



Phenotypic and Genotypic Characterization of Antimicrobial Resistance (AMR) Strains from the ATCC® Collection

ATCC, Manassas, VA 20110

Background

or other applications

to establish its phenotypic and genotypic profile.

- **Methylation profiles:** Most strains are provided with methylation data accessible through the ATCC® Genome

- Portal (genomes.atcc.org).

Methods

Susceptibility Testing

susceptibility was obtained using bioMérieux ETEST strips and MIC interpretations are based on CLSI M100-Ed35.

Genomic Analyses

available on the ATCC® Genome Portal.

Methylation Profiles

Methylation profiles were generated with ONT data using the Dorado basecaller version 0.8.0+accec121 (ONT). The methylation profile for many strains are available to customers through the ATCC® Genome Portal.

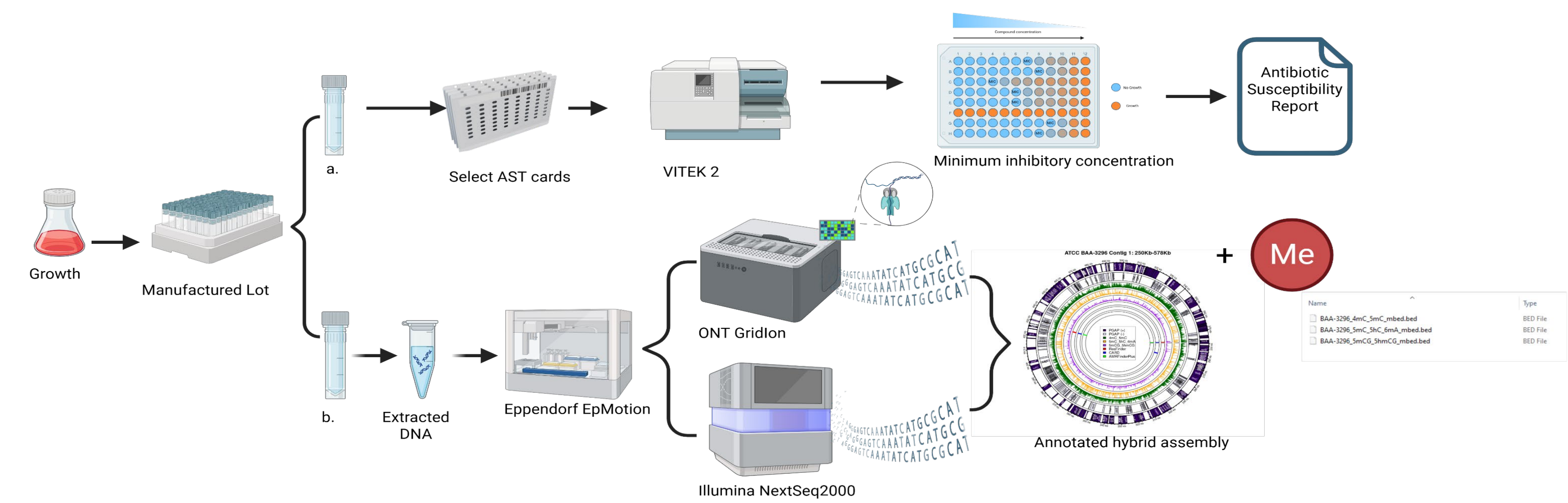


Figure 1: Bacterial strains are grown, viald, and QC'd according to ATCC® parameters. (A) The appropriate VITEK 2 AST cards (bioMérieux) are selected and loaded on to the VITEK 2 instrument to produce the MIC profile. Note: for *Neisseria gonorrhoeae*, antibiotic susceptibility was obtained using ETEST strips (bioMérieux). (B) DNA is extracted using a method best suited for organism type and next-generation sequencing (NGS) libraries are produced and loaded on the NextSeq 2000 (Illumina) and Gridlon (Oxford Nanopore Technologies) instruments. Data from both platforms were QC'd and a subsequent hybrid genome assembly was produced and annotated. ONT base called data was further analyzed for the presence of methylated nucleotides. Image created using BioRender.com

