

# Innovating Microbial QC

How MicroQuant™ Transforms Speed,  
Accuracy, and Resource Efficiencies

Nilay Chakraborty, PhD, MBA  
BioNexus Foundation Principal Scientist  
and Head of Cryobiology  
ATCC



# Introduction to today's speaker



## **Nilay Chakraborty, PhD, MBA**

BioNexus Foundation Principal Scientist and Head of Cryobiology

ATCC

Dr. Chakraborty specializes in biopreservation and innovative product development. With an engineering background and degrees from Indian Institute of Engineering Science and Technology and University of North Carolina, he pioneered biopreservation and cell-based technologies that focus on innovative delivery formats. During his tenure at the Center for Engineering in Medicine at Harvard Medical School, Massachusetts General Hospital, and Shriners Burns Hospital, he continued research in biopreservation, delivery formats, and cell-based technologies. Previously a tenured Associate Professor at the University of Michigan, Dearborn, he now leads ATCC's efforts in preservation sciences and advanced strategic biological product development.

# About ATCC

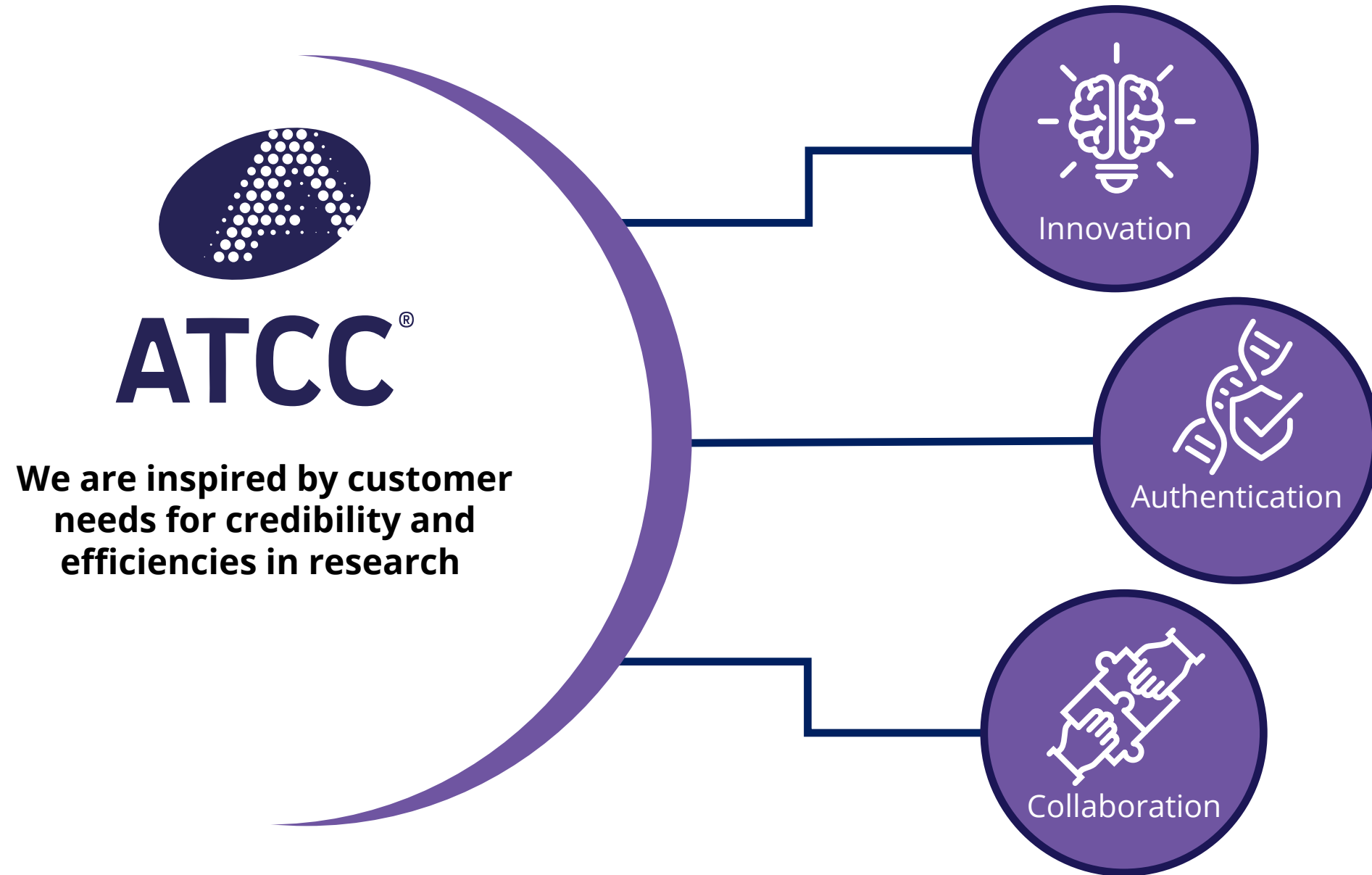


ATCC is a global leader in providing authenticated, high-quality biological resources and standards for industry, academia, and government.

- Founded in 1925
- Private, nonprofit, global biological resource center and standards organization
- Provides scientists with the biomaterials and resources:
  - 80,000+ microorganisms
  - 4,000+ cell lines
  - Standards & controls
  - Media, sera, and reagents
  - Advanced cell models
  - ATCC Genome Portal
    - 6,500 genomes
    - 3,000 transcriptomes
    - 500 exomes



