

Finding Your Perfect Match – Evolving Technologies for Bacterial Strain Typing

Brian Cantwell, Ph.D.
Scientist, ATCC



About ATCC

- Founded in 1925, ATCC is a non-profit organization with headquarters in Manassas, VA and an R&D and Services center in Gaithersburg, MD (225,000 sq. ft. total)
- World's largest and most diverse biological materials resource center
 - 5,000 cell biology products
 - 80,000 microorganisms
- ATCC collaborates with and supports the scientific community with industry-standard biological products and innovative solutions
- Talented team of 450+ employees; over one-third with advanced degrees



Established partner to global researchers and scientists

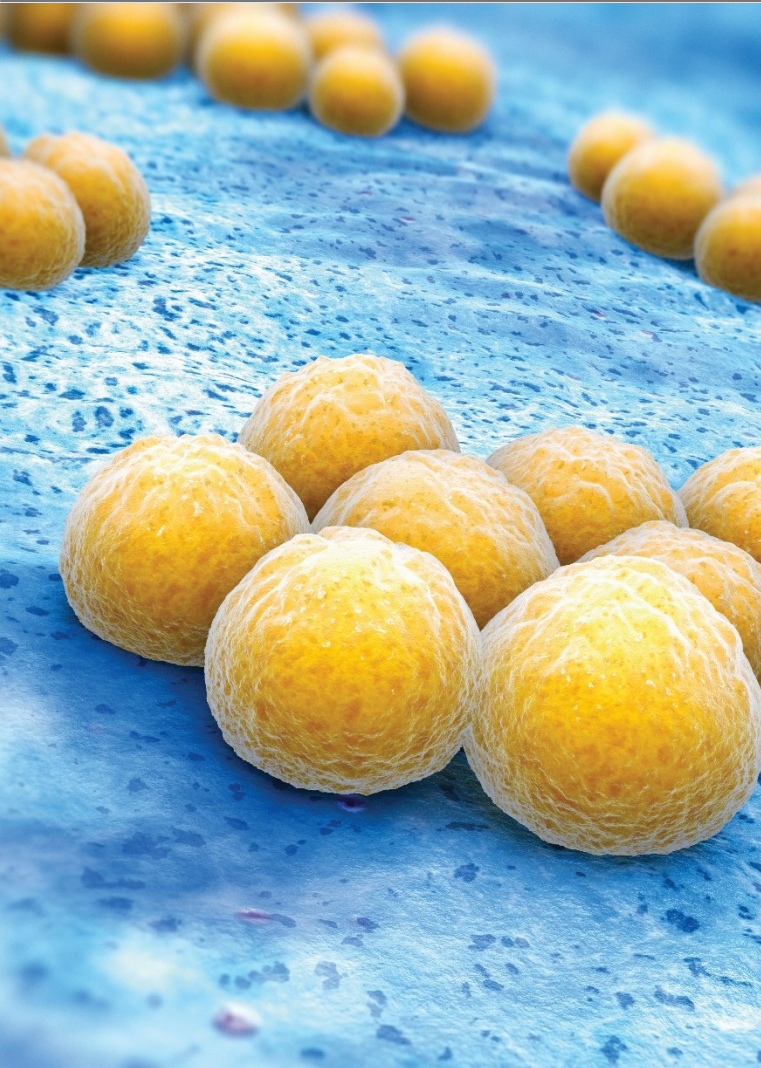


Outline



- Definition of a strain
- Why strain typing is important
- Phenotypic and genotypic methods for strain identification

Definition of a strain

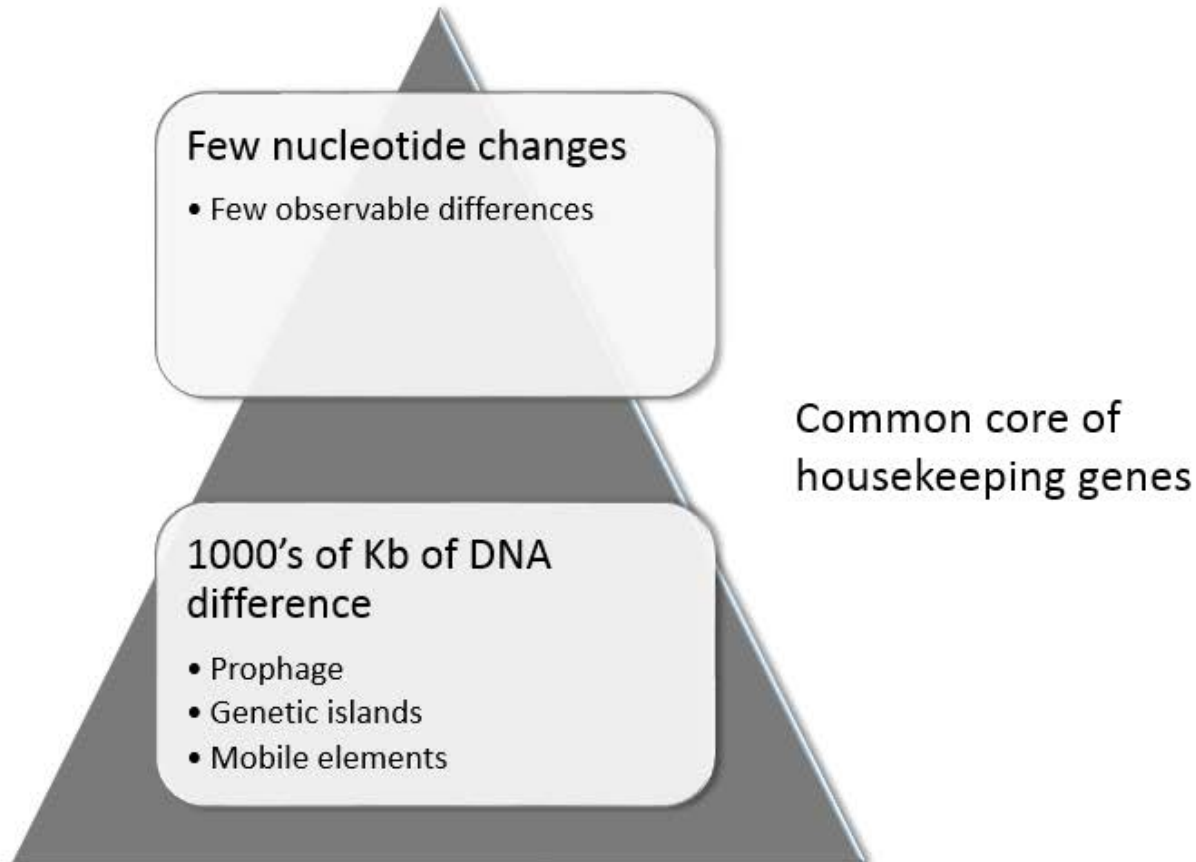


What is a “strain”?

- An individual isolate?
- A clonal population?
- A genetic variant?

A strain is a genetic variant or subtype of a microorganism below the level of species and subspecies

Strain variability



Strains of the same species can have a wide range of variation

Why is it important to identify strains?

