



# **Construction of an IGR IRES-mediated reporter gene expression** system based on RNA polymerase I for large scale screening of high-nucleic acid in Saccharomyces cerevisiae Duwen Zeng<sup>1</sup>, Chenxi Qiu<sup>2</sup>, Lili Xu<sup>1\*</sup>, Xiaoming Bao<sup>1,2,\*</sup>

1 State Key Laboratory of Biobased Material and Green Papermaking, Qilu University of Technology, Shandong Academy of Sciences, Jinan, 250353, China.

2 State Key Laboratory of Microbial Technology, Institute of Microbial Technology, Shandong University, Qingdao, 266237, China.

\* Corresponding author: Xiaoming Bao. E-mail address: baoxm@qlu.edu.cn

Introduction

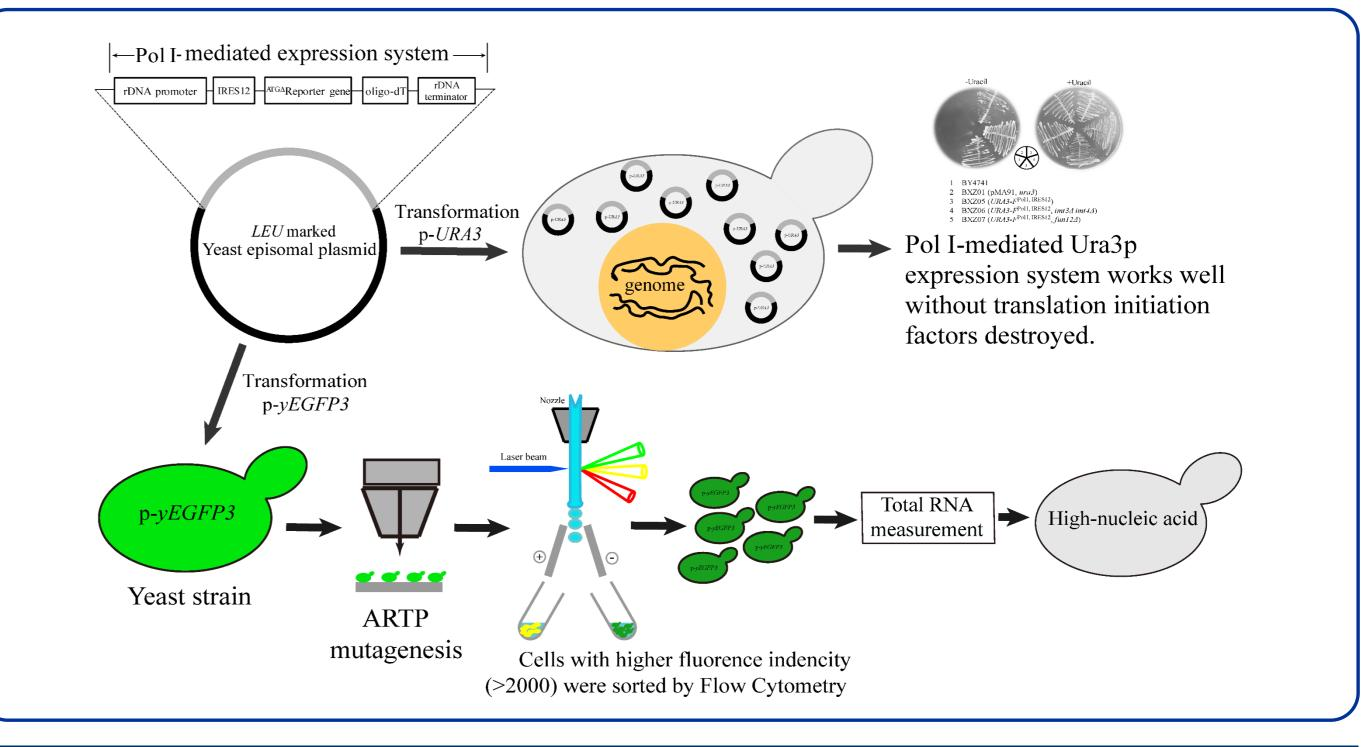
• Derivatives from Ribonucleic acid (RNA) are widely used in food and medicine homologous. rRNA, which is the most abundant RNAs, was synthesized by the tight regulation of RNA polymerase I (RNA Pol I) and other factors.