# Enabling scientific progress for over 100 years





# Innovative products for life science research



Biologics production



Cell & gene therapy development



Vaccine development

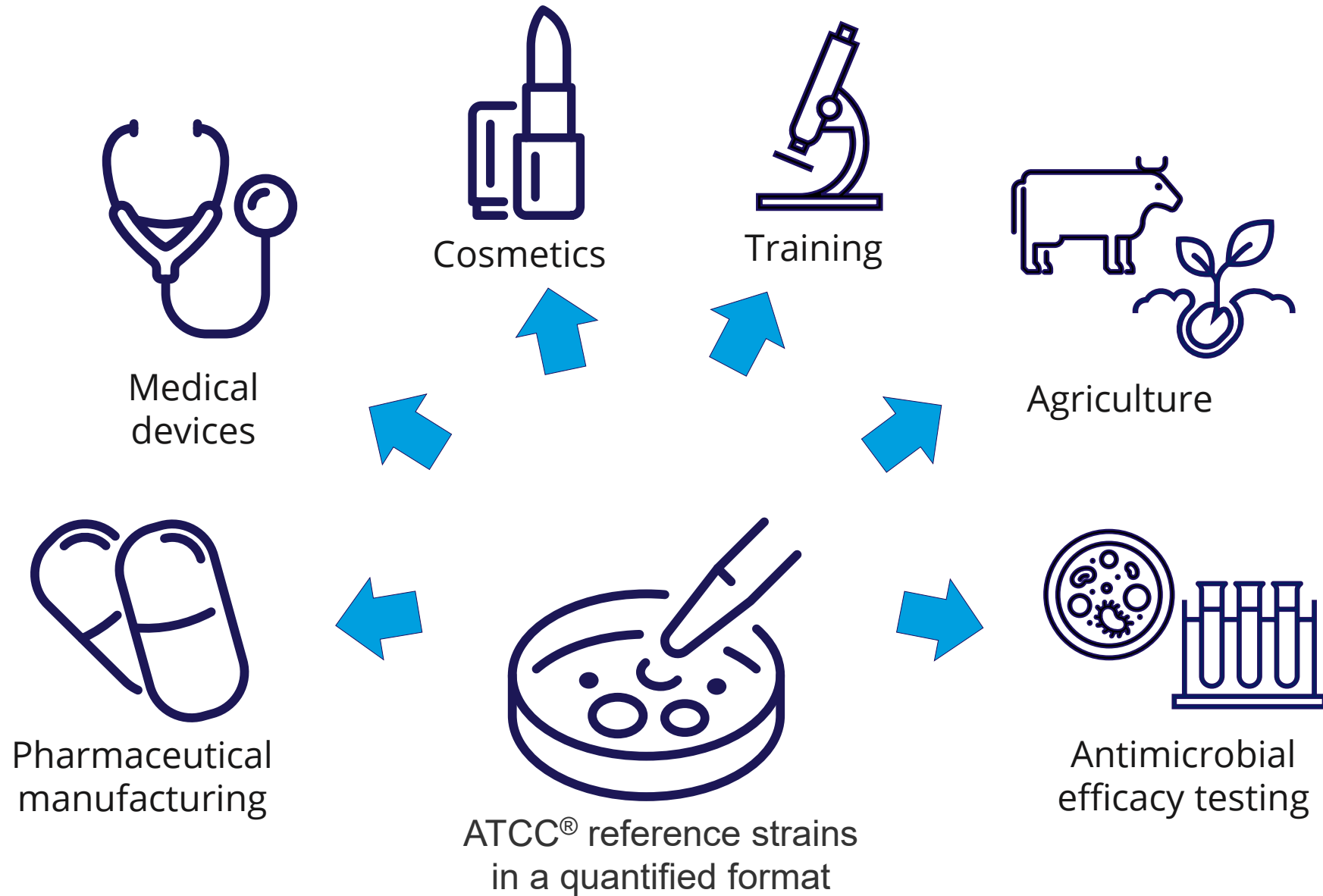


Quality & safety testing

## Comprehensive collection of advanced cell models and microbial reference materials that support biologics production

- Bioproduction cell lines and cell lines for enhanced virus production
- Analytical reference materials for residual host cell DNA testing
- Viral reference materials for evaluating the dose and potency of gene therapy products
- Purified polysaccharides and infectious disease strains for vaccine development
- Microbial quality control strains specified in standards and guidelines by organizations and regulatory agencies (USP, EP, ISO, FDA, CLSI, USDA, ASTM, AOAC, WHO)

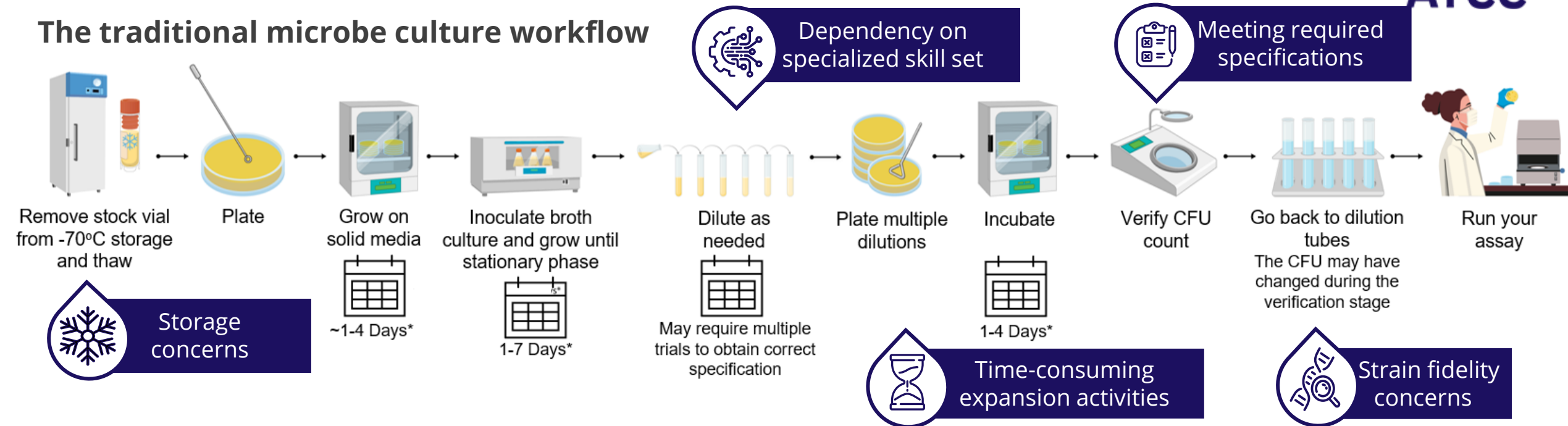
# Applications for quantified microbial control strains



# Challenges when using microbial reference strains



## The traditional microbe culture workflow





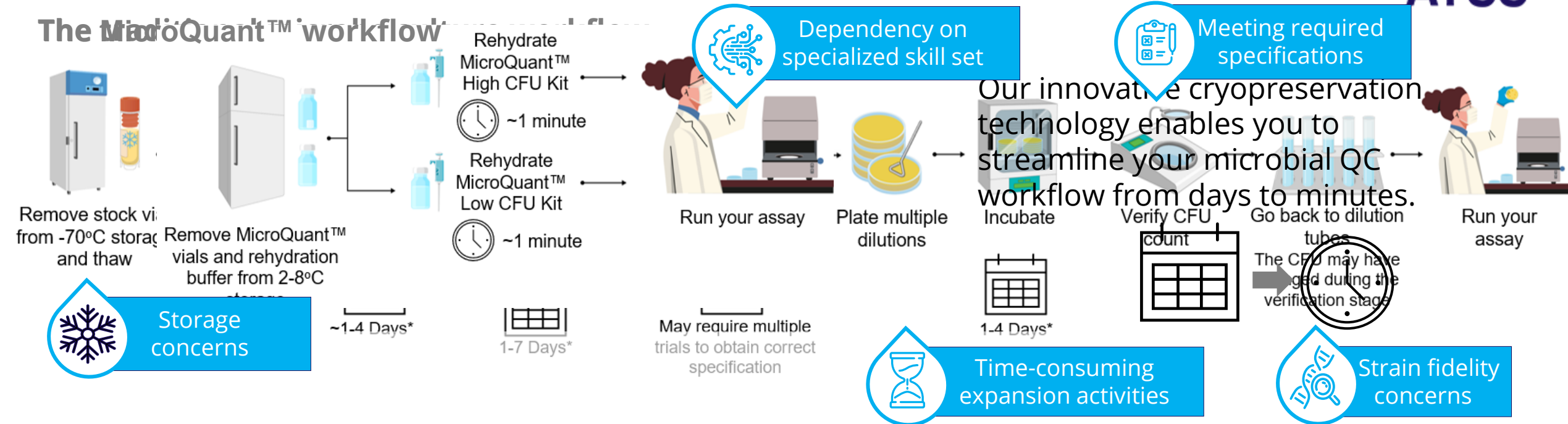
**MicroQuant™**  
by **ATCC®**



# Simplified workflow using MicroQuant™



## The MicroQuant™ workflow



# Meeting the challenge through cryobiology



ATCC developed an innovative preservation technique that delivers:

- A stable, pelleted format that rehydrates rapidly and uniformly
- Precise quantitation of biological materials
- Convenient refrigeration storage



# Introducing MicroQuant™ by ATCC®

Precision in every pellet, trust in every test



## A ready-to-use solution to help you streamline microbial QC testing

- Precisely quantitated in high-titer and low-titer pelleted formats
- Single-use format using an innovative, proprietary preservation technology
- Rapid, uniform rehydration in less than a minute
- Easy to store and ready to use anytime—no need to thaw
- Passage zero ATCC materials manufactured under ISO 17034

