

## Monolith Protocol MO-P-044

# Sc-tRNA – Elongator Subcomplex yElp456

tRNA is a class of RNA molecules that plays a role during the translation process – the synthesis of functional proteins. Recently, modifications of RNA molecules were shown to be linked with their biological activity. The elongator complex Elp is formed by two subcomplexes that consist of structurally arranged Elp1, Elp2, Elp3 (Elp123 subunit) and Elp4, ELp5, Elp6 (Elp456 subunit). The Elp456 subunit forms a heterohexameric ring-like structure and is involved in specific interactions with tRNA molecules.

protein – RNA interaction | elongator protein | tRNA binding | tRNA modifications

## A1. Target/Fluorescent Molecule

Sc-tRNA<sub>Ala</sub>

## A2. Molecule Class/Organism

RNA

*Saccharomyces cerevisiae*

## A3. Sequence/Formula

5' GGG CAC AUG GCG CAG UUG GUA GCG CGC UUC CCU UGC AAG GAA GAG GUC AUC GGU UCG AUU CCG  
GUU GCG UCC A 3'

## A4. Purification Strategy/Source

In vitro transcription of tRNA with T7 polymerase:

1. The template for transcription is prepared as double-stranded DNA in a standard PCR approach.<sup>1</sup>
2. Large scale purification of the template is achieved from 2L 2xYT-carbenicilin culture and extraction is performed with QIAfilter® plasmid MEGA or GIGA protocols (Qiagen™) – this process typically yields 3 – 6 mg of DNA.
3. The DNA is dissolved in water and linearized with BbsI or BsaI, using an enzyme concentration of 50 U/ml and overnight incubation at 37°C (BbsI) or 50°C (BsaI).
4. The large scale 10 – 40 mL transcription reaction contains 4 mM rNTPs, 1.3 mM CTP-Cy5, 70 µg/ml linearized DNA template, 1200 U/ml T7 RNA polymerase, 40 mM Tris-HCl (pH 8.1 @37°C), 1 mM spermidine, 5 mM DTT, 0.1% Triton-X, 1 U/ml inorganic pyrophosphatase and 40 mM MgCl<sub>2</sub>. The transcription reaction is incubated for 2 – 4 h at 37°C and stopped by addition of 50 mM EDTA. The transcripts could be stored at –20°C until RNA purification.

<sup>1</sup> Lukavsky and Puglisi, RNA (2004), 10:889–893

5. tRNAs are purified using an ÄKTA-FPLC system, equipped with a 50 mL superloop and three connected in tandem 5 mL HiTrap DEAE columns, using Buffer A (50 mM sodium phosphate, 150 mM sodium chloride and 0.2 mM EDTA, pH 6.5 at room temperature) and Buffer B (50 mM sodium phosphate, 2 M sodium chloride and 0.2 mM EDTA, pH 6.5 at room temperature). Columns are equilibrated with 3 CV of Buffer A. The stopped transcription reaction is loaded into the 50 mL superloop and weak anion-exchange chromatography is performed using the following gradient and collecting 10 mL fractions in sterile 15 mL tubes.
  - a. Gradient step 1: 0 – 70 mL, 0% B at 1 mL/min – loading sample on DAEA columns
  - b. Gradient step 2: 70 – 100 mL, 0% – 10% buffer B at 2 mL/min – wash free rNTPs
  - c. Gradient step 3: 100 – 380 mL, 10% – 30% buffer B at 2 mL/min – separate short abortive transcripts, RNA product, and plasmid DNA
  - d. Gradient step 4: 380 – 410 mL, 30% – 100% buffer B at 4 mL/min – columns wash
  - e. Gradient step 5: 410 – 455 mL, 100% buffer B at 4 mL/min – columns wash
  - f. Gradient step 6: 455 – 485 mL, 100% – 0% buffer B at 4 mL/min – columns re-equilibration
6. 5 µL of each fraction is analyzed by denaturing PAGE (8% acrylamide, 8 M urea)
7. Optional: In case RNA profiles display more than one peak along gradient suggesting alternative, conformations or multimerization of the RNA, each fraction should be analyzed by HPLC-SEC using e.g. a SEC-125 column and Buffer A.
8. Structurally homogeneous RNA is pooled, concentrated and equilibrated into appropriate buffers using 15 mL Centriprep centrifugal devices.

