

Monolith Protocol MO-P-067

# Bovine Serum Albumin – Zinc

Bovine serum albumin (also known as BSA) is a serum albumin protein derived from cows, which is often used as a protein concentration standard in lab experiments. Among its 583 amino acid residues, it contains 17 intrachain disulfide bonds as well as one free thiol group, which can be used for site-specific labeling with a thiol-reactive (maleimide) fluorescent dye. Apart from fatty acids and a variety of small molecules in the blood, BSA is also associated with the transport of metal ions such as zinc ( $Zn^{2+}$ ).

protein – ion interaction | Zn

---

## A1. Target/Fluorescent Molecule

Bovine serum albumin (BSA)

[uniprot.org/uniprot/P02769](https://uniprot.org/uniprot/P02769)

## A2. Molecule Class/Organism

Serum protein

*Bos taurus* (Bovine)

## A3. Sequence/Formula

DTHKSEIAHR FKDLGEEHFK GLVLIAFSQY LQQCPFDEHV KLVNELTEFA KTCVADESHA GCEKSLHTLF GDELCKVASL  
 RETYGDMA DC CEKQEPERNE CFLSHKDDSP DLPKLPDPN TLCDEFKADE KKF<sup>W</sup>GKLYE IARRHPYFYA PELLYYANKY  
 NGVFQECCQA EDKGACLLPK IETMREKVLA SSARQLRCA SIQKFGERAL KAW<sup>S</sup>VARLSQ KFPKAEFVEV TKLVTDLTKV  
 HKECCHGDL L ECADDRADLA KYICDNQDTI SSKLKECCDK PLLEKSHCIA EVEKDAIPEN LPPLTADFAE DKDVCKNYQE  
 AKDAFLGSFL YEYSRRHPEY AVSVLLRLAK EYEATLEEC AKDDPHACYS TVFDK<sup>L</sup>KHLV DEPQNLIKQN CDQFEKLG EY  
 GFQNALIVRY TRKVPQVSTP TLVEVSRSLG KVGTRCCTKP ESERPCTED YLSLILNRLC VLHEKTPVSE KVTKCCTESL  
 VNRRPCFSAL TPDETYVPKA FDEKLFTFHA DICTLPDTEK QIKKQTALVE LLKHKPKATE EQLKTVMENF VAFVDKCCAA  
 DDKEACFAVE GPKL<sup>V</sup>STQT ALA

## A4. Purification Strategy/Source

Carl Roth GmbH

[8076.2](#)

## A5. Stock Concentration/Stock Buffer

Powdered

## A6. Molecular Weight/Extinction Coefficient

66.5 kDa

$43,800 \text{ M}^{-1}\text{cm}^{-1}$  ( $\epsilon_{280}$ )

## A7. Dilution Buffer

Phosphate buffered saline (PBS, pH 7.4)<sup>1</sup>

## A8. Labeling Strategy

Monolith Protein Labeling Kit RED – MALEIMIDE 2nd Generation (MO-L014, NanoTemper Technologies GmbH)  
 1\* Labeling Buffer MALEIMIDE | 1\* Dye RED-MALEIMIDE 2nd Generation (10 µg) | 1\* B-Column

