

ATCC®

Credible leads to Incredible®

ATCC Genome Portal: Fungal Genome Assembly Improvements and Mitochondrial DNA Recovery

Emily A. White, MS; Joseph R. Petrone, PhD; Scott V. Nguyen, PhD, Nikhita P. Puthuveetil, MS; David A. Yarmosh, MS; Amy L. Reese, MS; John Bagnoli, BS; Briana Benton, BS; Jonathan L. Jacobs, PhD
ATCC, Manassas, VA 20110

Abstract

The utility of genomic data is constrained by the tools, assumptions, and reference quality available at the time the data is generated; consequently, many datasets contain biological signals that remain obscured until advances in analytical methods enable strategic reanalysis. Fungal mitochondrial genomes are a case in point: high-quality references now include 3,131 JGI-annotated and 892 complete RefSeq mitogenomes, yet 6,467 undeclared mitogenomes remain embedded in GenBank nuclear assemblies.¹ Mitogenomes further complicate analysis through extreme size variation (~2 kb to >200 kb), mixed architectures, absence of universal marker genes, and nuclear insertions (NUMTs).¹ Here, we compare 37 authenticated ATCC® Genome Portal (AGP) fungal assemblies against strain-matched RefSeq genomes using standard quality metrics. Because mitochondrial representation remains inconsistent across public databases, we assessed mitogenome integration in a subset of ATCC® assemblies using curated RefSeq references and explore the impacts of implementing a targeted workflow to separate mitochondrial and nuclear reads prior to assembly.

