

Prevent Analysis Variability by Using NGS-authenticated Microbial Genomes — Shift from Consensus to Discernible

Andrew Frank, Scientist Bioinformatician, ATCC

Credible Leads to Incredible™



About ATCC

- Founded in 1925, ATCC is a not-for-profit organization with HQ in Manassas, VA, and an R&D and Services center in Gaithersburg, MD
- World's largest, most diverse biological materials and information resource for microbes – the "gold standard"
- Innovative R&D company featuring gene editing, microbiome, NGS, and advanced models
- cGMP biorepository

- Partner with government, industry, and academia
- Leading global supplier of authenticated cell lines, microorganisms, and molecular standards
- Sales and distribution in 150 countries, 19 international distributors
- Talented team of 450+ employees, over onethird with advanced degrees



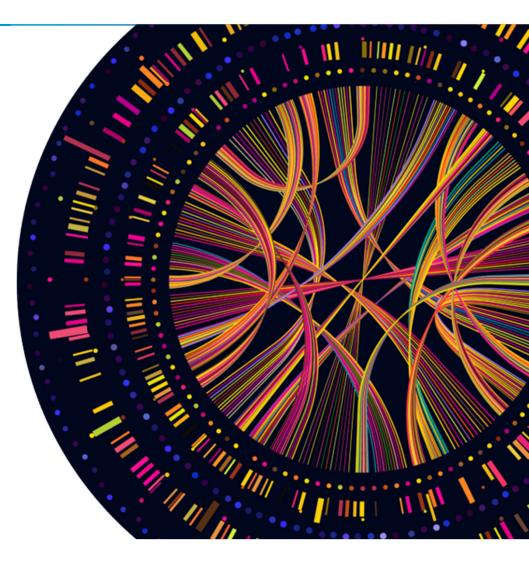
Overview

Enhanced Authentication Initiative

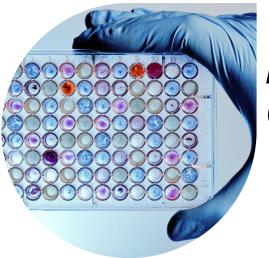
Enriching the characterization of ATCC's biological collections through whole-genome sequencing

ATCC Genome Portal

Providing reference-grade genomes matched to authenticated ATCC biological materials



Science is Changing



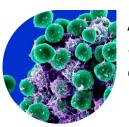
The Reproducibility Crisis



The Genomics Revolution



Modern Bacterial Strain Authentication Approaches



ATCC[®] 12228™ Staphylococcus epidermidis

ATCC	CERTIFICATE OF ANALYSIS			
	-			
Render	-			
		a presente proven 1981		
Name Price in Dates	-			
Aut Format	that sizes	a law		
	Ad applicable			
ap loadbas	11011			
Balled .		Specification .	-	
10,00		104411	1.4	
and a little		shape of		
and a state		11-694-48	1.404	
the of the strength		The reduced stage descended	No obtain and downed	
		(MA, N) 1084 (MA)	(No. 16 - 160 Mil	
			ter statig of take	
		Secondari 11 angli dan d	Accessible PCA angeliance of coloring genetic	
	(22) 10	canto polo	formale FS angliste of status parts formale all some spece	
I	Gent 10	International Sector Spectra Sector Spectra	hander of and an of	
l		international parts formated and some sparses and the source of the sour	An and a set of a set	
		International Sector Spectra Sector Spectra		
ĺ		International Sector Spectra Sector Spectra	An and a set of a set	

1. ATCC polyphasic testing

- Phenotypic testing
- Genotypic testing
- Functional analyses



2. Sequence the genome yourself

- Expensive but highly accurate
- Must be done as close to receipt of the item as possible, or else not reproducible



3. "Crowd-source" the genome

- Free but risky
- Requires bioinformatics expertise
- Issues with quality, traceability, etc.



