

Optimized IVT Platforms: Advancing RNA Quality and Scalability for Animal Biologics Innovation

Hélène Boyer¹, Laurence Delaurière¹, Dana Brecklin-Benassi¹, Daniela Torres Campana¹, Laura Alexander¹, Maria Dashek¹

¹Promega Corporation, 2800 Woods Hollow Rd, Madison, WI 53711



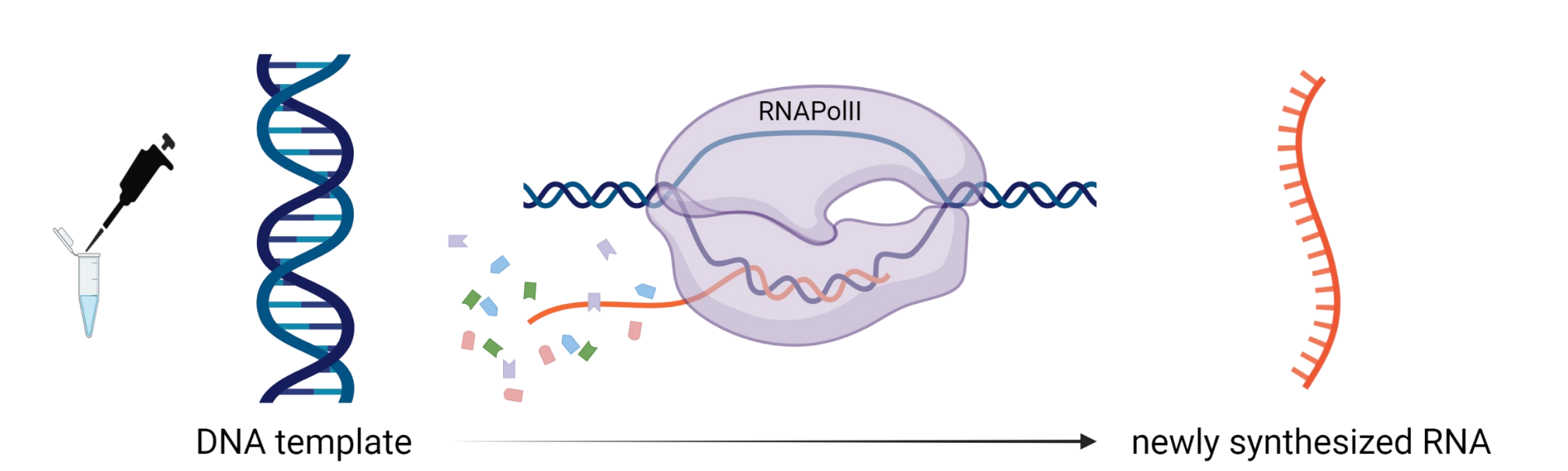
1. Introduction: Tackling the Challenges in Animal Biologics

In the rapidly evolving field of animal biologics development, producing high-quality RNA is essential for developing next-generation vaccines and therapeutic interventions. Advanced in vitro transcription (IVT) methodologies provide researchers and process engineers with scalable options to generate large quantities of high-integrity RNA. From efficient transcription systems to precise clean-up and automated workflows, these approaches address the rigorous demands of animal health applications by minimizing immunogenic risk and ensuring consistent results.

2. Advanced RNA Production Technologies

The RiboMAX™ Large Scale RNA Production Systems enable the production of **milligram amounts of RNA** from a linear DNA template. They are typically used to produce transcripts up to **5-6kb** in size, although much longer transcripts have been successfully generated.

Capped and uncapped RNA transcripts can be produced.



- Main Applications:**
- In vitro translation
 - In vivo translation
 - Synthesis of tRNA, rRNA, siRNA...
 - Synthesis of RNA virus genomes
 - Synthesis of ribozymes
 - Synthesis of guide RNA for CRISPR gene editing¹
 - cDNA library preparation²
 - Study of RNA:protein interaction^{3,4}
 - Study of RNA splicing
 - Drug discovery
 - mRNA therapeutics/vaccine
 - ...

3. Ensuring RNA Integrity and Purity: Comparative Clean-Up Solutions

ReliaPrep™ RNA Clean-Up and Concentration System
After in vitro transcription (20µl reaction volume), RNA was cleaned-up using either the standard protocol recommended in the technical manual (Phenol chloroform; N=3) or using the ReliaPrep™ RNA Clean-Up and Concentration System (N=4).
→ *Yields were lower with the ReliaPrep™ system, but RNA integrity remained comparable across methods.*

