



The  
Microbiology  
NETWORK

# Microbiological Quality Control as Described in the Compendia

Scott Sutton, Ph.D.

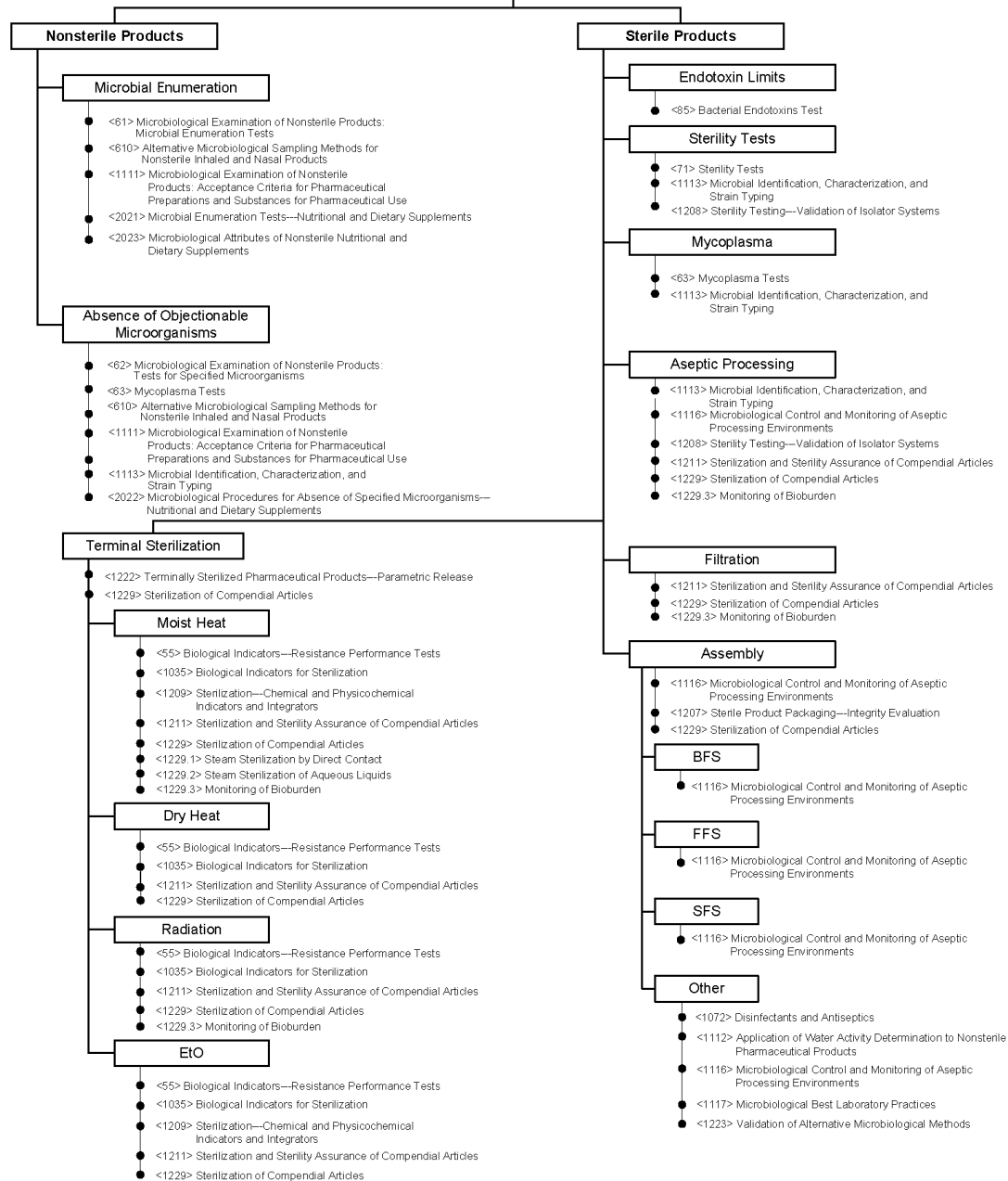
[www.microbiol.org](http://www.microbiol.org)

# Microbiological Quality Control as Described in the Compendia

- <51> Antimicrobial Effectiveness Test
- <61>, <62>, <1111> Microbial Limits Tests
- <71> Sterility Testing
- <1117> Best Microbiological Laboratory Practices



