

The Importance of Using Next-Generation Sequencing to Further Authenticate the ATCC Microbial Collections

Briana Benton Technical Manager, <u>ATCC</u>

Credible Leads to Incredible™



About ATCC

- Founded in 1925, ATCC is a not-for-profit organization with HQ in Manassas, VA, and an R&D and Services center in Gaithersburg, MD
- World's largest, most diverse biological materials and information resource for microbes – the "gold standard"
- Innovative R&D company featuring gene editing, microbiome, NGS, and advanced models
- cGMP biorepository

- Partner with government, industry, and academia
- Leading global supplier of authenticated cell lines, microorganisms, and molecular standards
- Sales and distribution in 150 countries, 19 international distributors
- Talented team of 450+ employees, over onethird with advanced degrees



Overview

Using next-generation sequencing to further authenticate the ATCC microbial collections

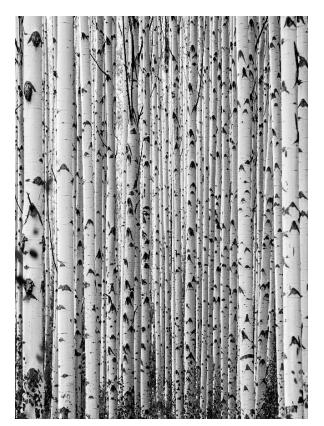
- Discuss why ATCC is committed to providing reference-quality genomes for items within the microbial collections
- Discuss some of the standardized processes and quality control criteria required for extracting, sequencing, and analyzing our reference-quality genomes
- Explore the ATCC Genome Portal





Providing reference-quality genomes

Why - Challenge # 1



- Public databases routinely host genomic data that is cited as "ATCC," but there often no traceability back to genuine ATCC curves an ATCC doesn't perform confirmation testing on public data.
 - How do researchers *know _____h data set to use?
 - Which is the "correct" he?
 - Close enough?
 - How do researchers have confidence in their selection?



Providing reference-quality genomes

Why - Challenge # 2



- How do we bring authentication into the genomics era while maintaining our commitment to our customers that we've fully and accurately authenticated our material?
- Typically, authentication* may refer to:
 - Morphology
 - Purity
 - Viability
 - Phenotypic testing
 - Genotypic testing
 - 16S ribosomal gene
 - ITS and D1D2

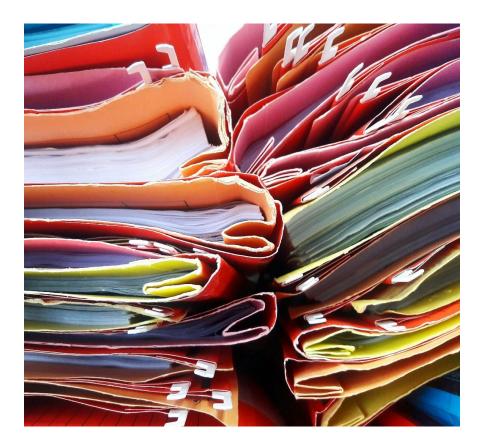


*not an inclusive list



Providing reference-quality genomes

Why - Challenge # 3



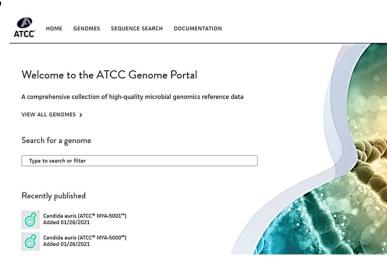
- Acknowledge there is a problem with reference genomes
- Work through a plan to address the problem
- How do we effectively and easily provide customers with genomic data while not diluting it or burying it in a public database?



The Enhanced Authentication Initiative

ATCC's solution to the authenticated reference genomes

- 2017-2018 Planning and proof-of-concept experiments
- 2018 Commitment
 - Laboratory and staffing resources
 - Instrumentation
 - Bioinformatics pipelines
- 2019 Launch of the Enhanced Authentication Initiative
 - June 2019 beta launch at ASM Microbe
 - Sept 2019 formal launch of the ATCC Genome Portal
 - Provide our customers with the whole-genome sequences of the specific, authenticated materials researchers need to generate credible data
 - o genomes.atcc.org





Overview

Using next-generation sequencing to further authenticate the ATCC microbial collections

Discuss why ATCC is committed to providing reference-quality genomes for items within the microbial collections