## A9. Labeling Procedure

1. Add 3 mL of dilution buffer to 20 mg of BSA to obtain a 100 µM solution. Mix carefully with a pipette to dissolve all protein and avoid creating air bubbles.
2. Mix 20 µL of 100 µM BSA with 180 µL of Labeling Buffer MALEIMIDE to obtain 200 µL of a 10 µM BSA solution.
3. Mix 2 µL of 500 mM TCEP with 198 µL of Labeling Buffer MALEIMIDE to obtain 200 µL of 5 mM TCEP.
4. Mix 3.2 µL of 5 mM TCEP with 196.8 µL of Labeling Buffer MALEIMIDE to obtain 200 µL of 80 µM TCEP.
5. Mix 50 µL of the 80 µM TCEP solution with 100 µL of the 10 µM BSA solution to obtain 150 µL of a 6.7 µM BSA, 26.7 µM TCEP solution (4x protein concentration).
6. Incubate for 30 minutes at room temperature.
7. Add 13.2 µL of DMSO to Dye RED-MALEIMIDE 2nd Generation (10 µg) to obtain a ~1 mM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
8. Mix 10 µL of the 1 mM dye solution with 40 µL of Labeling Buffer MALEIMIDE to obtain 50 µL of a 200 µM dye solution.
9. Add the 150 µL of the BSA-TCEP mix (step 5) to the 200 µM dye solution to obtain 200 µL of a 5 µM BSA, 20 µM TCEP, 50 µM dye solution (10x protein concentration, 2.5x TCEP concentration<sup>2</sup>, 5% DMSO).
10. Incubate for 30 minutes at room temperature in the dark.
11. 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.
12. 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.
13. Add 200 µL of the labeling reaction from step 9 to the center of the column and let sample enter the bed completely.
14. Add 400 µL of dilution buffer after the sample has entered and discard the flow through.
15. Place column in a new collection tube, add 500 µL of dilution buffer and collect the eluate.
16. Keep the labeled BSA (~2 µM) on ice in the dark.

## A10. Labeling Efficiency

Measurement of protein concentration and degree of labeling (DOL) using a NanoDrop™:

[nanotempertech.com/dol-calculator](https://nanotempertech.com/dol-calculator)

Absorbance A <sub>280</sub>	0.082	Protein concentration	1.76 µM
Absorbance A <sub>650</sub>	0.125	Degree-of-labeling (DOL)	0.36

<sup>1</sup> Detergents (e.g. TWEEN® 20) cannot be used as they bind to a similar binding site as Zn<sup>2+</sup> (Coverdale et al., *Metallomics*, 2019, 11, 1805).

<sup>2</sup> As maleimide can also react with TCEP, it is important to have an excess of maleimide over TCEP in the labeling reaction.

## B1. Ligand/Non-Fluorescent Binding Partner

Zn<sup>2+</sup>

## B2. Molecule Class/Organism

Ion

## B3. Sequence/Formula

ZnCl<sub>2</sub>

## B4. Purification Strategy/Source

Carl Roth GmbH

[T887.1](#)

## B5. Stock Concentration/Stock Buffer

Powdered

## B6. Molecular Weight/Extinction Coefficient

136.28 Da

## B7. Serial Dilution Preparation

1. Dissolve 0.34 g of ZnCl<sub>2</sub> in 10 mL of ddH<sub>2</sub>O to obtain a 250 mM ZnCl<sub>2</sub> stock solution. Add 4 drops of HCl (1.2 M) to acidify the solution and completely dissolve all ZnCl<sub>2</sub>.
2. Mix 2 μL of the 250 mM ZnCl<sub>2</sub> stock with 998 μL of ddH<sub>2</sub>O to obtain a 1 mL of a 500 μM ZnCl<sub>2</sub> solution<sup>3</sup>.
3. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of the 500 μM ZnCl<sub>2</sub> solution into tube **1**. Then, add 10 μL of ddH<sub>2</sub>O to tubes **1** 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. Mix 8 μL of labeled BSA (~2 μM) and 192 μL of dilution buffer containing 0.005% Pluronic® F-127 to obtain 200 μL of ~80 nM labeled BSA solution.
6. Add 10 μL of this solution to each tube from **16** to **1** and mix by pipetting.
7. Load capillaries immediately<sup>4</sup>.

