

Monolith Protocol MO-P-064

CCL19 – CCR7 Chemokine Receptor Peptide

Chemokines are small molecular weight proteins which bind to and activate G-protein-coupled receptors (GPCRs) and play a crucial role in cancer cell metastasis. CCR7 is a chemokine receptor which when expressed on the surface of a cell recognizes chemokine CCL19. Interaction of chemokine CCL19 and chemokine receptor CCR7 peptides is characterized as a model system for development of peptide-based PPI inhibitors.

protein – peptide interaction | chemokine | His-tag

A1. Target/Fluorescent Molecule

CCL19

uniprot.org/uniprot/Q99731

A2. Molecule Class/Organism

Chemokine

Homo sapiens (Human)

A3. Sequence/Formula

MALLLALSLL VLWTSAPPTL SGTNDAEDCC LSVTQKPIPG YIVRNFHYLL IKDGCRVPAV VFTTLRGRQL CAPPDQPWVE
RIIQRLQRTS AKMKRRSS

A4. Purification Strategy/Source

Recombinant, N-terminal His-tag, produced in E. coli

Creative Biomart, Shirley, USA

[CCL19-721H](#)

A5. Stock Concentration/Stock Buffer

20 µg lyophilized powder

Phosphate-buffered saline (PBS), pH 7.4, 0.05% TWEEN® 20

A6. Molecular Weight/Extinction Coefficient

10.3 kDa

14,230 M⁻¹cm⁻¹ (ε₂₈₀)

A7. Dilution Buffer

Phosphate-buffered saline (PBS), pH 7.4, 0.05% TWEEN® 20

A8. Labeling Strategy

Monolith His-Tag Labeling Kit RED-tris-NTA (MO-L008, NanoTemper Technologies GmbH)

1* 125 pmol RED-tris-NTA Dye

A9. Labeling Procedure

1. Suspend 20 µg CCL19 in 100 µL of PBS to obtain a 19.5 µM solution.
2. Suspend 125 pmol RED-tris-NTA Dye in 25 µL of dilution buffer to obtain a 5 µM dye solution.
3. Prepare a 100 nM dye solution by mixing 2 µL of dye (5 µM) and 98 µL of dilution buffer.
4. Prepare a 20 nM dye solution by mixing 20 µL of dye (100 nM) and 80 µL of dilution buffer.
5. Prepare a 200 nM CCL19 solution by mixing 1 µL of 19.5 µM CCL19 and 99 µL of dilution buffer.
6. Mix 100 µL of CCL19 (200 nM) with 100 µL of dye (20 nM).
7. Incubate for 30 minutes at room temperature in the dark.

A10. Labeling Efficiency

N/A

B1. Ligand/Non-Fluorescent Binding Partner

CCR7 peptide¹

B2. Molecule Class/Organism

Truncated chemokine receptor peptide

Homo sapiens (Human)

B3. Sequence/Formula

DDYIGDNTTV-NH₂

B4. Purification Strategy/Source

In-house synthesis²

B5. Stock Concentration/Stock Buffer

500 µM

Phosphate-buffered saline (PBS), pH 7.4, 0.05% TWEEN® 20

¹ Exemplified here at the truncated version 10.2.

² See [Fuchs et al., Angew. Chem. Int. Ed. 58 \(21\), 7138–7142 \(2019\)](#) for further information.

B6. Molecular Weight/Extinction Coefficient

1111 Da

1,490 M⁻¹cm⁻¹ (ε₂₈₀)

B7. Serial Dilution Preparation

1. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 µL of the 500 µM CCR7 peptide 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. Add 10 µL of labeled CCL19 (100 nM) to each tube from **16** to **1** and mix by pipetting.
 4. Incubate for 15 minutes at room temperature in the dark before loading capillaries.
-

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)

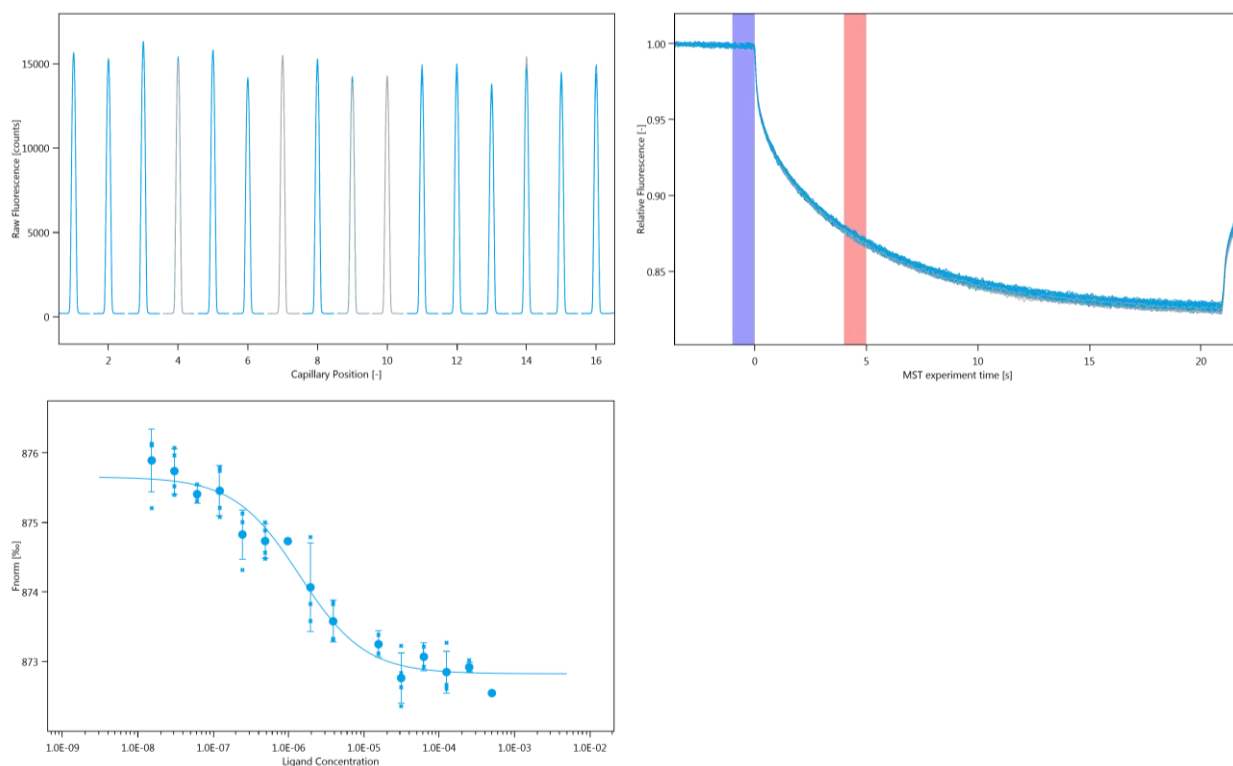
Phosphate-buffered saline (PBS), pH 7.4, 0.05% TWEEN® 20

50 nM CCL19 | 250 µM – 7.63 nM CCR7 | 25°C | medium MST power | 20% excitation power

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

$K_d = 1.4 \mu\text{M}$

Fuchs et al., *Angew. Chem. Int. Ed.* 58 (21), 7138–7142 (2019)³



D5. Reference Results/Supporting Results

N/A

E. Contributors

Cyrril Brunner⁴, Katarzyna Walkiewicz⁵

³ In the publication, final concentrations of 10 nM CCL19 and 5 nM RED-tris-NTA Dye have been used.

⁴ ETH Zurich, Zurich, Switzerland

⁵ NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](https://www.nanotempertech.com)