

## Monolith Protocol MO-P-058

# Beta-Amylase – Serpin-Z4

Both, amylases and serpins, are highly abundant enzymes in all kingdoms of life. Beta-Amylase is a glycohydrolase that is crucial for seed germination by breaking down starch into smaller carbohydrates that serve as an energy source. Serpins are known to inhibit various serine proteases, but have recently been shown to possess non-inhibitory functions as well.

protein – protein interaction | amylase | serpin | plant proteins | germination

---

### A1. Target/Fluorescent Molecule

Beta-Amylase

[uniprot.org/uniprot/P82993](https://uniprot.org/uniprot/P82993)

### A2. Molecule Class/Organism

Amylases

Barley (*Hordeum vulgare* cv. Harrington)

### A3. Sequence/Formula

MEVNVKGNVY QVYVMLPLDA VSVNNRFEKG DELRAQLRKL VEAGVDGVMV DVWVGLVEGK GPKAYDWSAY KQLFELVQKA  
 GLKLQAIMSF HQCGGNVGDA VNIPIQWVR DVGTRDPDIF YTDGHGTRNI EYLTLGVDNQ PLFHGRSAVQ MYADYMTSFR  
 ENMKEFLDAG VIVDIEVGLG PAGEMRYPSY PQSHGWSFPG IGEFICYDKY LQADFKA AAA AVGHPEW EFP NDAGQYNDTP  
 ERTQFFRDNG TYLTEKGRFF LAWYSNNLIK HGDRILDEAN KVFLGYKVQL AIKISGIHWW YKVPSHAAEL TAGYYNLHDR  
 DGYRTIARML KRHRASINFT CAEMRDSEQS SQAMSAPEEL VQQVLSAGWR EGLNVACENA LPRYDPTAYN TILRNARPHG  
 INQSGPPEHK LFGFTYLRLS NQLVEGQNYV NFKTFVDRMH ANLPRDPYVD PMAPLPRSGP EISIEMILQA AKPKLQPFPP  
 QEHTDLPVGP TGGMGGQAEQ PTCGMGGQVK GPTGGMGGQA EDPTSGMGGE LPATM

### A4. Purification Strategy/Source

Expressed as C-terminal His-tag fusion in *Escherichia coli* SHuffle cells

Purification via Ni-NTA affinity chromatography and gel filtration

### A5. Stock Concentration/Stock Buffer

2.95 mg/mL | 32  $\mu$ M

Phosphate-buffered saline (PBS, pH 7.4), 2 mM DTT, 5% glycerol

### A6. Molecular Weight/Extinction Coefficient

59.6 kDa

92,250  $M^{-1}cm^{-1}$  ( $\epsilon_{280}$ )

### A7. Dilution Buffer

Phosphate-buffered saline (PBS, pH 7.4), 0.1% TWEEN® 20

## A8. Labeling Strategy

Monolith Protein Labeling Kit RED – NHS (MO-L001, NanoTemper Technologies GmbH)

1\* Labeling Buffer NHS | 1\* NT-647-NHS (10 µg) | 1\* B-Column

## A9. Labeling Procedure

1. Add 69 µL of Labeling Buffer NHS to 31 µL of 32 µM Beta-Amylase to obtain 100 µL of a 10 µM solution.
2. Add 30 µL of DMSO to 10 µg NT-647-NHS to obtain a ~470 µM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
3. Mix 6.4 µL of the 470 µM dye solution with 93.6 µL of Labeling Buffer NHS to obtain 100 µL of a 30 µM dye solution (3x protein concentration).
4. Mix Beta-Amylase and dye in a 1:1 volume ratio (200 µL final volume, 3.2% 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 µL of the labeling reaction from step 4 to the center of the column and let sample enter the resin bed completely.
9. Add 400 µL of dilution buffer after the sample has entered and discard the flow through.
10. Place column in a new collection tube, add 500 µL of dilution buffer and collect the eluate.
11. Keep the labeled Beta-Amylase (~2 µM) on ice in the dark.

## A10. Labeling Efficiency

N/A

## B1. Ligand/Non-Fluorescent Binding Partner

Serpin-Z4

[uniprot.org/uniprot/P06293](https://uniprot.org/uniprot/P06293)

## B2. Molecule Class/Organism

Serpins

Barley (*Hordeum vulgare* cv. Harrington)

## B3. Sequence/Formula

MATTLATDVR LSIHQTRFA LRLRSAISSN PERAAGNVAF SPLSLHVALS LITAGAAATR DQLVAILGDG GAGDAKELNA  
 LAEQVVQFVL ANESSTGGPR IAFANGIFVD ASLSLKPSFE ELAVCQYKAK TQSVDFQHKT LEAVGQVNSW VEQVTTGLIK  
 QILPPGSVDN TTKLILGNAL YFKGAWDQKF DESNTKCDSF HLLDGSSIQT QFMSSTKKQY ISSSDNLKVL KLPYAKGHDK  
 RQFSMYILLP GAQDGLWLSLA KRLSTEPEFI ENHIPKQTV E VGRFQLPKFK ISYQFEASSL LRALGLQLPF SEEADLSEMV  
 DSSQGLEISH VFHKSFEVEN EEGTEAGAAT VAMGVAMSMP LKVDLVDFVA NHPFLFLIRE DIAGVVVFG HVTNPLISA

#### B4. Purification Strategy/Source

Expressed as C-terminal His-tag fusion in *Escherichia coli* SHuffle cells  
Purification via Ni-NTA affinity chromatography and gel filtration

