



Texas Children's Hospital

Rojelio Mejia<sup>1</sup>, Ashish Damania<sup>1</sup>, Rebecca Jeun<sup>1</sup>, Patricia Bryan<sup>1</sup>, Rubén Cimino<sup>2</sup>, Alejandro Krolewiecki<sup>2</sup>, Maria Elena Bottazzi<sup>1</sup>, Peter Hotez<sup>1</sup>, Ashley Luck<sup>3</sup>, Barton Slatko<sup>3</sup>

National School of Tropical Medicine, Baylor College of Medicine, Houston, TX<sup>1</sup>

Universidad Nacional de Salta, Salta Province, Argentina<sup>2</sup>

New England BioLabs, Inc. Ipswich, MA<sup>3</sup>



## Introduction

- >2 billion GI parasite infections worldwide
  - Poorest and resource-deprived communities
- Standard method of diagnosis: Stool microscopy
  - Sensitivity variable depending on prevalence, species, and concentration method
  - 50-90% sensitivity
  - Underestimates polyparasitism
- qPCR is rapid, quantitative, and high-throughput species-specific method
- GI parasites may disrupt normal intestinal microbiota
  - Decreased biodiversity is associated with disease
    - Malabsorption
    - Inflammatory bowel diseases

## Materials and methods

- Field site: Orán, Argentina
  - Peri-urban community
  - Temperate climate
- 99 patient samples
  - Asymptomatic children
  - Ages 2-10 years old
  - No recent antibiotics



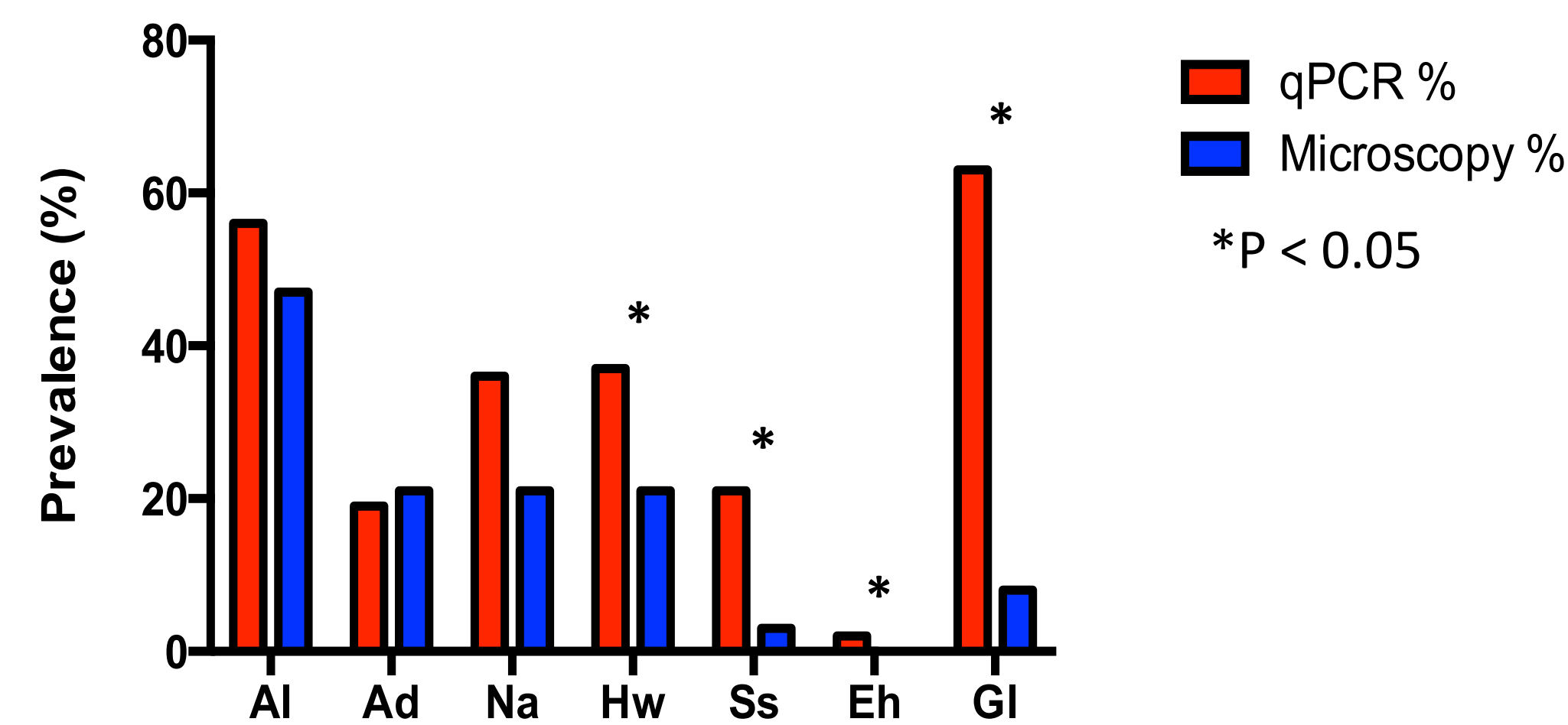
Stool samples evaluated by qPCR and microscopy for presence of :

