

Monolith X Protocol MOX-P-112

SecA – SecYEG DIBMA nanodisc

SecY is the main transmembrane subunit of the bacterial Sec pathway and a protein-secreting ATPase complex. Homologs of the SecYEG complex are found in eukaryotes and in archaea. SecA is a cell membrane associated subunit. Within this system the SecA ATPase forms a translocase complex with the SecYEG channel, thereby driving the movement of the protein substrate across the membrane.

protein – protein | membrane protein | nanodisc

A1. Target/Fluorescent Molecule

SecA

[uniprot.org/uniprot/P10408](https://www.uniprot.org/uniprot/P10408)

A2. Molecule Class/Organism

ATPase

E.coli

A3. Sequence/Formula

```
MLIKLLTKVF GSRNDRTLRR MRKVNIINA MEPEMEKLS D EELKGKTAEF RARLEKGEVL ENLIPEAFV VREASKRVFG
MRHFDVQLLG GMVLNERCIA EMRTGEGKTL TATLPAYLNA LTGKGVHVV V NDYLAQRDA ENNRPLFEFL GLTVGINLPG
MPAPAKREAY AADITYGTNN EYGFYLRDN MAFSPEERVQ RKLHYALVDE VDSILIDEAR TPLIISGPAE DSSEMYKRVN
KIIPHLIRQE KEDSETFQGE GHFSVDEKSR QVNLTGERGLV LIEELLVKEG IMDEGESLYS PANIMLMHHV TAALRAHALF
TRDVDYIVKD GEVIVDEHT GRMTQGRWWS DGLHQAVEAK EGVQIQNENQ TLASITFQNY FRLYEKLAGM TGTADTEAFE
FSSIYKLDTV VVPTNRPMIR KDLPLVYMT EAEKIQAIIE DIKERTAKGQ PVLVGTISIE KSELVSNET KAGIKHNVLN
AKFHANAAI VAQAGYPAV TIATNMAGRG TDIVLGGSWQ AEVAALNP AEQIEKIKAD WQVRHDAVLE AGGLHIIGTE
RHESRRIDNQ LRGRSGRQGD AGSSRFYLSM EDALMRIFAS DRVSGMMRKL GMKPGEAIEH PWVTKAIANA QRKVESRNF
IRKQLLEYDD VANDQRAIY SQRNELLDVS DVSETINSIR EDVFKATIDA YIPPPQSLEEM WDIPGLQERL KNDFDLPLI
AEWLDKEPEL HEETLRERIL AQSIQVYQK EEEVGAEMMR HFEGVMLQT LDSLWKEHLA AMDYLRQGIH LRGYAQKDPK
QEYKRESFSM FAAMLESKY EVISTLSKVQ VRMPREEVEL EQRRMEAR LAQMQQLSHQ DDDSA AAAAL AAQTGERKVG
RNDPCPCGSG KKYKQCHGRL Q
```

A4. Purification Strategy/Source

Institute of Biochemistry, Heinrich-Heine-Universität Düsseldorf

A5. Stock Concentration/Stock Buffer

1.6 mg/mL | 16 µM

20 mM Tris/HCl, pH 7.5, 50 mM potassium acetate, 20 % glycerol, 5 mM Magnesium acetate, 1 mM DTT

A6. Molecular Weight/Extinction Coefficient

102 kDa

76,000 M⁻¹cm⁻¹ (ε₂₈₀)

A7. Dilution Buffer

50 mM HEPES, pH 7.4, 150 mM KCl, 5 mM Magnesium chloride, 0.005% Pluronic F-127®

A8. Labeling Strategy

Monolith Protein Labeling Kit RED – NHS 2nd Generation (MO-L011, NanoTemper Technologies GmbH)
1* Dye RED-NHS 2nd Generation (10 µg) | 1* A-Column | 1* B-Column

A9. Labeling Procedure

1. Prepare 40 µL of 16 µM SecA.
2. Use the A-Column to perform a buffer exchange into dilution buffer.
 - a. Invert A-Column to suspend slurry and twist off bottom (twist slightly in both directions).
 - b. Loosen the cap of the column and place it in a 2 mL microcentrifuge collection tube.
 - c. Centrifuge at **1500 × g** for **1 min** to remove excess liquid.
 - d. Add 300 µL of dilution buffer and centrifuge at **1500 × g** for **1 min** (3x).
 - e. Place 40 µL of the 16 µM SecA solution in the center of the resin.
 - f. Place the sample in a **new** microcentrifuge collection tube and centrifuge at **1500 × g** for **2 min**.