• Saccharomyces cerevisiae is recognized as the ideal source of RNA. At present, RNA is mainly extracted from discarded beer yeast. Affected by the raw material supply, RNA production is far from meeting market demand in China. Therefore, breeding of yeast with high RNA content is the basic measure to solve this problem. However, RNA extraction determination procedure limited the implement of above strategy. There is an urgent need to develop effective methods for large scale screening of high nucleic acid yeast.

• Since RNA Pol I does not contain an extra C-terminal domain (CTD), its transcript (rRNA) cannot be added with 5' cap and 3' poly (A) tail, and therefore cannot be translated into proteins. The intergenic region (IGR) IRES (internal ribosome entry site) from *cricket* paralysis virus does not require any translation initiation factors to initiate protein translation.

• In this study, we constructed an IGR IRES-mediated GFP expression system based on RNA Pol I in S. cerevisiae to reflect the intracellular change of RNA content by the fluorescence intensity change of GFP, and established a high throughput screening system based on fluorescent-activated cell sorting (FACS).

#### Results +Uracil -Uracil 1) BXZ01 (pMA91, ura3)



Ideas and design

### • Conclusions

• We designed and constructed of an IGR IRES-mediated reporter gene expression system based on RNA Pol I for large scale screening of highnucleic acid in S. cerevisiae.

• In order to test this expression system working or not, we firstly designed a binary growth assay using the URA3 gene (encoding for 5-phosphate decarboxylase) as the reporter gene. The results indicated that the reporter gene URA3 was successfully expressed and the amount of Ura3p was sufficient to support the growth on an auxotrophic medium with no need to deletion any genes involved in protein translation initiation. • We used another reporter gene (GFP) to replace the original URA3 for high-throughout screening high rRNA synthesis cells. After a round of ARTP mutagenesis to disturb the rRNA synthesis, about 200,000 cells were analyzed using FACS, and 100 cells with fluorescence intensities higher than the threshold were large-scale sorted onto the selection medium. After single yeast separation, twelve pure mutants were finally obtained. Three mutants were randomly selected for further flow cytometry analysis and total RNA content measurement. The results showed that the mutants with higher fluorescence intensity than the threshold showed that RNA content improved by a maximum of 58% without changing its Pol I promoter sequence.

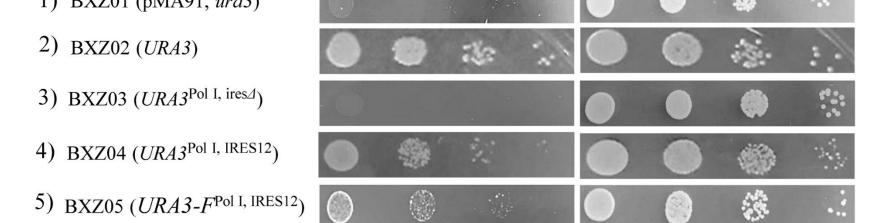


Fig.1. Cells containing Pol I-mediated URA3 expression system grew well on plates minus uracil

2)

amount of Ura3p

expressed by the expression

system based on RNA Pol I was

sufficient to support the growth on

----There is no need to destroy the

yeast cap-dependent translation

initiation mechanism, the reporter

gene URA3 can also expressed

an auxotrophic medium.

