



A Unified Product Structure for Innovative Product Development and Manufacturing of Reference Biomaterials in a Regulated Environment: The MicroQuant™ Case Study

Rahul Tevatia, PhD; Jyoti Jha, PhD; Quinn Osgood, BSE; Chad Madaloni, MBA; Nilay Chakraborty, PhD MBA
ATCC, 10801 University Blvd, Manassas, Virginia, 20110

TAP HERE TO
RETURN TO
KIOSK MENU

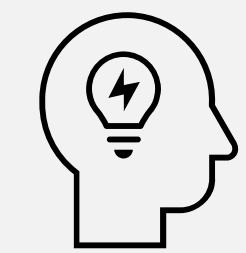
Introduction

Methods

Results

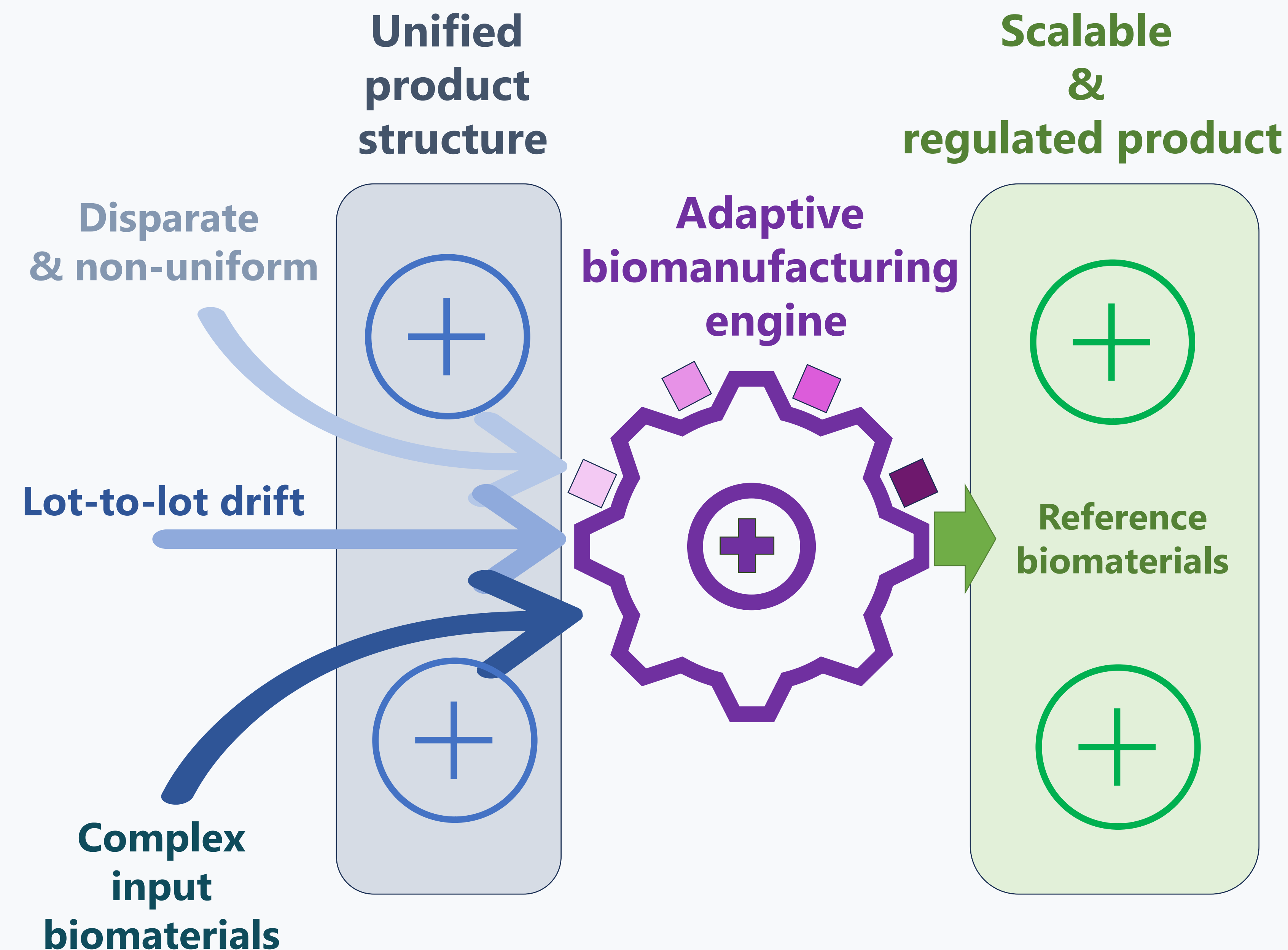
Discussion

Introduction



Problem:

- Translating novel biomaterials from R&D to regulated manufacturing is **slow and risky** 🕒
- Misalignment between innovation, development, manufacturing, and quality leads to:
 - Process inefficiencies
 - Variability and rework
 - Delayed timelines and regulatory risk



Rahul Tevatia, PhD

I am a Senior Scientist in the Cryobiology Department at ATCC. I have PhD in Chemical and Biomolecular Engineering and over 10 years of experience in solving complex problems in regulated biotech development and biopharmaceutical manufacturing.





A Unified Product Structure for Innovative Product Development and Manufacturing of Reference Biomaterials in a Regulated Environment: The MicroQuant™ Case Study

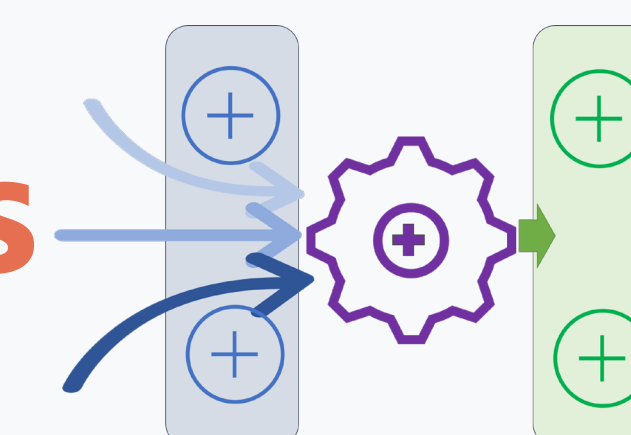