- Discuss some of the standardized processes and quality control criteria required for extracting, sequencing, and analyzing our reference-quality genomes
- Explore some of the features of the ATCC Genome Portal

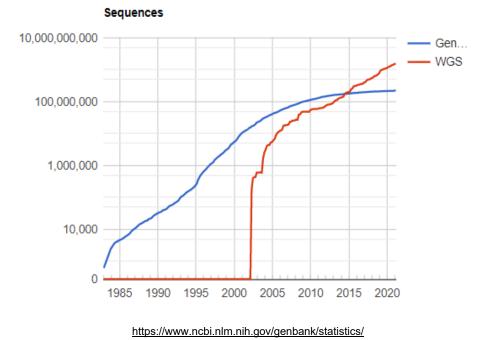




Reference genomes

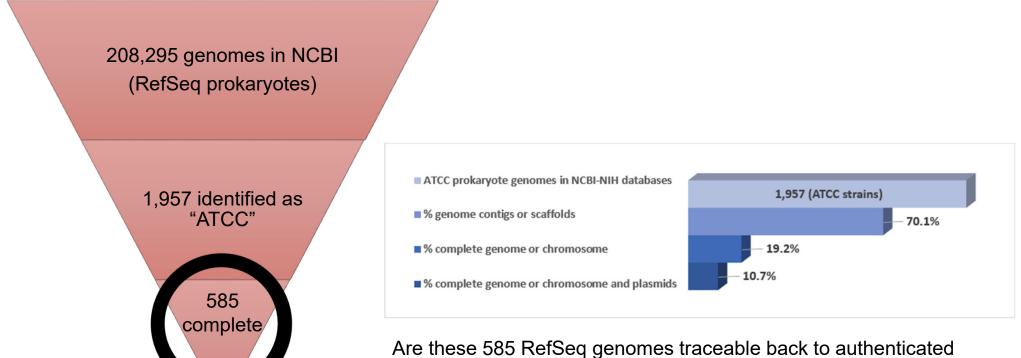
Where can researchers turn to for "reference" genomes?

- De facto standard
 - The sequence database for the entire public scientific community
 - Contains numerous genomes
 - Genomes submitted by a variety of labs
- Relatively little curation
- Highly variable quality
- NEVER authenticated by ATCC



ATCC

Reference genomes



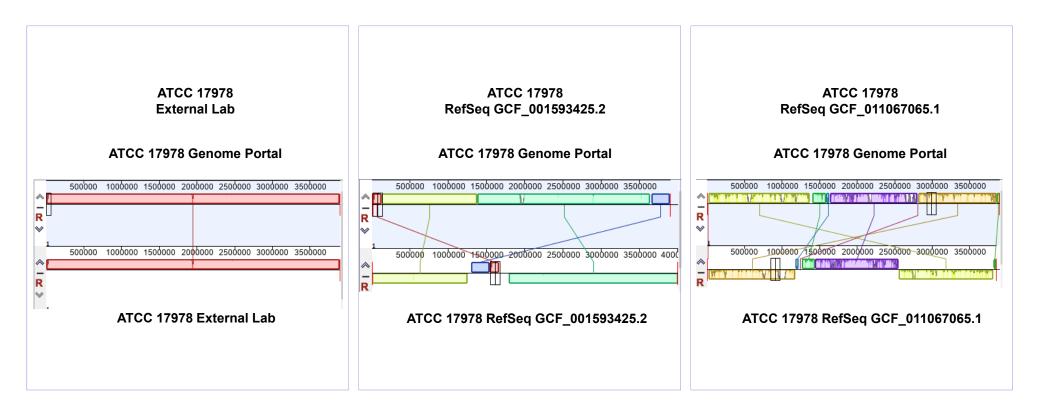
Are these 585 RefSeq genomes traceable back to authenticated ATCC cultures with well-documented growth and storage conditions?

