

DEVELOPMENT AND EVALUATION OF WHOLE CELL- AND GENOMIC DNA-BASED NEXT-GENERATION SEQUENCING (NGS) STANDARDS

Juan Lopera,* Monique Hunter,* Ray-Yuan Chuang,* Stephen King,* Megan Amselle,* Brian Chase,* Maria Mayda,* Kevin Zinn,* Samuel S. Minot,† Nicholas B. Greenfield,† and Dev Mittar.*

*ATCC, Manassas, VA, †One Codex, San Francisco, CA

ABSTRACT

ATCC has developed NGS Standards for use in a broad array of applications ranging from method optimization to data interpretation. These standards are fully sequenced, characterized, and authenticated mock microbial communities that mimic mixed metagenomic samples. They were developed as whole cell or nucleic acid preparations with even or staggered genomic DNA abundance, and medium or high levels of mock community complexity ranging from 10 to 20 strains per sample. Here, we explore the development and use of the NGS Standards.

INTRODUCTION

Advancement and accessibility of next-generation sequencing technologies have influenced microbiome analyses in tremendous ways, opening up applications in the areas of clinical, diagnostic, therapeutic, industrial, and environmental research. However, due to the complexity of 16S rRNA and metagenomic sequencing analysis, significant challenges can be posed by biases introduced during sample preparation, DNA extraction, PCR amplification, library preparation, sequencing, or data interpretation. Many researchers have published studies on these biases, and leaders in the microbiome field have highlighted the need for standardization.¹⁻⁴

One of the primary challenges in assay standardization is the limited availability of reference materials. To address these biases and provide a measure of standardization within microbiome research and applications, ATCC has developed a set of mock microbial communities, which includes lyophilized whole cells or genomic DNA, for use as NGS Standards in microbiome research. These standards mimic mixed metagenomics samples and comprise fully sequenced, characterized strains selected on the basis of select phenotypic and genotypic attributes. To further enhance the use of NGS Standards and eliminate the bias associated with data analysis, One Codex has developed a data analysis module in collaboration with ATCC that provides simple output in the form of true-positive, relative abundance, and false-negative scores for 16S rRNA community profiling and shotgun metagenomic sequencing.



DEVELOPMENT OF ATCC NGS STANDARDS

NGS Standards comprise fully sequenced, characterized strains selected on the basis of phenotypic and genotypic attributes, such as cell wall type (Gram stain classification), GC content, genome size, unique cell wall characteristics, and spore formation (Table 1). These standards were prepared as lyophilized whole cells or genomic DNA and were developed with even or staggered relative abundance and medium or high levels of mock community complexity (10 or 20 strains per sample) (Figure 1).