#### B5. Stock Concentration/Stock Buffer

2.06 mg/mL | 80  $\mu$ M  
Phosphate-buffered saline (PBS, pH 7.4), 2 mM DTT, 5 % glycerol

#### B6. Molecular Weight/Extinction Coefficient

42.3 kDa  
25,440  $M^{-1}cm^{-1}$  ( $\epsilon_{280}$ )

#### B7. Serial Dilution Preparation

1. Prepare a PCR-rack with 16 PCR tubes. Transfer 20  $\mu$ L of the 80  $\mu$ M Serpin-Z4 solution into tube **1**. Then, transfer 10  $\mu$ L of dilution buffer into tubes **2** to **16**.
2. Prepare a 1:1 serial dilution by transferring 10  $\mu$ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10  $\mu$ L from tube **16** to get an equal volume of 10  $\mu$ L for all samples.
3. Mix 7  $\mu$ L of labeled Beta-Amylase with 193  $\mu$ L of dilution buffer to obtain 200  $\mu$ L of  $\sim$ 70 nM Beta-Amylase.
4. Add 10  $\mu$ L of 70 nM labeled Beta-Amylase to each tube from **16** to **1** and mix by pipetting.
5. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

#### D1. MST System/Capillaries

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

#### D2. MST Software

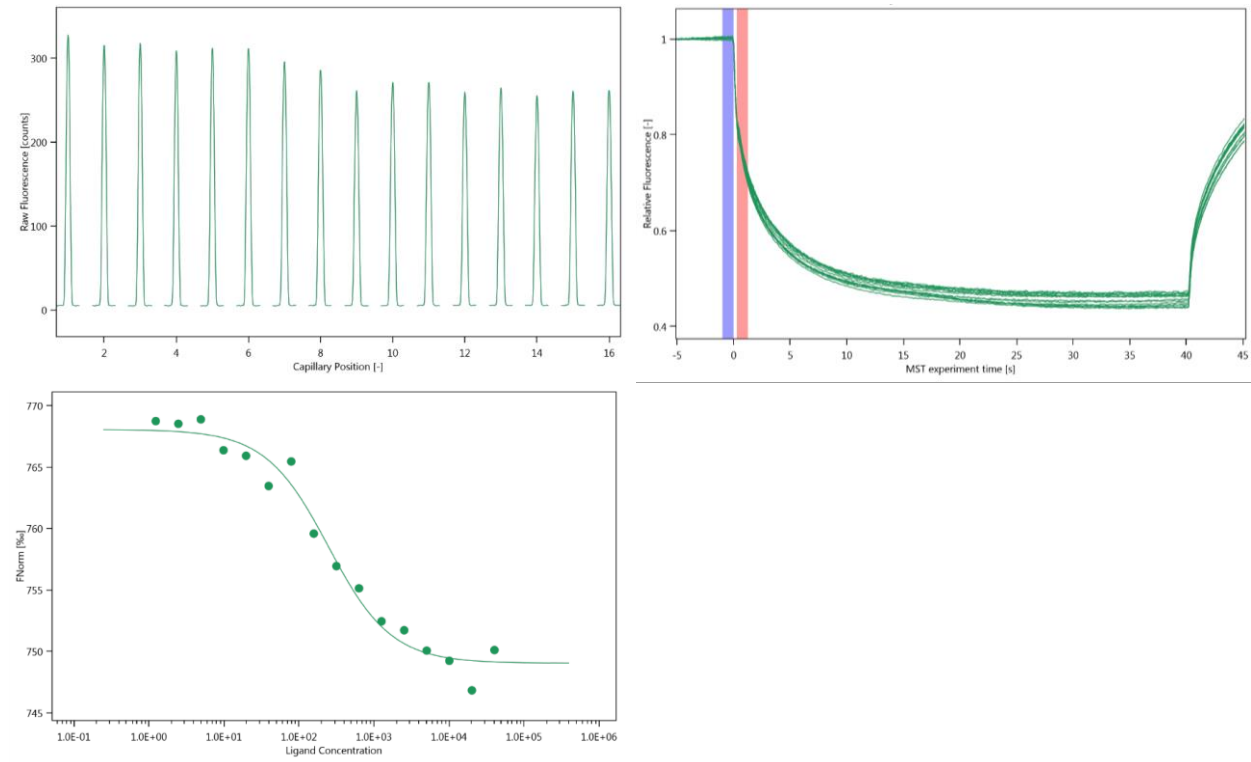
MO.Control v1.6 (NanoTemper Technologies GmbH)  
[nanotempertech.com/monolith-mo-control-software](https://nanotempertech.com/monolith-mo-control-software)

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

Phosphate-buffered saline (PBS, pH 7.4), 0.1% TWEEN<sup>®</sup> 20  
35 nM Beta-Amylase | 40.5  $\mu$ M – 1.24 nM Serpin-Z4 | 25°C | low MST Power | 50% excitation power

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

$K_d = 215 \pm 14$  nM



#### D5. Reference Results/Supporting Results

$K_d = 105$  nM      MicroScale Thermophoresis (MST) – inverted assay

$K_d = 316$  nM      Surface Plasmon Resonance (SPR)

$K_d = 142$  nM      Surface Plasmon Resonance (SPR) – inverted assay

[Cohen and Fluhr, Plant Direct \(2018\) 1–12](#)

#### E. Contributors

Maja Cohen<sup>1</sup>, Robert Fluhr<sup>1</sup>, Timm Hassemer<sup>2</sup>

<sup>1</sup> Department of Plant Sciences, The Weizmann Institute of Science, Rehovot, Israel

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