MBP-FRIDAY-651 June 14, 2024



Does Long-Read Sequencing Technology ATCC[®] Produce Superior Viral Genome Assemblies? Credible leads to Incredible®

Corina L. Tabron, MS; Nikhita P. Puthuveetil, MS; Jade L. Kirkland, MS; Kaitlyn Gaffney, MS; Noah Wax, MS; James Duncan, BS; Robert Marlow, BS; Stephen King, MS; Ana Fernandes, BS; John Bagnoli, BS; Briana Benton, BS; Jonathan L. Jacobs, PhD | ATCC, Manassas, VA 20110

Background

Authenticated and traceable genomic data is vital for reproducible science, whether for preclinical studies, drug discovery, host-virus interaction, therapeutic development, or countless other applications. While the genomes for many of ATCC[®]'s viruses are available in public databases, as previously shown, these reference genomes published by third parties are often error-prone, incomplete, or generated using a variety of methods with a lack of supporting metadata, making downstream analysis challenging.¹ To address this problem, we developed reproducible next-generation sequencing and genome assembly workflows to produce viral assemblies for over three hundred viruses within our diverse collection. Our viral assemblies are generally produced with only short-read sequencing technology. Here, we set out to determine if we could produce higher-quality assemblies using long-read sequencing technology, starting with DNA viruses.

Table 2: Genome assembly length and contiguous DNA fragment counts were recorded for 11 DNA viruses produced by each assembly method.

| ATCC® | Virus | Assembly Length | | | Contigs | | |
|----------|---------------------|-----------------------|---------|---------|-----------|------|--------|
| | | Illumina [®] | ONT® | Hybrid | Illumina® | ONT® | Hybrid |
| VR-1890™ | Human adenovirus 1 | 11,076 | 33,987 | 35,783 | 15 | 1 | 1 |
| VR-5™ | Human adenovirus 5 | 7,481 | 35,872 | 35,725 | 5 | 1 | 1 |
| VR-1504™ | Human adenovirus 10 | 1,163 | 33,937 | 34,783 | 2 | 1 | 1 |
| VR-1097™ | Human adenovirus 20 | 18,385 | 34,753 | 35,729 | 23 | 1 | 2 |
| VR-1952™ | Human adenovirus 33 | 34,196 | 34,006 | 34,811 | 29 | 1 | 1 |
| VR-1407™ | Human adenovirus 49 | 34,623 | 35,122 | 34,915 | 26 | 1 | 1 |
| VR-1090™ | Human adenovirus D | 34,477 | 34,395 | 35,020 | 23 | 1 | 1 |
| VR-260™ | Human herpesvirus 1 | 111,863 | 115,246 | 134,147 | 102 | 2 | 6 |
| VR-733™ | Human herpesvirus 1 | 129,692 | 132,133 | 134,672 | 78 | 3 | 4 |
| VR-3393™ | Human herpesvirus 2 | 137,564 | 168,060 | 138,296 | 7 | 1 | 7 |
| VR-1781™ | Human herpesvirus 2 | 125,735 | 136,351 | 138,133 | 124 | 1 | 4 |



Figure 1: Standard ATCC[®] Genome Portal publication workflow.



Results

Figure 3: CheckV genome completeness comparison across three different assembly methods for 11 viral strains.



Methods

Table 1: Preliminary extraction kit selection. We currently generate viral assemblies from starting material extracted with the QIAGEN[®] QIAamp[®] Viral RNA Mini Kit (catalog no. 52904, QIAGEN[®], MD, USA). This method supports our minimum DNA concentration requirement of 1 ng for library preparation on Illumina[®] platforms. However, to integrate long-read sequencing into our viral pipeline, a method that yields concentrations ≥5 ng was required to be in line with our internal Oxford Nanopore Technologies[®] library preparation protocols. This was achieved with the QIAGEN[®] MinElute[®] Media Kit (catalog no. 57414, QIAGEN[®], MD, USA).

| | | Qubit [®] (ng/µL) | | | |
|----------|--------------------------|---------------------------------------|--|--|--|
| ATCC® | Organism | QIAamp [®] Viral Mini Kit | QIAamp [®] MinElute [®] Media Kit | | |
| VR-1090™ | Human adenovirus D | 3.67 | 20.7 | | |
| VR-197™ | Simian adenovirus | 1.30 | 9.05 | | |
| VR-3393™ | Human herpesvirus 2 | 1.19 | 7.29 | | |
| VR-1491™ | Human gammaherpesvirus 4 | 3.22 | 25.2 | | |





Figure 4: Graphical Fragment Assembly B (GFA) files were created by each assembler during the assembly process. The GFAs were then visualized using Bandage.⁴ (A) The GFAs for ATCC[®] VR-1890[™] and ATCC[®] VR-260[™] for each assembly method are shown above. Overall, the graphs created using only Illumina[®] sequencing reads indicate poor assemblies. The assemblies improve with the inclusion of ONT[®] reads or with only assembling with ONT[®] reads, as these graphs are much less complex and have

Α



Figure 2: Workflow used to produce viral assemblies. Isolated nucleic acids from viruses (see Table 1) were divided and processed for short-read sequencing on the Illumina[®] MiSeq[®] and for long-read sequencing on the Oxford Nanopore[®] GridION[®]. Unicycler² and Flye³ were used to generate individual de novo assemblies from the ONT[®] and Illumina[®] reads, as well as hybrid de novo assemblies from data produced on both platforms. Final assemblies were then evaluated for quality.

References

- 1.Yarmosh DA, et al. Comparative Analysis and Data Provenance for 1,113 Bacterial Genome Assemblies. 7(3): e0007722, 2022.
- 2.Wick RR, et al. Unicycler: Resolving bacterial genome assemblies from short and long sequencing reads. PLoS Comput Biol 13(6): e1005595, 2017.
- 3.Kolmogorov M, et al. Assembly of long, error-prone reads using repeat graphs. Nature Biotechnol 37: 540-546, 2019.
- 4. Wick RR, et al. Bandage: interactive visualisation of de novo genome assemblies. Bioinformatics, 31(20): 3350-3352, 2015.

fewer edges. (B) Close-up image of the tangled Illumina[®] -only assembly for ATCC[®] VR-260TM.

Conclusions

- Our findings show that the inclusion of long-read sequencing technology can improve genome assembly quality for DNA viruses as several strains assembled with long-reads exhibited a higher genome completeness and a total assembly length closer to what is expected for the strain.
- With the inclusion of long reads, assemblies had a drastic decrease in contig counts compared to their short-read only counterparts.
- Based on these results, utilizing both short-read and long-read technology in our viral assembly pipelines generates higher-quality assemblies.
- Further work will be conducted to optimize workflows dedicated to processing RNA viruses using long-read sequencing technology.

ATCC 10801 University Boulevard, Manassas, Virginia 20110-2209 Phone: 800.638.6597 Email: sales@atcc.org Web: www.atcc.org

© 2024 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. MiSeq and Illumina is a registered trademark of Illumina, Inc. Oxford Nanopore Technologies and GridION are registered trademarks of Oxford Nanopore Technologies plc. QIAGEN, MinElute, and QIAamp are registered trademarks of the QIAGEN Group. Qubit and Thermo Fisher Scientific are registered trademarks of Thermo Fisher Scientific, Inc.