

## Monolith Protocol MO-P-056

# ALDH21 – NADP<sup>+</sup>

Aldehyde dehydrogenases constitute a superfamily of NAD(P)<sup>+</sup>-dependent enzymes, which detoxify aldehydes produced in various metabolic pathways. Plants comprise 13 ALDH families, five of them, ALDH11, ALDH19, ALDH21, ALDH22 and ALDH23, unique to plants. ALDH21 from the moss *P. patens* codes for a tetrameric NADP<sup>+</sup>-dependent succinic semialdehyde dehydrogenase (SSALDH), which converts succinic semialdehyde, an intermediate of the c-aminobutyric acid (GABA) shunt pathway, into succinate in the cytosol. Lower plant species including some green algae, non-vascular plants (bryophytes) as well as the oldest vascular plants (lycophods) and ferns (monilophytes) possess a unique aldehyde dehydrogenase (ALDH) gene named ALDH21. However, the gene is absent in flowering plants. The expression of ALDH21 gene is strongly induced by a desiccation and turns back upon rehydration in the twisted mosses *S. ruralis* and *S. caninervis*. Also, the overexpression of ALDH21 in cotton and tobacco improves their drought and salt tolerance.

protein – small molecule interaction | plant biology | ALDH

### A1. Target/Fluorescent Molecule

ALDH21

[uniprot.org/uniprot/A9SS48](http://uniprot.org/uniprot/A9SS48)

### A2. Molecule Class/Organism

Cytosolic enzyme

*Physcomitrella patens*

### A3. Sequence/Formula

MTLGHMVQKA KESSGDVTPK KYNIFLASKP VDGRKWLVD TNKYTNDVAA KVPQATHKDI DDAIDAAVAA APAMAAMGAY  
 ERKAVLEKVV AELKNRFEEI AQLTMESGK PIKDARGEVT RTIDTFQVAA EESVRIYGEH IPLDISARNK GLQGIVKKFP  
 IGPVSMVSPW NFPLNLVAHK VAPAIAVGCP FVLKPARSTP LSALILGEIL HKIEELPLGA FSILPVSRED ADMFTVDERF  
 KLLTFTGSGP IGWDMKARAG KKKVVMELGG NAPCIVDDYV PDLDYTIQRL INGGFYQGGQ SCIHMQRLYV HERLYDEVKE  
 GFVAAVKKLK MGNPFEEPTY LGPMISESAA KGIEDWVKEA VAKGGKLLTG GNRKGAFIEP TVIEDVPIEA NARKEEIFGP  
 VVLLYKYSDF KEAVKECNNT HYGLQSGIFT KDLNKAFYAF EHMEVGGVIL NDSPALRVDS QPYGGLKDSG IQREGVKYAM  
 DDMLETKVLV MRNVGTL

### A4. Purification Strategy/Source

Total RNA was isolated from *Physcomitrella* ('Gransden 2004' strain) at the protonema stage using the RNAqueous kit and plant RNA isolation aid and treated twice with the Turbo DNase-free kit (Thermo Fisher Scientific). The cDNA was synthesized using Superscript III RT (Thermo Fisher Scientific) and oligo dT primers. The sequence coding for PpALDH21 (1494 bp) was amplified using gene-specific primers with SacI and KpnI sites (5'-CAAGAGCTCAATGACGCTCGGCCATAT-3', 5'-ATCGGTACCTACAAGGTCCAACATTCC-3') and the Accuprime Pfx polymerase (Thermo Fisher Scientific). The gene was further cloned into a pCDFDuet vector from Merck and then transformed into T7 express *E. coli* cells (New England Biolabs). Protein production was carried out at 20°C overnight. ALDH21 was purified on a HisPur Cobalt Spin Column (Thermo Fisher Scientific) and further on a HiLoad 26/60 Superdex 200 column (GE Healthcare scientific) into 20 mM Tris-HCl buffer, pH 8.0, 100 mM NaCl and 2% (w/v) glycerol. Protein content was measured using the Coomassie plus (Bradford) protein assay kit with bovine serum albumin (BSA) as a standard (Thermo Fisher Scientific).

