

The ATCC Genome Portal: The Genomics Database of a 100-year-old Culture Collection



Scott V Nguyen, PhD; Ana Fernandes, BS; John Bagnoli, BS; Briana Benton, BS; Jonathan L Jacobs, PhD | ATCC, Manassas, VA 20110

Introduction

The ATCC® Genome Portal (AGP, <https://genomes.atcc.org>) is a reference database of authenticated and traceable whole-genome sequence assemblies generated directly from strain vials from one of the largest and most diverse culture collections in the world. As a non-profit biorepository, thousands of scientists and institutions have deposited numerous microbes, fungi, protists, viruses, cell lines, plasmids, and patented strains into ATCC over the past century.

Purpose:
The AGP provides a resource to the scientific community as many genomes that are in public genomic databases may represent sequences from strains that have been passed between laboratories, are mislabeled, or have accumulated genetic drift arising from laboratory domestication. Thousands of foodborne pathogens and food related microbes have been deposited into the collection, with isolates representing over a century of research. Currently, there are over 5,750 reference quality genomes of prokaryotes, fungi, viruses, and protists available in the AGP with new additional assemblies uploaded each quarter.

Methods

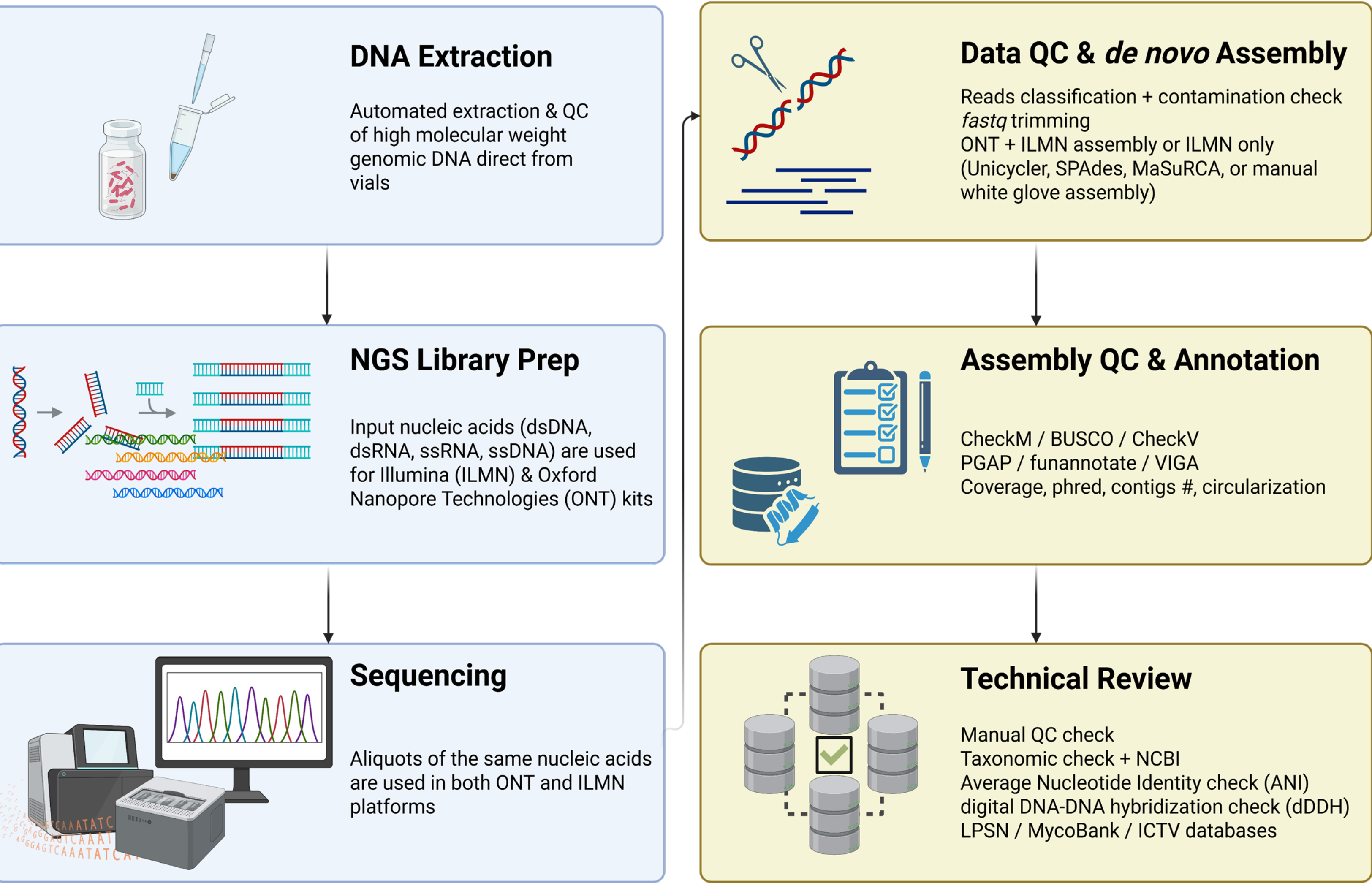


Figure 1: Standard pipeline for data provenance and sequencing of strains at ATCC. Created with BioRender.com.

Results

An Authenticated Reference Database for a Non-profit Culture Collection

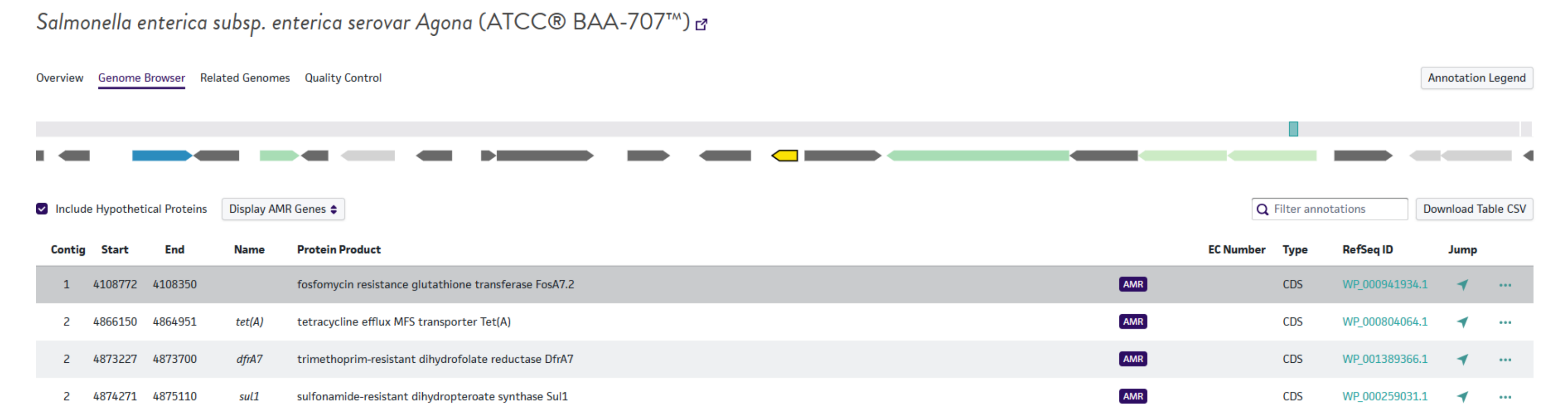


Figure 2: The AGP provides an easy-to-use Genome Browser for strains in the database.

What Reference Sequence To Use?