Methods

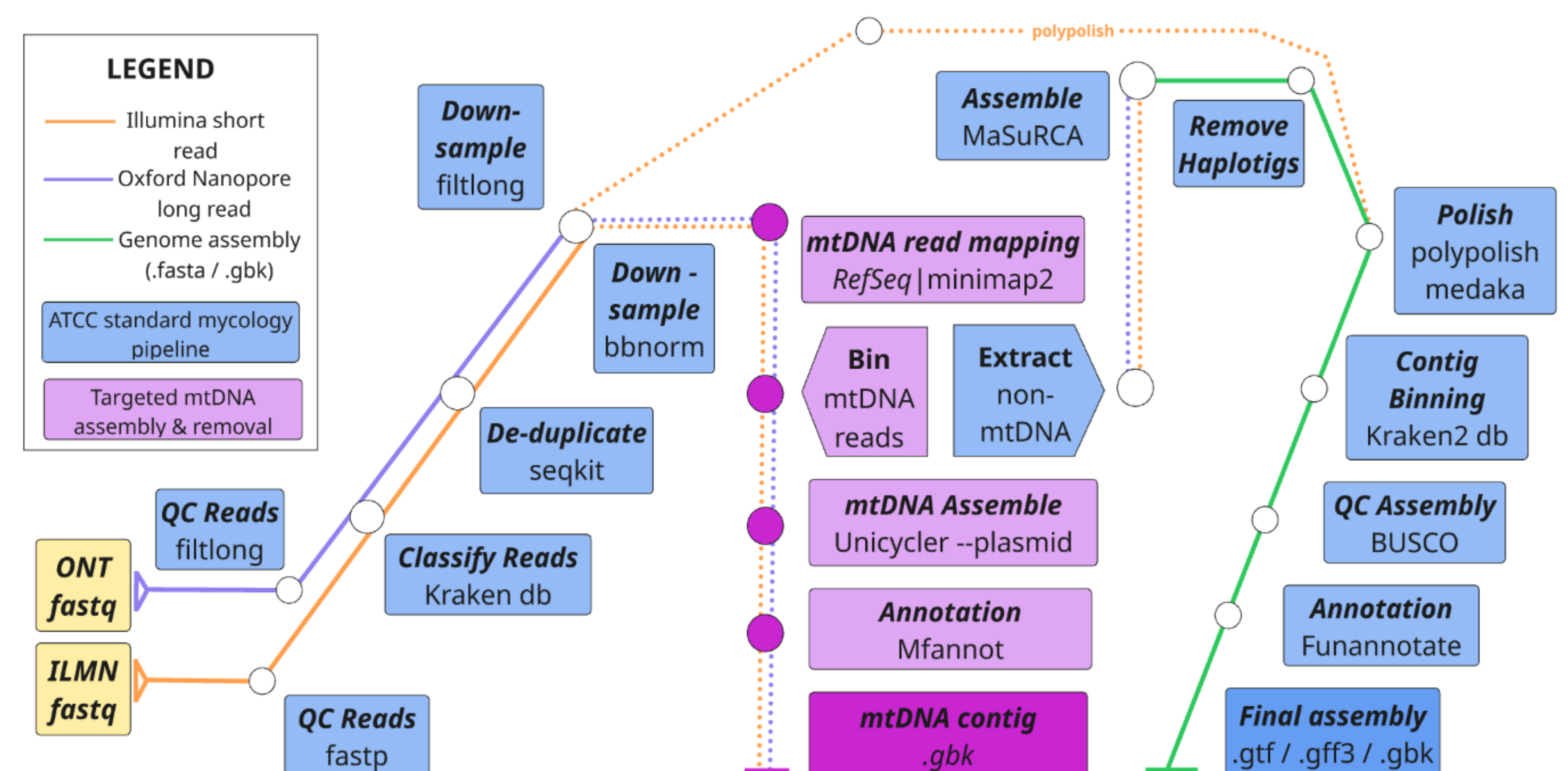


Figure 1: AGP fungal genome assembly pipeline highlighting targeted mitochondrial extraction (non-standard workflow). A de novo hybrid assembly workflow (blue) with a derivative mitochondrial extraction pipeline (pink). After downsampling, Illumina and ONT reads (dotted lines) branch to support parallel processing: mitochondrial reads are identified via alignment to strain-matched RefSeq mitogenomes, extracted, and assembled independently, while mitochondrial-mapped reads are removed from nuclear read sets. Following mtDNA extraction, Illumina and ONT reads are combined in a hybrid MaSuRCA assembly. A secondary dotted orange branch from the initial pre-processed Illumina reads represents downstream polishing (e.g., Polypolish).

Results: Mitogenome

Table 1: *Rhodotorula mucilaginosa* (ATCC® 2503™) assembly contiguity and BUSCO completeness metrics before and after mtDNA removal.

Assembly Version	Contigs (n)	N50 (nt)	Genome Size	Completeness
ATCC® 2503™ [standard]	24	1,614,595	20.4Mb	92.70%
ATCC® 2503™ [targeted]	28	1,270,344	20.6 Mb	96.30%

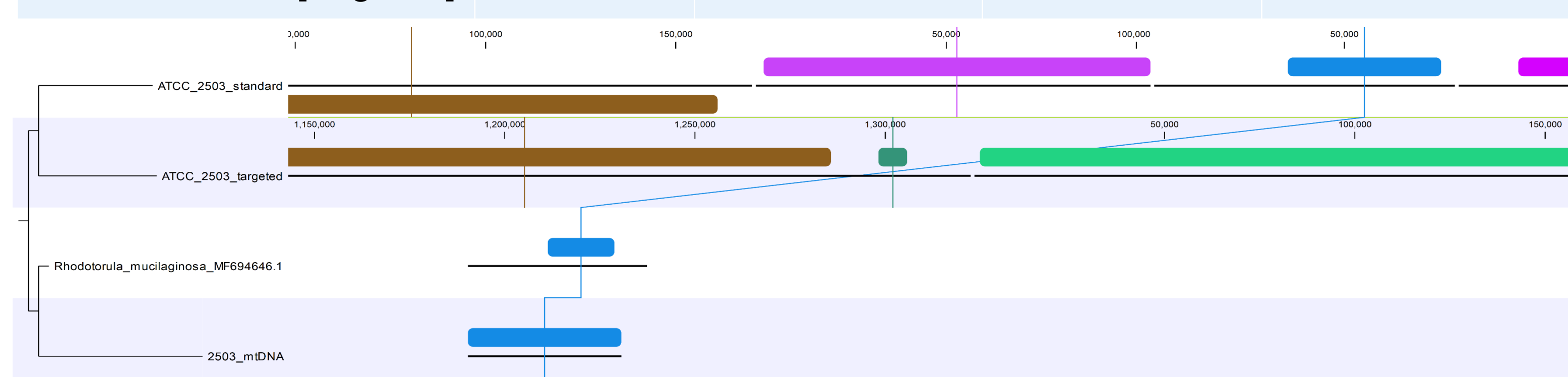


Figure 2: CLC whole-genome alignment of *Rhodotorula mucilaginosa* (ATCC_2503) assemblies to the RefSeq mitogenome (MF69464.1; 47,023 bp). Homology is color-coded. The targeted assembly excludes the mitochondrial contig (blue), and the de novo mtDNA assembly (bottom track; 40,289 bp) shares 99.74% ANI with the reference.