ATCC[°]

Evaluation of genome sequences from public databases

Product	NCBI existing reference genomes	NCBI assembly level (plasmids)	Sequencing technology and coverage	# of SNPs	# of indels	Average coverage (variants)
5	GCA_001593425.2	Complete Genome	Illumina (300.0x)	14	5	210.1
	GCA_000015425.1*	Complete Genome (2)	Not available	118	656	152.7
	GCA_014672775.1	Complete Genome (1)	PacBio (399.24x)	15	87	170.4
Acinetobacter baumannii (ATCC [®] 17978™)	GCA_013372085.1	Complete Genome (2)	Illumina, Nanopore (80x)	14	2	210.2
	GCA_004797155.2	Complete Genome (2)	PacBio (247.19x)	28	62	162.1
	GCA_001077675.1	Complete Genome (1)	Illumina, PacBio (153x)	15	6	135.9
	GCA_011067065.1	Complete Genome (2)	PacBio (231.08x)	60227	2486	165.6
Candida albicans	GCA_015227795.1	3, 081 Contigs	NovaSeq (16x)	10174	1573	265.6
(ATCC [®] 10231™)	GCA_002276455.1	2,219 Scaffolds	HiSeq (95x)	13408	2390	274.6
Meyerozyma guilliermondii (ATCC® 6260™)	GCF_000149425.1	9 RefSeq Scaffolds	Not available	505	1973	278.2
	GCA_006942155.1	9 Contigs	ONT+MiSeq (240x)	74	386	223.3
Clavispora lusitaniae (ATCC [®] 42720™)	GCF_000003835.1	9 RefSeq Scaffolds	Not available	587	2336	265.6
	GCA_003675505.1	109 Scaffolds	NextSeq (182x)	102	5142	236.9

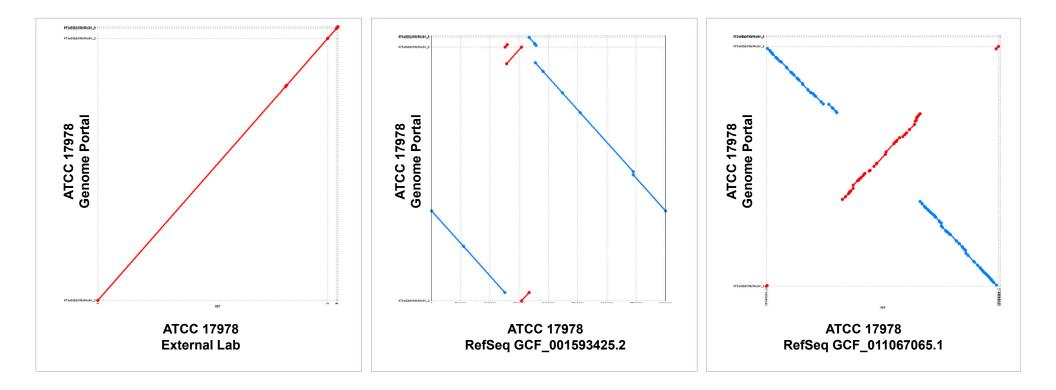
Evaluation of public sequences for ATCC 17978





Evaluation of public sequences for ATCC 17978

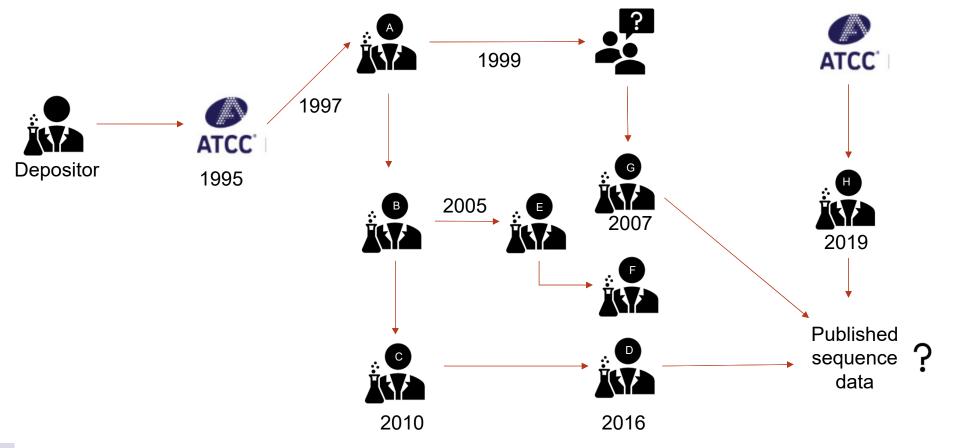
MUMmer alignment with the de novo ATCC 17978 versus GenBank RefSeq genome assemblies GCF_001593425.2 and GCF_011067065.1



ATCC°

13

Genomics data and a traceability and reproducibility crisis



ATCC[®]