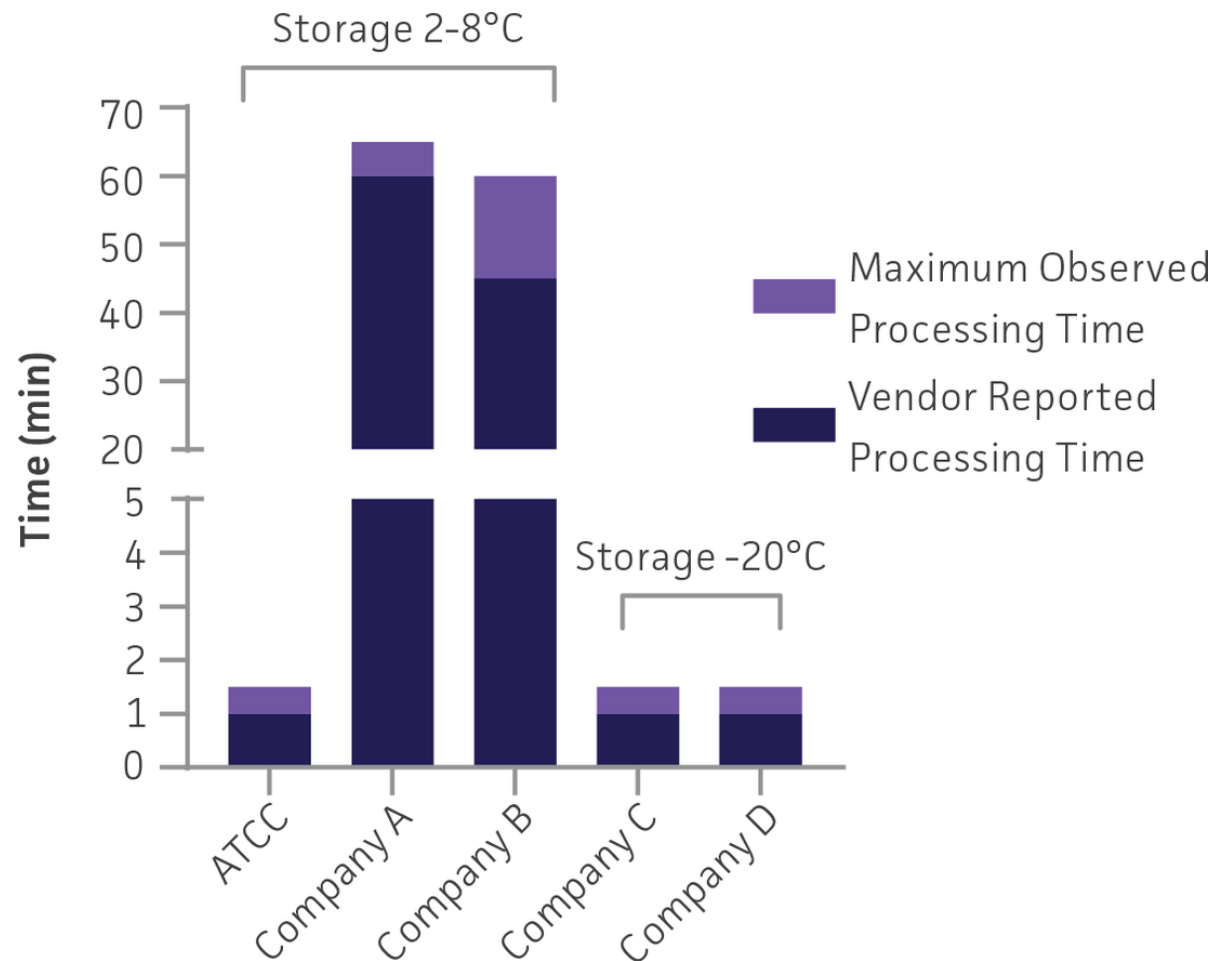






# MicroQuant™ solves the 'limited resources' problem

An easy-to-use control product with storage at convenient temperatures

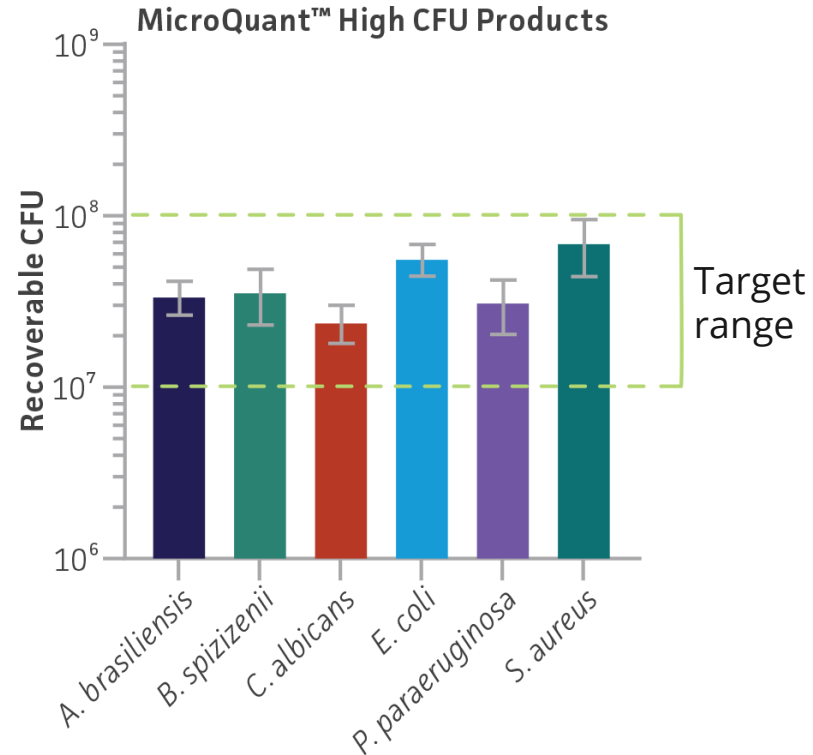
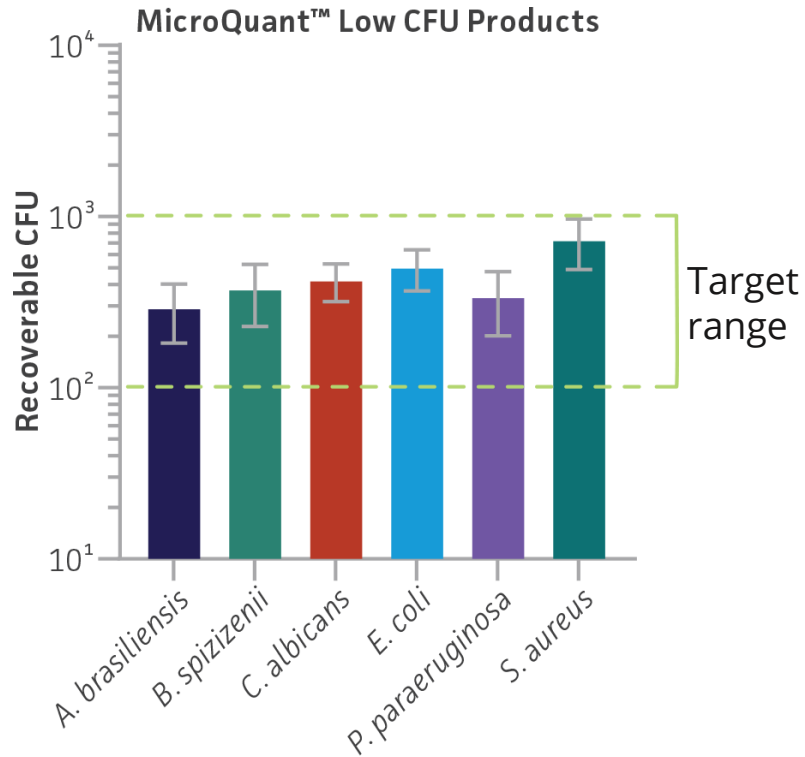


MicroQuant™ only requires **refrigerated storage.**

**It outperforms competitive products** stored by refrigeration and performs at least as well as products stored at -20°C.

