

Monolith Protocol MO-P-027

30S Ribosomal Subunit – YjeQ

The ribosome is responsible for translating the genetic code into functional proteins and is often a target for antimicrobial research. Specifically, the 30S ribosomal subunit decodes mRNA to ensure high fidelity during protein synthesis. Although much is known about protein translation, the details surrounding ribosome biogenesis remain largely elusive. Ribosome biogenesis is a complex process that requires the assembly of multiple proteins and rRNA molecules to form a functional 70S ribosome in bacteria. This process is aided by multiple assembly factors that transiently interact with assembling ribosome particles to aid in their maturation. A central protein involved in this process is the GTPase YjeQ, which has been associated with aiding in the later stages of 30S maturation.

protein – protein interaction | ribosome assembly | GTPase | biogenesis

A1. Target/Fluorescent Molecule

Prokaryotic small ribosomal subunit (30S ribosomal subunit)

A2. Molecule Class/Organism

Ribonucleoprotein
Escherichia coli

A3. Sequence/Formula

N/A¹

A4. Purification Strategy/Source

Isolated from wild type *E. coli* (BW25113)
Purified using ultracentrifugation over sucrose cushions and gradients²

A5. Stock Concentration/Stock Buffer

4.25 mg/mL | 5 μ M
10 mM Tris-HCl, pH 7.5, 1.1 mM magnesium acetate, 60 mM NH₄Cl, 0.5 mM EDTA, 2 mM 2-mercaptoethanol

A6. Molecular Weight/Extinction Coefficient

850 kDa
13,700,000 M⁻¹cm⁻¹ (ϵ_{260})

¹ The 30S ribosomal subunit is a complex of the 16S ribosomal RNA (rRNA) and 19 proteins.

² See [Leong et al., RNA, 19, 789-802 \(2013\)](#) for further details.

A7. Dilution Buffer

20 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM MgCl₂, 1 mM DTT, 0.05% TWEEN® 20, 0.4 mg/mL BSA, 1 mM GMP-PNP³

A8. Labeling Strategy

Monolith Protein Labeling Kit RED – Maleimide (MO-L004, NanoTemper Technologies GmbH)

1x Labeling Buffer NHS | 1x A-Column | 1x 10 µg NT-647-MALEIMIDE dye | 1x B-Column

A9. Labeling Procedure

1. Prepare 100 µL of a 5 µM 30S ribosomal subunit solution.
2. Use Spin Column A to perform a buffer exchange into Labeling Buffer NHS **supplemented** with 10 mM MgCl₂.
 - a. Invert Spin Column A to suspend slurry and twist off bottom (twist slightly in both directions).
 - b. Loosen the cap of the column and place it in a 1.5 mL microcentrifuge collection tube.
 - c. Centrifuge at **1500 × g** for **1 min** to remove excess liquid.
 - d. Add 300 µL of Labeling Buffer NHS with MgCl₂ and centrifuge at **1500 × g** for **1 min** (3x).
 - e. Place 100 µL of the 5 µM 30S ribosomal subunit 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 100 µL of ~4 µM 30S ribosomal subunit (~80 % recovery).
3. Add 30 µL of DMSO to 10 µg NT-647-MALEIMIDE dye to obtain a ~470 µM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
4. Mix 3.2 µL of the 470 µM dye solution with 96.8 µL of Labeling Buffer NHS with MgCl₂ to obtain 100 µL of a 15 µM dye solution (~4x protein concentration).
5. Mix 30S ribosomal subunit and dye in a 1:1 volume ratio (200 µL final volume).
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 200 µL of the labeling reaction from step 5 to the center of the column and let sample enter the bed completely.
10. Add 400 µL of dilution buffer after the sample has entered and discard the flow through.
11. Place column in a new collection tube, add 500 µL of dilution buffer and collect the eluate.
12. Keep labeled 30S ribosomal subunit (~0.8 µM) on ice in the dark.

A10. Labeling Efficiency

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

nanotempertech.com/dol-calculator

Absorbance A ₂₆₀	7.0	Protein concentration	0.51 µM
Absorbance A ₆₅₀	0.06	Degree-of-labeling (DOL)	0.47

³ GMP-PNP was substituted with GDP for the negative control.

