

Monolith Protocol MO-P-079

Antibody – TNT

The illegal use of explosives by terrorists and other criminals is an increasing issue in public spaces, such as airports, railway stations, highways, sports venues, theaters, and other large buildings. The fast and extremely sensitive detection of explosives is one of the most relevant tasks to guarantee security in areas of public access. 2,4,6-trinitrotoluene (TNT) is a small chemical compound, which is best known as an explosive material with convenient handling properties. In this study the binding between an antibody and TNT was analyzed.

protein – small molecule interaction | TNT | Explosives

A1. Target/Fluorescent Molecule

Monoclonal antibody Clone A.1.1.1

A2. Molecule Class/Organism

Immunoglobulin IgG1

mouse

A3. Sequence/Formula

N/A

A4. Purification Strategy/Source/Batch-No.

Hydrophobic Interaction Chromatography | SDIX (Newark,USA) | 107415

A5. Stock Concentration/Stock Buffer

10.9 g/L | 73 μ M

PBS with 0.05 % NaN_3 ,

A6. Molecular Weight/Extinction Coefficient

150 kDa

A7. Assay Buffer

8.1 mM disodium phosphate, 1.5 mM potassium dihydrogen phosphate, 2.7 mM potassium chloride, 137 mM sodium chloride, pH 7.4, 0.1% BSA

A8. Labeling Strategy

Fluorescent labeling of the monoclonal antibody A.1.1.1 for MST measurements was done using the Dy-654-NHS labeling kit (#654-01) according to the manufacturer's protocol (Dyomics GmbH, Jena, Germany).

A9. Labeling Procedure

1. Prepare 100 mL coupling buffer containing 137 mM NaCl and 100 mM phosphate (pH 7.8).
 2. Prepare 100 mL PBS (8.1 mM disodium phosphate, 1.5 mM potassium dihydrogen phosphate, 2.7 mM potassium chloride, 137 mM sodium chloride, pH 7.4)
 3. Prepare 100 μ L of a 13 μ M A.1.1.1 solution in PBS.
 4. Dissolve 200 μ g of Dy-654-NHS in 40 μ L of dry amine-free DMF to obtain a 4.53 mM solution (5g/L).
 5. Equilibrate a SpinTrap G-25 column (Cytiva) four times with 140 μ L coupling buffer and centrifuge at 800 g for one minute and 4°C.
 6. Add the 100 μ L A.1.1.1 solution (2 g/L, 13 μ M) to the SpinTrap, subsequently add 40 μ L of coupling buffer, centrifuge and collect the 140 μ L eluate.
 7. Add 1.68 μ L of the 4.53 mM Dy-654-NHS solution (7.6 nmol) to the 140 μ L antibody solution and mix immediately by careful resuspending with the pipette. Incubate for two hours at RT in a shaker at 800 rpm.
 8. Equilibrate a SpinTrap four times with 140 μ L of PBS as described above. Add the labeling reaction on the SpinTrap and centrifuge at 800 g for one minute at 4°C. As eluate 140 μ L of light blue solution of approx. 10 μ M A1.1.1-Dy654 are obtained.
 9. Mix 1.5 μ L of the labeled A.1.1.1 with 7 498 μ L assay buffer to obtain 7 500 mL of ~2 nM A.1.1.1-Dy-654
-

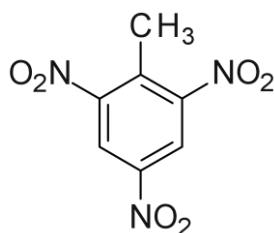
B1. Ligand/Non-Fluorescent Binding Partner

2,4,6-trinitrotoluene (TNT)

B2. Molecule Class/Organism

Small molecule (explosives)

B3. Sequence/Formula



$C_6H_2(NO_2)_3CH_3$

B4. Purification Strategy/Source/Batch-No.

Supplied by BAM division 2.3

B5. Stock Concentration/Stock Buffer

4.54 g/L | 20 mM

Absolute Ethanol

B6. Molecular Weight

227 g/mol

B7. Serial Dilution Preparation

1. Dilute 5 μ L of 20 mM TNT stock in 995 μ L absolute ethanol to obtain a 100 μ M TNT solution
2. Dilute 2 μ L of 100 μ M TNT solution with 398 μ L assay buffer to obtain a 500 nM TNT solution
3. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μ L of the 500 nM TNT solution into tube **1**. Then, transfer 10 μ L of assay 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. Add 10 μ L of 2 nM labeled A.1.1.1-Dy-654 to each tube from **16** to **1** and mix by pipetting.
6. Load capillaries directly.

