

# rRNA removal and optimization of host-bacterial expression data

Jonathan Shaffer, M.B.A., Ph.D.  
Associate Director of Research & Development

## Legal disclaimer

QIAGEN products shown here are intended for molecular biology applications. These products are not intended for the diagnosis, prevention or treatment of a disease.

For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN kit handbook or user manual. QIAGEN kit handbooks and user manuals are available at [www.qiagen.com](http://www.qiagen.com) or can be requested from QIAGEN Technical Services or your local distributor.

# Agenda

Background

---

FastSelect overview

---

FastSelect results: Mammalian, bacterial, mammalian + bacterial

---

Summary

---



# Agenda

Background

---

FastSelect overview

---

FastSelect results: Mammalian, bacterial, mammalian + bacterial

---

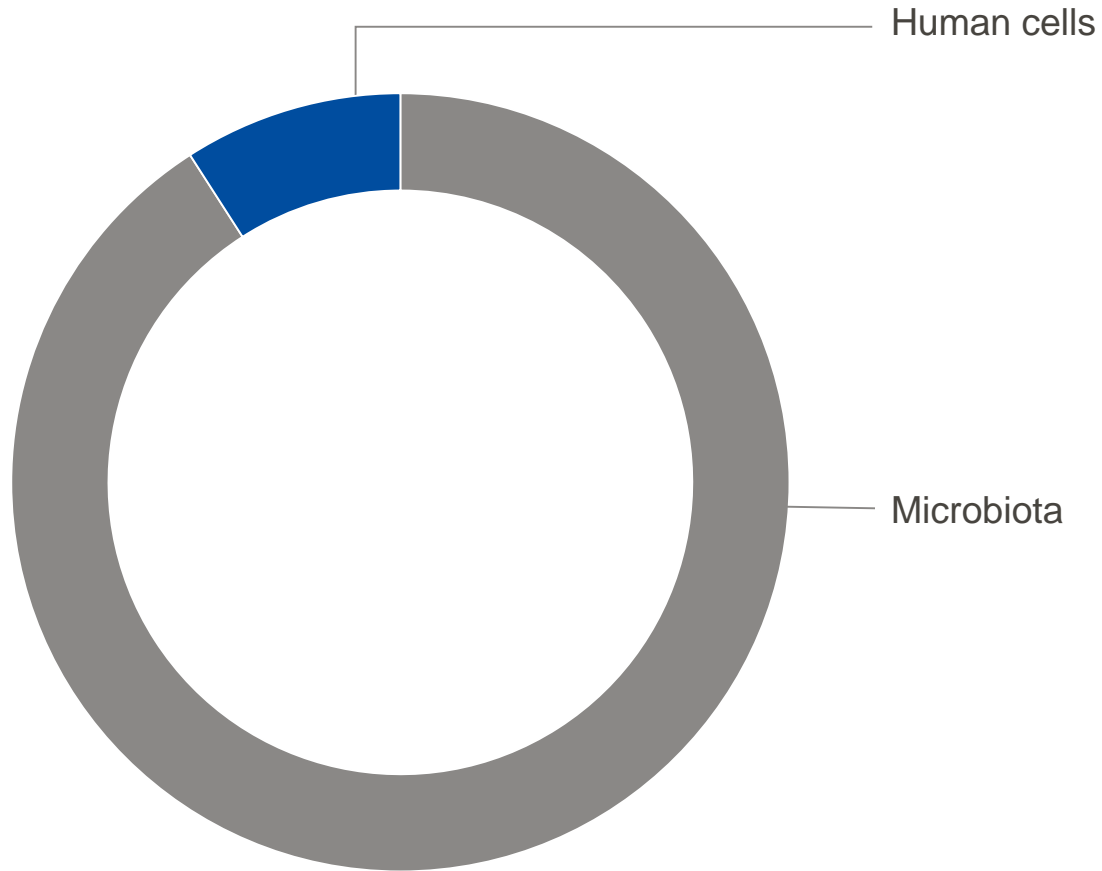
Summary

---



# Humans or 'superorganisms'?

## Total number of cells: Human cells versus microbiota



## Cellular composition of the 'superorganism'

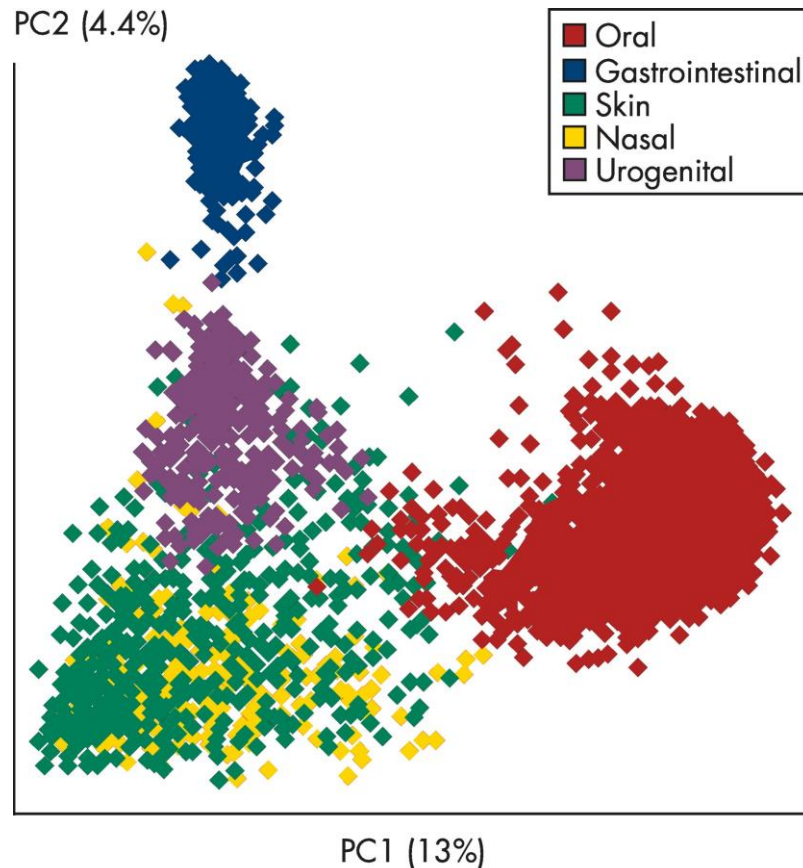
Estimation of the number of microbial cells that live in and on the human body; human cells are outnumbered by a factor of 10

Nomenclature:

- **Microbiota** are the microbes that live in a specific location, e.g., the human body, the gut, soil, etc.
- **Metagenomics** is the study of the collection of genomes derived from a specific sample or community
- **Metatranscriptomics** is the study of the RNA expression of genes from a community sample to interpret the physiological state of that community at that time

# Microbiota composition

## Microorganisms cluster by body site



Cataloguing efforts by the NIH Human Microbiome Project suggest:

- Around 10,000 organisms live with us
- Around  $8 \times 10^6$  genes constitute this 'second genome'

Identifying microbiota in healthy individuals revealed:

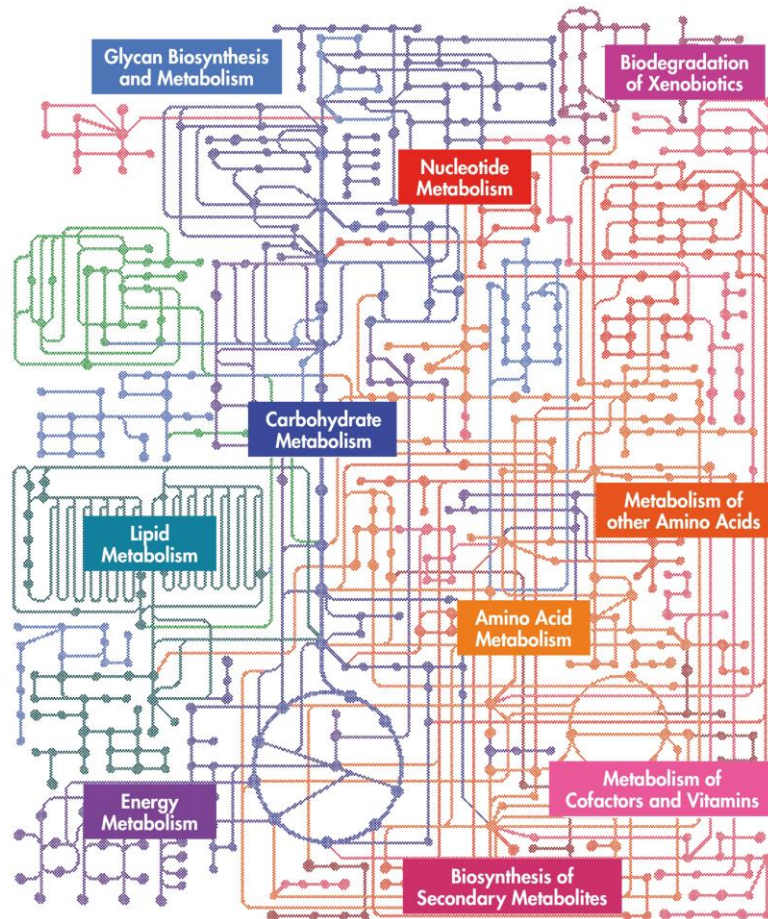
- Different body sites have unique communities
- Race, age, gender, weight or ethnicity can affect microbiota composition

Human Microbiome Project Consortium (2012). Structure, function and diversity of the healthy human microbiome. *Nature*, 486(7402), 207–214.