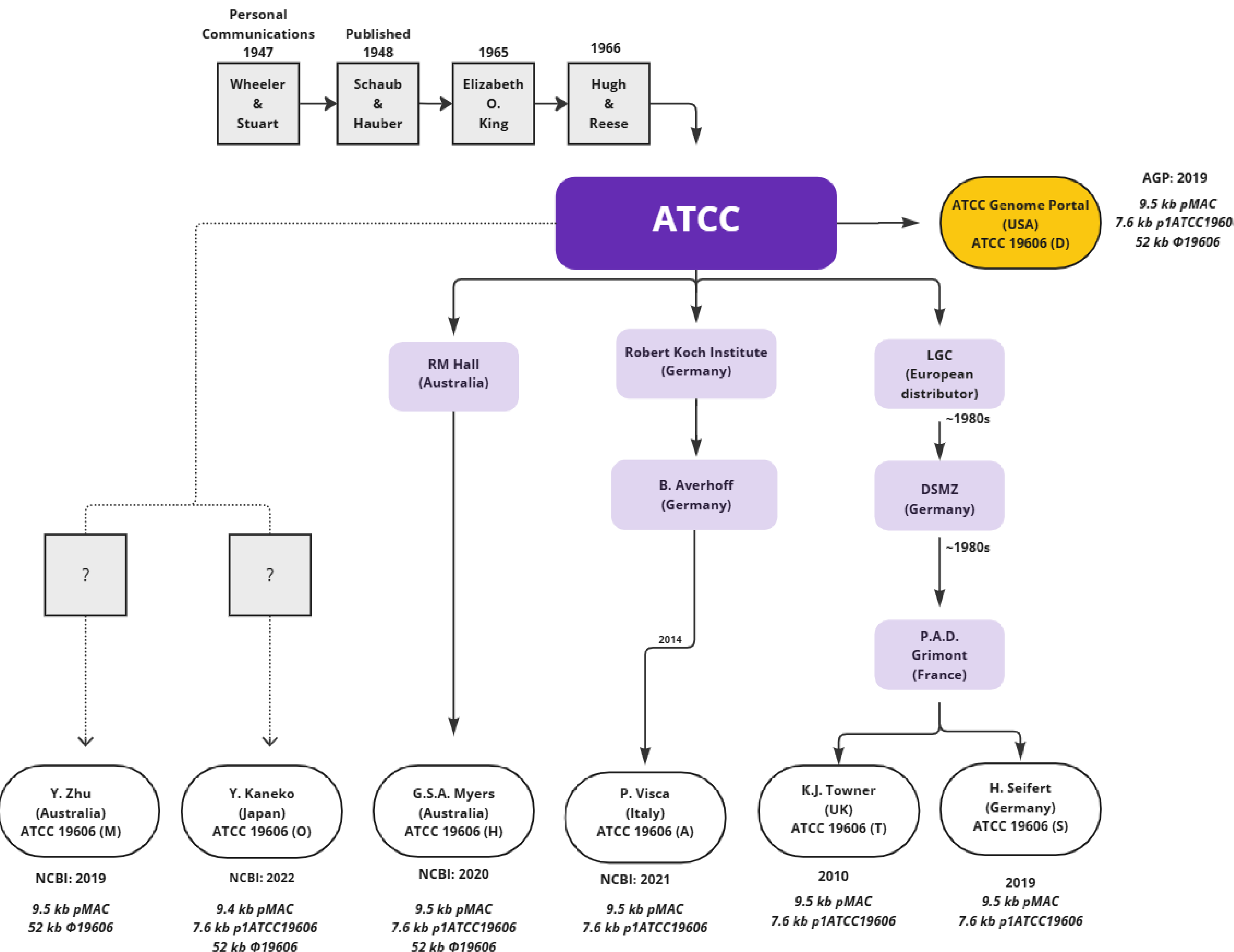


Figure 3: Multiple assemblies of *Acinetobacter baumannii* strain 2208.

Artuso *et al.* 2022 investigated the genomes assemblies of the type strain of *Acinetobacter baumannii* in NCBI (ATCC 19606). RefSeq genomes varied by thousands of SNPs and indels with missing plasmids and prophages.

Comparisons of AGP vs NCBI RefSeq Assemblies

Yarmosh *et al.* 2022 investigated direct comparisons between ATCC genome assemblies that are in NCBI RefSeq and the AGP. Over 98% of AGP assemblies were more complete and of higher quality than RefSeq representatives.

