

## Monolith Protocol MO-P-008

# Maltose Binding Protein – Maltose (label-free)

Maltose binding protein (MBP) is part of the periplasmic transport system of *Escherichia coli* and involved in the transport of maltose into the bacterium. It binds the disaccharide once it has crossed the outer membrane, and then assists its translocation across the inner membrane. Additionally, it is often used as a fusion tag for protein purification or solubilization.

protein – small molecule interaction | carbohydrate | conformational change | label-free

## A1. Target/Fluorescent Molecule

Maltose/maltodextrin-binding periplasmic protein (MBP)

[uniprot.org/uniprot/PDAEX9](https://uniprot.org/uniprot/PDAEX9)

## A2. Molecule Class/Organism

Periplasmic protein

*Escherichia coli*

## A3. Sequence/Formula

KIEEGKLVIV INGDKGYNGL AEVGKKFEKD TGIKVTVEHP DKLEEKFPQV AATGDGPDII FWAHDRFGGY AQSGLLAEIT  
 PDKAFQDKLY PFTWDAVRYN GKLIAYPIAV EALSIIYNKD LLPNPPKTWE EIPALDKELK AKGKSALMFN LQEPYFTWPL  
 IAADGGYAFK YENGKYDIKD VGVDNAGAKA GLTFLVDLIK NKHMNADTDY SIAEAAFNKG ETAMTINGPW AWSNIDTSKV  
 NYGVTVLPTF KGQPSKPFVG VLSAGINAAS PNKELAKEFL ENYLLTDEGL EAVNKDKPLG AVALKSYEEE LAKDPRIAAT  
 MENAQKGEIM PNIPQMSAFW YAVRTAVINA ASGRQTVDEA LKDAQTRITK

## A4. Purification Strategy/Source

N/A

## A5. Stock Concentration/Stock Buffer

0.5 mg/mL | 12 µM

Phosphate-buffered saline, 10% glycerol, 0.1% Pluronic® F-127

## A6. Molecular Weight/Extinction Coefficient

42 kDa

66,350 M<sup>-1</sup>cm<sup>-1</sup> (ε<sub>280</sub>)

## A7. Dilution Buffer

Phosphate-buffered saline (PBS, pH 7.4)

## A8. Labeling Strategy

Trp and Tyr fluorescence

## A9. Labeling Procedure

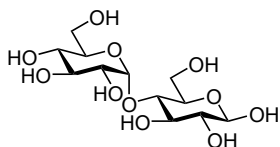
1. Mix 4  $\mu\text{L}$  of Pluronic® F-127 (5%) with 187  $\mu\text{L}$  of PBS.
2. Add 8.3  $\mu\text{L}$  of 12  $\mu\text{M}$  MBP to obtain 200  $\mu\text{L}$  of a 500 nM MBP solution in PBS with 0.1% Pluronic® F-127.

## A10. Labeling Efficiency

N/A

## B1. Ligand/Non-Fluorescent Binding Partner

D-(+)-Maltose monohydrate (maltose)



