

Efficient rRNA depletion using the new RiboMinus Bacteria 2.0 Kit to uncover the complexity of Bacterial transcriptomes

Charmaine San Jose Hinahon, Madhu Jasti, Varvara Dubovskaja, Angie Cheng and Xingwang Fang

Thermo Fisher Scientific, 2130 Woodward St, Austin, TX 78744

ABSTRACT

Global transcriptomic analysis is often hindered by the high abundance of ribosomal RNA (rRNA) content in bacterial cells, comprising up to 98% of RNA in cells. To remove unwanted rRNA and enrich for valuable protein coding transcripts of interest, rRNA depletion by capture probe hybridization has been the method of choice for RNA-seq, because it maintains relative transcriptome levels. The existing RiboMinus™ Bacteria design has had limited success with partially degraded total RNA, due to an abundance of fragmented rRNA contamination in compromised RNA that the current design of eight locked nucleic acid (LNA) probes does not address. Presently, we evaluate the efficiency of a new RiboMinus™ Bacteria 2.0 Kit in depleting rRNA from various inputs (100ng-5ug) of *E. coli* total RNA, which utilizes a new probe mix design. The RiboMinus™ procedure entails hybridization of a new set of 3'-biotin labeled ssDNA oligonucleotide probes (called RiboMinus™ Pan-Prokaryote Probe Mix) targeting sequence-specific regions of 16S, 23S rRNA and 5S rRNA. This RNA/probe hybrid is then removed from the sample using streptavidin-coated magnetic beads, leaving behind rRNA-depleted RNA. Sequencing results indicate that the newly developed RiboMinus™ Pan-Prokaryote Probe Mix in the RiboMinus™ Bacteria 2.0 workflow exhibits superior rRNA depletion efficiency compared to the existing kit, with less than 10% of total reads mapping to 16S and 23S rRNA vs. 90-95% of total reads mapping to rRNA using the existing kit. Removal of rRNA also results in a 10-fold increase of protein coding transcripts, from 5% in the existing kit to 55% in RiboMinus™ Bacteria 2.0. In addition, rRNA removal increases sensitivity for detection of noncoding transcripts. Finally, we also demonstrate that the newly developed RiboMinus™ Bacteria 2.0 Kit utilizing the new RiboMinus™ Pan-Prokaryote Probe Mix exhibits high ribosomal removal efficiency when *Salmonella enterica* RNA samples were tested.

MATERIALS AND METHODS

rRNA-depleted RNA was obtained from *E. coli* DH5α total RNA using the new RiboMinus™ Bacteria 2.0 Kit. Input amounts of 100ng-5μg total RNA were hybridized with the RiboMinus™ Pan-Prokaryote Probe Mix. The probe mix is a mixture of biotinylated DNA oligos designed against conserved regions of rRNA across a wide variety of pan-prokaryote species (see Table 1). Next, the rRNA-probe complexes are removed from the total RNA by capture with streptavidin-conjugated RiboMinus™ Magnetic Beads. The resulting rRNA depleted RNA is concentrated and purified with magnetic Nucleic Acid Binding Beads (see Figure 2). Approximately, 50ng of rRNA-depleted RNA was used in the whole transcriptome library workflow as described in the Ion Total RNA-Seq Kit v2 manual. Libraries were sequenced using the Ion GeneStudio™ S5 Systems.

RNA-Seq data analysis. FastQC (Babraham Bioinformatics, [http://www.bioinformatics.bbsrc.ac.uk/projects/fastqc/]) used to evaluate the quality of raw reads. All samples were aligned to reference genomes using universal RNA-seq aligner STAR and alignment metrics collected (alignment rate, chimeric re [https://www.ncbi.nlm.nih.gov/pubmed/23104886] ads fraction, duplication rates, etc). Biotype distribution was calculated based on QORTS read counts ([https://bmcbioinformatics.biomedcentral.com/articles/10.1186/s12859-015-0670-5]).