# Chart 10. Microbiology



# Antimicrobial Efficacy Test

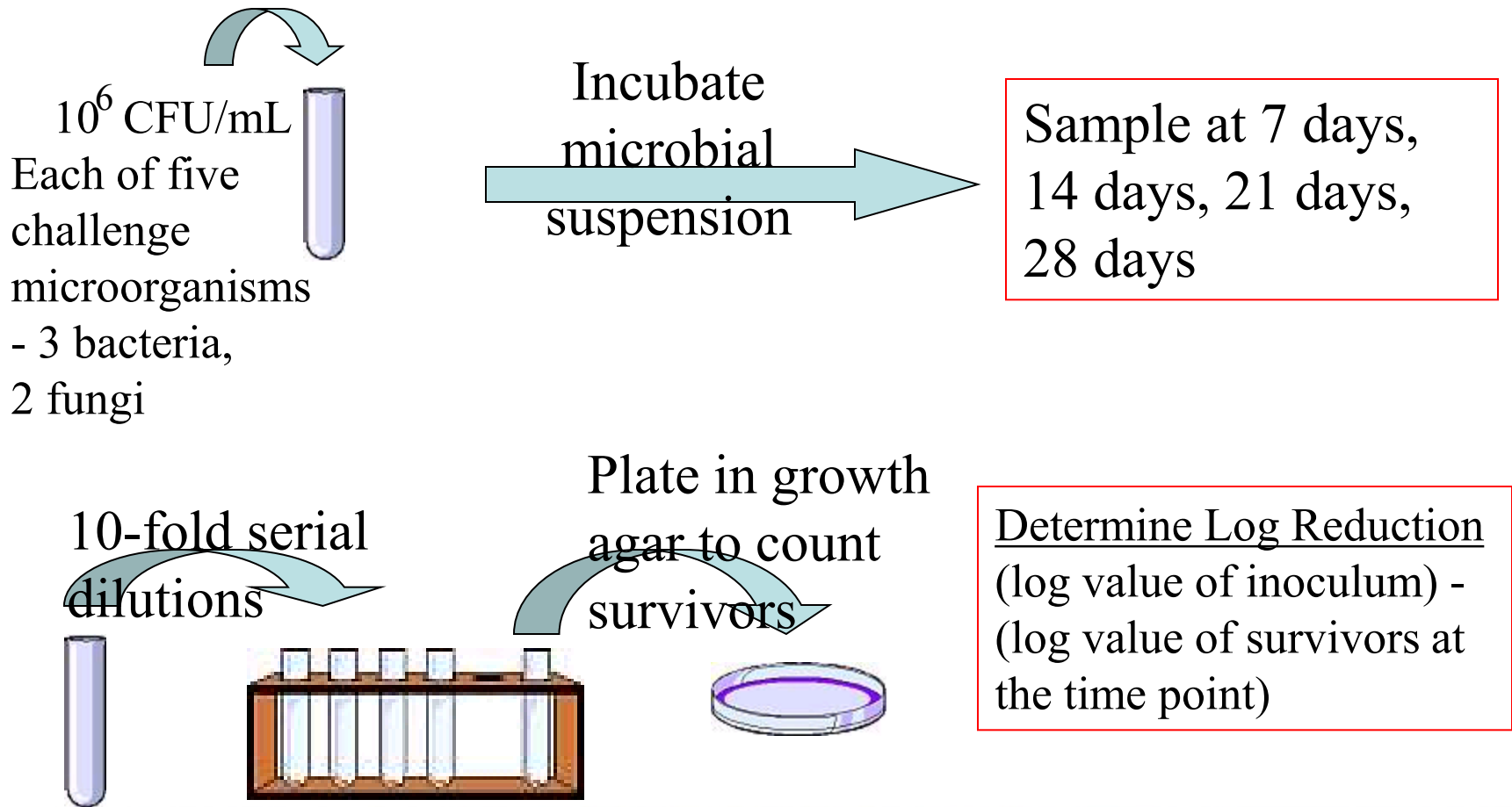
## USP <51>

- Designed to demonstrate the ability of a multidose product to withstand microbial challenge.
- High-level challenge, frequent sampling of the challenge suspension for survivors
- Completely rewritten for clarity in 2015
- Method Suitability included in test





# Antimicrobial Efficacy Test



# USP Criteria for Passage

	Log <sub>10</sub> Reduction			
	<u>7 Day</u>	<u>14 Day</u>	<u>21 Day</u>	<u>28 Day</u>
1A: Bacteria	1.0	3.0	--	NI
1A: Fungi	--	NI	--	NI
1B: Bacteria	--	2.0	--	NI
1B: Fungi	--	NI	--	NI
1C: Bacteria	--	1.0	--	NI
1C: Fungi	--	NI	--	NI
2: Bacteria	--	NI	--	NI
2: Fungi	--	NI	--	NI



# AET Categories

<u>Category</u>	<u>Product Description</u>
1	Injections and other parenterals, otic, sterile nasal products, and ophthalmics
2	Topical Products
3	Oral Products
4	Liquid Antacids



# Scope of AET

- It can
  - Provide relative estimates of the biological activity of a preservative system in a particular formulation at a particular time.
- It cannot
  - Predict the preservative efficacy of the multidose finished product in all patients hands under all conditions.