BXZ08

Line 1, cells with empty plasmid pMA91 as a negative control. Line 2, cells with Pol Fig.2. Pol I-mediated Ura3p expression was confirmed by western blot. The II-mediated expression cassette (TEF1p-URA3-PGK1t) as a positive control. Line 3, proteins produced by BXZ01 (pMA91, ura3), and BXZ05 (URA3-FPol I, cells containing the Pol I-mediated URA3 expression system with no IRES element <sup>IRES12</sup>) were prepared and incubated with the anti-FLAG tag antibody to showed no growth on the minus uracil plate. Line 4, cells containing the Pol Idetect expression of Ura3p with a C-terminal FLAG tag. mediated URA3 expression system worked well. Line 5, FLAG labeling on the Cterminal end did not impact the Ura3p function.

----The reporter gene URA3 was successfully expressed using the expression system based on RNA Pol I by the growth assay and western blot analysis.

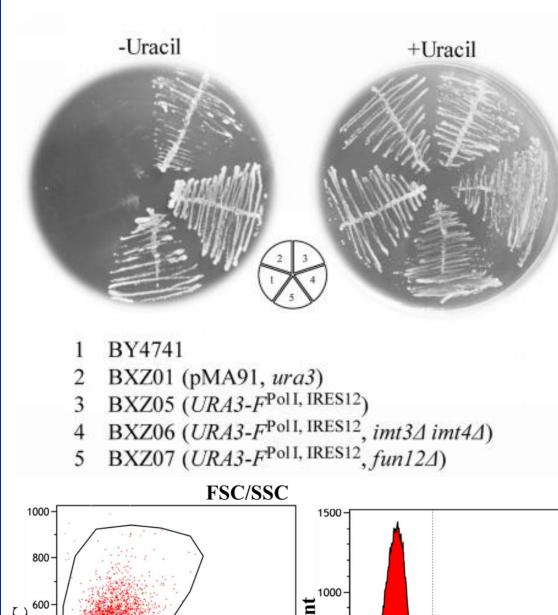


Fig.3. Destroying the cap-dependent factors was translation initiation unnecessary for Ura3p expressing strains. Regions 1 and 2, negative control. Region 3, the Pol I-mediated expression system worked well without destroying translation initiation factors related genes. Region 4, the Pol I-mediated expression system worked well with genes IMT3 and IMT4 deleted. Region 5, the Pol I-mediated expression system worked well with gene