FIGURE 1. WHOLE TRANSCRIPTOME RNA-SEQ WORKFLOW



TABLE 1 – LIST OF PHyla & SUB-PHyla REPRESENTATIVE SPECIES IN PROBE DESIGN

ARCHAEA	Proteobacteria	Gram-positives	Cyanobacteria	Spirochaetes	Green Sulphur Bacteria	Bacteroidetes	Planctomycetes	Chlamydiae	Radioresistant Micrococci	Green Non-Sulphur Bacteria	Thermotogae	Aquificae	Acidobacteria	Fibrobacteres
Thaumarchaeota Crenarchaeota	α β γ δ ε	Actinobacteria Firmicutes Photosynthetic Bacteria	Spirochaetes Leptospiras	Chlorobi	Bacteroides Flavobacterium	Thermophilus Deinococcus	Chloroflexi	Aquifex						

Table 1. RiboMinus™ Pan-Prokaryote Probe Mix is designed against 76 representative species of bacteria and archaeobacteria from the common phyla and sub-phyla listed in Table 1. The Probe Mix consists of 3' biotinylated ssDNA oligos designed to target 5S, 16S, and 23S rRNA. The probe mix contains over 100 designs and has been tested against *E. coli*, *Bifidobacterium longum subsp. infantis*, *Staphylococcus aureus*, *Bacteroides thetaiotaomicron* and *Salmonella enterica*.