# Major Sources of Variability

- Preparation and handling of inocula
  - Source of culture
  - Growth conditions
    - Solid vs liquid
    - Length of time/Temperature incubated
  - Harvesting, buffer composition, storage conditions
- Recovery Conditions
  - Determined in Method Suitability Study
- Plating conditions
  - Colony counting rules
  - Math

**Proactive  
Documentation**



# Unique Microorganisms

FDA regulatory focus in recent years –

- *Bacillus cereus*  
Gram-positive, facultatively aerobic sporeformer.  $\beta$ -hemolytic and may be an emerging pathogen.
- *Burkholderia cepacia*  
Pseudomonad, clearly pathogenic to cystic fibrosis patients who can develop pneumonia.  
**THE CONCERN IS THAT CF PATIENTS ARE EXPOSED TO PRODUCTS THAT MAY LEAD TO PNEUMONIA.**



# Microbiological Quality Control as Described in the Compendia

- <51> Antimicrobial Effectiveness Test
- <61>, <62>, <1111> Microbial Limits Tests
- <71> Sterility Testing
- <1117> Best Microbiological Laboratory Practices



# Compendial Microbial Limits Tests

## Internationally Harmonized

- USP <61> Enumeration, EP 2.6.12
- USP <62> Specified Organisms, EP Chapter 2.6.13
  - Test for *Staphylococcus aureus*
  - Test for *P. aeruginosa*
  - Test for *Salmonella spp*
  - Test for *Escherichia coli*
  - Test for Bile-tolerant Gram-negative Bacteria
  - Test for *Clostridia*
  - Test for *Candida albicans*
- USP <1111> Guidance on Microbial Quality, EP 5.1.4



# USP <61> - Enumeration

- Sampling Plan
- Categories
- Methodology
  - Membrane Filtration
  - Plate Count: Pour Plate
  - Plate Count: Spread Plate
  - Most Probable Number (MPN)
- Method Suitability and Growth Promotion Requirements





# <62> - Specific Tests

- Test for *Staphylococcus aureus*
- Test for *P. aeruginosa*
- Test for *Salmonella spp*
- Test for *Escherichia coli*
- Test for Bile-tolerant Gram-negative Bacteria
- Test for *Clostridia*
- Test for *Candida albicans*



# “Microbial Attributes”

## USP <1111> EP 5.1.4

- One page in USP, two tables:
  - Table 1 – Acceptance criteria for Microbial Quality of Nonsterile Dosage Forms
  - Table 2 – Acceptance Criteria of Microbiological Quality of Nonsterile Substances for Pharmaceutical Use (TAMC- $10^3$  CFU/g or mL, TYMC- $10^2$ )
- Significance of other organisms should be evaluated



# Evaluation of Other Organisms

- Route of Administration
- Nature of the Product
- Method of Application
- Intended Recipient
- Use of Immunosuppressive Drugs
- Presence of Disease, Wounds, Organ Damage



# Microbiological Quality Control as Described in the Compendia

- <51> Antimicrobial Effectiveness Test
- <61>, <62>, <1111> Microbial Limits Tests
- <71> Sterility Testing
- <1117> Best Microbiological Laboratory Practices



# Sterility Testing

- Two separate tests
  - Membrane Filtration
  - Direct Transfer
- 20 Units, 2 media & temperatures
- Requires Growth
  - Incubation period - 14 days





# Membrane Filtration

- Filter required amount of product through two filters
- Neutralize/Rinse
  - 3 100 mL volumes suggested
  - Formulations for dilution fluids suggested
- One filter into Soybean Casein Digest Broth (SCDB or TSB) – incubate at 20-25°C for 14 days
- One filter into Fluid Thioglycollate Medium (FTM) – incubate at 30-35°C for 14 days



# Direct Inoculation

- Place required amount of product into sufficient recovery medium (with neutralizers?)
  - Soybean Casein Digest Broth (SCDB or TSB)
    - incubate at 20-25°C for 14 days
  - Fluid Thioglycollate Medium (FTM) – incubate at 30-35°C for 14 days



# Product Requirements

- Minimum Quantity per Unit for Each Medium detailed in Table 2 of chapter
- Minimum Number of Units to be tested detailed in Table 3 of chapter



# Method Suitability Test

Can we neutralize any antimicrobial properties of the medication?

Use specified challenge organisms

Use specified total amounts of products



# Method Suitability Test for Each Challenge Organism

- Filter maximal amount of medication to be tested
- Filter 2 volumes (100 mL?) of diluting fluid
- Add third volume, inoculate with <100 CFU challenge organism
- Filter
- Show microbial growth from filter in relevant medium at relevant temperature in 5 days





# Scope of Sterility Tests

- They can provide a recognized, standardized test.
- They cannot prove product sterility.
  - Limited sample size
  - Can only show what can grow



# Microbiological Quality Control as Described in the Compendia

- <51> Antimicrobial Effectiveness Test
- <61>, <62>, <1111> Microbial Limits Tests
- <71> Sterility Testing
- <1117> Best Microbiological Laboratory Practices



# USP <1117> Importance Aspects of Control

- Control of Media
- Control of Test Strains
- Control of Equipment
- Lab Lay-out and Operations
- Sample Handling
- Microbiological Media Incubation Times
- Training of Staff
- Laboratory Resources
- Control of Data and Documentation
- Interpretation of Results



# Maintenance of Microbiological Cultures

- Must be handled carefully
- Confirm ID of culture from culture collection before use
- Resuscitate cultures as per manufacturer's instructions
- Use a seed lot technique, do not enter a master vial more than once or refreeze stock
- Track number of passages (“Any form of subculturing is considered to be a transfer/passage”).