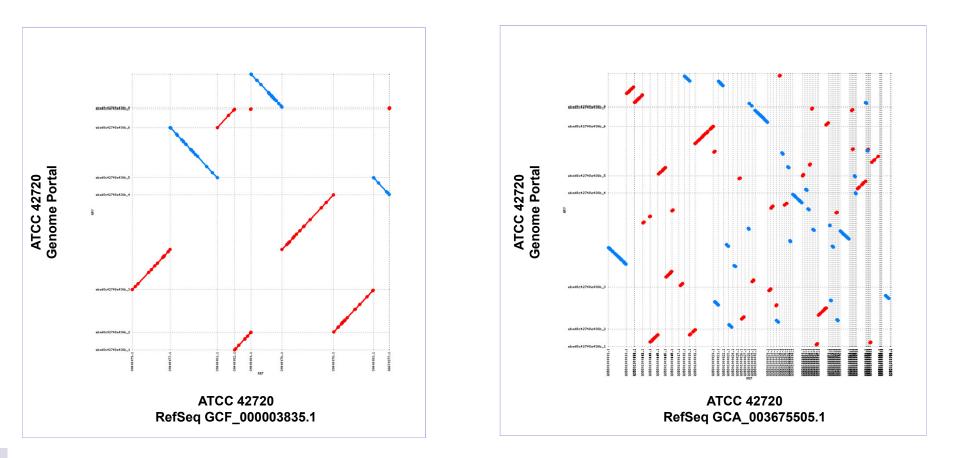
Evaluation of genome sequences from public databases

Product	NCBI existing reference genomes	NCBI assembly level (plasmids)	Sequencing technology and coverage	# of SNPs	# of indels	Average coverage (variants)
Acinetobacter baumannii (ATCC® 17978™)	GCA_001593425.2	Complete Genome	Illumina (300.0x)	14	5	210.1
	GCA_000015425.1*	Complete Genome (2)	Not available	118	656	152.7
	GCA_014672775.1	Complete Genome (1)	PacBio (399.24x)	15	87	170.4
	GCA_013372085.1	Complete Genome (2)	Illumina, Nanopore (80x)	14	2	210.2
	GCA_004797155.2	Complete Genome (2)	PacBio (247.19x)	28	62	162.1
	GCA_001077675.1 GCA_011067065.1	Complete Genome (1) Complete Genome (2)	Illumina, PacBio (153x) PacBio (231.08x)	15 60227	6	135.9 165.6
Candida albicans	GCA 015227795.1	3, 081 Contigs	NovaSeq (16x)	10174	1573	265.6
(ATCC [®] 10231™)	GCA 002276455.1	2,219 Scaffolds	HiSeq (95x)	13408	2390	274.6
Meyerozyma	GCF 000149425.1	9 RefSeq Scaffolds	Not available	505	1973	278.2
guilliermondii			ONT+MiSeq			
(ATCC [®] 6260™)	<u>CCA_000942155.1</u>	9 Contigs	(240X)	74	386	223.3
Clavispora lusitaniae (ATCC [®] 42720™)	GCF_000003835.1	9 RefSeq Scaffolds	Not available	587	2336	265.6
	GCA_003675505.1	109 Scaffolds	NextSeq (182x)	102	5142	236.9

ATCC[°]

Evaluation of public sequences for ATCC 42720

MUMmer whole genome alignments of ATCC de-novo genome assembly of ATCC 42720 versus GenBank RefSeq genome assemblies GCF_000003835.1 and GCA_003675505.1



Overview

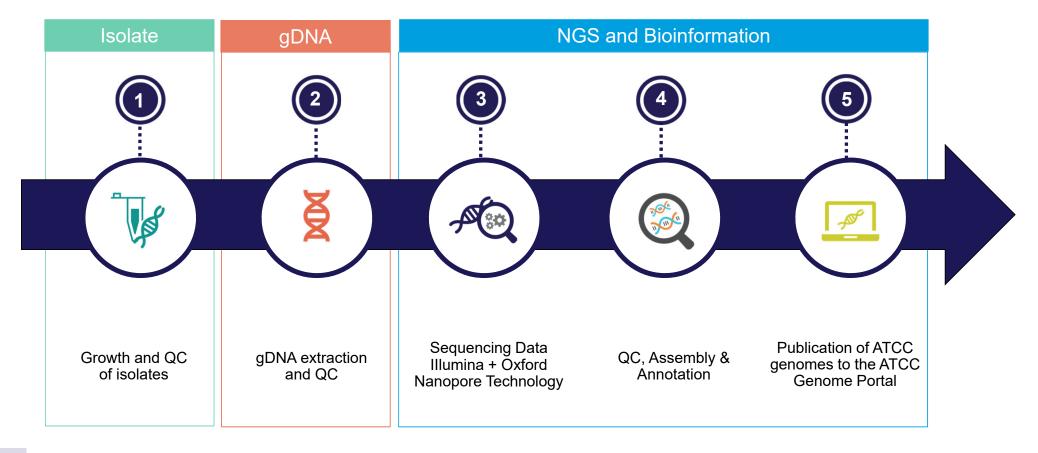
Using next-generation sequencing to further authenticate the ATCC microbial collections