The Enhanced Authentication Initiative

- 1. Enrich the characterization of ATCC biological collections
- Provide our customers with the wholegenome sequences of the specific, authenticated materials researchers need to generate credible data

ATCC[®]

ATCC Genome Portal

The ATCC Genome Portal is a cloud-based platform that enables users to easily browse genomic data and metadata by simply logging into the portal.



Download whole-genome sequences and annotations of ATCC materials



Search for nucleotide sequences or genes within genomes

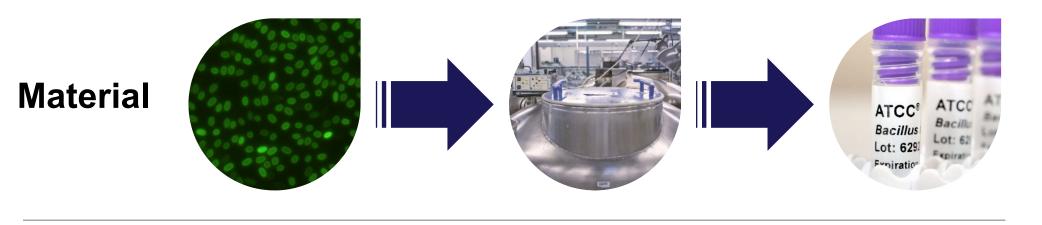
.1.1

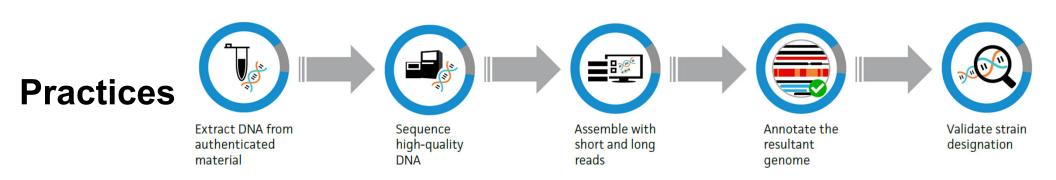
View genome assembly metadata and quality metrics

genomes.atcc.org



Differentiating ATCC by Material and Practices







ATCC Bioinformatics Best Practices

87

Extract DNA with optimized, proprietary protocols to get the best quality input material Avoid "garbage in, garbage out" data and results

Sequence on both Illumina[®] and Oxford Nanopore Technologies[®] Impose strict quality control thresholds to save only highest quality reads per instrument

Combine data from both technologies to achieve high-quality, complete genomes We perform hybrid genome assembly, leveraging the strengths of both platforms while avoiding biases and weaknesses of them individually

Annotate genomes to enable gene-level analyses Provide users with a reliable annotation, so they can identify genes of interest rapidly to enable gene-specific research

Confirm strain identity against highly curated databases We've partnered with One Codex, an industry leader in bioinformatic bacterial identification, to verify our strain designations by using *k*-mer and genomic distance approaches

ΔΤϹϹ



ATCC Extraction Quality Control

ATCC [®] no.	Species	PicoGreen [®] (ng/µL)	A _{260/} A ₂₈₀	DNA fragment size (range)**
8739DX™*	Escherichia coli	101.9	1.92	49.5 kb (1.5 – >60 kb)
13048DX™*	Klebsiella aerogenes	98.1	1.86	49.5 kb (1.6 – >60 kb)
11828DX™*	Cutibacterium acnes	197.7	1.84	29.8 kb (0.8 – >60 kb)
6538DX™*	Staphylococcus aureus	97.8	1.85	32.9 kb (2.7 – >60 kb)
BAA-2797DX™*	Pseudomonas aeruginosa	153.3	1.99	44.1 kb (1.1 – >60 kb)
824D-5™	Clostridium acetobutylicum	73.8	2.05	12.5 kb (4.6 – 57.8 kb)
6538D-5™	Staphylococcus aureus	37.1	2.00	26.2 kb (6.9 – >60 kb)
27774D-5™	Desulfovibrio desulfuricans	69.2	1.99	58.5 kb (13.3 – >60 kb)
11842D-5™	Lactobacillus delbrueckii	64.8	2.02	41.9 kb (6.1 – >60 kb)
15697D-5™	Bifidobacterium longum	76.2	1.95	51.3 kb (10.5 – >60 kb)