Identify novel isolates

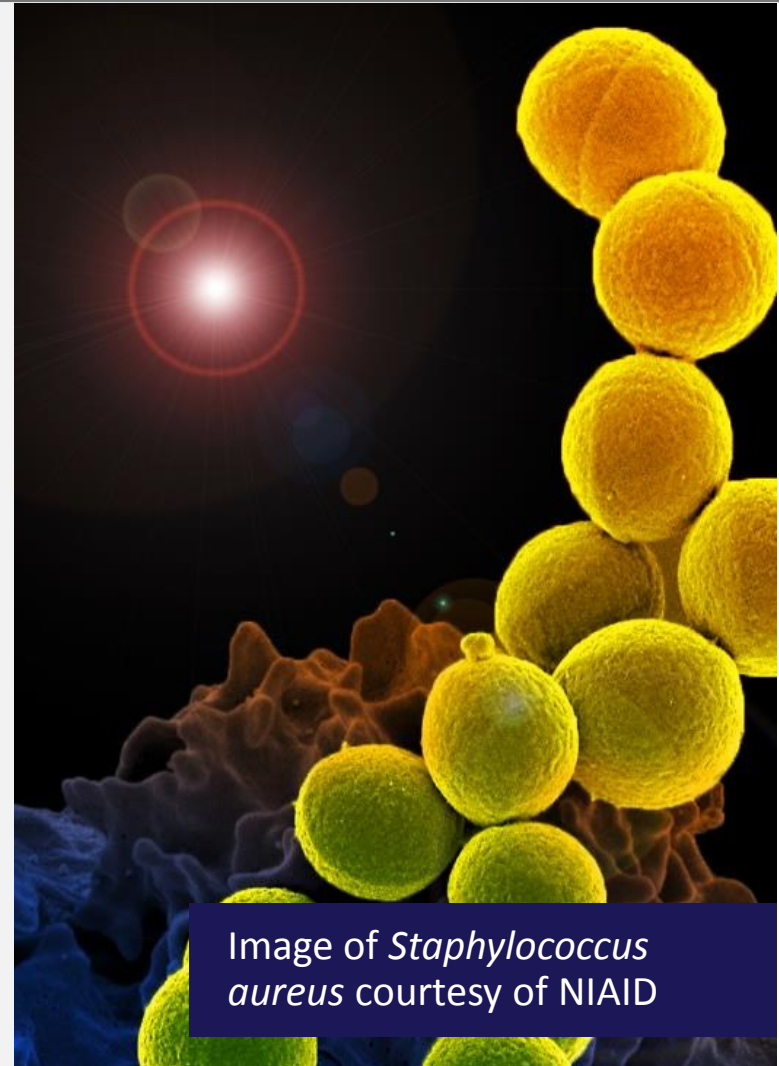
- Defining type strains and variants

Clinical impact

- Identifying antimicrobial-resistant strains and virulent strains

Epidemiology

- Tracking outbreak sources and modes of transmission



Identifying strains



Phenotypic Methods

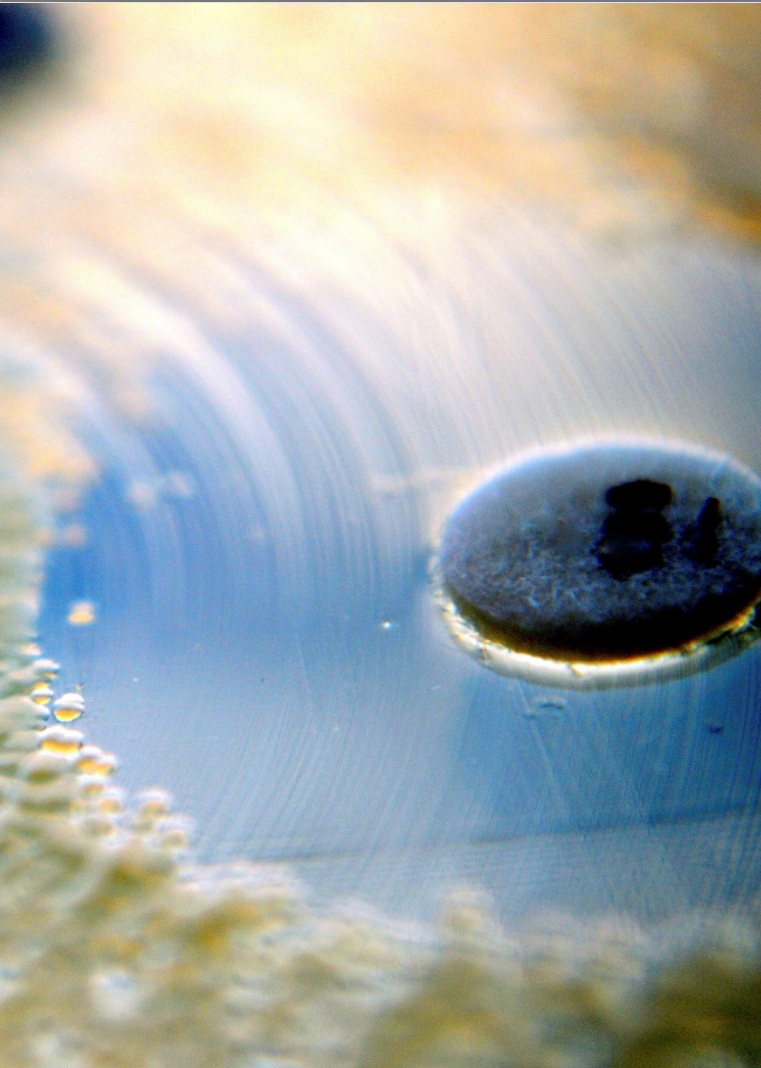
- Biotyping
- Phage typing
- Serotyping



Genotypic Methods

- Electrophoretic analysis
- PCR analysis
- Sequence analysis

Biotyping

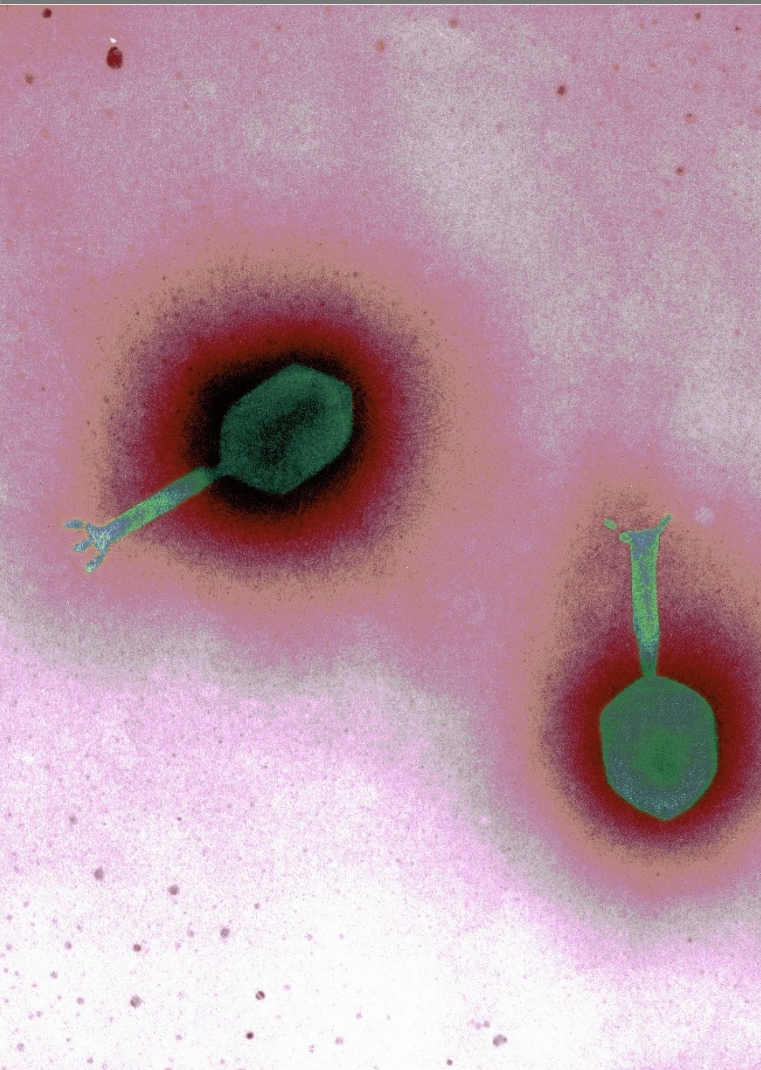


The use of differential tests, in addition to typical morphological and functional analyses used to determine the genus and species