FIGURE 2 – RIBOMINUS™ BACTERIA 2.0 DEPLETION WORKFLOW

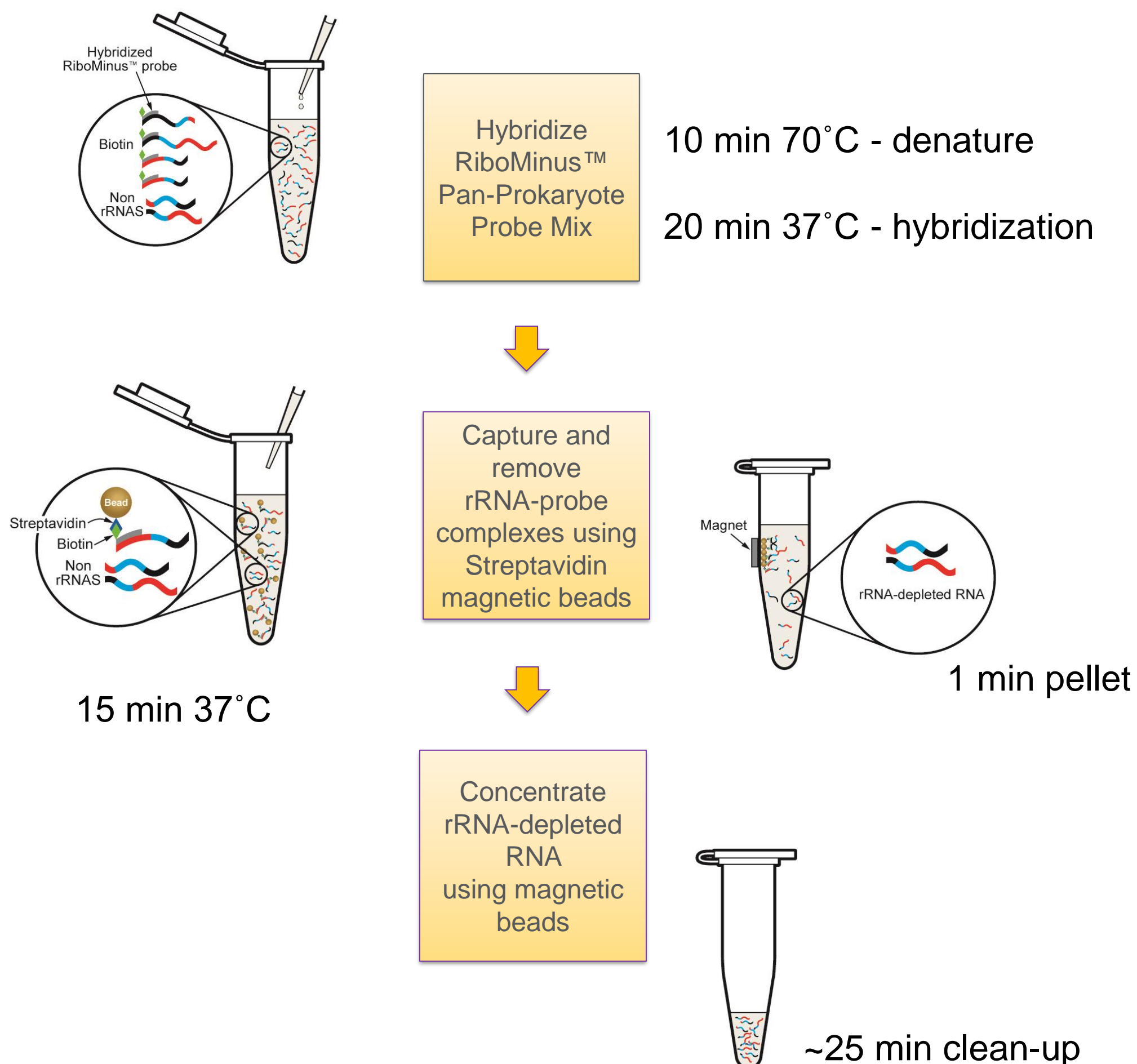