* New NGS-ready DNA product ** Main peak reported



NGS-Ready DNA Products Now Available



NGS-Ready DNA products (-DX) now available to customers

- ≥ 10 µg of DNA
- DNA fragment sizes ≥ 20 kb
- Launched with <u>50</u> of the most popular ATCC bacterial strains
- Check them out at atcc.org/NGS-ready



ATCC Bioinformatics Best Practices



Extract DNA with optimized, proprietary protocols to get the best quality input material Avoid "garbage in, garbage out" data and results

Sequence on both Illumina[®] and Oxford Nanopore Technologies[®] Impose strict quality control thresholds to save only highest quality reads per instrument

Combine data from both technologies to achieve high-quality, complete genomes We perform hybrid genome assembly, leveraging the strengths of both platforms while avoiding biases and weaknesses of them individually

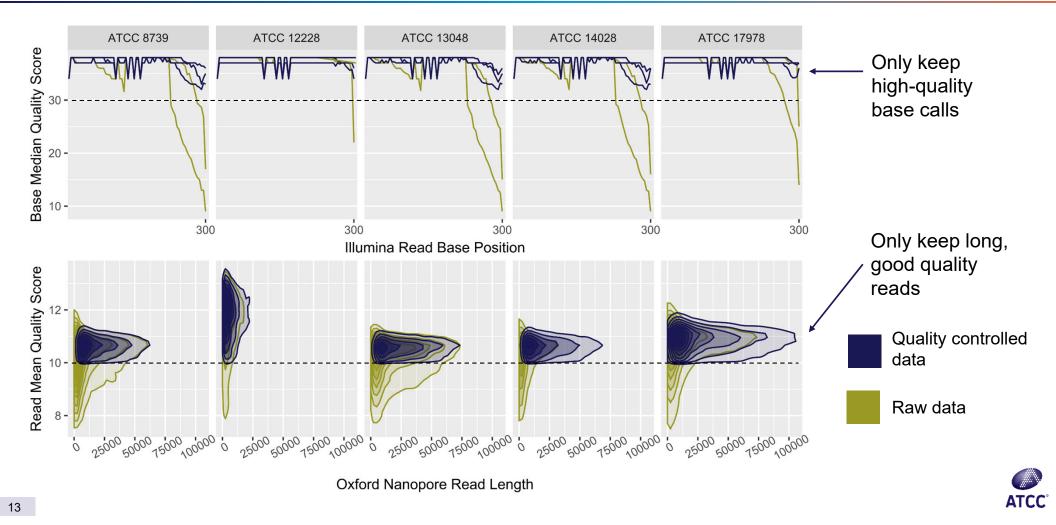
Annotate genomes to enable gene-level analyses Provide users with a reliable annotation, so they can identify genes of interest rapidly to enable gene-specific research

Confirm strain identity against highly curated databases We've partnered with One Codex, an industry leader in bioinformatic bacterial identification, to verify our strain designations by using *k*-mer and genomic distance approaches