### A5. Stock Concentration/Stock Buffer

33 mg/mL | 583  $\mu$ M (calculated for monomer) or 146  $\mu$ M (calculated for the whole tetramer)  
 20 mM Tris-HCl buffer, pH 8.0, 100 mM NaCl, 4% (w/v) glycerol

### A6. Molecular Weight/Extinction Coefficient

55.56 kDa (for 6xHis-tagged monomer)  
 46,090  $\text{M}^{-1}\text{cm}^{-1}$  ( $\epsilon_{280}$ )

### A7. Dilution Buffer

50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM  $\text{MgCl}_2$ , 0.05% TWEEN<sup>®</sup> 20, supplemented with 1 mg/mL BSA

### A8. Labeling Strategy

Monolith Protein Labeling Kit RED – NHS (MO-L001, NanoTemper Technologies GmbH)  
 1\* Labeling Buffer NHS | 1\* 10  $\mu$ g Dye RED-NHS | 1\* B-Column

### A9. Labeling Procedure

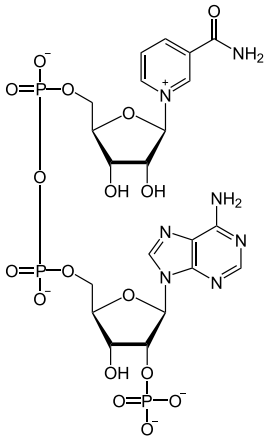
1. Add 95  $\mu$ L of Labeling Buffer NHS to 5  $\mu$ L of 200  $\mu$ M calculated as monomer PpALDH21 to obtain 100  $\mu$ L of a 20  $\mu$ M solution.
2. Add 30  $\mu$ L of DMSO to 10  $\mu$ g Dye RED-NHS to obtain a ~470  $\mu$ M solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
3. Mix 12.5  $\mu$ L of the 470  $\mu$ M dye solution with 87.5  $\mu$ L of Labeling Buffer NHS to obtain 100  $\mu$ L of a 60  $\mu$ M dye solution (3x protein concentration).
4. Mix PpALDH21 and dye in a 1:1 volume ratio (200  $\mu$ L final volume, 6.25% final DMSO concentration).
5. Incubate for 30 minutes at room temperature in the dark.
6. In the meantime, remove the top cap of the B-Column and pour off the storage solution. Remove the bottom cap and place with adapter in a 15 mL tube.
7. Fill the column with dilution buffer and allow it to enter the packed resin bed completely by gravity flow. Discard the flow through collected. Repeat this step 3 more times.
8. Add 200  $\mu$ L of the labeling reaction from step 4 to the center of the column and let sample enter the resin bed completely.
9. Add 300  $\mu$ L of dilution buffer after the sample has entered and discard the flow through.
10. Place column in a new collection tube, add 600  $\mu$ L of dilution buffer and collect the eluate.
11. Keep the labeled PpALDH21 (~3.3  $\mu$ M) on ice in the dark.

### A10. Labeling Efficiency

N/A

## B1. Ligand/Non-Fluorescent Binding Partner

NADP<sup>+</sup> (β-Nicotinamide Adenine Dinucleotide)



## B2. Molecule Class/Organism

Small organic coenzyme

## B3. Sequence/Formula

C<sub>21</sub>H<sub>27</sub>N<sub>7</sub>NaO<sub>17</sub>P<sub>3</sub> (sodium salt)

## B4. Purification Strategy/Source

Applichem Pancreac

[A1394,0001](#)

## B5. Stock Concentration/Stock Buffer

30.6 mg/mL | 40 mM

50 mM Tris-HCl buffer substance at pH 7.5, 150 mM NaCl, 10 mM MgCl<sub>2</sub>

## B6. Molecular Weight/Extinction Coefficient

765.39 Da

## B7. Serial Dilution Preparation

1. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of the 40 mM NADP<sup>+</sup> solution 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. Mix 2.5 μL of labeled PpALDH21 (~3.3 μM) with 197.5 μL of dilution buffer to obtain 200 μL of ~40 nM PpALDH21. Add 10 μL of this solution to each tube from **16** to **1** and mix by pipetting.
4. Incubate tubes for 10 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 v1.5 (NanoTemper Technologies GmbH)

MO.AffinityAnalysis v2.1 (NanoTemper Technologies GmbH)