Table 1: Selection attributes for strains included in NGS Standards

Senus species	ATCC Number	Gram Status	Genome Size (Mb)	% GC	16S Copies	GenBank ID	Special Features	Microbiome
Bacillus cereus	10987™	Positive	5.42	35.2	12	NC 003909.8	Endospores former	Soil
Rifidobacterium dolescentis	<u>15703</u> ™	Positive	2.09	59.2	5	NC_008618.1	Anaerobe	Gut
Clostridium beijerinckii	35702™	Positive	6.49	30	14	NC_009617.1	Spores former	Gut/soil
Peinococcus radiodurans	BAA-816 [™]	Negative	3.29	66.7	7	NC_001263.1	Thick cell wall	Gut/environme
nterococcus faecalis	47077 tm	Positive	3.36	37.5	4	NC_017316.1	Biofilm producer	Gut
scherichia coli	<u>700926</u> ™	Negative	4.64	50.8	7	NC_000913.3	Facultative anaerobe	Gut
actobacillus gasseri	<u>33323</u> ™	Positive	1.89	35.3	6	NC_008530.1	Nuclease producer	Vaginal/gut
hodobacter sphaeroides	<u>17029</u> ™	Negative	4.60	68.8	3	NZ_AKVW01000001.1	Metabolically diverse	Aquatic
aphylococcus epidermidis	<u>12228</u> ™	Positive	2.56	31.9	5	NC_004461.1	Thick cell wall	Skin/mucosa
reptococcus mutans	<u>700610</u> ™	Positive	2.03	36.8	5	NC_004350.2	Facultative anaerobe	Oral
cinetobacter baumannii	<u>17978</u> ™	Negative	4.34	39	6	NZ_CP009257.1	Filaments, capsule	Environment
ctinomyces odontolyticus	<u>17982</u> ™	Positive	2.39	65.5	2	NZ_DS264586.1	Type 1 fimbriae	Oral
acteroides vulgatus	<u>8482</u> ™	Negative	5.16	42.2	7	NC_009614.1	Anaerobe	Gut
elicobacter pylori	<u>700392</u> ™	Negative	1.67	38.9	2	NC_000915.1	Helix shaped	Stomach/gut
eisseria meningitidis	BAA-335 [™]	Negative	2.27	51.5	4	NC_003112.2	Diplococcus	Respiratory tra
orphyromonas gingivalis	<u>33277</u> ™	Negative	2.35	48.4	4	NC_010729.1	Anaerobe, collagenase	Oral
opionibacterium acnes	<u>11828</u> ™	Positive	2.56	60	4	NC_006085.1	Aerotolerant anaerobe	Skin
seudomonas araeruginosa	<u>9027</u> ™	Negative	6.26	66.6	4	NC_009656.1	Facultative anaerobe	Skin
taphylococcus aureus	BAA-1556 [™]	Positive	2.82	32.8	6	NC_007795.1	Thick cell wall	Skin/repiratory
reptococcus agalactiae	BAA-611 [™]	Positive	2.16	35.6	7	NC_004116.1	Serogroup B	Vagina
Even Distrib	oution	Staggere	d Distribu	ıtion	■ Ac	inetobacter baumannii	■ Actinomyce:	s odontolyticus
Bacter.a	ATCC® MSA-1003™			■ Ra	ocillus cereus	■ Bacteroides		
					fidobacterium adolescen		■ Clostridium beijerinckii	
20				-	einococcus radiodurans	■ Enterococcu	•	
					■ Es	cherichia coli	■ Helicobacte	r pylori
					■La	ctobacillus gasseri	■ Neisseria me	eningitidis
, a				■ Pc	orphyromonas gingivalis	■ Propionibac	terium acnes	
10 Bacteria 2-ASW @201V	1000™	ATOS	LAGA 4001	TM	■ Ps	eudomonas aeruginosa	■ Rhodobacte	r sphaeroides
σ <u> </u>		AICC	MSA-1001		■ St.	aphylococcus aureus	■ Staphylococ	cus epidermidis
ATCC® MSA-2	2003				0.0	apriyes social dan sus	1.1.	

Figure 1: ATCC NGS Standards. $\underline{ATCC\ MSA-1000,^{TM}\ MSA-1001,^{TM}\ MSA-1002,^{TM}\ }$ and $\underline{MSA-1003,^{TM}\ }$ are genomic DNA standards, and ATCC $\underline{MSA-2002,^{TM}\ }$ and $\underline{MSA-2003,^{TM}\ }$ are lyophilized whole cell standards.

USING ATCC NGS STANDARDS TO EVALUATE PCR AMPLIFICATION, LIBRARY PREPARATION, AND SEQUENCING

To evaluate factors that contribute to biases associated with PCR amplification, library preparation, and sequencing, we performed an inter-laboratory comparison following the earth microbiome protocol, which targets the V4 region for 16S community profiling. The data revealed significant inter-laboratory variability in the number of true positives (70–95%) and false positives (83–100%) as well as in the relative abundances (Figure 2A). We also compared three different regions of the 16S rRNA gene (V1V2, V3V4, and V4) using the genomic DNA NGS Standards. The results revealed that only the V1/V2 region of the 16S rRNA gene was able to profile the bacteria to the species level (true positives = V1V2: 90–100%, V3V4: 90–95%, V4: 90–95%, along with significant differences in the expected verses observed relative abundances) (Figure 2B). Overall, the data clearly reveals the need for reference standards to standardize critical methods used in microbiome analyses.