# Microbiological Quality Control as Described in the Compendia

- <51> Antimicrobial Effectiveness Test
- <61>, <62>, <1111> Microbial Limits Tests
- <71> Sterility Testing
- <1117> Best Microbiological Laboratory Practices







The  
Microbiology  
NETWORK

# Thank you for your attention

Scott Sutton, Ph.D.

[scott.sutton@microbiol.org](mailto:scott.sutton@microbiol.org)

+1 585-298-0767

<http://www.linkedin.com/in/scottvwsutton>

Twitter - [@microbiologynet](https://twitter.com/microbiologynet)

[www.microbiol.org](http://www.microbiol.org)





# ALWAYS IMPROVING QC ATCC SOLUTIONS YOU TRUST FOR THE QUALITY OF YOUR BRAND

Liz Kerrigan

Director of Standards, Sales & Marketing, ATCC

March 19, 2015



THE ESSENTIALS OF LIFE SCIENCE RESEARCH  
GLOBALLY DELIVERED™

# About ATCC

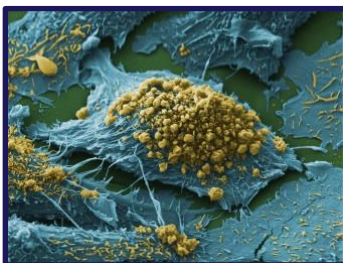
- Founded in 1925, ATCC is a non-profit organization with headquarters in Manassas, VA
- World's premiere biological materials resource and standards development organization
- ATCC collaborates with and supports the scientific community with industry-standard products and innovative solutions
- Broad range of biomaterials
  - Continuous cell lines, iPSCs, primary cells, and hTERT immortalized cells
  - Bacteria, fungi, yeasts, protists, and viruses
  - Microbial and tumor cell panels
  - Genomic and synthetic nucleic acids
  - Media, sera, and reagents



# ATCC complete solutions for pharmaceutical QC



- Proficiency testing programs
  - PHARMASSURE – Test materials for chemical and microbiological analyses as well as sterility testing



- Mycoplasma quality control
  - Titered strains
  - Quantitative nucleic acids



- ATCC<sup>®</sup> Minis
  - ATCC Genuine Cultures<sup>®</sup> packaged in a convenient, single-use glycerol stock



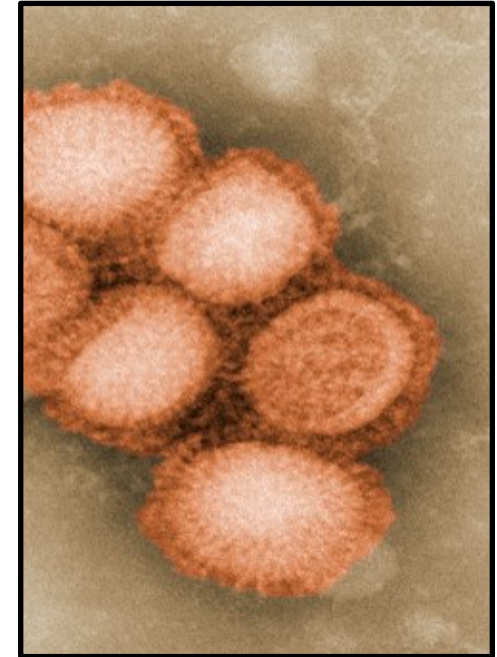
# ATCC Genuine Cultures®

- **Microbial Portfolio**

- 18,000 bacterial strains
- 3,000 human and animal viruses
- 50,000 yeast and fungal strains
- 2,000 parasites
- Nearly 1,000 ATCC Genuine Nucleics® from the collection

- **Brand Recognition**

- Standard/reference cultures
- Organizations and regulatory agencies specify ATCC cultures (USP, ISO, FDA, CLSI, USDA, ASTM, AOAC, and more)
- Over 475 microbial cultures recommended as reference strains in microbial collection
- Cultures are always authenticated



# Passage matters

USP clearly states that the working cultures used for testing should not be more than five passages from the ATCC reference culture.

The USP 37-NF32 <51> states:

“The viable microorganisms used in the test must not be more than five passages removed from the original ATCC culture.”



