

Monolith Protocol MO-P-018

Bovine Serum Albumin – Ibuprofen

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. Ibuprofen is a nonsteroidal anti-inflammatory drug that is used for treating pain, fever, and inflammation. It binds to a hydrophobic pocket of BSA (Sudlow site II) which is the principal binding site for drugs.

protein – small molecule interaction | maleimide

A1. Target/Fluorescent Molecule

Bovine serum albumin (BSA)

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^WGKYLIE IARRHPYFYA PELLYYANKY
 NGVFQECCQA EDKGACLLPK IETMREKVL A SSARQLRCA SIQKFGERAL KAW^SVARLSQ KFPKAEFVEV TKLVTDLTKV
 HKECCHGDL L ECADDRADLA KYICDNQDTI SSKLKECCDK PLLEKSHCIA EVEKDAIPEN LPPLTADFAE DKDVCKNYQE
 AKDAFLGSFL YEYSRRHPEY AVSVLLRLAK EYEATLEECC AKDDPHACYS TVFDK^LKHLV DEPQNLIKQN CDQFEKLG EY
 GFQNALIVRY TRKVPQVSTP TLVEVSRSLG KVGTRCCTK P ESERMPCTED YLSLILNRLC VLHEKTPVSE KVTKCCTESL
 VNRRPCFSAL TPDETYVPKA FDEKLFTFHA DICTLPDTEK QIKKQTALVE LLKHKPKATE EQLKTVMENF VAFVDKCCAA
 DDKEACFAVE GPKL^VSTQT 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 M⁻¹cm⁻¹ (ε₂₈₀)

A7. Dilution Buffer

Phosphate buffered saline (PBS, pH 7.4)¹

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 concentration², 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

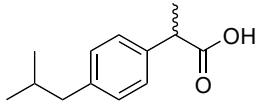
Absorbance A ₂₈₀	0.082	Protein concentration	1.76 µM
Absorbance A ₆₅₀	0.125	Degree-of-labeling (DOL)	0.36

¹ Detergents (e.g. TWEEN® 20) cannot be used, as they also bind to BSA and would compete with ibuprofen binding.

² 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

Ibuprofen



B2. Molecule Class/Organism

Non-steroidal anti-inflammatory drug

B3. Sequence/Formula

$C_{13}H_{18}O_2$

B4. Purification Strategy/Source

Sigma-Aldrich GmbH

1110

B5. Stock Concentration/Stock Buffer

Powdered

B6. Molecular Weight/Extinction Coefficient

206.28 Da

$346 \text{ M}^{-1}\text{cm}^{-1}$ (ϵ_{272})

