

Enable Your Drug Development with cGMP Cell Banking and Storage Services



Shalmica Jackson, PhD Marketing Segment Manager, ATCC

Credible Leads to Incredible™



About ATCC



- World's largest, most diverse biological materials and information resource for microbes – the "gold standard"
- Innovative R&D company featuring gene editing, microbiome, NGS, advanced models
- cGMP cell banking and biorepository services

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



Overview

- The origin of GMP
- What is cGMP and why is it important?
- cGMP regulations and drug manufacturing
- Biologics manufacturing and cell banking
- ATCC's cGMP cell banking and storage services



ATCC



The Origin of Good Manufacturing Practices (GMP)





What Are Current Good Manufacturing Practices (cGMP)?

- Requirements for methods, facilities, and controls used in manufacturing
- Ensures that products meet specifications and Quality/Regulatory standards
- Covers requirements for Quality Control and Production



ATCC°



5

Why Is cGMP important?

- Public and consumer safety
- Minimizes risks
- Ensures safety, efficacy, and quality
- Reduces health hazards
- Improves health care economics







cGMP Regulations

US FDA Title 21 Code of Federal Regulations (CFR) Parts:

- 11 Establishes the US FDA regulations on electronic records and electronic signatures (ERES)
- 210 /211 cGMP for pharmaceuticals
- 600/610 Covers biological products (*eg*, mAbs, vaccines, and blood)
- 820 Quality System Regulation (QSR) for medical devices





Drug Manufacturing

Drug Type	Small Molecule Drug	Large Molecule/ Biologic Drug
Example	Aspirin	IgG Antibody
Size	180 Da	150 kDa
Complexity Analogy	Bicycle ~ 20 lbs 150 parts	Private jet ~ 25,000 lbs 6,000,000 parts
Process Flow Schematic	$ \begin{array}{c} \bullet \\ \bullet $	
Manufacturing Process	For recovery unit Feed tanks Solvent Acetic anhydride Solvent Crystallizer Crystallizer Crystallizer Solvent Crystallizer Solvent Solvent Filter Filter Filter Solvent Filter Solvent Filter Solvent Filter Solvent Filter Solvent Filter To recovery unit	Cell expansion (for each batch) Cuese the cell ine (the use initial cell all (the use initial cell

Challenges with Biologics Manufacturing



ATCC[°]

Importance of Cell Banking & Storage





Cell Banking Market Drivers & Restraints

Global cell banking market is growing rapidly, but is challenged by regulatory climate and high costs

Market Restraint #1 Increasing regulatory stringency Market Driver #1 regulatory requirements Limited in-house storage and Clonality documentation requirements testing capabilities Host cell line history traceability Market Restraint #2 MCB & WCB guality control Increasing cell line testing demands Long-term continuity of 03 production processes 03 Market Restraint #3 High barrier to entry due to operational costs

- Significant investment in infrastructure and QMS
- Complex multi-step processes



Full-service outsourcing

Market Driver #2

- Improved operating efficiency
- Increased pipeline productivity

Market Driver #3

- Increased expenditures on healthcare (biologics)
- Advancements in regenerative medicine, cell & gene therapy, and tissue engineering
- Novel cryopreservation technologies and surging research in cell line development

11

ATCC – The Partner of Choice

- A pioneer in cell banking and storage
- Use of the most reliable and standardized procedures and equipment
- Provide critical starting materials for pharmaceuticals and biologics
- Trusted global source for cell authentication and uncompromised quality
- Avoid time and labor-intensive in-house production and eliminate associated risks and costs





ATCC Quality Overview

ATCC Quality Management Systems



- Maintain ISO 9001 certification across the broad organization
- ISO 13485 certification for synthetic standards, IVD test kits, and selected standards, controls and reagents
- Accredited for ISO 17025 and ISO 17034



Gaithersburg MD cGMP Facility

- Newly commissioned in 2021
- FDA 21 CFR part 11 compliant automated record keeping
- Our new purpose-built facility is the first to hold the distinction of being 21 CFR 600 and 610 compliant
- Examples of cGMP projects:
 - Expansion of mesenchymal stem cells, and production of exosomes (Pharma application)
 - MCB/WCB cell banks genetically modified CHO-K1 cell (Pharma application)
 - MCB/WCB KG-1A cells (Pharma application)