# MicroQuant™ solves 'process complexity'

The value to teams who make internal control banks



## What does this mean for you?

- No culture management
- Reduced QC testing
- Reduced inventory management
- Enhanced consistency
- Faster workflows

*Each bar represents a minimum of 12 datapoints, error bars show standard deviation.*

# MicroQuant™: Precise and consistent quantitation

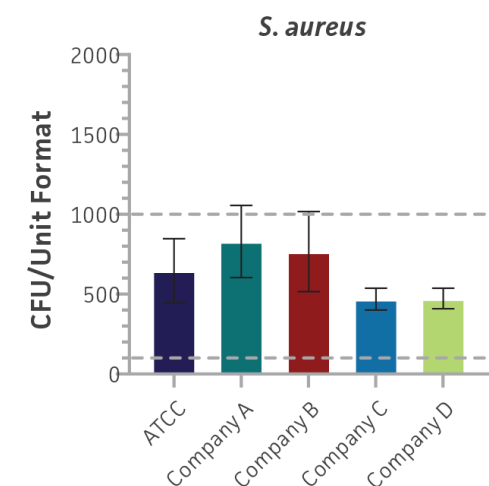
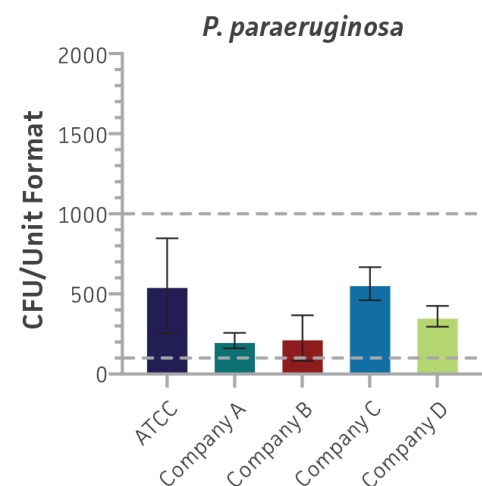
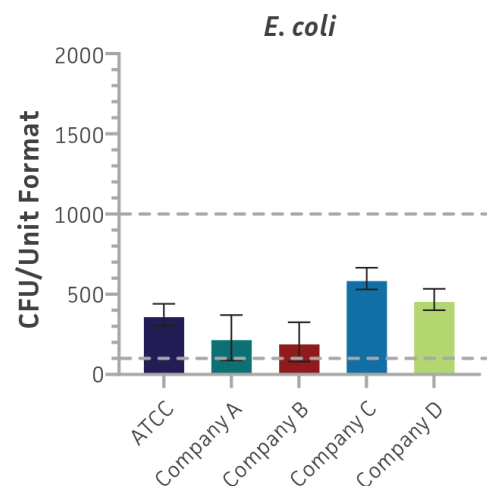
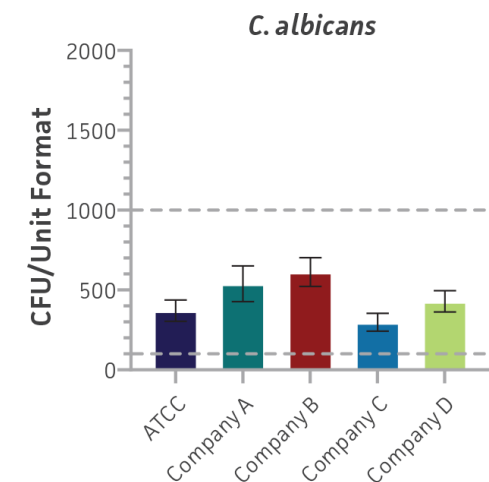
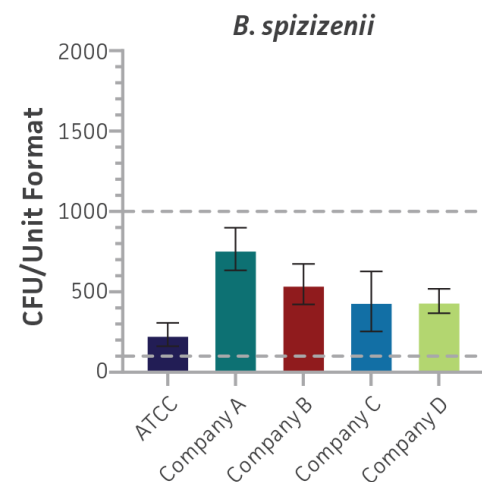
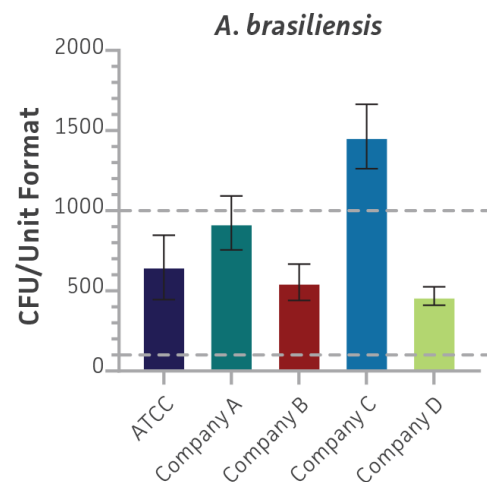
Precise and accurate low CFU pellets for bioburden testing



In benchmark tests of comparable low CFU products, MicroQuant™ performs as well as or better than other products on the market.

## Benchmarking for to 1,000 CFU

Each bar represents a minimum of 12 data points, error bars show standard deviation.