B7. Serial Dilution Preparation

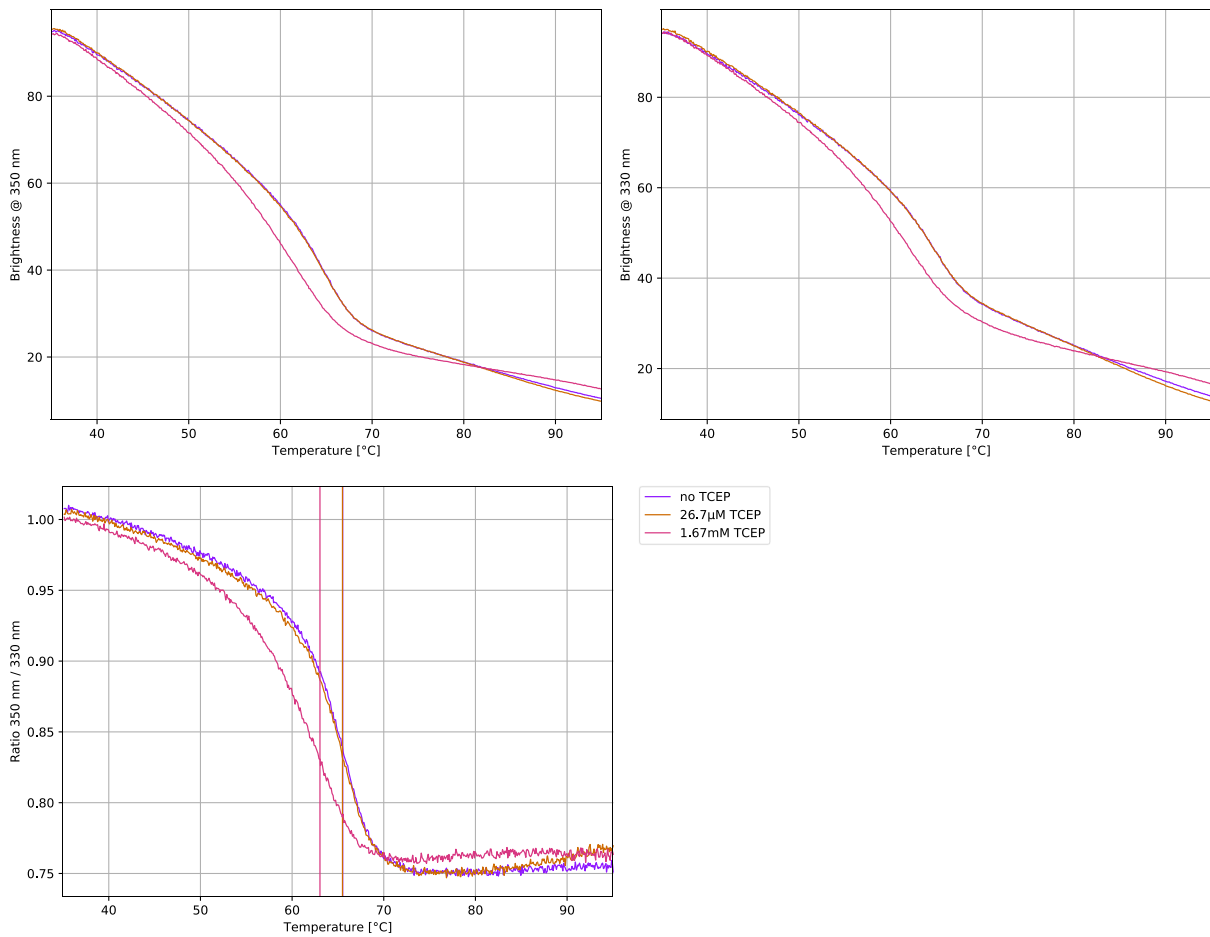
1. Dissolve 20.6 mg of ibuprofen in 1 mL of DMSO to obtain a 100 mM ibuprofen stock solution.
2. Mix 2 μL of the 100 mM ibuprofen stock with 98 μL of dilution buffer to obtain 100 μL of a 2 mM ibuprofen solution.
3. Mix 4 μL of DMSO with 196 μL of dilution buffer to obtain 200 μL of an 2% DMSO solution.
4. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μL of the 2 mM ibuprofen solution into tube **1**. Then, transfer 10 μL of the 2% DMSO solution into tubes **2** to **16**.
5. 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.
6. Mix 2 μL of unlabeled BSA (100 μM) with 2 μL of labeled BSA ($\sim 2 \mu\text{M}$) and 196 μL of dilution buffer to obtain 200 μL of a 1 μM unlabeled BSA, $\sim 20 \text{ nM}$ labeled BSA solution.
7. Add 10 μL of this solution to each tube from **16** to **1** and mix by pipetting.
8. Incubate for 5 minutes at room temperature in the dark before loading capillaries.

C. Applied Quality Checks

Validation of TCEP activity using Tycho NT.6:

nanotempertech.com/tycho

no TCEP	10 μ L of 10 μ M BSA + 5 μ L of Labeling Buffer MALEIMIDE	$T_i = 65.5^\circ\text{C}$
26.7 μ M TCEP	10 μ L of 10 μ M BSA + 5 μ L of 80 μ M TCEP	$T_i = 65.5^\circ\text{C}$
1.67mM TCEP	10 μ L of 10 μ M BSA + 5 μ L of 5 mM TCEP	$T_i = 63.0^\circ\text{C}^3$

