

## Monolith Protocol MO-P-045

# ATP Binding Cassette Transporter – Phosphite

The periplasmic binding proteins from bacterial phosphite ATP-binding cassette transporters are involved in the uptake of phosphorus compounds, which are essential for cell structure, metabolism and signaling. In many natural habitats the availability of inorganic phosphate is low enough to limit microbial growth, and under these phosphate-depleted conditions some bacteria utilize phosphite as an alternative source of phosphorus.

protein – ion interaction | ABC transporters | phosphite | His<sub>6</sub>-tag

---

### A1. Target/Fluorescent Molecule

ABC transporter periplasmic phosphonate-binding protein (Te\_PtxB)

[uniprot.org/uniprot/Q11919](https://uniprot.org/uniprot/Q11919)

### A2. Molecule Class/Organism

ATP binding cassette transporter

*Trichodesmium erythraeum* (strain IMS101)

### A3. Sequence/Formula

MLGLILKKNL FTVLFLACLS LVSCSNSNIQ KSEKANKPQK LVVALLPDES AATVIQNNKG LEMYLENKLN  
 KDIELFVSTD YSSMIEVASK GRDLAYFGP LSYVLAKTKS NIEPFAALEK DGKNTYQALV IGNAEAGINS  
 YEKIEGKIMA YGDQASTSSH LIPKSMKQK QLKAGENYEE VFVGAHDAVA IAVANGKAQA GGLSKPIFTA  
 LIERGTIDKN KVIIIAESKP FPQYPWTMRS DLDSELKTI QQAFLELEDK AILKPFKADA FTLVTDQDYD  
 VVRNLGEVLE LNFEQLNK

### A4. Purification Strategy/Source

Expressed in *E. coli* BL21 (DE3) with N-terminal His<sub>6</sub>-tag, purified by immobilized Ni-affinity chromatography and gel filtration chromatography

Produced at the University of Sheffield<sup>1</sup>

### A5. Stock Concentration/Stock Buffer

250 μM

25 mM Tris-HCl, pH 7.4, 200 mM NaCl

### A6. Molecular Weight/Extinction Coefficient

33 kDa

20,400 M<sup>-1</sup>cm<sup>-1</sup> (ε<sub>280</sub>)

---

<sup>1</sup> See Bisson et al., *Nature Communications* 2017, 8(1), 1–12 for full details.

**A7. Dilution Buffer**

50 mM HEPES, pH 7.4, 250 mM NaCl, 0.05% TWEEN® 20

**A8. Labeling Strategy**

Monolith His-Tag Labeling Kit RED-tris-NTA (MO-L008, NanoTemper Technologies GmbH)

1\* 125 pmol RED-tris-NTA Dye

**A9. Labeling Procedure**

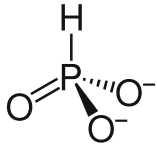
1. Suspend 125 pmol RED-tris-NTA Dye in 25  $\mu$ L of dilution buffer to obtain a 5  $\mu$ M dye solution.
2. Prepare a 100 nM dye solution by mixing 2  $\mu$ L of dye (5  $\mu$ M) and 98  $\mu$ L of dilution buffer.
3. Adjust the concentration of Te\_PtxB to 200 nM in dilution buffer in a volume of 100  $\mu$ L.
4. Mix 100  $\mu$ L of Te\_PtxB protein (200 nM) with 100  $\mu$ L of dye (100 nM).
5. Incubate for 30 minutes at room temperature in the dark.
6. Centrifuge the labeled Te\_PtxB for 10 min at 4°C and 15,000  $\times$  g.