- Growth in differential media
- Colony morphology
- Cell morphology
- Motility
- Antimicrobial resistance

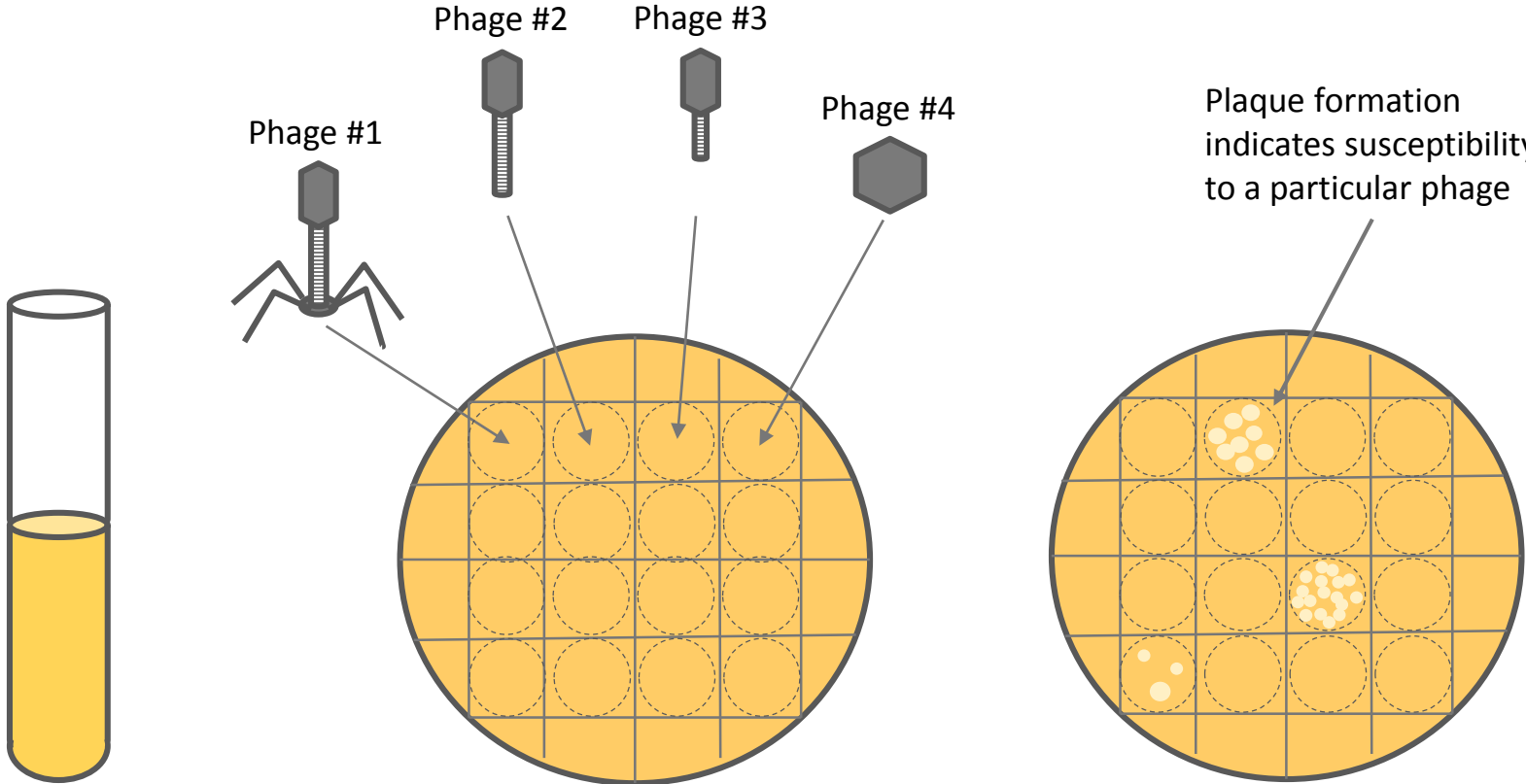
These assays have limited discriminatory power

Phage typing



- Divides species into strains on the basis of susceptibility to bacteriophage lysis
- Uses common lab equipment and techniques
- Good discriminating power, but limited by panels of available phages and susceptibility to mutation

Phage typing



Prepare a lawn of the bacterial test strain

Spot with aliquots of each phage in the panel

Evaluate the pattern of phage susceptibility to identify the strain

Serotyping



Image of *Salmonella* serotype Typhi courtesy of the CDC

- Groups strains on the basis of antigenic variation on the cell surface
- Convenient and widely used
 - Rapid and easy to perform
 - Requires the use of basic lab equipment
- Limited by the availability of antibodies against specific antigens

Serotyping

Bacterial species	Antigens
<i>Escherichia coli</i>	LPS oligosaccharide Flagellar subunit
<i>Salmonella</i> sp.	LPS oligosaccharide Flagellar subunit Capsule
<i>Pseudomonas</i> sp.	LPS oligosaccharide
<i>Streptococcus pneumoniae</i>	Capsule
<i>Haemophilus influenzae</i>	Capsule

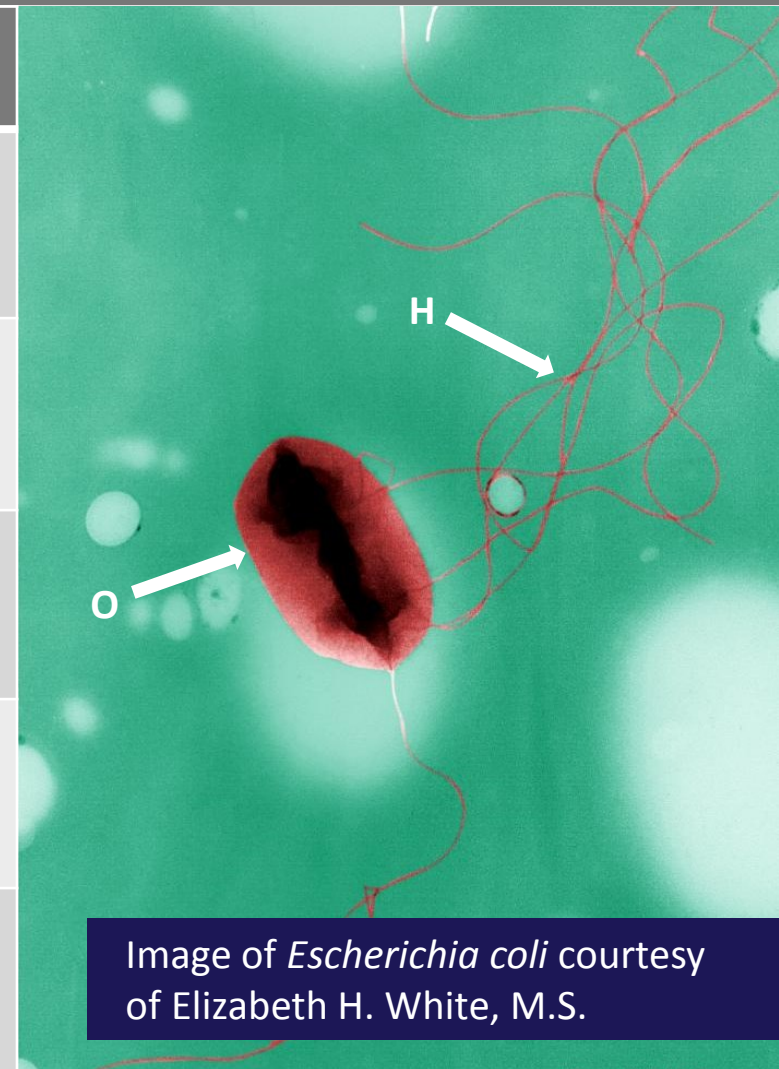


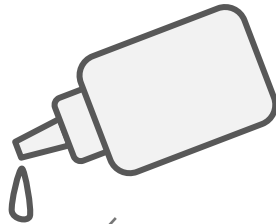
Image of *Escherichia coli* courtesy of Elizabeth H. White, M.S.