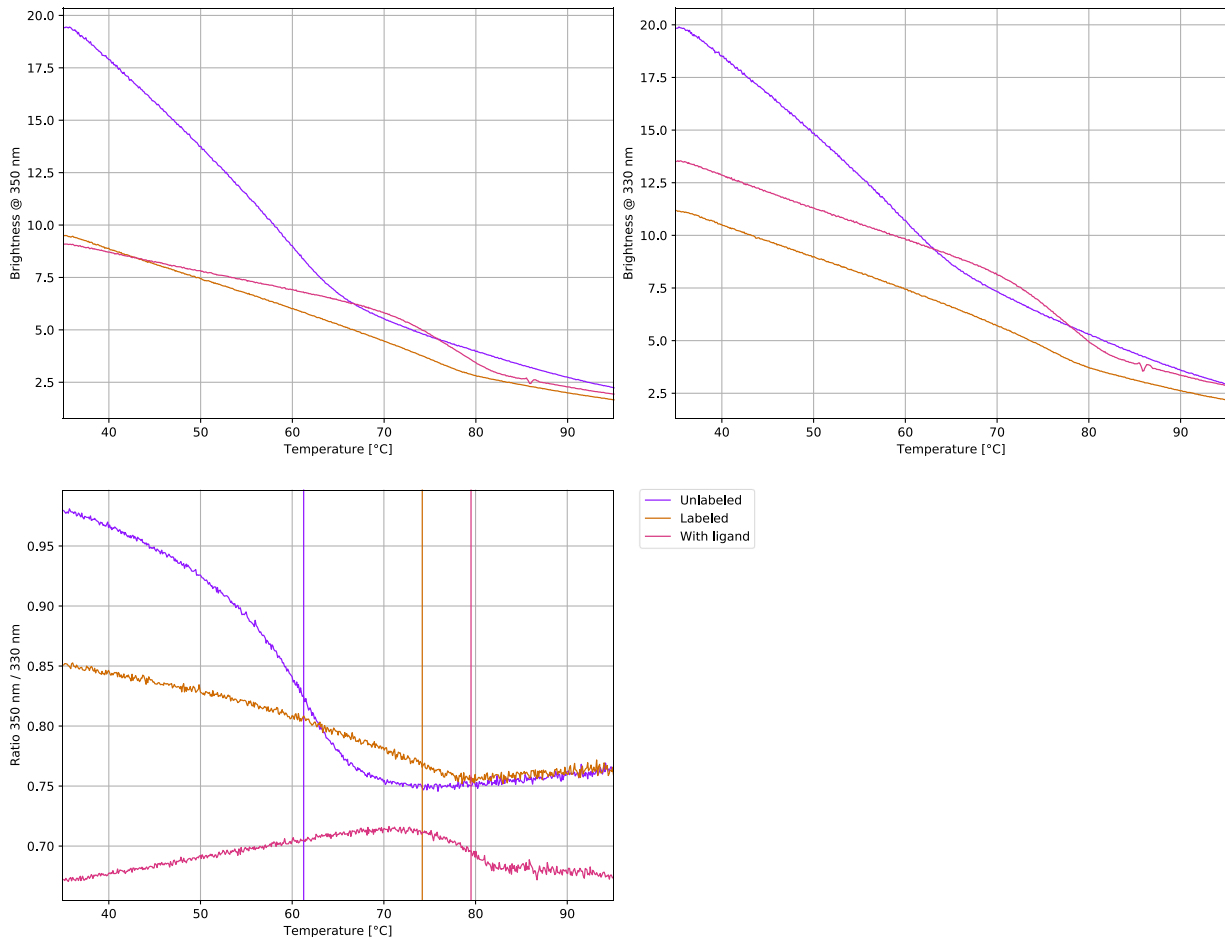


³ While 26.7 μ M TCEP has no effect on the structural integrity of BSA, 1.67 mM TCEP significantly destabilizes the protein, presumably by the reduction of intrachain disulfide bonds.