- |                                   |                                       |
|-----------------------------------|---------------------------------------|
| <i>Ascaris lumbricoides</i> (Al)  | <i>Strongyloides stercoralis</i> (Ss) |
| <i>Ancylostoma duodenale</i> (Ad) | <i>Giardia lamblia</i> (GI)           |
| <i>Necator americanus</i> (Na)    | <i>Cryptosporidium</i> species (C)    |
| <i>Trichuris trichiura</i> (Tt)   | <i>Entamoeba histolytica</i> (Eh)     |

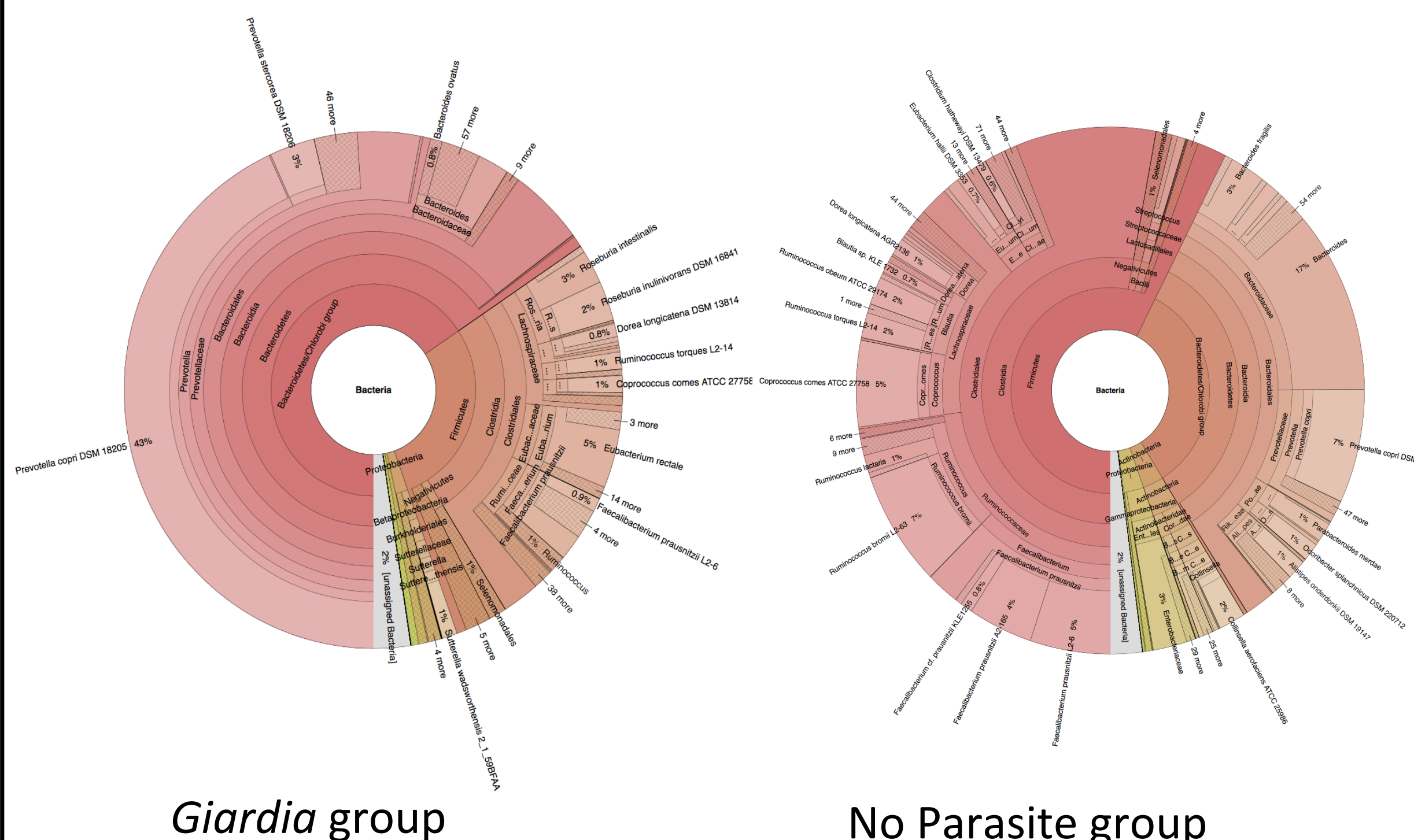
- qPCR required 50 mg samples
- Microscopy required 2 g samples
  - McMaster technique (semi-quantitative)
- NEBNext<sup>®</sup> Microbiome DNA Enrichment Kit
- NEBNext<sup>®</sup> Ultra<sup>™</sup> DNA Library Prep Kit for Illumina<sup>®</sup>
- Illumina MiSeq<sup>®</sup> "shotgun" sequencing

