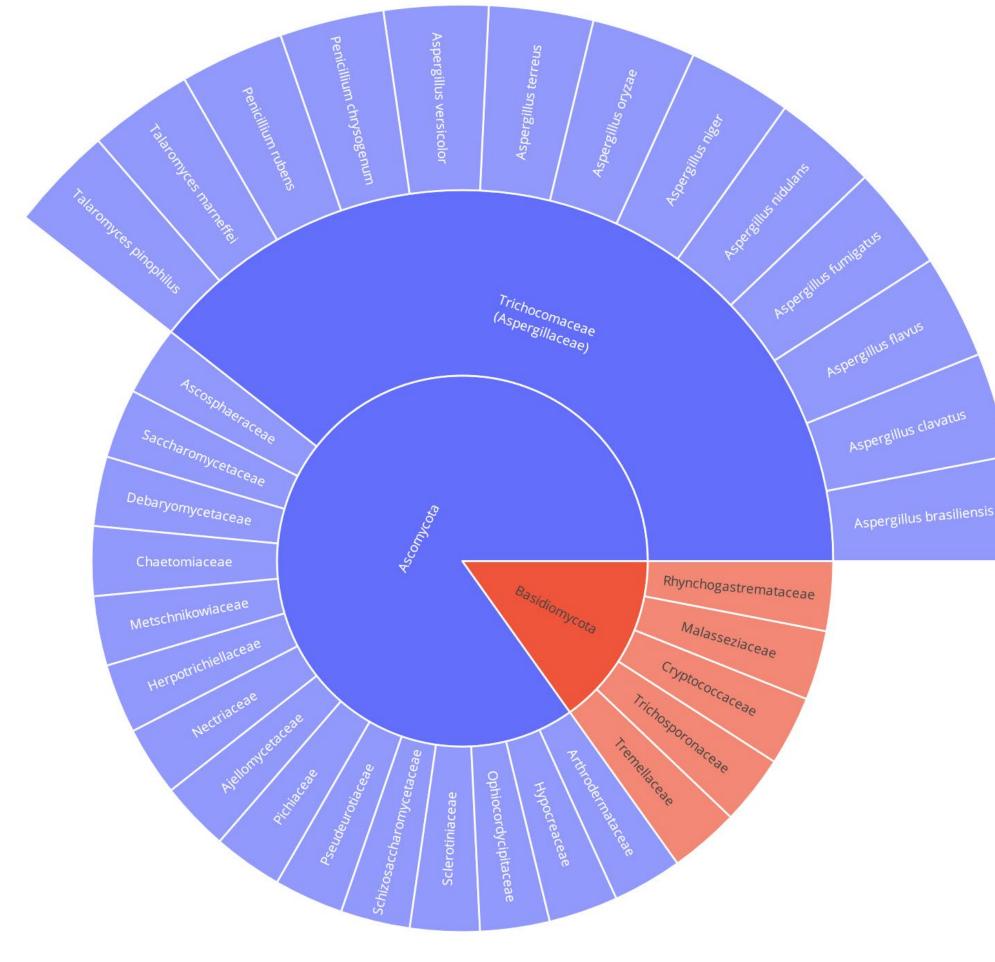
High-quality genome assemblies and biosynthetic gene clusters annotation from laboratory reference fungal strains

Ford Combs, MS; Juan Lopera, PhD; Nikhita Puthuveetil, MS; David Yarmosh, MS; Stephen King, MS; Marco Riojas, PhD; Amanda Pierola, BS; John Bagnoli, BS; Briana Benton, BS; Jonathan Jacobs, PhD ATCC, Manassas, VA 20110

Background

Fungi produce a wide variety of secondary metabolites that play an important role in human health and disease and have great potential as drug candidates and therapeutics. A comparative analysis of public assemblies and ATCC's internally produced assemblies showed that ATCC's assemblies typically had higher quality. Further analysis using genome annotation and mining tools indicated high levels of biosynthetic potential in the ATCC mycology collection.



