

Biosynthesis of P(3HB-co-LA) from Renewable Carbon Source by Metabolically Engineered *Escherichia coli*

Xiangju Wei, JingXian Lu , Ju Wu , Zhimin Li, Qin Ye, Hui Wu *

State Key Laboratory of Bioreactor Engineering, East China University of Science and Technology, Shanghai 200237

E. mail: hwu@ecust.edu.cn

Abstract

P(LA-co-3HB) has been considered as a good alternative to petroleum-based plastic as it possesses several desirable properties such as biodegradability, biocompatibility, compostability, and low toxicity to humans. P(LA-co-3HB) can be produced in engineered *Escherichia coli* harboring the genes encoding LA-polymerizing enzyme (LPE) and monomer-supplying enzymes. In this study, the LA content in the polymer was increased through the strategies of weakened respiratory chain, promoter engineering and carbon source optimization. The results showed that weakened respiratory chain level can increase the component of LA in P(3HB-co-LA) from 5.1 mol% to 14.1 mol%; After optimization the expression of pct_{th} via promoter engineering, the LA component in the polymer can be further increased to 19.5 mol%. When the xylose was used instead of glucose as carbon, the LA content in wild strain was increased significantly, the highest LA content reached 30.6 moL%. Then, mixed sugar (glucose: xylose = 7:3) and lignocellulosic hydrolysate were used as carbon resource, respectively, the LA component of the polymer is improved. This study suggested that lignocellulose as a abundant and low-cost source can be used to produce the Lactic acid-based polymer.

Introduction

Strategy

Utilization of promoter engineering of *pct_{th}* to improve P(LA-co-3HB) biosynthesis



Resource saving and environmentfriendly.



- ✓ Xylose as carbon can increase the LA content of P(LA-co 3HB).
- lignocellulosic hydrolysate from the maize straw has the potential to be anabundant and inexpensive source of xylose for the production of value-added products.



Weaken respiratory chain:

deletion of UbiX to decrease the co-enzyme Q8 level, weakened respiratory chain, increase lactic acid content

> **Promoter engineering:** strengthen the transcription of the key gene pct_{th}

Carbon source optimization:

lignocellulosic hydrolysate, an agricultural waste, as a anabundant and inexpensive source advantageous for lactic acid-based polymer

✓ Effect of *trc* promoter





When the *trc* promoter was was used to express pct_{th} , the LA content of JX041-03 reached **19.5mol%**.

Biosynthesis of P(LA-co-3HB) through different carbon resources

✓ Effect of xylose

Strain	Carbon	Polymer content	LA fraction
	source	(wt%)	(mol%)
MG-01	Glucose	76.3	5.1
	xylose	76.1	24.5
JX04-01	Glucose	79.2	9.1
	xylose	38.0	2.7
JX041-01	Glucose	81.7	14.1
	xylose	28.6	7.5
MG-02	Glucose	82.2	(3.6
	xylose	70.1	6.3
JX04-02	Glucose	62.9	6.9
	xylose	42.7	2.5
JX041-02	Glucose	64.0	13.4
	xylose	33.6	7.9
MG-03	Glucose	83.1	(10.0)
	xylose	72.0	30.6
JX04-03	Glucose	70.0	5.9
	xylose	18.8	5.5
JX041-03	Glucose	69.8	19.5
	xylose	22.7	11.3

 ✓ Effect of glucose and xylose as co-carbon resource (G:X=7:3)

Strain	Carbon	Polymer	LA fraction
	source	content (wt%)	(mol%)
	Glucose	76.3	5.1
MG-01	mixture	72.0	12.0
IV 04 01	Glucose	79.2	9.1
JX04-01	mixture	38.4	10.3
IV0/1 01	Glucose	81.7	14.1
JA041-01	mixture	35.5	13.0
	Glucose	82.2	3.6
WIG-02	mixture	90.5	1.5
JX04-02	Glucose	62.9	6.9
	mixture	42.0	8.3
JX041-02	Glucose	64.0	13.4
	mixture	26.8	15.3
MG-03	Glucose	83.1	10.0
	mixture	88.0	5.4
JX04-03	Glucose	70.0	5.9
	mixture	53.7	8.8
JX041-03	Glucose	69.8	19.5
	mixture	51.5	13.3

Strains and plasmids

lS
opionicun
lium
ım

When xylose was used as carbon resource, the wild type strains' LA content of copolymer has been improved obviously, especially MG-03 reached **30.6mol%**. When the mixture(glucose:xylose=7:3) was used, most strains' LA content in the copolymer were increased.

Biosynthesis of P(LA-co-3HB) through lignocellulosic hydrolysate

Strain	Carbon	Polymer	LA fraction
	source	content(wt%)	(mol%)
MG-01	Glucose	76.3	5.1
	Hydrolysate	91.2	6.9
JX04-01	Glucose	79.2	9.1
	Hydrolysate	37.4	24.6
JX041-01	Glucose	81.7	14.1
	Hydrolysate	31.5	25.0



Results

Effect of weakening the respiratory chain on P(LA-co-3HB) biosynthesis





Deletion of *ubiX* and *dld* showed a positive effect on lactate content, the JX041-01 produced P(3HBco-14.1mol%LA), the lactate content was 2.76-fold of that produced by MG-01.

Acknowlegments This study was supported by the National Natural Science Foundation of China (Grant No. 21776083), and the Fok Ying-Tong Education Foundation, China (Grant No. 161017). The lignocellulosic hydrolysate is mainly composed with gluose(69%) and xylose(31%), results shows it is advantageous for producing the lactic acid-based polymer.

Conclusions

- ✓ Construct the heterologous synthetic pathway of P(3HB-co-5.1mol%LA)
- ✓ Weaken the respiratory chain, the production of P(3HB-co-14.1mol%LA)
- ✓ Strengthen the promoter of the key enzyme, the production of P(3HB-co-19.5mol%LA)
- ✓ After the carbon optimization, the production of P(3HB-co-30.6mol%LA)
- \checkmark Innovatively used the lignocellulosic hydrolysate, an agricultural waste, we synthesize degradable,

biocompatible copolymer

References

[1] Yang T H , et al. Biotechnology & Bioengineering, 2010, 105(1):150-160.
[2] Wu H , et al. Metabolic Engineering, 2015, 28:159-168.
[3] Choi S Y , et al. Nature Biotechnology, 2016, 34:435-441.