

Monolith X Protocol MOX-P-113

SecYEG DIBMA nanodisc – SecA

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

SecYEG DIBMA nanodisc

SecY

uniprot.org/uniprot/P0AGA2

SecE

uniprot.org/uniprot/P0AG96

SecG

uniprot.org/uniprot/P0AG99

A2. Molecule Class/Organism

Translocon

E.coli

A3. Sequence/Formula

SecY

MAKQPGDLDFQ SAKGGLGELK RRLLFVIGAL IVFRIGSFIP IPGIDAAVLA KLEEQQRGTI IEMFNMFSGG ALSRASIFAL
GIMPYISASI IIQLLTVVHP TLAEIKKEGE SGRRKISQYT RYGTLVLAIF QSIGIATGLP NMPGMQGCVI NPGFAFYFTA
VVS LVTGTMF LMWLGEQITE RGIGNGISII IFAGIVAGLP PAIAHTIEQA RQGD LHF LVL LLVAVLVFAV TFFVVFVERG
QRRIVVNYAK RQQR RRVYAA QSTHLPLKVN MAGVIPAIFA SSIILFPATI ASWFGGGGTGW NWLTTISLYL QPGQPLYVLL
YASAIFFAF FYTALVFNPR ETADNLKKS G AFVPGIRPGE QTAKYIDKVM TRLT LVGALY ITFIALIPEF MRDAMKVPFY
FGGTSL L I V V VIMDFMAQV QTLMMSSQYE SALKKANLKG YGR

SecE

MSANTEAQGS GRGLEAMKVV VVVALLLVAI VGNYLYRDIM LPLRALAVVI LIAAAGGVAL LTTKGKATVA FAREARTEVR
KVIWPT RQET LHTTLIVAAV TAVMSLILWG LDGILVRLVS FITGLRF

SecG

MYEALLVFL IVAIGLVGLI MLQQGKGADM GASFGAGASA TLFGSSGSGN FMTRMTALLA TLFFIISLVL GNINSNKTNK
GSEWENLSAP AKTEQTQPAA PAKPTSDIPN

A4. Purification Strategy/Source

L148C single-cysteine mutant

Institute of Biochemistry, Heinrich-Heine-Universität Düsseldorf

A5. Stock Concentration/Stock Buffer

0.54 mg/mL | 7.5 μ M
50 mM HEPES/KOH, pH 7.4, 150 mM KCl, 5 % Glycerol

A6. Molecular Weight/Extinction Coefficient

72 kDa
72,000 M⁻¹cm⁻¹ (ϵ_{280})

A7. Dilution Buffer

50 mM HEPES, pH 7.4, 150 mM KCl, 5 mM Magnesium acetate¹

A8. Labeling Strategy

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

A9. Labeling Procedure

1. Prepare 40 μ L of 7.5 μ M SecYEG.
2. Add 22 μ L of DMSO to Dye RED-MALEIMIDE 2nd Generation (10 μ g) to obtain a ~600 μ M solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
3. Mix 2 μ L of the 600 μ M dye solution with 38 μ L of dilution buffer to obtain 40 μ L of a 30 μ M dye solution (4x protein concentration).
4. Mix SecYEG and dye in a 1:1 volume ratio (80 μ L final volume, 2.5% final DMSO concentration).
5. Incubate for 60 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 80 μ L of the labeling reaction from step 4 to the center of the column and let sample enter the resin bed completely.
9. Add 600 μ L of dilution buffer after the sample has entered and discard the flow through.
10. Place column in a new collection tube, add 400 μ L of dilution buffer and collect the eluate.
11. Prepare a 5 mg/mL BSA solution in dilution buffer. Then, add 8 μ L of the 5 mg/mL BSA solution to the 400 μ L labeled SecYEG (~1 μ M) to obtain a final BSA concentration of 0.1 mg/mL.
12. Prepare 10 μ L aliquots and store at -80°C.

A10. Labeling Efficiency

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

nanotempertech.com/dol-calculator

Absorbance A ₂₈₀	0.072	Protein concentration	~0.96 μ M
Absorbance A ₆₅₀	0.094	Degree-of-labeling (DOL)	~0.51