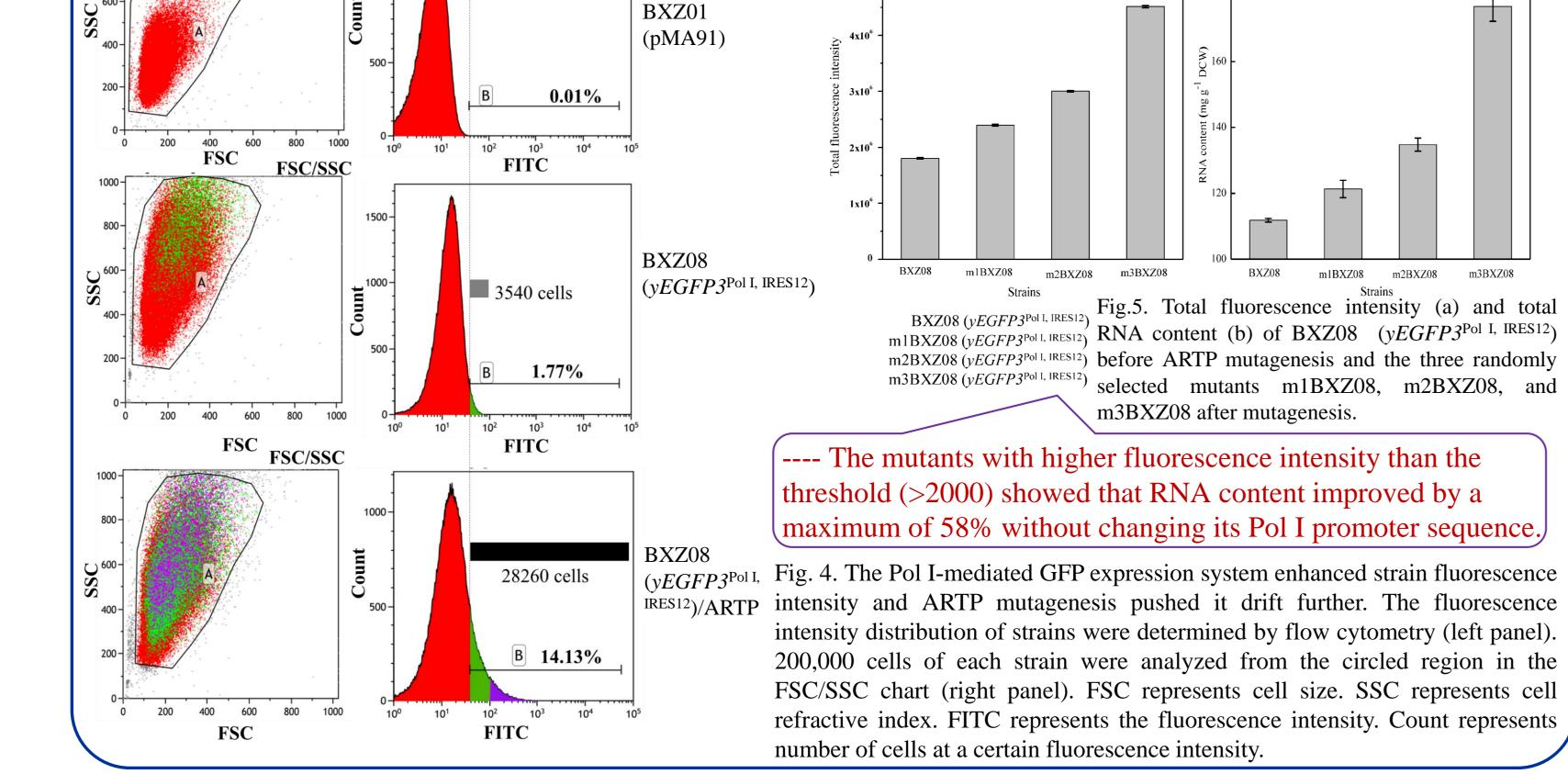
FUN12 deleted. ----Large scale screening of the yeast cells with the fluorescence intensities higher than the threshold using FACS. The histogram of fluorescence intensity and counts for the ARTP mutants changed, which indicated that that ARTP can induce mutation of transcription factors related to rRNA synthesis.

well.

----The

#### • References

- 1. Hodgman CE &, Jewett MC. Characterizing IGR IRES-mediated translation initiation for use in yeast cell-free protein synthesis. N Biotechnol. 2014, 31(5):499-505.
- 2. Thompson SR, Gulyas initiation in Sarnow P. KD, Internal Saccharomyces cerevisiae mediated by an initiator tRNA/eIF2independent internal ribosome entry site element. Proc Natl Acad Sci U S A. 2001, 98(23):12972-7.



- 3. Sasano Y, Kariya T, Usugi S, Sugiyama M, Harashima S. Molecular breeding of Saccharomyces cerevisiae with high RNA content by harnessing essential ribosomal RNA transcription regulator. AMB Express. 2017, 7(1):32.
- 4. Li B, Liu Y, Wang L, Hong J, Chen Y, Ying H. RNA accumulation in Candida tropicalis based on cofactor engineering. FEMS Yeast Res. 2019, 19(3).
- 5. Thompson SR. Tricks an IRES uses to enslave ribosomes. Trends Microbiol. 2012, 20(11):558-66.

## Acknowledgements

This work was supported by the National Natural Science Foundation of China (31801515), Shandong Provincial Natural Science Foundation (ZR2019PC010), the Key R&D Program of Shandong Province (2017CXGC1105).