# Microbial strain authentication



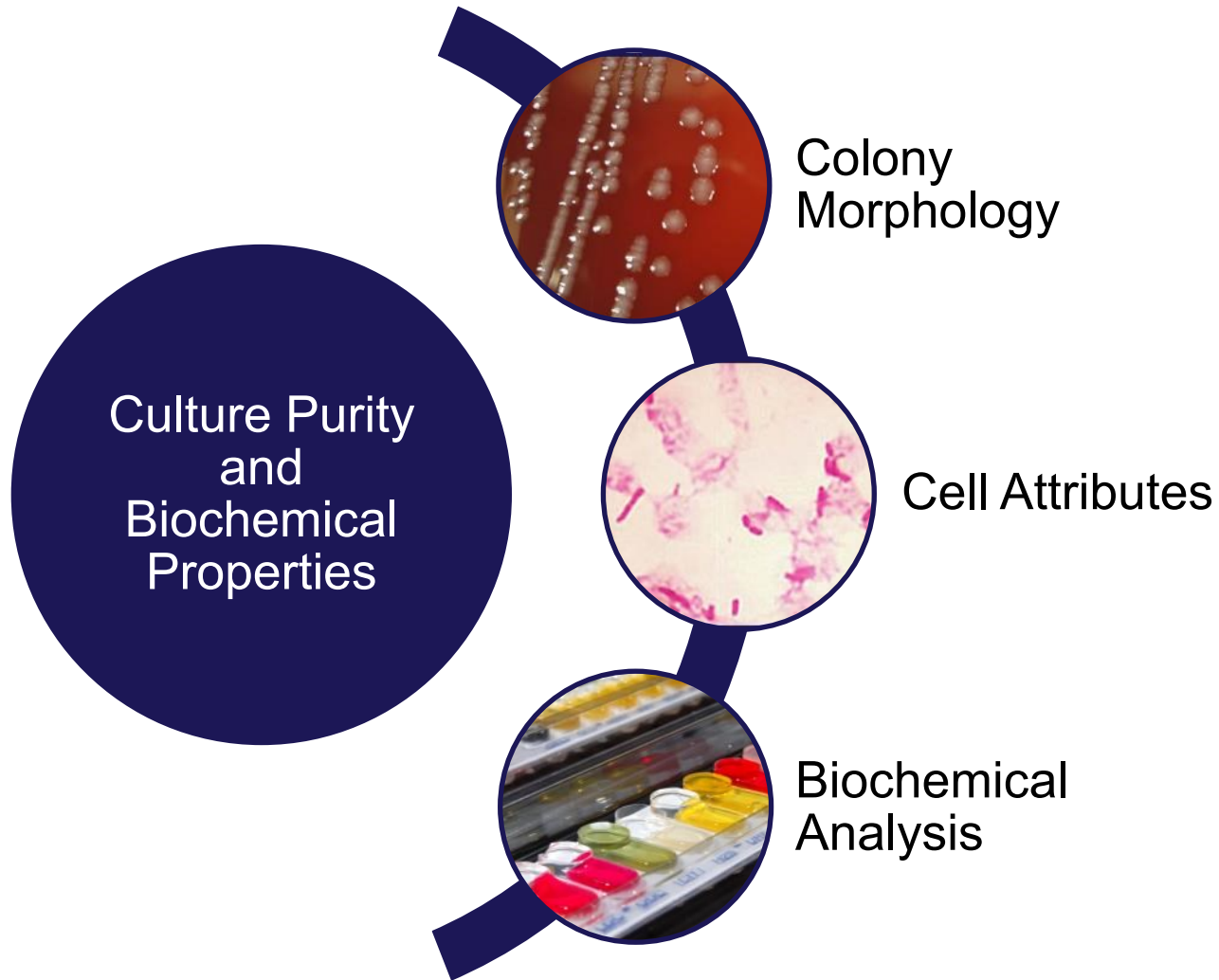
ATCC utilizes both classical and modern techniques

