



Optimization of cfRNA detection via sequencing: Addressing workflow challenges

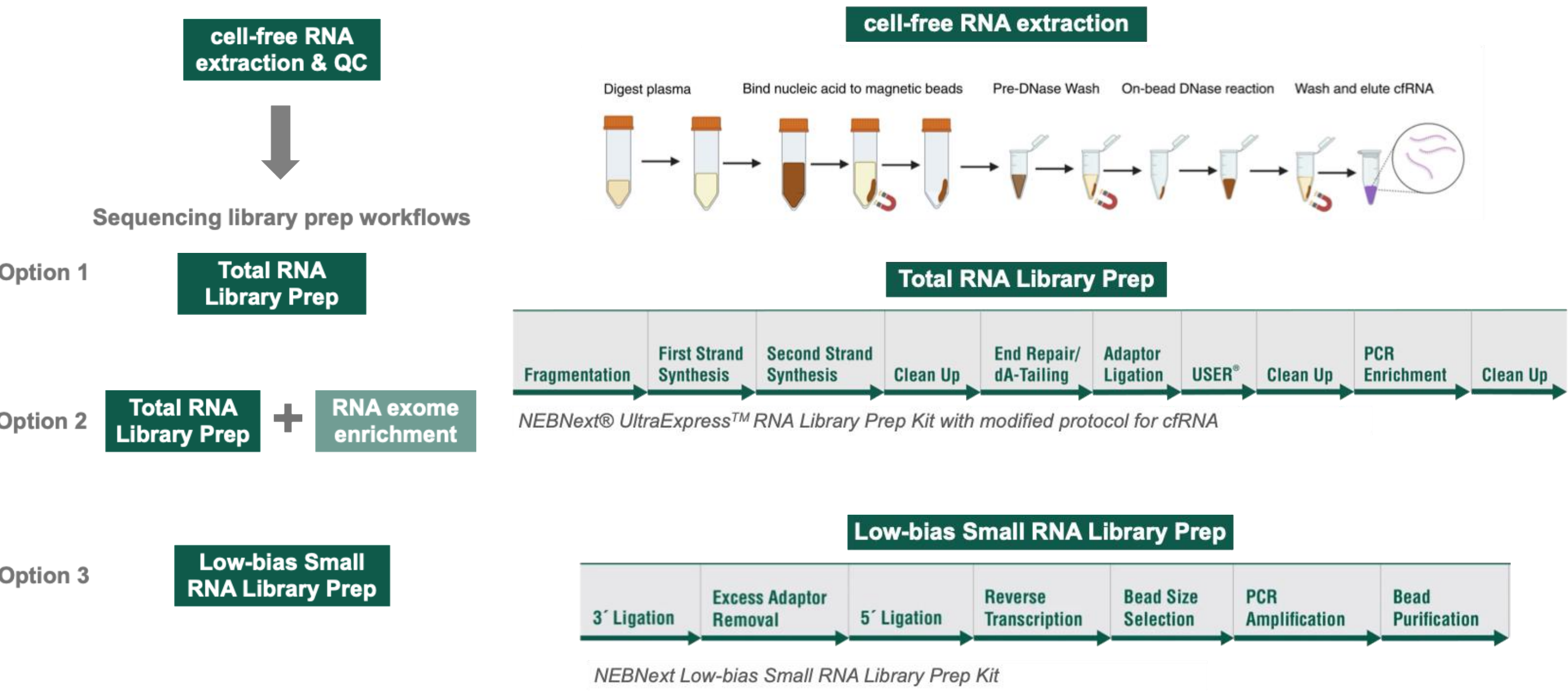
Chen Song, Chelsea Pinegar, Anagha Kadam, Heather Raimer Young, Gautam Naishadham, Keerthana Krishnan
New England Biolabs, Ipswich MA, USA

Background

Circulating cell-free RNA (cfRNA) has emerged as a promising biomarker for non-invasive diagnostics, offering dynamic insights into gene expression and disease states. Dynamic changes in cfRNA signatures can reflect both pathological and physiological processes, making it valuable for early disease detection, treatment monitoring, and prognostic assessments. However, clinical adoption of cfRNA-based diagnostics remains limited, largely due to technical and biological challenges throughout the workflow, including plasma processing, cfRNA extraction, quality control, and library preparation for sequencing. Here we show results from an optimized workflow for cfRNA detection, from extraction to sequencing.

