

Monolith Protocol MO-P-073

Antibody – Morphine (competitive assay)

Heroin is a drug with one of the highest mortality rates. Therefore, there is an urgent need to develop alternative heroin abuse treatments. Recently, vaccines have been explored as a potential treatment modality for substances of abuse because they do not produce unwanted neurological side effects and have the potential to be utilized as preventive therapeutics against drug overdose. MicroScale Thermophoresis (MST) is a reliable method for measuring antibody binding affinities with a potential for use in vaccine development. Here, we show how the affinity of antibodies to heroin and its major metabolites such as morphine can be detected in a competitive MST assay using a fluorescent heroin analog. Moreover, the assay can even be used to determine affinities of polyclonal antibodies in sera.

antibody – small molecule interaction | heroin hapten | vaccines | competitive assay

A1. Target/Fluorescent Molecule

Anti-morphine antibody (ab1060)

A2. Molecule Class/Organism

Monoclonal antibody
Mouse (Mus musculus)

A3. Sequence/Formula

N/A

A4. Purification Strategy/Source

Protein G purified
Abcam (Cambridge, MA, USA)
[ab1060](#)

A5. Stock Concentration/Stock Buffer

1 mg/mL | ~6.7 μ M
pH 7.4

A6. Molecular Weight/Extinction Coefficient

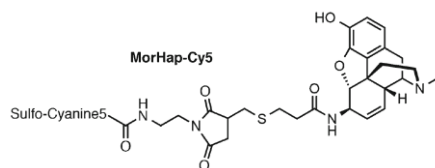
~150 kDa

A7. Dilution Buffer

10 mM Na₂HPO₄, 1.8 mM KH₂PO₄, 2.7 mM KCl, 137 mM NaCl, pH 7.4, 0.05% Tween® 20
(Dulbecco's phosphate buffered saline, DPBS-Tween)

A8. Labeling Strategy

Competitive binding assay with MorHap-Cy5¹, a fluorescent heroin analog.



A9. Labeling Procedure

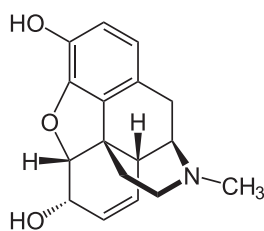
1. Mix 100 μ L of MorHap-Cy5 (1 nM) and 100 μ L of ab1060 (40 nM) to obtain 200 μ L of a 0.5 nM MorHap-Cy5, 20 nM ab1060 solution.²
2. Incubate for 20 minutes at room temperature in the dark.

A10. Labeling Efficiency

N/A

B1. Ligand/Non-Fluorescent Binding Partner

Morphine



B2. Molecule Class/Organism

Opiate

B3. Sequence/Formula

C₁₇H₁₉NO₃

B4. Purification Strategy/Source

Lipomed (Cambridge, MA, USA)

¹ For the synthesis of MorHap-Cy5, see [Torres et al., Analytical and Bioanalytical Chemistry \(2018\) 410:3885–3903](#).

² As the K_d between MorHap-Cy5 and ab1060 is ~5 nM (c.f. section D4), a final concentration of 10 nM ab1060 is sufficient for complex formation, but still low enough for accurate K_d determination of the morphine – ab1060 interaction.

B5. Stock Concentration/Stock Buffer

1 mg/mL | 3.5 mM
Methanole

B6. Molecular Weight/Extinction Coefficient

285.34 Da

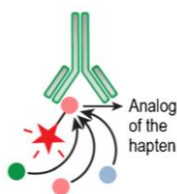
B7. Serial Dilution Preparation

Direct binding assay

1. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of a 3.2 μM solution of ab1060 into tube **1**. Then, transfer 10 μL of dilution buffer into tubes **2** to **16**.
2. Prepare a 1:1 serial dilution by transferring 10 μL from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10 μL from tube **16** to get an equal volume of 10 μL for all samples.
3. Add 10 μL of MorHap-Cy5 (0.5 nM) to each tube from **16** to **1** and mix by pipetting.
4. Incubate for 20 minutes at room temperature in the dark before loading capillaries.

