

Monolith Protocol MO-P-049

B. burgdorferi CspA – C9

Borrelia burgdorferi is a Gram-negative bacterium that causes Lyme disease. These spirochetes are spread by ticks and if the infection is left untreated, severe and chronic symptoms develop, including neuroborreliosis, carditis and arthritis. To be able to survive in the human host, this bacterium has evolved several sophisticated mechanisms to evade the human innate immune system. One strategy is to bind complement proteins and thus inhibit the complement cascade directly. CspA is a surface exposed protein that binds the complement protein C9 and thereby blocks the complement attack and contributes to bacterial survival in the human host.

protein – protein interaction | complement evasion

A1. Target/Fluorescent Molecule

CspA

uniprot.org/uniprot/A0A0H3BZN5

A2. Molecule Class/Organism

Bacterial surface protein/immune evasion protein

Borrelia burgdorferi

A3. Sequence/Formula

MTKAKLNIIK LNIIAMILTL ICTSCAPFSK IDPKANANTK PPKITNPGEN TQNFEDKSGD LSASDEKIME TIASELKAIG
 KELEDQKKEE NIQIAKIAKE KDFDLSTFKV GPYDLIDEDI QMKIKRTLVS SLDYKKNIE KLKEILEILK KNSKHNYIIG
 RLIYHISW^SSI QFQIEQNLEL IQNGVENLSQ EESKSLMQI KSNLEIKQRL KKTLNETLKV YNQNTQDNEK ILAEHFNKYY
 KDFDTLKPAF Y

A4. Purification Strategy/Source

Construction and expression of plasmids harboring the *cspA* gene encoding for CspA and purification of the proteins were previously described.¹

A5. Stock Concentration/Stock Buffer

8.3 mg/mL | 320 μM

A6. Molecular Weight/Extinction Coefficient

25.9 kDa

¹ Kraiczky et al., J Biol Chem 279 (2004) 2421-2429

A7. Dilution Buffer

50 mM Tris-HCl, pH 7.6, 150 mM NaCl, 10 mM MgCl₂, 0.05% TWEEN® 20

A8. Labeling Strategy

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

1* Labeling Buffer NHS | 1* 10 µg RED-NHS dye | 1* B-Column

A9. Labeling Procedure

1. Add 96.9 µL of Labeling Buffer NHS to 3.1 µL of 320 µM CspA to obtain 100 µL of a 10 µM solution.
2. Add 30 µL of DMSO to 10 µg RED-NHS dye 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 CspA 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 6 to the center of the column and let sample enter the bed completely.
9. Add 300 µL of dilution buffer after the sample has entered and discard the flow through.
10. Place column in a new collection tube, add 600 µL of dilution buffer and collect the eluate.
11. Keep the labeled CspA (~1.6 µM) on ice in the dark.

A10. Labeling Efficiency

N/A

B1. Ligand/Non-Fluorescent Binding Partner

Complement component 9 (C9)

uniprot.org/uniprot/P02748

B2. Molecule Class/Organism

Human complement protein

Homo sapiens (Human)

B3. Sequence/Formula

MSACRSFAVA ICILEISILT AQYTTSYDPE LTESSGSASH IDCRCMPWSE WSQCDPCLRQ MFRSRSIEVF GQFNGKRCTD
 AVGDRRQCVP TEPCEDAEDD CGNDFQCSTG RCIKMRLRCN GDNDCGDFSD EDDCESEPRP PCRDRVVEES ELARTAGYGI
 NILGMDPLST PFDNEFYNGL CNRDRDGNTL TYYRRPWVVA SLIYETKGEK NFRTEHYEEQ IEAFKSIIQE KTSNFNAAIS
 LKFTPTETNK AEQCCEETAS SISLHGKGSF RFSYSKNETY QLFLSYSSKK EKMFLHVKGE IHLGRFVMRN RDVVLTTTFV
 DDIKALPTY EKGEYFAFLE TYGTHYSSSG SLGGLYELIY VLDKASMKRK GVELKDIKRC LGYHLDVSLA FSEISVGAEF
 NKDDCVKRG EGRAVNITSEN LIDDVVSLIR GGTRKYAFEL KEKLLRGTVI DVTDFVNWAS SINDAPVLIS QKLSPiYNLV
 PVKMKNHLK KQNLERAIED YINEFSVRKC HTCQNGGTVI LMDGKCLCAC PFKFEGIACE ISKQKISEGL PALEFPNEK

B4. Purification Strategy/Source

Purified from human serum (see manufacturer website)

Complement Technology

[A126](#)

B5. Stock Concentration/Stock Buffer

1.1 mg/mL | 15 μ M

10 mM sodium phosphate, pH 7.3, 145 mM NaCl

B6. Molecular Weight/Extinction Coefficient

71 kDa

B7. Serial Dilution Preparation

1. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μ L of the 15 μ M C9 solution into tube **1**. Then, transfer 10 μ L of dilution buffer into tubes **2** to **16**.
2. Prepare a 1:1 serial dilution by transferring 10 μ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10 μ L from tube **16** to get an equal volume of 10 μ L for all samples.
3. Mix 4 μ L of labeled CspA with 196 μ L of dilution buffer to obtain 200 μ L of ~32 nM CspA.
4. Add 10 μ L of labeled CspA (~32 nM) 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 Red (NanoTemper Technologies GmbH)

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

D2. MST Software

MO.Affinity Analysis v2.3 (NanoTemper Technologies GmbH)

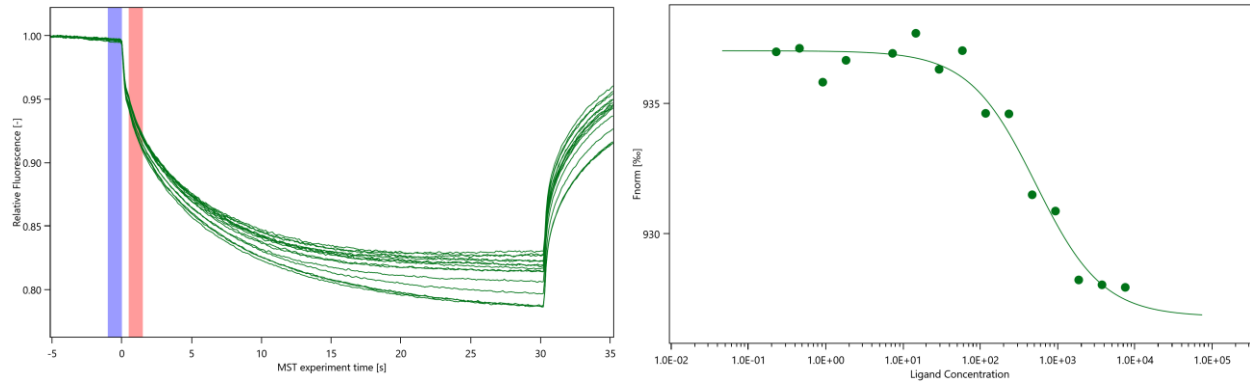
nanotempertech.com/monolith/#monolith-software

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

50 mM Tris-HCl, pH 7.6, 150 mM NaCl, 10 mM MgCl₂, 0.05% TWEEN® 20
 16 nM CspA | 7.5 μM – 229 pM C9 | 25°C | medium MST power | 50% excitation power

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

$K_d = 559$ nM



D5. Reference Results/Supporting Results

N/A

E. Contributors

Teresia Hallström²

² NanoTemper Technologies GmbH, München, Germany | nanotempertech.com