- Discuss why ATCC is committed to providing reference-quality genomes for items within the microbial collections
- Discuss some of the standardized processes and quality control criteria required for extracting, sequencing, and analyzing our reference-quality genomes
- Explore some of the features of the ATCC Genome Portal



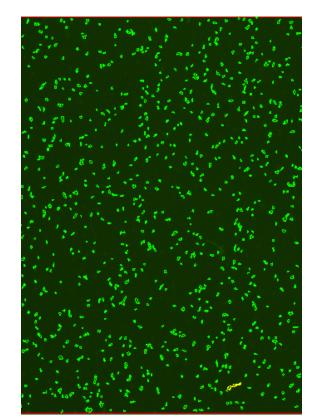


Authenticated physical material coupled with reference-quality genome sequences



Extraction of gDNA

- Start with a fresh culture grown according to ATCC's item-specific manufacturing process
- Determine the cell count
 - Typically start with ≥10⁹ cells/mL
- The "best" extraction method depends on the organism
- Simply recovering DNA is not good enough
 - -Concentration
 - Measured by Qubit[™] or Picogreen[®]
 - -Purity
 - o Measured with NanoDrop™
 - A_{260/280} ≥1.7 to ≤ 2.1
 - -Quality and Integrity
 - Fragment size is measured by Fragment Analyzer[™]



Fusobacterium nucleatum ATCC[®] 25586™ 6.58 x 10⁸ cells /mL



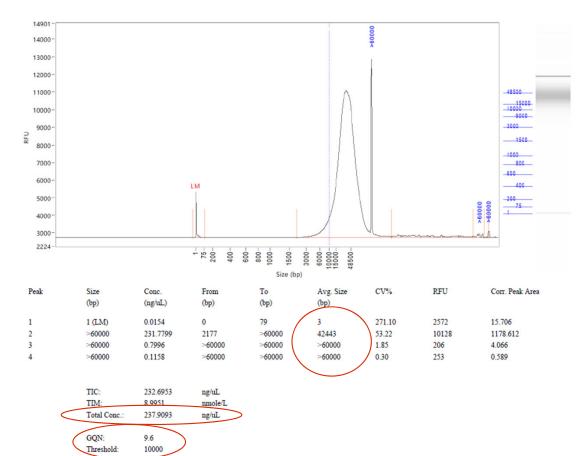
ATCC extraction quality control

ATCC [®] no.	Species	Qubit (ng/µL)	A _{260/} A ₂₈₀	DNA fragment size (range)**
8739™	Escherichia coli	101.9	1.92	49.5 kb (1.5 – >60 kb)
13048™	Klebsiella aerogenes	98.1	1.86	49.5 kb (1.6 – >60 kb)
11828™	Cutibacterium acnes	197.7	1.84	29.8 kb (0.8 – >60 kb)
6538™	Staphylococcus aureus	97.8	1.85	32.9 kb (2.7 – >60 kb)
BAA-2797™	Pseudomonas aeruginosa	153.3	1.99	44.1 kb (1.1 – >60 kb)
824™	Clostridium acetobutylicum	73.8	2.05	12.5 kb (4.6 – 57.8 kb)
6538™	Staphylococcus aureus	37.1	2.00	26.2 kb (6.9 – >60 kb)
27774™	Desulfovibrio desulfuricans	69.2	1.99	58.5 kb (13.3 – >60 kb)
11842™	Lactobacillus delbrueckii	64.8	2.02	41.9 kb (6.1 – >60 kb)
15697™	Bifidobacterium longum	76.2	1.95	51.3 kb (10.5 – >60 kb)