Results

Table 1: Species diversity of antimicrobial-resistant strains

Species	Strains	VITEK 2 Card(s) or ETEST strips	Gram stain
<i>Acinetobacter baumannii</i>	13	AST-XN09 and AST-GN69	Negative
<i>Citrobacter braakii</i>	1	AST-XN09 and AST-GN69	Negative
<i>Citrobacter freundii</i>	4	AST-XN09 and AST-GN69	Negative
<i>Enterobacter asburiae</i>	1	AST-XN09 and AST-GN69	Negative
<i>Enterobacter cloacae</i>	3	AST-XN09 and AST-GN69	Negative
<i>Enterobacter hormaechei</i>	2	AST-XN09 and AST-GN69	Negative
<i>Escherichia coli</i>	17	AST-XN09 and AST-GN69	Negative
<i>Klebsiella oxytoca</i>	1	AST-XN09 and AST-GN69	Negative
<i>Klebsiella pneumoniae</i>	14	AST-XN09 and AST-GN69	Negative
<i>Klebsiella quasipneumoniae</i>	2	AST-XN09 and AST-GN69	Negative
<i>Neisseria gonorrhoeae</i>	5	BioMerieux Etest strips	Negative
<i>Proteus mirabilis</i>	3	AST-XN09 and AST-GN69	Negative
<i>Pseudomonas aeruginosa</i>	15	AST-GN69 and AST-XN06	Negative
<i>Pseudomonas paraaeruginosa</i>	1	AST-GN69 and AST-XN06	Negative
<i>Serratia surfactantifaciens</i>	1	AST-GN69 and AST-XN06	Negative
<i>Staphylococcus aureus</i>	2	AST-GP72 and AST-GP75	Positive
<i>Streptococcus pneumoniae</i>	15	AST-ST02	Positive

Enterobacter hormcheae subsp. *xiangfangensis*

(Gu et al.) Sutton et al.

[Download genome](#)

LEARN ABOUT THE ATCC GENOME PORTAL >

This drug-resistant bacterium was isolated in 2012 from the urine of a 57-year-old female in Turkey. This product can be used for all stages of the discovery and development process for novel antimicrobial and therapeutic, molecular-based detection assays, and updated efficacy protocols.

BACTERIA

2 Product Classes

Product category	Bacteria
Product type	Drug-resistant bacterium
Classification	Enterobacteriaceae
Strain designation	E8960
Type strain	No
Genome deposited strain	Yes
Former name	Enterobacter cloacae (Clostrid) Hormachee and Edwards
Isolation source	Urine
Geographical isolation	Turkey
Applications	Antibiotic-resistance research Bioinformatics Drug development
Product format	Frozen
Storage conditions	-80°C or colder

Documentation

- Product sheet
- Certificate of analysis
- Safety data sheet

Antimicrobial resistance report

Figure 2: Example of product page with metadata and available documentation.