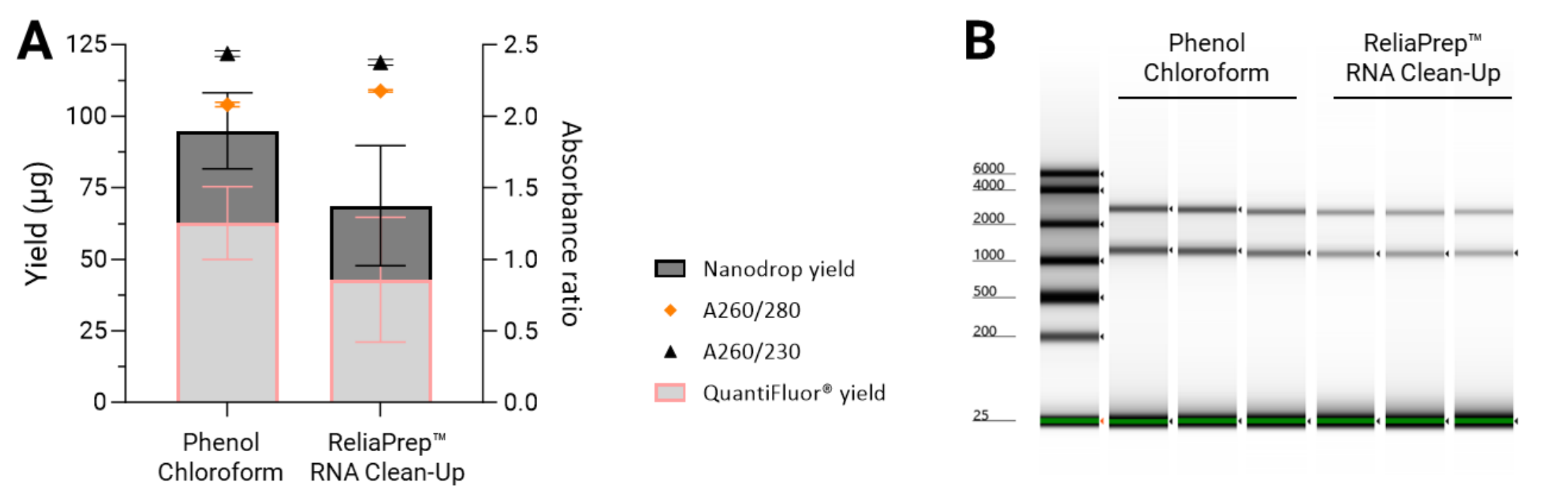


Figure 1. Comparison of phenol chloroform and ReliaPrep™ RNA Clean-Up methods. **A.** Yields (in µg) expressed as mean ± standard deviation, determined using Nanodrop™ or Quantifluor® RNA System (left y axis). Absorbance ratios (mean ± standard deviation) are represented on the right y axis. **B.** RNA integrity was assessed using an RNA ScreenTape.

Maxwell® RSC miRNA Tissue Kit
After in-vitro transcription (100µl reaction volume), RNA was cleaned-up using the Maxwell® RSC miRNA Tissue Kit, from an input volume of 20µl, 50µl or 100µl (N=3 for each).
→ *Under comparable conditions, yields matched those from the phenol chloroform protocol, with linear scalability from 20 to 100µl.*

