Increasing Oxford Nanopore Technologies Throughput for the ATCC Genome Portal



Expanding Genome Publication Rates through an Improved Understanding of Bacterial Genome Assemblies

Credible leads to Incredible®

Samuel R. Greenfield, BS; Nikhita Puthuveetil, MS; James Duncan, BS; Amanda Pierola, BS; Corina Tabron, MS; David Yarmosh, MS; Amy Reese, MS; Stephen King, MS; Robert Marlow, BS; Ford Combs, PhD; Ana Fernandes, BS; John Bagnoli, BS; Briana Benton, BS; Jonathan Jacobs PhD ATCC, Manassas, VA 20110

Background

ATCC has made a commitment to the scientific community to provide reference-quality, whole-genome sequencing data for items within our collection through the ATCC Genome Portal. Our Enhanced Authentication Initiative enriches the characterization of our biological collections by using next-generation sequencing to provide reference-quality genomes to the scientific community; however, given the fact that we have only published 2,158 of our 78,700 items, a higher throughput method must be implemented to increase our publication rate. We sought to evaluate our current DNA requirements to determine if lower thresholds could generate similar quality bacterial assemblies, allowing for more rapid and automated DNA extractions and therefore an increased number of genomes being published to the ATCC Genome Portal.

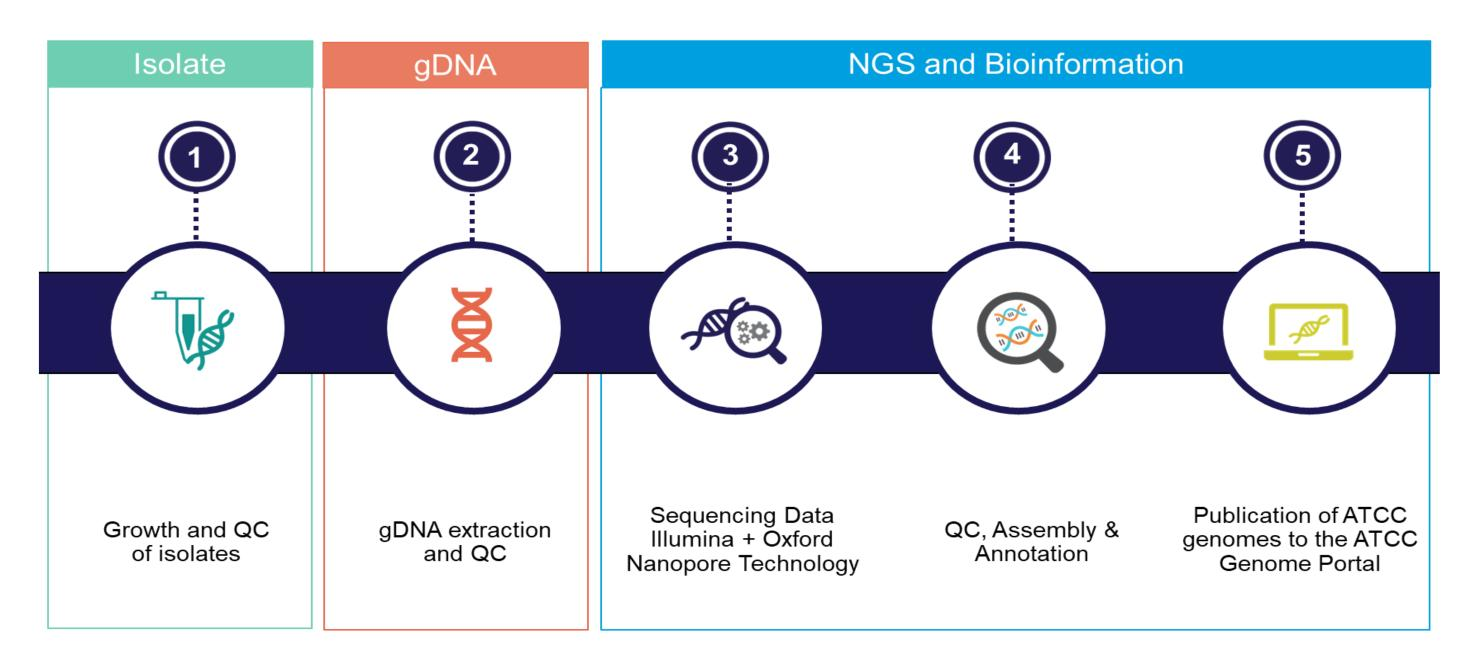


Figure 1. A hybrid assembly method combining Illumina® and Oxford Nanopore Technologies® (ONT) data to assemble the highest quality bacterial genomes is used.