Serotyping

Culture



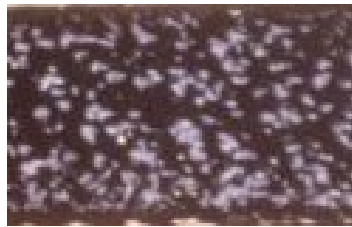
Antibody solution



Mix

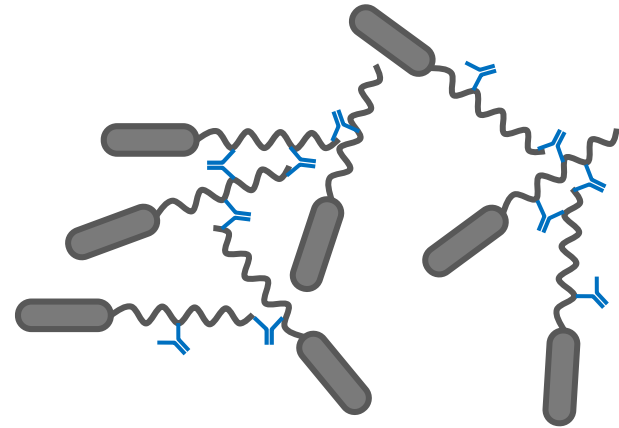


(-)

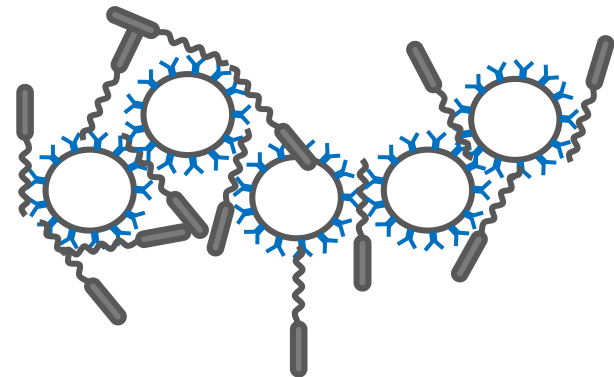


(+)

Direct agglutination



Indirect agglutination
(Latex beads)



Genotypic methods

Genome fragment size

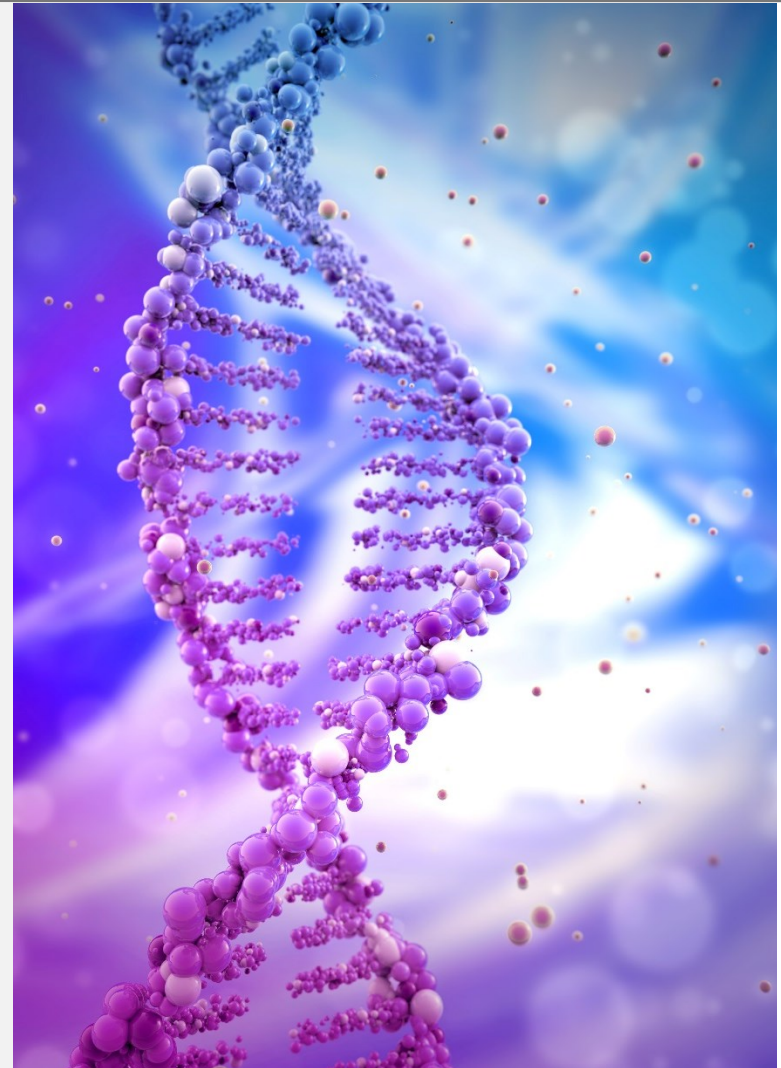
- Electrophoresis-based methods

Gene presence

- PCR-based detection of genes

Gene sequence

- Nucleotide sequence of one or more genes



Bacterial genome

Core Genome

- Set of housekeeping genes in all strains
- Evolves slowly
- Conserved positions

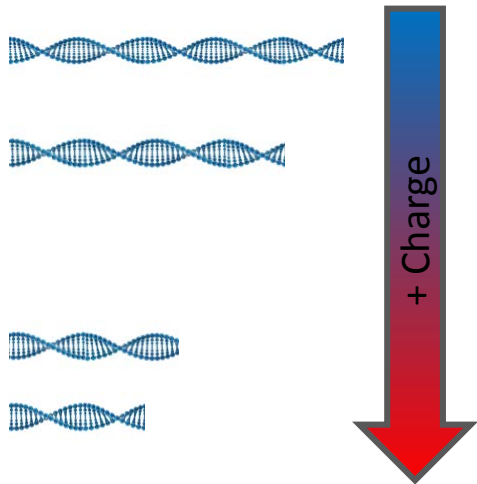
Accessory Genes

- Additional functions in a subset of strains
- Includes virulence genes, resistance genes, prophage, etc.
- Evolves quickly
- Frequent rearrangement, insertion/deletion

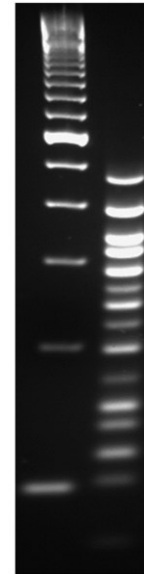
Electrophoretic analysis



Fragment genomic DNA using restriction enzymes



Separate by size using electrophoresis



Visualize fragments and compare pattern to database

Restriction fragment length polymorphism



DNA cut into small fragments

- Generates large number of fragments
- Visualized by hybridization probe

Difference in banding pattern

- Lost or gained restriction enzyme sites
- Insertions or deletions

Examples

- Ribotyping – 16S rDNA probe
- *Clostridium* toxinotyping – PCR of highly variable toxin locus