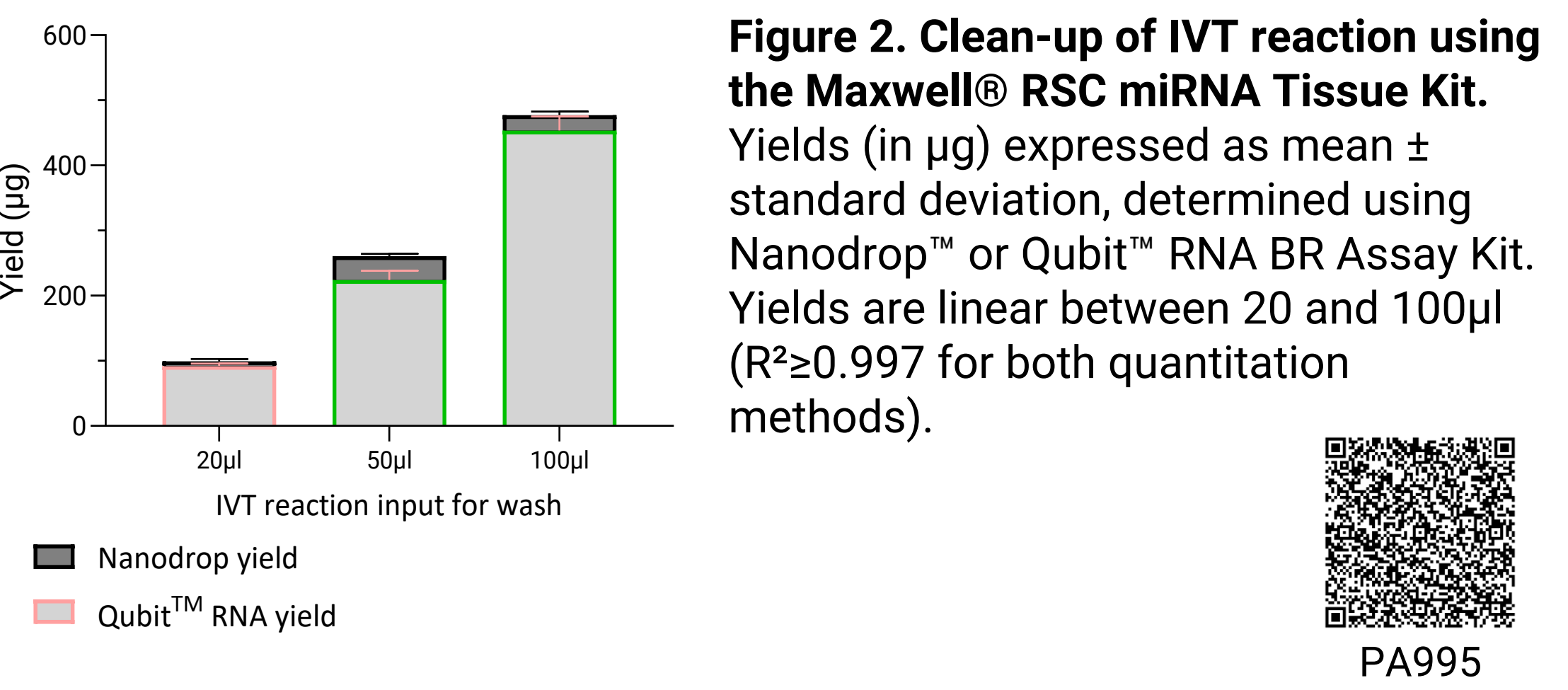
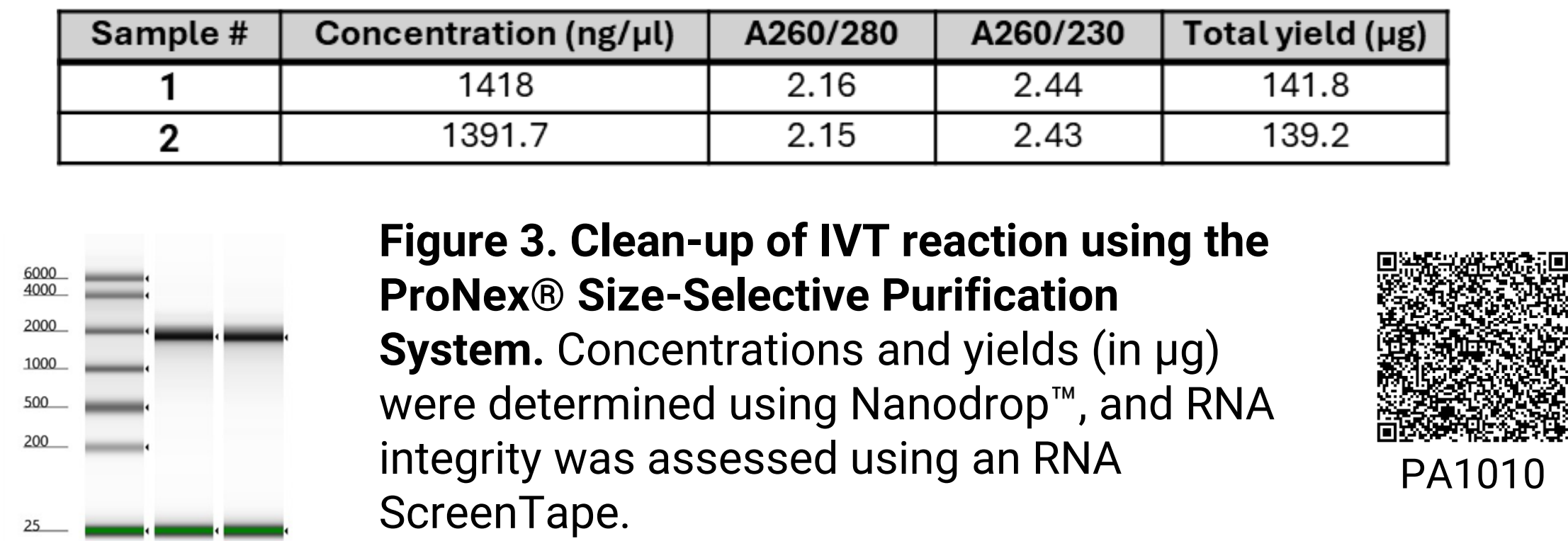