ΔΤϹϹ



Sequencing QC – Read Trimming/Filtering



Sequencing QC – Changing Conditions

Test	Species	ATCC [®] no.	Reference Genome	Analysis	# of SNPs	# of indels	Number of variants	Variant coverage
A h	Mycoplasma hominis	23114™	GCA_000085865.1	Preparation 1	14	10	24	1042.1
				Preparation 2	14	10	24	900.0
	Cutibacterium acnes	11828™	GCA_000231215.1	Preparation 1	28	37	65	121.1
				Preparation 2	28	39	67	128.6
В	Clostridium acetobutylicum	824D-5™	GCA_000008765.1	Kit 1	171	55	226	95.9
				Kit 2	170	55	225	202.0
	Aeromonas hydrophila	7966D-5™	GCA_000014805.1	Kit 1	1	1	2	216.8
				Kit 2	1	1	2	203.0
С	Escherichia coli	700926™	GCA_000005845.2	Extraction 1	0	1	1	137.0
				Extraction 2	0	1	1	186.3
	Streptococcus pyogenes	19615™	GCA_000743015.1	Extraction 1	2	44	46	314.2
				Extraction 2	2	41	43	460.2

Test A: Same DNA sequenced using two different DNA preparations

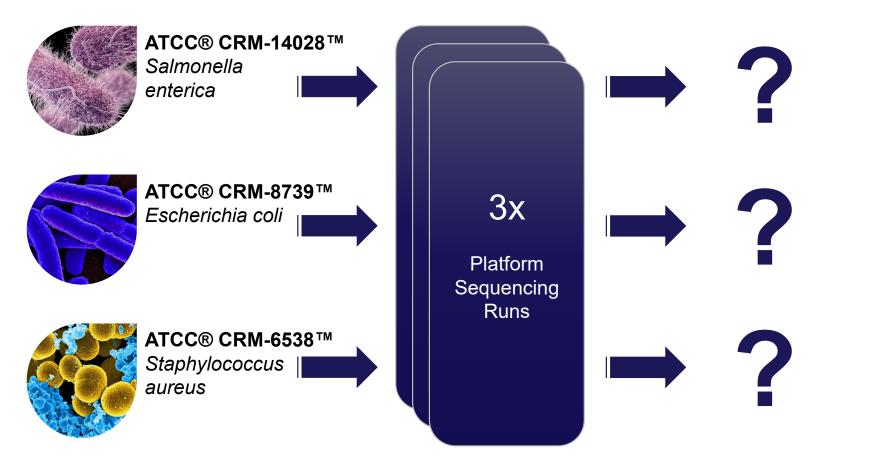
Test B: Same DNA sequenced with two different library kits

Test C: Same strain extracted with two different methods



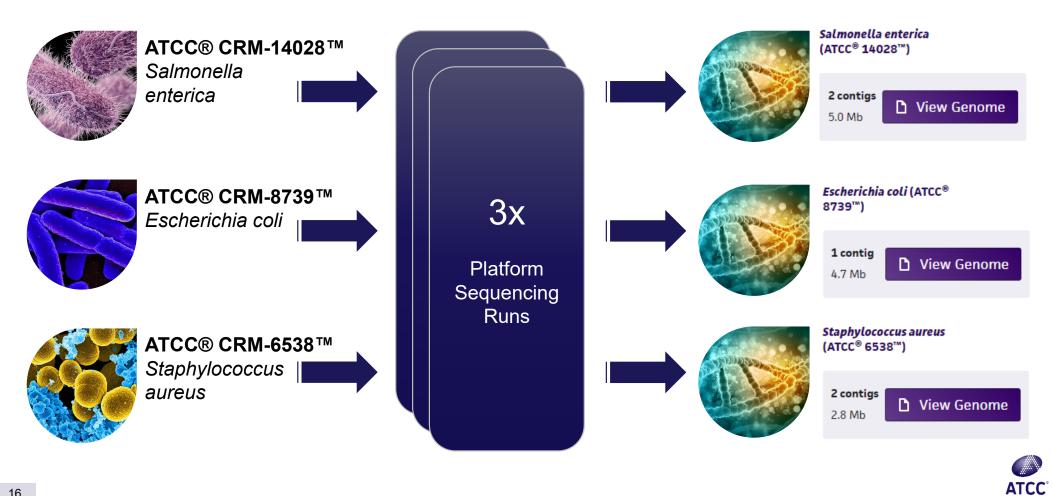
14

Application – NGS Proficiency Testing





Application – NGS Proficiency Testing



ATCC Bioinformatics Best Practices

Extract DNA with optimized, proprietary protocols to get the best quality input material Avoid "garbage in, garbage out" data and results