## Results

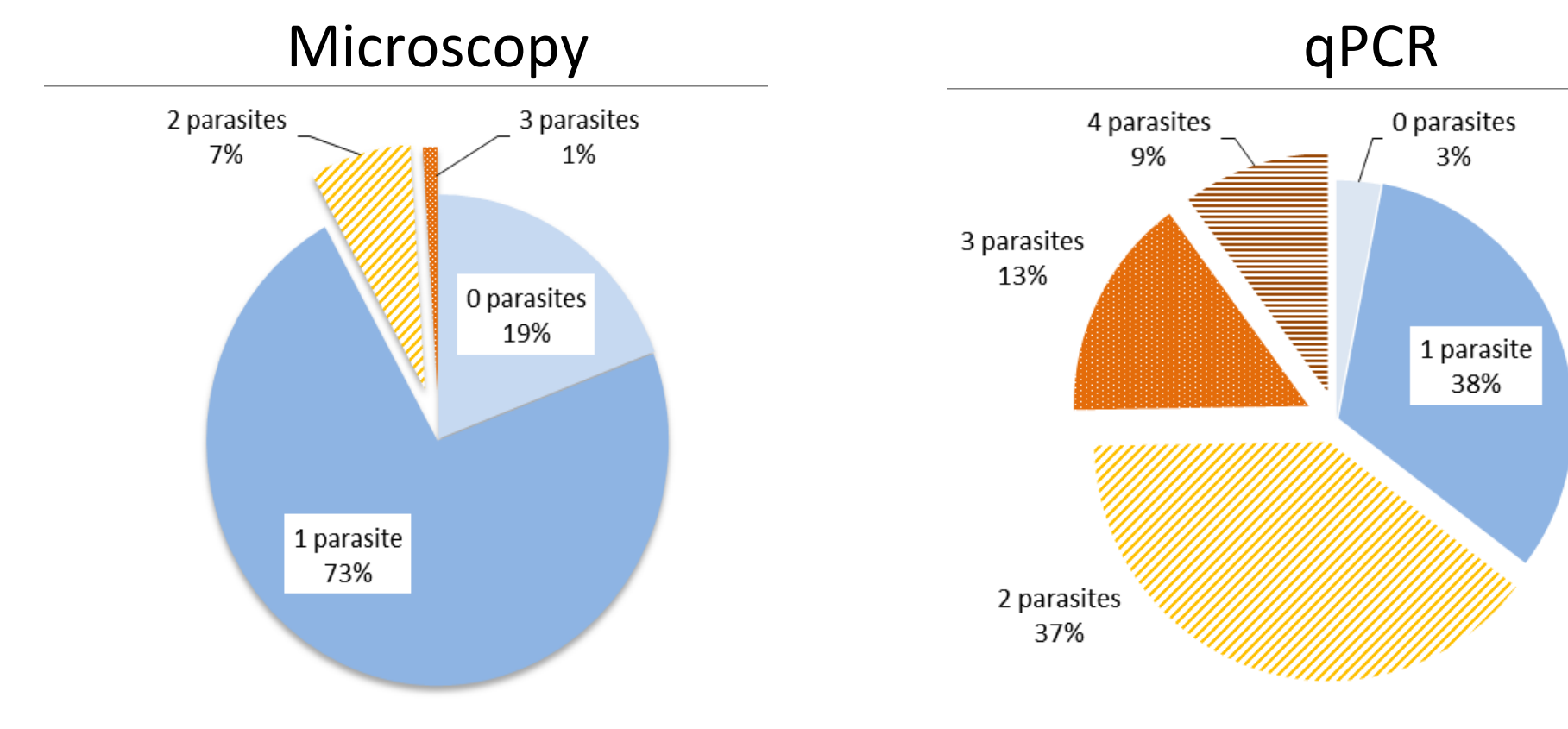
- qPCR (ITS region) (ABI 7500) identified more cases of *Ascaris*, hookworm (Hw), *Strongyloides*, *Entamoeba histolytica* and *Giardia* infection than microscopy. (Tt, C) no positives