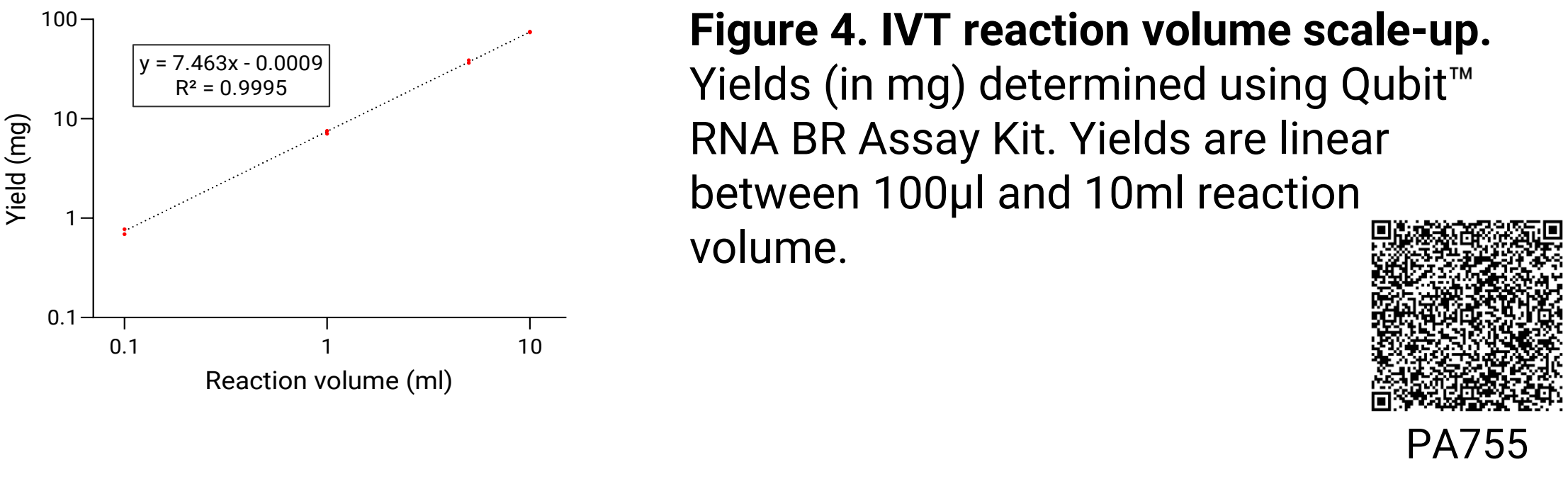
Figure 2. Clean-up of IVT reaction using the Maxwell® RSC miRNA Tissue Kit. Yields (in µg) expressed as mean ± standard deviation, determined using Nanodrop™ or Qubit™ RNA BR Assay Kit. Yields are linear between 20 and 100µl (R²≥0.997 for both quantitation methods).

ProNex® Size-Selective Purification System
After in vitro transcription (20µl reaction volume), RNA was cleaned-up using the ProNex® Size-Selective Purification System (N=2).
→ *The ProNex® Size-Selective Purification System produced higher yields than phenol chloroform, with good RNA integrity.*



4. Innovative Workflows: From Circular RNA to Long ssDNA

In vitro transcription reaction volume scale-up
RNAs were generated by in vitro transcription in a 100µl, 1ml, 5ml or 10ml reaction volume (N=2 for each), and dye-based quantified.
→ *In vitro transcription can be scaled to 10 ml while maintaining consistent RNA yields.*



Synthesis of circular RNA using the RiboMAX™ System
Circular RNAs can be synthesized in vitro as shown by sequencing of the circRNA junction and ScreenTape analysis.
→ *Reaction products are a mix of circRNA and linear precursors, which can be isolated by gel purification or RNase R digestion.*