# Complexity and function of genomic content

## Function of microbiome enables individual survival

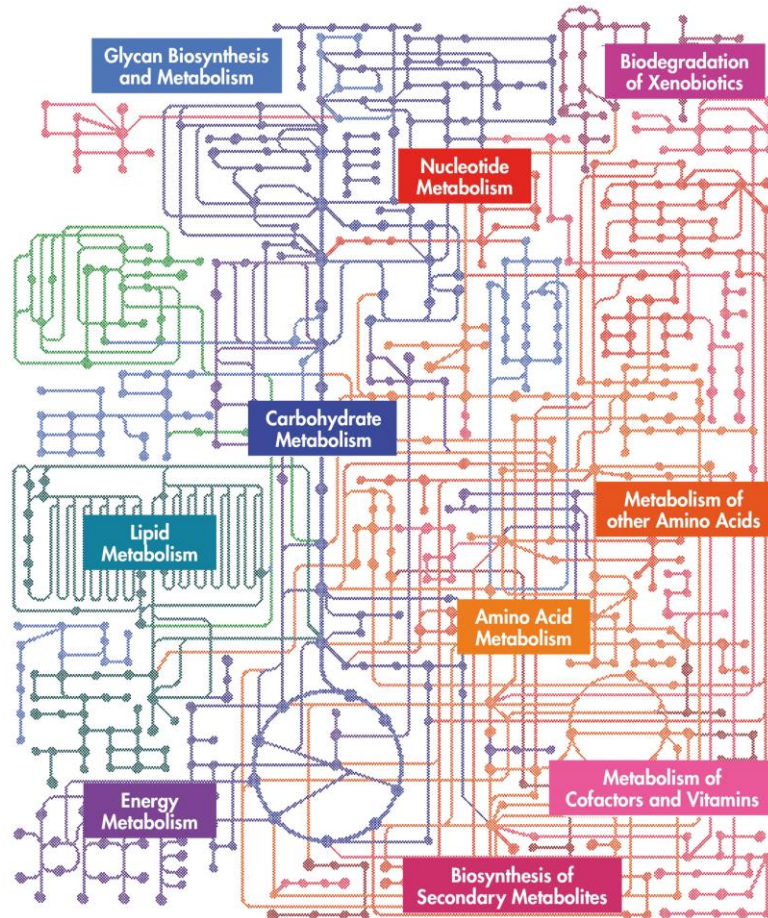


Each organism has developed its:

- Genetic content for its own survival in a specific environment
- Metabolism tuned to local nutrient sources
- Virulence factors for stable colonization
- Antibiotic resistance genes to metabolize toxins

# Complexity and function of genomic content (continued)

## Function of microbiome enables individual survival



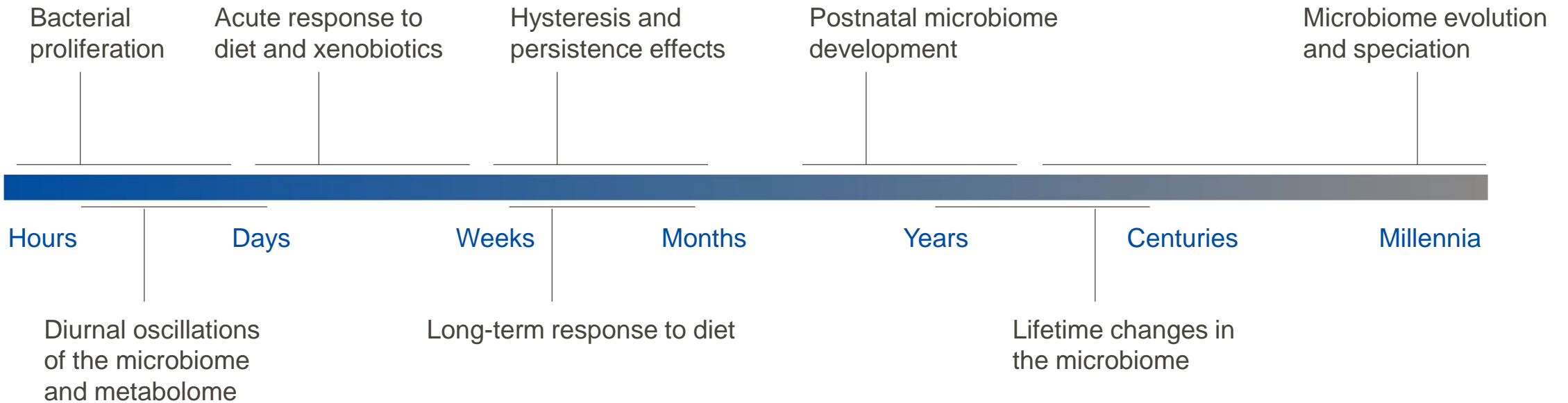
## Understanding the microbiota

- DNA – who is or was there
- RNA – who is alive and what they are doing
- RNA can be used to identify:
  - Metabolic status
  - Expression of virulence factors
  - Expression of antibiotic resistance
  - Both host and microbiota