Pulsed-field gel electrophoresis

Bacterial cells imbedded in gel plugs

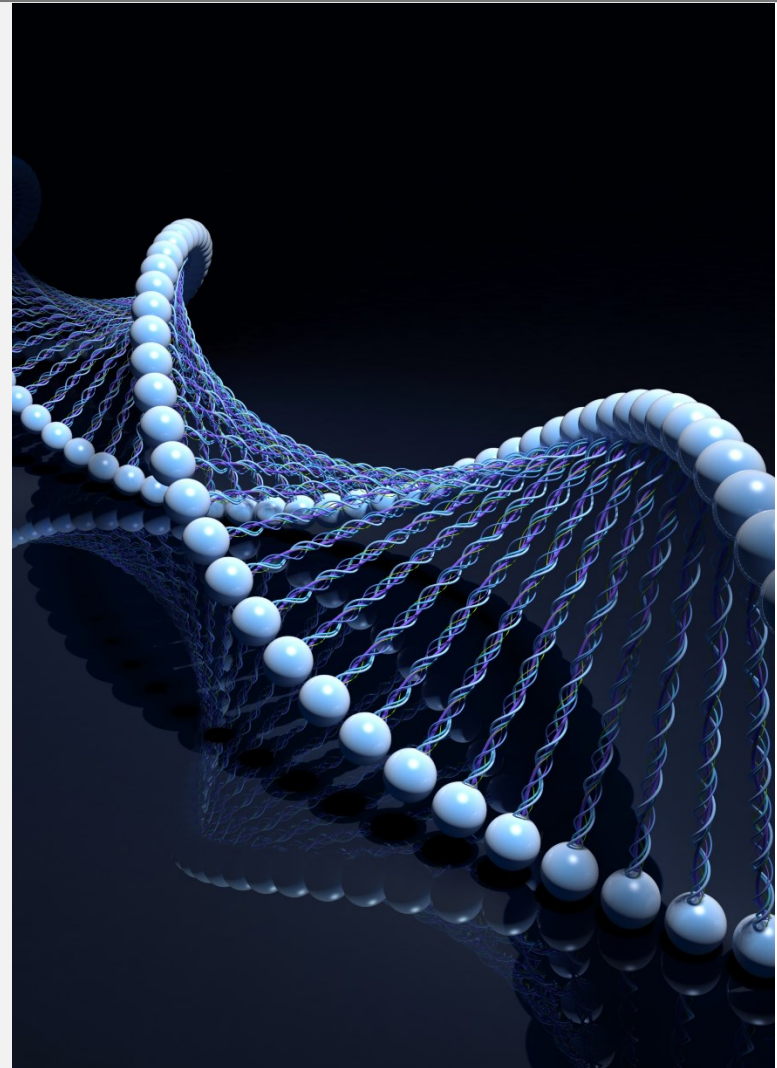
- Cells lysed and DNA digested in plugs
- Generate 10-20 large fragments

Separate DNA by using a pulsed-field gel apparatus

- Directly visualize fragments

Compare pattern to those within an established database

- Software stores and compares imaged gels



Pulsed-field gel electrophoresis databases

PulseNet

- Established by CDC in 1996
- Searchable database of PFGE patterns
- Patterns provided by local health labs
- CDC cluster analysis looks for patterns in new uploads

PulseNet International

- Success of PulseNet led to similar efforts in other regions
- Advanced networks in Europe, Canada, etc.
- Challenges in developing world include access to PFGE equipment and data

Multiple Locus Variable-number Tandem Repeat Analysis

Variable-number tandem repeats

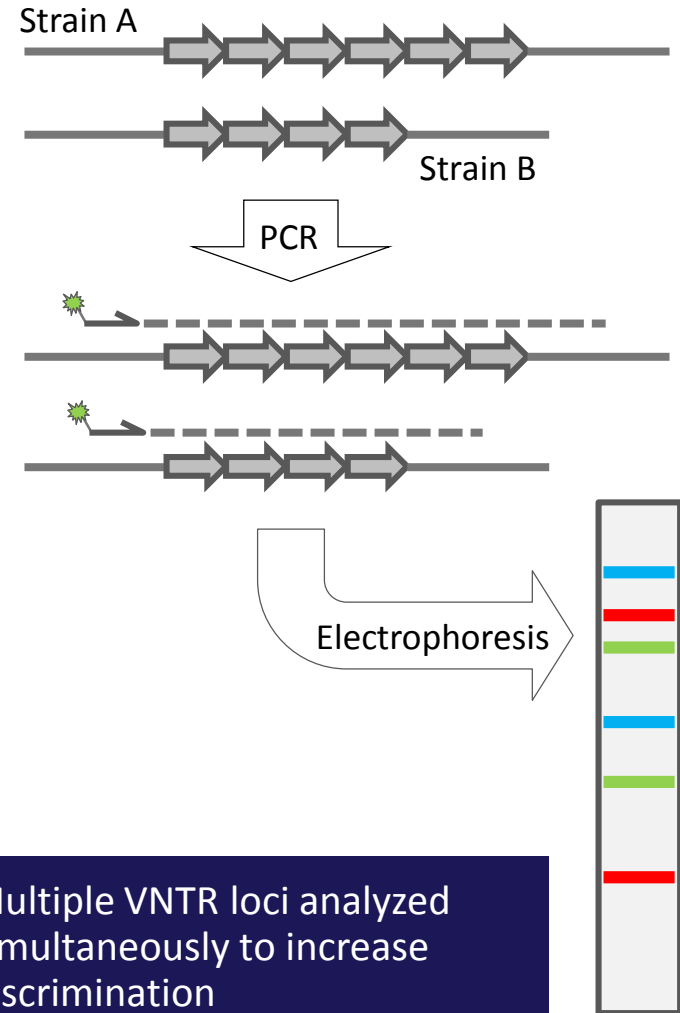
- Short repeated nucleotide sequences
- Found in most organisms
- The number of repeats varies between strains

MLVA

- Amplify tandem repeat loci by PCR and separate by electrophoresis
- Visualize using fluorescently labeled primers

MLVA resources

- PulseNet, MLVABank
- Strain databases and protocols for different MLVA equipment
- PulseNet Databases include *Salmonella* and *E. coli* O157:H7
- MLVABank has databases for ~20 species



Multiple VNTR loci analyzed simultaneously to increase discrimination

Amplified Fragment Length Polymorphism

...GATTA CCGCTATCTCA... Digest
...C TAATGCCGATAGAGT...