Rahul Tevatia, PhD; Jyoti Jha, PhD; Quinn Osgood, BSE; Chad Madaloni, MBA; Nilay Chakraborty, PhD MBA
ATCC, 10801 University Blvd, Manassas, Virginia, 20110

TAP HERE TO
RETURN TO
KIOSK MENU

Introduction

Methods

Materials & Methods



Results

Discussion

Quality-by-design control strategy

Critical quality attribute (CQA) and critical process parameter (CPP) mapping
Aligned with ICH Q8(R2), Q9, Q10, and Q11 as well as ISO 17034, ISO/IEC 17025, and ISO 9001

CONTROL STRATEGY

In-process monitoring · Specification limits · Risk-based decision gates · Deviation management · Change control

CQAs

Stage 1

- Organism identity (ATCC® Seed-stock: Passage zero)
- Supplier-certified traceability
- Inventory management

Stage 2

- Sterility of growth-medium and buffers
- Target colony forming units (CFU)
- No contamination

Stage 3

- Pellet integrity index
- Water activity ($AW < 0.2$)
- CFU/pellet is in intended use

Stage 4

- CFU/vial release criterion
- Strain identity confirmed
- Pellet specification met

Stage 5

- Certified property value
- Shelf-life statement
- Identity statement

Stage 6

- Temperature-controlled transit
- Customer feedback
- Customer complaints assessment

Material procurement

Culture & preparation

Lyo-casting & packaging

Quality control

Value assignment (CoA)

Distribution
Customer relation management (CRM)