- Phenotypic analysis
- Genotypic & proteotypic analyses
- Functional analysis

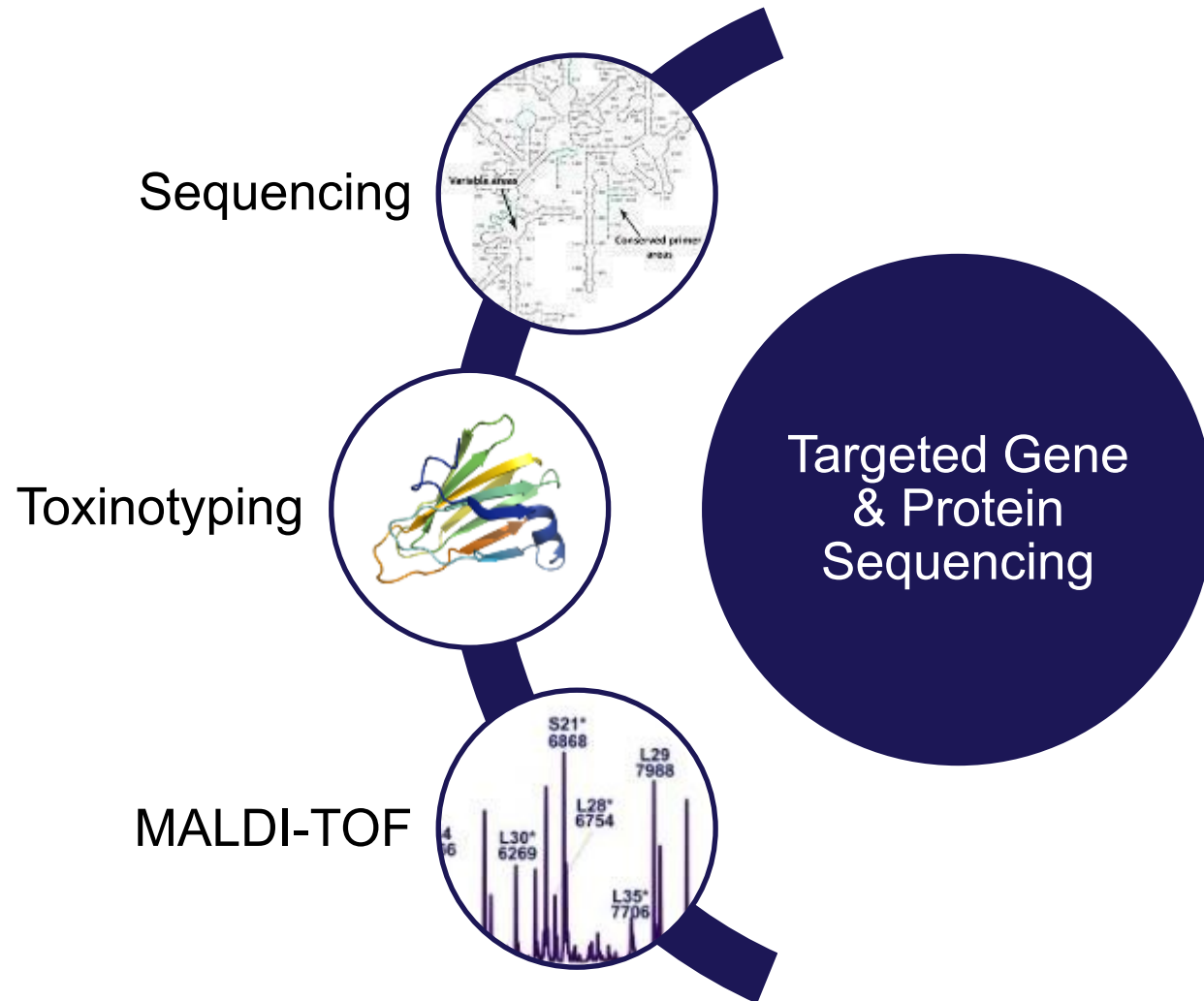
No single method of identification is sufficient