# MicroQuant™: Precise and consistent quantitation

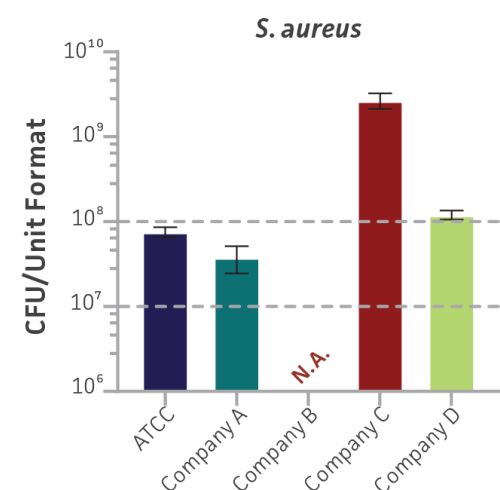
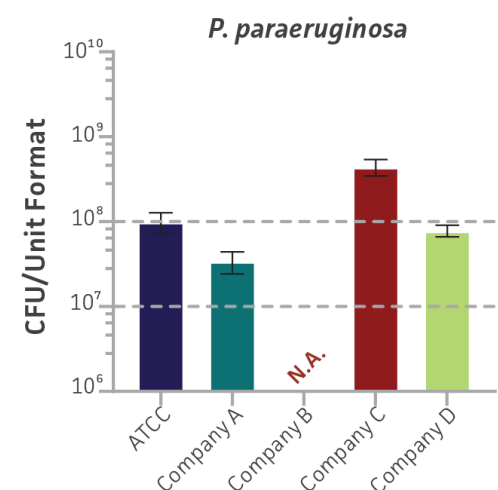
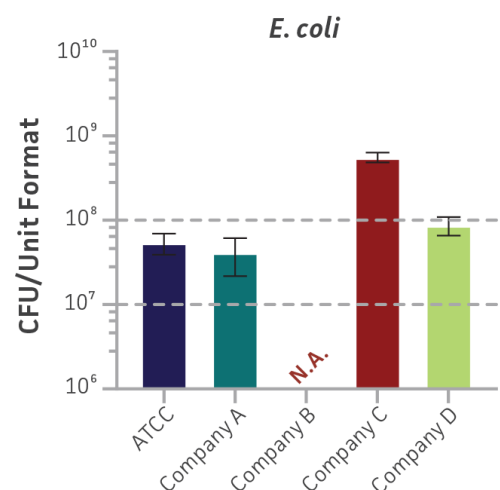
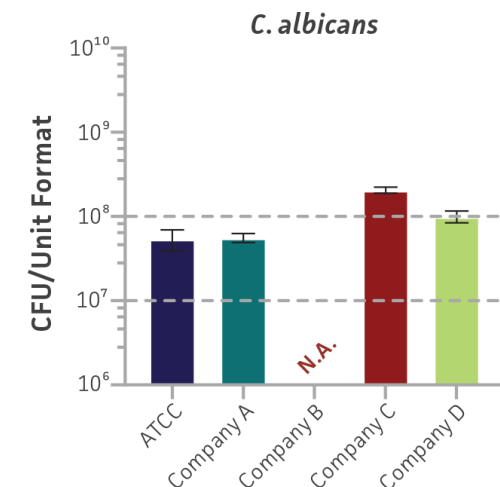
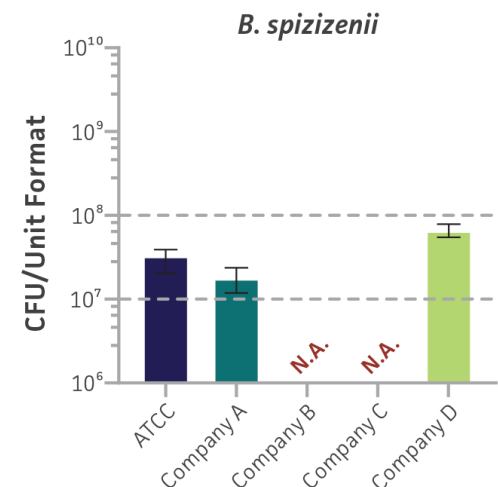
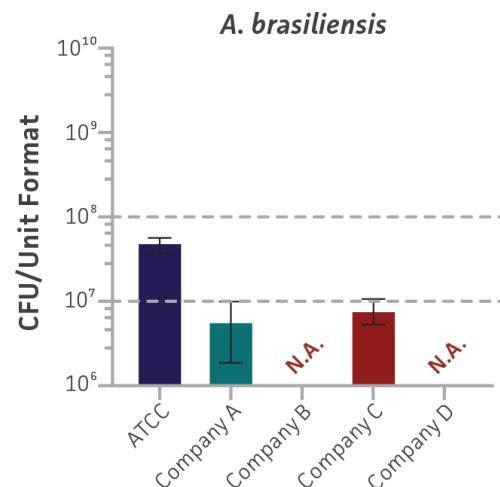
## Precise & accurate high CFU pellets for antimicrobial effectiveness testing



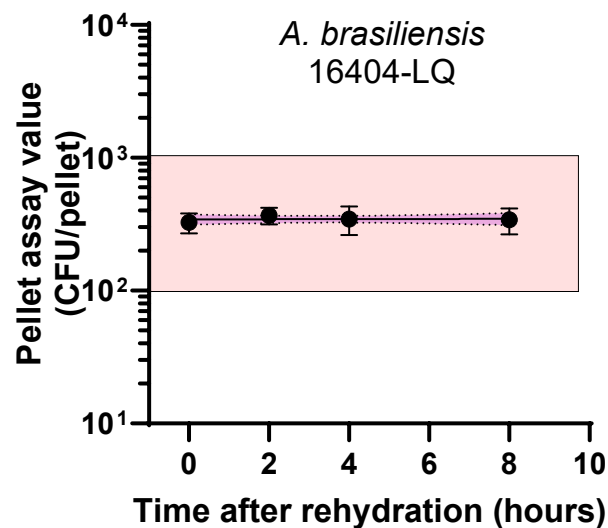
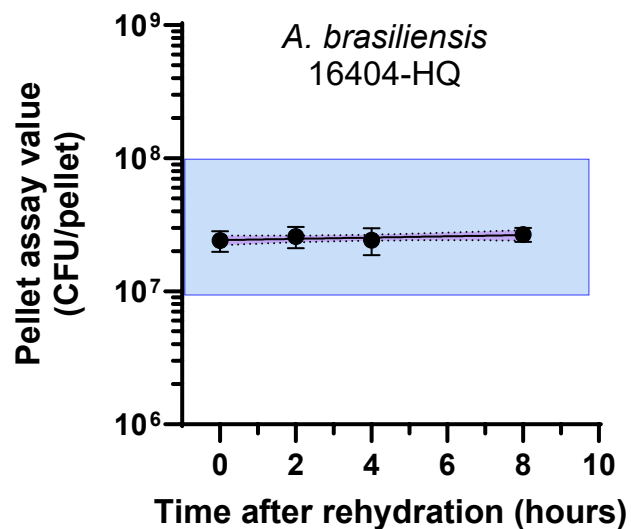
In benchmark tests of comparable high CFU products, MicroQuant™ performs as well as or better than other products on the market.

### Benchmarking for $10^7$ to $10^8$ CFU products

Each bar represents a minimum of 12 data points, error bars show standard deviation. N.A. = Not available at time of testing.



# Usability of MicroQuant™ products after rehydration



Organism	ATCC catalog	Usability (hours)	
		4°C	Room temp 22.3 ± 1.5°C
<i>A. brasiliensis</i>	16404-HQ-PACK™	8	8
	16404-LQ-PACK™	8	8
<i>B. spizizenii</i>	6633-HQ-PACK™	8	7
	6633-LQ-PACK™	8	8
<i>C. albicans</i>	10231-HQ-PACK™	8	8
	10231-LQ-PACK™	8	8
<i>E. coli</i>	8739-HQ-PACK™	8	8
	8739-LQ-PACK™	8	8
<i>P. paraeruginosa</i>	9027-HQ-PACK™	8	7
	9027-LQ-PACK™	8	8
<i>S. aureus</i>	6538-HQ-PACK™	8	8
	6538-LQ-PACK™	8	8
<i>C. sporogenes</i>	19404-HQ-PACK™	8	8
	19404-LQ-PACK™	8	8
<i>C. acnes</i>	11827-HQ-PACK™	8	8
	11827-LQ-PACK™	8	8

# Recovery of low CFU MicroQuant with selective media

Based on USP<62> guidelines for bioburden testing on selective media



Selective Medium for USP<62>	Test Strains / Target Organism	Recovery compared to TSA (%)
Violet Red Bile Glucose Agar	<i>E. coli</i> (ATCC <sup>®</sup> 8739-LQ-PACK <sup>™</sup> )	71 ± 19
	<i>P. paraeruginosa</i> (ATCC <sup>®</sup> 9027-LQ-PACK <sup>™</sup> )	68 ± 18
MacConkey Agar	<i>E. coli</i> (ATCC <sup>®</sup> 8739-LQ-PACK <sup>™</sup> )	88 ± 16
Xylose Lysine Deoxycholate Agar	<i>S. enterica</i> (ATCC <sup>®</sup> 14028-LQ-PACK <sup>™</sup> )	70 ± 14
Cetrimide Agar	<i>P. paraeruginosa</i> (ATCC <sup>®</sup> 9027-LQ-PACK <sup>™</sup> )	82 ± 11
	<i>E. coli</i> (ATCC <sup>®</sup> 8739-LQ-PACK <sup>™</sup> )	Negative Control
Mannitol Salt Agar	<i>S. aureus</i> (ATCC <sup>®</sup> 6538-LQ-PACK <sup>™</sup> )	83 ± 11
	<i>E. coli</i> (ATCC <sup>®</sup> 8739-LQ-PACK <sup>™</sup> )	Negative Control
Reinforced Medium for Clostridia	<i>C. sporogenes</i> (ATCC <sup>®</sup> 11437-LQ-PACK <sup>™</sup> )	100 ± 14
	<i>C. sporogenes</i> (ATCC <sup>®</sup> 19404-LQ-PACK <sup>™</sup> )	114 ± 13
Sabouraud Dextrose Agar	<i>C. albicans</i> (ATCC <sup>®</sup> 10231-LQ-PACK <sup>™</sup> )	107 ± 29