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

nanotempertech.com/tycho

Unlabeled	1.6 μ L of 100 μ M BSA + 98.4 μ L of dilution buffer	$T_i = 61.3^\circ\text{C}$
Labeled	8 μ L of B-column eluate (~ 2 μ M BSA) + 2 μ L of 2% DMSO	$T_i = 74.1^\circ\text{C}$
With ligand	8 μ L of B-column eluate (~ 2 μ M BSA) + 2 μ L of 2 mM ibuprofen	$T_i = 79.5^\circ\text{C}$



D1. MST System/Capillaries

Monolith NT.115^{Pico} 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

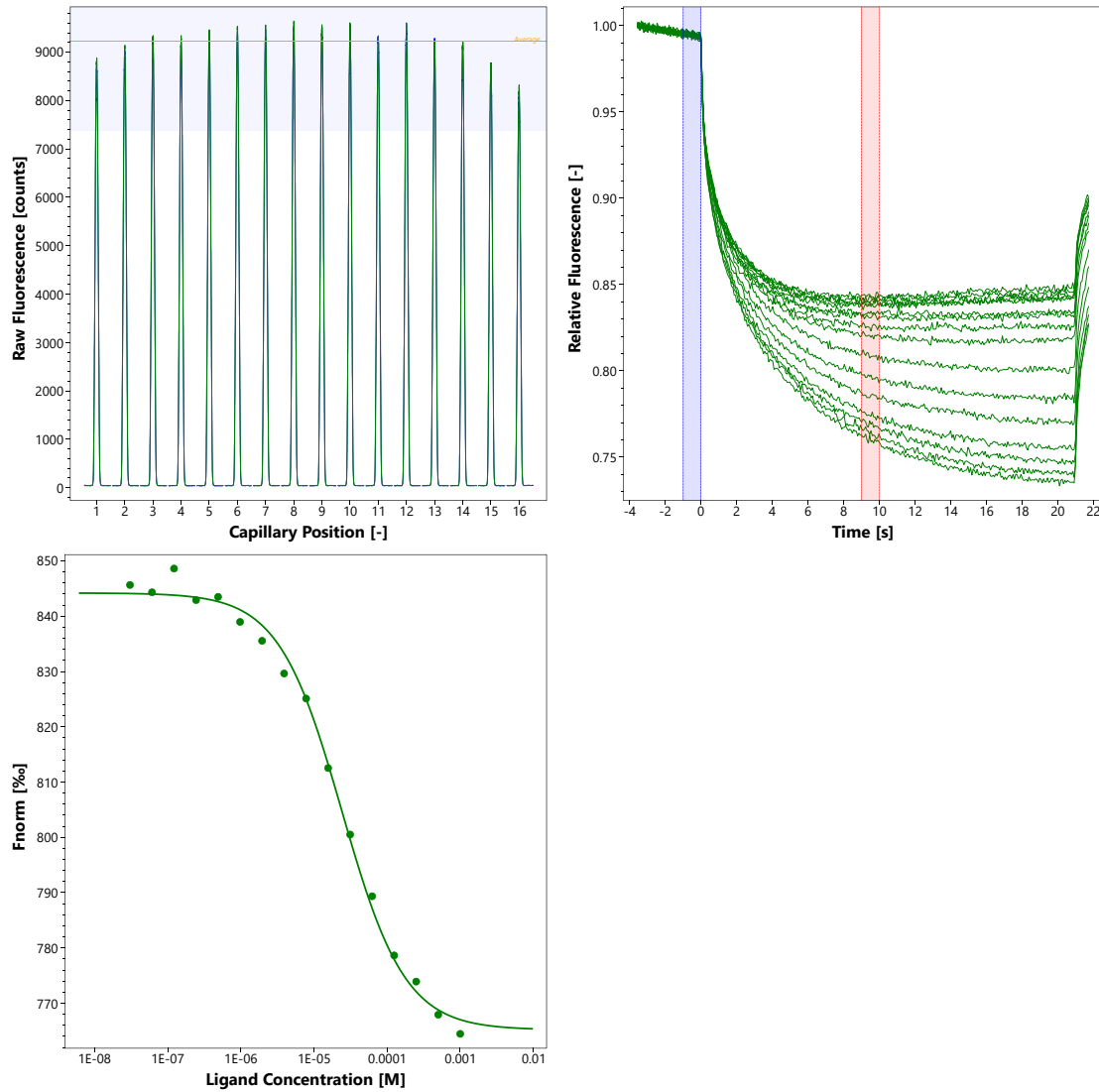
D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