#### A5. Stock Concentration/Stock Buffer

280 µg/ml | 12 µM

20 mM HEPES pH7.5, 50 mM NaCl, 50 mM KCl, 1 mM MgCl<sub>2</sub>, 2 mM DTT

#### A6. Molecular Weight/Extinction Coefficient

24.9 kDa

940,500 M<sup>-1</sup>cm<sup>-1</sup> (ε<sub>260</sub>)

#### A7. Dilution Buffer

20 mM HEPES, pH 7.5, 50 mM NaCl, 0.05% TWEEN® 20, 2 mM DTT

#### A8. Labeling Strategy

Labeled during in-vitro transcription by internal covalent attachment of Cy5

#### A9. Labeling Procedure

Labeling protocol available from New England Biolabs® for RNA synthesis with modified nucleotides (E2050)™. The molar ratio of modified NTP to standard NTP was 1:3.

#### A10. Labeling Efficiency

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

[nanotempertech.com/dol-calculator](https://nanotempertech.com/dol-calculator)

Absorbance A <sub>260</sub>	17.424	Protein concentration	18.5 µM
Absorbance A <sub>647</sub>	0.134	Degree-of-labeling (DOL)	0.0289

## B1. Ligand/Non-Fluorescent Binding Partner

yElp456 wt

## B2. Molecule Class/Organism

tRNA binding protein  
*Saccharomyces cerevisiae*

## B3. Sequence/Formula

Elp4

MSFRKRGEIL NDRGSGLRGP LLRGPPRTSS TPLRTGNRRA PGNVPLSDTT ARLKKLNIAD ESKTKMGLDS SHVGVRPSA  
TSQPTTSTGS ADLDSILGHM GLPLGNSVLV EEQSTTEFHS ILGKLFAAQG IVHNRISDSS ADKTRNGDTH VIVLSLNQMF  
AKELPGIYKG SRKQMKKNLI SEEEKSVTVQ NLNETQRSTP SRYKDLKIAW KYKLADEKRL GSPDRDDIQQ NSEYKDYNHQ  
FDITTRLMPA PIASELTFIA PTQPVSTILS QIEQTIKRNQ KKLIRIVIPS LLHPAMYPPK MFESSEIIGL MHGVRSLVKK  
YYERVVLFAS ISIDIITPPL LVLLRNMFDV VINLEPFNQE MTEFLERVYK SQPGKIQHGL VHILKLPVFT DRGEMRVLKS  
EWAFKNGRKK FEIEQWIGIPV DDAEGSAASE QSHSHSHSDE ISHNIPAKKT KISLDY

Elp5

MASSSHNPVI LLKRILSLTE SSPFILCLDS IAQTSYKLIQ EFVHQSKSKG NEYPIVYISF ETVNKPSYCT QFIDATQMDF  
VHLVKQIISY LPAATATQAK KHMVIIDSLN YISTEYITRF LSEIASPHCT MVATYHKDIK DENRTVIPDW NNNYPDKLTL  
LQFMATTIVD IDVVLGTGLD TEEVSELLNE FRIPRGLNND IFQLRLVNKR KSGRSLEYDF IVNSNTHEYE LLSTTKQEEE  
SSSNGLETPE MLQGLTTFNL GTSNKQKLAK DQVALPFLEA QSFQGGGAIW YEYKDDDDYD EEDPYEDPF