...NNNGATTACGGCTATCTCA... Add
...NNNCTAATGCCGATAGAGT... Adapter

...NNNGATTACGGC → PCR w/
| | | | | | | | | | matching
...NNNCTAATGCCGATAGAGT... primer

...NNNGATTAC TAG ⓧ PCR w/
| | | | | | | | | | mismatch
...NNNCTAATGCCGATAGAGT... primer

Base pairs at end of primer match to subset of fragments

3 base extension will match ~1/64 fragments

Process DNA with RFLP

- Frequent restriction sites
- Many smaller fragments (~50 to ~1,000 bp)

Visualize subset of fragments

- Non-specific adapter annealed to the restriction site
- PCR primers bind adapter and restriction sites
- Separate fragments by electrophoresis and visualize

No prior sequence knowledge required

- Environmental samples
- Eukaryotes
- Commercial kits

Comparative genomic fingerprinting

Targets the presence of accessory genes

- Target selection enabled by the availability of genome sequences

Multiplex PCR and electrophoresis to detect ~40 genes

Isolates groups into subtypes based on patterns of genes present

- Assays available for *Campylobacter*, *Arcobacter*, and *Escherichia*



Gene sequence

DNA template and primer

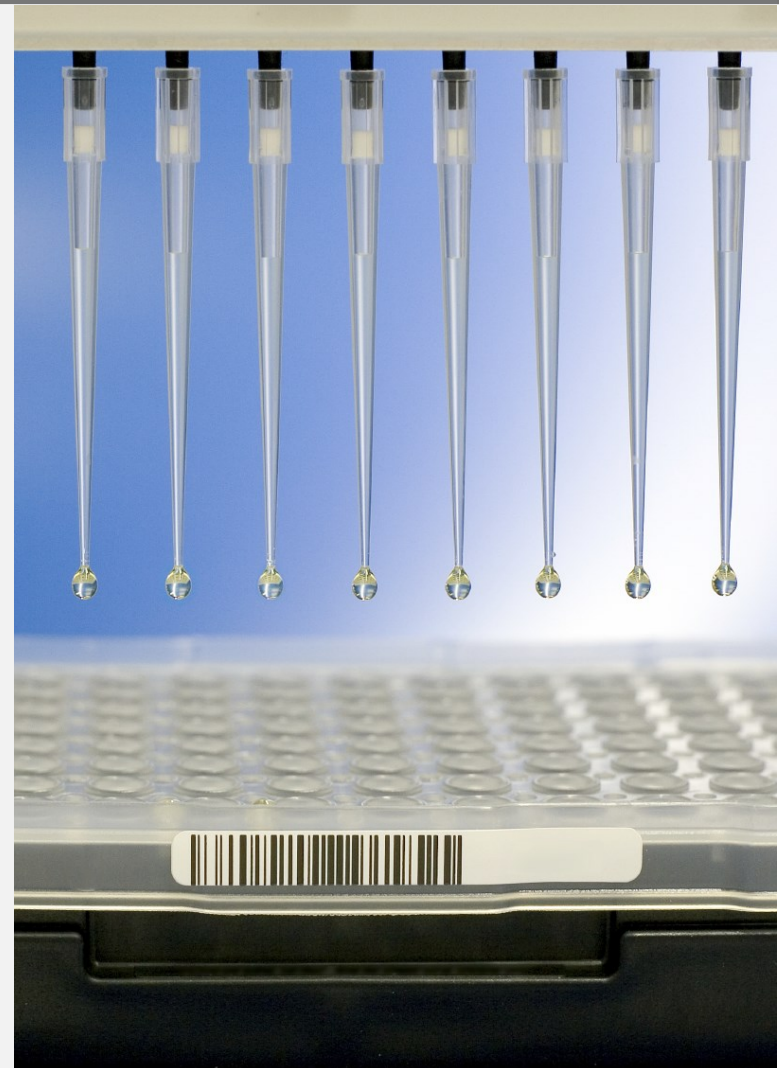
- Plasmid DNA
- Amplified PCR product
- Specific oligonucleotide primer

Amplification

- Incorporates nucleotides labeled with fluorescent dyes
- Sequence length up to ~700 bp

Analysis

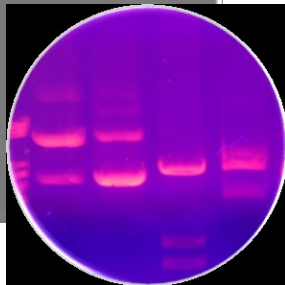
- Fluorescent dyes detected to determine nucleotide sequence
- Sequence results compared to sequence databases – NCBI, etc.



Gene sequence

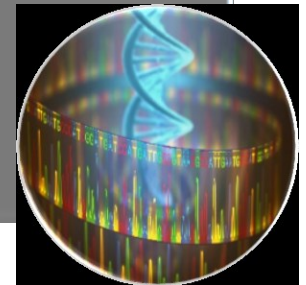
- Detects sequence changes that produce:
 - Altered restriction sites
 - Changes in size

Electrophoresis



- Detects all sequence changes within the area examined

Sequencing



Gene sequence

Methicillin-resistant *Staphylococcus aureus*

- *spa*
- *SCCmec*

Clostridium difficile

- *slpA* (molecular serotyping)

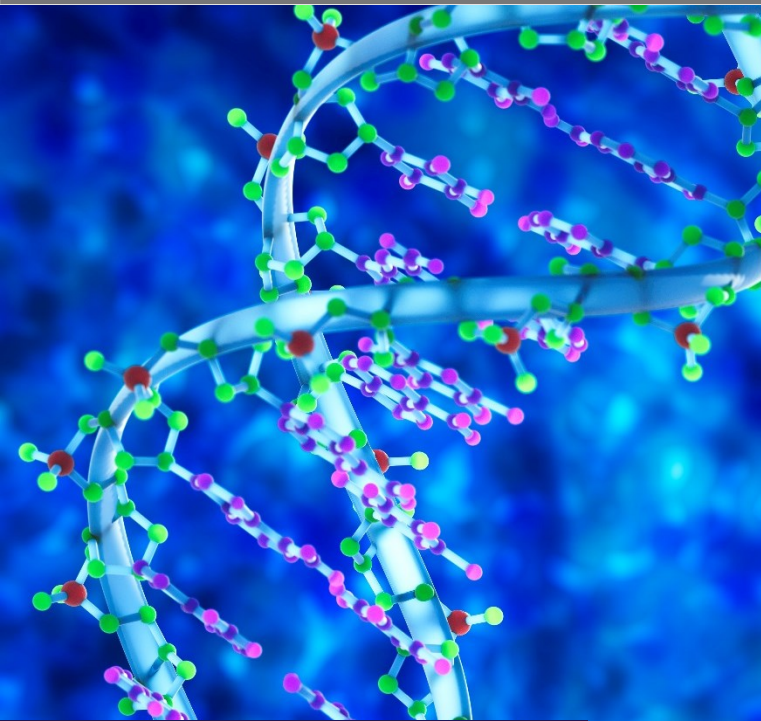
Rickettsia

- *ompA*



Genes need to have sufficient variability to discriminate and sufficient conservation to be in many strains

Multilocus sequence typing



Similar methodology to single gene sequence

Leverages greater number of genes for better discrimination

PCR amplify and sequence the defined regions of housekeeping genes – usually 7



Assign the allele number to each unique DNA sequence of each gene



Compare to database to assign sequence type number to each unique combination of alleles

Multilocus sequence typing

MLST databases

- Establish primer sets to define sequenced regions of each species
- Collect allele sequences and combinations from researchers
- Define sequence types for allele combinations
- Gather databases from multiple species into general MLST

MLST information

- Easy to store sequence text versus gel images
- Easy and fast to search text files

Multilocus sequence typing database

PubMLST Databases Downloads BIGSdb Contact Site map

Google™ Custom Search

Search

PubMLST

Welcome to PubMLST - Public databases for molecular typing and microbial genome diversity.