The collected flow-through should yield around 40 µL of ~12 µM SecA (~75% recovery).
3. Add 25 µL of DMSO to Dye RED-NHS 2nd Generation (10 µg) to obtain a ~600 µM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
4. Mix 2 µL of the 600 µM dye solution with 38 µL of dilution buffer to obtain 40 µL of a 30 µM dye solution (2.5x protein concentration).
5. Mix SecA and dye in a 1:1 volume ratio (80 µL final volume, 2.5% final DMSO concentration).
6. Incubate for 30 minutes at room temperature in the dark.
7. 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.
8. 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.
9. Add 80 µL of the labeling reaction from step 5 to the center of the column and let sample enter the resin bed completely.
10. Add 520 µL of dilution buffer after the sample has entered and discard the flow through.
11. Place column in a new collection tube, add 480 µL of dilution buffer and collect the eluate.
12. Keep the labeled SecA (~1 µM) on ice in the dark.

A10. Labeling Efficiency

N/A

B1. Ligand/Non-Fluorescent Binding Partner

SecYEG DIBMA nanodisc

SecY

uniprot.org/uniprot/P0AGA2

SecE

uniprot.org/uniprot/P0AG96

SecG

uniprot.org/uniprot/P0AG99

B2. Molecule Class/Organism

Translocon

E.coli

B3. Sequence/Formula

SecY

MAKQPGLDLFDQ SAKGGLGELK RRLLFVIGAL IVFRIGSFIP IPGIDAAVLA KLLEQQRGTI IEMFNMFSGG ALSRASIFAL
GIMPYISASI IIQLLTVVHP TLAEIKKEGE SGRRKISQYT RYGTLLVLAIF QSIGIATGLP NMPGMQGCVI NPGFAFYFTA
VVSLVTGTMF LMWLGQEITE RGIGNGISII IFAGIVAGLP PAIAHTIEQA RQGDHLFLVL LLVAVLVFAV TFFVVFVERG
QRRIVVNYAK RQQGRRVYAA QSTHLPLKVN MAGVIPAIFA SSIILFPATI ASWFGGGTGW NWTTLTISLYL QPGQPLYVLL
YASAIIFFAF FYTALVFNPR ETADNLKKSQ AFVPGIRPGE QTAKYIDKVM TRLTIVGALY ITFIALIPEF MRDAMKVPFY
FGGTSLLIVV VVIMDFMAQV QTLMMSSQYE SALKKANLKG YGR

SecE

MSANTEAQGS GRGLEAMKWV VVALLLVIAI VGNLYLRDIM LPLRALAVVI LIAAAGGVAL LTTKGKATVA FAREARTEVR
KVIWPTREQET LHTTLIVAAV TAVMSLILWG LDGILVRLVS FITGLRF

SecG

MYEALLVVFL IVAIGLVGLI MLQQGKGADM GASFGAGASA TLFSSSGSGN FMTRMTALLA TLFFIISLVL GNINSNKTNK
GSEWENLSAP AKTEQTQPAA PAKPTSDIPN

B4. Purification Strategy/Source

L148C single-cysteine mutant

Institute of Biochemistry, Heinrich-Heine-Universität Düsseldorf

B5. Stock Concentration/Stock Buffer

0.54 mg/mL | 7.5 μ M

50 mM HEPES/KOH, pH 7.4, 150 mM KCl, 5 % Glycerol

B6. Molecular Weight/Extinction Coefficient

72 kDa

72,000 $\text{M}^{-1}\text{cm}^{-1}$ (ϵ_{280})