** Main peak reported



Fragment analysis of gDNA



- Corynebacterium tuberculostearicum (ATCC[®] 35692[™])
- Total concentration: 234 ng/µL
- Average fragment size: ≥42,000bp
- GQN: 9.6 with a threshold of 10,000bp
 - "Genomic Quality Number"
 - 96% of the sample contains fragments larger than 10,000 bp



Library preps for both Illumina[®] and Oxford Nanopore Technologies[®]

Illumina

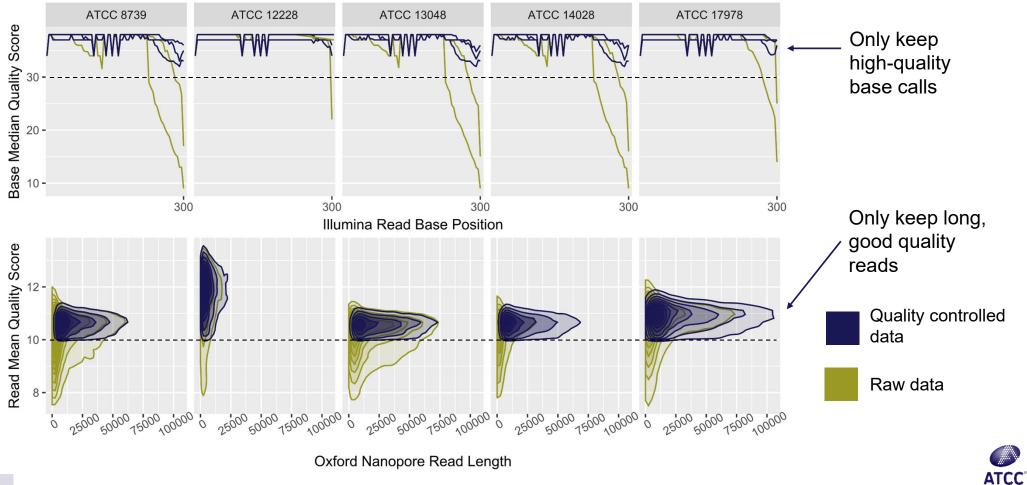
- DNA libraries are prepared using Illumina's DNA Prep kit and unique dual indexes (Cat. # 20018705)
- RNA libraries are prepared using NEBNext Ultra II RNA Library Prep Kit (Cat # E7770S)
- Sequenced on the MiSeq[®] or NextSeq[®] instrument
 - Paired-end read set per sample
 - Multiplexing is based on the estimated genome size
 - Data necessary to generate at least 100X coverage of the genome
- Reads are adapter trimmed using the adapter trimming option on the Illumina instrument

Oxford Nanopore Technologies

- Libraries are prepared using ONT's Ligation Sequencing Kit (SQK-LSK109) with the Native Barcoding Expansion kit (EXP-NBD104 or EXP-NBD114)
- Sequenced on the GridION using the version 9.4.1 flow cell
- The quantity of samples typically multiplexed is based on the estimated genome size of the given organism.
- Flow cells are run for 48-72 hours
- Barcode detection, demultiplexing, and barcode trimming are completed on the instrument, parallel to the run



Sequencing QC – Read trimming/filtering



QC Metrics for both Illumina and Oxford Nanopore Technologies

Illumina

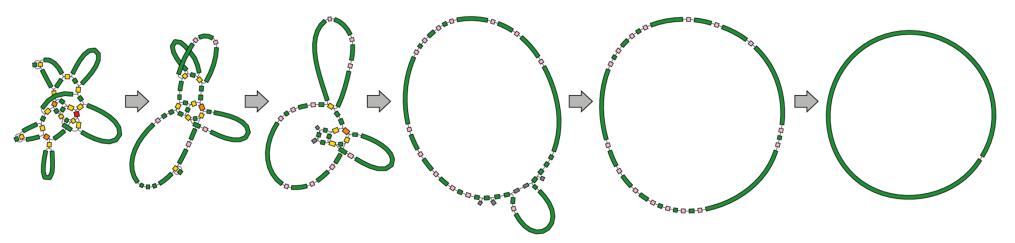
- Remove low-quality regions and adapter sequences
- This also ensures removal of adapter sequences otherwise missed by Illumina software
- Assess the quality of the read set by using FastQC
- Illumina reads must pass the following quality control:
 - Median Q score, all bases > 30
 - Median Q score, per base > 25
 - Ambiguous content (% N bases) < 5%