Expert cGMP Manufacture

Highlights

- Focus on Mammalian cells and stem cells
- Over 5,000 cell lines available from ATCC's fully authenticated collection -or- use your own proprietary cells
- Healthy cells, as well as cells derived from diseased tissues
- cGMP master and working cell banks (MCB and WCB)



ATCC°

Our Capabilities

	cGMP		
Compliance Standard	21 CFR 600, 610		
Customers	Biologics manufacturers Assay developers (potency & toxicology)	Human cells, tissues, and cellular and tissue-based product manufacturers (HCT/Ps)	
Cell Types	Mammalian cells Stem cells Primary cells	Expansion of cells (CDU)	
Cell Bank Types	 Production banks – cells will be used to make a "product" – produced in campaign "mode" Non-production banks - typically, bioassay banks – non-campaign – strict acceptance criteria 		
QC Release Testing	 Mycoplasma testing (Direct/indirect & PCR) Endotoxin testing Adventitious agents testing CO1 assay (interspecies testing) Sterility testing Virus panel Other tests as requested by customer and/or recommended for cell type 		



16

cGMP Cell Banking at ATCC



ATCC[®]

- Over 5K ATCC cell lines
- Customer's proprietary cells

cGMP Biorepository

- Dedicated and validated units, with access restricted to cGMP-trained staff with QC and QA oversight.
- Tracking labels are supplied to ensure the proper segregation immediately upon receipt at ATCC.
- Separate storage boxes for each cell bank or multiple banks in one box (customer's option). Only one customer per storage box.
- In addition to cGMP-compliant storage we also offer ISO 9001:2008
- We have storage at LN₂, -80°C, -20°C, and 2-8°C available
- Cell, microbe, protein, and nucleic acid storage options too
- Safe segregation of biological materials
- Cold-chain support and management
- On-call, after-hours personnel; 24/7 temperature monitoring
- Direct quality oversight of all unit entry and retrieval activities





Summary

- GMP is important for consumer safety
- Ensures that products meet specifications and quality/regulatory standards
- Biologics manufacturing is complex and requires additional considerations
- Use authenticated cell lines with established pedigrees under validated cGMP cell banking procedures
- Preserve the integrity of biologics cell banks long-term with ATCC cGMP biorepository services
- For more information please visit:

www.atcc.org/cGMP





References

- https://commons.wikimedia.org/wiki/File:Harvey_W._Wiley_%28masthead_picture%29.jpg
- https://www.crf-usa.org/bill-of-rights-in-action/bria-24-1-b-upton-sinclairs-the-jungle-muckraking-the-meatpacking-industry.html
- https://www.sciencephoto.com/media/1057609/view/sample-from-1937-sulfanilamide-poisoning-incident
- <u>https://www.history.com/news/7-things-you-may-not-know-about-the-jungle</u>
- <u>https://nhd-purefoodanddrugact.weebly.com/thesis.html</u>
- https://www.goodreads.com/book/show/13825242-federal-food-drug-and-cosmetic-act
- Alf van Beem, CC0, via Wikimedia Commons, <u>https://seita.com.co/Descargas/Aplicaciones/APN_Aspirin_Crystallization.pdf</u>
- ICH Quality Guidelines (Q5A Q5E Quality of Biotechnological Products)





ATCC[°]

Thank you and questions

Coming attractions:

hTERT-immortalized Primary Epidermal Cells: Key Components in Complex Toxicological Models July 28 at 12:00 PM EST Kevin Grady, BS

Skin Microbiome: Considerations, Applications, and Future Directions August 11 at 12:00 PM EST Tasha Santiago-Rodriguez, PhD

Proteomic Studies Using Tumor-derived Organoids (Title TBA) September 22 at 12:00 PM EST Chia-Fang Tsai, PhD

For more information about ATCC cGMP services:

www.atcc.org/cGMP

©2022 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.

