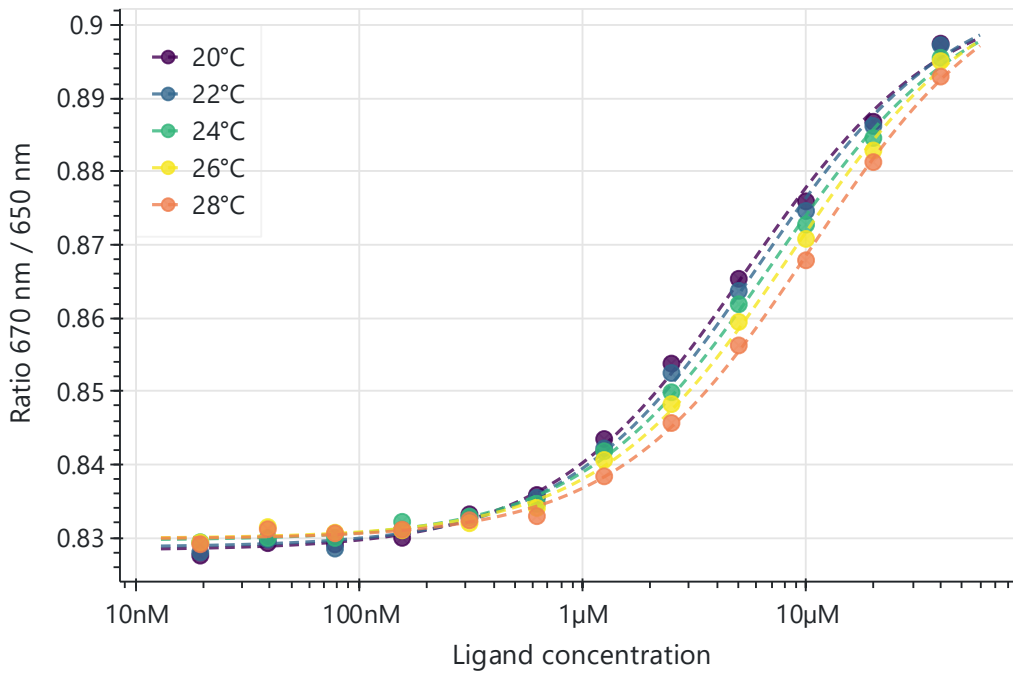
20 mM HEPES, pH 7.4, 150 mM NaCl, 2 mM MgCl<sub>2</sub>, 0.5 mM TCEP, 0.01 % TWEEN® 20, 2% DMSO  
 10 nM KRAS<sup>G12D</sup> (2 nM dye) | 40 μM -19.5 nM BI-2852 | 20°C | 100% Excitation Power

## D4. Monolith Results (Dose Response)

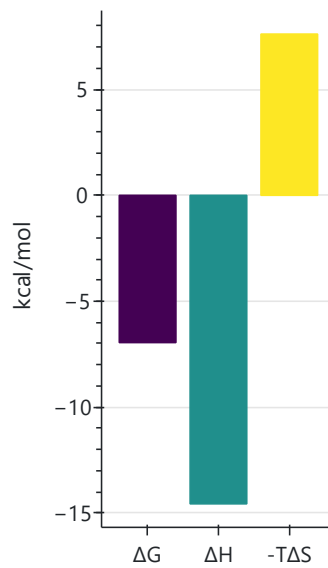
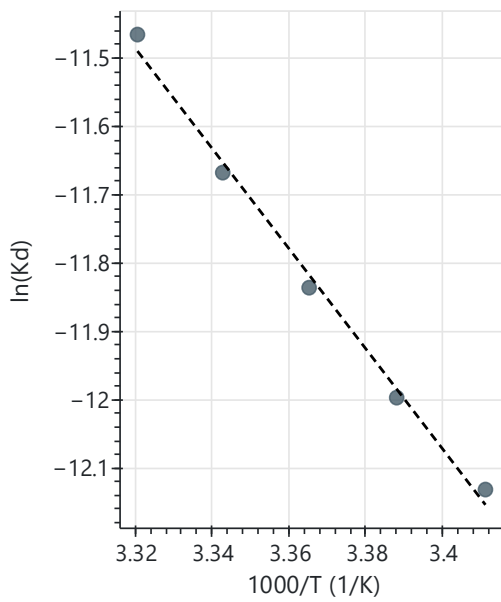
KRAS<sup>G12D</sup> - BI-2852 |  $K_d = 5.21 \pm 0.22 \mu\text{M}$  (S/N = 103.3)