TSA-Tryptic soy agar

# How MicroQuant™ compares to similar products



Features	MicroQuant™ by ATCC®	Company A	Company B	Company C	Company D
Processing time ~1 minute	✓			✓	✓
2-8°C refrigeration storage	✓	✓	✓		
Manufactured under ISO 17034	✓	✓	✓	✓	✓
Includes all strains for USP <51> (High CFU format)	✓	✓*		✓*	✓*
Includes all strains for USP <61> (Low CFU format)	✓	✓	✓	✓	✓
Product kit includes rehydration buffer	✓	✓	✓	✓	
Sourced from ATCC®	✓	✓	✓		

\*Some strains were not available for purchase at the time of testing



# MicroQuant™ saves time and cost

## Versus preparing internal controls



Based on USP<62> guidelines for bioburden testing on selective media

**86%**

of labor time is saved when using MicroQuant™ versus preparing internal controls manually

**83%**

cost reduction is achieved using MicroQuant™ versus using internally prepared controls



# Pharmacopeias supported by MicroQuant™

## For microbial quality control testing



- Antimicrobial Effectiveness Testing (USP <51>, EP 5.1.3, JP 4.05 I)
- Test for *Burkholderia cepacia* Complex (USP <60>)
- Microbial Enumeration Tests (USP <61>, EP 2.6.12, JP 4.05 I)
- Tests for Specified Microorganisms (USP <62>, EP 2.6.13, JP 4.05 II)
- Sterility Test (USP <71>, EP 2.6.1, JP 4.06)
- Respiration-Based Microbiological Methods for Short-Life Products (USP <72>)
- Disinfectants and Antiseptics (USP <1072>)



# Summary for MicroQuant™



- Streamline microbial quality control testing
- Precisely quantitated
  - High-titer (HQ;  $10^7$  to  $10^8$  CFU per vial)
  - Low-titer (LQ; 100 to 1,000 CFU per vial)
- Single-use format for fast assay setup & minimal handling.
- Developed from traceable, original source materials provided at Passage zero
- Immediate rehydration at room temperature with an 8-hour usability window
- Stable storage at 2-8°C
- Manufactured under an ISO 17034–accredited process



# What MicroQuant™ products are available?

# Explore the MicroQuant™ portfolio

## Low CFU = 100 to 1,000 CFU/vial



ATCC® No.	Description
10231-LQ-PACK™	MicroQuant™ <i>Candida albicans</i> , low CFU
9027-LQ-PACK™	MicroQuant™ <i>Pseudomonas paraeruginosa</i> , low CFU
6538-LQ-PACK™	MicroQuant™ <i>Staphylococcus aureus</i> subsp. <i>aureus</i> , low CFU
16404-LQ-PACK™	MicroQuant™ <i>Aspergillus brasiliensis</i> , low CFU
8739-LQ-PACK™	MicroQuant™ <i>Escherichia coli</i> , low CFU
6633-LQ-PACK™	MicroQuant™ <i>Bacillus spizizenii</i> , low CFU
25416-LQ-PACK™	MicroQuant™ <i>Burkholderia cepacia</i> , Low CFU, Pack of 5
BAA-245-LQ-PACK™	MicroQuant™ <i>Burkholderia cenocepacia</i> , Low CFU, Pack of 5
BAA-247-LQ-PACK™	MicroQuant™ <i>Burkholderia multivorans</i> , Low CFU, Pack of 5
14028-LQ-PACK™	MicroQuant™ <i>Salmonella enterica</i> , Low CFU, Pack of 5
19404-LQ-PACK™	MicroQuant™ <i>Clostridium sporogenes</i> , Low CFU, Pack of 5
11437-LQ-PACK™	MicroQuant™ <i>Clostridium sporogenes</i> , Low CFU, Pack of 5
11827-LQ-PACK™	MicroQuant™ <i>Cutibacterium acnes</i> , Low CFU, Pack of 5
10106-LQ-PACK™	MicroQuant™ <i>Penicillium chrysogenum</i> , Low CFU, Pack of 5
11229-LQ-PACK	MicroQuant™ <i>Escherichia coli</i> , low CFU (Pack of 5)
15442-LQ-PACK	MicroQuant™ <i>Pseudomonas aeruginosa</i> , low CFU (Pack of 5)
19659-LQ-PACK	MicroQuant™ <i>Bacillus subtilis</i> , low CFU (Pack of 5)
11709-LQ-PACK	MicroQuant™ <i>Penicillium chrysogenum</i> , low CFU (Pack of 5)



# Explore the MicroQuant™ portfolio

## High CFU = $10^7$ to $10^8$ CFU/vial



ATCC® No.	Description
10231-HQ-PACK™	MicroQuant™ <i>Candida albicans</i> , high CFU
9027-HQ-PACK™	MicroQuant™ <i>Pseudomonas paraeruginosa</i> , high CFU
6538-HQ-PACK™	MicroQuant™ <i>Staphylococcus aureus</i> subsp. <i>aureus</i> , high CFU
16404-HQ-PACK™	MicroQuant™ <i>Aspergillus brasiliensis</i> , high CFU
8739-HQ-PACK™	MicroQuant™ <i>Escherichia coli</i> , high CFU
6633-HQ-PACK™	MicroQuant™ <i>Bacillus spizizenii</i> , high CFU
25416-HQ-PACK™	MicroQuant™ <i>Burkholderia cepacia</i> , High CFU, Pack of 5
BAA-245-HQ-PACK™	MicroQuant™ <i>Burkholderia cenocepacia</i> , High CFU, Pack of 5
BAA-247-HQ-PACK™	MicroQuant™ <i>Burkholderia multivorans</i> , High CFU, Pack of 5
14028-HQ-PACK™	MicroQuant™ <i>Salmonella enterica</i> , High CFU, Pack of 5
19404-HQ-PACK™	MicroQuant™ <i>Clostridium sporogenes</i> , High CFU, Pack of 5
10106-HQ-PACK™	MicroQuant™ <i>Penicillium chrysogenum</i> , High CFU, Pack of 5
11229-HQ-PACK	MicroQuant™ <i>Escherichia coli</i> , high CFU (Pack of 5)
15442-HQ-PACK	MicroQuant™ <i>Pseudomonas aeruginosa</i> , high CFU (Pack of 5)
19659-HQ-PACK	MicroQuant™ <i>Bacillus subtilis</i> , high CFU (Pack of 5)
11709-HQ-PACK	MicroQuant™ <i>Penicillium chrysogenum</i> , high CFU (Pack of 5)





# Explore the MicroQuant™ portfolio

- For U.S., 20% discount pricing built into the panels below
- For non-U.S., check with your ATCC Distributor
- [www.atcc.org/support/determine-your-distributor](http://www.atcc.org/support/determine-your-distributor)

Find the  
“Find Distributor”  
link in the  
webinar console



ATCC® No.	Product Name	Description
MQ-51™	MicroQuant™ Antimicrobial Effectiveness Panel	A panel comprising the five boxes of microbial challenge organisms cited in USP <51>.
MQ-61™	MicroQuant™ Microbial Examination of Nonsterile Products Panel	A panel comprising the five boxes of microbial challenge organisms cited in USP <61>.
MQ-60™	MicroQuant™ Microbial Examination of Nonsterile Products: Tests for Burkholderia Cepacia Complex Panel	A panel comprising the five boxes microbial challenge organisms cited in USP <60>.
MQ-62™	MicroQuant™ Microbial Examination of Nonsterile Products: Tests for Specified Microorganisms	A panel comprising the seven boxes of microbial challenge organisms cited in USP <62>.
MQ-71™	MicroQuant™ Sterility Tests Panel	A panel comprising the six boxes of microbial challenge organisms cited in USP <71>.