Oxford Nanopore Technologies

- ONT ultra-long reads are critical for scaffolding over the low-complexity regions of bacterial or fungal genomes during hybrid assembly, but they have limited influence in determining base identity given enough Illumina coverage.
- All data is trimmed and filtered for low-quality regions
- The quality control metrics used across all ONT read sets produced are:
 - Minimum mean Q score, per read > 10
 - Minimum read length > 5000
- To perform this quality control step, we employ NanoFilt on demultiplexed ONT read sets in addition to barcode sequence removal during demultiplexing



Hybrid genome assembly

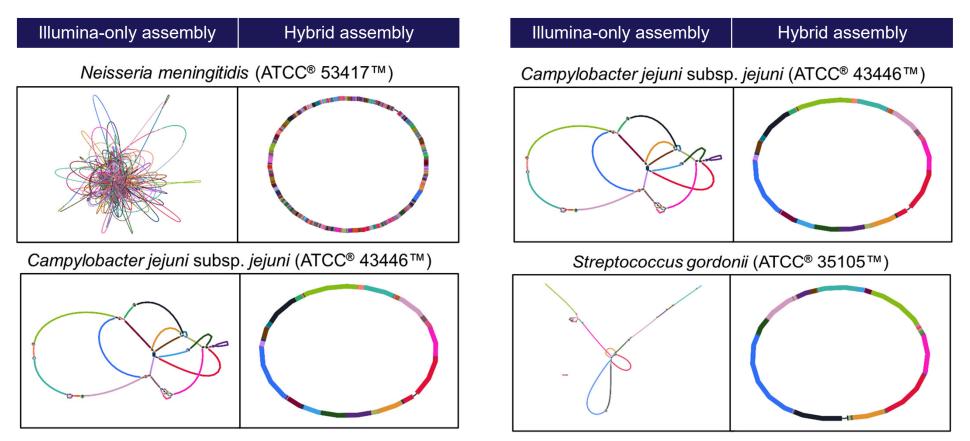


Illumina-only

genome	Long reads mapped to a tangled region creates a resolved bridge	Completed
assembly	Successively applying bridges resolves the structure of the genome	hybrid assembly
150 bp reads		