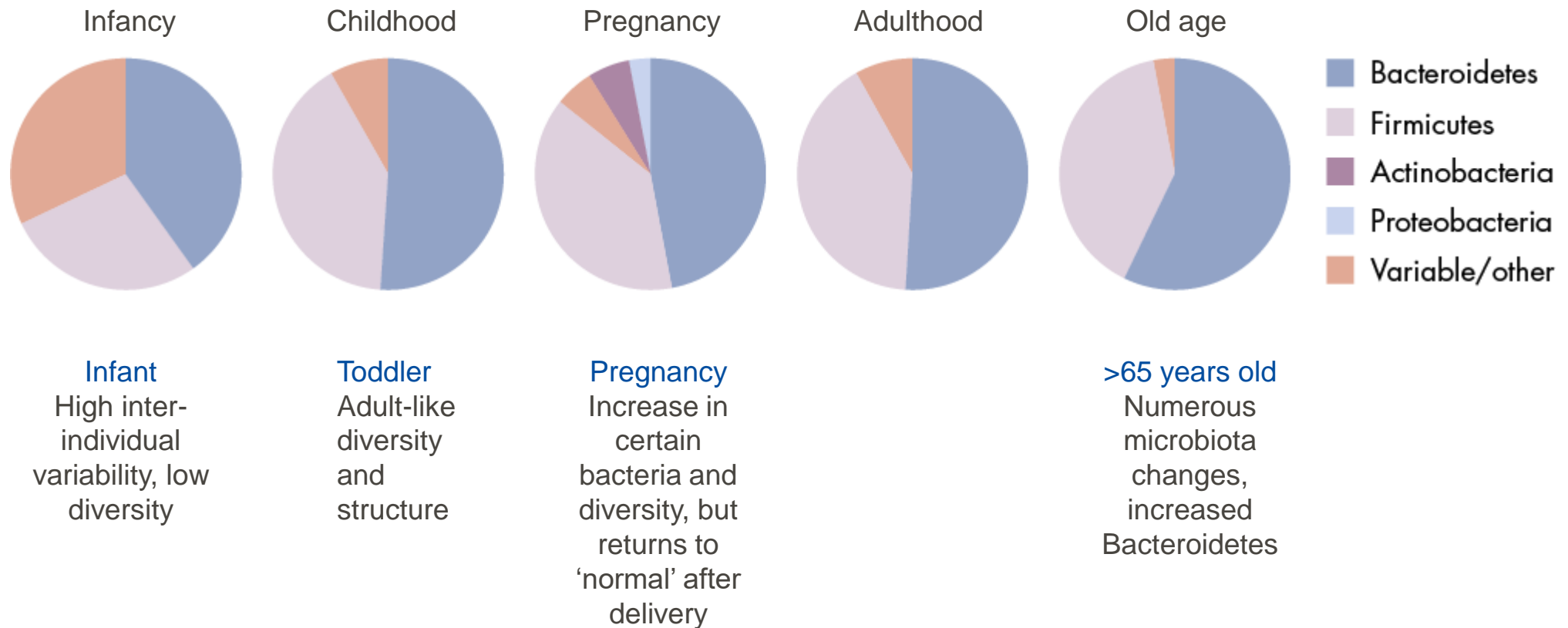


# Microbiota dynamics important for host physiology

## Function of microbiome enables individual survival



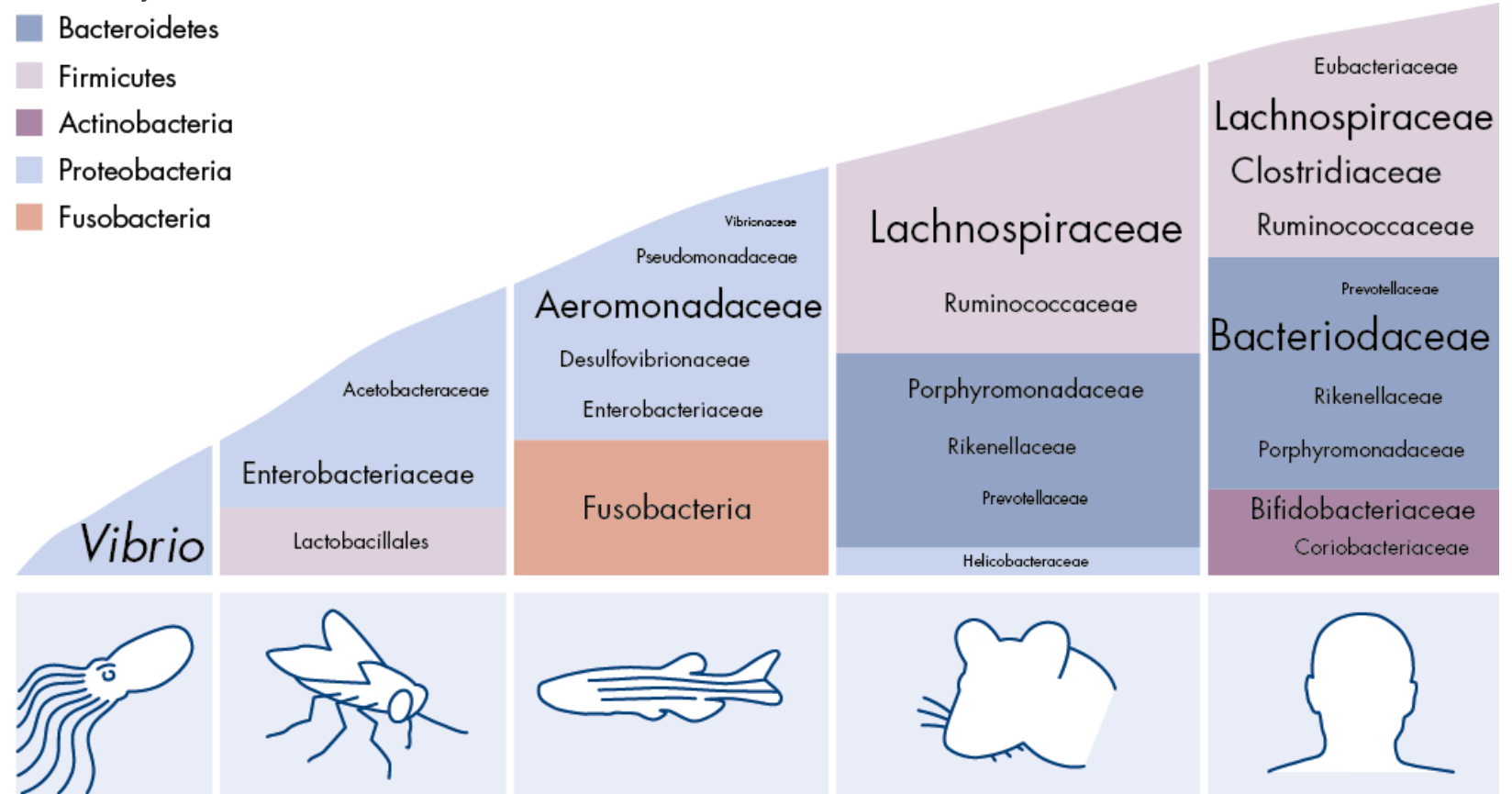
# Intestinal microbiota changes throughout the course of life



Source: Kostic, A.D., Howitt, M.R. and Garrett, W.S. (2013) Exploring host-microbiota interactions in animal models and humans. *Genes Dev.*27:701-718

# Broad microbiota trends exist across species, especially at the phylum level

- **Phyla:** Represented by color
- **Abundance:** Lower taxonomic levels is indicated by font size.

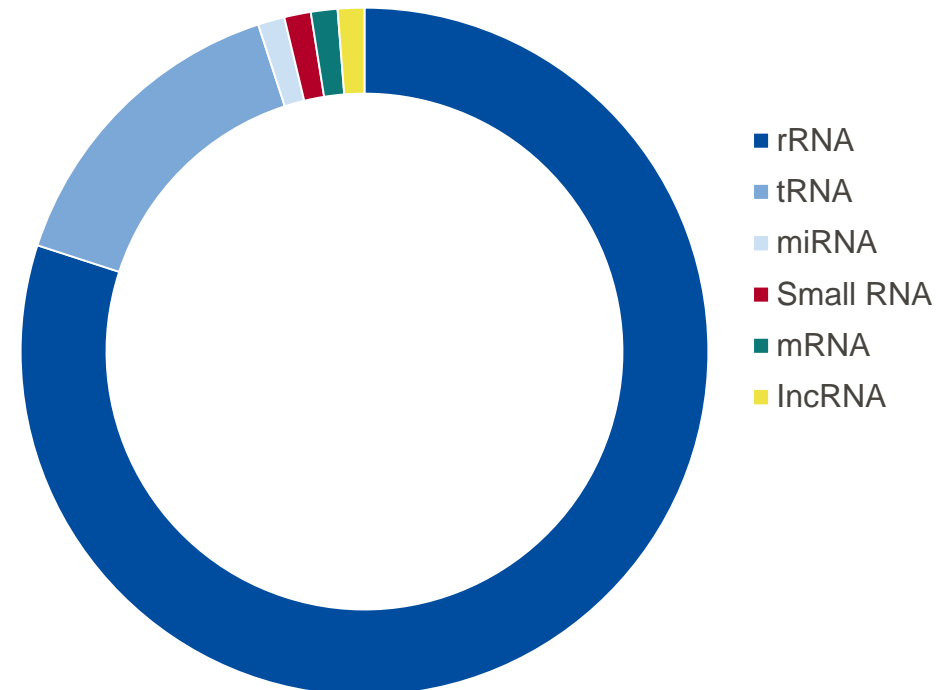


Source: Kostic, A.D., Howitt, M.R. and Garrett, W.S. (2013) Exploring host-microbiota interactions in animal models and humans. Genes Dev.27:701-718

# Library enrichment/depletion strategies: Whole transcriptome NGS

Enrichment or depletion is necessary to maximize reads from the RNAs of interest

- Typical RNA composition in a cell: >80% ribosomal RNA
- Highly abundant transcripts consume a lot of reads
- Enrichment or depletion is used to obtain more reads from the RNAs of interest, such as:
  - mRNA
  - lncRNA