---

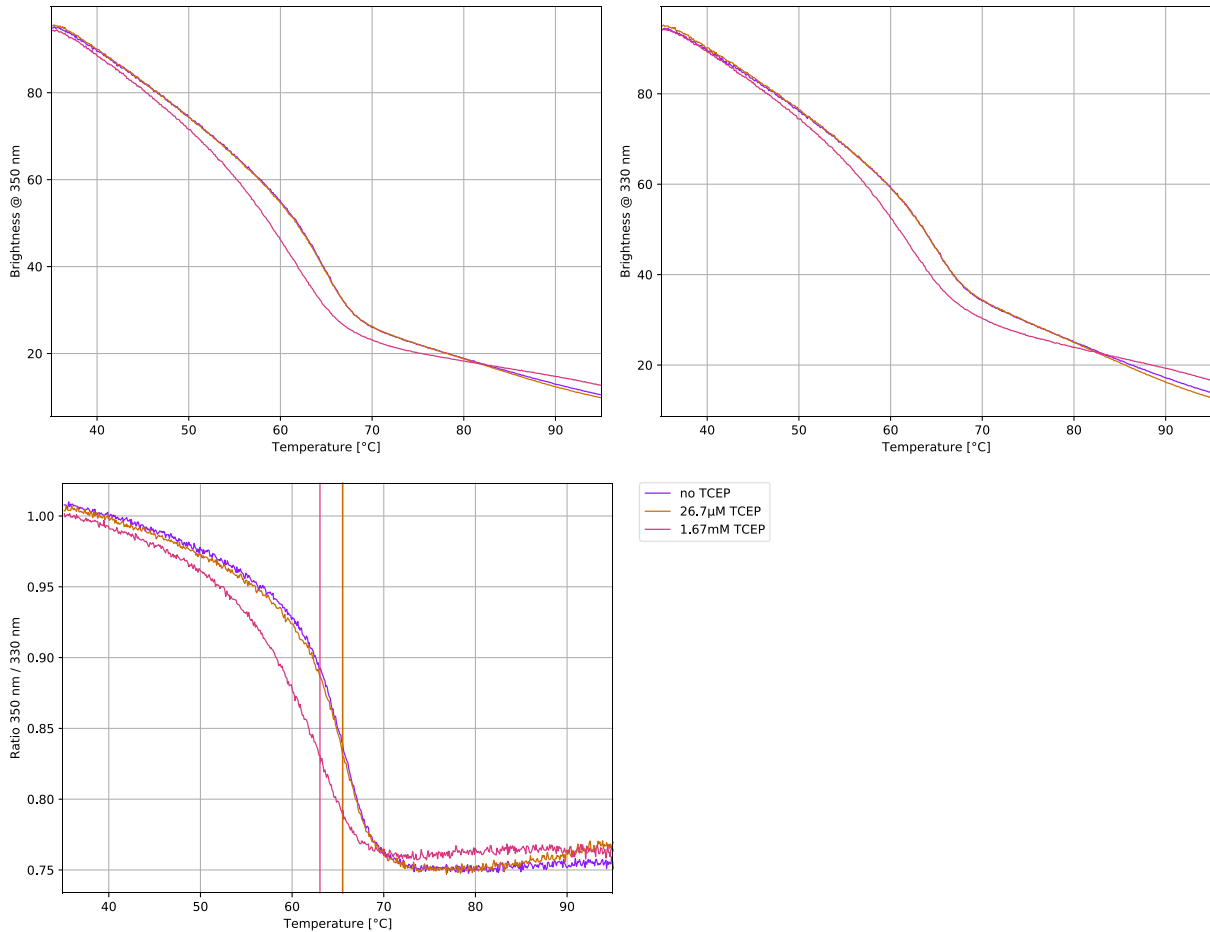
<sup>3</sup> Due to the low solubility of ZnCl<sub>2</sub> at non-acidic pH and higher ionic strength, as well as its tendency to form insoluble Zn(OH)<sub>2</sub> at pH > 7, the dilution series is prepared in ddH<sub>2</sub>O (see also [Krezel et al., Archives of Biochemistry and Biophysics 611 \(2016\) 3-19](#)).