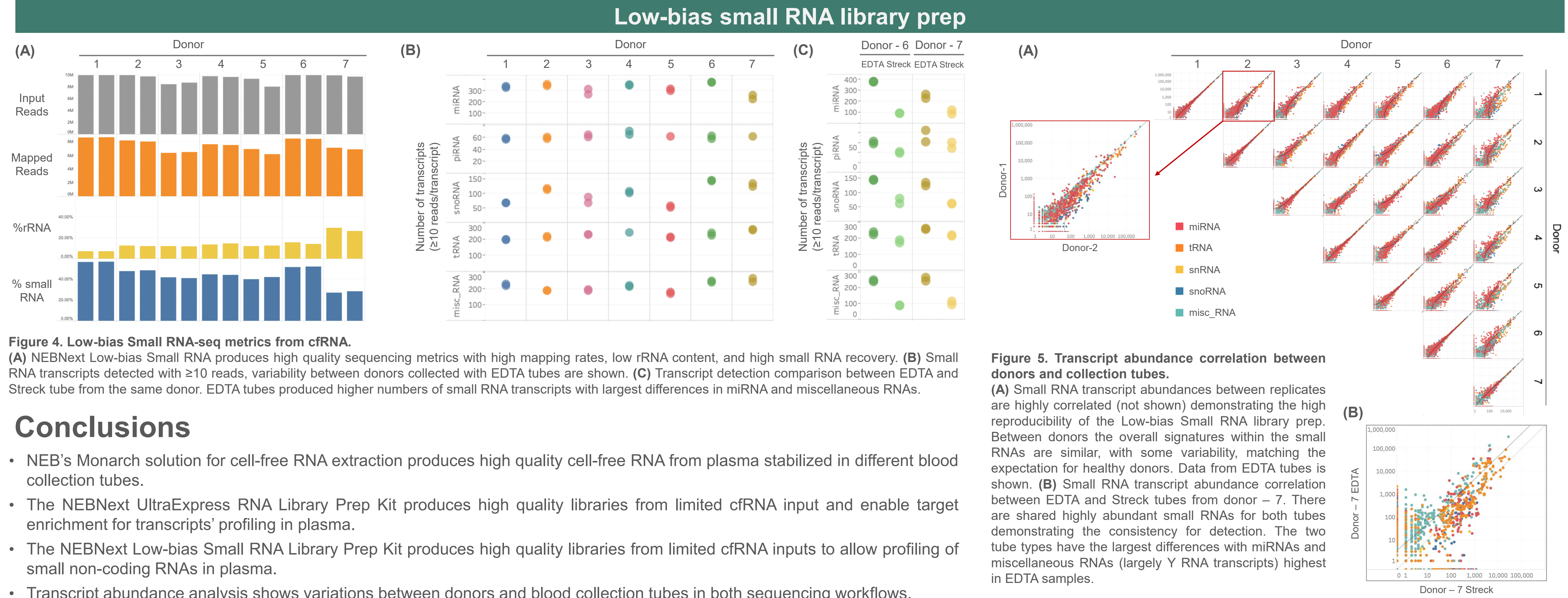
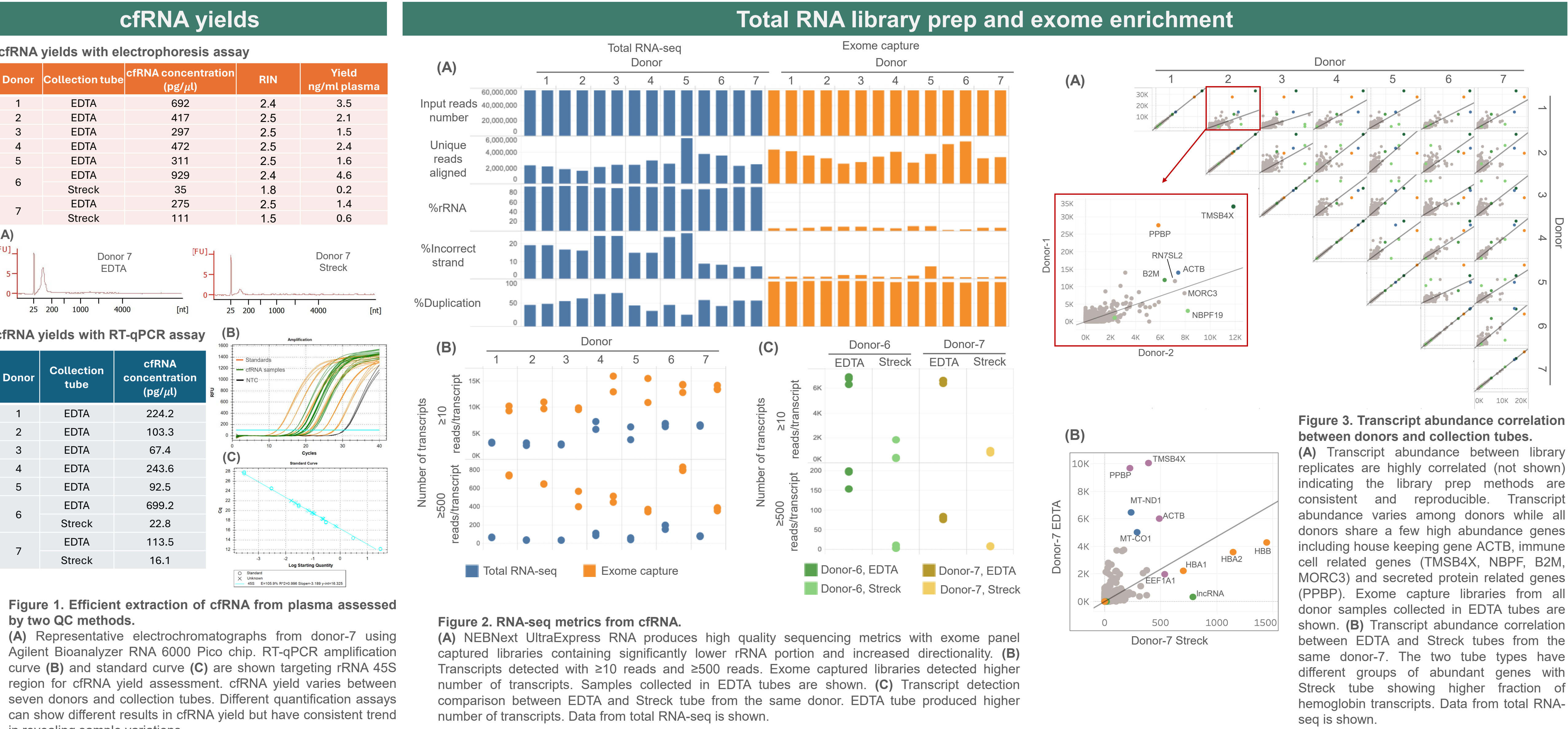
Methods

Following standard plasma collection and storage methods, we used a magnetic bead-based cfRNA extraction protocol developed to maximize RNA yield and minimize DNA contamination. cfRNA quality and quantity were measured by automated electrophoresis and qPCR assays. Sequencing library prep assays were performed on these samples, to generate basic quality metrics, including total RNA-seq, target enrichment with RNA exome panel and small RNA library prep. The assays optimized for mRNA or small RNA were used for assessing the respective expected performances.



Please reach out to NEB technical support if you are interested in applying these in-development protocols.

Results



Conclusions

- NEB's Monarch solution for cell-free RNA extraction produces high quality cell-free RNA from plasma stabilized in different blood collection tubes.
- The NEBNext UltraExpress RNA Library Prep Kit produces high quality libraries from limited cfRNA input and enable target enrichment for transcripts' profiling in plasma.
- The NEBNext Low-bias Small RNA Library Prep Kit produces high quality libraries from limited cfRNA inputs to allow profiling of small non-coding RNAs in plasma.
- Transcript abundance analysis shows variations between donors and blood collection tubes in both sequencing workflows.