B7. Serial Dilution Preparation

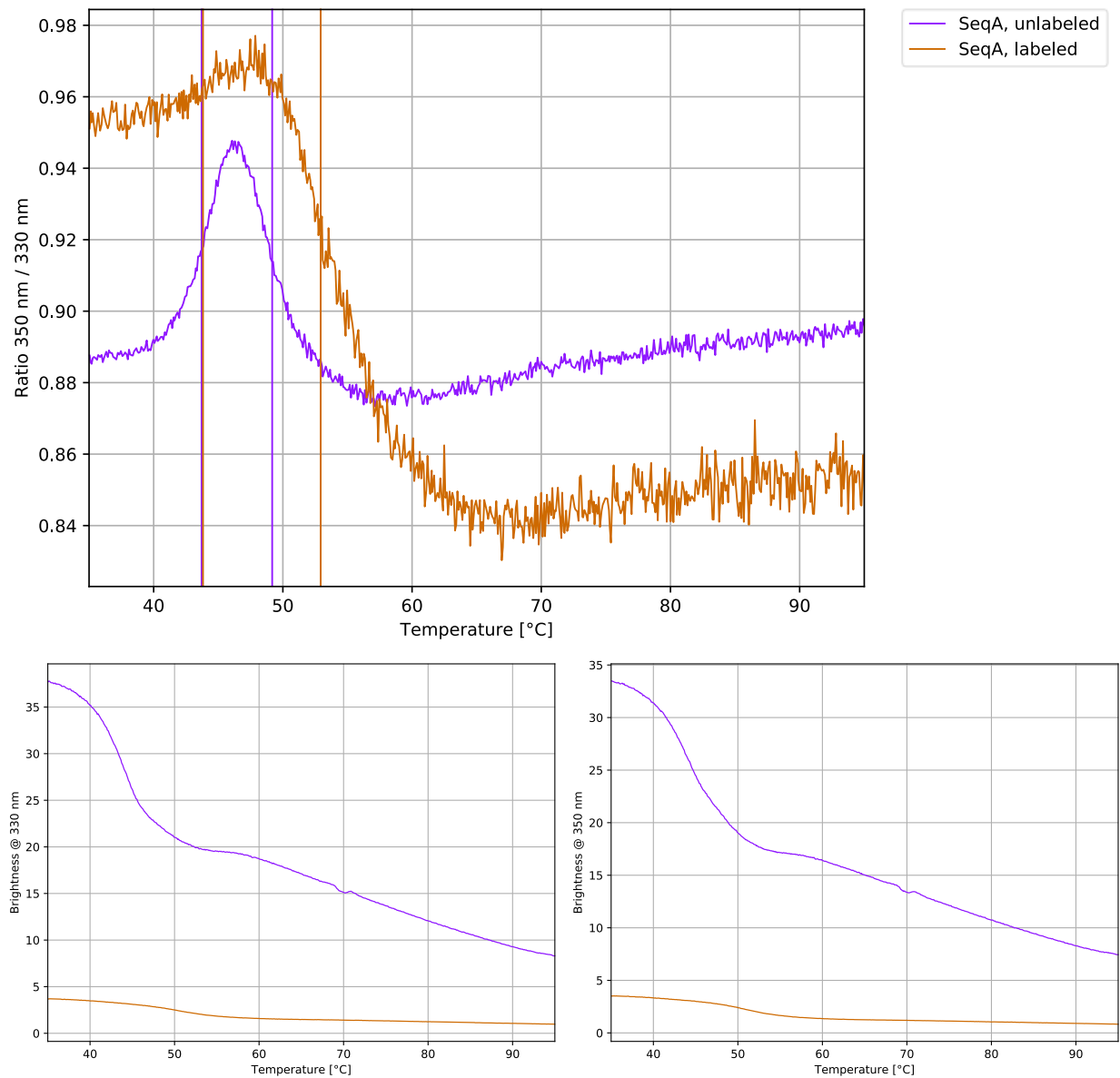
1. Prepare a PCR-rack with 16 PCR tubes. Mix 5.3 μL of 7.5 μM SecYEG DIBMA nanodisc with 14.7 μL of dilution buffer in tube **1**. Then, transfer 10 μL of dilution buffer into tubes **2** to **16**.
2. Prepare a 1:1 serial dilution by transferring 10 μL from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10 μL from tube **16** to get an equal volume of 10 μL for all samples.
3. Mix 8 μL of the labeled SecA ($\sim 1 \mu\text{M}$) with 192 μL of dilution buffer to obtain 200 μL of a $\sim 40 \text{ nM}$ SecA solution.
4. Add 10 μL of labeled SecA ($\sim 40 \text{ nM}$) to each tube from **16** to **1** and mix by pipetting.
5. Incubate for 30 minutes on ice in the dark before loading capillaries.