# More information about MicroQuant™

# Access more information



App. Note & ASM Microbe Conference poster sharing results shown within this webinar

ATCC

APPLICATION NOTE

### DEVELOPMENT AND PERFORMANCE EVALUATION OF MICROQUANT™ – QUANTITATIVE, RAPIDLY REHYDRATING REFERENCE MATERIALS FOR STREAMLINED MICROBIAL QUALITY CONTROL TESTING

Quinn Osgood, BSE; Jyoti K Jha, PhD; Rahul Tevatia, PhD; Cara Wilder, PhD; Nilay Chakraborty, PhD, MBA  
ATCC, Manassas VA

#### ABSTRACT

Microbial quality control testing is a cornerstone of pharmaceutical product development and essential for the safety, efficacy, and purity of biological products. To ensure the accuracy of these tests, pharmaceutical manufacturers need full confidence that their microbial controls are authenticated, low passage, and precisely quantitated. Leveraging innovative preservation technologies, ATCC has developed our best-in-class microbial quality control strains in a rapidly rehydrating pelleted format that delivers consistent quantitation. In this study, we demonstrate the performance of this novel suite of products in streamlining microbial quality control testing.

#### INTRODUCTION

Microbial contamination of pharmaceutical products is one of the leading causes of product recalls in the pharmaceutical industry.<sup>1,2</sup> These recalls can lead to extensive financial losses, regulatory consequences, brand erosion, and reputational damage.<sup>3,4</sup> More importantly, microbial contamination can significantly affect patient health and safety, potentially resulting in permanent disabilities or life-threatening conditions. To ensure product safety and integrity, pharmaceutical manufacturers are required by their respective government agencies to perform robust quality control procedures that, when properly followed, can help identify microbial contamination prior to product release.

Pharmaceutical manufacturers in the United States perform tests such as antimicrobial effectiveness testing, environmental monitoring, growth promotion testing, bioburden testing, sterility testing, and suitability testing as recommended by the United States Pharmacopeia (USP).<sup>5-8</sup> To ensure the accuracy of these tests, the regulatory guidance recommends the use of test strains that are maintained by reputable culture collections like ATCC. While ATCC has consistently met the needs for microbial quality control testing by providing authenticated reference materials, the laboratories performing microbial quality control testing often face infrastructure and resource limitation challenges when preparing control strains for testing.

For instance, many testing laboratories choose to develop and maintain their own internal control banks of reference materials. Creation and maintenance of the control banks requires complex expansion protocols and extensive laboratory space and equipment. These processes can be error prone. Precise quantitation of the reference materials is also a challenge. Several quality control tests require reference materials to be prepared in either a low-titer or high-titer format. Meeting these specifications is traditionally obtained by preparing a culture of the required strain, which can require a significant investment in time and resources. Additionally, accurate quantification and passage of the microbial strains require highly skilled microbiologists to be involved. All these factors can make the process of microbial quality control testing a costly, time-intensive, and often inconvenient affair.

AES-SUNDAY-5128  
June 22, 2025

## Precisely Quantitated Reference Materials for Microbial Quality Control Testing

Jyoti K Jha, PhD; Emmanuel Bofo-Asare, MS; Rahul Tevatia, PhD; Quinn Osgood, BSE; Fadima Samboal, MS; Joshua Yeroshsky, MS; Aqliah Kermani, MS; Viraj Patel, MS; Nilay Chakraborty, PhD, MBA  
ATCC, Manassas, VA 20110

#### Abstract

MicroQuant™ is ATCC's new product line of quantitated microbial reference materials designed for quality control in biopharmaceutical, cosmetic, food safety, and wastewater testing. This novel product suite was developed using an innovative cryopreservation technology to create pre-quantitated microbial strains in a stable pelleted format that delivers consistent and reproducible results, easy workflows, and quick turnaround times. MicroQuant™ meets ISO 17034 standards and supports guidelines from organizations like USP, EP, and JP. Unlike other products, MicroQuant™ controls are made from original ATCC® strains, ensuring better strain identity and authentication. They provide consistent quantitation, instant rehydration at room temperature, and stable storage at 2-8°C. In this study, we highlight the stability of MicroQuant™ products during transport and rehydration, and we showcase how they compare to similar products with regard to processing time, storage, and stability.

#### Schematic use of MicroQuant™

**MicroQuant™ HQ/LQ workflow**

**Figure 1: MicroQuant™ assay workflow.** MicroQuant™ products are supplied as a kit, each kit contains five pellets of a specific organism (either HQ or LQ titer) and five vials of the respective rehydration buffer.

#### Application of MicroQuant™ high (HQ) and low (LQ) titer pellets

**Table 1: Application of MicroQuant™ products**

Species	ATCC® No.	Format	Commercial usage	Other industrial uses
<i>Aspergillus brasiliensis</i>	16404-HQ-PACK™	Available	USP <61> (107) [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Bacillus subtilis</i>	16810-HQ-PACK™	Available	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Candida albicans</i>	10231-HQ-PACK™	Available	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Enterobacter coli</i>	12221-HQ-PACK™	Available	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Escherichia coli</i>	8739-HQ-PACK™	Available	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Pseudomonas aeruginosa</i>	9027-HQ-PACK™	Available	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Staphylococcus aureus</i>	6358-HQ-PACK™	Available	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Burkholderia cepacia</i>	11427-HQ-PACK™	Q1-2025	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Burkholderia multivorans</i>	11428-HQ-PACK™	Q1-2025	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Clostridium sporogenes</i>	11429-HQ-PACK™	Q1-2025	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Clostridium sporogenes</i>	11430-HQ-PACK™	Q1-2025	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing
<i>Serratia enteritidis</i>	14028-HQ-PACK™	Q1-2025	USP <61> (71) [EP 2.2] [EP 2.6.13] [JP 10]	Food, media, QC, and pharma testing

#### Quantitation of MicroQuant™ production batches

**Figure 2: Analysis of MicroQuant™ colony forming units (CFU) per pellet from production batches.** Assay property values for (A) MicroQuant™ HQ and (B) MicroQuant™ LQ products. The data were obtained from nine random pellets per product with two technical replicates; the assay was performed by three independent analysts. Mean CFU/pellet and minimum and maximum CFU/pellet values are shown here. The blue and red shaded regions indicate the product specification for the MicroQuant™ HQ and LQ, respectively. CFU of the products was determined as described in Figure 1.

#### Transportation stability of ATCC® MicroQuant™

**Figure 3: Comparing the stability of MicroQuant™ products before and after transportation.** Assay property values for the pellets of (A) MicroQuant™ HQ and (B) MicroQuant™ LQ products. To assess transportation stability, MicroQuant™ kits were shipped at 2-8°C temperatures. Products were assayed to determine CFU/pellet before and after shipping. The average time during transportation was about 4 days. Control: CFU/pellet of MicroQuant™ HQ or LQ pellets before shipping. RT: CFU/pellet of pellets retrieved after room temperature transportation (Average temperature obtained from logger was 24.5°C). 4°C: CFU of pellets retrieved after 4°C transportation (Average temperature obtained from logger was 4.2°C). ns: not significant. The Y-axis indicates the CFU/pellet values of the product, and the X-axis indicates shipping conditions. A total of 9 random pellets and two technical replicates were used for CFU/pellet determination. The blue and red boxes indicate the product specification for the MicroQuant™ HQ and LQ, respectively.

#### Short-term stability (usability) of ATCC® MicroQuant™

**Figure 4: Short-term stability of MicroQuant™ after rehydration.** Assay property values for the pellets of (A) MicroQuant™ HQ and (B) MicroQuant™ LQ products were resuspended in rehydration buffer and then immediately diluted and plated (HQ) or plated (LQ) for "0 hour" reading. The remaining amounts of resuspended samples were stored at 4°C. Samples were removed from storage at specified intervals and were then diluted and plated (HQ) or directly plated (LQ). Plates were incubated and CFUs were counted following ATCC's guidelines. The blue and red boxes indicate product specification for the MicroQuant™ HQ and LQ, respectively. Above each panel, product number, and the microorganism names are included.