## B2. Molecule Class/Organism

Carbohydrate

## B3. Sequence/Formula

$\text{C}_{12}\text{H}_{22}\text{O}_{11}$

## B4. Purification Strategy/Source

Sigma-Aldrich GmbH

M9171

## B5. Stock Concentration/Stock Buffer

Powdered

## B6. Molecular Weight/Extinction Coefficient

360.31 Da

## B7. Serial Dilution Preparation

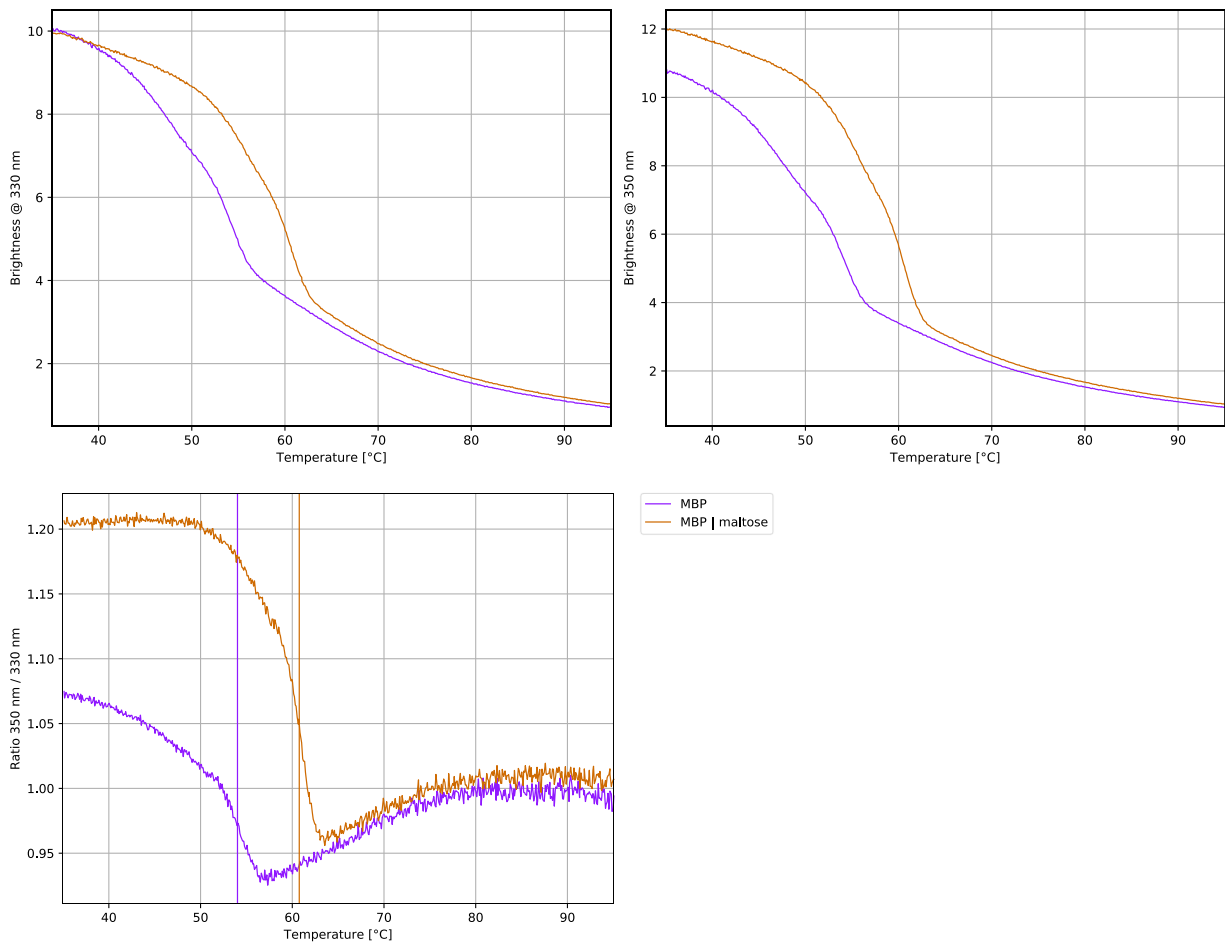
1. Dissolve 10 mg of maltose monohydrate in 55.5  $\mu\text{L}$  of ddH<sub>2</sub>O to obtain a 500 mM maltose solution.
2. Mix 4  $\mu\text{L}$  of 500 mM maltose with 196  $\mu\text{L}$  of dilution buffer to obtain 200  $\mu\text{L}$  of a 10 mM maltose solution.
3. Mix 8  $\mu\text{L}$  of 10 mM maltose with 192  $\mu\text{L}$  of dilution buffer to obtain 200  $\mu\text{L}$  of a 400  $\mu\text{M}$  maltose solution.
4. Prepare a PCR-rack with 16 PCR tubes. Transfer 20  $\mu\text{L}$  of the 400  $\mu\text{M}$  maltose solution into tube **1**. Then, transfer 10  $\mu\text{L}$  of dilution buffer into tubes **2** to **16**.
5. Prepare a 1:1 serial dilution by transferring 10  $\mu\text{L}$  from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10  $\mu\text{L}$  from tube **16** to get an equal volume of 10  $\mu\text{L}$  for all samples.
6. Add 10  $\mu\text{L}$  of 500 nM MBP to each tube from **16** to **1** and mix by pipetting.
7. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

## C. Applied Quality Checks

Validation of structural integrity and functionality of MBP using Tycho NT.6:

[nanotempertech.com/tycho](https://nanotempertech.com/tycho)

MBP	5 $\mu\text{L}$ of 500 nM MBP + 5 $\mu\text{L}$ of dilution buffer (tube <b>16</b> )	$T_i = 54.0^\circ\text{C}$
MBP   maltose	5 $\mu\text{L}$ of 500 nM MBP + 5 $\mu\text{L}$ of 400 $\mu\text{M}$ maltose (tube <b>1</b> )	$T_i = 60.8^\circ\text{C}$



## D1. MST System/Capillaries

Monolith NT.LabelFree (NanoTemper Technologies GmbH)

Capillaries Monolith NT.LabelFree (MO-Z022, NanoTemper Technologies GmbH)

