

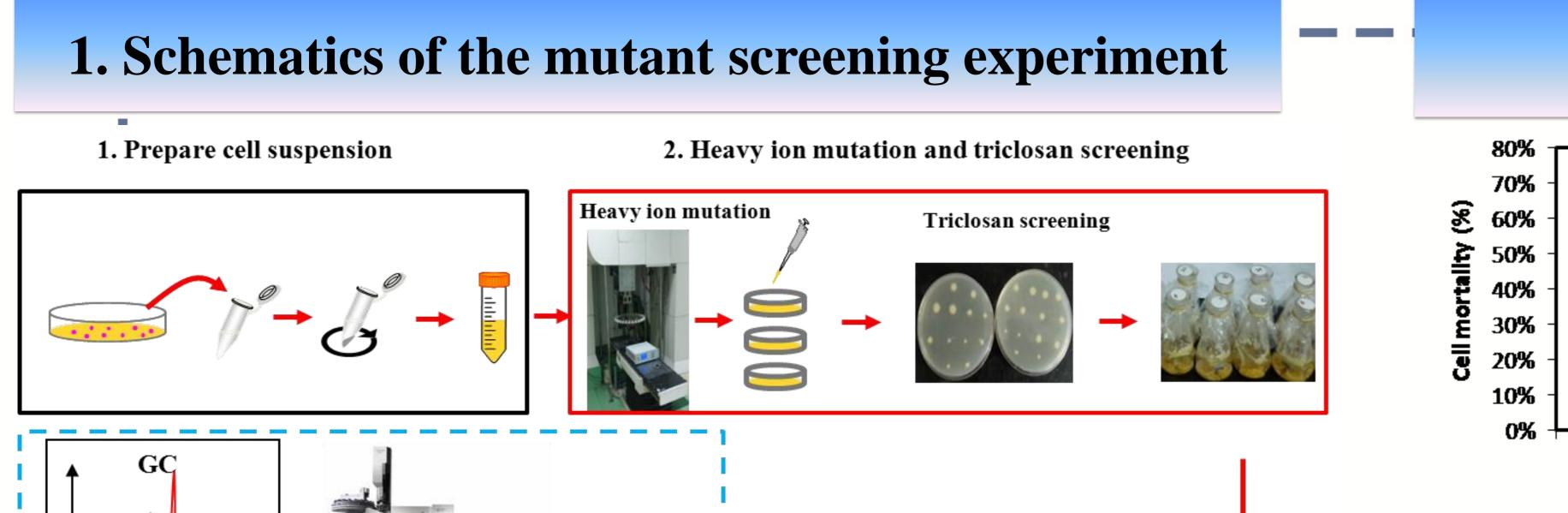
Heavy ion mutagenesis combined with triclosan screening provides a new strategy for improving the arachidonic acid yield in Mortierella alpina

Huidan Zhang, Dong Lu, Xin Li, Yingang Feng, Qiu Cui, Xiaojin Song*

CAS Key Laboratory of Biofuels and Shandong Key Laboratory of Synthetic Biology, Qingdao Institute of Bioenergy and Bioprocess Technology Chinese Academy of Sciences

Abstract

Arachidonic acid (ARA), which is a ω -6 polyunsaturated fatty acid, has a wide range of biological activities and is an essential component of cellular membranes in some human tissues. Mortierella alpina is the best strain for industrial production of ARA. To increase its yield of arachidonic acid, heavy ion beam irradiation mutagenesis of Mortierella alpina was carried out in combination with triclosan and octyl gallate treatment. The obtained mutant strain F-23 ultimately achieved an ARA yield of 5.26 g L⁻¹, which is 3.24 times higher than that of the wild-type strain. In addition, quantitative real-time PCR confirmed that the expression levels of fatty acid synthase (FAS), $\Delta 5$ -desaturase, $\Delta 6$ -desaturase, and $\Delta 9$ -desaturase were all significantly up-regulated in the mutant F-23 strain, especially $\Delta 6$ - and $\Delta 9$ -desaturase, which were up-regulated 3- and 2-fold, respectively. This study confirmed a feasible mutagenesis breeding strategy for improving ARA production and provided a mutant of Mortierella alpina with high ARA yield.