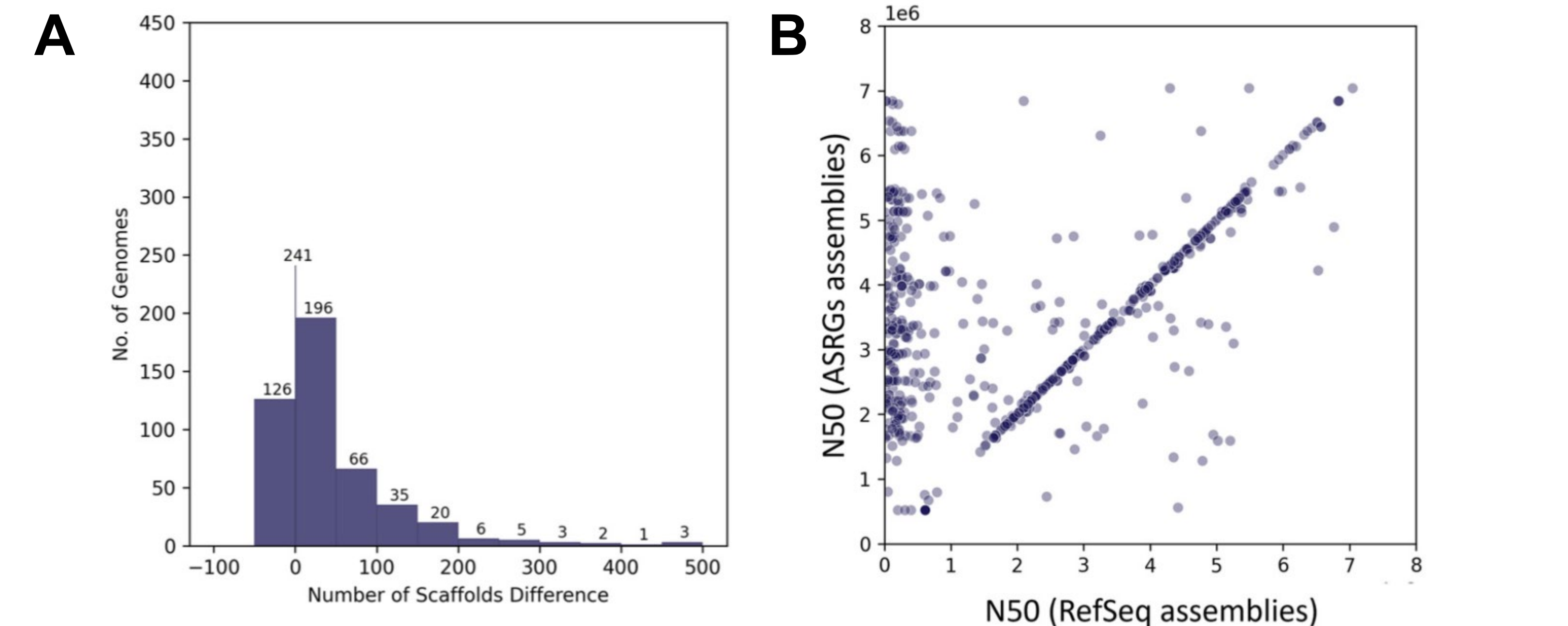


Figure 4: Comparison of AGP and NCBI RefSeq Assemblies. (A) NCBI RefSeq assemblies had more scaffolds when compared to ATCC Standard Reference Genomes (ASRGs). (B) Most ASRGs have better N50s when compared to RefSeq counterparts.

Conclusions

A majority of assemblies in AGP are of higher quality than equivalent assemblies in public genomic databases. Additionally, some of these assemblies represent strains and/or species that have no representation in public genomic databases. The AGP is a unique and authenticated resource for reference genomes.

References

- Benton B, et al. The ATCC Genome Portal: Microbial Genome Reference Standards with Data Provenance. Microbiol Resour Anounc 10(47): e0081821, 2021. PubMed: 34817215
- Yarmosh et al. Comparative Analysis and Data Provenance for 1,113 Bacterial Genome Assemblies. mSphere 7(3): e0007722, 2022. PubMed: 35491842
- Nguyen et al. The ATCC genome portal: 3,938 authenticated microbial reference genomes. Microbiol Resour Anounc 13(2): e0104523, 2024. PubMed: 38289057
- Nguyen et al. Reclassification of atypical *Moraxella catarrhalis* ATCC 23246 as *Moraxella veridica* sp. nov. Int J Syst Evol Microbiol 75(5): ePub. PubMed: 40397495
- Artuso I, Lucidi M, Visaggio D, Capecchi G, Lugli GA, Ventura M, Visca P. Genome diversity of domesticated *Acinetobacter baumannii* ATCC 19606T strains. Microb Genom. 2022 Jan;8(1):000749. PubMed: 8914354.

Learn more about the ATCC® Genome Portal



Visit and explore the ATCC® Genome Portal