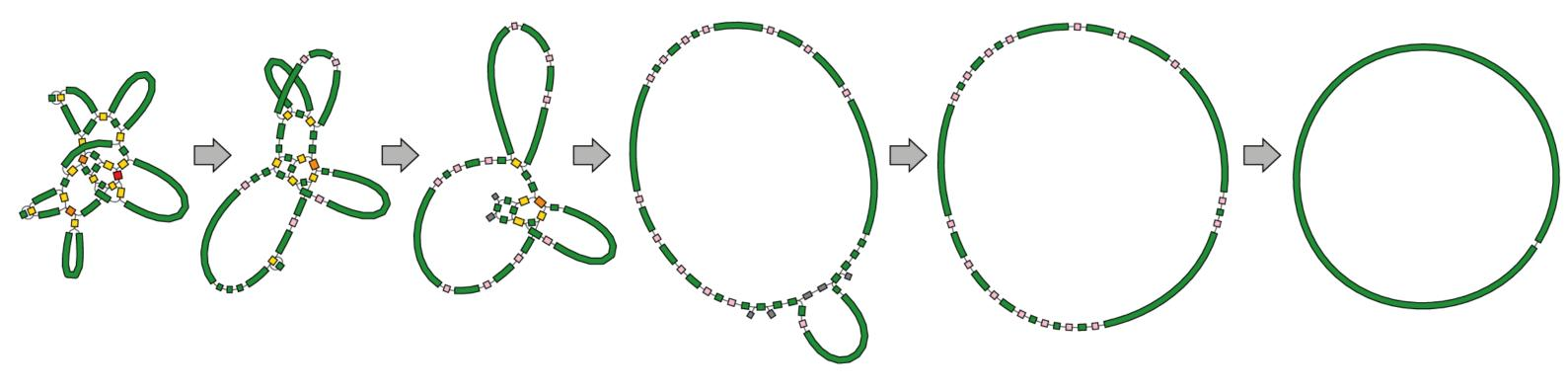
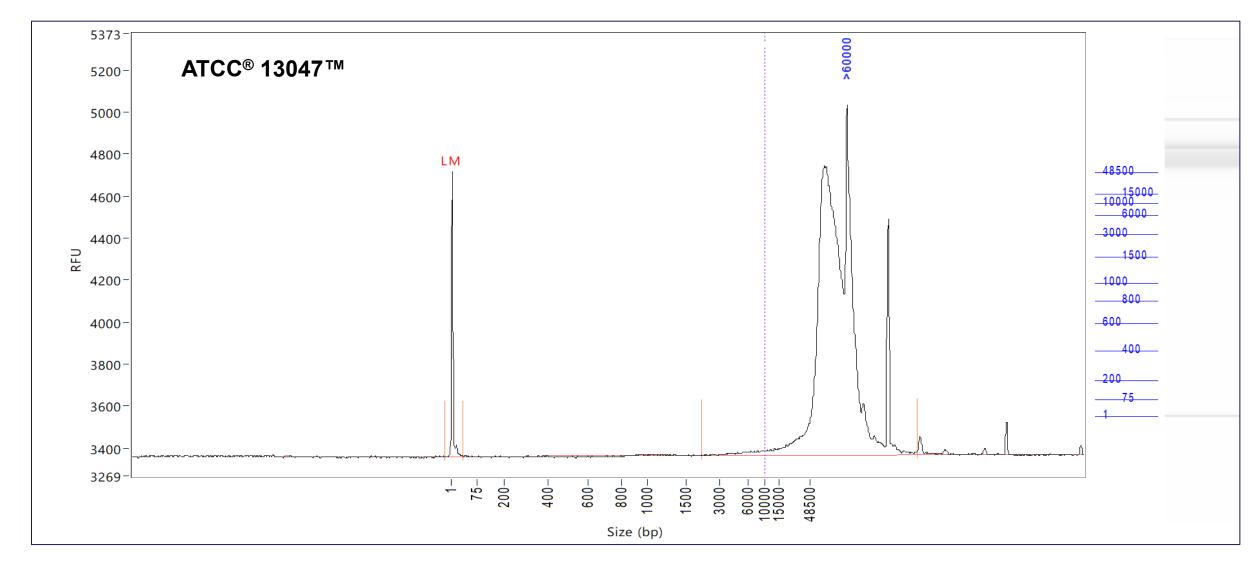


Figure 2. ONT long-reads are used to resolve tangled Illumina-only assemblies. To be published to the ATCC Genome Portal, bacterial assemblies are required to have <30 contigs, >100X Illumina depth, and both >95% completeness and <5% contamination analyzed by checkM.² Image reproduced from https://github.com/rrwick/Unicycler.