Elp6

MGSVQRQDLV LFSDQSVLPA HFFQDSNSHN LFFITHQSCT QPLWMINALV ETHVLGSPSS LNESSSSMLP SSTRSHAVLA  
SFIHEQNYFT NSLNKLKIPS NNYNVLDFLS DFIVNNIHNK PRDKILSDVL AKFSAAIQNN PTDTIVIIIEQ PELLLSLVSG  
LTCSELNNKF ITPLLRQCKV LIIVSNSDIF NIDEYDASVH SSNLQNFYKS SFIKSMINLN LNPLKTGFAK DVTGSLHVCR  
GGAPIATSNT SLHVVENEYL YLNEKESTKL FYR

## B4. Purification Strategy/Source

The purification strategy included trial-and-error process of expression and purification of different combinations of full-length Elp4, Elp5, Elp6 (including single 6xHis-tagged protein), followed by mass spectrometry analysis to define stable, interacting and well-expressing hexameric Elp456 protein construct.

The final construct, pBS3576 encodes 6xHis-tag-Elp4<sub>1-273</sub>, Elp5<sub>1-270</sub>, Elp6<sub>66-426</sub> in tandem for co-expression using pET24d derived vector. Constructs were expressed in *E. coli* (BL21 Gold (DE3) incubated at 18°C overnight for 12 – 15 h). Bacterial cells were lysed using a French press in 50 mM HEPES (pH 7.5), 300 mM NaCl, 10 mM imidazole, 1 mM DTT, 5% (v/v) glycerol and protease inhibitors. The soluble fraction was cleared by centrifugation (20,000× g for 45 min at 4°C). 6-His-tagged complexes were purified using Ni-NTA affinity chromatography followed by gel filtration with a Superdex 200 column (GE Healthcare™) in 20 mM Hepes-NaOH (pH 7.5), 150 mM NaCl, 2 mM DTT.

## B5. Stock Concentration/Stock Buffer

9 mg/mL | 43.7 μM  
20 mM HEPES pH 7.5, 50 mM NaCl, 2 mM DTT

## B6. Molecular Weight/Extinction Coefficient

206 kDa (theoretical: 136.2 kDa)