ATCC

Figure 1: Taxonomic diversity of fungi on the ATCC Genome Portal. The innermost partition includes two phyla: Ascomycota and Basidiomycota; the middle partition includes each of the families of the phyla found in the collection; and the outermost partition displays each of the species of the Trichocomaceae family found in the collection. While many biosynthetic gene clusters (BGCs) were found in the mycology collection, a particularly large and diverse set was found in *Trichocomaceae*.

ATCC Enhanced Authentication Initiative: Mycology Collection

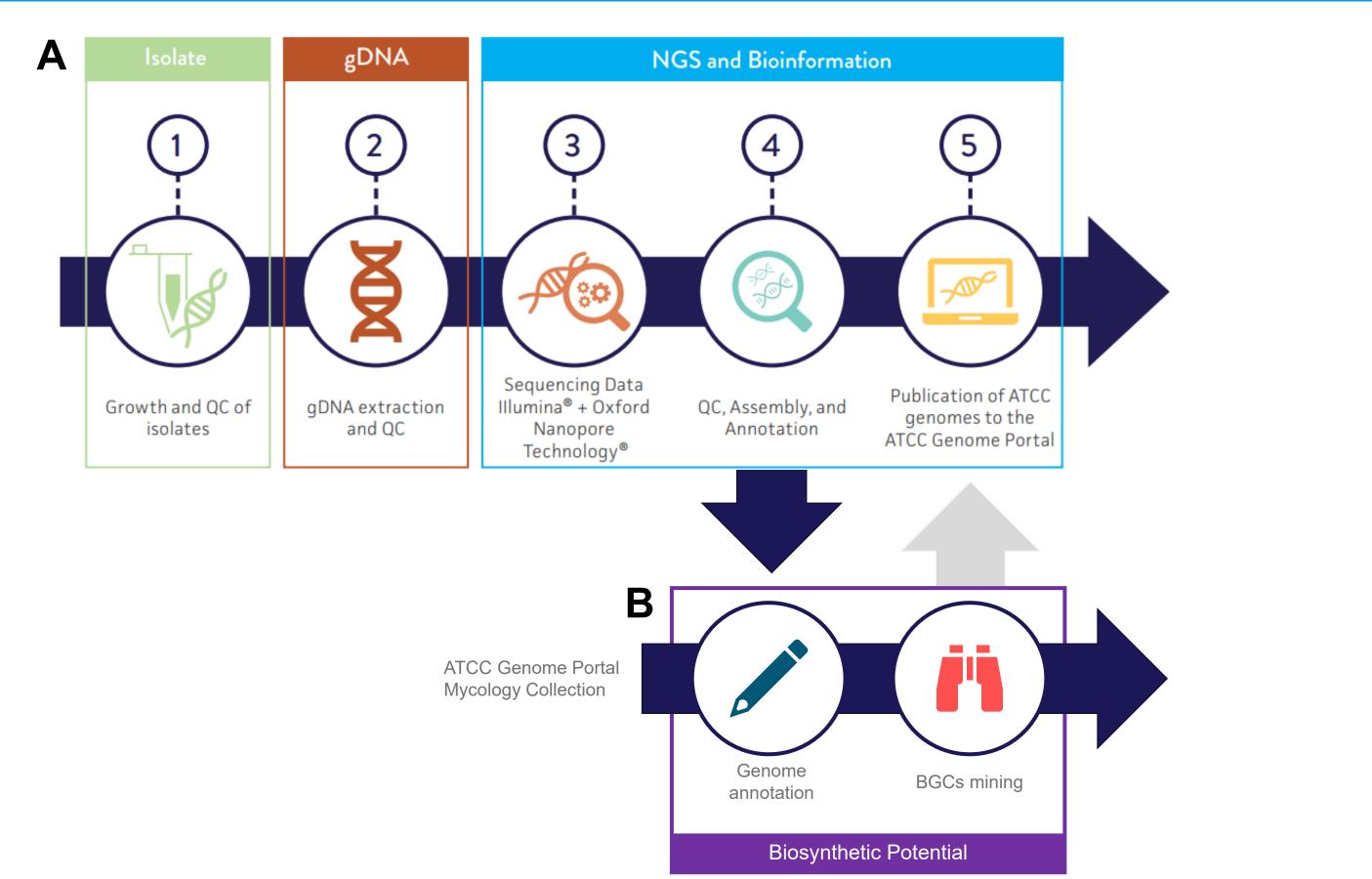


Figure 2: Workflows related to this study. (A) The ATCC genome portal workflow used to produce high-quality, authenticated genomes. (B) The biosynthetic potential workflow used in this study. ATCC Mycology Collection genomes were annotated using BRAKER and then mined using antiSMASH to detect BGCs.