**A10. Labeling Efficiency**

N/A

### B1. Ligand/Non-Fluorescent Binding Partner

Phosphite (as sodium phosphite dibasic pentahydrate)



### B2. Molecule Class/Organism

Inorganic compound

### B3. Sequence/Formula

$\text{Na}_2\text{HPO}_3 \cdot 5\text{H}_2\text{O}$

### B4. Purification Strategy/Source

Sigma-Aldrich, UK

04283

### B5. Stock Concentration/Stock Buffer

Powdered

### B6. Molecular Weight/Extinction Coefficient

216.04 Da

### B7. Serial Dilution Preparation

1. Dissolve 5.4 g of sodium phosphite dibasic pentahydrate in 20 mL of ddH<sub>2</sub>O, titrate the phosphite to pH 7.4 and make up to 25 mL in a volumetric flask to obtain a 1 M sodium phosphite solution.
2. Prepare 30 μL of an 80 μM solution of phosphite in dilution buffer.
3. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of the 80 μM phosphite solution into tube **1**. Then, transfer 10 μL of dilution buffer tubes **2** to **16**.
4. 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.
5. Add 10 μL of 100 nM labeled Te\_PtxB to each tube from **16** to **1** and mix by pipetting.
6. Incubate for 15 minutes at room temperature in the dark before loading capillaries.

## D1. MST System/Capillaries

Monolith NT.115 Red (NanoTemper Technologies GmbH)  
 Premium Capillaries Monolith NT.115 (MO-K025, NanoTemper Technologies GmbH)

## D2. MST Software

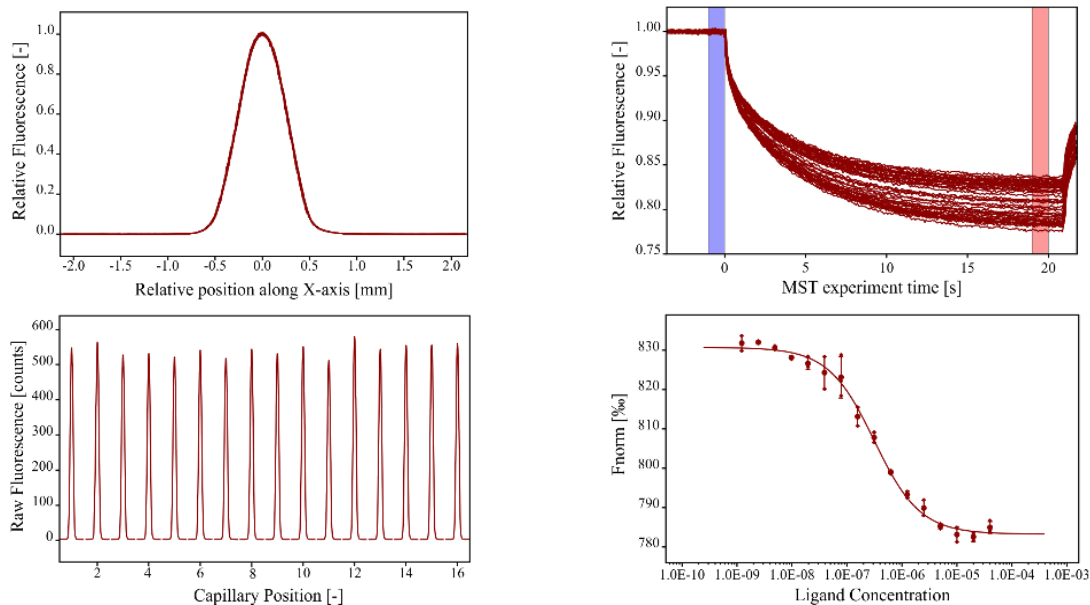
MO.Control v1.6 (NanoTemper Technologies GmbH)  
[nanotempertech.com/monolith-mo-control-software](http://nanotempertech.com/monolith-mo-control-software)

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

50 mM HEPES, pH 7.4, 250 mM NaCl, 0.05% TWEEN® 20  
 50 nM Te\_PtxB | 40 μM – 1.2 nM phosphite | 22°C | medium MST power | 20% excitation power

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

$K_d = 170 \pm 70$  nM [Bisson et al., Nature Communications 2017, 8\(1\), 1–12](#)



## D5. Reference Results/Supporting Results

$K_d = 289 \pm 64$  nM Isothermal Titration Calorimetry (ITC)  
[Bisson et al., Nature Communications 2017, 8\(1\), 1–12](#)

## E. Contributors

Nathan B.P. Adams<sup>2</sup>, Claire Hatty<sup>3</sup>

<sup>2</sup> Department of Molecular Biology and Biotechnology, University of Sheffield, UK

<sup>3</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](http://nanotempertech.com)