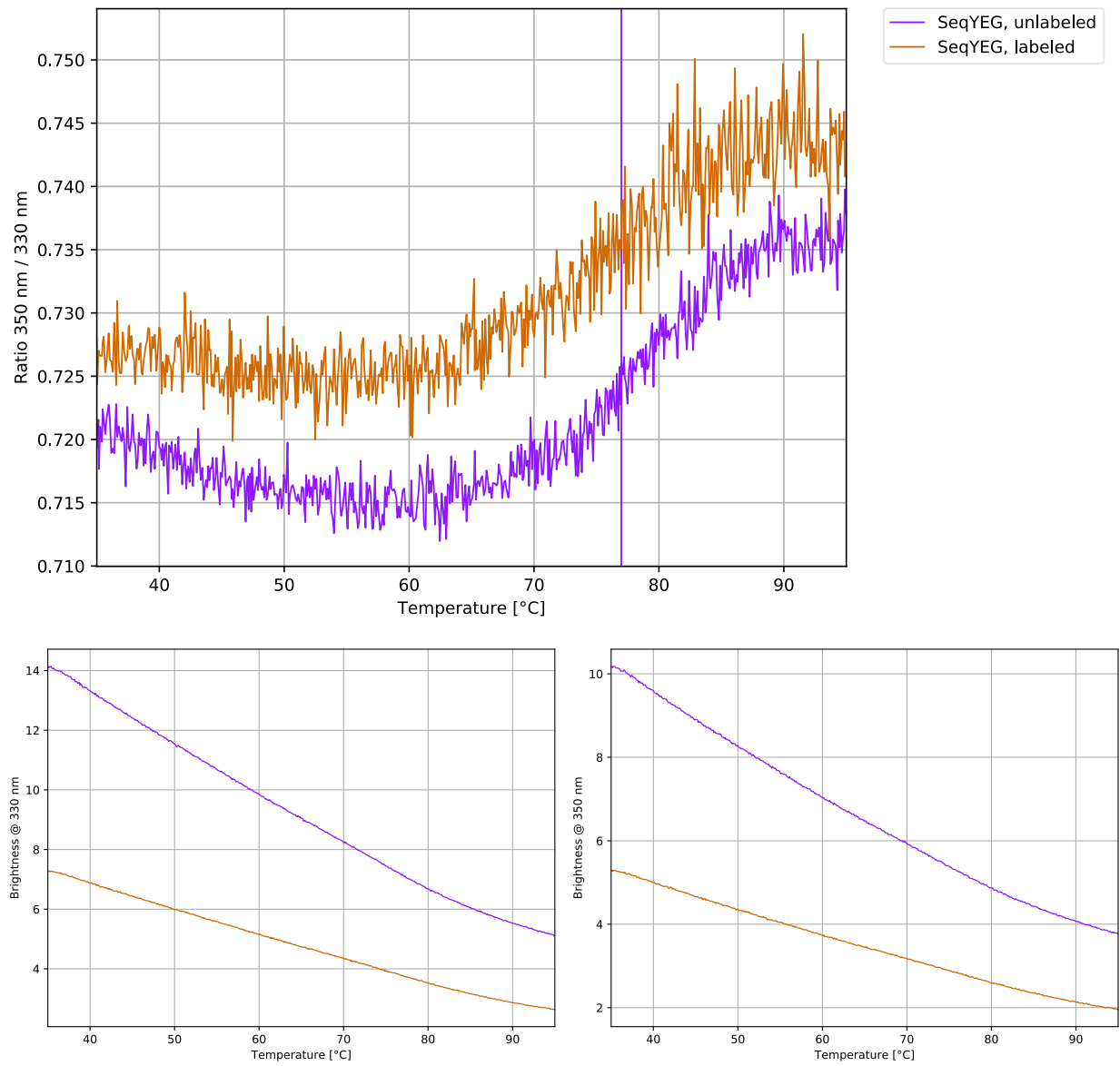
¹ Note that no detergent can be used due to the nanodisc. Instead, BSA is added after the labeling reaction is completed.