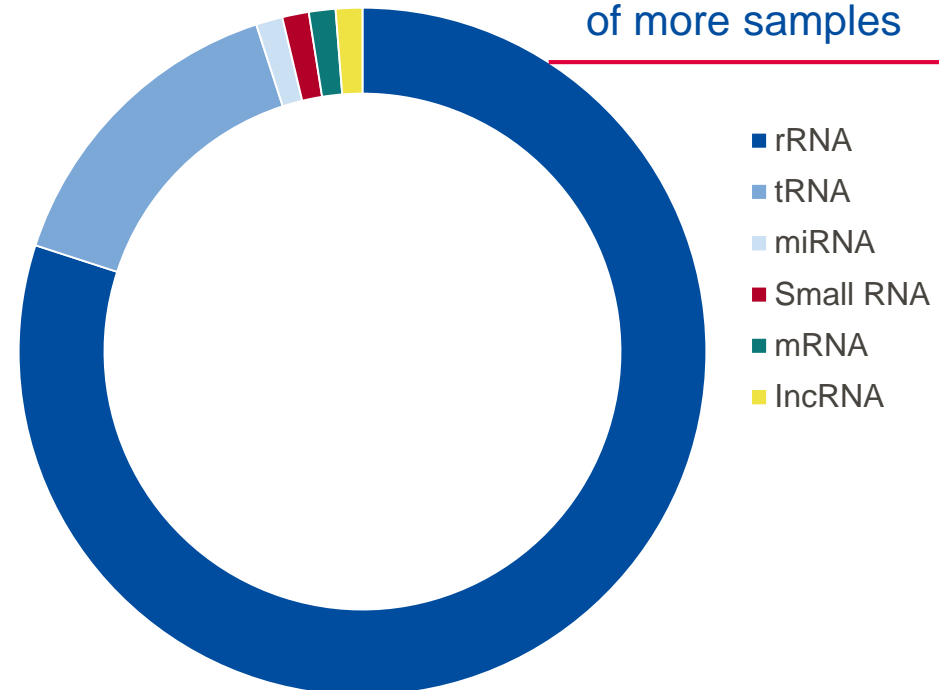


# Library enrichment/depletion strategies: Whole transcriptome NGS

Enrichment or depletion is necessary to maximize reads from the RNAs of interest

- Typical RNA composition in a cell: >80% ribosomal RNA
- Highly abundant transcripts consume a lot of reads
- Enrichment or depletion is used to obtain more reads from the RNAs of interest, such as:
  - mRNA
  - lncRNA

Removing rRNA will increase sensitivity and decrease the cost of sequencing, allowing analysis of more samples





# Library enrichment/depletion strategies: Whole transcriptome NGS

## Enrichment/depletion strategies

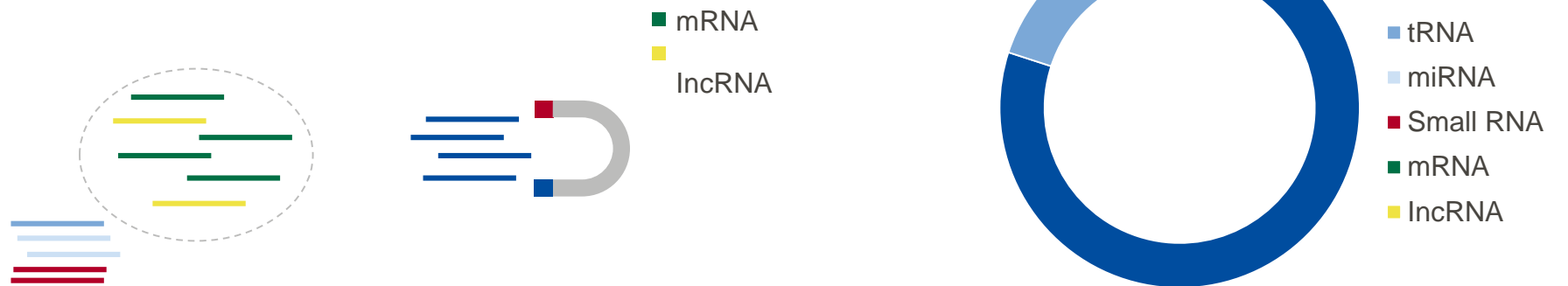
### Poly(A) enrichment

- Not useful for fragmented samples
- Not useful for prokaryotic samples



### rRNA depletion

- The 'catch all' method



- rRNA depletion is the only choice for metatranscriptomics.

# Agenda

Background

---

FastSelect overview

---

---

FastSelect results: Mammalian, bacterial, mammalian + bacterial

---

Summary

---



## Remove bacterial rRNA only: FastSelect –5S/16S/23S overview

### What is FastSelect –5S/16S/23S?

- Fragmentation and pan-bacterial (5S/16S/23S) rRNA depletion module

### Number of reactions:

- 24, 96 and 384

### How does it work?

- Inhibits reverse transcription of its specific targets

### Coverage:

- Designed to block **community level** cDNA synthesis of 5S, 16S and 23S rRNA
- Designed against SILVA 16S sequences (nearly 600,000 unique entries), SILVA 23S sequences (nearly unique 170,000 entries) and 5S rRNA Database (over 7,200 unique entries)
- Theoretically blocks >95% cDNA synthesis of all 5S, 16S and 23S rRNA sequences
  - In practice, results will vary, based on the exact composition of the sample

### Total RNA input:

- 20 ng to 1 µg

### Tested RNA library prep kit compatibility:

- QIAseq Stranded Total RNA Lib Kit (QIAGEN Group), TruSeq® Stranded (Illumina®, Inc), NEBNext® Ultra II Directional (New England Biolabs, Inc)
- FastSelect is compatible with most RNA library prep kits

# Remove mammalian rRNA and globin mRNA only: FastSelect –rRNA HMR and –Globin

## How does FastSelect –rRNA HMR and –Globin work?

- Inhibits reverse transcription of specific targets
- Removes cytoplasmic and mitochondrial rRNA and/or globin mRNA

## Species covered in a single-tube:

- Human, mouse, rat (HMR) and other mammalian species
- HMR removes 95–99% rRNA from cow, horse, sheep and hamster samples
- HMR removes 80–90% rRNA from dog, chicken, rabbit, pig and monkey samples

## RNA compatibility:

- Total RNA: Use FastSelect –rRNA HMR (include –Globin if working with whole blood)
- Poly(A) enriched RNA: Use FastSelect –Globin if working with whole blood

## Sample compatibility:

- Cell lines, tissues (fresh/frozen), FFPE tissues, blood and biofluids

## Total RNA input:

- 1 ng – 1 µg

## Tested RNA library prep kit compatibility:

- QIAseq Stranded Total RNA Lib Kit (QIAGEN), TruSeq Stranded (Illumina), NEBNext Ultra II Directional RNA Library Prep Kit (NEB), KAPA RNA HyperPrep Kit (Roche Group)
- FastSelect is compatible with most RNA library prep kits

## Remove both bacterial and host rRNA: FastSelect –5S/16S/23S plus –rRNA HMR

### Catalog numbers

- QIAseq FastSelect –5S/16S/23S: 335925, 335927 or 335929
- QIAseq FastSelect –rRNA HMR: 334386, 334378 or 334388
- QIAseq FastSelect –Globin: 334376, 334377 or 334378
- QIAseq FastSelect –rRNA/Globin: 335376, 335377 or 335378

### Where do I find the protocol?

- QIAseq FastSelect –5S/16S/23S Handbook

### How long does the protocol take?

- Only 1 hour, which includes fragmentation time!

### How does it work?

- Just like the individual products through inhibition of rRNA reverse transcription

### Total RNA input:

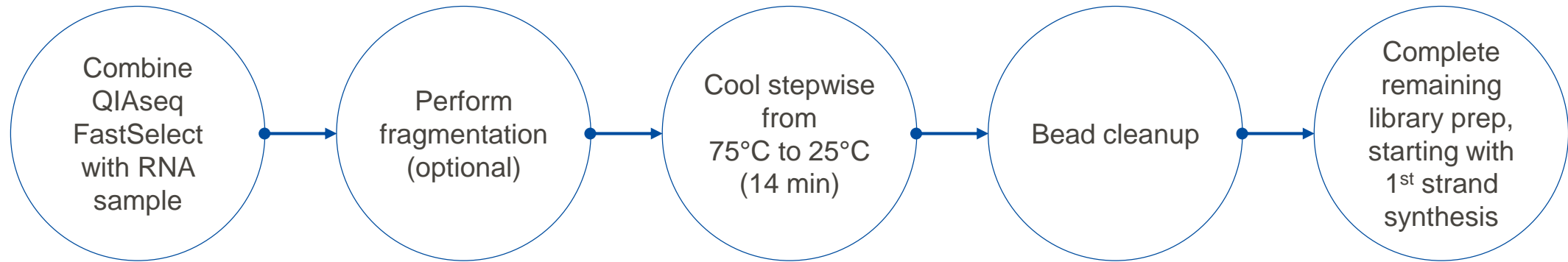
- 20 ng to 1 µg

### Tested RNA library prep kit compatibility:

- QIAseq Stranded Total RNA Lib Kit (QIAGEN Group), TruSeq Stranded (Illumina, Inc), NEBNext Ultra II Directional (New England Biolabs, Inc)
- FastSelect is compatible with most RNA library prep kits