FIGURE 3 – ANALYSIS OF rRNA-DEPLETED RNA

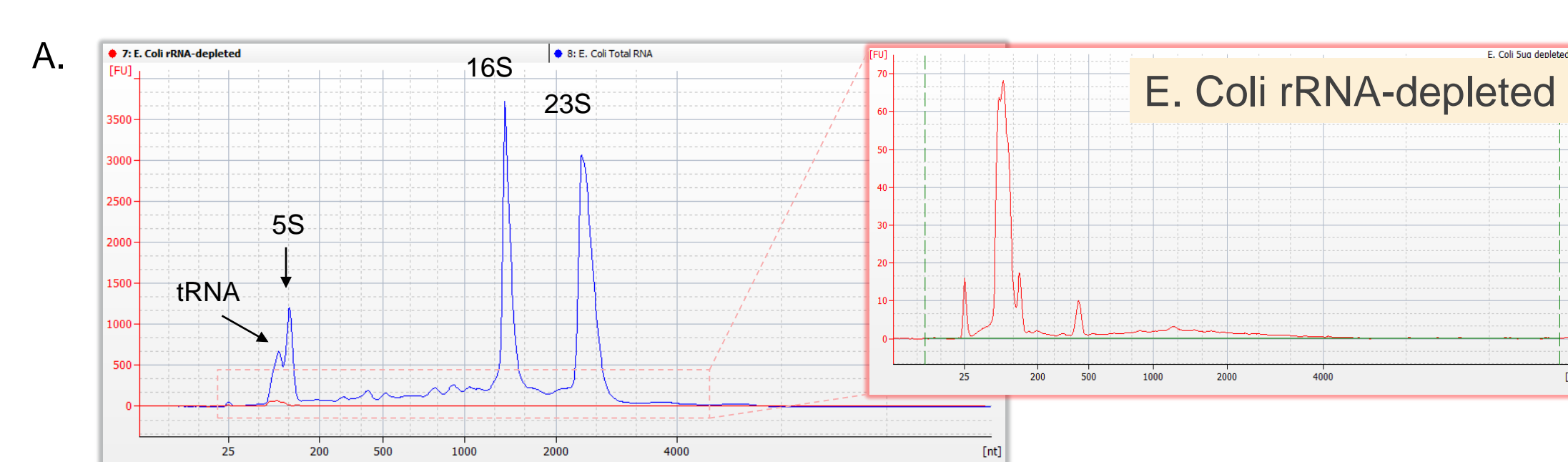
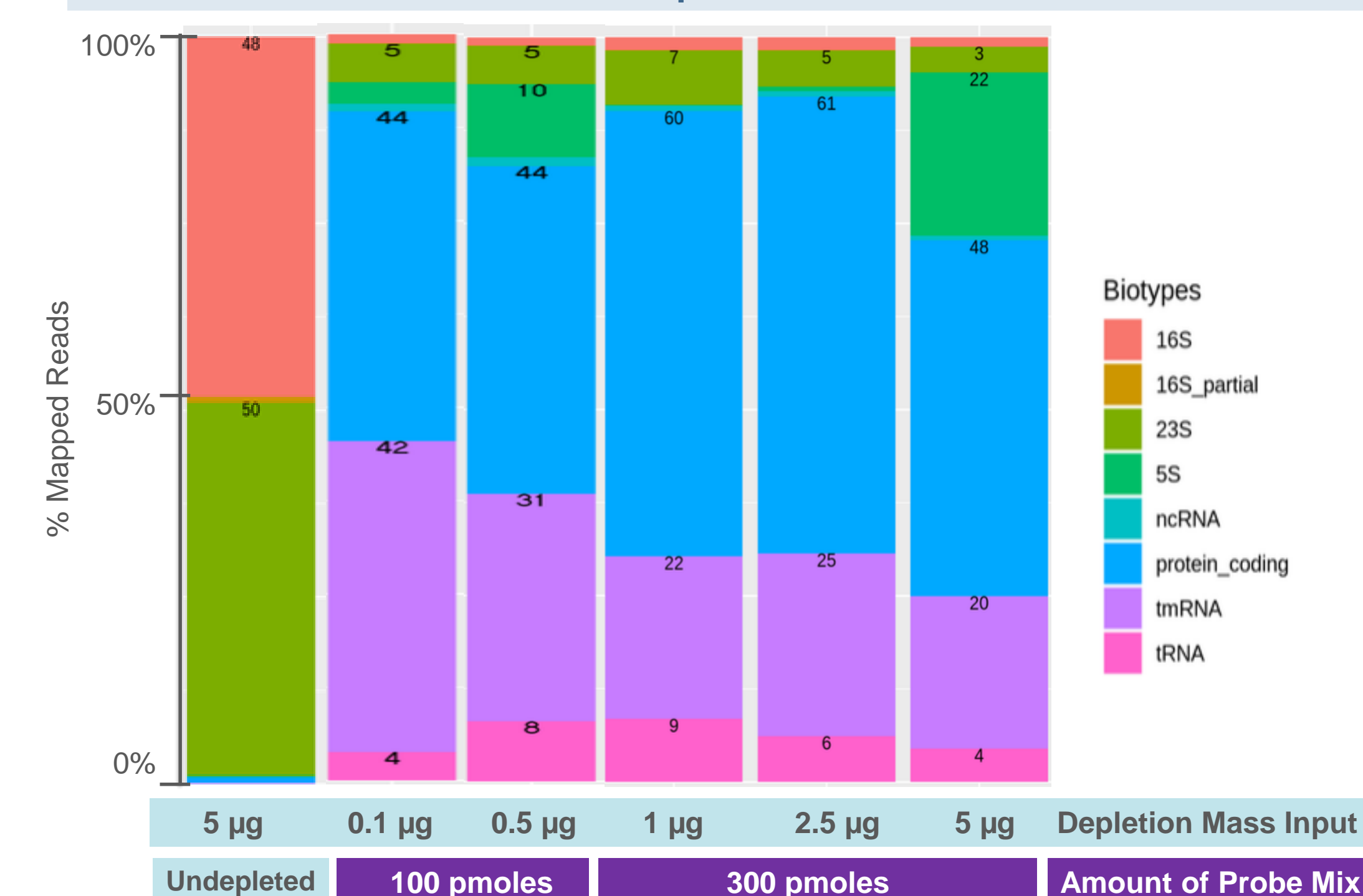