- [Log in or register for a PubMLST account](#)
- [Databases](#)
- [Download MLST definitions](#) - allelic profiles and sequences from all publicly accessible databases
- [BIGSdb](#) - software that runs most of the databases on this site
- [Documentation](#) - user and curator guides for BIGSdb
- [MLST schemes hosted on other sites](#)
- [Information about all species MLST databases and published schemes](#)
- [Recent publications that use or mention MLST](#)
- [News and updates](#)
- [Site map](#)
- [Policy document](#)
- [RESTful Application Programming Interface \(API\)](#)
- [International mirrors of this site: Primary | NL1 | NO1 | UK4 | US1](#)

Please [contact us](#) if you would like us to host a MLST database for a particular organism, or have a request for new functionality.

The primary PubMLST site is hosted at [The Department of Zoology](#), University of Oxford, UK and is funded by The Wellcome Trust.

News

2017-01-17: MLST schemes for *Macrocooccus canis* and *M. caseolyticus* have been developed by Vincent Perreten and Christian Strauss, University of Berne, Switzerland, and are now available.

2016-11-11: MLST schemes for *Brucella spp.* have been developed by Adrian Whatmore and colleagues, Animal and Plant Health Agency, UK, and are now available.

2016-11-09: MLST schemes for *Mycoplasmma iowae* and *M. synoviae* have been developed by Mohamed El-Gazzar and Mostafa Ghanem, The Ohio State University, USA, and are now available.

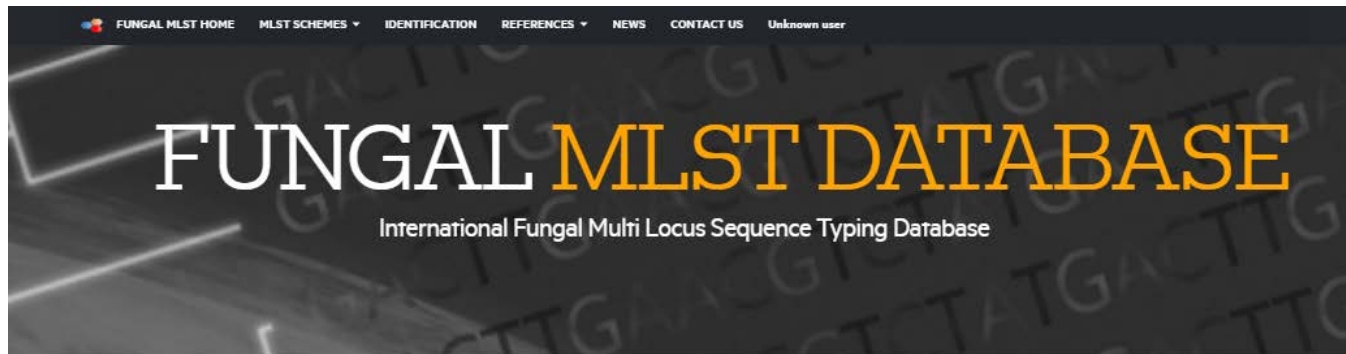
[Latest updates](#)

<http://pubmlst.org/>

University of Oxford, Wellcome Trust



Multilocus sequence typing database



<p>CG MLST</p> <p><i>Cryptococcus gattii</i></p> <p>Go</p>	<p>CN MLST</p> <p><i>Cryptococcus neoformans</i></p> <p>Go</p>	<p>PJ MLST</p> <p><i>Pneumocystis jirovecii</i></p> <p>Go</p>
<p>SAP MLST</p> <p><i>Scedosporium apiospermum</i></p> <p>Go</p>	<p>SAU MLST</p> <p><i>Scedosporium aurantiacum</i></p> <p>Go</p>	<p>PB MLST</p> <p><i>Scedosporium boydii</i></p> <p>Go</p>
<p>BA MLST</p> <p><i>Bipolaris australiensis</i></p> <p>Go</p>	<p>BH MLST</p> <p><i>Bipolaris hawaiiensis</i></p> <p>Go</p>	<p>BS MLST</p> <p><i>Bipolaris spicifera</i></p> <p>Go</p>

Multilocus sequence typing

Example species available in PubMLST databases

Bacteria			Eukaryotes
<i>Escherichia coli</i> #1	<i>Acinetobacter baumannii</i>	<i>Bacillus cereus</i>	<i>Candida albicans</i>
<i>Enterobacter cloacae</i>	<i>Borrelia</i> spp.	<i>Staphylococcus aureus</i>	<i>Candida tropicalis</i>
<i>Salmonella enterica</i>	<i>Burkholderia pseudomallei</i>	<i>Streptococcus pneumoniae</i>	<i>Aspergillus fumigatus</i>
<i>Klebsiella pneumoniae</i>	<i>Campylobacter jejuni</i>	<i>Arcobacter</i> spp.	<i>Penicillium marneffeii</i>
<i>Yersinia pseudotuberculosis</i>	<i>Helicobacter pylori</i>	<i>Wolbachia</i> spp.	<i>Kudoa septempunctata</i>
<i>Vibrio parahaemolyticus</i>	<i>Neisseria</i> spp.	<i>Xylella fastidiosa</i>	<i>Trichomonas vaginalis</i>
92 MLST schemes			9 MLST schemes

Multilocus sequence typing

PubMLST Database home Contents

Log in

Download allele sequences

Select loci by scheme | Alphabetical list | All loci by scheme

MLST

Locus	Download	Type	Alleles	Length (setting)	Min length	Max length	Full name/product	Curator(s)	Last updated
arcC		DNA	436	Fixed: 456 bp	435	457		K. Jolley	2017-01-23
aroE		DNA	588	Fixed: 456 bp	453	456		K. Jolley	2017-01-23
glpF		DNA	523	Fixed: 465 bp	450	466		K. Jolley	2017-01-23
gmk		DNA	297	Fixed: 417 bp	417	420		K. Jolley	2017-01-30
pta		DNA	468	Fixed: 474 bp	474	474		K. Jolley	2017-01-31
tpi		DNA	429	Fixed: 402 bp	402	402		K. Jolley	2017-01-30
yqiL		DNA	511	Fixed: 516 bp	516	516		K. Jolley	2017-01-23

Staphylococcus aureus

31,885 isolates

3,775 sequence types

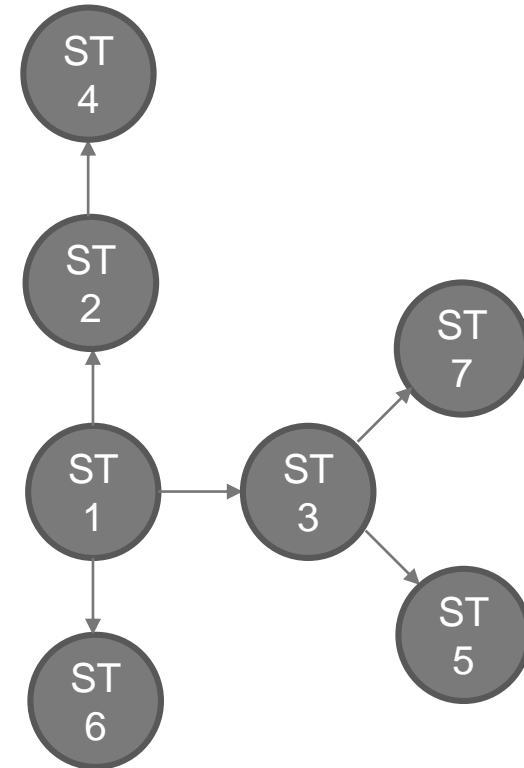
glpF_8	AGCTGGCGCGAAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	ST	arcC	aroE	glpF	gmk	pta	tpi	yqiL
glpF_22	AGCTGGCGCGAAATTAGGTGTTTTCTCTACAGCA ^T CGGCTATTAAGAATTACTTTGCCAACTTTT	1	1	1	1	1	1	1	1
glpF_309	AGCTGGCGCG ^A AAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	2	2	2	2	2	2	2	26
glpF_237	AGCTGGCGCGAAATTAGGTGTTTTCTCTACAGCACCG ^A GCTATTAAGAATTACTTTGCCAA ^T TTTT	3	1	1	1	9	1	1	12
glpF_324	AGCTGGCGCGAAATTAGGTGTTTTCTCTACAGCACCG ^A GCTATTAAGAATTACTTTGCCAA ^T TTTT	4	10	10	8	6	10	3	2
glpF_205	AGCTGGCGCG ^A AAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	5	1	4	1	4	12	1	10
glpF_82	AGCTGGCGCGAAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	6	12	4	1	4	12	1	3
glpF_267	AGCTGGCGCGAAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	7	5	4	1	4	4	6	3
glpF_358	AGCTGGCGCGAAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	8	3	3	1	1	4	4	3
glpF_298	AGCTGGCGCGAAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	9	3	3	1	1	1	1	10
glpF_212	AGCTGGCGCG ^A AAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	10	8	7	6	2	9	9	7
glpF_49	AGCTGGCGCG ^A AAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	11	1	24	1	4	12	1	10
glpF_207	AGCTGGCGCG ^A AAATTAGGTGTTTTCTCTACAGCACCGGCTATTAAGAATTACTTTGCCAACTTTT	12	1	3	1	8	11	5	11
		13	1	3	1	10	11	5	11
		14	1	13	1	1	12	11	13