CPPs

Stage 1

- Seed-stock storage (Liquid N₂ vapor phase)
- Growth-media composition
- Lyo-formulation
- Expiry / usable window

Stage 2

- Incubation temperature & time
- Aseptic handling / dispensing

Stage 3

- Lyophilization profile
- Inert gas purge & fill
- Cold-chain storage (4°C)

Stage 4

- Incubation conditions
- MALDI-TOF / 16s rRNA / ITS
- Plate count enumeration method

Stage 5

- Statistical analysis (GUM)
- Values reported: Mean, St dev., 95% CV
- Identity confirmation
- CoA issuance

Stage 6

- ISTA-aligned pack configuration
- Domestic: 4°C packaging overnight shipping
- International: dry ice shipment

QMS DOCUMENTATION

MasterControl SOPs / Work Instructions · Batch Records · Training & competence records · Calibration (IOQ) · CAPA / deviations · ISO 17034 compliance

CQA – A physical, chemical, biological, or microbiological property or characteristic that should be within an appropriate limit, range, or distribution to ensure the desired product quality. (ICH Q8(R2))

CPP – A process parameter whose variability has an impact on a CQA and therefore should be monitored or controlled to ensure the process produces the desired product quality. (ICH Q8(R2))

Process bands represents the validated process of MicroQuant™ production (ATCC®, Germantown, MD) and QC-value assignment-distribution (ATCC®, Manassas, VA)



A Unified Product Structure for Innovative Product Development and Manufacturing of Reference Biomaterials in a Regulated Environment: The MicroQuant™ Case Study

Rahul Tevatia, PhD; Jyoti Jha, PhD; Quinn Osgood, BSE; Chad Madaloni, MBA; Nilay Chakraborty, PhD MBA
ATCC, 10801 University Blvd, Manassas, Virginia, 20110