Sequence on both Illumina[®] and Oxford Nanopore Technologies[®] Impose strict quality control thresholds to save only highest quality reads per instrument

Combine data from both technologies to achieve high-quality, complete genomes We perform hybrid genome assembly, leveraging the strengths of both platforms while avoiding biases and weaknesses of them individually

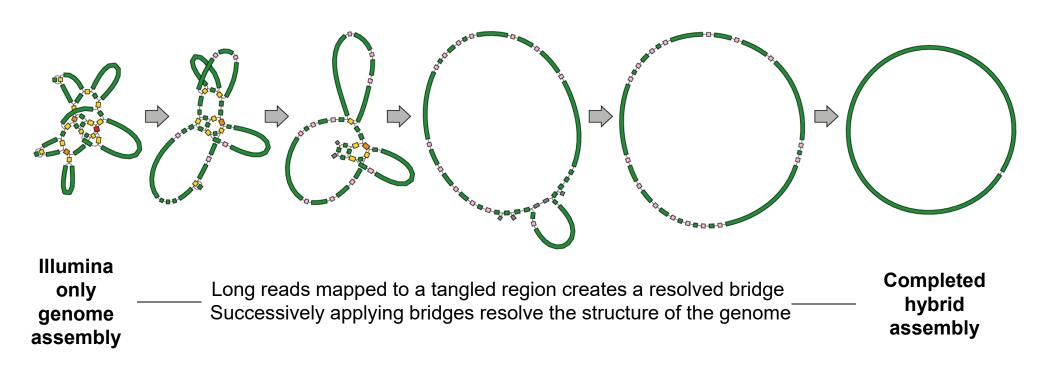
Annotate genomes to enable gene-level analyses Provide users with a reliable annotation, so they can identify genes of interest rapidly to enable gene-specific research

Confirm strain identity against highly curated databases We've partnered with One Codex, an industry leader in bioinformatic bacterial identification, to verify our strain designations by using *k*-mer and genomic distance approaches