B1. Ligand/Non-Fluorescent Binding Partner

YjeQ (RsgA)

uniprot.org/uniprot/P39286

B2. Molecule Class/Organism

GTPase

Escherichia coli

B3. Sequence/Formula

```
MSKNKLSKGQ QRRVNANHQR RLKTSKEKPD YDDNLFGEPE EGVISRFGM HADVESADGD VHRCNIRRTI RSLVTGDRVV
WRPGKPAEEG VNVKGIVEAV HERTSVLTRP DFYDGVKPIA ANIDQIVIVS AILPELSLNI IDRYLVACET LQIEPIIVLN
KIDLLDDEGM AFVNEQMDIY RNIGYRVLMV SSHTQDGLKP LEEALTGRIS IFAGQSGVGK SLLNALLGL QKEILTNDIS
DNSGLGQHTT TAARLYHFPH GGDVIDSPGV REFGLWHLEP EQITQGFVEF HDYLGGLCKYR DCKHDTDPGC AIREAVEEGK
IAETRFENYH RILEMAQVK TRKNFSDTDD
```

B4. Purification Strategy/Source

YjeQ was overexpressed in BL21-AI competent cells using the pDEST17-yjeQ plasmid with an N-terminal His₆-tag cleavable by tobacco etch virus (TEV) protease⁴. One liter of LB medium was inoculated with 10 mL of saturated overnight culture and cells were grown to OD₆₀₀ = 0.6 by incubation at 37 °C and shaking at 225 rpm in an Excella E24 incubator (New Brunswick). Expression was induced with 0.2% L-arabinose. Cells were then induced for 3 hours at 37 °C and harvested by centrifugation at 3700 × g for 15 min. Cell pellets were washed with 1X PBS buffer (137 mM NaCl, 2.7 mM KCl, 8.1 mM Na₂HPO₄ at pH 7.4) and resuspended in 20 mL of lysis buffer (50 mM Tris-HCl at pH 8.0, 10% [w/v] sucrose, 100 mM NaCl) containing a protease inhibitor cocktail (cCompleteTM Mini Protease Inhibitor Cocktail Tablets, Roche).

The cell suspension was passed through a French pressure cell at 1400 kg/cm² three consecutive times to lyse the cells. The lysate was spun at 39,200 × g for 45 min to clear cell debris and the supernatant was collected. All the following steps were performed at 4 °C. NaCl was added to the supernatant of the YjeQ overexpression cells to a concentration of 0.5 M. Clarified cell lysates of YjeQ were passed through a 0.45-µm syringe filter (Millipore) and loaded to a HiTrap Metal Chelatin2 Column (GE Healthcare Life Sciences) previously equilibrated with 50 mM Tris-HCl at pH 8.0, 0.5 M NaCl and 5% [v/v] glycerol. Nonspecifically bound proteins were washed with incremental step-wise increases in the concentration of imidazole from 45 mM to 90 mM. YjeQ was eluted with 240 mM imidazole. Purity of the fractions was monitored by SDS-PAGE and fractions containing pure (>95%) YjeQ protein were collected and pooled together.

The N-terminal His₆-tag was removed by digestion with TEV protease at a ratio of 10:1 (YjeQ:TEV) during overnight dialysis against buffer containing 50 mM Tris-HCl at pH 8.0, 60 mM imidazole and 0.2 M NaCl. Any overnight precipitate was removed by spinning at 12,000 × g for 10 min and the supernatant was collected and loaded onto a HiTrap Metal Chelating Column previously equilibrated with 50 mM Tris-HCl at pH 8.0, 0.2 M NaCl and 60 mM imidazole. Fractions were collected, and their purity evaluated by SDS-PAGE and Coomassie Brilliant Blue staining. Fractions containing pure untagged YjeQ were pooled and dialyzed against 50 mM Tris-HCl at pH 8.0, 5% [v/v] glycerol overnight. A 10 kDa-cutoff filter (Amicon) was used to concentrate the protein and the purified YjeQ was frozen in liquid nitrogen and stored at -80 °C.

⁴ See [Jomaa et al., RNA, 17, 2026-2038 \(2011\)](#) for further details.

B5. Stock Concentration/Stock Buffer

2.0 mg/mL | 50 μ M
 50 mM Tris-HCl, pH 8.0, 5% glycerol

B6. Molecular Weight/Extinction Coefficient

39.1 kDa
 24,410 $M^{-1}cm^{-1}$ (ϵ_{280})