TAP HERE TO RETURN TO KIOSK MENU

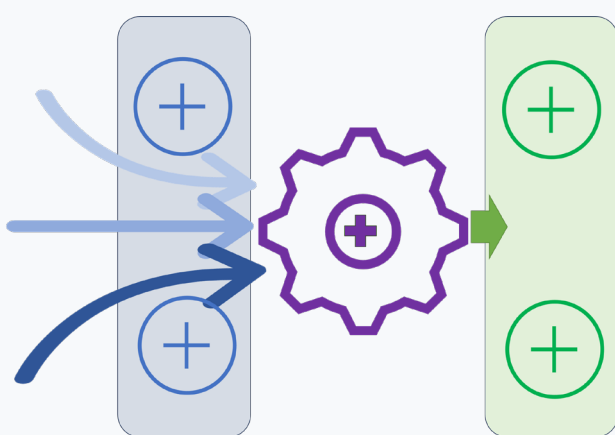
Introduction

Methods

Results

Results

Discussion

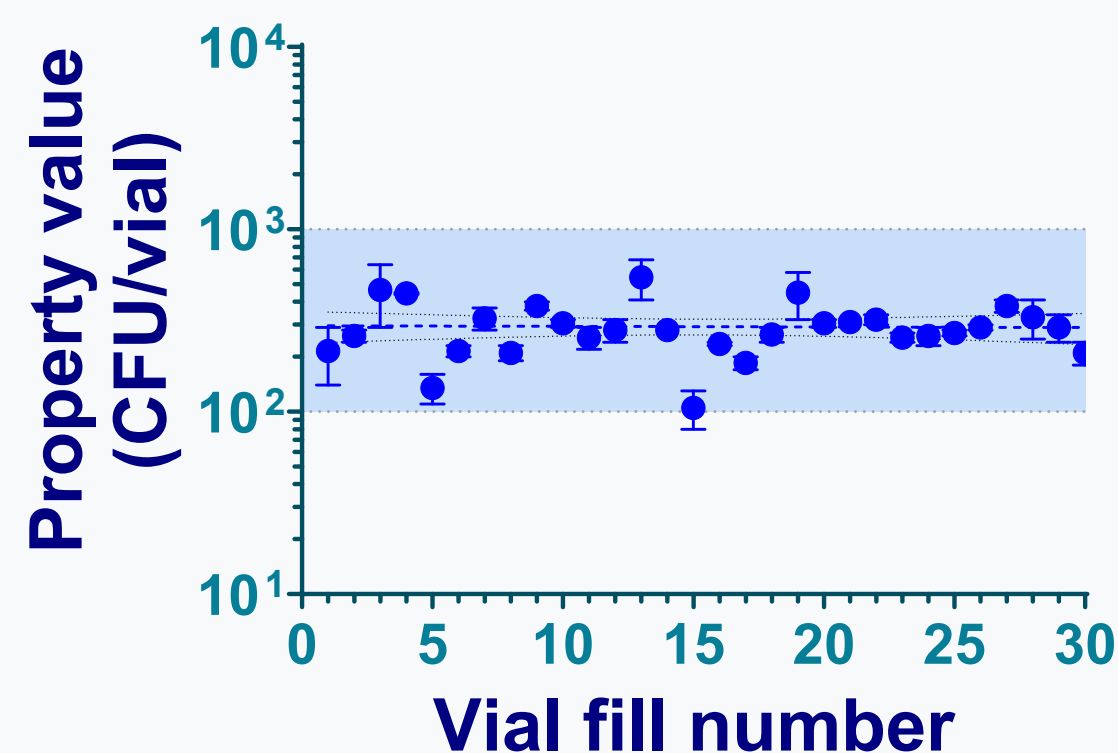


ATCC® MicroQuant™ (ATCC® 16404-LQ-PACK™)