10801 University Boulevard, Manassas, Virginia 20110-2209

© 2021 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks owned by the American Type Culture Collection unless indicated otherwise. Illumina is a registered trademark of Illumina, Inc. Oxford Nanopore is a registered trademark of Oxford Nanopore Technologies Limited.

Comparison of ATCC and Public Assembly Quality

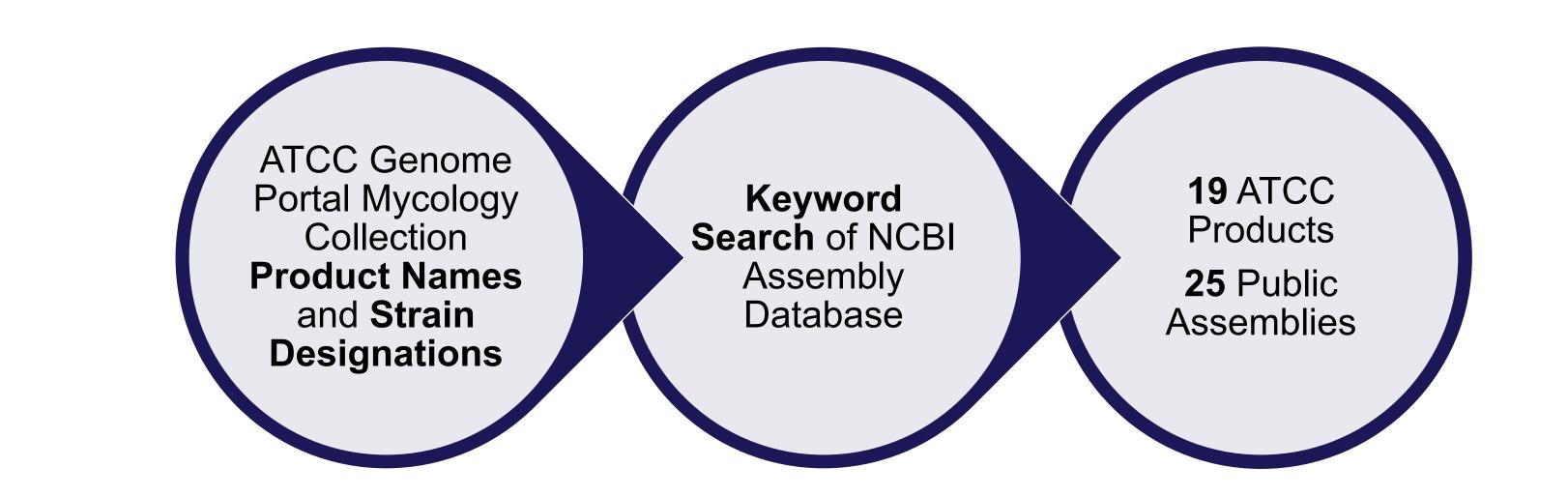


Figure 3: Process description for finding public assemblies associated with ATCC mycology collection products. The product name and strain designations for each ATCC mycology collection product were used as keywords in a search of the NCBI assembly database.