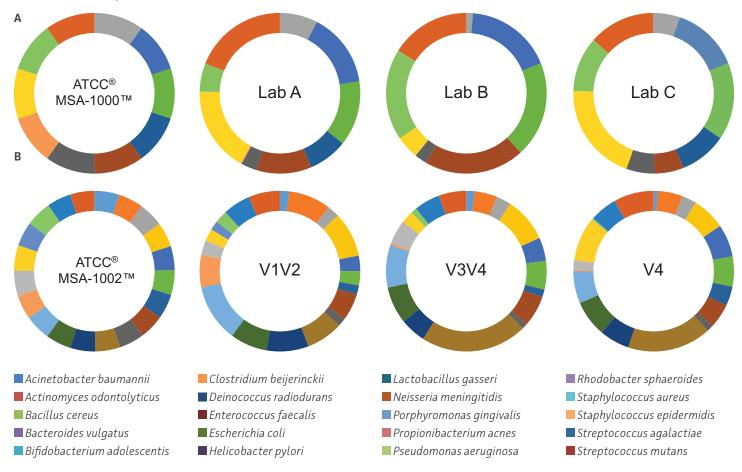
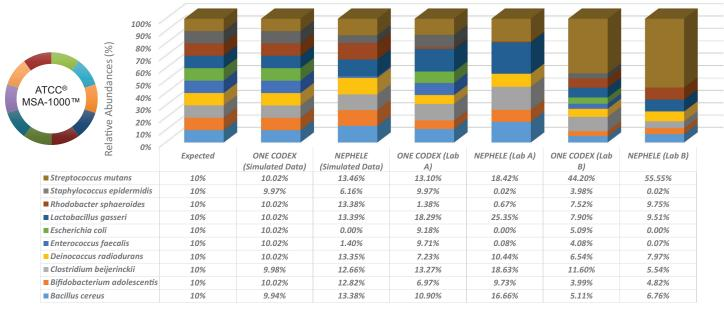


Figure 2: The use of standards during PCR amplification, library preparation, and sequencing. A) Inter-laboratory variations in identity and relative abundances. 16S rRNA V4 sequence data from different laboratories. Percent ratios of expected and observed organisms in the even genomic mock community comprising 10 organisms (ATCC MSA-1000). The blinded samples were sent to commercial vendors where they used their standard 16S protocol (Earth Microbiome Project) on the Illumina® platform. B) Choice of 16S rRNA primer regions affects identity and relative abundances. 16S rRNA community profiling results from the ATCC MSA-1002 standard using primer sets covering the V3V4, V1V2, and V4 regions on the Illumina® platform. The FASTQ files were analyzed using One Codex. For details on how to calculate the true-positive, false-positive, and relative abundance score, visit http://app.onecodex.com/atcc.

USING ATCC NGS STANDARDS TO EVALUATE DATA ANALYSIS

To evaluate biases associated with variations between data analysis platforms, we compared simulated data sets and two laboratory data sets that were produced using the ATCC MSA-1000 genomic DNA standard on the One Codex and NEPHELE⁵ data analysis platforms (Figure 3). The simulated data were generated by using a next-generation sequencing read simulator (ART)⁶ and the GenBank ID data and 16S rRNA copy number from each individual bacterial strain (Table 1). Here, all data were generated using primers against the V1/V2 region of the 16S rRNA gene. The results indicate that the One Codex platform identified all bacteria at the species level, while the NEPHELE platform identified bacteria at the genus level with wide variations in the relative abundances.



Analysis Platform

Figure 3: Data analysis platform affects identification and relative abundances. Simulated data sets and two laboratory data sets generated using the <u>ATCC MSA-1000</u> genomic DNA standard using the 16S rRNA (V1/V2) primer set were evaluated on the One Codex and NEPHELE (https://nephele.niaid.nih.gov) platforms.

COMBINING THE ATCC NGS STANDARDS WITH THE POWER OF ONE CODEX

To further enhance the use of ATCC NGS Standards and eliminate the bias associated with data analysis, we developed a data analysis module in collaboration with One Codex (https://app.onecodex.com/atcc) to provide simple output in the form of true-positive, relative abundance, and false-positive scores for 16S community profiling and shotgun metagenomic sequencing methods. Here, we compared the One Codex data analysis module side-by-side with three other commonly used analysis platforms. The results demonstrated significant variations among the number of true positives, the relative abundances, and the inability to identify all organisms at the species level (Figure 4). In contrast, the One Codex analysis tool, which was specifically customized for the ATCC NGS Standards, generated relative abundances close to the expected ratio.