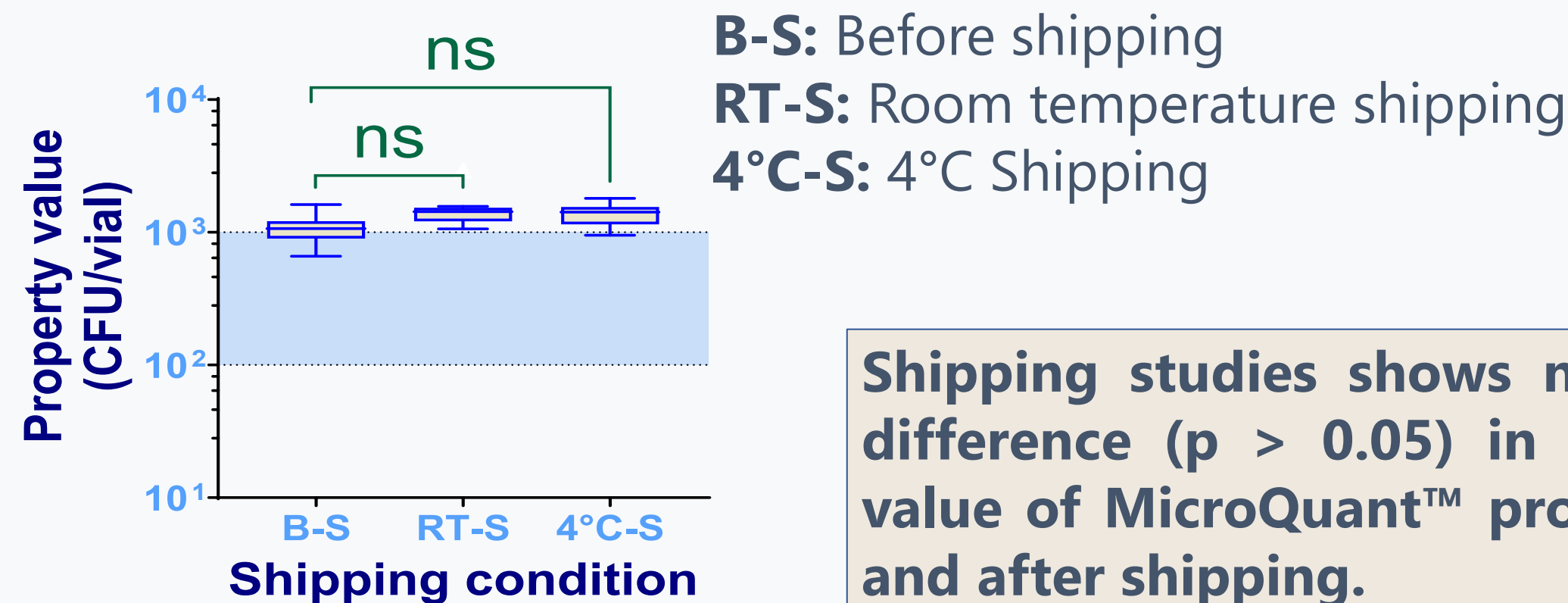
(A) Homogeneity

Random Vial Filling trend

Parameter	Best fit values	95% CI
Slope	-5.077	-11.82 to 1.66
Y-intercept (CFU/vial)	756	656 to 856
R ² value	0.0456	
Deviation from zero	Not significant (p > 0.05)	