<sup>4</sup> Longer incubation times can lead to BSA aggregation at high ZnCl<sub>2</sub> concentrations, as well as precipitation of Zn<sub>2+</sub> in the phosphate buffer.

Validation of TCEP activity using Tycho NT.6:

[nonotempertech.com/tycho](http://nonotempertech.com/tycho)

no TCEP	10 $\mu$ L of 10 $\mu$ M BSA + 5 $\mu$ L of Labeling Buffer MALEIMIDE	$T_i = 65.5^\circ\text{C}$
26.7 $\mu$ M TCEP	10 $\mu$ L of 10 $\mu$ M BSA + 5 $\mu$ L of 80 $\mu$ M TCEP	$T_i = 65.5^\circ\text{C}$
1.67mM TCEP	10 $\mu$ L of 10 $\mu$ M BSA + 5 $\mu$ L of 5 mM TCEP	$T_i = 63.0^\circ\text{C}$ <sup>5</sup>

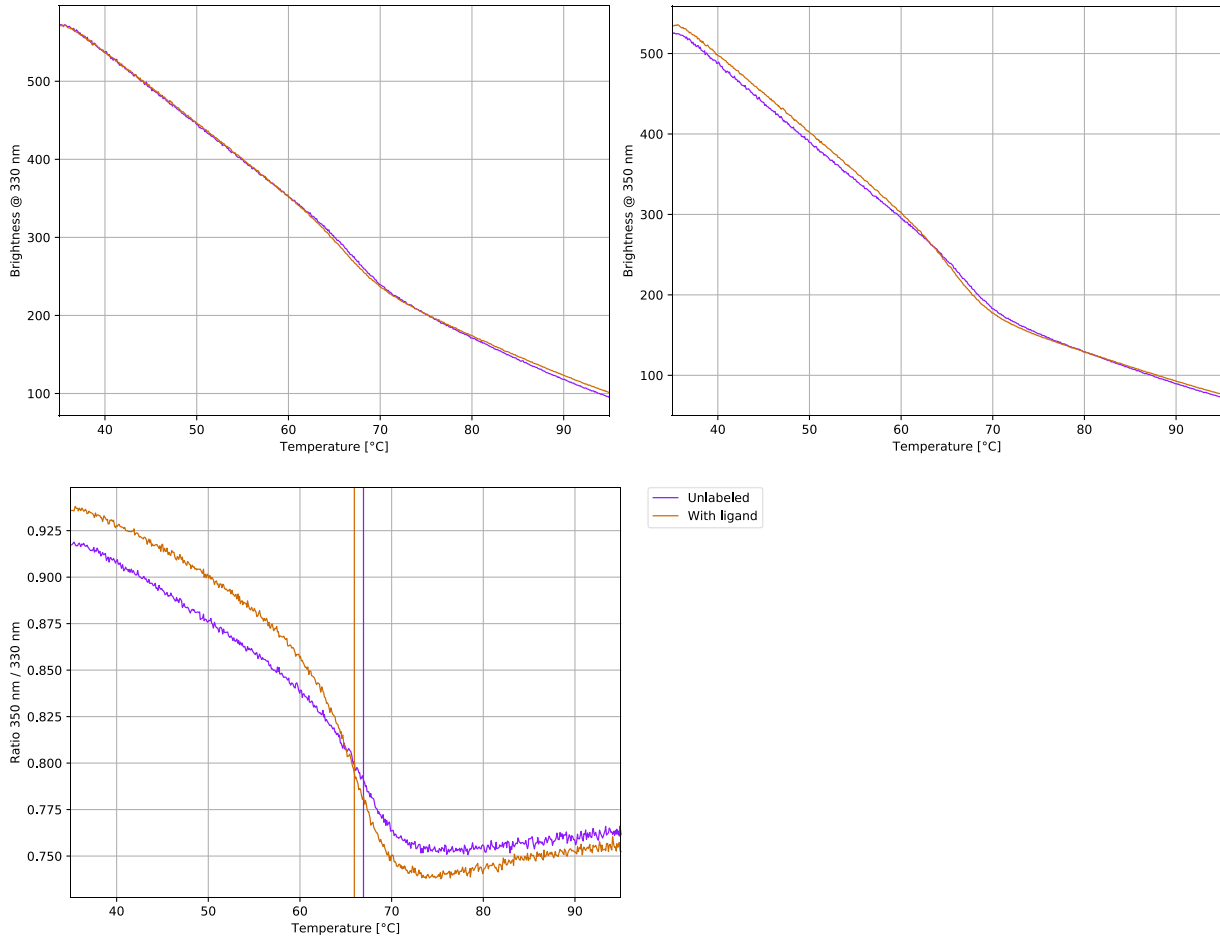