Results: Strain Comparisons

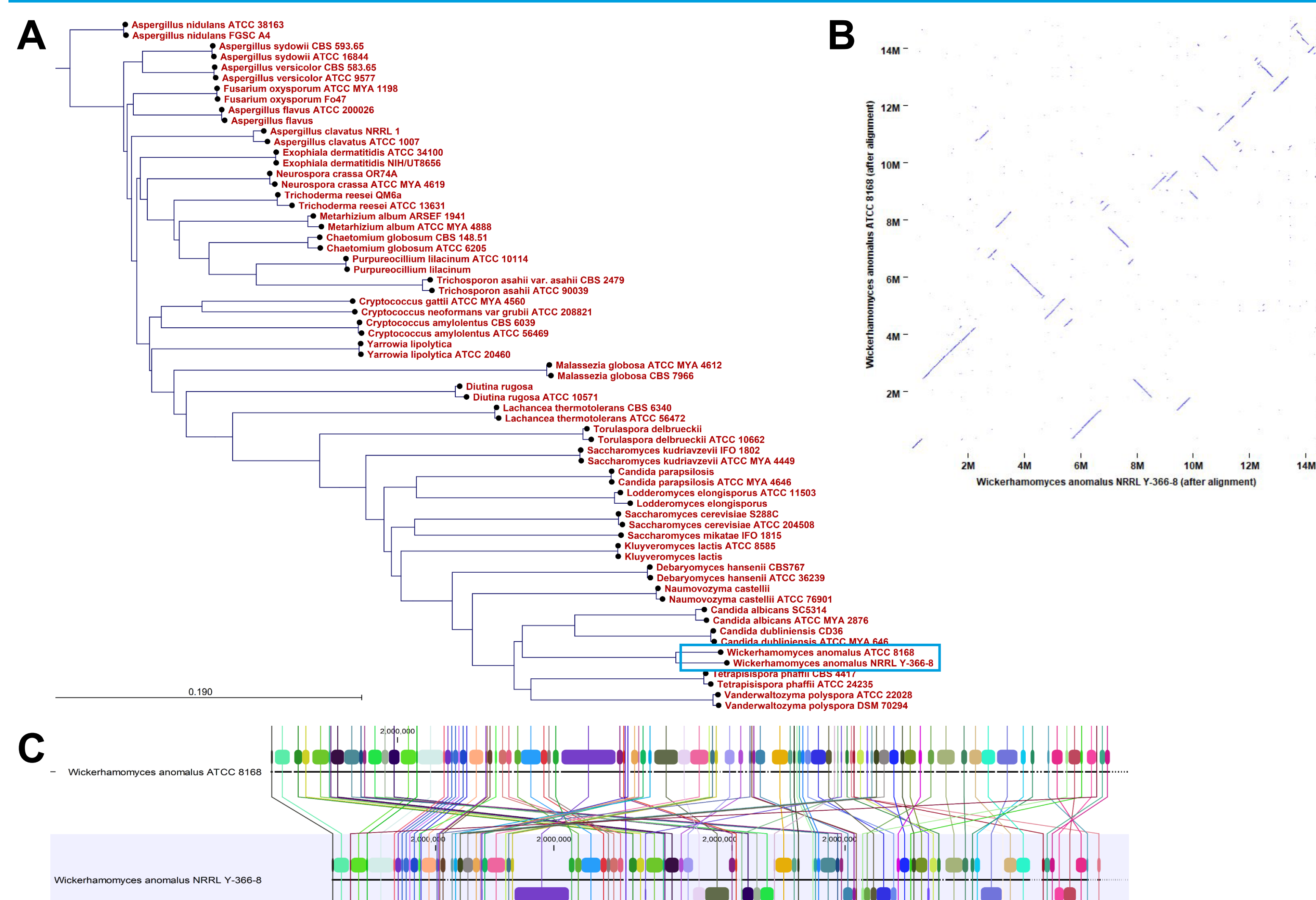


Figure 3: Comparison of strain-matched AGP and RefSeq fungal assemblies. (A) K-mer tree showing close clustering for most pairs; outliers suggest contamination, misassembly, or quality differences. (B) CLC Dot plot and (C) whole-genome alignment for *W. anomalus* (blue box in [A]) show structural differences despite >97% ANI and ~91% gene-level identity.