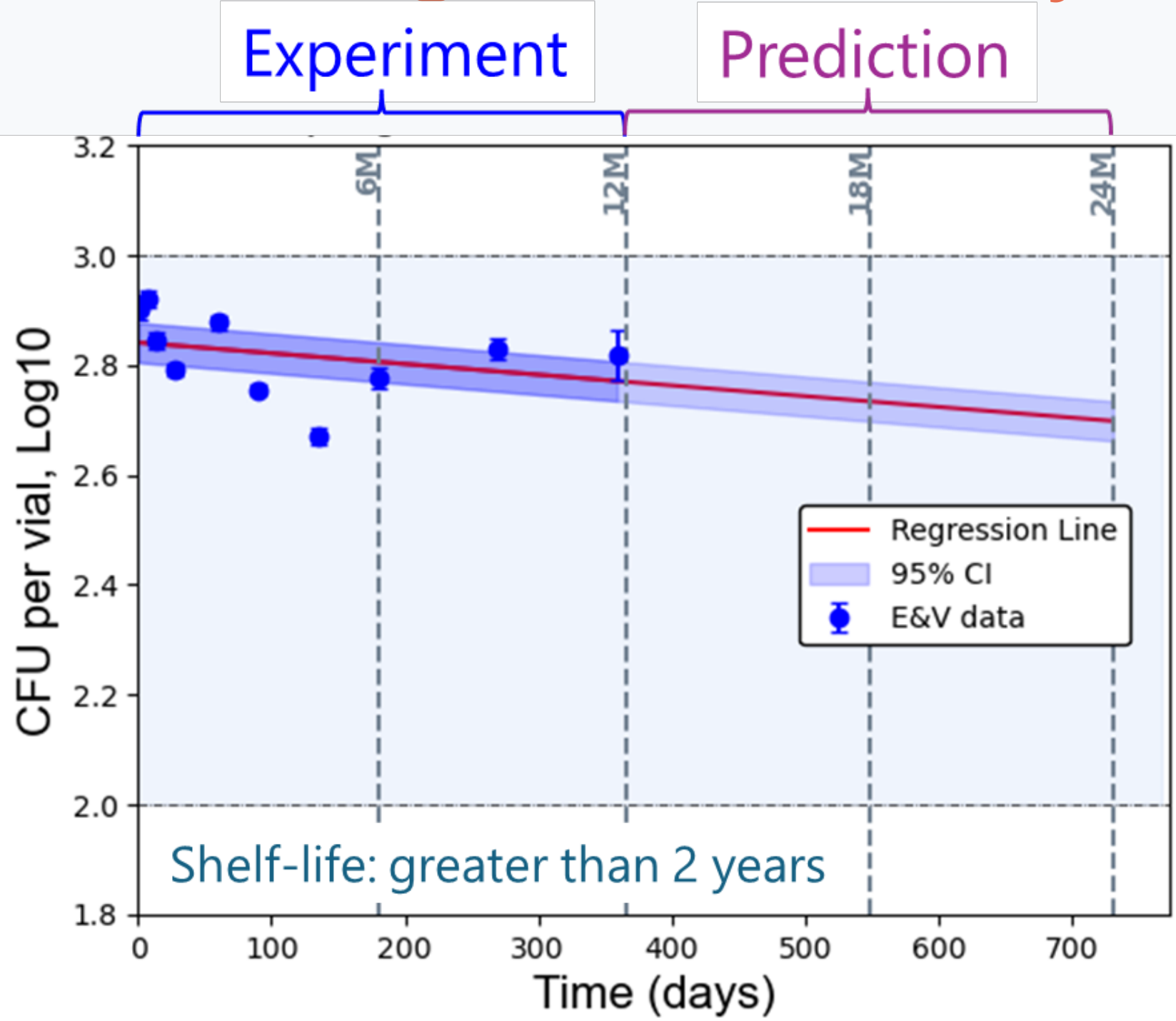


(C) Transport stability



Shipping studies shows no significant difference (p > 0.05) in the property value of MicroQuant™ products before and after shipping.

(B) Long-term stability



Storage temperature: 4±2°C
Testing temperature: 22.5±2.5°C (room temp.)

ISTA aligned pack configuration



(D) Scalable Model

Replicable blueprint for ISO 17034-compliant RM development and manufacturing

	1. Feasibility Scope & Design	2. Execute & Verify Characterize batch	3. Production Scale & Release	4. Monitor Post release
Homogeneity Between / within units	Not Applicable	✓	✓	✓
Stability Long-term / shelf-life	Not Applicable	✓	✓	✓
Transport Domestic & International	Not Applicable	✓	Not Applicable	✓
Use stability End-user stability	Not Applicable	✓	✓	Not Applicable
Characterization Property value & traceability	✓	✓	✓	✓
Uncertainty GUM propagation	Not Applicable	✓	✓	✓



A Unified Product Structure for Innovative Product Development and Manufacturing of Reference Biomaterials in a Regulated Environment: The MicroQuant™ Case Study

Rahul Tevatia, PhD; Jyoti Jha, PhD; Quinn Osgood, BSE; Chad Madaloni, MBA; Nilay Chakraborty, PhD MBA
ATCC, 10801 University Blvd, Manassas, Virginia, 20110