## Workflow for removal of both bacterial and host rRNA



Mix and match 'FastSelects' (you can add all three or even more)

- **FastSelect –5S/16S/23S:** Bacterial rRNA
- **QIAseq FastSelect –rRNA HMR:** Mammalian rRNA
- **QIAseq FastSelect –Globin:** Mammalian Globin

# Agenda

Background

---

FastSelect overview

---

FastSelect results: Mammalian, bacterial, mammalian + bacterial

---

Summary

---

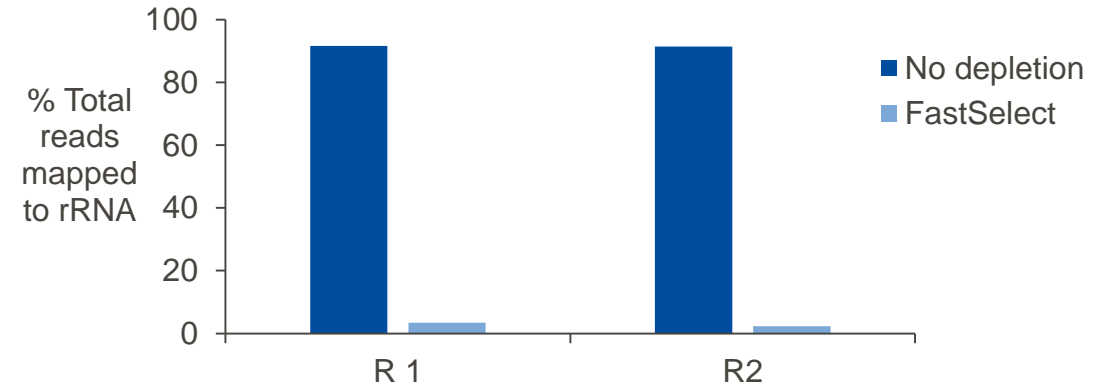


# FastSelect –rRNA HMR robustly removes rRNA, freeing up reads for gene detection

## Experimental overview

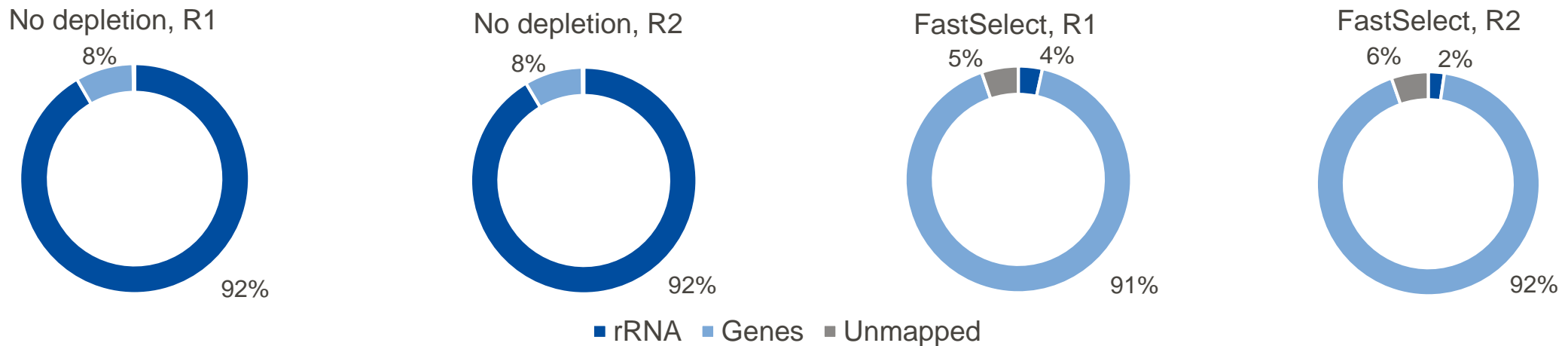
- Sample: 100 ng, Universal Human Reference RNA (UHRR)
- Depletion: No depletion, FastSelect –rRNA HMR
- Library prep: QIAseq Stranded Total RNA Lib Kit
- Sequencing: NextSeq 550 (2 x 75 bp)
- Mapping: CLC Genomics Workbench

## FastSelect substantially removes rRNA



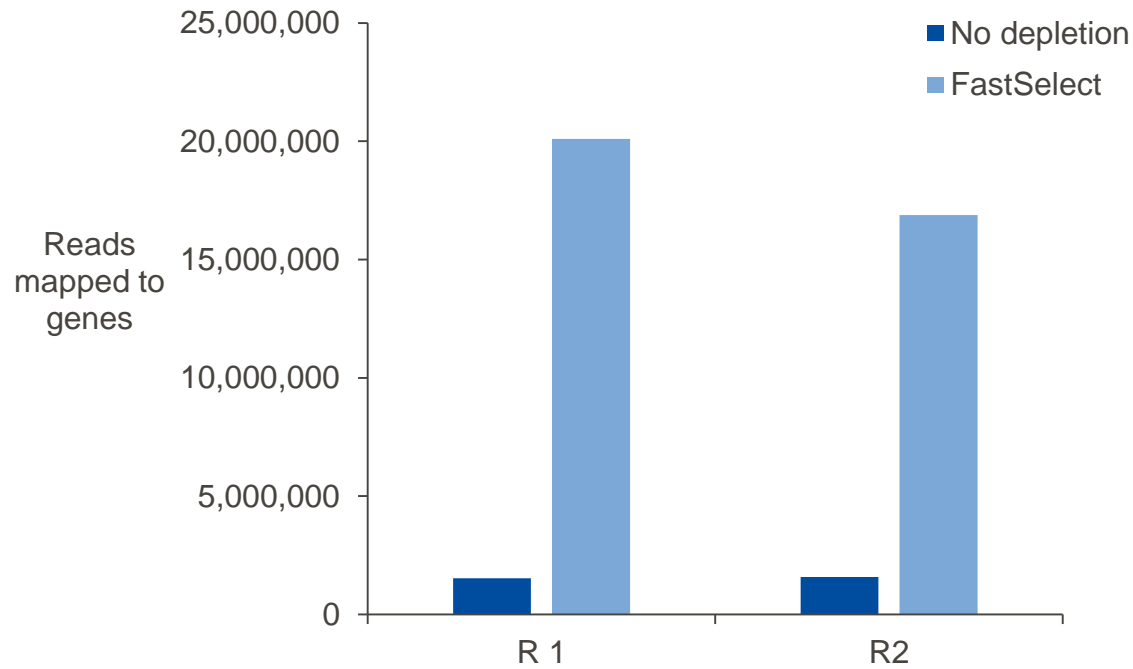
## FastSelect improves read mapping metrics, removing contaminating rRNAs

Breakdown of mapped reads

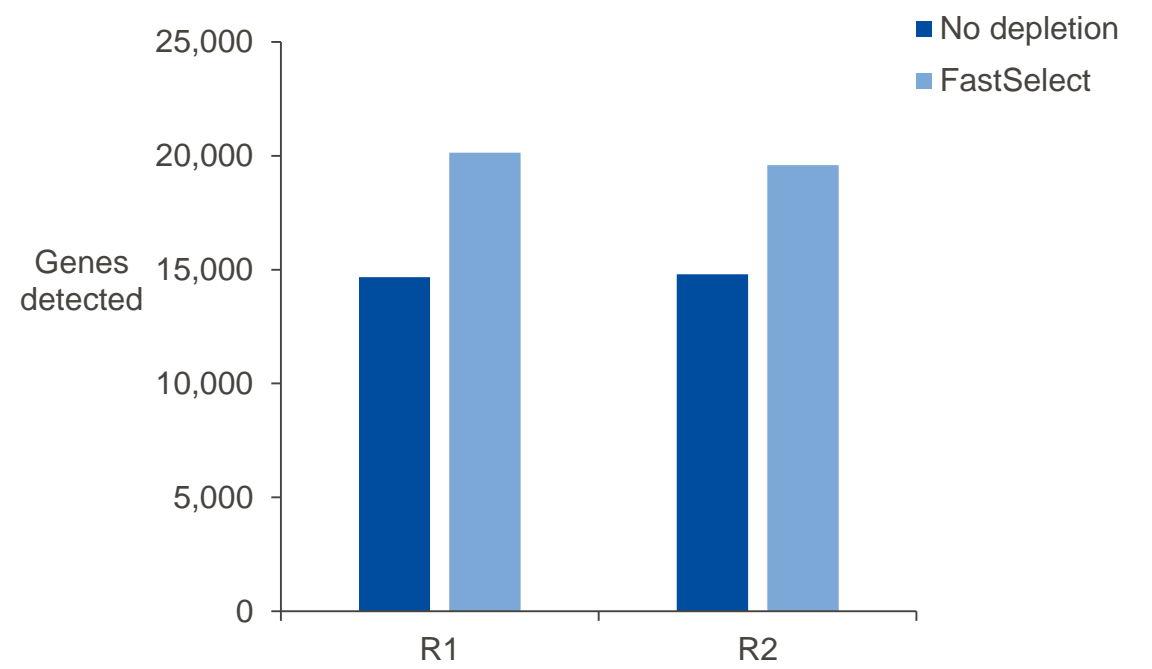


# FastSelect –rRNA HMR robustly removes rRNA, freeing up reads for gene detection

FastSelect dramatically increases reads mapped to genes



FastSelect increases the number of genes detected



- FastSelect removes nearly all human rRNA, freeing up reads for robust gene detection.