C. Tycho

Validation of structural integrity of labeled SecA using Tycho NT.6:

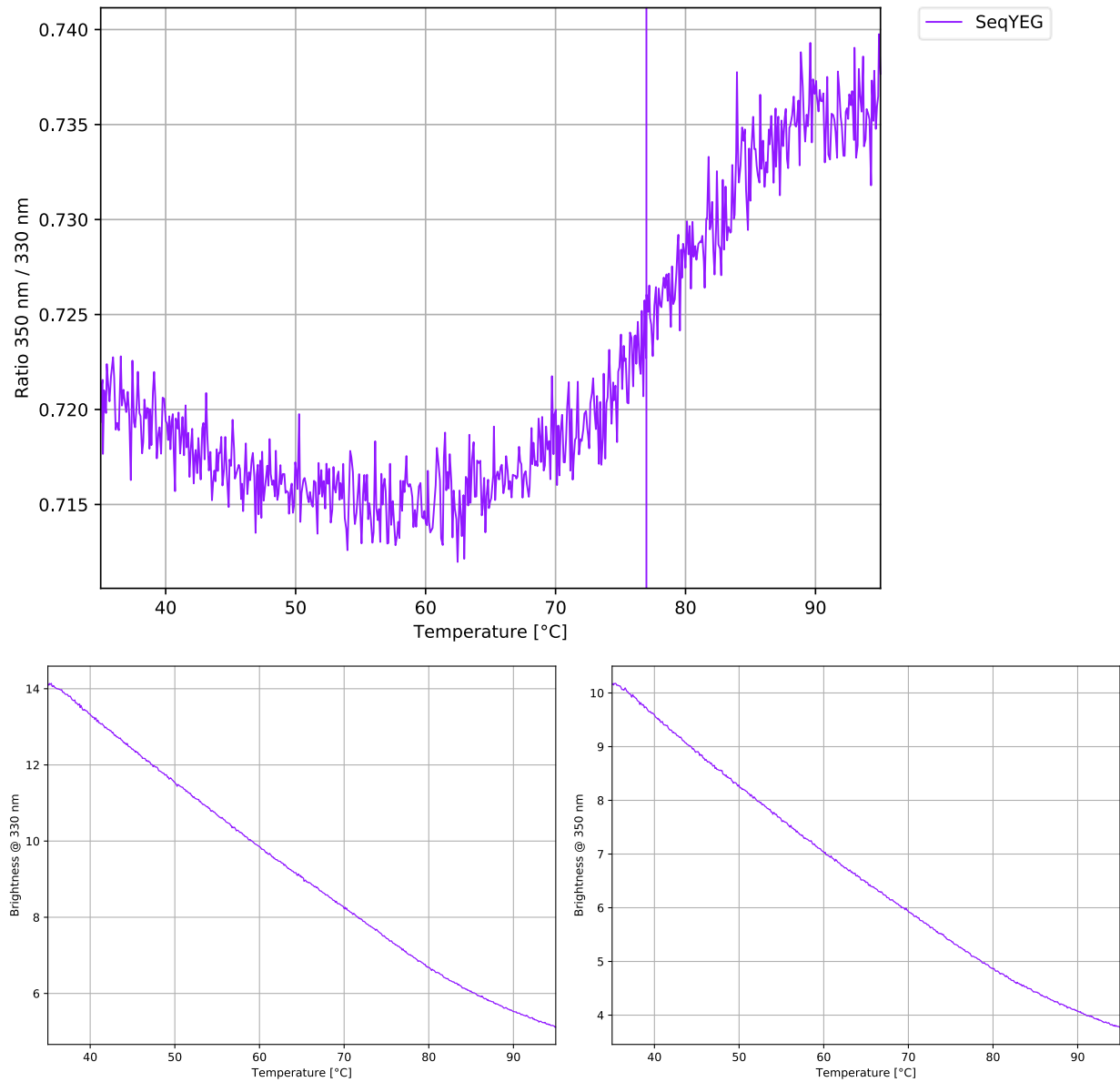
nanotempertech.com/tycho

SecA, unlabeled	0.6 µL of 16 µM SecA + 9.4 µL of dilution buffer	T _{i,1} = 43.7°C T _{i,2} = 49.2°C
SecA, labeled	10 µL of B-Column eluate (~1 µM)	T _{i,1} = 43.8°C T _{i,2} = 52.9°C



Validation of structural integrity of SecYEG using Tycho NT.6:
nonotempertech.com/tycho

SecYEG	0.7 µL of 7.5 µM SecYEG DIBMA nanodisc + 9.3 µL of dilution buffer	T _i = 77.0°C
--------	--	-------------------------



D1. Monolith System/Capillaries

Monolith X (NanoTemper Technologies GmbH)

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

D2. Monolith Software

MO.Control v2.4.2 (NanoTemper Technologies GmbH)