TAP HERE TO
RETURN TO
KIOSK MENU

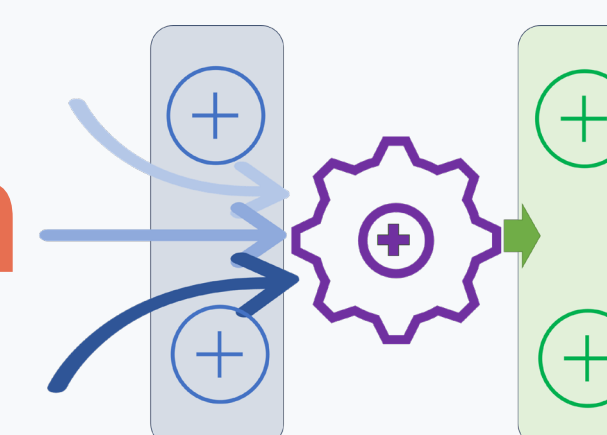
Introduction

Methods

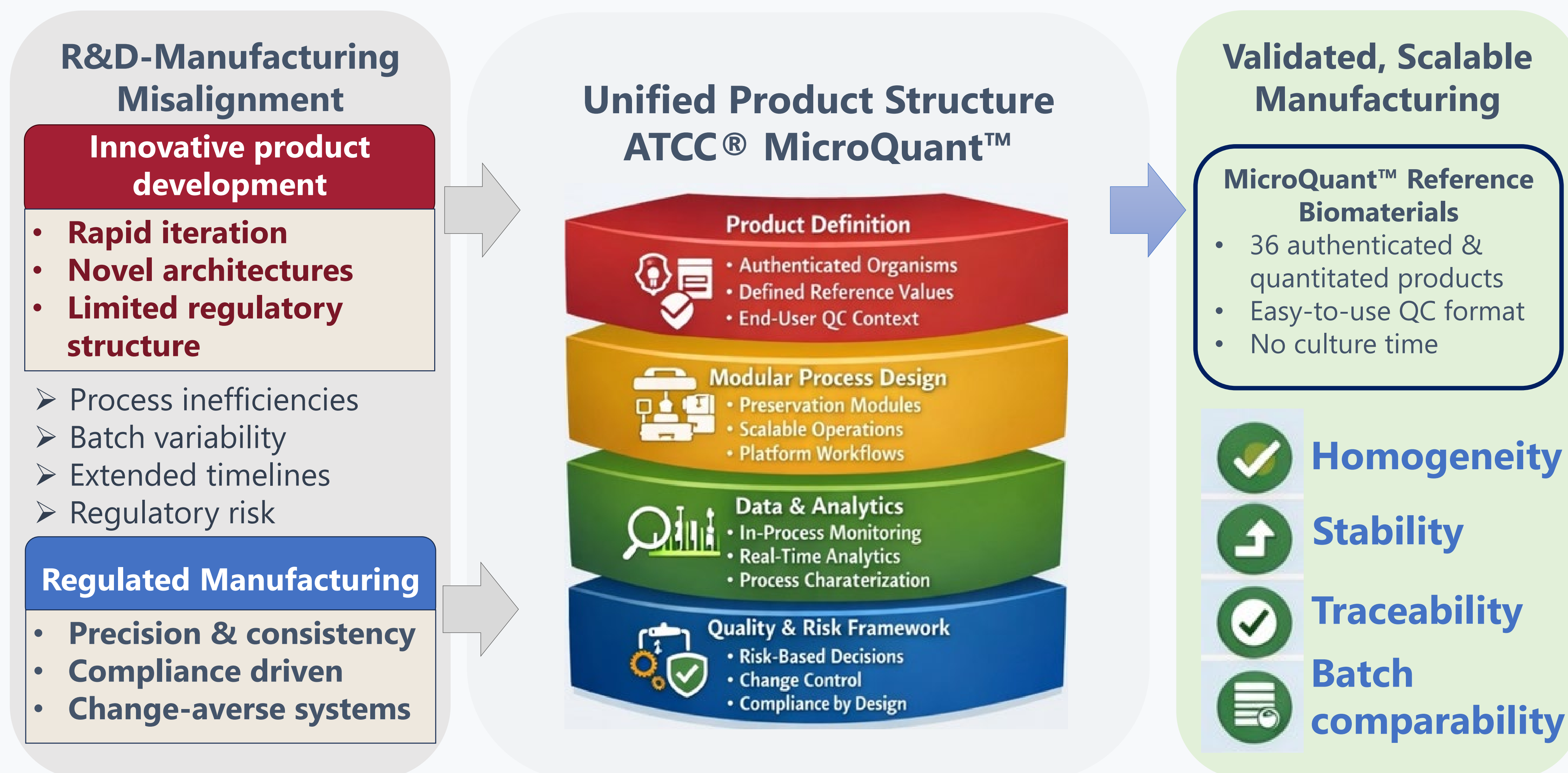
Discussion

Results

Discussion



(A) Unified product structure enabling translation of novel biomaterials into regulated manufacturing



A cohesive product structure bridges innovation and compliance, enabling scalable, accredited production of next-generation reference biomaterials.

(B) R&D to Manufacturing: Process and Quality Maturation