# FastSelect –5S/16S/23S: Robust rRNA removal from bacterial communities

## Experimental overview

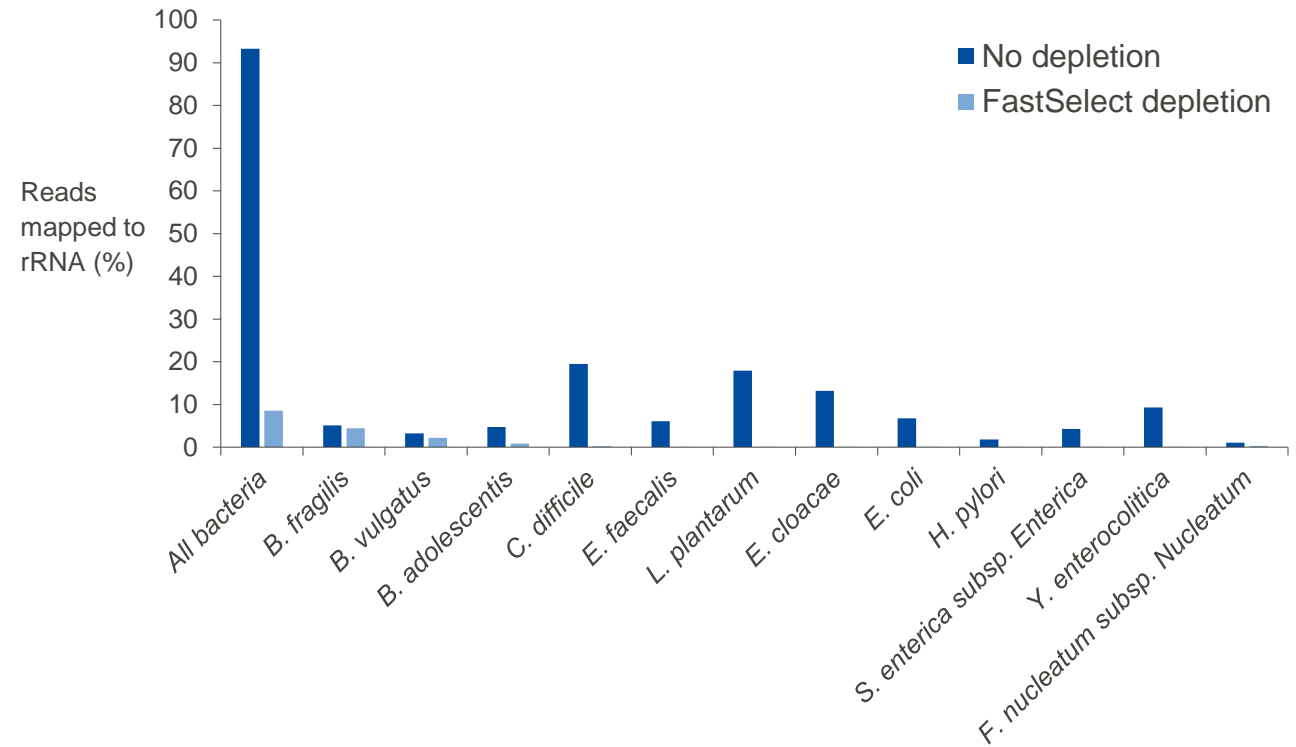
- Sample: 1 µg, Gut Microbiome Whole Cell Mix (ATCC)
- Depletion: No depletion, FastSelect –5S/16S/23S
- Library prep: QIAseq Stranded Total RNA Lib Kit
- Sequencing: NextSeq 550 (2 x 75 bp)
- Mapping: CLC Genomics Workbench

## FastSelect substantially removes rRNA

Sample	Percentage of reads mapped to bacterial rRNA (total)	
	No treatment	FastSelect –5S/16S/23S
ATCC Gut Microbiome (12 bacteria)	93.28	8.54

- FastSelect removes nearly 90% of all rRNA
- When individual species are mapped, FastSelect removes rRNA from a broad range of species

## Robust depletion of rRNA from individual species

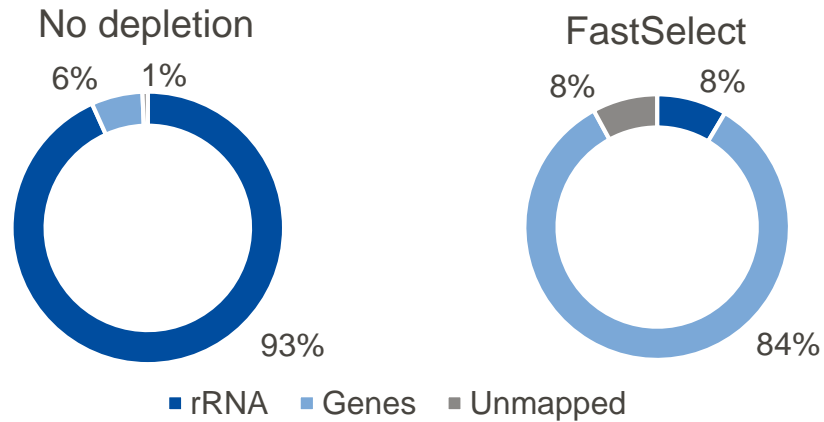




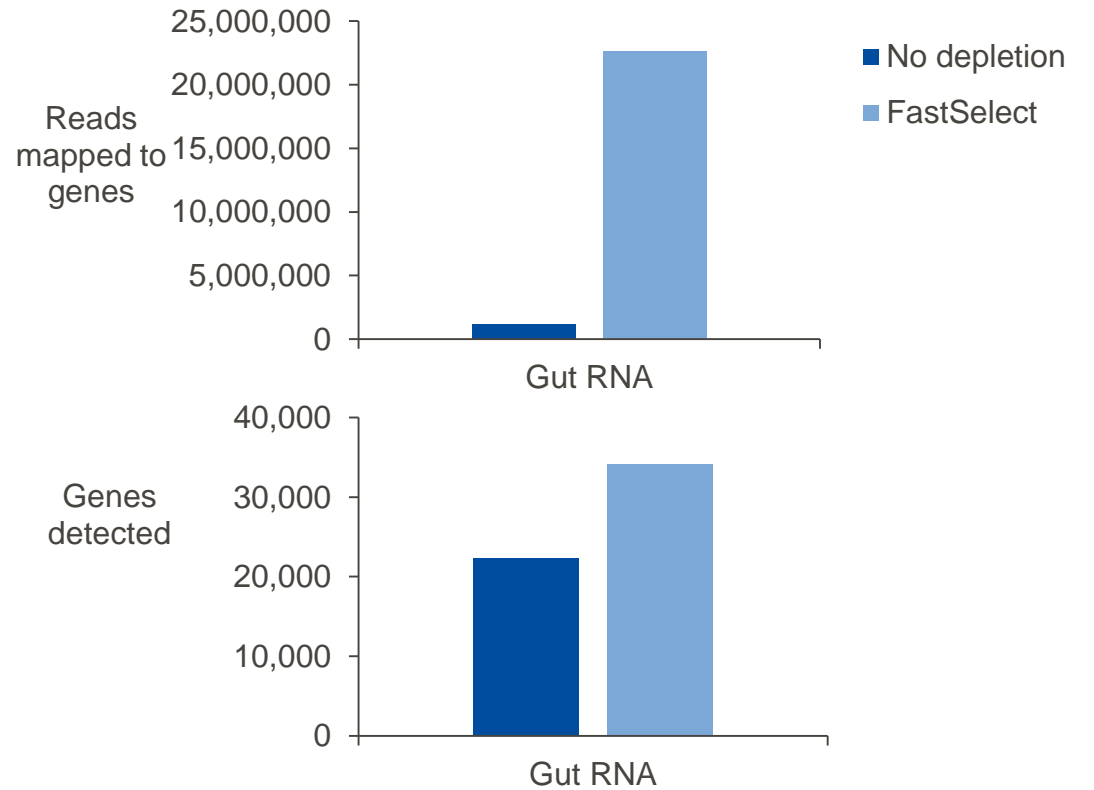
# FastSelect –5S/16S/23S: Robust rRNA removal from bacterial communities