C. Tycho

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

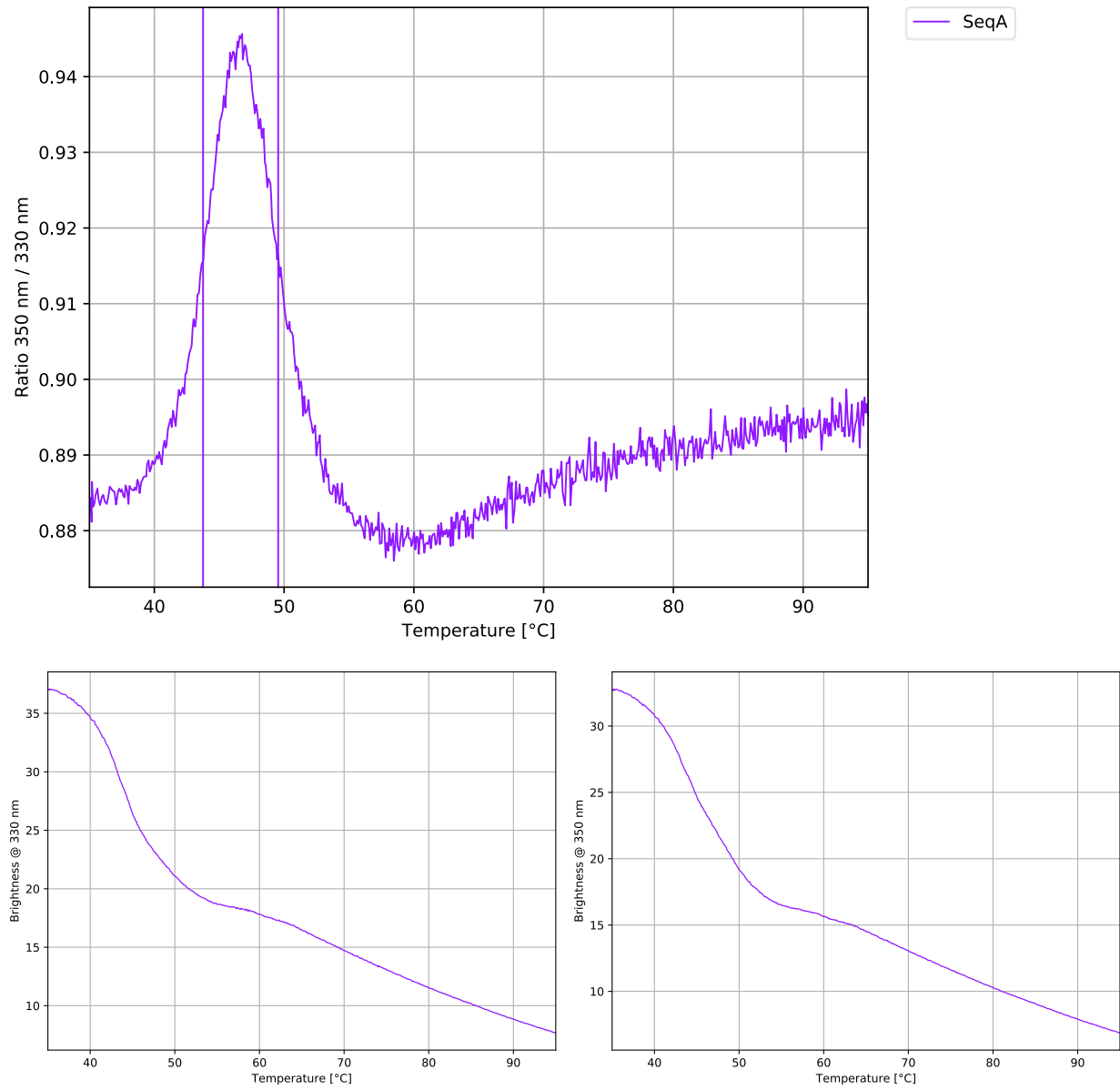
nanotempertech.com/tycho

SecYEG, unlabeled	0.35 μ L of 7.5 μ M SecYEG + 9.7 μ L of dilution buffer	$T_i = 77.0^{\circ}\text{C}$
SecYEG, labeled	5 μ L of B-Column eluate ($\sim 1 \mu\text{M}$) + 5 μ L of dilution buffer	$T_i = 77.0^{\circ}\text{C}$



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

SecA	0.6 µL of 16 µM SecA + 9.4 µL of dilution buffer	$T_{i,1} = 43.8^{\circ}\text{C}$ $T_{i,2} = 49.5^{\circ}\text{C}$
------	--	---



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

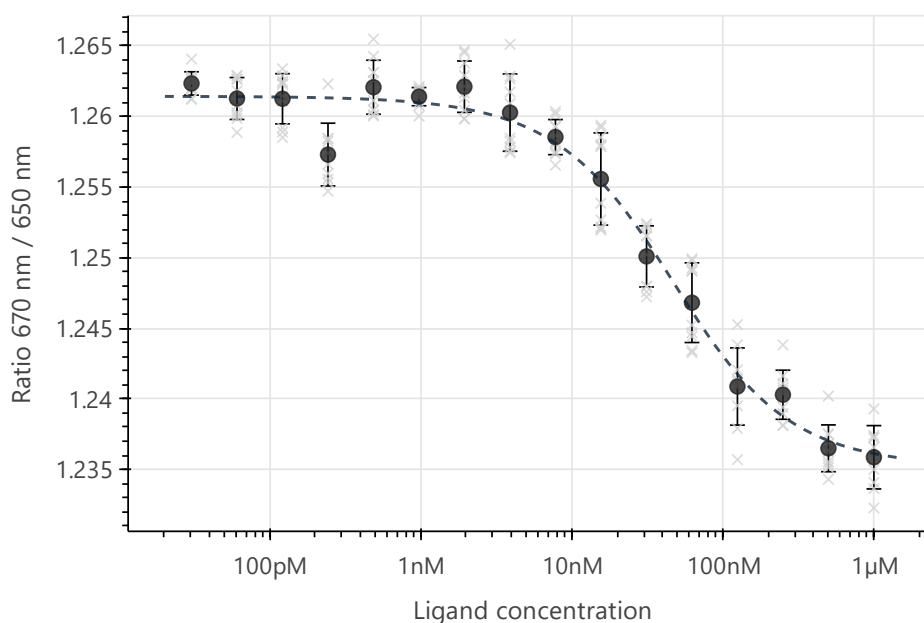
D3. Monolith Experiment (Assay Buffer/Concentrations/Temperature/ Excitation Power)

50 mM HEPES, pH 7.4, 150 mM KCl, 5 mM Magnesium acetate, 0.05 mg/mL BSA

20 nM SecYEG | 1 μ M – 30.6 pM SecA | 20°C | 100% excitation power

D4. Monolith Results (Dose Response)

$K_d = 36.8 \pm 7.9$ nM (S/N = 19.2)



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 Langer², Alexej Kedrov³

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

³ Institute of Biochemistry, Heinrich-Heine-Universität, Düsseldorf, Germany