## D2. MST Software

MO.Control v1.6 (NanoTemper Technologies GmbH)

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

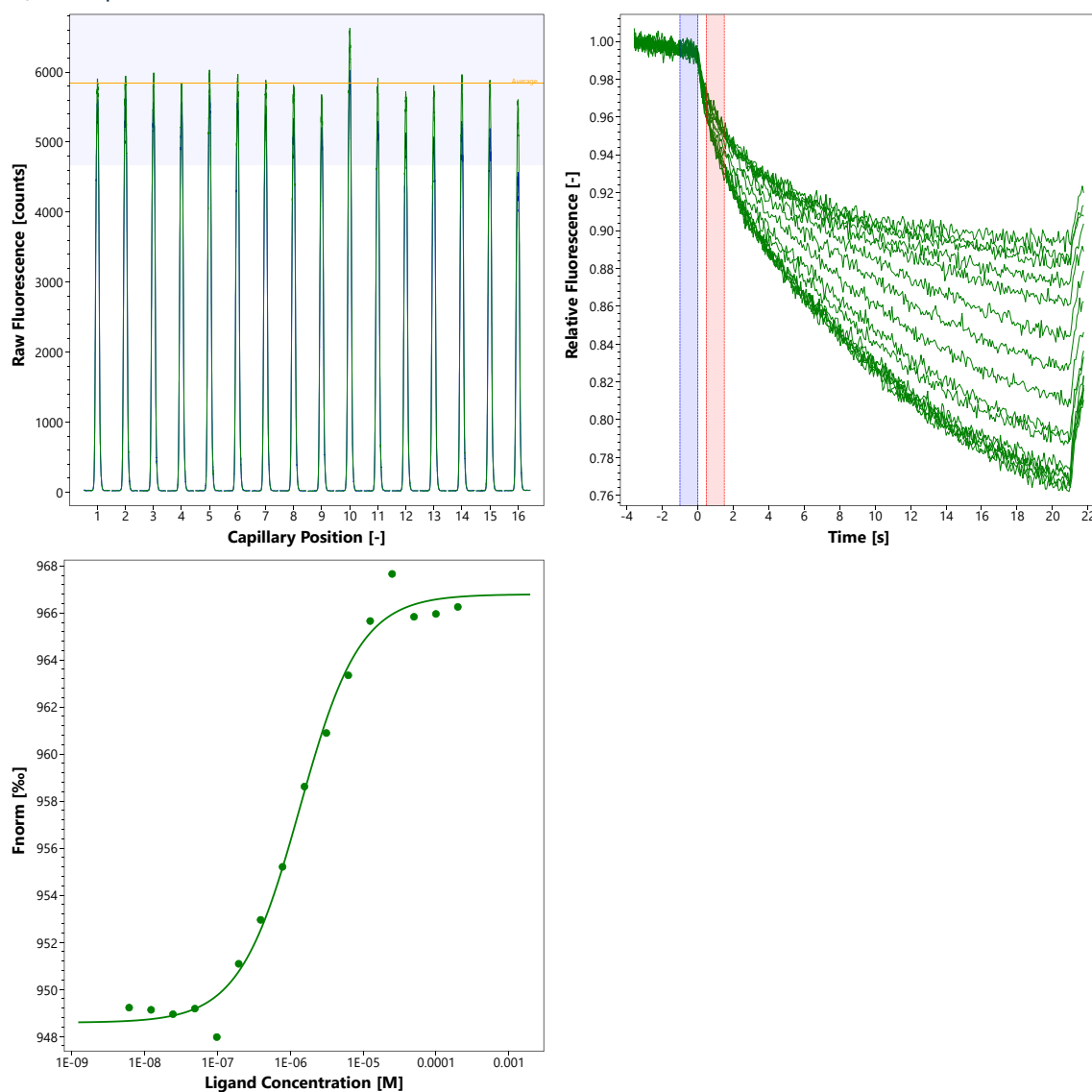
## D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

Phosphate-buffered saline (PBS, pH 7.4), 0.05% Pluronic® F-127

250 nM MBP | 200  $\mu$ M – 6.1 nM maltose | 25°C | medium MST power | 10% excitation power

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

$K_d = 1.22 \mu$ M



## D5. Reference Results/Supporting Results

$K_d = 1.2 \mu\text{M}$       Intrinsic fluorescence changes  
[Telmer and Shilton, J Biol Chem 278 \(2003\) 34555-34567](#)

---

## E. Contributors

Andreas Langer<sup>1</sup>

---

<sup>1</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](https://www.nanotempertech.com)