## FastSelect dramatically improves read mapping metrics

Breakdown of mapped reads



## FastSelect dramatically increases reads mapped to genes, resulting in a substantial increase in genes detected



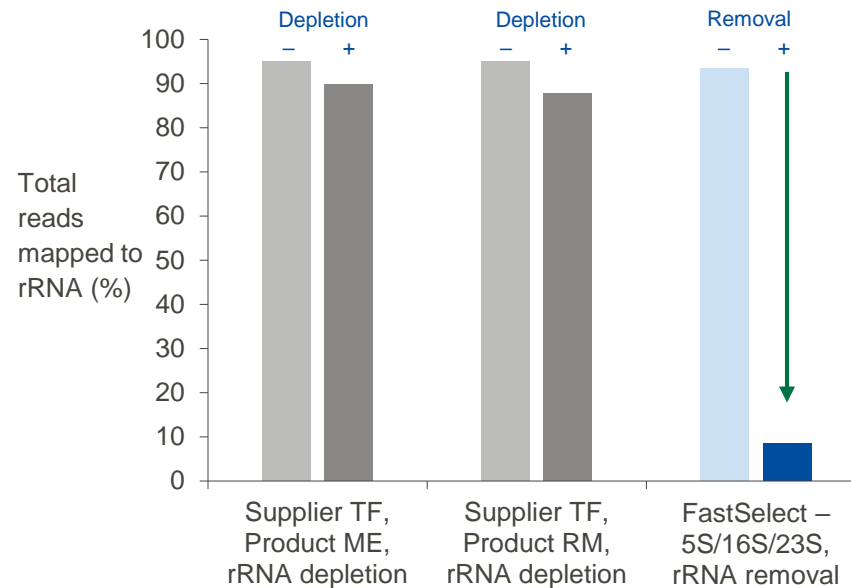
- FastSelect removes substantial amounts of bacterial rRNA, even from communities, freeing up reads for gene detection.

# FastSelect –5S/16S/23S dramatically outperforms the other supplier's kits

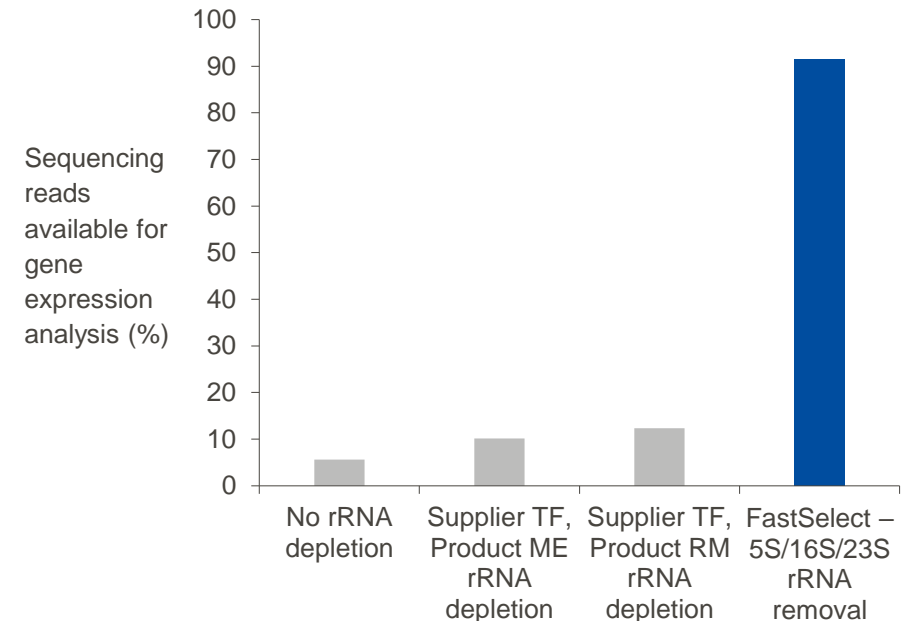
## Experimental overview

- Sample: 1 µg, Gut Microbiome Whole Cell Mix (ATCC)
- Depletion: No depletion, Supplier TF – Product ME, Supplier TF – Product RM, FastSelect –5S/16S/23S
- Library prep: QIAseq Stranded
- Sequencing: NextSeq 550
- Mapping: CLC Genomics Workbench

## FastSelect robustly removes rRNA, while the other supplier's kits do not



## FastSelect frees up reads for gene detection



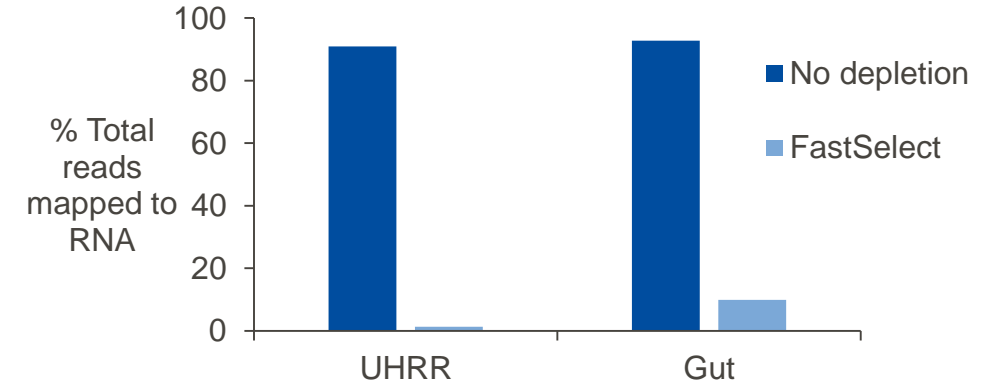
- FastSelect efficiently removes rRNA, while Supplier TF's rRNA depletion products do not
- The robust rRNA removal frees up a substantial amount of sequencing reads (9x compared to the other supplier's kits) for gene expression analysis

# FastSelect –5S/16S/23S plus –rRNA HMR remove bacterial and host rRNA

## Experimental overview

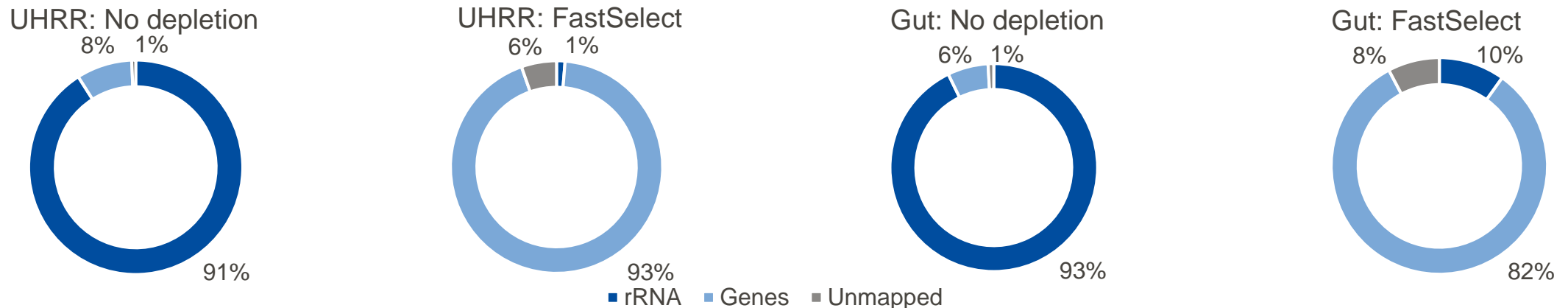
- Sample: 100 ng, different ratios of Universal Human (H) + Gut (G) RNA
- Depletion: No depletion, FastSelect –5S/16S/23S + FastSelect –rRNA HMR
- Library prep: QIAseq Stranded Total RNA Lib Kit
- Sequencing: NextSeq 550 (2 x 75 bp)
- Mapping: CLC Genomics Workbench