ATCC [®]	Organism	Qubit [®] (ng/μL)	A ₂₆₀ /A ₂₈₀ Ratio	Mean fragment size (kb)		
13047™	Enterobacter cloacae subsp. cloacae	32.7	1.85	>60		

Figure 3. We currently require DNA fragments for ONT sequencing to have an average fragment length of 20 kb, 50% of fragments to be larger than 10 kb, and at least 1000 ng of starting material. These standards are only consistently achieved using QIAGEN[®] Genomic-Tip (catalog no. 10223, QIAGEN, MD, USA), a time-consuming kit but in line with current ONT standards for the native barcoding and ligation sequencing with a R9.4.1 flow cell.

Methods

A hybrid assembly was performed using Unicyler³ with default parameters for each read set (see table 1) using filtered ONT and unfiltered Illumina reads. This constituted 63 separate assemblies. All ONT data was generated by native barcoding gDNA using LSK109 and NBD-104/114 on an R9.4.1 flow cell (Oxford Nanopore Technologies, UK). Illumina short read data was generated on a MiSeqTM from libraries created with Illumina's DNA Prep kit and unique dual indexes (catalog no. 20018705 and 20027213, Illumina, CA, USA).

Table 1. Illumina and Nanopore fastq files for nine bacterial samples published to the ATCC Genome Portal were gathered, evaluated using NanoPlot,¹ and filtered to maximum length thresholds of 5, 7.5, 10, 15, 20, 25-200 kb, and no threshold using Filtlong.⁴ Cells highlighted red failed to meet our previously stated publication standards while cells highlighted green met these standards.

ATCC®	Genome Size (bp)	5kb		7.5kb		10kb		15kb		20kb		25-200kb		No Max Filter	
		ONT Reads	N50	ONT Reads	N50	ONT Reads	N50	ONT Reads	N50	ONT Reads	N50	ONT Reads	N50	ONT Reads	N50
23114™	669,676	6,341	4,348	4,159	6,642	3,122	8,836	2,105	13,171	1,586	17,519	554	50,961	127,862	25,386
33453™	793,473	8,173	4,110	5,236	6,352	3,869	8,550	2,565	12,973	1,976	17,007	786	41,796	101,107	24,675
49145™	1,716,585	9,635	3,398	13,630	4,848	11,605	6,869	6,592	11,333	4,912	15,124	1,668	42,433	33,015	24,708
14799™	2,207,181	33,589	3,092	20,362	4,740	15,362	6,265	10,209	9,599	7,665	12,783	2,032	45,144	61,016	25,593
33236™	2,281,739	4,575	3,340	6,336	4,757	7,581	6,013	9,261	8,398	10,341	10,519	2,435	38,019	13,756	25,088
13047™	5,608,020	15,128	3,202	19,433	4,415	22,629	5,680	26,817	8,059	29,732	10,376	5,527	41,574	37,777	25,076
700441™	6,570,388	83,574	3,356	55,942	4,981	43,089	6,515	29,975	9,496	23,048	12,568	5,418	51,859	239,555	25,250
700084™	6,947,810	21,809	3,232	29,470	4,662	36,047	6,283	31,033	10,627	18,933	15,799	7,821	35,689	91,493	25,225
BAA-477™	7,070,224	82,821	2,955	78,957	4,215	49,360	6,352	29,131	10,465	21,281	14,370	6,274	47,351	158,833	25,417