#### Comparison of processing time and storage condition

**Figure 5: Comparison of processing time and storage temperature of ATCC's MicroQuant™ with other products.** The processing time for MicroQuant™ was compared with the processing time for the formats available from Company A-D in the market. The storage temperature of ATCC, Company A, and Company B products is 4°C, whereas the storage temperature of Company C and Company D products is -20°C.

#### MicroQuant™ stability compared to similar products

**Table 2: Shelf-life comparison of MicroQuant™ HQ with products from other companies**

Organism	ATCC® No.	ATCC	Company A	Company B	Company C	Company D
<i>A. brasiliensis</i>	16404-HQ-PACK™	>1	<2	NA	2	NA
<i>B. subtilis</i>	16810-HQ-PACK™	>1	<2	NA	NA	<2
<i>C. albicans</i>	10231-HQ-PACK™	>1	<2	NA	2	<2
<i>E. coli</i>	8739-HQ-PACK™	>1	2	NA	2	<2
<i>P. aeruginosa</i>	9027-HQ-PACK™	>1	<2	NA	<2	<2
<i>S. aureus</i>	6358-HQ-PACK™	>1	<2	NA	<2	<2
Product storage		4°C				-20°C

**Table 3: Shelf-life comparison of MicroQuant™ LQ with products from other companies**

Organism	ATCC® No.	ATCC	Company A	Company B	Company C	Company D
<i>A. brasiliensis</i>	16404-LQ-PACK™	>1	<2	>1	<2	<1
<i>B. subtilis</i>	16810-LQ-PACK™	>1	<2	>1	2	<2
<i>C. albicans</i>	10231-LQ-PACK™	>1	<2	>1	<2	<2
<i>E. coli</i>	8739-LQ-PACK™	>1	<2	>1	<2	<2
<i>P. aeruginosa</i>	9027-LQ-PACK™	>1	<2	>1	2	<2
<i>S. aureus</i>	6358-LQ-PACK™	>1	<2	>1	<2	<2
Product storage		4°C				-20°C

NA - Products were not available from the companies at the time the study was conducted.  
\*Equivalent products available from Company C and D, respectively.  
Company A-D are products from other companies currently available in the market.  
ATCC's MicroQuant™ shelf-life is based on a combination of real-time stability studies, accelerated stability studies, and stability modeling.

#### Conclusions

ATCC® MicroQuant™ is an innovative product suite designed to streamline microbial quality control testing.

- Precisely quantitated in high-titer (HQ, 10<sup>10</sup> to 10<sup>11</sup> CFU per pellet) and low-titer (LQ, 100 to 1,000 CFU per pellet) formats.
- Single-use format enables fast assay setup and minimal handling.
- Immediate rehydration at room temperature with 8 hours of usability window.
- Stable storage at 2-8°C.
- Manufactured under an ISO 17034-accredited process.
- Developed from traceable, original source materials (Passage zero).

ATCC | Credible leads to Incredible®

Page 1

ATCC 10801 University Boulevard, Manassas, Virginia 20110-2209 Phone: 800.638.6597 Email: sales@atcc.org Web: www.atcc.org

© 2025 American Type Culture Collection. ATCC product identifiers marked with the TM symbol are trademarks owned by the American Type Culture Collection.

29

© 2026 American Type Culture Collection. ATCC product identifiers marked with the TM symbol are trademarks owned by the American Type Culture Collection.



# Access more information

Whitepaper with end-user feedback



YouTube videos

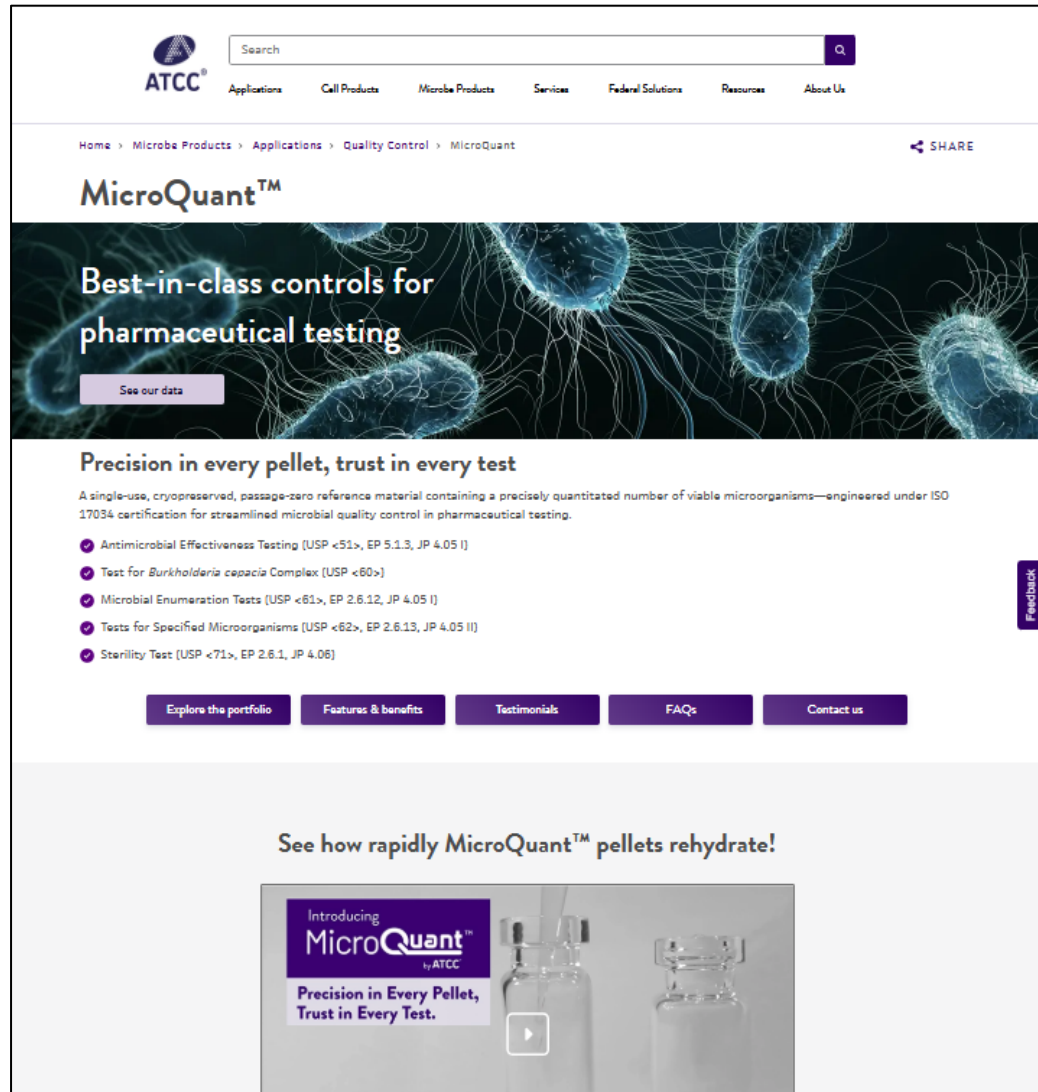


## MicroQuant: A Framework of Innovation

330K views • 9 months ago



# Access more information about MicroQuant™



Visit [www.atcc.org/microquant](http://www.atcc.org/microquant)

to view more information including:

- Application Note
- Whitepaper
- Conference posters
- Conference presentations
- Collaboration poster
- Flyers
- Brochures
- Instructions for use
- And more

For non-US, find your distributor:

[www.atcc.org/support/determine-your-distributor](http://www.atcc.org/support/determine-your-distributor)