ΔΤϹϹ

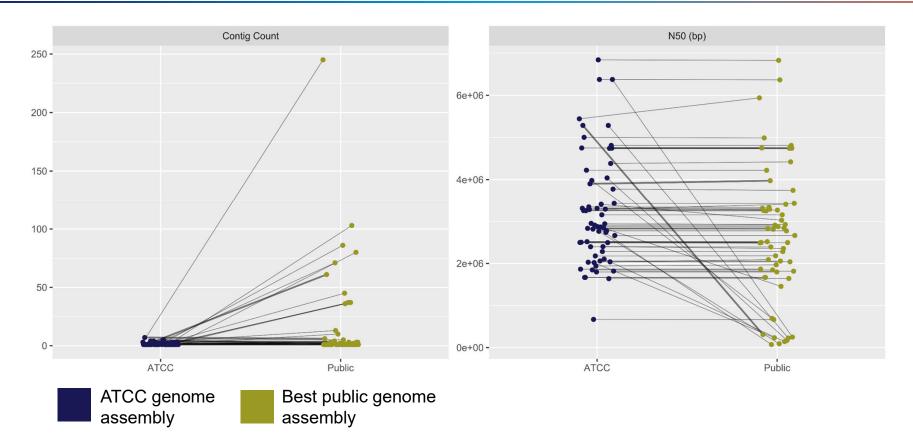


Hybrid Genome Assembly





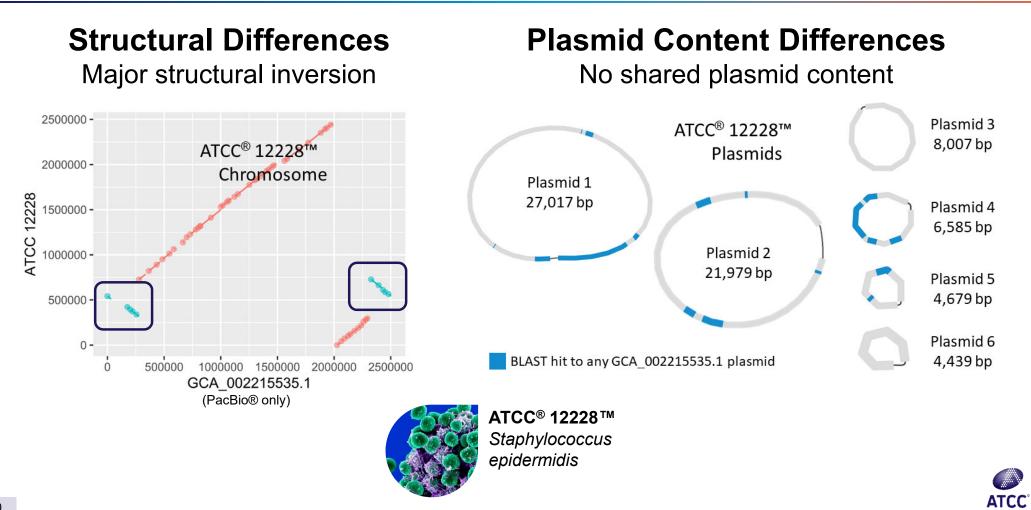
ATCC's Assemblies Substantially Improve Upon Public Assemblies



Downward trend in contig count and **upward** trend in N50 indicate ATCC is producing more complete genomes than the best publicly available alternative genome



ATCC's Assemblies Substantially Improve Upon Public Assemblies



ATCC Bioinformatics Best Practices



Extract DNA with optimized, proprietary protocols to get the best quality input material Avoid "garbage in, garbage out" data and results

Sequence on both Illumina[®] and Oxford Nanopore Technologies[®] Impose strict quality control thresholds to save only highest quality reads per instrument

Combine data from both technologies to achieve high-quality, complete genomes We perform hybrid genome assembly, leveraging the strengths of both platforms while avoiding biases and weaknesses of them individually

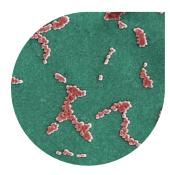
Annotate genomes to enable gene-level analyses

Provide users with a reliable annotation, so they can identify genes of interest rapidly to enable gene-specific research

Confirm strain identity against highly curated databases We've partnered with One Codex, an industry leader in bioinformatic bacterial identification, to verify our strain designations by using *k*-mer and genomic distance approaches

ΔΤϹϹ

Application – Variation in AMR Genes



ATCC[®] BAA-2803™

Acinetobacter baumannii Global Priority Superbug

- ATCC hosts 57 strains of WHO priority pathogens the Global Priority Superbug collection
 - Genomes being made available on the ATCC Genome Portal
 - Rich source data
 - Susceptibility data including MIC values and susceptibility profiles for targeted drugs
- More info at atcc.org/globalprioritysuperbugs
- In A. baumannii, mutations of gyrA can confer high levels of resistance to fluoroquinolone antibiotics (i.e. Ciprofloxacin)
- The ATCC Genome Portal can rapidly search for sequences of interest within our genome collection, up to 4000X faster than BLAST



ATCC Bioinformatics Best Practices

87

Extract DNA with optimized, proprietary protocols to get the best quality input material Avoid "garbage in, garbage out" data and results

Sequence on both Illumina[®] and Oxford Nanopore Technologies[®] Impose strict quality control thresholds to save only highest quality reads per instrument

Combine data from both technologies to achieve high-quality, complete genomes We perform hybrid genome assembly, leveraging the strengths of both platforms while avoiding biases and weaknesses of them individually

