

Enhanced Rolling Circle Amplification Performance with a Newly Engineered phi29 DNA Polymerase

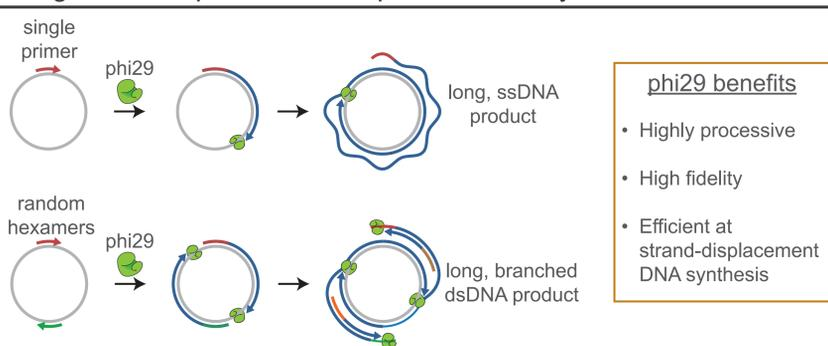
Kyle B. Vrtis, Yanxia Bei, Sean Lund, Léa Chuzel, Yinhua Zhang, Nathan A. Tanner, Jennifer L. Ong, Gregory C. Patton, Nicole M. Nichols

New England Biolabs, Inc., Ipswich, MA 01938, USA

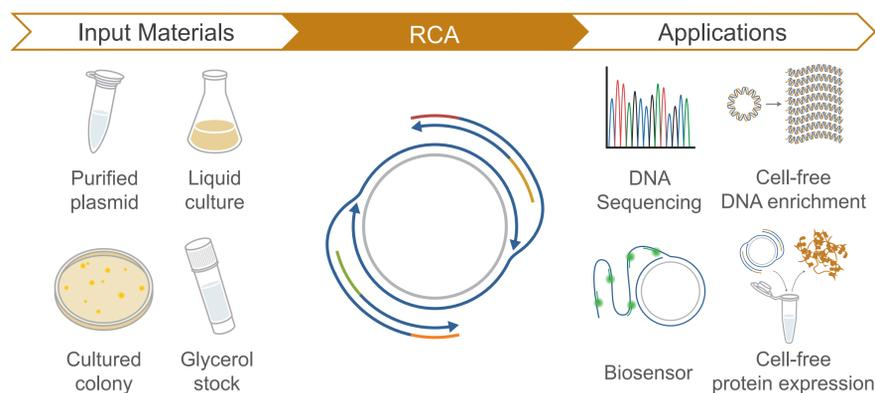
Introduction

Isothermal DNA amplification techniques exponentially increase the amount of DNA in a sample without the constraint of thermocycling. New England Biolabs (NEB) offers a wide range of innovative solutions in isothermal amplification. One of them, Rolling Circle Amplification (RCA) has become an invaluable tool for many biotechnology applications including: DNA sequencing, cell-free DNA synthesis, cell-free protein expression, biomolecular detection, and DNA hydrogel generation. The most widely used DNA polymerase for RCA is phi29, which possesses high processivity and fidelity, and strong strand-displacement activity. In the recently launched NEB phi29-XT RCA kit, we have incorporated an engineered phi29 DNA polymerase, phi29-XT, that shares the positive characteristics of the wild-type phi29, but also generates higher yield in less time than the wild-type enzyme. Additionally, while wild-type phi29 works optimally at 30°C, phi29-XT performs best at 42°C. We have used this kit downstream of DNA assembly technologies, such as GoldenGate and NEBuilder HiFi Assembly, to perform high-throughput cell- and vector-free protein expression. Furthermore, the kit also enabled high-throughput fosmid amplification directly from bacterial cells, followed by ONT sequencing for *de novo* fosmid DNA assembly. Taken together, these applications demonstrate the utility of the phi29-XT RCA kit for isothermal amplification of circular DNA templates.

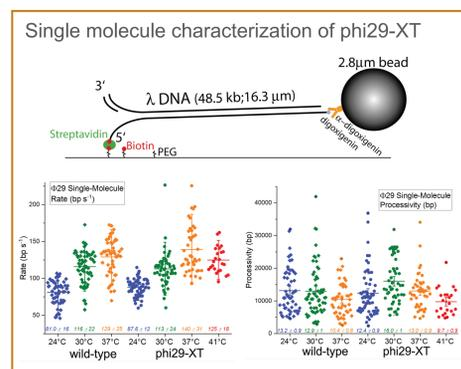
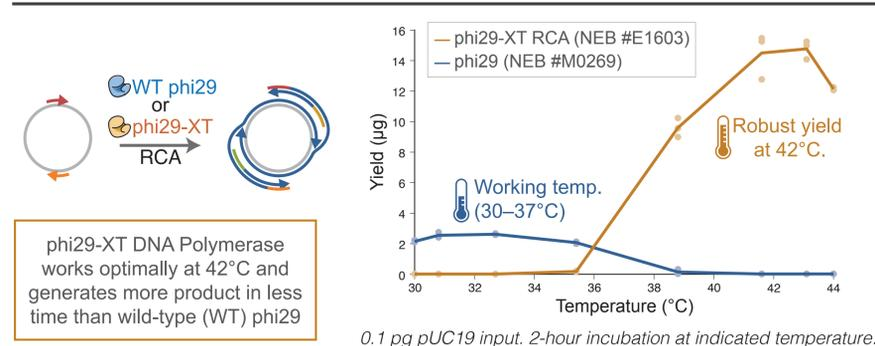
Rolling Circle Amplification with phi29 DNA Polymerase