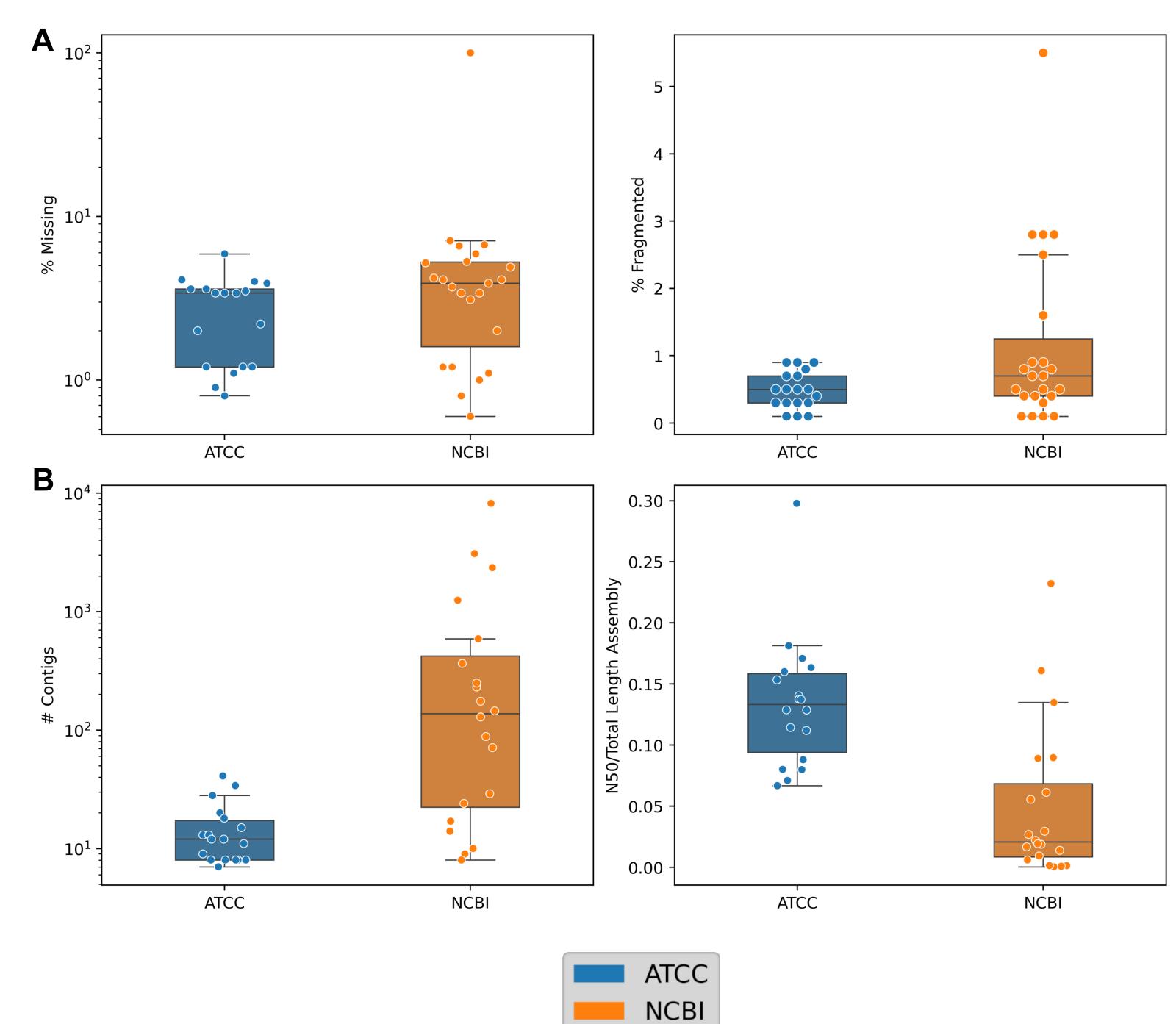


Figure 4: Comparison results of ATCC Genome Portal assemblies and their associated public assemblies. (A) Scatterplots showing BUSCO percent missing and percent fragmented of each assembly. (B) Scatterplots showing contig count and N50/Total Length of Assembly of each assembly.