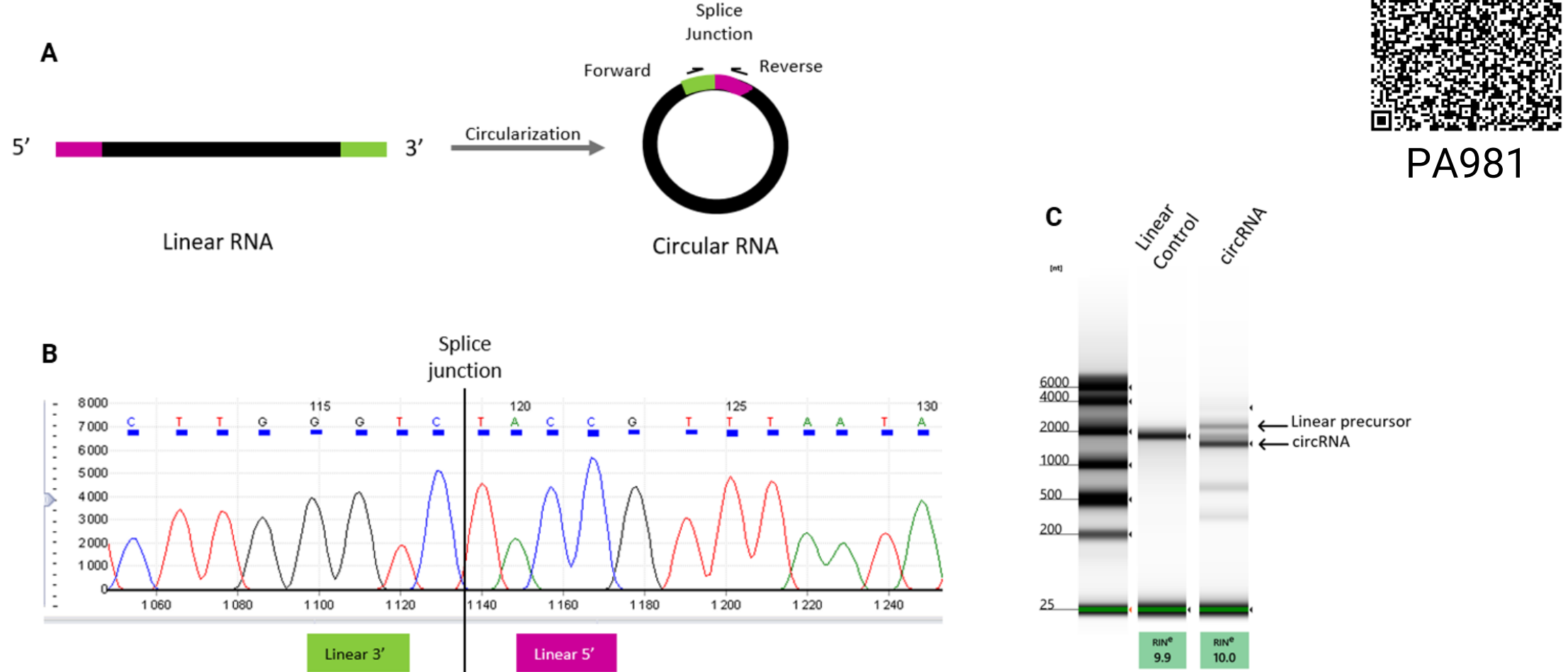


Figure 5. circRNA synthesis. **A.** Schematic Representation of the junction and primers used for sequencing. **B.** Electropherogram obtained by Sanger sequencing confirming the RNA circularization. **C.** RNA circularization was visualized using an RNA ScreenTape.

Workflow: long ssDNA template synthesis for CRISPR insertion
Long ssDNA can be synthesized using a multistep workflow that includes in vitro transcription, followed by appropriate purification and reverse transcription steps.