Octyl gallate sreening

gallate sreening

3. 5-Fluorouracil mutagenesis and octyl

5-FU mutagenesis

2. Screening of high-yield ARA mutants with triclosan and octyl gallate

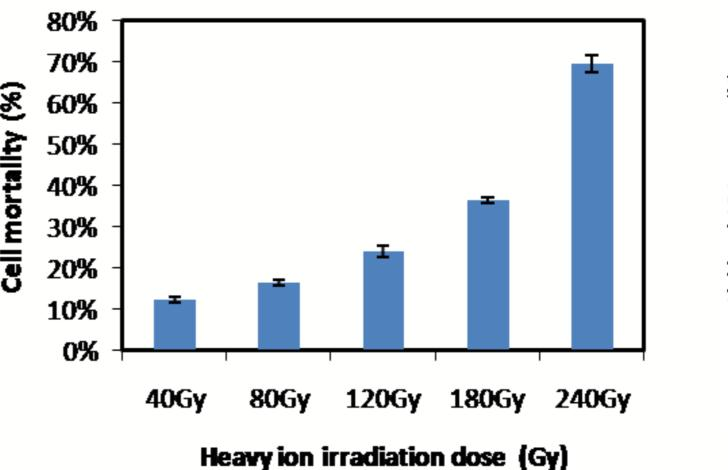


Fig.2 Mortality of *M.alpina*

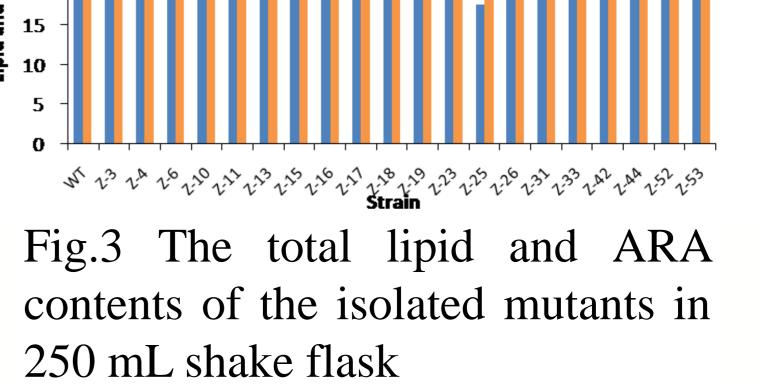
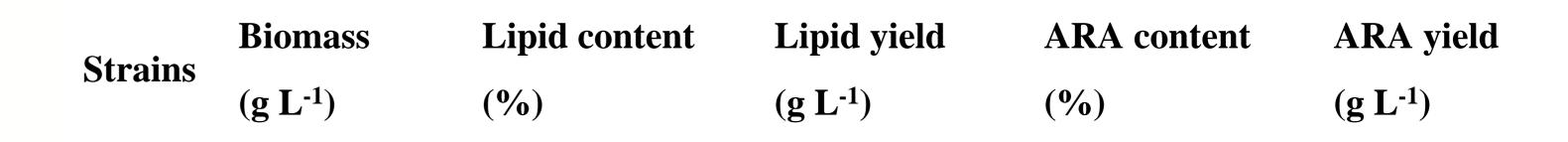


Table 1.Productivity of wild-type strain, mutant strain Z-44 and F-23





qRT-PCR

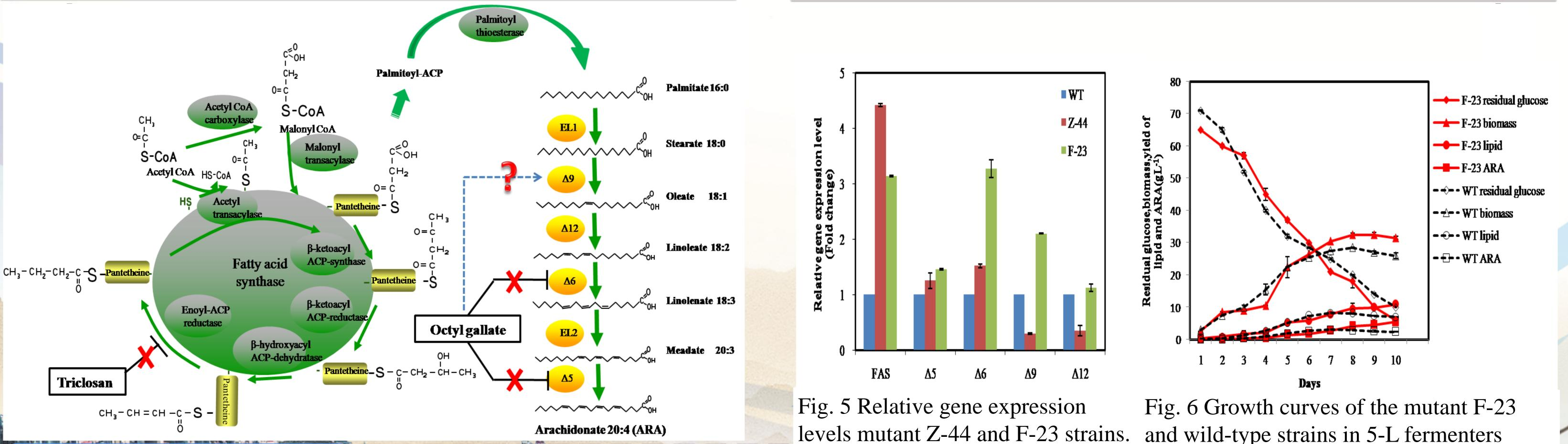
4. Verification and analysis of fermentor

Fig.1 Shematics of the mutant screening experiment

3.The ARA biosynthesis pathway and the action sites of inhibitors in Mortierella alpina

1.24 ± 0.18 WT 19.8 ± 0.36 4.60 ± 0.56 27.04 ± 0.63 23.24 ± 0.46 25.2 ± 0.93 38.3 ± 0.97 3.80 ± 0.65 **Z-44** 9.65 ± 0.26 39.37 ± 0.28 28.2 ± 0.36 38.0 ± 0.25 10.72 ± 0.41 49.08 ± 0.17 5.26 ± 0.48 **F-23**

4.Genetic stability and batch fermentation of mutant F-23



and wild-type strains in 5-L fermenters

5.Conclusions

 \bullet In this study, mutant F-23 was selected after heavy ion beam irradiation combined with triclosan and octyl gallate treatment. Compared with the wild-type strain, the total lipid and ARA yields were increased by 1.33 and 3.24 times, respectively.

•These results provide a good method and strategy for screening microorganisms with high yield of unsaturated fatty acids.

Authors: Name: Huidan Zhang; Xiaojin Song E-mail:zhanghd@qibebt.ac.cn; songxj@qibebt.ac.cn **Tel:**+86 532 80662705; **Fax:** +86 532 80662707; Post Address: No.189 Songling Road, Laoshan District, Qingdao 266101, Shandong Province, China