Phone: 800.638.6597

Comparison of ATCC and NCBI Assembly Quality

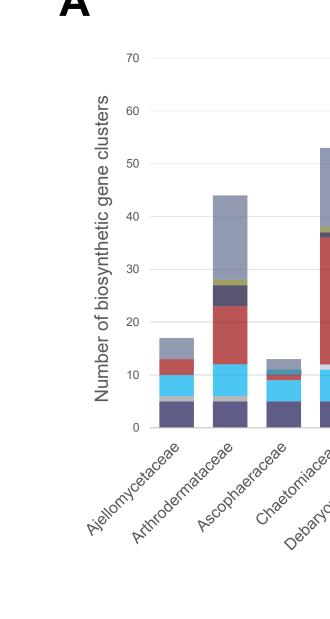


Figure 5: Charts depicting the total number of BGCs of different types found in the fungal families from the ATCC Genome Portal Mycology Collection. (A) All fungal families except for Trichocomaceae and (B) Trichocomaceae only, separated because of its significantly greater quantity of BGCs.

Aspergillus niger ATCC 1015

- Aspergillus_nidulans_ATCC_38163

Aspergillus_flavus_ATCC_204304

- Aspergillus oryzae ATCC 42149

Aspergillus terreus ATCC 20542

- Aspergillus_clavatus_ATCC 1007

Penicillium rubens ATCC 28089

- Aspergillus_flavus_ATCC_9643

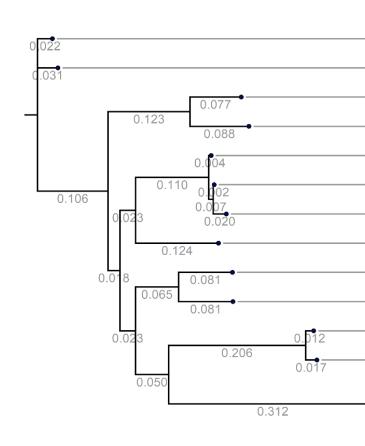


Figure 6: Phylogenetic tree (ML) of *Trichocomaceae* family and associated number and types of BGCs found in each species.