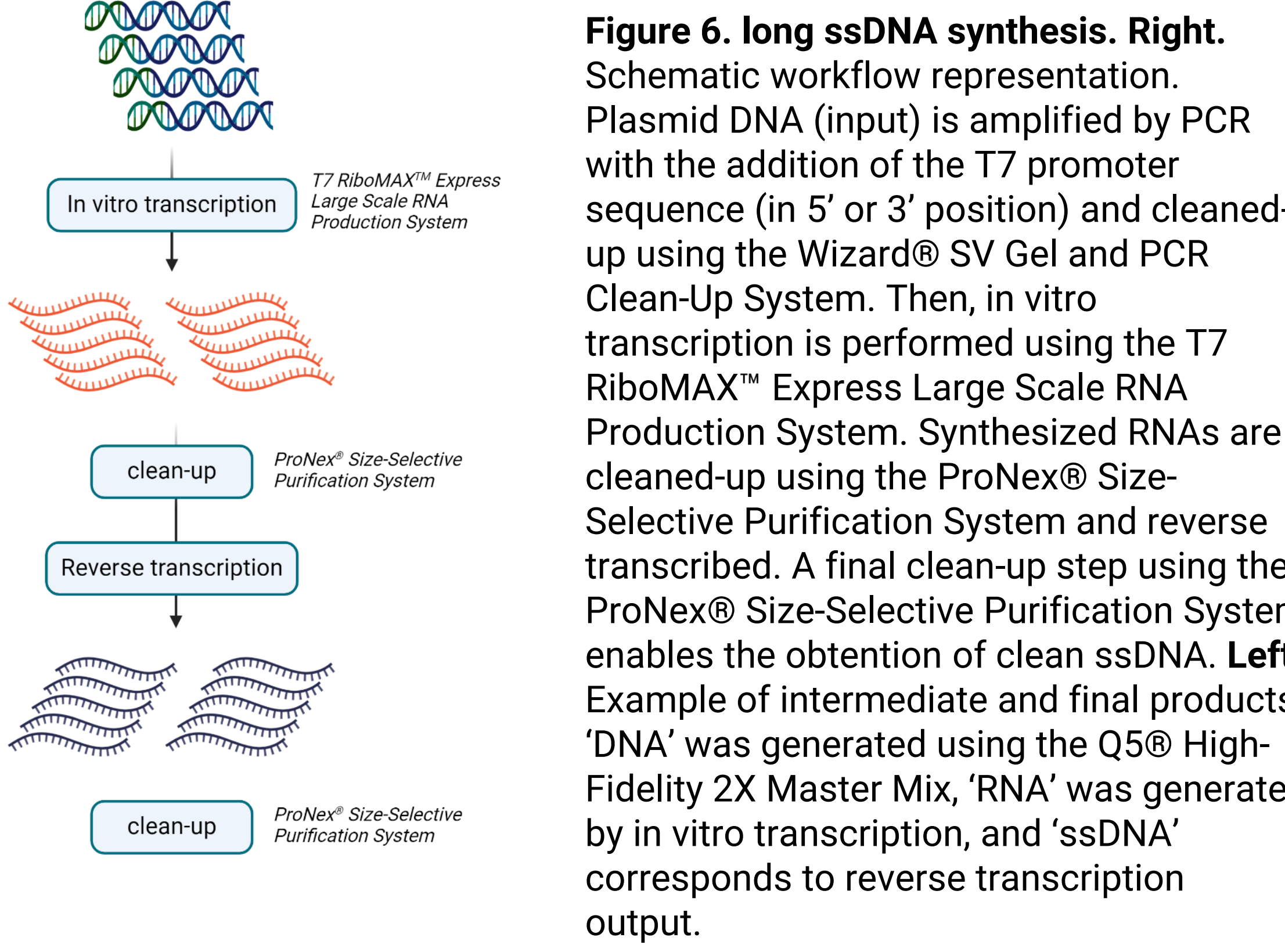
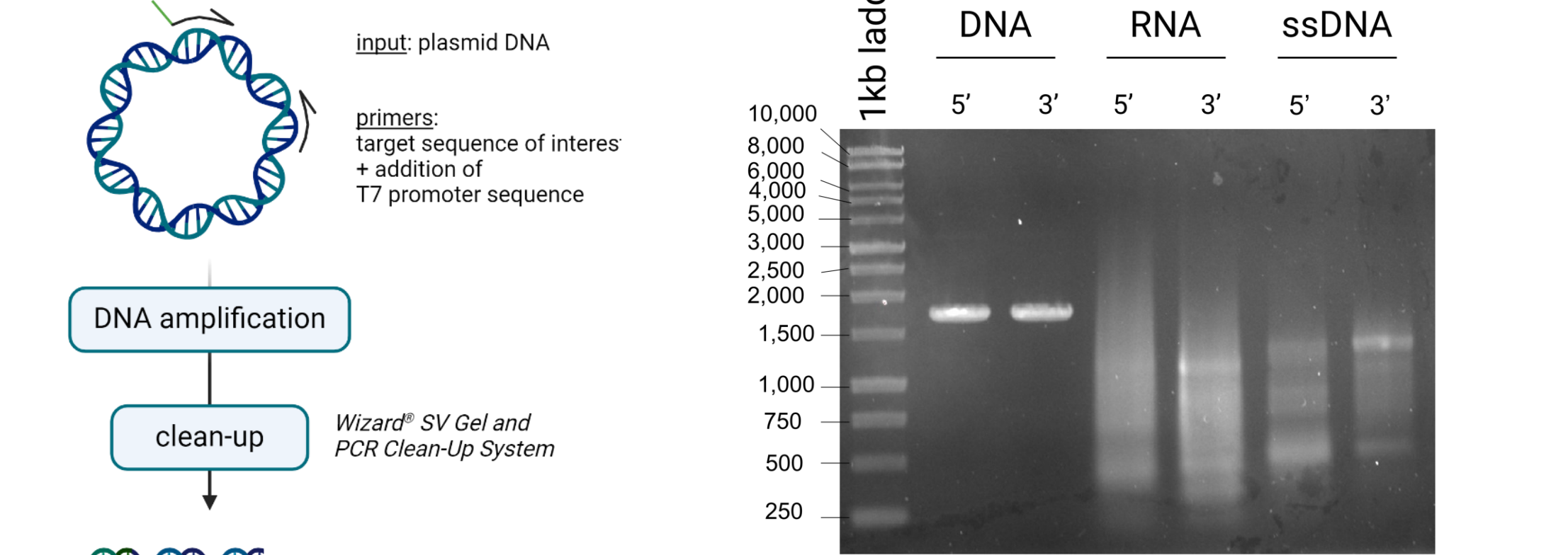
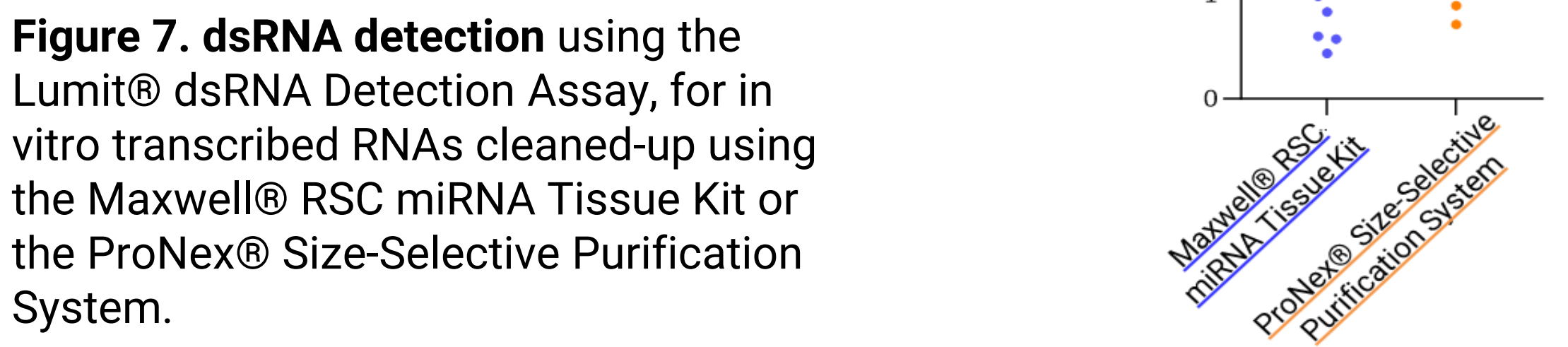