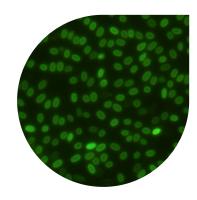
Annotate genomes to enable gene-level analyses Provide users with a reliable annotation, so they can identify genes of interest rapidly to enable gene-specific research

Confirm strain identity against highly curated databases

We've partnered with One Codex, an industry leader in bioinformatic bacterial identification, to verify our strain designations by using *k*-mer and genomic distance approaches

ΔΤϹϹ

Enhanced Authentication Reveals Hidden Biodiversity





- Taxonomy of *Bacillus* continues to be rapidly updated this decade
- Sequenced and assembled genomes for 61 of ATCC's Bacillus cereus and Bacillus subtilis group strains, some of our oldest strains
- Measured genomic distance to type strains of all taxonomically valid *Bacillus*
 - -12 items are being moved into *Bacillus* species described only in the last decade
 - -1 item is being moved into an older existing *Bacillus* species
 - -9 items are being moved to taxonomically undescribed species



Application – Ascertain Strain Identity



ATCC® 33397™ Streptococcus anginosus



ATCC® **27335™** Streptococcus intermedius

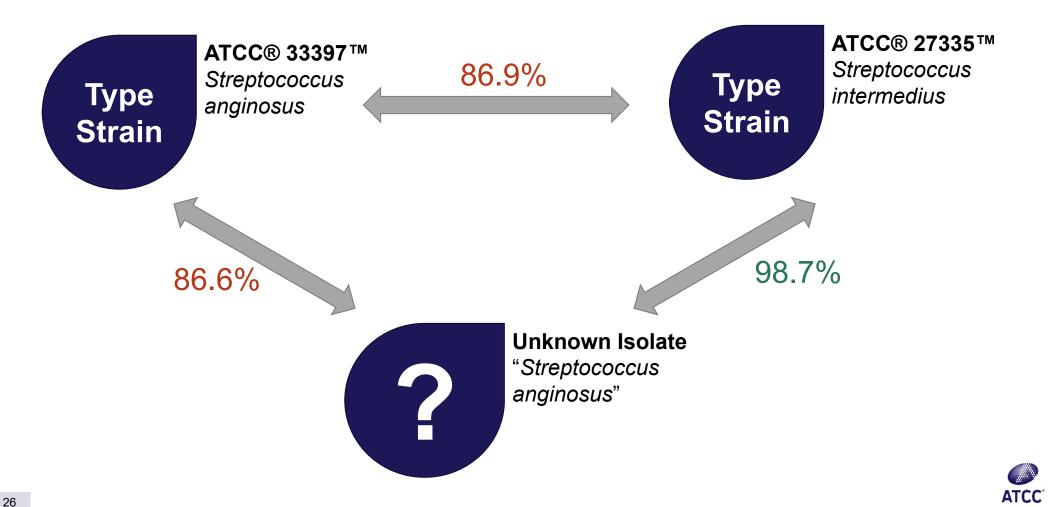


Unknown Isolate "Streptococcus anginosus"



25

Application – Ascertain Strain Identity



ATCC Genome Portal

The ATCC Genome Portal is a cloud-based platform that enables users to easily browse genomic data and metadata by simply logging into the portal.



Download whole-genome sequences and annotations of ATCC materials



Search for nucleotide sequences or genes within genomes

.1.1

View genome assembly metadata and quality metrics

genomes.atcc.org





Thank you



ATCC Genome Portal

The ATCC Genome Portal is a cloud-based platform that enables users to easily browse genomic data and metadata by simply logging into the portal.



Download whole-genome sequences and annotations of ATCC materials



Search for nucleotide sequences or genes within genomes

.1.1

29

View genome assembly metadata and quality metrics

genomes.atcc.org

© 2019 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.