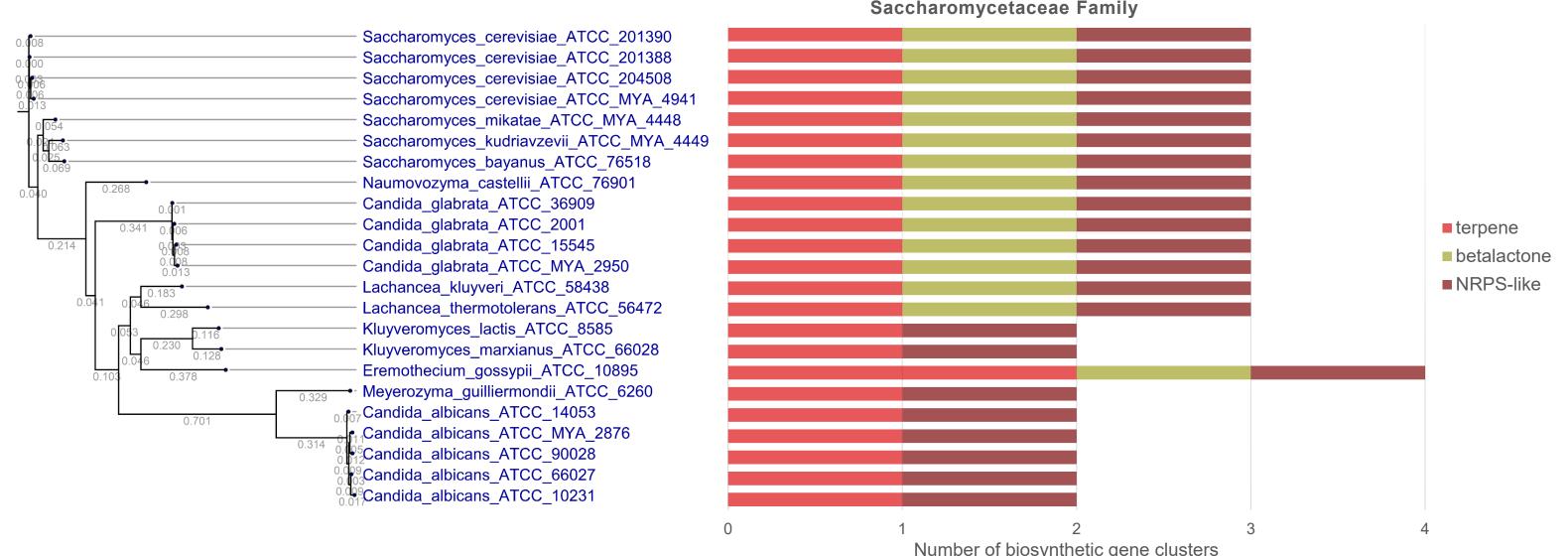


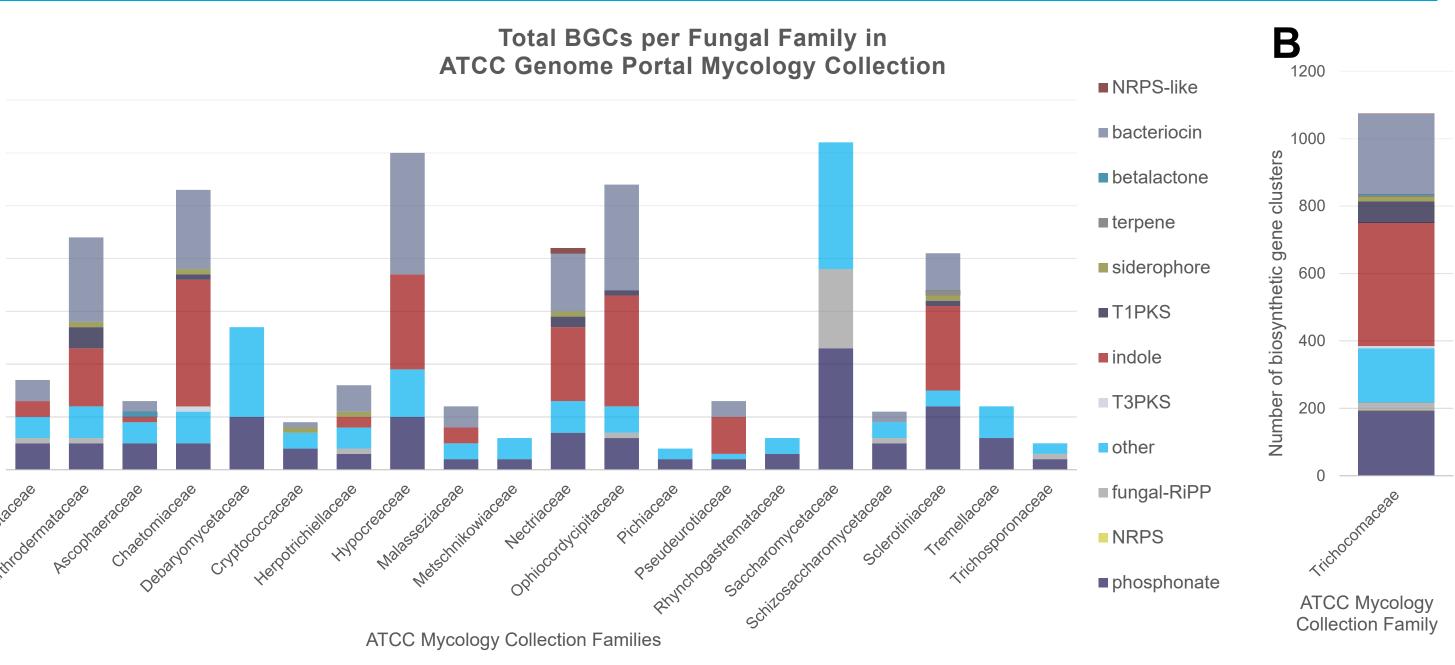
Figure 7: Phylogenetic tree (ML) of Saccharomycetaceae family and associated number and types of BGCs found in each species.

Conclusions

The ATCC Genome Portal mycology collection contains dozens of high-quality, authenticated assemblies that are linked to physical materials. Annotation and genome mining revealed a large presence of BGCs in these genomes, particularly in the *Trichocomaceae* family.



BGC Mining and Phylogenetic Analysis



BGCs found in ATCC's Genome Portal Mycology Collection Trichocomaceae Family

BGCs found in ATCC's Genome Portal Mycology Collection