FIGURE 3A. Agilent® 2100 Bioanalyzer® profiles using the RNA 6000 Pico Assay showing an overlay of *E. coli* rRNA-depleted RNA (left) and an enlarged detailed trace of rRNA-depleted RNA (aka RiboMinus RNA) (right). rRNA-depleted RNA is shown in red and undepleted total RNA in blue. 5 μg of *E. coli* (RIN 8.7) and 300 pmoles of RiboMinus™ Pan-Prokaryote Probe Mix was used to perform the RiboMinus™ Bacteria 2.0 workflow.

B. RNA-Seq detected transcripts by biotypes for *E. coli* rRNA-depleted libraries



C. RNA-Seq detected transcripts by biotypes for *E. coli* Manual vs Automated KingFisher rRNA depletion workflow

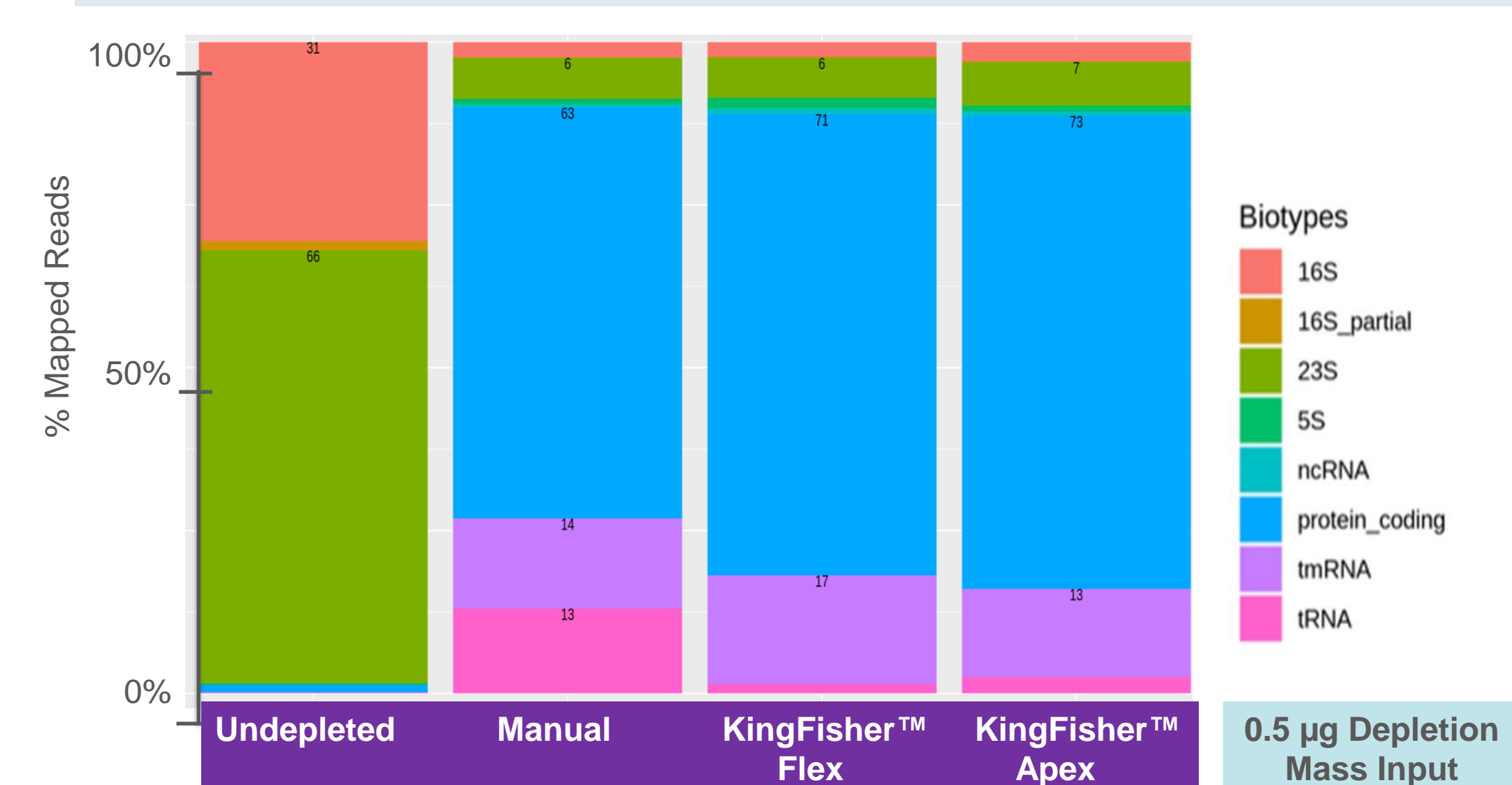


FIGURE 3B. Mapping Statistics Performance of RiboMinus Bacteria 2.0 Kit with *E. coli* using 100 ng to 5 μg mass input total RNA into depletion. Compared to undepleted, samples treated with the new Pan-Prokaryote Probe Mix show a much greater percentage of reads mapping to protein coding genes, from <2% in undepleted samples to >44% in rRNA-depleted samples. **3C.** Mapping Statistics comparing manual and automated depletion performance using KingFisher™ instruments on *E. coli* total RNA.