Figure 6. long ssDNA synthesis. **Right.** Schematic workflow representation. Plasmid DNA (input) is amplified by PCR with the addition of the T7 promoter sequence (in 5' or 3' position) and cleaned-up using the Wizard® SV Gel and PCR Clean-Up System. Then, in vitro transcription is performed using the T7 RiboMAX™ Express Large Scale RNA Production System. Synthesized RNAs are cleaned-up using the ProNex® Size-Selective Purification System and reverse transcribed. A final clean-up step using the ProNex® Size-Selective Purification System enables the obtention of clean ssDNA. **Left.** Example of intermediate and final products. 'DNA' was generated using the Q5® High-Fidelity 2X Master Mix, 'RNA' was generated by in vitro transcription, and 'ssDNA' corresponds to reverse transcription output.

5. Quality Control: Reducing Immunogenicity with dsRNA Detection

dsRNA is a common by-product of in vitro transcription responsible for immunogenicity and decreased protein translation efficiency⁵. The Lumit® dsRNA Detection Assay was developed to quantitate dsRNA in biological samples.



6. Automation for Consistency and Scale

Automation streamlines production, reduces hands-on time, and ensures reproducibility in process development. An automated method has been developed on the CyBio Felix Liquid Handler from Analytik Jena, enabling the automation of the synthesis of RNA from a linear DNA input.

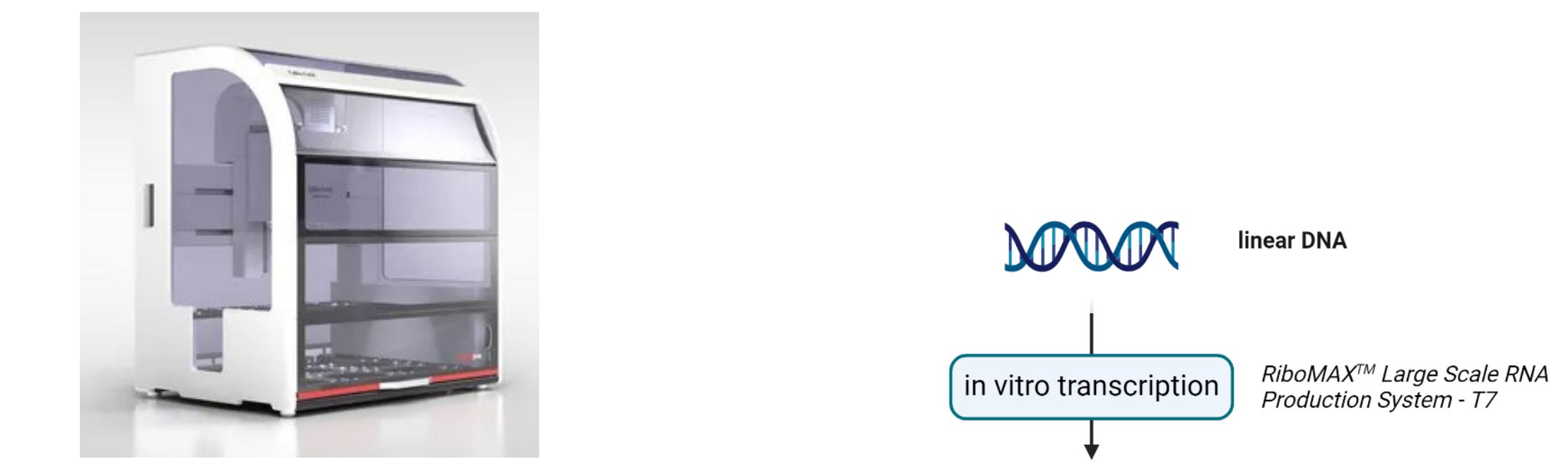


Figure 8. Automated synthesis of RNA from a linear DNA input. Schematic representation of the automated workflow. Linear DNA is used as input. In vitro transcription is carried out using the RiboMAX™ Large Scale RNA Production System T7 (automated dispensing of the reaction mix onto the linear input plate). The synthesized RNA is cleaned-up using the ProNex® Size-Selective Purification System.