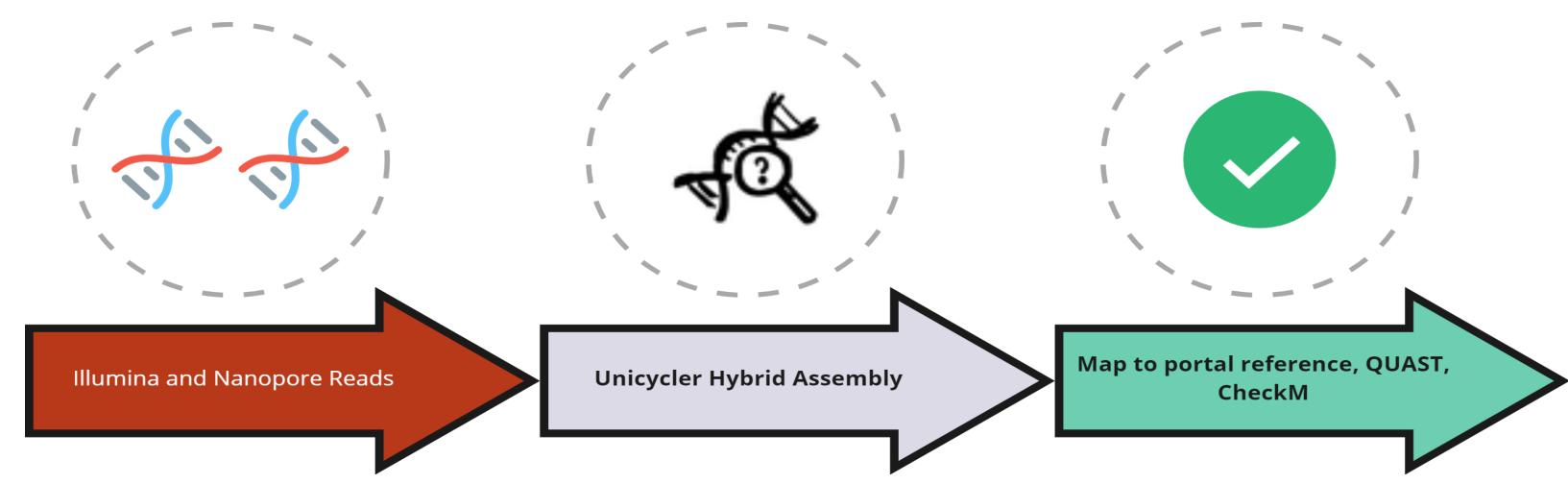


Figure 4. The resulting assemblies were evaluated by mapping assemblies to our own ATCC Genome Portal reference, running CheckM² for completeness and contamination metrics and QUAST⁵ for assembly metadata. Each was evaluated based on current ATCC Genome Portal publication requirements.

Results

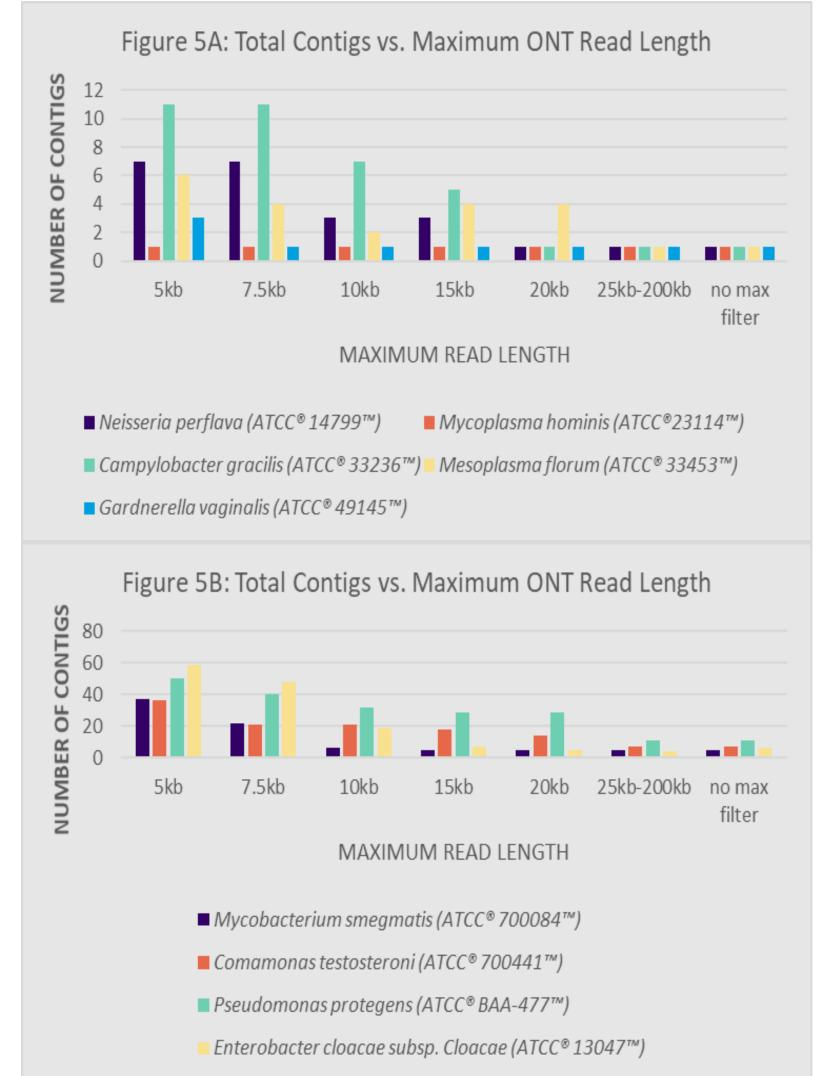


Figure 5(A&B). Read length with corresponding number of contigs for each bacterial genome assembly.