D. RNA-Seq detected transcripts by biotypes for *Salmonella enterica* rRNA-depleted libraries

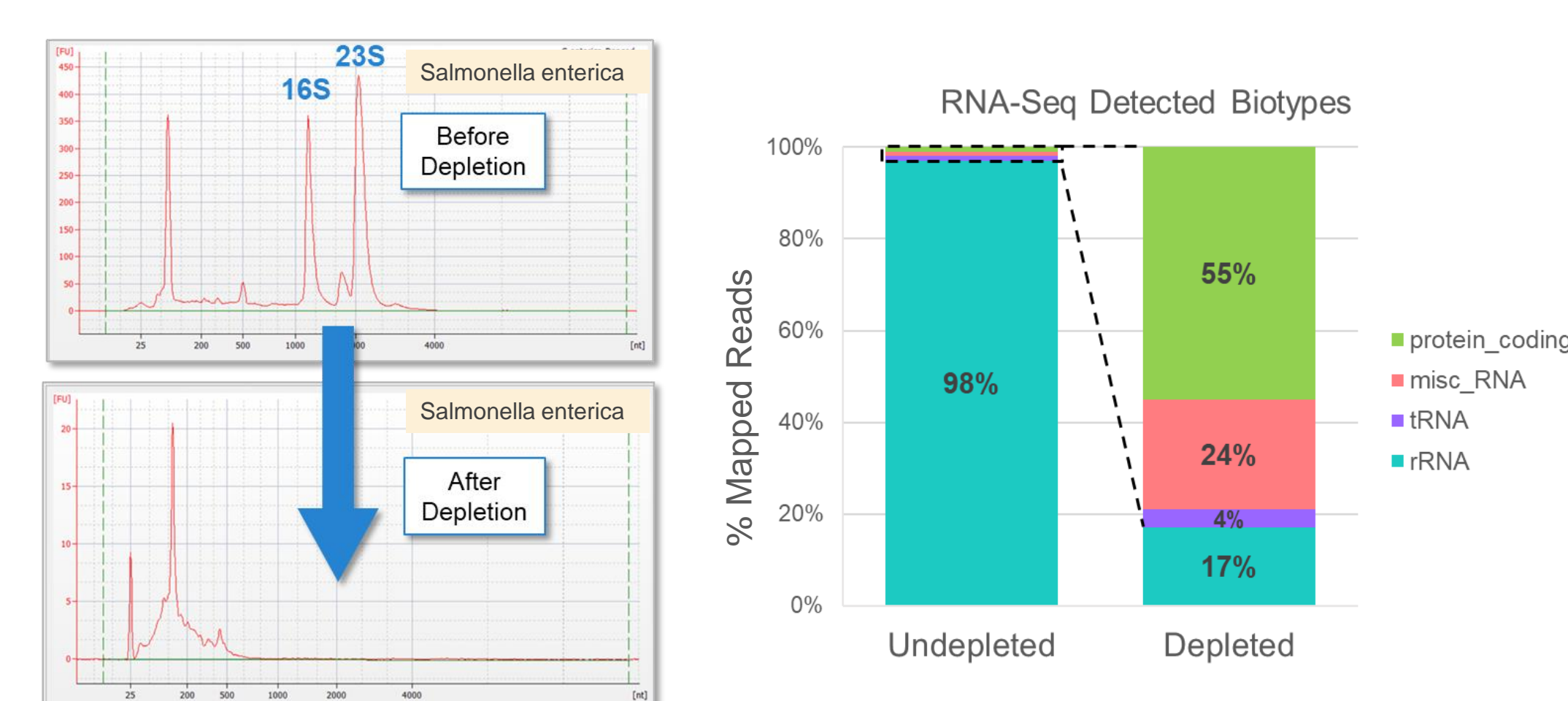


FIGURE 3D. Mapping Statistics Performance of RiboMinus Bacteria 2.0 Kit with *Salmonella enterica* using 1 μg mass input total RNA into depletion. Compared to undepleted, samples treated with the new Pan-Prokaryote Probe Mix show a much greater percentage of reads mapping to protein coding genes, from <2% in undepleted samples to >55% in rRNA-depleted samples.