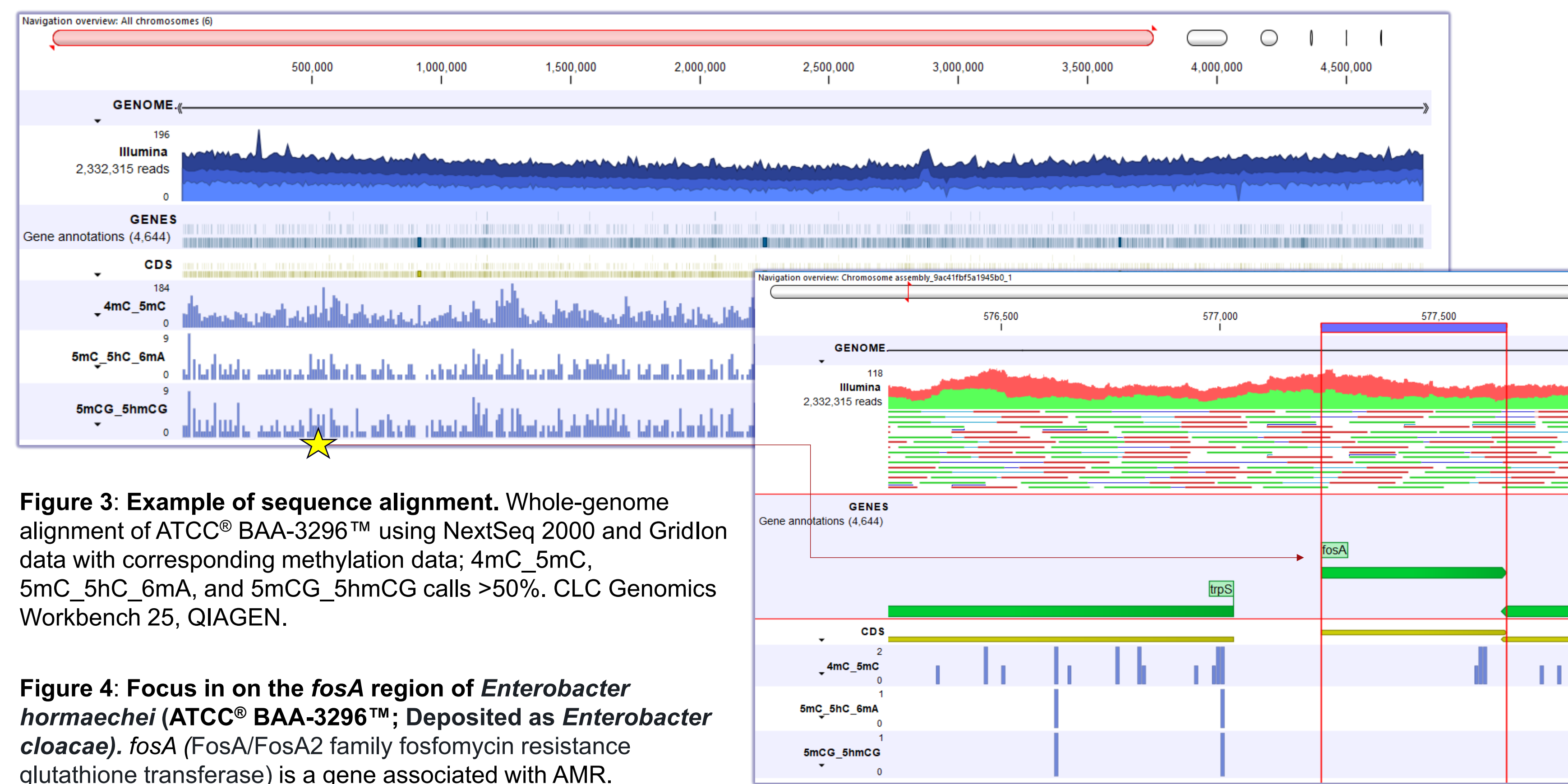


Figure 3: Example of sequence alignment. Whole-genome alignment of ATCC® BAA-3296™ using NextSeq 2000 and Gridlo data with corresponding methylation data; 4mC_5mC, 5mC_5hC_6mA, and 5mCG_5hmCG calls >50%. CLC Genomics Workbench 25, QIAGEN.

Figure 4: Focus in on the *fosA* region of *Enterobacter hormaechei* (ATCC® BAA-3296™; Deposited as *Enterobacter cloacae*). *fosA* (FosA/FosA2 family fosfomycin resistance glutathione transferase) is a gene associated with AMR.

→ **Table 2:** MIC values and interpretations of 16 *Klebsiella* spp.

[illegible]

(R) resistant, (I) intermediate, and (S) susceptible. A report outlining the MIC values and interpretation of susceptibility accompanies each strain. (Supplemental handout available)

► **Table 3:** MIC values and interpretations of 15 *Streptococcus pneumoniae* strains.

[illegible]

(R) resistant, (I) intermediate, and (S) susceptible. A report outlining the MIC values and interpretation of susceptibility accompanies each strain. (Supplemental handout available)

Conclusion

Here, we highlight one hundred clinically relevant and extensively characterized antimicrobial-resistant isolates that are available to the research community. Each strain comes with:

- **Genetic data** – The complete and assembled genome sequence, annotated with antibiotic resistance genes and methylation data. Data is available on genomes.atcc.org

Overall, this diverse collection of highly characterized AMR strains provides a valuable resource for diagnostics and therapeutic development.

Explore our AMR clinical isolates at
www.atcc.org/AMR



Find the whole-genome sequences on the ATCC Genome Portal at genomes.atcc.org