D1. MST System/Capillaries

Monolith Pico (NanoTemper Technologies GmbH)

Capillaries Monolith NT.115 (MO-K022, NanoTemper Technologies GmbH)

D2. MST Software

MO.Control v1.6 (NanoTemper Technologies GmbH)

nanotempertech.com/monolith-mo-control-software

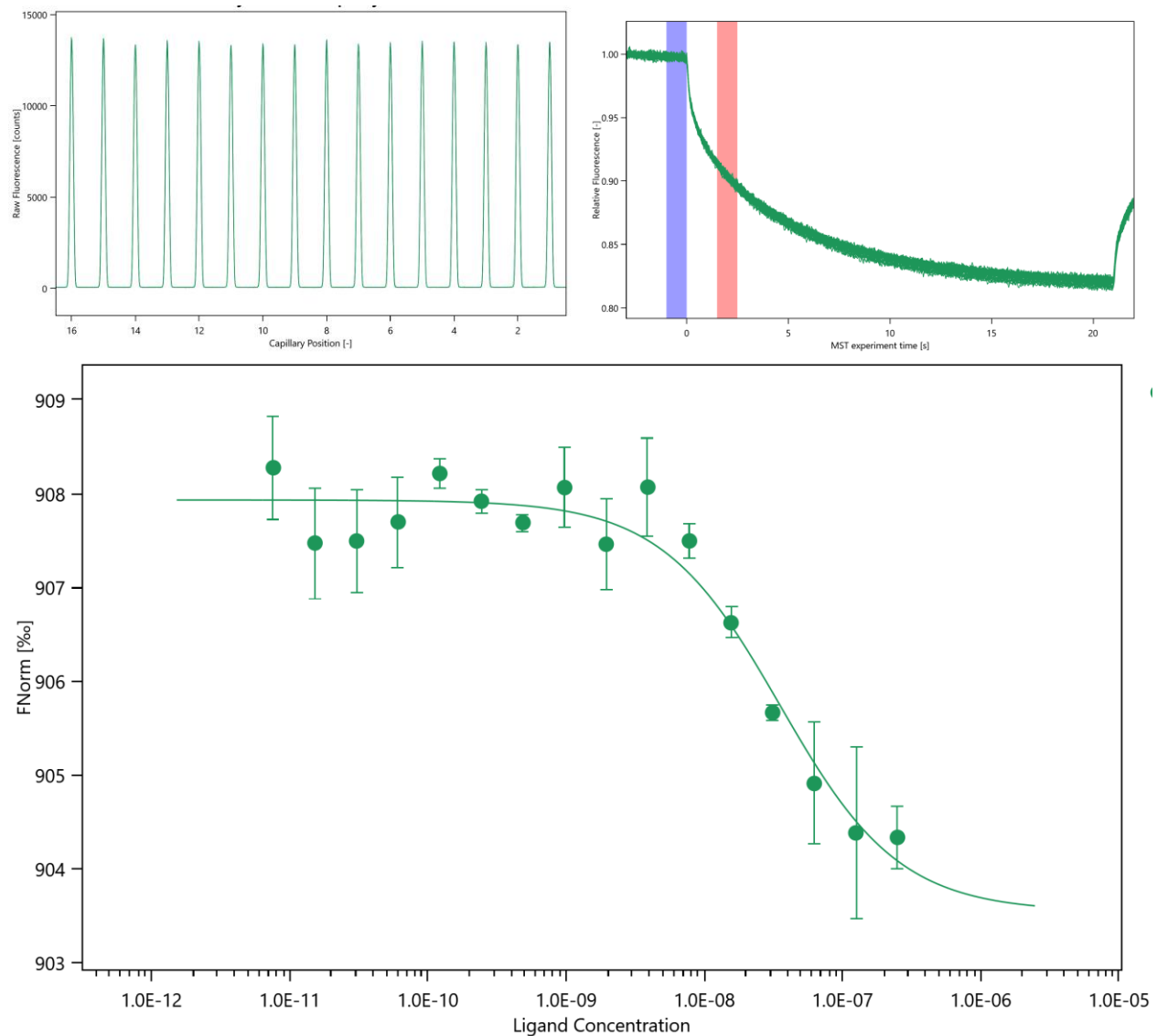
D3. MST Experiment (Buffer/Concentrations/Temperature/MST Power/Excitation Power)

137 mM NaCl, 9.6 mM phosphate, 0.1 % BSA, pH 7.4

1 nM A.1.1.1-Dy654 | 7.6 pM – 500 nM TNT | 22°C | Medium MST Power | 30 % Excitation Power

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

$K_D = 37 \pm 7$ nM



D5. Supporting Results

Paul et al., *Biosensors*, 10(8): 89 (2020)

E. Contributors

Martin Paul | Division Protein Analysis, Federal Institute for Materials Research and Testing (BAM)¹

Michael G. Weller | Division Protein Analysis, Federal Institute for Materials Research and (BAM)¹

Dr. Teresia Hallström | NanoTemper Technologies GmbH² | nanotempertech.com

¹ Richard-Willstätter-Straße 11, 12489 Berlin, Germany

² Flößergasse 4, 81369 München, Germany