FIGURE 4 – COMPARISON TO RIBOZERO

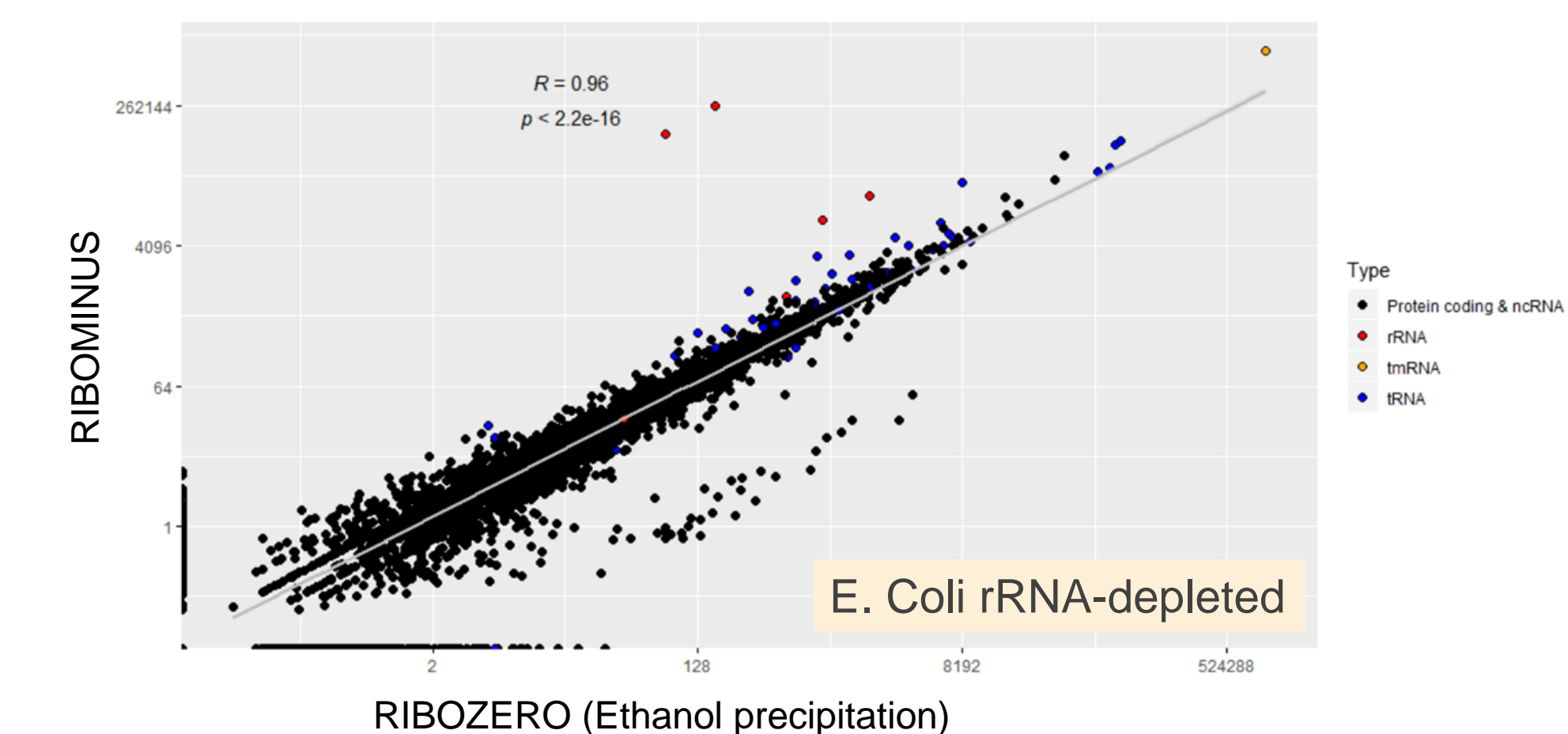


FIGURE 4A. Scatter plot showing pairwise count comparisons with high correlation between RiboZero and RiboMinus Bacteria 2.0. $R=0.96$.

BENEFITS OF NEW RIBOMINUS BACTERIA 2.0 KIT

- Efficient depletion of ribosomal RNA: 5S, 16S, and 23S targets
- Compatible with a wide range of pan-prokaryote species
- Broad input range: 100ng - 5μg total RNA
- < 2 hours manual workflow for 12 depletion samples
- Workflow is automatable on KingFisher instruments
- Compatible with meta-transcriptomics studies

CONCLUSIONS

- We developed a fast and efficient workflow for depletion of ribosomal RNA from Bacterial total RNA samples towards transcriptome analysis.
- Ribosomal RNA is consistently and efficiently removed.
- RNA-Seq data indicates that recovered rRNA-depleted RNA (aka RiboMinus™ RNA) is enriched for protein coding genes.
- This new kit is also applicable to microbiome preps

REFERENCES

- Petrova, O.E.; Garcia-Alcalde F.; Zampaloni, C. and Sauer, K.: Comparative evaluation of rRNA depletion procedures for the improved analysis of bacterial biofilm and mixed pathogen culture transcriptomes. Nature Scientific Reports (7). doi: 10.1038/srep41114.
- Hartley SW, Mullikin JC (2015) QoRTs: A comprehensive toolset for quality control and data processing of RNA-Seq experiments. BMC Bioinformatics 16:224.

ACKNOWLEDGEMENTS

We are grateful to the Austin Sample Prep teams and Vilnius Bioinformatics and Production teams for their assistance during the development of this workflow.

Also, we would like to thank our alpha and early access customers for their time and feedback.

TRADEMARKS/LICENSING

For Research Use Only. Not for use in diagnostic procedures. © 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.