<sup>5</sup> While 26.7  $\mu$ M TCEP has no effect on the structural integrity of BSA, 1.67 mM TCEP significantly destabilizes the protein, presumably by the reduction of intrachain disulfide bonds.

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

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

Unlabeled	10 $\mu$ L of ddH <sub>2</sub> O + 10 $\mu$ L of 100 $\mu$ M BSA	T <sub>i</sub> = 66.9°C
With ligand	8 $\mu$ L of ddH <sub>2</sub> O + 2 $\mu$ L of 500 $\mu$ M ZnCl <sub>2</sub> + 10 $\mu$ L of 100 $\mu$ M BSA	T <sub>i</sub> = 65.9°C



**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)

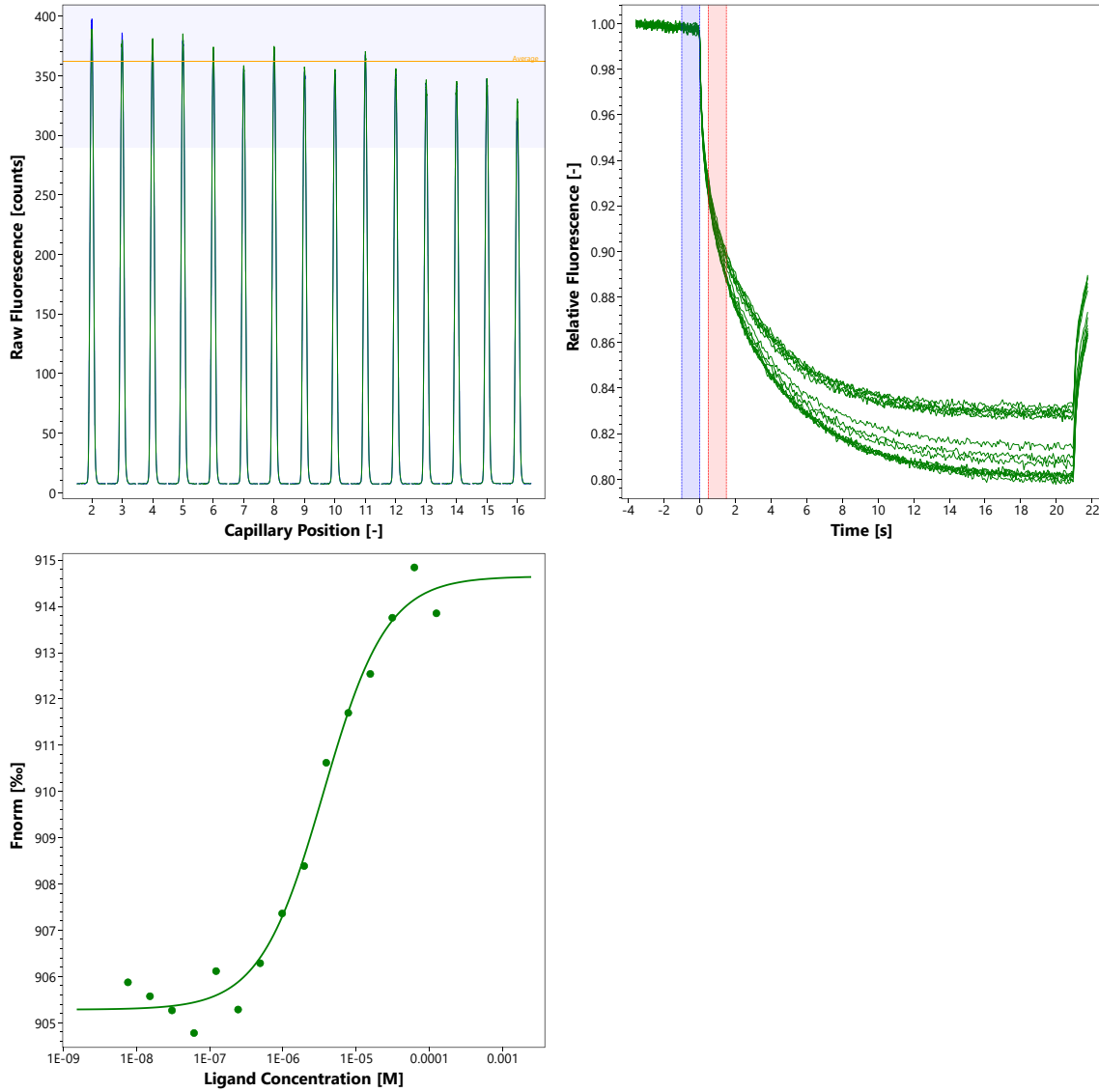
[nanotempertech.com/monolith-mo-control-software](http://nanotempertech.com/monolith-mo-control-software)