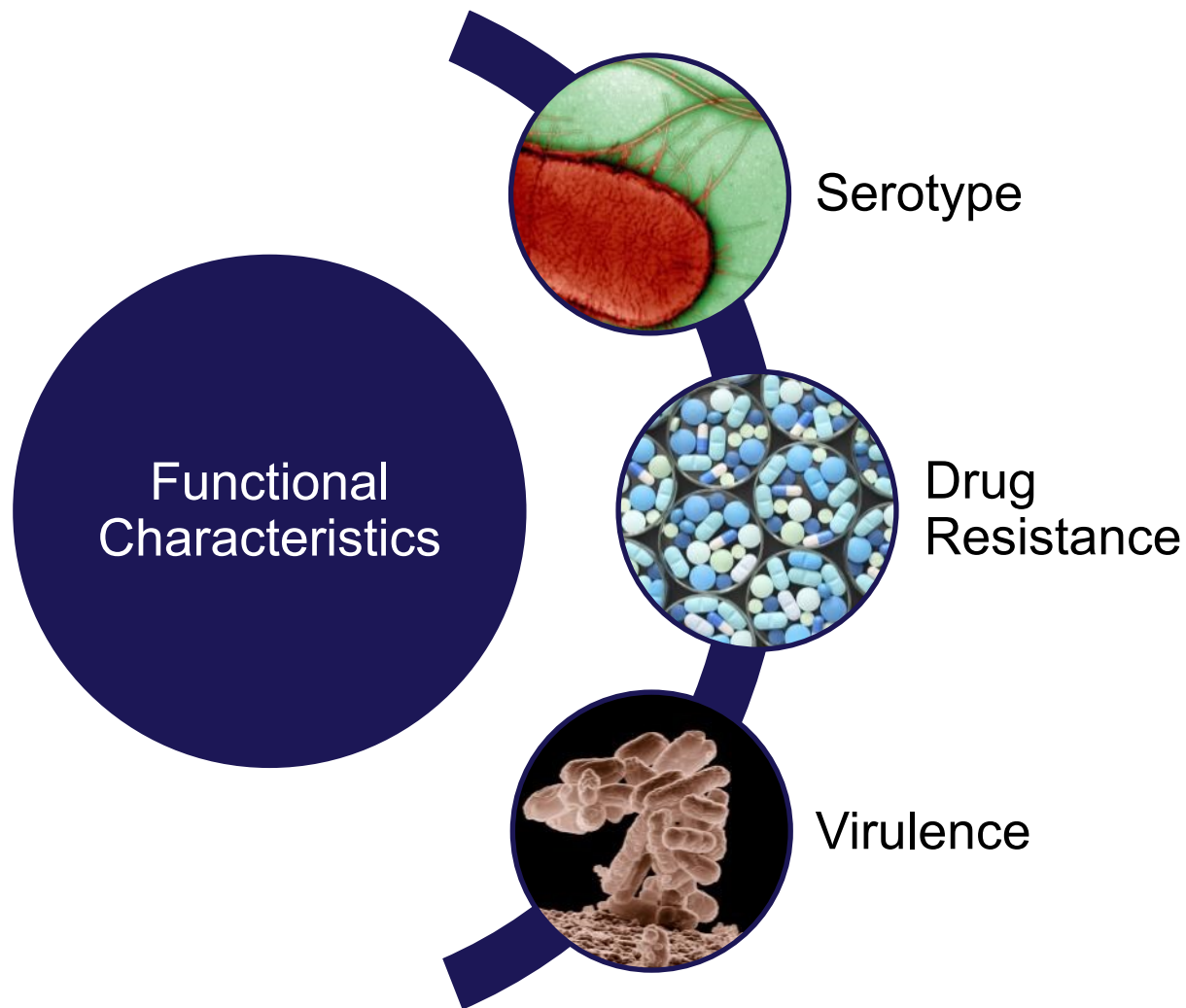
# Phenotypic testing



# Genotypic & proteotypic testing

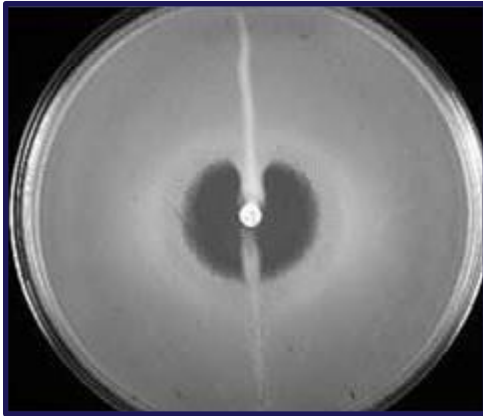


# Functional testing



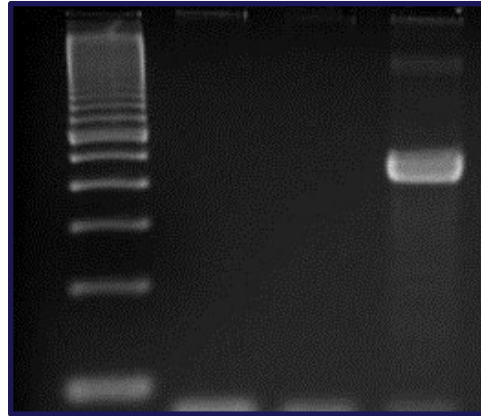
# Verification of drug resistance

## Modified Hodge Test



Recommended by CLSI and the CDC for the detection of carbapenemase production.

## Endpoint PCR



Endpoint PCR used to detect the presence or absence of genes required for antibiotic production.

## Antibiotic Profiling



VITEK used to analyze resistance to various antibiotic classes

Penicillins

Cephalosporins

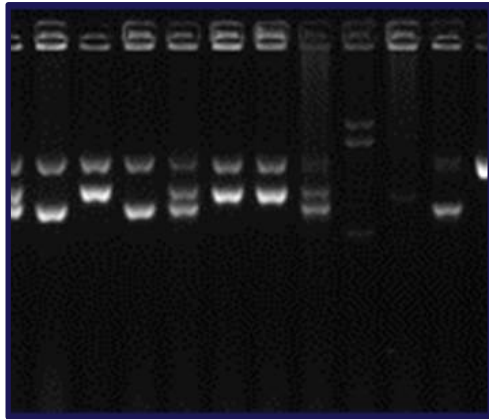
Carbapenems

Quinolones

Aminoglycosides

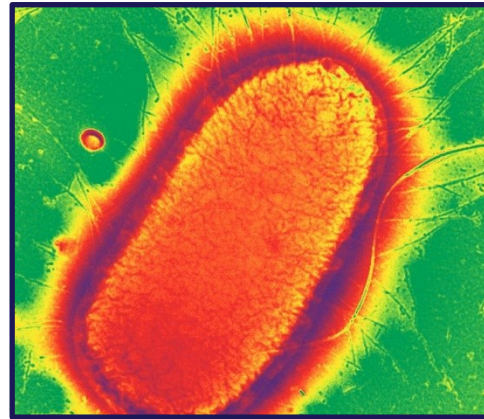
# Verification of virulence

## Endpoint PCR



Endpoint PCR used to detect the presence or absence of genes required for the production of toxins and other virulence factors.

## Toxin Production

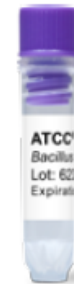


Detection of toxin production using an Enzyme Immunoassay (EIA)



# ATCC® Minis

BIG QUALITY  
in a **TINY TUBE**



actual size



Trusted ATCC Genuine Cultures® are offered in a convenient single-use “mini” format that saves you precious time and resources.

- Six pack of ready-to-use strains in glycerol stock
- Glass-free mini-cryovials with 2D barcode for easy storage
- Peel-off labels for fast and reliable recordkeeping

ATCC® Minis are authenticated and backed by ATCC polyphasic testing – ensuring the consistency and reliability that you have come to trust from ATCC Genuine Cultures®.