Multilocus sequence typing

Sequence types are grouped together on the basis of similarity

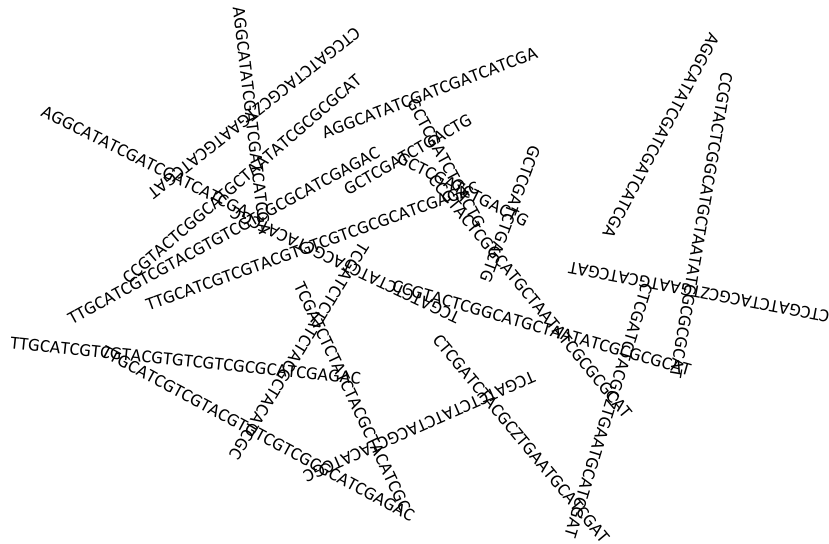
- Identical at number of alleles

Can be used to track changes to a population



Each arrow represents a change at one locus relative to the original sequence type

Whole genome sequence



Fragments are aligned and assembled into larger contigs

Pyrosequencing generates a very large number of small (~30-50 bp) fragments

- Automated process

Contigs are assembled to produce finished chromosomes

- Manual process

CCGTACTCGGCATGCTAATATCTCGATCTA
CTCGATCTACGCZTGAATGCATCGAT
TCGATTATCGATCGATCATCGA

...CCGTACTCGGCATGCTAATATCTCGATCTACGCZTGAATGCATCGATTATCGATCGATCATCGA...

Whole genome sequence

Multi-locus analysis

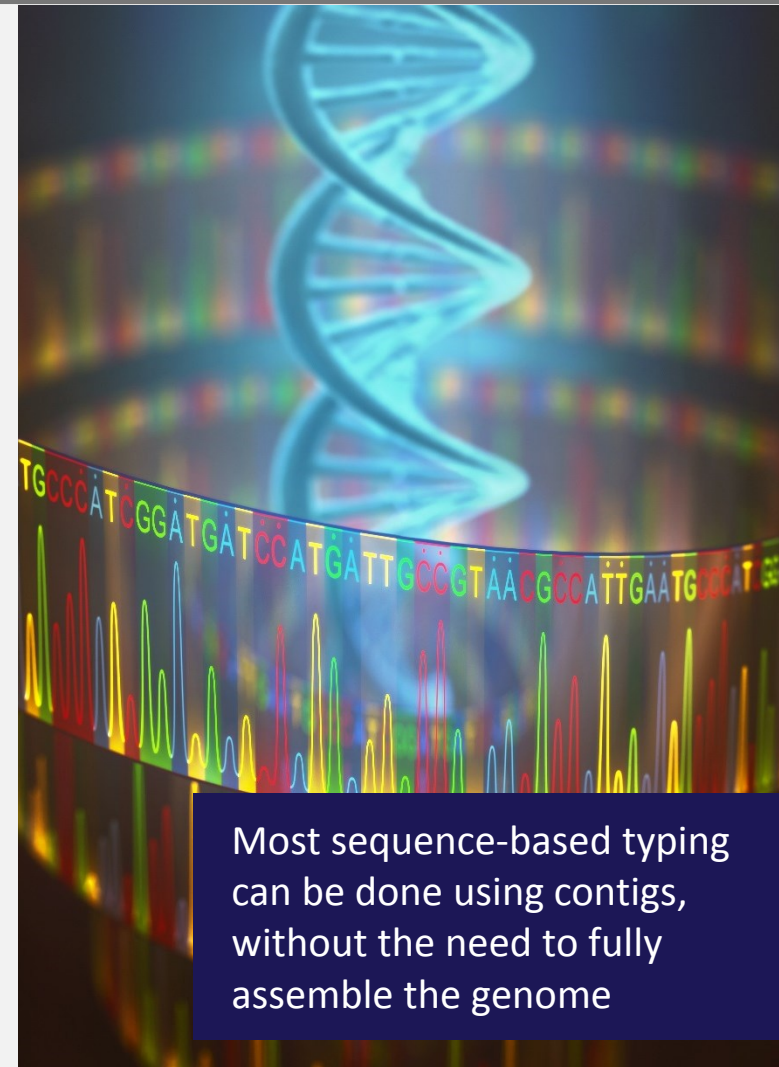
- MLST
- Pairwise SNP comparison
- Phylogenetic analysis

Single locus analysis

- Single gene typing (*e.g.* MRSA *spa*)
- Molecular typing
 - Identify serotype by specific alleles

Molecular detection of phenotypic typing markers

- Virulence genes, antimicrobial resistance genes, biochemical pathways, phage receptors



Whole genome sequence



Limitations to whole genome sequencing

- Cost
- Bioinformatics expertise

Currently used by national labs

- Centers for Disease Control and Prevention
- United States Department of Agriculture

Summary



- Strain typing identifies characteristics that differ between isolates of the same species
- Phenotypic typing methods, such as serotyping, offer speed and ease of use
- Genetic typing methods assess strain differences at the DNA level
 - Electrophoretic techniques look for changes in DNA fragment patterns
 - Sequencing techniques directly analyze the DNA sequence of one or more genes
- Genetic typing offers increasing levels of discrimination but requires more specialized equipment and training

Disclaimers



© 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.

Thank you for joining today!

Register for more ATCC webinars at
www.atcc.org/webinars

View our collection of antimicrobial-resistant reference materials at
www.atcc.org/superbugs