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

0.5X Phosphate buffered saline (PBS, pH 7.4), 0.0025% Pluronic® F-127<sup>6</sup>  
 40 nM BSA | 250 μM – 7.6 nM ZnCl<sub>2</sub> | 23°C | high MST power | 40% excitation power

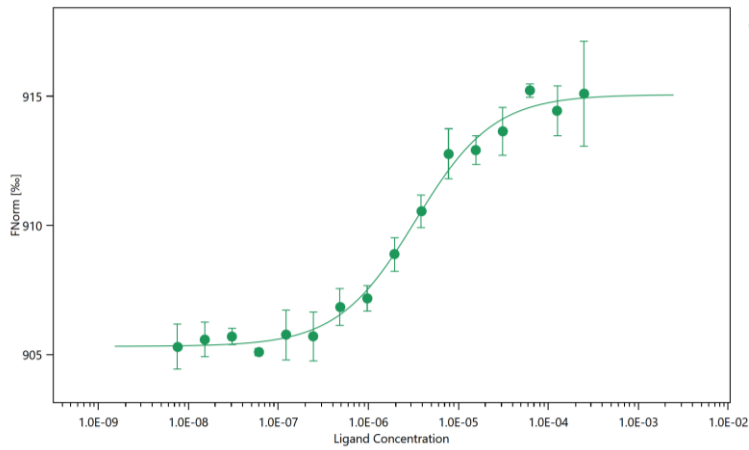
### D4. MST Results (Capillary Scan/Time Traces/Dose Response)

$K_d = 3.5 \mu\text{M}$



<sup>6</sup> Higher concentrations of detergent interfere with Zn<sup>2+</sup> binding due to the shared binding sites.

$K_d = 3.3 \pm 0.4 \mu\text{M}$  (n = 3)



### D5. Reference Results/Supporting Results

$K_d = 0.3 \mu\text{M}$  Competitive spectrophotometry  
 Oyashi et al., *Journal of Inorganic Biochemistry* 75 (1999) 213–218

### E. Contributors

Andreas Langer<sup>7</sup>

<sup>7</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](http://nanotempertech.com)