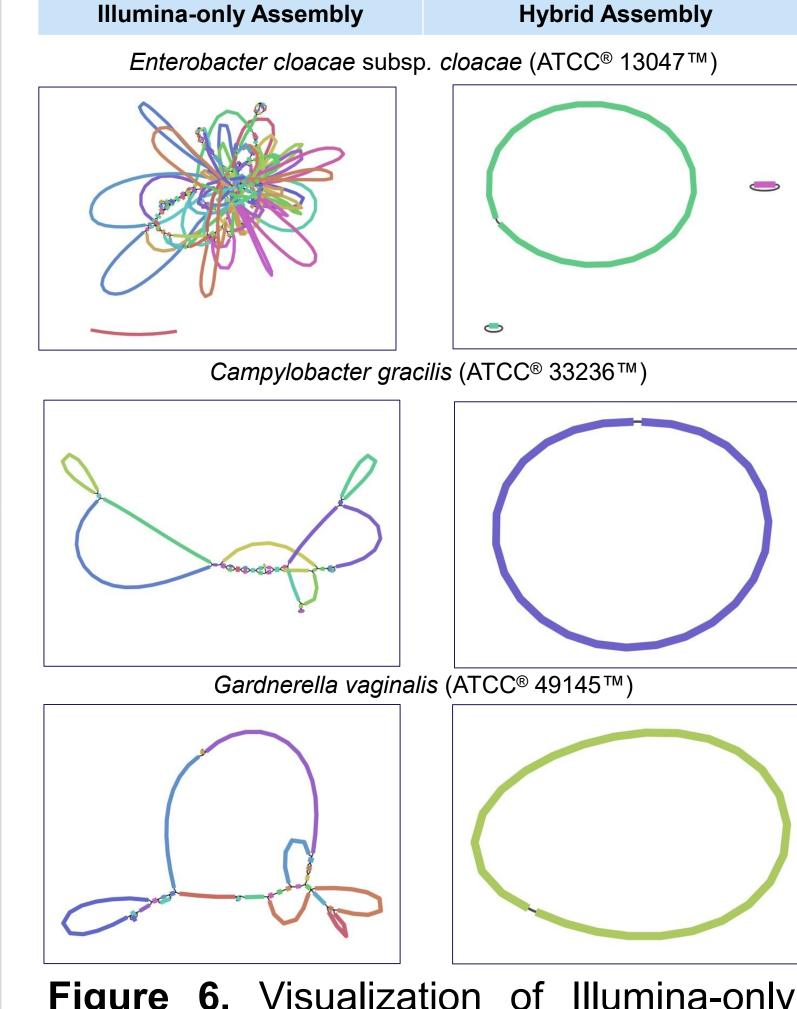


Figure 6. Visualization of Illumina-only and hybrid de novo assemblies. From our analysis, at ONT read lengths >8-12 kb, the hybrid assemblies meet ATCC's requirements for publication.

Our study provides in silico evidence to credit hybrid assemblies derived from smaller fragments. As seen by the filtered maximum 15 kb read N50s, we are confident that an average fragment length of >8-12 kb will provide successful high quality bacterial genome assemblies for ATCC's Genome Portal. With the decrease in average fragment length, we have been able to implement QIAGEN EZ1® extraction robots (QIAGEN, MD, USA), which have led to a 4X increase in extraction efficiency compared to the QIAGEN genomic-tip.

References

- 1. De Coster, Wouter (2016) NanoPlot [Source code]
- https://github.com/wdecoster/NanoPlot
- 2. Parks, D, Skennerton, C, Imelfort, M (2014) CheckM [Source code] https://github.com/Ecogenomics/CheckM
- 3. Wick, R (2016) Unicycler [Source code] https://github.com/rrwick/Unicycler
 4. Wick, R (2017) Filtlong [Source code] https://github.com/rrwick/Filtlong
- 5. Gurevich, A, et al. (2018) QUAST [Source code] https://github.com/ablab/quast



Learn more about our Enhanced Authentication Initiative