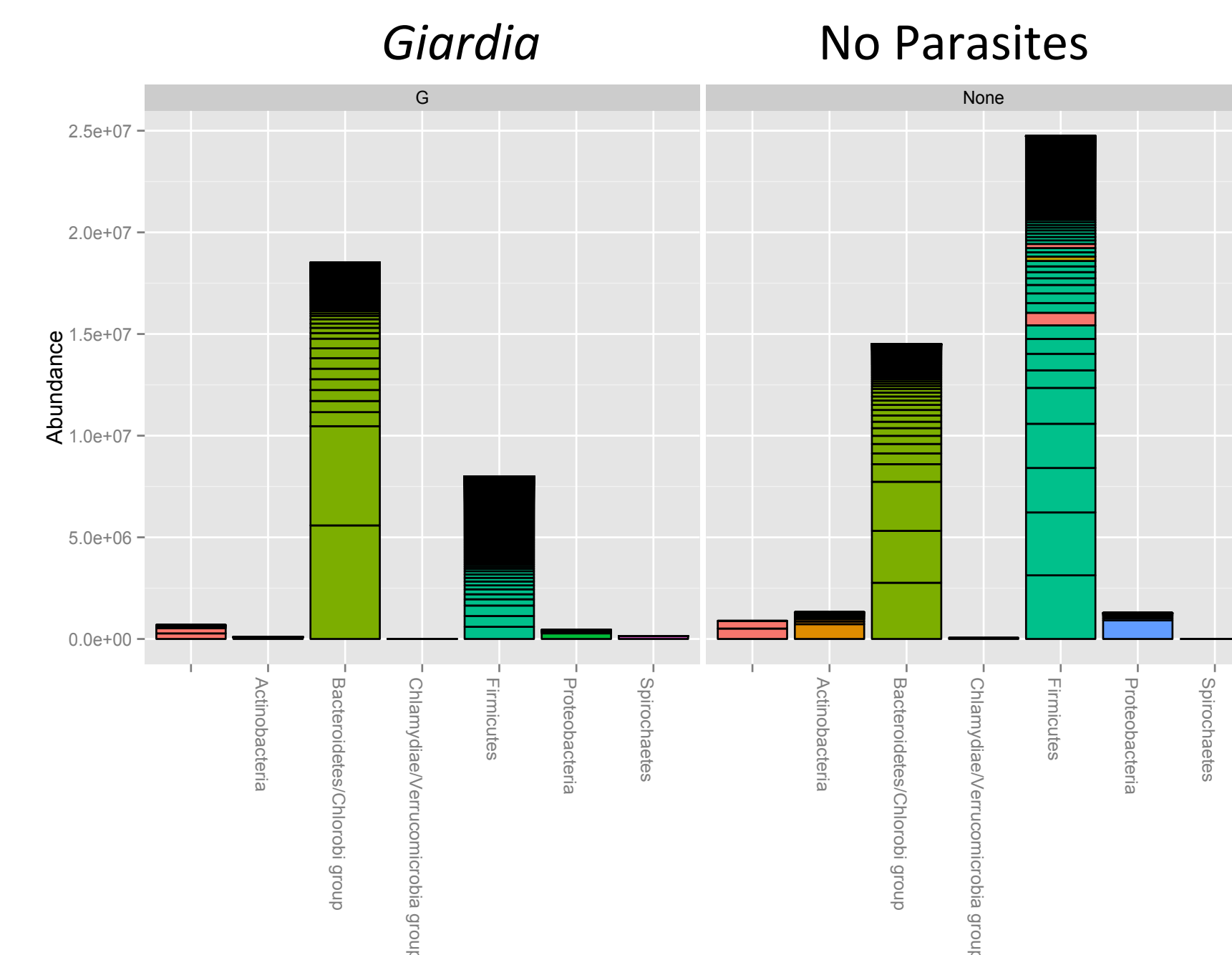
- Decreased intestinal bacterial biodiversity in combined *Giardia* group versus No parasite group