Temperature dependence of BI-2852 binding to KRAS<sup>G12D</sup>



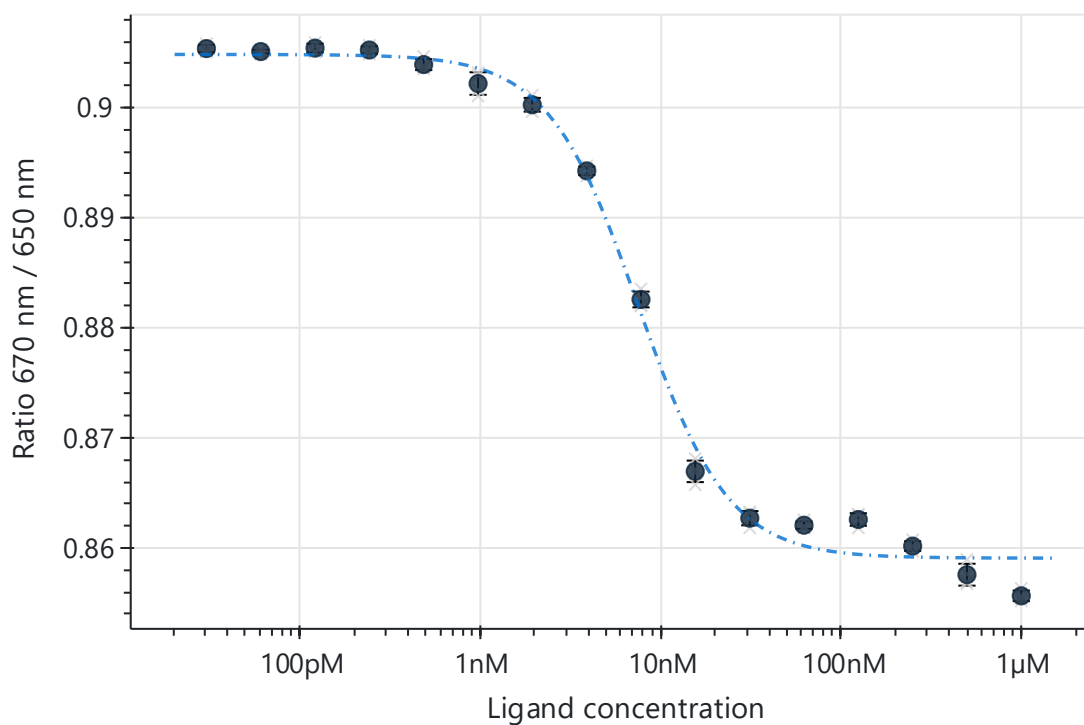
Van't Hoff analysis<sup>2</sup>



$\Delta H = -14.6 \pm 0.7$  kcal/mol  
 $\Delta S = -25.5 \pm 2.3$  cal/mol/K

<sup>2</sup> 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  $\Delta G$  calculation.

KRAS<sup>G12D</sup> (in complex with 40 μM BI-2852) – MRTX1133 | EC<sub>50</sub> = 7.56 ± 0.55 nM (S/N = 28.4) | K<sub>i</sub> = 0.3 nM<sup>3</sup>



## D5. Reference Results/Supporting Results

BI-2852	K <sub>D</sub> = 2.0 μM (GDP KRAS <sup>G12D</sup> )	Isothermal Titration Calorimetry (ITC) <a href="#">Kessler et al, PNAS 116, 15823–15829 (2019)</a>
MRTX1133	K <sub>D</sub> = 0.2 pM	Surface Plasmon Resonance (SPR) <a href="#">Wang et al, J. Med. Chem. 2022, 65, 3123–3133 (2022)</a>

## E. Contributors

Andreas Langer<sup>4</sup>

<sup>3</sup> Calculated with Nanotemper Technologies' K<sub>i</sub> calculator (input: 40.000 nM BI-2852 for 2. concentration of fluorescent competitor and 10 nM labelled KRAS<sup>G12D</sup> for 3. concentration of unlabeled target | [nanotempertech.com/user-tools/ki-calculator](https://nanotempertech.com/user-tools/ki-calculator))

<sup>4</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](https://nanotempertech.com)