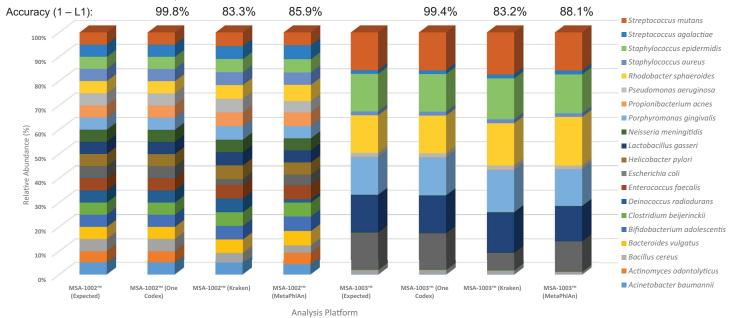


Figure 4: Data analysis platform affects identification and relative abundances. ATCC MSA-1002 and ATCC MSA-1003 were used to compare the performance of the One Codex, Kraken, and MetaPhlAn data analysis platforms. The percentages located above the bars represent the overall accuracy between platforms, as compared using L1-distance. Only One Codex demonstrated the accuracy necessary to robustly quantify microbiome sequencing errors.

CONCLUSIONS

Our data clearly reveals the need for standardization in microbiome analyses. Here, we demonstrate that bacterial identification and the evaluation of relative abundances in mixed samples can be affected by the 16S rRNA region chosen for amplification, general inter-laboratory differences, and variations between data analysis platforms. ATCC NGS Standards combined with the One Codex data analysis module provide a comprehensive solution for standardizing data from a wide range of sources, and generating consensus among microbiome applications and analyses.

REFERENCES

- 1 Wesolowska-Andersen A, et al. Choice of bacterial DNA extraction method from fecal material influences community structure as evaluated by metagenomic analysis. Microbiome 2: 19, 2014.
- 2 Brooks JP, et al. The truth about metagenomics: quantifying and counteracting bias in 16S rRNA studies. BMC Microbiol 15: 66, 2015.
- 3 Yuan S, Cohen DB, Ravel J, Abdo Z, Forney L J. Evaluation of methods for the extraction and purification of DNA from the human microbiome. PLoS One 7: e33865, 2012.
- 4 Wagner Mackenzie B, Waite D, Taylor M. Evaluating variation in human gut microbiota profiles due to DNA extraction method and inter-subject differences. Front Microbiol 6:130, 2015.
- 5 Office of Cyber Infrastructure and Computational Biology (OCICB), National Institute of Allergy and Infectious Diseases (NIAID). Nephele. http://nephele.niaid.nih.gov, 2016.
- 6 Huang W, Li L, Myers JR, Marth GT. ART: a next-generation sequencing read simulator. Bioinformatics 28(4): 593-594, 2012.

ACKNOWLEDGEMENTS

We would like to thank Dr. Stefan J. Green, Director, DNA Services Facility, University of Illinois at Chicago, for the generation of sequencing data and for his valuable input.

ABOUT ONE CODEX

One Codex is the leading bioinformatics platform for microbial genomics, supporting taxonomic and functional analysis of metagenomic (WGS), 16S, and other sequencing data. We specialize in creating robust, scalable, and secure bioinformatics solutions for metagenomics and microbial genomics, with a strong focus on ease of use. Founded in 2014, the One Codex platform counts thousands of users across leading academic institutions, biotechnology companies, and public sector organizations. One Codex is built on top of Amazon Web Services and is the only microbial genomics offering providing HIPAA-level security, as well as other strong compliance and audit guarantees. To learn more, visit http://www.onecodex.com.













©2023 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise. Droplet digital PCR and Bio-Rad are trademarks of Bio-Rad Laboratories, Inc.