## B7. Serial Dilution Preparation

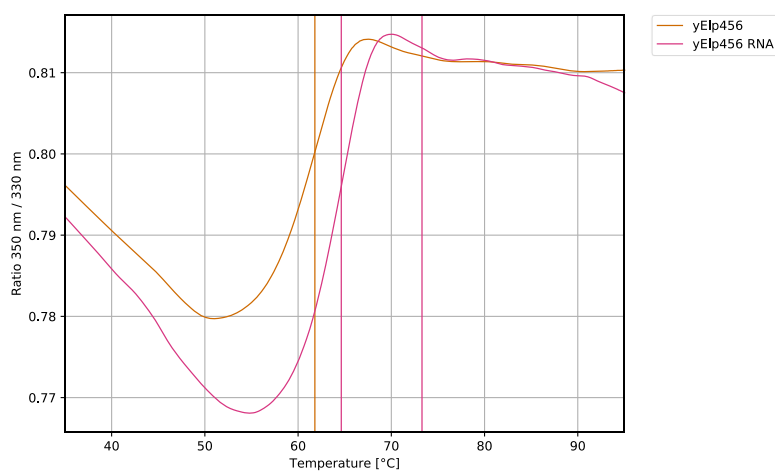
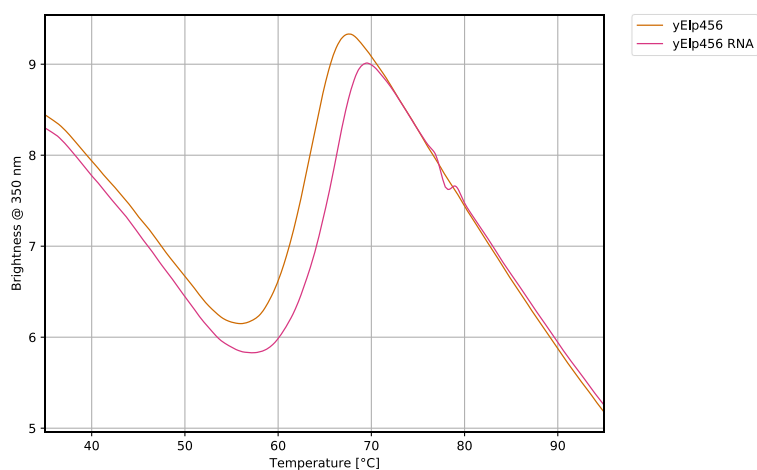
1. Add 193  $\mu\text{L}$  of dilution buffer to 7  $\mu\text{L}$  of labeled Sc-tRNA<sub>Ala</sub> to obtain 200  $\mu\text{L}$  of a 420 nM solution.
2. Add 16  $\mu\text{L}$  of dilution buffer to 14  $\mu\text{L}$  of 43.7  $\mu\text{M}$  yElp456 to obtain 30  $\mu\text{L}$  of a 20  $\mu\text{M}$  solution.
3. Prepare a PCR-rack with 16 PCR tubes. Transfer 20  $\mu\text{L}$  of the 20  $\mu\text{M}$  yElp456 solution into tube **1**. Then, transfer 10  $\mu\text{L}$  of dilution buffer into tubes **2** to **16**.
4. Prepare a 1:1 serial dilution by transferring 10  $\mu\text{L}$  from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10  $\mu\text{L}$  from tube **16** to get an equal volume of 10  $\mu\text{L}$  for all samples.
5. Add 10  $\mu\text{L}$  of 420 nM labeled Sc-tRNA<sub>Ala</sub> to each tube from **16** to **1** and mix by pipetting.
6. Incubate for 20 minutes at room temperature in the dark before loading capillaries.

## C. Applied Quality Checks

Validation of structural integrity and functionality of yElp456 using Tycho NT.6:

[nanotempertech.com/tycho](https://nanotempertech.com/tycho)

yElp456	yElp456	$T_i = 61.8^\circ\text{C}$
yElp456 RNA	yElp456 + Sc-tRNA <sub>Ala</sub>	$T_i = 64.6^\circ\text{C}$



## 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.AffinityAnalysis v2.3 (NanoTemper Technologies GmbH)

[nanotempertech.com/monolith/#monolith-software](https://nanotempertech.com/monolith/#monolith-software)

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

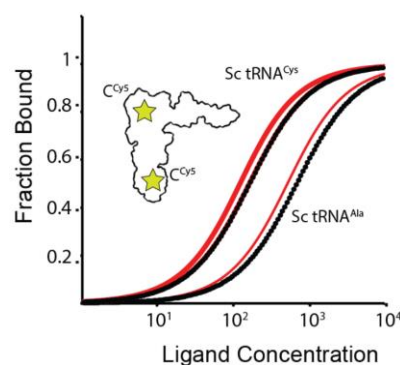
20 mM HEPES pH 7.5, 50 mM NaCl, 0.05% TWEEN® 20, 2 mM DTT

210 nM Sc-tRNA<sub>Ala</sub> | 10 µM – 31 nM yElp456 | 20°C | low MST power | 20% excitation power

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

$K_d = 605.7 \pm 17$  nM

[Kojic et al., Nature Communications 9:3195 \(2018\)](#)



## D5. Reference Results/Supporting Results

N/A

## E. Contributors

Monika Gaik<sup>2</sup>, Sebastian Glatt<sup>2</sup>, Ting-Yu Lin<sup>2</sup>, Jakub Nowak<sup>3</sup>

<sup>2</sup> Malopolska Centre of Biotechnology, Jagiellonian University, Krakow, Poland

<sup>3</sup> NanoTemper Technologies GmbH, München, Germany | [nanotempertech.com](https://nanotempertech.com)