nanotempertech.com/monolith-mo-control-software

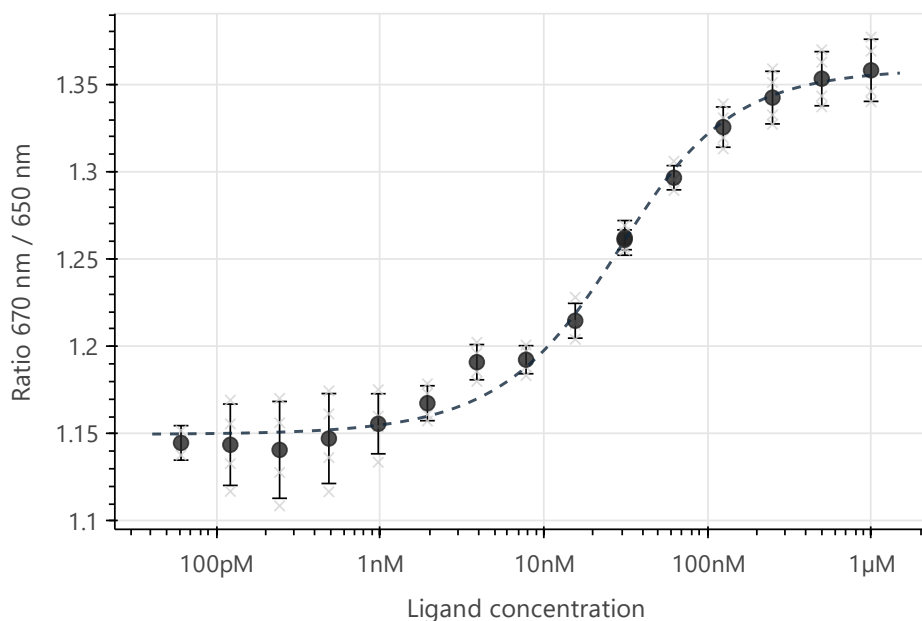
D3. Monolith Experiment (Assay Buffer/Concentrations/Temperature/ Excitation Power)

50 mM HEPES, pH 7.4, 150 mM KCl, 5 mM Magnesium chloride, 0.005% Pluronic F-127®

20 nM SecA | 1 μ M – 30.5 pM SecYEG DIBMA nanodisc | 20°C | 100% excitation power

D4. Monolith Results (Dose Response)

$K_d = 18.3 \pm 2.6$ nM (S/N = 28.5)



D5. Reference Results/Supporting Results

$K_d = 50$ nM MicroScale Thermophoresis (MST)

[unpublished results](#)

$K_d = 4.5$ nM Surface Plasmon Resonance (SPR)

[unpublished results](#)

E. Contributors

Andreas Langer¹, Alexej Kedrov²

¹ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com

² Institute of Biochemistry, Heinrich-Heine-Universität, Düsseldorf, Germany