<https://nanotempertech.com/monolith-mo-control-software/>

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

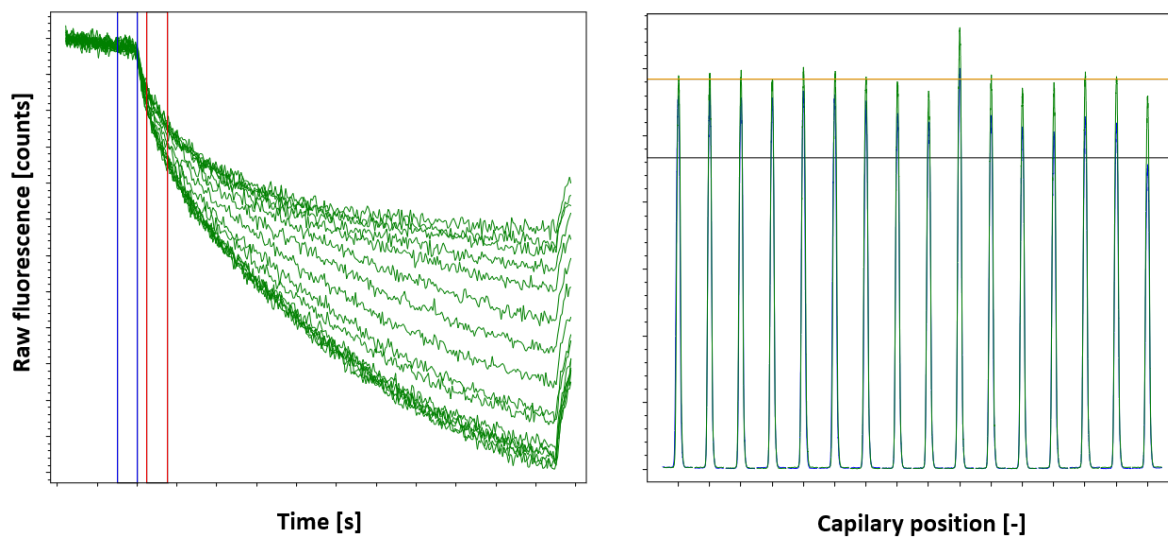
50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM MgCl<sub>2</sub>, 0.05% TWEEN® 20, supplemented with 1 mg/mL BSA  
 20 nM ALDH21 | 20 mM – 0.61 μM NADP<sup>+</sup> | 25°C | high MST power | 20% excitation power

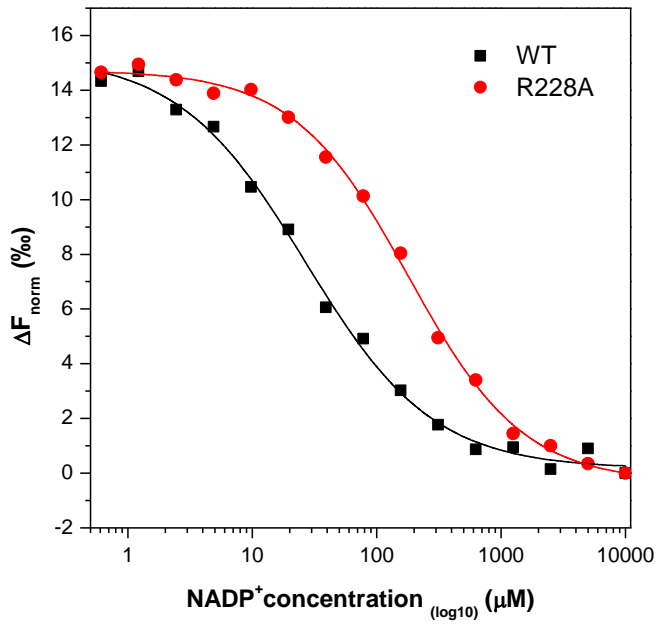
## D4. MST Results (Capillary Scan/Time Traces/Dose Response)

$K_d = 31 \pm 8 \mu\text{M}$  (wt PpALDH21)

$K_d = 282 \pm 50 \mu\text{M}$  (R228A mutant of PpALDH21)

[Kopečný et al. \(2017\) Plant J. 92, 229-243](#)





## D5. Reference Results/Supporting Results

$K_m = 24 \pm 1 \mu\text{M}$  (*wt* PpALDH21)

$K_m = 375 \pm 12 \mu\text{M}$  (R228A mutant of PpALDH21)

Kinetically measured  $K_m$  values for the coenzyme NADP<sup>+</sup>

[Kopečný et al. \(2017\) Plant J. 92, 229-243](#)

## E. Contributors

Dr. David Kopečný<sup>1</sup>, Dr. Radka Končítková<sup>1</sup>, Dr. Jakub Nowak<sup>2</sup>

<sup>1</sup> Palacky University Olomouc, Olomouc, Czech Republic

<sup>2</sup> NanoTemper Technologies GmbH, Krakow, Poland | [nanotempertech.com](http://nanotempertech.com)