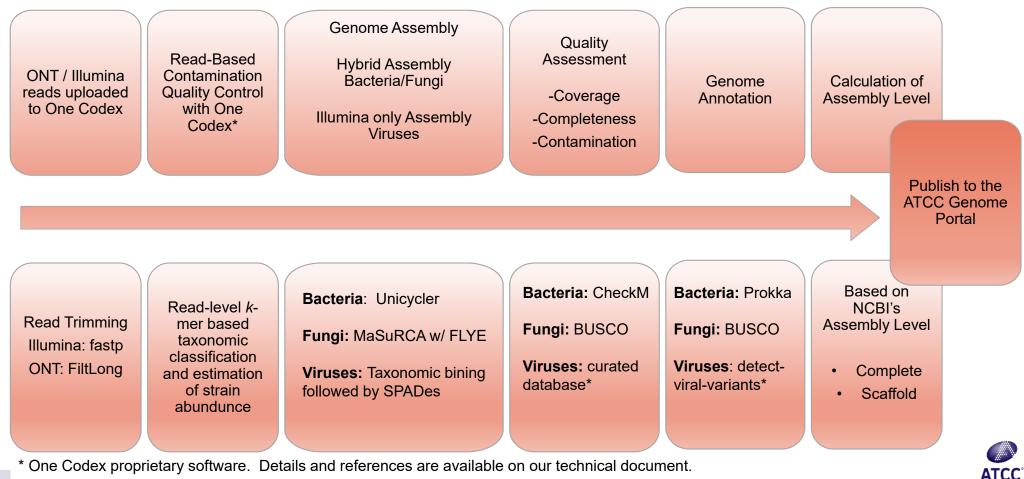


Advantage of hybrid assemblies



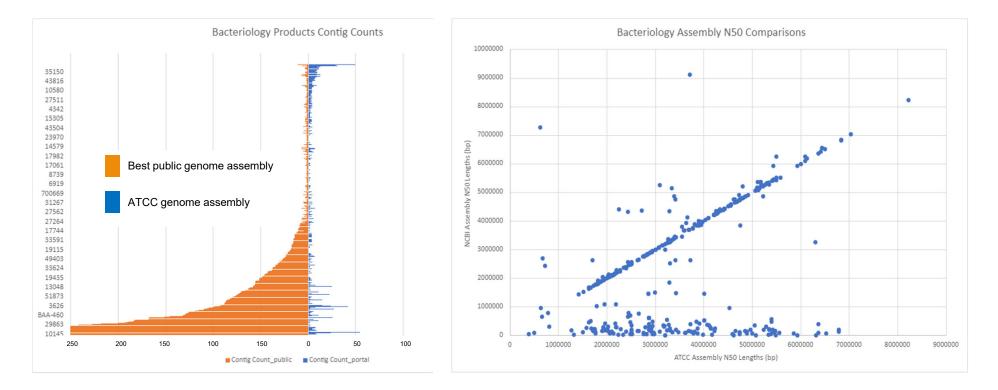
ATCC

ATCC genome assembly process



27

ATCC assemblies improve upon public assemblies



The **downward** trend in contig count and the **upward** trend in N50 indicate the ATCC produced genomes are of higher quality

Overview

Using next-generation sequencing to further authenticate the ATCC microbial collections

- Discuss why ATCC is committed to providing reference-quality genomes for items within the microbial collections
- Discuss some of the standardized processes and quality control criteria required for extracting, sequencing, and analyzing our reference-quality genomes
- Explore some of the features of the ATCC Genome Portal





ATCC Genome Portal

The ATCC Genome Portal is a cloud-based platform that enables users to easily browse genomic data and metadata by simply logging into the portal



Download whole-genome sequences and annotations of ATCC materials



Search for nucleotide sequences or genes within genomes

.....

View genome assembly metadata and quality metrics

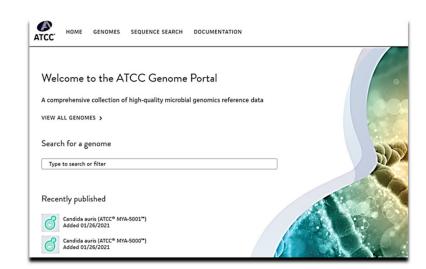
genomes.atcc.org



Summary

Using next-generation sequencing to further authenticate the ATCC microbial collections

- Discussed why ATCC is committed to providing reference-quality genomes for items within the microbial collections
 - traceability and reproducibility crisis
 - authentication in the genomics era
 - provide customers with easily accessible genomic data
- Discussed some of the standardized processes and quality control criteria required for extracting, sequencing, and analyzing our reference-quality genomes
 - gDNA extraction and QC
 - NGS library preps
 - Data QC
 - genome assembly process
- Explored the ATCC Genome Portal

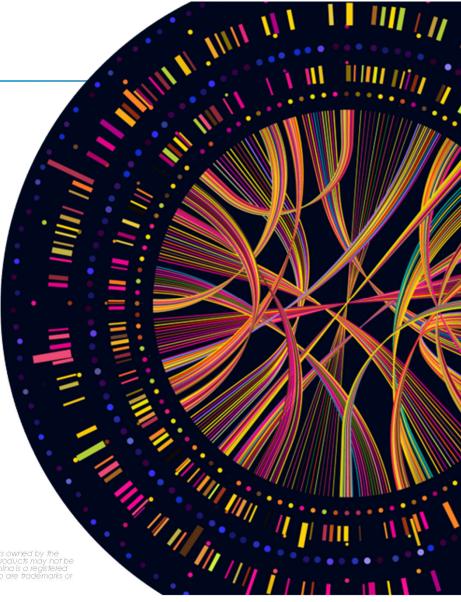




Acknowledgements

- Amanda Pierola, BS
- Brian Shapiro, PhD
- David Yarmosh, MS
- John Bagnoli, BS
- Juan Lopera, PhD
- Jonathan Jacobs, PhD
- P. Ford Combs, MS
- Nikhita Putheveetil, MS
- Samuel Greenfield, BS
- Stephen King, MS
- Our partners at One Codex

© 2021 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. These products are for laboratory use only. Not for human or diagnostic use. ATCC products may not be resold, modified for resale, used to provide commercial services, or to manufacture commercial products without prior ACC written approval. Illumina is a registered trademark of Illumina, Inc. Oxford Nanopore is a registered trademark of Oxford Nanopore Technologies Limited. PicoGreen, Qubit, and NanoDrop are trademarks or registered trademarks of Thermo Fisher Scientific. Fragment Analyzer is a trademark of Agilent Technologies, Inc.





Thank you