Highly versatile RCA assay



phi29-XT DNA Polymerase performance

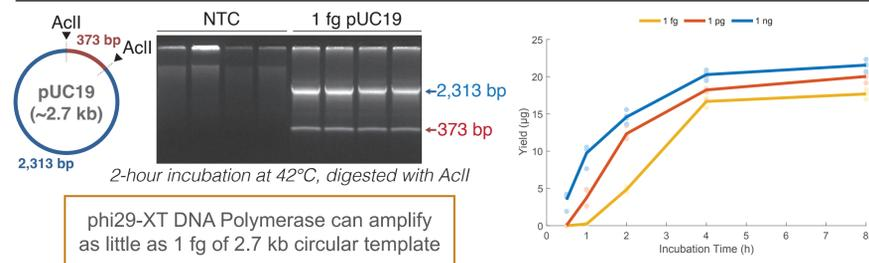


DNA Polymerase	Error Rate	Fidelity, Relative to Taq
Taq	1.6×10^{-4}	1
Wild-type phi29	1.5×10^{-5}	10.7
phi29-XT	9.0×10^{-6}	17.8

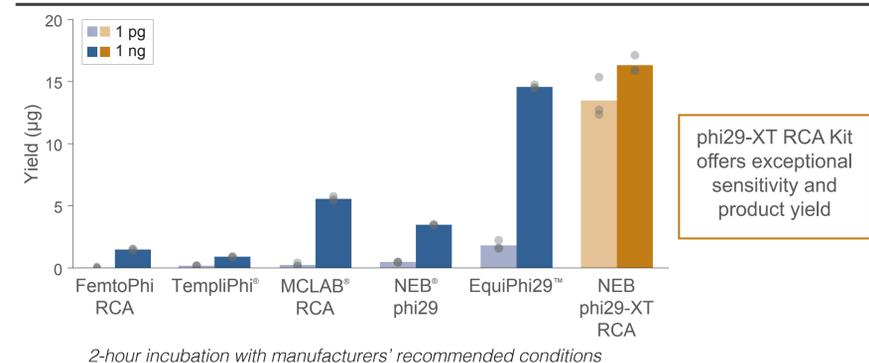
phi29-XT synthesizes with a rate, processivity, and fidelity comparable to wild-type phi29 DNA polymerase

Note: Single molecule and fidelity experiments were not carried out with phi29-XT reaction buffer

phi29-XT DNA Polymerase sensitivity

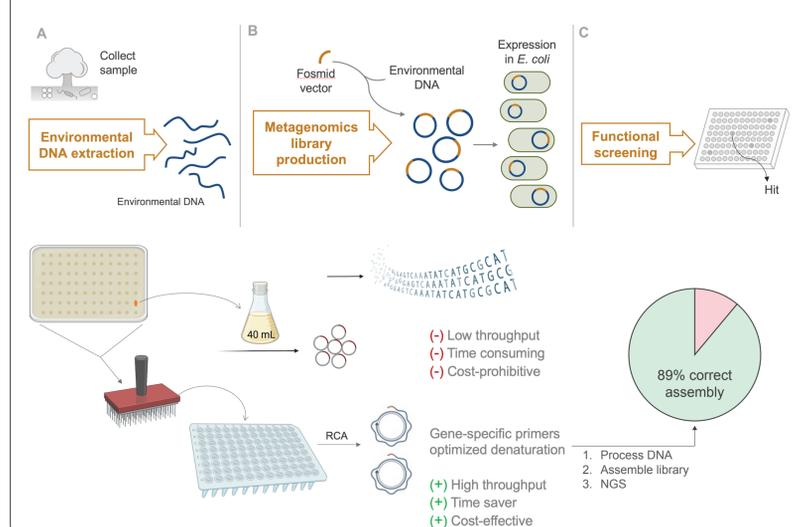


Comparison of phi29-XT to other commercially available phi29s

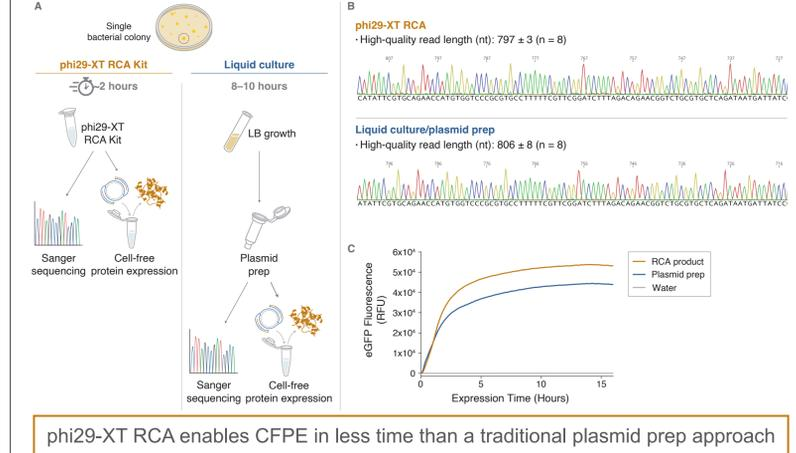


phi29-XT DNA Polymerase-mediated applications

phi29-XT RCA enables high-throughput metagenomics screening



Direct colony RCA enables cell-free protein expression (CFPE)



Conclusions

- The newly engineered phi29 DNA polymerase, phi29-XT, has improved sensitivity and RCA yield over the wild-type phi29 DNA polymerase
- The NEB phi29-XT RCA Kit (E1603) is highly versatile, fast, and simple-to-use

References

- phi29-XT RCA Kit. www.neb.com/E1603
- Potapov V, Ong JL (2017) PLoS ONE 12(1): e0169774.
- Tanner, N, van Oijen, A (2010) Methods in Enzymology 475: 259-278
- Chuzel, L., et al., (2018) J Biol Chem 293:18138-18150