Phosphate buffered saline (PBS, pH 7.4), 1% DMSO

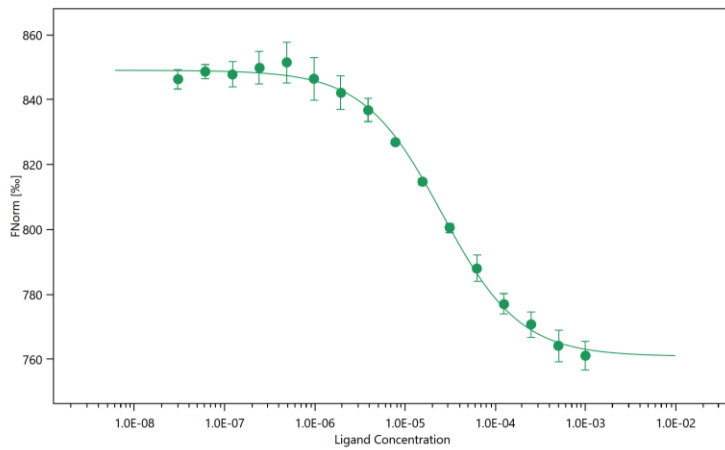
0.5 μ M (unlabeled), 10 nM (labeled) BSA | 1 mM – 30 nM ibuprofen | 25°C | high MST power | 20% excitation power

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

$K_d = 23.9 \mu\text{M}$



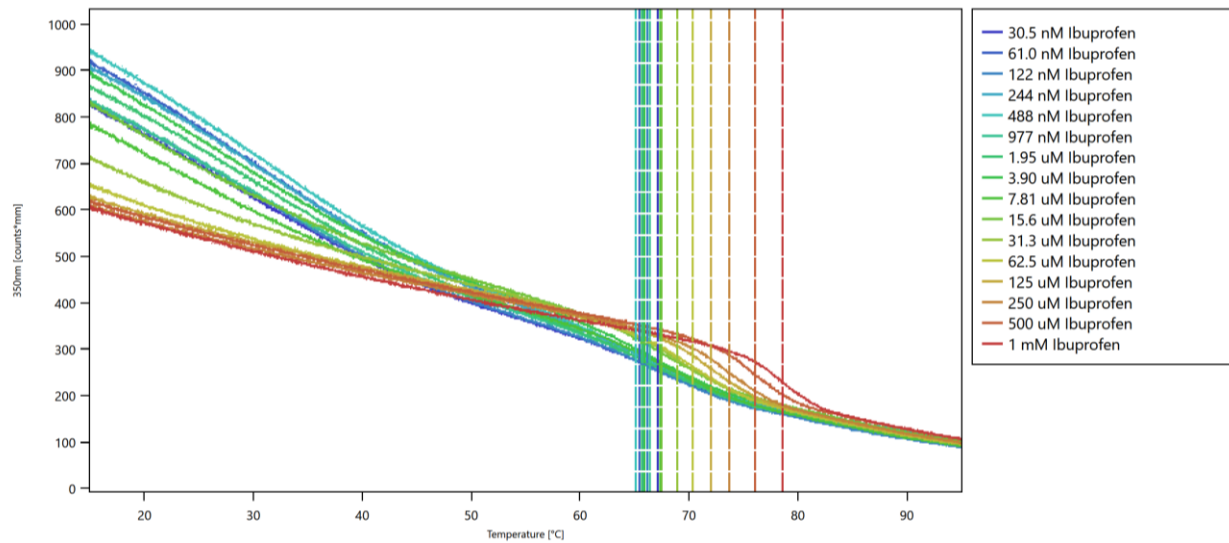
$K_d = 25.3 \pm 1.7 \mu\text{M}$ ($n = 4$)



D5. Reference Results/Supporting Results

Confirmation of ibuprofen binding to BSA by a Trp fluorescence quenching & thermal shift measurement:

nanotempertech.com/prometheus



E. Contributors

Andreas Langer⁴

⁴ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com