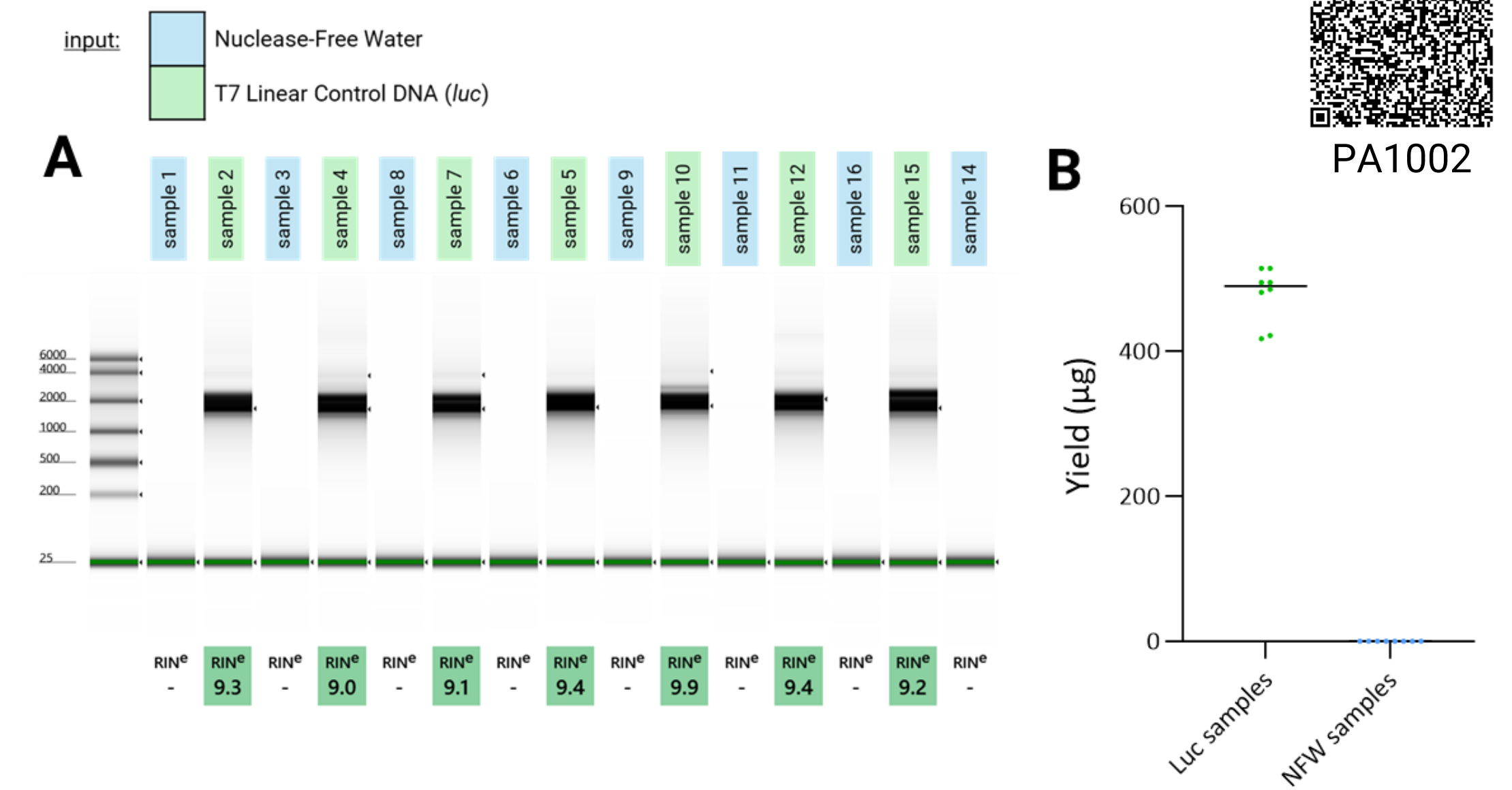


Figure 9. Automated transcription and clean-up on the CyBio Felix. Control RNA was generated and cleaned-up according to the workflow represented above in an automated fashion on the CyBio Felix, according to a checkerboard pattern (linear control RNA and Nuclease-Free Water were used as input). Eluates were analyzed using the TapeStation 4200 and quantified using the Qubit™ RNA HS Assay Kit on the Quantus™ Fluorometer.

→ *RNA was successfully produced in an automated fashion on the CyBio Felix, with no cross contamination detected.*

7. Conclusions

- Unlock scalable, high-quality RNA production tailored for animal biologics development.
- Leverage advanced clean-up methods, dsRNA detection, and automation to streamline workflows and boost reproducibility.
- Drive innovation in animal biologics development through these integrated technologies.

8. References

1. Bai, H. *et al.* (2020) CRISPR/Cas9-mediated precise genome modification by a long ssDNA template in zebrafish. *BMC Genomics* **21**, 67.
2. Fujimori, S. *et al.* (2012) Next-generation sequencing coupled with a cell-free display technology for high-throughput production of reliable interactome data. *Sci Rep* **2**, 691.
3. Tang, S.J. *et al.* (2020) Cis- and trans-regulations of pre-mRNA splicing by RNA editing enzymes influence cancer development. *Nat Commun* **11**, 799.
4. Wang, S. *et al.* (2019) PbTTG1 forms a ribonucleoprotein complex with polypyrimidine tract-binding protein PbPTB3 to facilitate the long-distance trafficking of PbWoxT1 mRNA. *Plant Sci.* **280**, 424–32.
5. Xin M, *et al.* (2018). An origin of the immunogenicity of in vitro transcribed RNA. *Nucleic Acids Research*, **14**, issue 10, 5239-5249