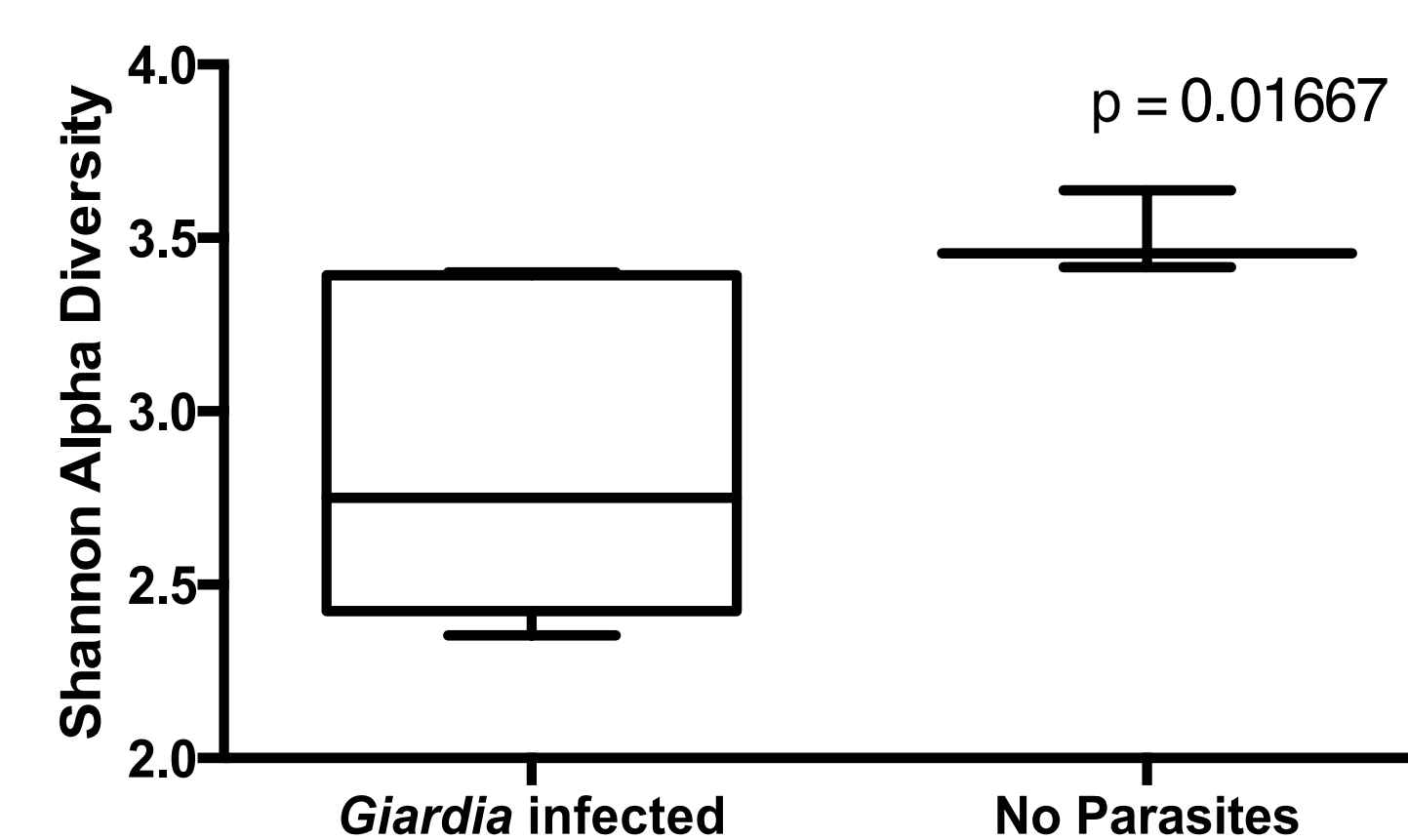
- qPCR identifies more polyparasitism than microscopy



- Giardia* infected group had higher abundance of *Bacteroidetes* compared to No Parasites group with higher *Firmicutes* (p < 0.05)



- Giardia* infection decreases intestinal bacterial biodiversity



## Conclusions

- qPCR can detect more parasites than microscopy
  - Ascaris* 91.3% Sens, 90.5% NPV
  - Hookworm 95.5% Sens, 98.4% NPV
  - Strongyloides* 100% Sens, NPV
  - Giardia* 87.5% Sens, 97.2% NPV
- qPCR can identify polyparasitism better than microscopy
  - Important for treatment selection
- GI parasitic infections at high prevalence
  - qPCR detected *Giardia* 6 x more than microscopy
- Giardia* infected group had decreased intestinal microbiota biodiversity (p = 0.01667)
  - Giardia* infected group (2.7)
  - No Parasite group (3.45)
- Giardia* infected group had significant increases in *Bacteroidetes* specifically *Prevotella* species
- Useful for epidemiology and morbidity studies
  - Surveillance after mass drug administration and vaccine programs
  - Expand understanding of morbidity and malnutrition
  - Cost is less than \$1.00 US per patient to screen for these parasites
- Future directions
  - Correlate quantity of parasite DNA with clinical outcomes
  - Associate morbidity to changes in microbiome
  - Treat children with anti-parasitics and evaluate changes in microbiome

## Acknowledgements

Funding for this project was provided by the National School of Tropical Medicine and New England BioLabs, Inc.