B7. Serial Dilution Preparation

1. Centrifuge YjeQ and 30S ribosomal subunits for $14\,000 \times g$ for 10 min and carefully remove supernatant.
 2. Add 2 μ l of unlabeled YjeQ to 48 μ l dilution buffer to obtain 50 μ l of a 2 μ M solution.
 3. Prepare a PCR-rack with 16 PCR tubes. Transfer 20 μ l of the 2 μ M YjeQ solution into tube **1**. Then, transfer 10 μ l of dilution buffer into tubes **2** to **16**.
 4. 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.
 5. Mix 20 μ l of labeled 30S ribosomal subunit with 180 μ l of dilution buffer to obtain 200 μ l of ~80 nM 30S.
 6. Add 10 μ l of 80 nM labeled 30S ribosomal subunit to each tube from **16** to **1** and mix by pipetting.
 7. Incubate for 10 minutes at room temperature in the dark before loading capillaries.
-

D1. MST System/Capillaries

Monolith NT.115 Red (NanoTemper Technologies GmbH)
 Hydrophobic Capillaries Monolith NT.115 (NanoTemper Technologies GmbH)

D2. MST Software

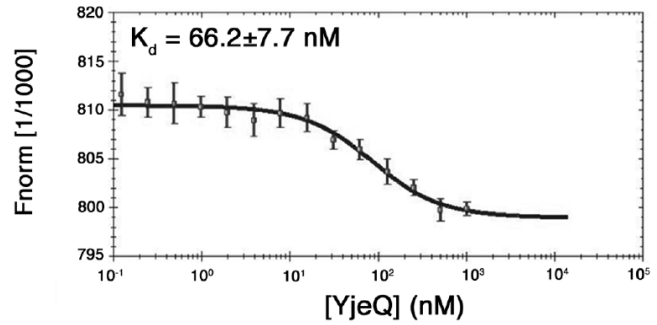
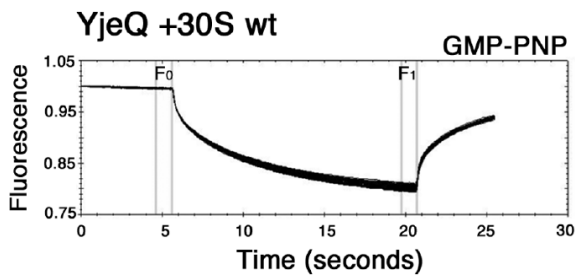
MO.Control v1.6 (NanoTemper Technologies GmbH)
nanotempertech.com/monolith-mo-control-software

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

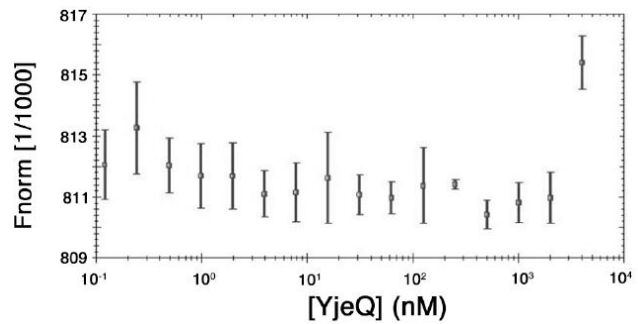
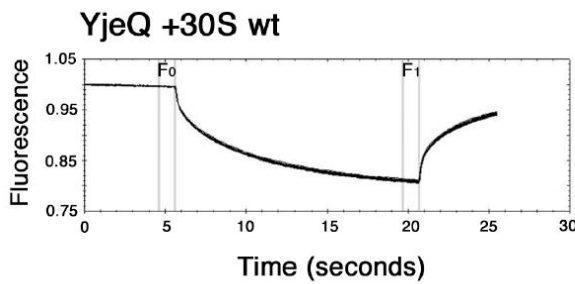
20 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM $MgCl_2$, 1 mM DTT, 0.05% TWEEN® 20, 0.4 mg/mL BSA, 1 mM GMP-PNP
 40 nM 30S ribosomal subunit | 0.03 nM – 1 μ M YjeQ | 25°C | high MST power | 70% excitation power

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

$K_d = 66.2 \pm 7.7$ nM (with GMP-PNP) [Thurlow et al. \(2016\), Nucleic Acids Research, 44\(20\), 9918-9932](#)



$K_d = \text{N/A}$ (with GDP, negative control)



D5. Reference Results/Supporting Results

Interaction of YjeQ with the 30S subunit

[Daigle et al., Journal of Bacteriology, 186, 1381-1387 \(2004\)](#) | [Jegonathan et al., RNA, 21, 1203-1216 \(2015\)](#) | [Razi et al., PNAS, 114, 3396-3403 \(2017\)](#)

E. Contributors

Brett Thurlow⁵

⁵ NanoTemper Technologies, Inc., South San Francisco, United States | nanotempertech.com