# ATCC® Minis USP QC organisms

2009: International USP <61> & US <62> Harmonization of US, Japan, EU Pharmacopeia<sup>1</sup>

Organism	ATCC® No.	USP 51 Antimicrobial	USP 61 Media	USP 62 Media
<i>Pseudomonas aeruginosa</i>	9027-MINI-PACK™	X	X	X
<i>Staphylococcus aureus</i>	6538-MINI-PACK™	X	X	X
<i>Candida albicans</i>	10231-MINI-PACK™	X	X	X
<i>Aspergillus brasiliensis</i>	16404-MINI-PACK™	X	X	
<i>Escherichia coli</i>	8739-MINI-PACK™	X		X
<i>Bacillus subtilis</i>	6633-MINI-PACK™		X	
<i>Salmonella enterica</i>	14028-MINI-PACK™			X
<i>Clostridium sporogenes</i>	11437-MINI-PACK™		X	

[www.atcc.org/Minis](http://www.atcc.org/Minis)

# ATCC® Minis general QC organisms

We are adding to the ATCC® Minis QC menu monthly  
Visit our website regularly for updates: [www.atcc.org](http://www.atcc.org)

Organism	ATCC® No.	Organism	ATCC® No.
<i>Escherichia coli</i>	25922-MINI-PACK™	<i>Escherichia coli</i>	11775-MINI-PACK™
<i>Staphylococcus aureus</i>	25923-MINI-PACK™	<i>Staphylococcus epidermidis</i>	12228-MINI-PACK™
<i>Pseudomonas aeruginosa</i>	27853-MINI-PACK™	<i>Clostridium sporogenes</i>	19404-MINI-PACK™
<i>Enterobacter aerogenes</i>	13048-MINI-PACK™	<i>Streptococcus pyogenes</i>	19615-MINI-PACK™
<i>Enterococcus faecalis</i>	29212-MINI-PACK™	<i>Burkholderia cepacia</i>	25416-MINI-PACK™
<i>Klebsiella pneumoniae</i>	10031-MINI-PACK™	<i>Streptococcus pneumoniae</i>	49619-MINI-PACK™
<i>Escherichia coli</i>	11229-MINI-PACK™		

[www.atcc.org/Minis](http://www.atcc.org/Minis)

# Features & benefits

## ATCC® Minis

- Glass-free, convenient, easy-to-use “Mini” format that saves you precious time and resources

## QC Strains

- The same trusted ATCC Genuine Cultures® you’ve come to know and trust

## Glycerol Stock

- The same ready-to-plate format many labs create in-house from ATCC strains

## 500 µL Tube

- Mini-cryovial takes less room than a standard cryovial
- ATCC® Minis Working Rack helps you keep them upright

## 96-well Format

- Smaller storage space needed
- ATCC® Minis Storage Box securely locks for safe keeping

## 6 Pack

- Convenient pack size allows you to plan ahead

## Peel-off Label

- Apply to lab notebooks, plates, and flasks for proper labelling and tracking

## 2D Barcode

- Easily ties into LIS/ LIMS for sample tracking

## -80°C Storage

- The same storage many labs use for their lab-created glycerol stocks

## Labeled Exp.

- Required by many labs, and helps manage inventory

# Mini conveniences

## ATCC® Minis

Authenticated

Gold standard

Ready-to-plate

Single-use

Glass-free

## ATCC Mini Accessories & Pack

Storage Box	A lock tight storage solution for your ATCC® Minis. The ATCC® Minis Storage box with locking lid allows you to store ATCC® Minis in a convenient 96-well format that saves you precious freezer space.
Rack	A working rack to keep all your ATCC® Minis upright and ready-to-use. The 96-well format and low profile design allows you to easily work with and identify each of your ATCC® Minis.
Cap Tool	ATCC® Minis' little helper. The Cap Tool allows for aseptic opening of ATCC® Minis by fitting into the cap. Give it a twist and you're ready-to-plate.
QC Pack	ATCC has conveniently bundled the USP recommended QC organism ATCC® Minis 6 packs, with an overall cost savings of over 10% versus buying each individually.



# Thank you!

Register for more webinars in the ATCC “*Excellence in Research*” webinar series at [www.atcc.org/webinars](http://www.atcc.org/webinars).

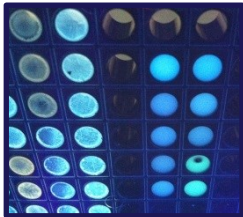


**April 2, 2015**

**10:00 AM, 3:00 PM EST**

James Clinton, Ph.D.

ATCC Transfection Reagents – Powerful Tools to Enable Genetic Manipulation



**May 21, 2015**

**10:00 AM, 3:00 PM EST**

Jodie Lee, M.S.

Seeing is Believing – Reporter-labeled Microbial Control Strains

**Thank you for joining today!**  
**Please send additional questions to [tech@atcc.org](mailto:tech@atcc.org)**