Competitive binding assay

1. Mix 2.3 μL of the 3.5 mM morphine stock solution with 998 μL of dilution buffer to obtain 1 mL of a 8 μM morphine solution.
2. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of the 8 μM morphine solution into tube **1**. Then, transfer 10 μL of dilution buffer into tubes **2** to **16**.
3. Prepare a 1:1 serial dilution by transferring 10 μL from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10 μL from tube **16** to get an equal volume of 10 μL for all samples.
4. Add 10 μL of the 0.5 nM MorHap-Cy5, 20 nM ab1060 solution to each tube from **16** to **1** and mix by pipetting.
5. Incubate for 20 minutes at room temperature in the dark before loading capillaries.



D1. MST System/Capillaries

Monolith NT.115^{Pico} Red (NanoTemper Technologies GmbH)
 Capillaries Monolith NT.115 (MO-K022, NanoTemper Technologies GmbH)

D2. MST Software

MO.Control v1.6 | MO.Affinity Analysis 2.2.4 (NanoTemper Technologies GmbH)
nanotempertech.com/monolith-mo-control-software

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

Direct binding assay

10 mM Na₂HPO₄, 1.8 mM KH₂PO₄, 2.7 mM KCl, 137 mM NaCl, pH 7.4, 0.05% Tween® 20
 0.25 nM MorHap-Cy5 | 3.2 μM – 97 pM ab1060 | 25°C | low MST power | 20% excitation power

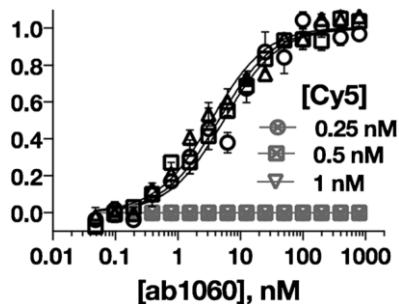
Competitive binding assay

10 mM Na₂HPO₄, 1.8 mM KH₂PO₄, 2.7 mM KCl, 137 mM NaCl, pH 7.4, 0.05% Tween® 20
 0.25 nM MorHap-Cy5, 10 nM ab1060 | 4 μM – 122 pM morphine | 25°C | low MST power | 20% excitation power

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

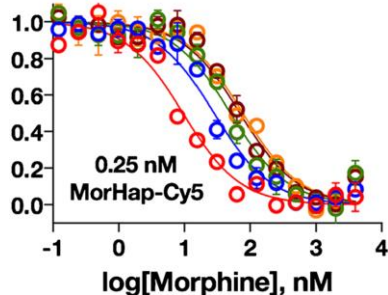
Direct binding assay

$K_d = 5.29 \pm 1.67$ nM (MorHap-Cy5 – ab1060)



Competitive binding assay

$EC_{50} = 28.2 \pm 1.00$ nM | $K_i = 6.87 \pm 2.99^3$ (ab1060 – morphine)



³ For calculation of K_i , see also the NanoTemper 'FAQ Competitive Binding Assay'.

D5. Reference Results/Supporting Results

$K_d = 1.97 \pm 0.17$ nM	Equilibrium dialysis with ultra-performance liquid chromatography/ tandem mass spectrometry (ED-UPLC/MS/MS) Torres et al., Analytical and Bioanalytical Chemistry (2018) 410:3885–3903
$K_d = 2$ nM	UV-VIS Abcam

E. Contributors

Oscar B. Torres⁴, Wyatt Strutz⁵, Andreas Langer⁶

⁴ Laboratory of Adjuvant and Antigen Research, Walter Reed Army Institute of Research, Silver Spring, MD, USA

⁵ NanoTemper Technologies, Inc., Cambridge, MA, USA

⁶ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com