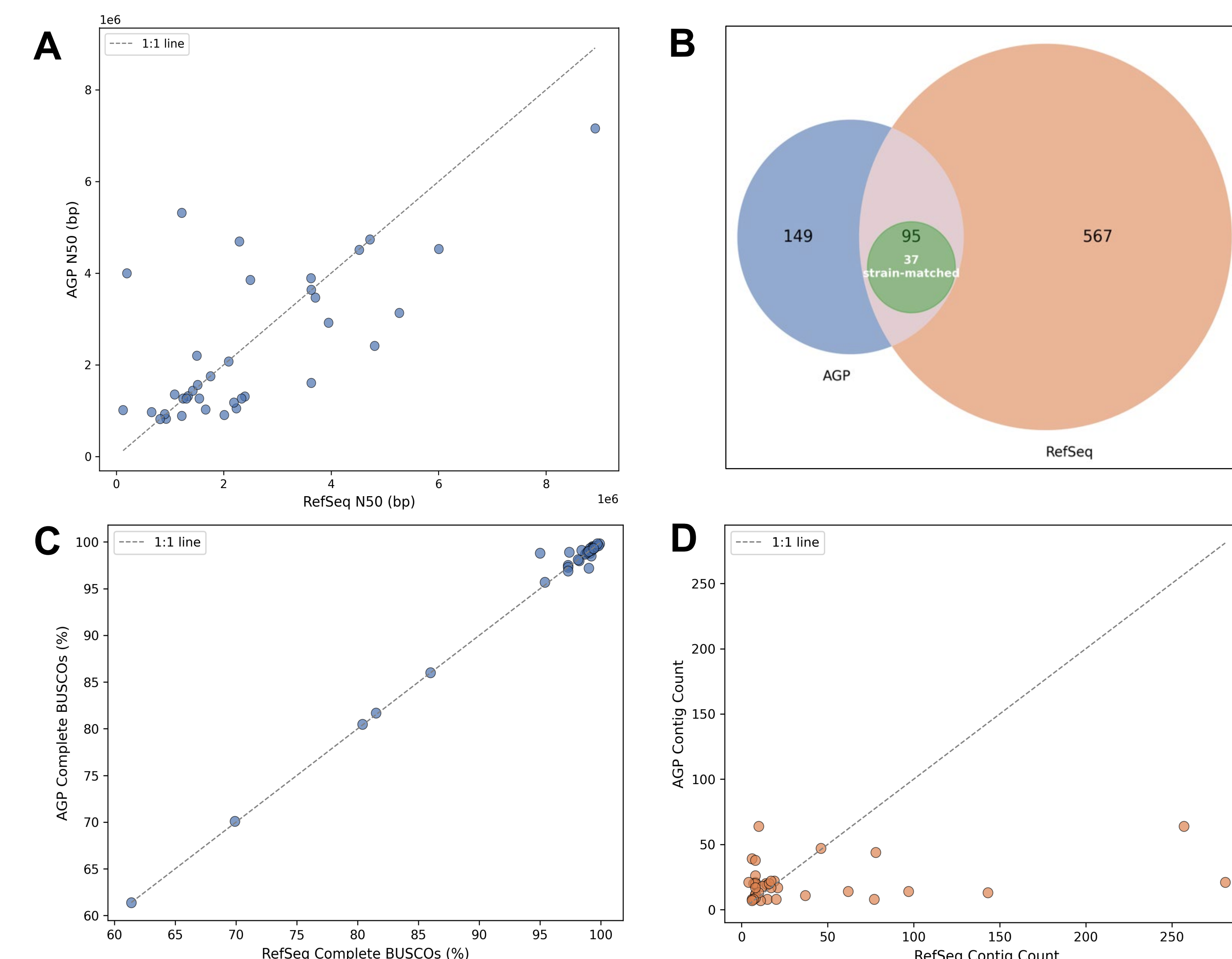


Figure 4: Comparison of AGP and strain-matched RefSeq fungal assembly (n = 37) quality metrics. (A) N50 values. (B) Species representation Venn diagram. (C) BUSCO completeness (%; eukaryota_odb12). (D) Contig count.

Conclusions

- ATCC® assemblies exhibit reduced fragmentation relative to strain-matched RefSeq genomes while maintaining comparable BUSCO completeness.
- Read-recruitment against taxa-specific mitochondrial references, followed by binned reassembly, further improves nuclear genome quality by removing embedded mitochondrial sequences.
- These findings support continued refinement of fungal assembly workflows via provenance-aware, reference-guided approaches.



Learn more about the ATCC® Genome Portal

References

1. Ahrendt SR, et al. Comparative mitogenomics of kingdom Fungi - evolutionary insights and metagenomic applications. *Nucleic Acids Res* 54(2): gkaf1419, 2026. PubMed: 41533582