## FastSelect substantially removes rRNA



## FastSelect improves read mapping metrics, removing contaminating rRNAs

Breakdown of mapped reads

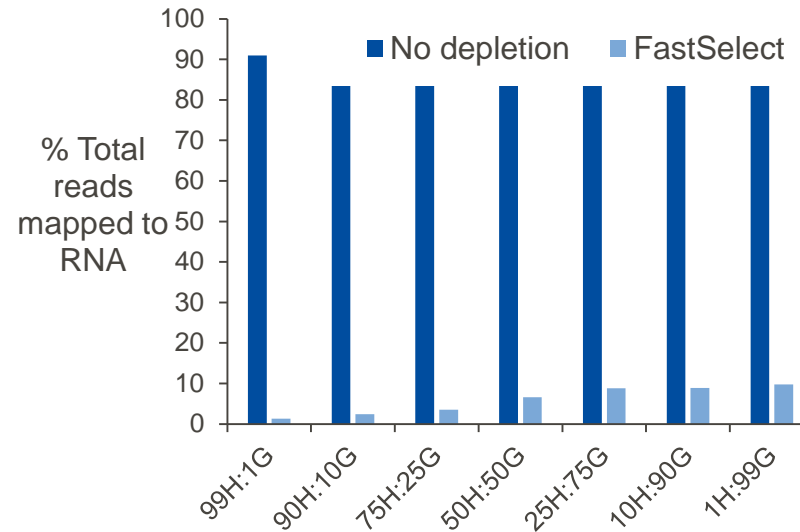


# FastSelect –5S/16S/23S plus –rRNA HMR remove bacterial and host rRNA

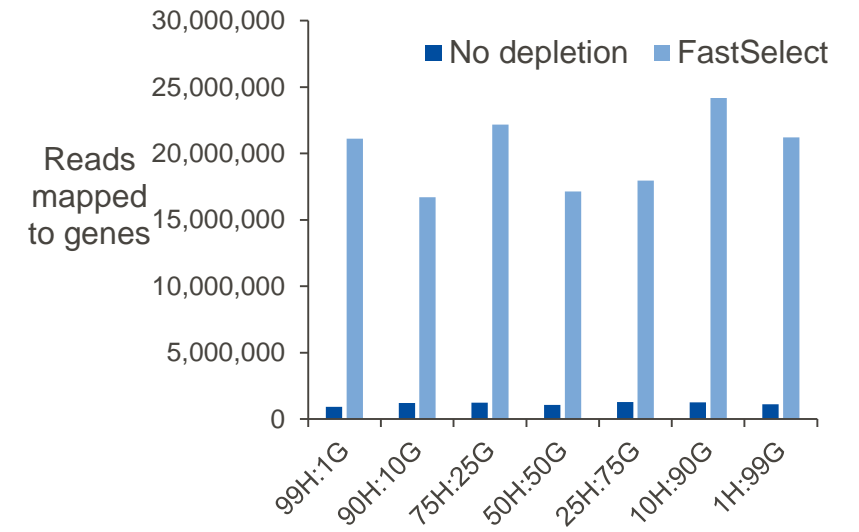
## Mixed samples

- 99% Human, 1% Gut (99H:1G)
- 90% Human, 10% Gut (90H:10G)
- 75% Human, 25% Gut (75H:25G)
- 50% Human, 50% Gut (50H:50G)
- 25% Human, 75% Gut (25H:75G)
- 10% Human, 90% Gut (10H:90G)
- 1% Human, 99% Gut (1H:99G)

## FastSelect –5S/16S/23S plus –rRNA HMR robustly removes rRNA



## FastSelect dramatically increases reads mapped to human + bacterial genes



- FastSelect –rRNA HMR + –5S/16S/23S efficiently removes rRNA from mixed human and bacteria samples
- This dramatically increases the reads mapped to human and bacterial genes (13- to 23-fold increase)

# Agenda

Background

---

FastSelect overview

---

FastSelect results: Mammalian, bacterial, mammalian + bacterial

---

Summary

---



## FastSelect is compatible with the QIAseq Stranded Total RNA Lib Kit

### Other compatible kits:

- TruSeq Stranded (Illumina)
- NEBNext Ultra II Directional (NEB)
- KAPA RNA HyperPrep (Roche)



- FastSelect is compatible with most RNA library prep kits.

# Summary

## FastSelect –rRNA HMR and –Globin

1

Removes cytoplasmic and mitochondrial rRNA and/or globin mRNA by inhibiting reverse transcription of specific targets

- Thirty percent faster than FastSelect V1, now in a single-tube
- Covers human, mouse, rat (HMR) and other mammalian species
- Compatible with several sample types, RNA types and input amounts and RNA library prep kits

## FastSelect –5S/16S/23S

2

Fragmentation and pan-bacterial (5S/16S/23S) rRNA depletion module which also works by inhibiting reverse transcription of specific targets

- Blocks community level cDNA synthesis of 5S, 16S and 23S rRNA
- Compatible with a range of RNA input amounts and most RNA library prep kits

## FastSelect –rRNA HMR plus –5S/16S/23S

3

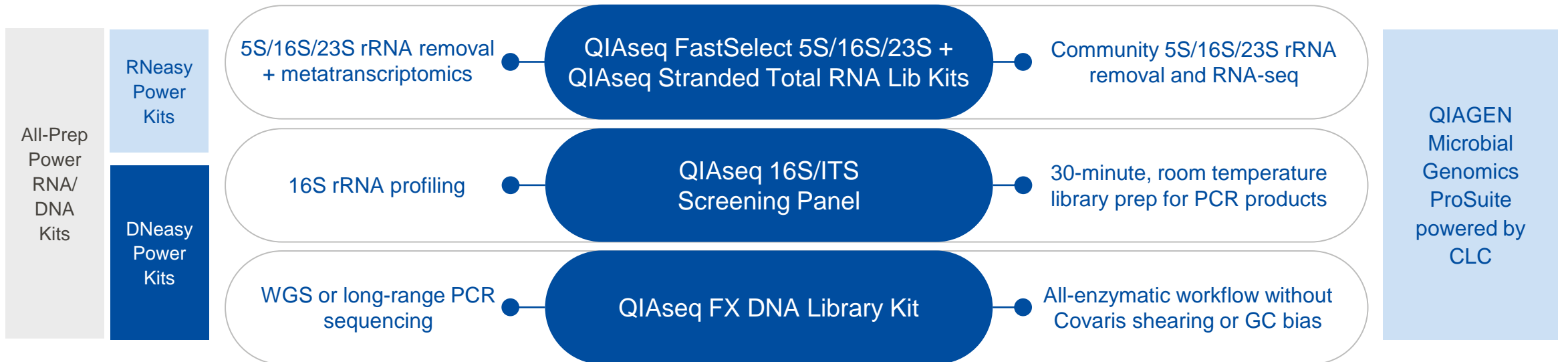
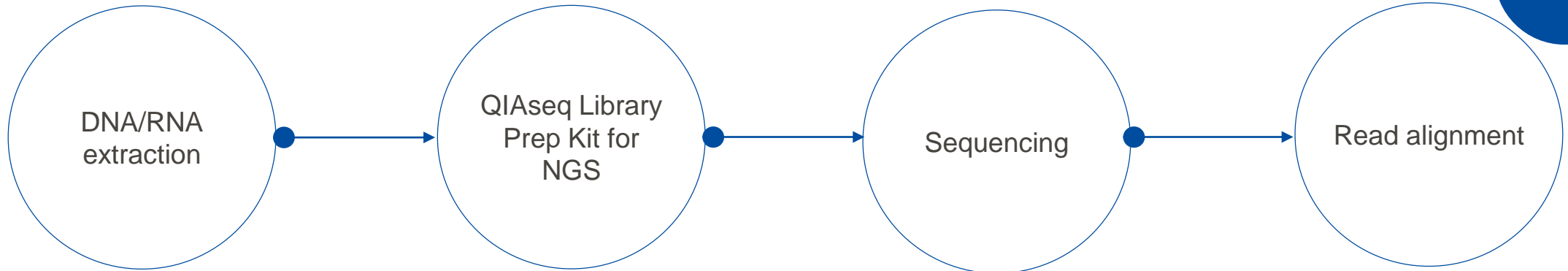
The FastSelect –rRNA HMR plus –5S/16S/23S protocol removes both human and bacterial rRNA


- One-hour protocol
- Expandable to include any species
- Dramatically improves sequencing read budgets by robustly removing rRNA



# Full QIAseq portfolio for metagenomics and metatranscriptomics

For  
Illumina





Thank you for attending.  
Questions?

Jonathan Shaffer, M.B.A., Ph.D.  
[Jonathan.Shaffer@qiagen.com](mailto:Jonathan.Shaffer@qiagen.com)